

Custom and Pre-Selected Kinase Profiling
to fit your Budget and Needs!

Kinase Profiling Book

As of October 1, 2023

128

Tyrosine
Kinases

197

Serine/Threonine
Kinases

27

Lipid
Kinases

Carna Biosciences, Inc.

Profiling Assays available from Carna Biosciences, Inc.

As of October 1, 2023

Page	Kinase Name	Assay Platform
5	ABL(ABL1)	MSA
5	ABL(ABL1)[E255K]	MSA
5	ABL(ABL1)[T315I]	MSA
5	ACK(TNK2)	MSA
5	AKT1	MSA
6	AKT2	MSA
6	AKT3	MSA
6	ALK	MSA
6	ALK[C1156Y]	MSA
6	ALK[F1174L]	MSA
7	ALK[G1202R]	MSA
7	ALK[G1269A]	MSA
7	ALK[L1196M]	MSA
7	ALK[R1275Q]	MSA
7	ALK[T1151_L1152insT]	MSA
8	EML4-ALK	MSA
8	NPM1-ALK	MSA
8	AMPKα1/β1/γ1(PRKAA1/B1/G1)	MSA
8	AMPKα2/β1/γ1(PRKAA2/B1/G1)	MSA
8	ARG(ABL2)	MSA
9	AurA(AURKA)	MSA
9	AurA(AURKA)/TPX2	MSA
9	AurB(AURKB)/INCENP	MSA
9	AurC(AURKC)	MSA
9	AXL	MSA
10	BLK	MSA
10	BMX	MSA
10	BRK(PTK6)	MSA
10	BRSK1	MSA
10	BRSK2	MSA
11	BTK	MSA
11	BTK[C481S]	MSA
11	BUB1/BUB3	MSA
11	CaMK1α(CAMK1)	MSA
11	CaMK1δ(CAMK1D)	MSA
12	CaMK2α(CAMK2A)	MSA
12	CaMK2β(CAMK2B)	MSA
12	CaMK2γ(CAMK2G)	MSA
12	CaMK2δ(CAMK2D)	MSA
12	CaMK4	MSA
13	CDC7/ASK	MSA
13	CDK1(CDC2)/CycB1	MSA
13	CDK2/CycA2	MSA
13	CDK2/CycE1	MSA
13	CDK3/CycE1	MSA
14	CDK4/CycD3	MSA
14	CDK5/p25	MSA
14	CDK6/CycD3	MSA
14	CDK7/CycH/MAT1	MSA
14	CDK9/CycT1	MSA
15	CGK2(PRKG2)	MSA
15	CHK1(CHEK1)	MSA
15	CHK2(CHEK2)	MSA
15	CK1α(CSNK1A1)	MSA
15	CK1γ1(CSNK1G1)	MSA
16	CK1γ2(CSNK1G2)	MSA
16	CK1γ3(CSNK1G3)	MSA
16	CK1δ(CSNK1D)	MSA
16	CK1ε(CSNK1E)	MSA
16	CK2α1/β(CSNK2A1/B)	MSA
17	CK2α2/β(CSNK2A2/B)	MSA
17	CLK1	MSA
17	CLK2	MSA
17	CLK3	MSA
17	CRIK(CIT)	MSA
18	CSK	MSA
18	DAPK1	MSA
18	DCAMKL2	MSA
18	DDR1	MSA
18	DDR2	MSA
19	DGKα(DGKA)	ADP-Glo
19	DGKβ(DGKB)	ADP-Glo
19	DGKγ(DGKG)	ADP-Glo
19	DGKδ(DGKD)	ADP-Glo
19	DGKε(DGKE)	ADP-Glo
20	DGKζ(DGKZ)	ADP-Glo
20	DGKη(DGKH)	ADP-Glo
20	DGKθ(DGKQ)	ADP-Glo
20	DGKi(DGKI)	ADP-Glo
20	DGKk(DGKK)	ADP-Glo
21	DYRK1A	MSA
21	DYRK1B	MSA
21	DYRK2	MSA
21	DYRK3	MSA
21	EEF2K	MSA

Page	Kinase Name	Assay Platform
22	EGFR	MSA
22	EGFR[C797S/L858R]	MSA
22	EGFR[d746-750]	MSA
22	EGFR[d746-750/C797S]	MSA
22	EGFR[d746-750/T790M]	MSA
23	EGFR[d746-750/T790M/C797S]	MSA
23	EGFR[D770_N771insNPG]	MSA
23	EGFR[L858R]	MSA
23	EGFR[L861Q]	MSA
23	EGFR[T790M]	MSA
24	EGFR[T790M/C797S/L858R]	MSA
24	EGFR[T790M/L858R]	MSA
24	EPHA1	MSA
24	EPHA2	MSA
24	EPHA3	MSA
25	EPHA4	MSA
25	EPHA5	MSA
25	EPHA6	MSA
25	EPHA7	MSA
25	EPHA8	MSA
26	EPHB1	MSA
26	EPHB2	MSA
26	EPHB3	MSA
26	EPHB4	MSA
26	Erk1(MAPK3)	MSA
27	Erk2(MAPK1)	MSA
27	Erk5(MAPK7)	MSA
27	FAK(PTK2)	MSA
27	FER	MSA
27	FES	MSA
28	FGFR1	MSA
28	FGFR1[V561M]	MSA
28	FGFR2	MSA
28	FGFR2[V564I]	MSA
28	FGFR3	MSA
29	FGFR3[K650E]	MSA
29	FGFR3[K650M]	MSA
29	FGFR3[V555L]	MSA
29	FGFR3[V555M]	MSA
29	FGFR4	MSA
30	FGFR4[N535K]	MSA
30	FGFR4[V550E]	MSA
30	FGFR4[V550L]	MSA
30	FGR	MSA
30	FLT1	MSA
31	FLT3	MSA
31	FLT4	MSA
31	FMS(CSF1R)	MSA
31	FRK	MSA
31	FYN[isoform a]	MSA
32	FYN[isoform b]	MSA
32	GSK3α(GSK3A)	MSA
32	GSK3β(GSK3B)	MSA
32	Haspin(GSG2)	MSA
32	HCK	MSA
33	HER2(ERBB2)	MSA
33	HER4(ERBB4)	MSA
33	HGK(MAP4K4)	MSA
33	HIPK1	MSA
33	HIPK2	MSA
34	HIPK3	MSA
34	HIPK4	MSA
34	HPK1(MAP4K1)	MSA
34	IGF1R	MSA
34	IKKα(CHUK)	IMAP
35	IKKβ(IKBKB)	MSA
35	IKKε(IKBKE)	MSA
35	INSR	MSA
35	IRAK1	IMAP
35	IRAK4	MSA
36	IRR(INSRR)	MSA
36	ITK	MSA
36	JAK1	MSA
36	JAK2	MSA
36	JAK3	MSA
37	JNK1(MAPK8)	MSA
37	JNK2(MAPK9)	MSA
37	JNK3(MAPK10)	MSA
37	KDR	MSA
37	KIT	MSA
38	KIT[D816E]	MSA
38	KIT[D816V]	MSA
38	KIT[D816Y]	MSA
38	KIT[T670I]	MSA
38	KIT[V560G]	MSA

Page	Kinase Name	Assay Platform
39	KIT[V654A]	MSA
39	LATS1/MOBKL1A	MSA
39	LATS2/MOBKL1A	MSA
39	LCK	MSA
39	LOK(STK10)	MSA
40	LTk	MSA
40	LYNa	MSA
40	LYNb	MSA
40	MAP4K2	MSA
40	MAPKAPK2	MSA
41	MAPKAPK3	MSA
41	MAPKAPK5	MSA
41	MARK1	MSA
41	MARK2	MSA
41	MARK3	MSA
42	MARK4	MSA
42	MELK	MSA
42	MER(MERTK)	MSA
42	MET	MSA
42	MET[D1228H]	MSA
43	MET[M1250T]	MSA
43	MET[Y1235D]	MSA
43	MINK(MINK1)	MSA
43	MNK1(MKNK1)	MSA
43	MNK2(MKNK2)	MSA
44	MRCKα(CDC42BPA)	MSA
44	MRCKβ(CDC42BPB)	MSA
44	MSK1(RPS6KA5)	MSA
44	MSK2(RPS6KA4)	MSA
44	MSSK1(STK23)	MSA
45	MST1(STK4)	MSA
45	MST2(STK3)	MSA
45	MST3(STK24)	MSA
45	MST4	MSA
45	MUSK	MSA
46	NDR1(STK38)	MSA
46	NDR2(STK38L)	MSA
46	NEK1	MSA
46	NEK2	MSA
46	NEK4	MSA
47	NEK6	MSA
47	NEK7	MSA
47	NEK9	MSA
47	NIM1K(MGC42105)	MSA
47	NuaK1	MSA
48	NuaK2	MSA
48	p38α(MAPK14)	MSA
48	p38β(MAPK11)	MSA
48	p38γ(MAPK12)	MSA
48	p38δ(MAPK13)	MSA
49	p70S6K(RPS6KB1)	MSA
49	p70S6KB2(RPS6KB2)	MSA
49	PAK1	MSA
49	PAK2	MSA
49	PAK4	MSA
50	PAK5(PAK7)	MSA
50	PAK6	MSA
50	PASK	MSA
50	PBK	MSA
50	PDGFRα(PDGFRα)	MSA
51	PDGFRα(PDGFRα)[D842V]	MSA
51	PDGFRα(PDGFRα)[T674I]	MSA
51	PDGFRα(PDGFRα)[V561D]	MSA
51	PDGFRβ(PDGFRβ)	MSA
51	PDHK2(PDK2)	MSA
52	PDHK4(PDK4)	MSA
52	PDK1(PDPK1)	MSA
52	PEK(EIF2AK3)	IMAP
52	PGK(PRKG1)	MSA
52	PHKG1	MSA
53	PHKG2	MSA
53	PIK3CA/PIK3R1	ADP-Glo
53	PIK3CA[E542K]/PIK3R1	ADP-Glo
53	PIK3CA[E545K]/PIK3R1	ADP-Glo
53	PIK3CA[H1047R]/PIK3R1	ADP-Glo
54	PIK3CA[P539R]/PIK3R1	ADP-Glo
54	PIK3CA[R88Q]/PIK3R1	ADP-Glo
54	PIK3CB/PIK3R1	ADP-Glo
54	PIK3CD/PIK3R1	ADP-Glo
54	PIKFYVE(PIP5K3)	ADP-Glo
55	PIM1	MSA
55	PIM2	MSA
55	PIM3	MSA
55	PIP4K2A	ADP-Glo
55	PIP4K2B	ADP-Glo
56	PIP5K1A	ADP-Glo
56	PIP5K1B	ADP-Glo
56	PIP5K1C	ADP-Glo
56	PIP5KL1	ADP-Glo
56	PKACα(PRKACA)	MSA

Page	Kinase Name	Assay Platform
57	PKACβ(PRKACB)	MSA
57	PKACγ(PRKACG)	MSA
57	PKCα(PRKCA)	MSA
57	PKCβ1(PRKCB1)	MSA
57	PKCβ2(PRKCB2)	MSA
58	PKCγ(PRKCG)	MSA
58	PKCδ(PRKCD)	MSA
58	PKCε(PRKCE)	MSA
58	PKCζ(PRKCZ)	MSA
58	PKCη(PRKCH)	MSA
59	PKCθ(PRKCQ)	MSA
59	PKCi(PRKCI)	MSA
59	PKD1(PRKD1)	MSA
59	PKD2(PRKD2)	MSA
59	PKD3(PRKD3)	MSA
60	PKN1	IMAP
60	PKR(EIF2AK2)	IMAP
60	PLK1	MSA
60	PLK2	IMAP
60	PLK3	MSA
61	PRKX	MSA
61	PYK2(PTK2B)	MSA
61	QIK(SNF1LK2)	MSA
61	RET	MSA
61	RET[G691S]	MSA
62	RET[M918T]	MSA
62	RET[S891A]	MSA
62	RET[Y791F]	MSA
62	ROCK1	MSA
62	ROCK2	MSA
63	RON(MST1R)	MSA
63	ROS(ROS1)	MSA
63	RSK1(RPS6KA1)	MSA
63	RSK2(RPS6KA3)	MSA
63	RSK3(RPS6KA2)	MSA
64	RSK4(RPS6KA6)	MSA
64	SGK	MSA
64	SGK2	MSA
64	SGK3(SGKL)	MSA
64	SIK(SNF1LK)	MSA
65	skMLCK(MYLK2)	MSA
65	SLK	MSA
65	SPHK1	MSA
65	SPHK2	MSA
65	SRC	MSA
66	SRM(SRMS)	MSA
66	SRPK1	IMAP
66	SRPK2	MSA
66	SYK	MSA
66	TAK1-TAB1(MAP3K7)	MSA
67	TAOK2	MSA
67	TBK1	MSA
67	TEC	MSA
67	TIE2(TEK)	MSA
67	TNIK	MSA
68	TNK1	MSA
68	TRKA(NTRK1)	MSA
68	TRKB(NTRK2)	MSA
68	TRKC(NTRK3)	MSA
68	TSSK1	MSA
69	TSSK2	MSA
69	TSSK3	MSA
69	TXK	MSA
69	TYK2	MSA
69	TYRO3	MSA
70	WNK1	MSA
70	WNK2	MSA
70	WNK3	MSA
70	YES(YES1)	MSA
70	YES(YES1)[T348I]	MSA
71	ZAP70	MSA

<< Cascade Assay >>

Page	Kinase Name	Assay Platform
72	BRAF	MSA
72	BRAF[V600E]	MSA
72	COT(MAP3K8)	MSA
72	DLK(MAP3K12)	MSA
72	MAP2K1	MSA
73	MAP2K2	MSA
73	MAP2K3	MSA
73	MAP2K4	MSA
73	MAP2K5	MSA
73	MAP2K6	MSA
74	MAP2K7	MSA
74	MAP3K1	MSA
74	MAP3K2	MSA
74	MAP3K3	MSA
74	MAP3K4	MSA
75	MAP3K5	MSA
75	MLK1(MAP3K9)	MSA
75	MLK2(MAP3K10)	MSA
75	MLK3(MAP3K11)	MSA
75	MOS	MSA
76	RAF1	MSA

- The Kinase Company -

Carna Biosciences, Inc.

TEL : +81-78-302-7091 / FAX: +81-78-302-7086

Email: info@carnabio.com

www.carnabio.com

ABL(ABL1)

Product code 08-001

Full-length human ABL [2-1130(end) amino acids of accession number NP_005148.2] was expressed as N-terminal His-tagged protein (126 kDa) using baculovirus expression system. His-tagged ABL was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: ABLtide
ATP (μM) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 75
IC50 at 1 mM ATP (nM)	: 1300

ABL(ABL1)[E255K]

Product code 08-094

Full-length human ABL [2-1130(end) amino acids and E255K of accession number NP_005148.2] was expressed as N-terminal His-tagged protein (126 kDa) using baculovirus expression system. His-tagged ABL[E255K] was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: ABLtide
ATP (μM) Km app / Bin	: 17 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 140
IC50 at 1 mM ATP (nM)	: 4500

ABL(ABL1)[T315I]

Product code 08-093

Full-length human ABL [2-1130(end) amino acids and T315I of accession number NP_005148.2] was expressed as N-terminal His-tagged protein (126 kDa) using baculovirus expression system. His-tagged ABL[T315I] was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: ABLtide
ATP (μM) Km app / Bin	: 4 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 6.4
IC50 at 1 mM ATP (nM)	: 890

ACK(TNK2)

Product code 08-196

Human ACK, catalytic domain [110-476 amino acids of accession number NP_005772.3] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-ACK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: WASP peptide
ATP (μM) Km app / Bin	: 97 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.2
IC50 at 1 mM ATP (nM)	: 3.8

AKT1

Product code 01-101

Human AKT1, catalytic domain [104-480(end) amino acids of accession number NP_005154.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-AKT1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Crosstide
ATP (μM) Km app / Bin	: 31 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.7
IC50 at 1 mM ATP (nM)	: 22

AKT2

Product code 01-102

Human AKT2, catalytic domain [120-481(end) amino acids of accession number NP_001617.1] was co-expressed as N-terminal GST-fusion protein (69 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-AKT2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Crosstide
ATP (μM) Km app / Bin	: 110 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 5.2
IC50 at 1 mM ATP (nM)	: n.a.

AKT3

Product code 01-103

Human AKT3, catalytic domain [108-479(end) amino acids of accession number NP_005456.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-AKT3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Crosstide
ATP (μM) Km app / Bin	: 54 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.2
IC50 at 1 mM ATP (nM)	: n.a.

ALK

Product code 08-518

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 57 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.5
IC50 at 1 mM ATP (nM)	: 15

ALK[C1156Y]

Product code 08-530

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and C1156Y, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[C1156Y] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 64 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.9
IC50 at 1 mM ATP (nM)	: 11

ALK[F1174L]

Product code 08-519

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and F1174L, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[F1174L] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 49 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.4
IC50 at 1 mM ATP (nM)	: 21

ALK[G1202R]

Product code 08-544

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and G1202R, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[G1202R] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 31 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 7.3
IC50 at 1 mM ATP (nM)	: 69

ALK[G1269A]

Product code 08-537

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and G1269A, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[G1269A] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.36
IC50 at 1 mM ATP (nM)	: 1.6

ALK[L1196M]

Product code 08-529

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and L1196M, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[L1196M] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 57 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.66
IC50 at 1 mM ATP (nM)	: 4.3

ALK[R1275Q]

Product code 08-520

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and R1275Q, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[R1275Q] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 84 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.3
IC50 at 1 mM ATP (nM)	: 16

ALK[T1151_L1152insT]

Product code 08-539

Human ALK , cytoplasmic domain [1058-1620(end) amino acids and T1151_L1152insT, I1461V, K1491R, D1529E of accession number NP_004295.2] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-ALK[T1151_L1152insT] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 110 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 6.5
IC50 at 1 mM ATP (nM)	: 16

EML4-ALK

Product code 08-516

Fused gene of human fusion EML4-ALK [1-1059 amino acids of accession number BAF73611.1] was expressed as N-terminal GST-fusion protein (145 kDa) using baculovirus expression system. GST-EML4-ALK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 43 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.9
IC50 at 1 mM ATP (nM)	: 16

NPM1-ALK

Product code 08-517

Fused gene of human fusion NPM1-ALK [1-680 amino acids of accession number BAA08343.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-NPM1-ALK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 57 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.4
IC50 at 1 mM ATP (nM)	: 14

AMPK α 1/ β 1/ γ 1(PRKAA1/B1/G1)

Product code 02-113

Full-length human AMPK α 1 [1-550(end) amino acids of accession number NP_006242.4] was co-expressed as N-terminal GST-fusion protein (90 kDa) with GST-PRKAB1 [1-270(end) amino acids of accession number NP_006244.2] and PRKAG1 [1-331(end) amino acids of accession number NP_002724.1] using baculovirus expression system. GST-AMPK α 1/ β 1/ γ 1 was purified by using glutathione sepharose chromatography and activated with His-tagged CaMKK1. Activated GST-AMPK α 1/ β 1/ γ 1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SAMS peptide
ATP (μ M) Km app / Bin	: 130 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.41
IC50 at 1 mM ATP (nM)	: 0.87

AMPK α 2/ β 1/ γ 1(PRKAA2/B1/G1)

Product code 02-114

Full-length human AMPK α 2 [1-552(end) amino acids of accession number NP_006243.2] was co-expressed as N-terminal GST-fusion protein (89 kDa) with GST-PRKAB1 [1-270(end) amino acids of accession number NP_006244.2] and PRKAG1 [1-331(end) amino acids of accession number NP_002724.1] using baculovirus expression system. GST-AMPK α 2/ β 1/ γ 1 was purified by using glutathione sepharose chromatography and activated with His-tagged CaMKK1. Activated GST-AMPK α 2/ β 1/ γ 1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SAMS peptide
ATP (μ M) Km app / Bin	: 100 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.79
IC50 at 1 mM ATP (nM)	: n.a.

ARG(ABL2)

Product code 08-102

Full-length human ARG [2-1161(end) amino acids of accession number NP_001161708.1] was expressed as N-terminal GST-fusion protein (153 kDa) using baculovirus expression system. GST-ARG was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: ABL tide
ATP (μ M) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 27
IC50 at 1 mM ATP (nM)	: 400

AurA(AURKA)

Product code 05-101

Full-length human AurA [1-403(end) amino acids of accession number NP_940835.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-AurA was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.8
IC50 at 1 mM ATP (nM)	: 17

AurA(AURKA)/TPX2

Product code 05-186

Full-length human AurA [1-403(end) amino acids of accession number NP_940835.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-AurA was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 1.7 / 2
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 6.1
IC50 at 1 mM ATP (nM)	: n.a.

AurB(AURKB)/INCENP

Product code 05-102

Full-length human AurB [1-344(end) amino acids and M298T of accession number NP_004208.2] was co-expressed as N-terminal GST-fusion protein (66 kDa) with His-tagged INCENP(INBOX) [803-918(end) amino acids of accession number NP_001035784.1] using baculovirus expression system. GST-AurB was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 7.1
IC50 at 1 mM ATP (nM)	: 62

AurC(AURKC)

Product code 05-103

Full-length human AurC [1-275(end) amino acids of accession number NP_003151.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system. GST-AurC was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.1
IC50 at 1 mM ATP (nM)	: 18

AXL

Product code 08-107

Human AXL, cytoplasmic domain [464-885(end) amino acids of accession number NP_001690.2] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-AXL was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 32 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 7.9

BLK

Product code 08-164

Full-length human BLK [1-505(end) amino acids of accession number NP_001706.2] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-BLK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 62 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.6
IC50 at 1 mM ATP (nM)	: 17

BMX

Product code 08-179

Full-length human BMX [1-675(end) amino acids of accession number NP_001712.1] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-BMX was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 75 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 45

BRK(PTK6)

Product code 08-165

Full-length human BRK [2-451(end) amino acids of accession number NP_005966.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-BRK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 250 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 260
IC50 at 1 mM ATP (nM)	: 390

BRSK1

Product code 02-115

Full-length human BRSK1 [1-778(end) amino acids of accession number NP_115806.1] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system. GST-BRSK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHK tide
ATP (μ M) Km app / Bin	: 30 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.27
IC50 at 1 mM ATP (nM)	: 0.57

BRSK2

Product code 02-116

Full-length human BRSK2 [1-674(end) amino acids of accession number ABA17261.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system. GST-BRSK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHK tide
ATP (μ M) Km app / Bin	: 31 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.31
IC50 at 1 mM ATP (nM)	: n.a.

BTK

Product code 08-180

Full-length human BTK [2-659(end) amino acids of accession number NP_000052] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-BTK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 22 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 24
IC50 at 1 mM ATP (nM)	: 93

BTK[C481S]

Product code 08-547

Full-length human BTK [2-659(end) amino acids and C481S of accession number NP_000052] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-BTK[C481S] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 37
IC50 at 1 mM ATP (nM)	: 170

BUB1/BUB3

Product code 05-187

Full-length human BUB1 [1-1085 (end) amino acids of accession number NP_004327] was co-expressed as N-terminal GST-fusion protein (149 kDa) with DYKDDDDK tagged BUB3 [1-328 (end) amino acids of accession number NP_004716] using baculovirus expression system. GST-BUB1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: H2A peptide
ATP (μM) Km app / Bin	: 2.9 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 300
IC50 at 1 mM ATP (nM)	: n.a.

CaMK1α(CAMK1)

Product code 02-104

Full-length human CaMK1α [1-370(end) amino acids of accession number NP_003647.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-CaMK1α was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 750 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 16
IC50 at 1 mM ATP (nM)	: 16

CaMK1δ(CAMK1D)

Product code 02-106

Full-length human CaMK1δ[1-357(end) amino acid of accession number NP_065130.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-CaMK1δ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Synapsin peptide
ATP (μM) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.5
IC50 at 1 mM ATP (nM)	: n.a.

CaMK2α(CAMK2A)

Product code 02-109

Full-length human CaMK2α [1-478(end) amino acids of accession number NP_741960.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-CaMK2α was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 33 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.75
IC50 at 1 mM ATP (nM)	: 3.2

CaMK2β(CAMK2B)

Product code 02-110

Full-length human CaMK2β [1-503 amino acids of accession number NP_742078.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-CaMK2β was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 19 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.54
IC50 at 1 mM ATP (nM)	: 5.8

CaMK2γ(CAMK2G)

Product code 02-112

Full-length human CaMK2γ[1-518(end) amino acids of accession number NP_751910.1] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-CaMK2γ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 23 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.39
IC50 at 1 mM ATP (nM)	: 1.1

CaMK2δ(CAMK2D)

Product code 02-111

Full-length human CaMK2δ [1-478 amino acids of accession number NP_742113.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-CaMK2δ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 6.3 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.26
IC50 at 1 mM ATP (nM)	: 0.64

CaMK4

Product code 02-108

Full-length human CaMK4 [1-473(end) amino acids of accession number NP_001735.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-CaMK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 140
IC50 at 1 mM ATP (nM)	: 1000

CDC7/ASK

Product code 05-109

Full-length human CDC7 [1-574(end) amino acids of accession number NP_003494.1] was co-expressed as N-terminal GST-fusion protein (92 kDa) with Dbf4(ASK) [1-674(end) amino acids of accession number NP_006707.1] using baculovirus expression system. GST-CDC7 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MCM2 peptide
ATP (μM) Km app / Bin	: 2.8 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 16
IC50 at 1 mM ATP (nM)	: 1600

CDK1(CDC2)/CycB1

Product code 04-102

Full-length human CDK1 [1-297(end) amino acids of accession number NP_001777.1] was co-expressed as N-terminal GST-fusion protein (61 kDa) with CyclinB1 [1-433(end) amino acids of accession number NP_114172.1] using baculovirus expression system. GST-CDK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Histone H1
ATP (μM) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.3
IC50 at 1 mM ATP (nM)	: 32

CDK2/CycA2

Product code 04-103

Full-length human CDK2 [1-298(end) amino acids of accession number NP_001789.2] was co-expressed as N-terminal GST-tagged protein (61 kDa) with GST-CyclinA2 [1-432(end) amino acids of accession number NP_001228.1] using baculovirus expression system. GST-CDK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Histone H1
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 7.1

CDK2/CycE1

Product code 04-165

Full-length human CDK2 [1-298(end) amino acids of accession number NP_001789.2] was co-expressed as N-terminal GST-tagged protein (61 kDa) with CyclinE1 [1-410(end) amino acids of accession number NP_001229.1] using baculovirus expression system. GST-CDK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Histone H1
ATP (μM) Km app / Bin	: 130 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.8
IC50 at 1 mM ATP (nM)	: 10

CDK3/CycE1

Product code 04-104

Full-length human CDK3 [1-305(end) amino acids of accession number NP_001249.1] was co-expressed as N-terminal GST-fusion protein (62kDa) with CyclinE1 [1-410(end) amino acids of accession number NP_001229.1] using baculovirus expression system. GST-CDK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Histone H1
ATP (μM) Km app / Bin	: 1000 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.4
IC50 at 1 mM ATP (nM)	: 3.4

CDK4/CycD3

Product code 04-105

Full-length human CDK4 [1-303(end) amino acids of accession number NP_000066.1] was co-expressed as N-terminal GST-fusion protein (61 kDa) with human GST-CyclinD3 [1-292(end) amino acids and S259A of accession number NP_001751.1] using baculovirus expression system. GST-CDK4/CycD3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 200 / 200
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 13
IC50 at 1 mM ATP (nM)	: 52

CDK5/p25

Product code 04-106

Full-length human CDK5 [1-292(end) amino acids of accession number NP_004926.1] was co-expressed as N-terminal GST-fusion protein (60 kDa) with p25 [99-307(end) amino acids of accession number NP_003876.1] using baculovirus expression system. GST-CDK5 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Histone H1
ATP (μM) Km app / Bin	: 10 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.5
IC50 at 1 mM ATP (nM)	: 86

CDK6/CycD3

Product code 04-107

Full-length human CDK6 [1-326(end) amino acids of accession number NP_001250.1] was co-expressed as N-terminal GST-fusion protein (64 kDa) with human GST-CyclinD3 [1-292(end) amino acids and S259A of accession number NP_001751.1] using baculovirus expression system. GST-CDK6/CycD3 was purified by using glutathione sepharose chromatography and activated with His-CDK7/CycH/MAT1. Activated GST-CDK6/CycD3 was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 330 / 300
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 58
IC50 at 1 mM ATP (nM)	: 110

CDK7/CycH/MAT1

Product code 04-108

Full-length human CDK7 [1-346(end) amino acids of accession number NP_001790.1] was co-expressed as N-terminal GST-fusion protein (66 kDa) with CyclinH [1-323(end) amino acids of accession number NP_001230.1] and MAT1 [1-309(end) amino acids of accession number NP_002422.1] using baculovirus expression system. GST-CDK7 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CTD3 peptide
ATP (μM) Km app / Bin	: 32 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 17
IC50 at 1 mM ATP (nM)	: 120

CDK9/CycT1

Product code 04-110

Full-length human CDK9 [1-372(end) amino acids of accession number NP_001252.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-CyclinT1 [1-726(end) amino acids of accession number NP_001231.2] using baculovirus expression system. GST-CDK9 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDK9 substrate
ATP (μM) Km app / Bin	: 9.4 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 5.2
IC50 at 1 mM ATP (nM)	: 130

CGK2(PRKG2)

Product code 01-143

Full-length human CGK2 [1-762(end) amino acids of accession number NP_006250.1] was expressed as N-terminal GST-fusion protein (114 kDa) using baculovirus expression system. GST-CGK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μM) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.88
IC50 at 1 mM ATP (nM)	: n.a.

CHK1(CHEK1)

Product code 02-117

Full-length human CHK1 [1-476(end) amino acids of accession number NP_001265.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-CHK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μM) Km app / Bin	: 50 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.24
IC50 at 1 mM ATP (nM)	: 1.1

CHK2(CHEK2)

Product code 02-162

Full-length human CHK2 [1-543(end) amino acids of accession number NP_009125.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-CHK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μM) Km app / Bin	: 51 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 11
IC50 at 1 mM ATP (nM)	: 25

CK1α(CSNK1A1)

Product code 03-101

Full-length human CK1α [1-337(end) amino acids of accession number NP_001883.4] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system. GST-CK1α was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μM) Km app / Bin	: 4.1 / 5
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 150
IC50 at 1 mM ATP (nM)	: >10000

CK1γ1(CSNK1G1)

Product code 03-105

Full-length human CK1γ1 [1-422(end) amino acids of accession number NP_071331.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-CK1γ1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μM) Km app / Bin	: 6.3 / 5
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 1300
IC50 at 1 mM ATP (nM)	: n.a.

CK1 γ 2(CSNK1G2)

Product code 03-106

Full-length human CK1 γ 2 [1-415(end) amino acids of accession number NP_001310.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-CK1 γ 2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μ M) Km app / Bin	: 10 / 10
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 510
IC50 at 1 mM ATP (nM)	: n.a.

CK1 γ 3(CSNK1G3)

Product code 03-107

Full-length human CK1 γ 3 [1-447(end) amino acids of accession number NP_004375.2] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system. GST-CK1 γ 3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μ M) Km app / Bin	: 3.2 / 5
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 920
IC50 at 1 mM ATP (nM)	: n.a.

CK1 δ (CSNK1D)

Product code 03-103

Human CK1 δ , catalytic domain [1-294 amino acids of accession number NP_001884.2] was expressed as N-terminal GST-fusion protein (61 kDa) using E. coli expression system. GST-CK1 δ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μ M) Km app / Bin	: 7.7 / 10
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 25
IC50 at 1 mM ATP (nM)	: 570

CK1 ϵ (CSNK1E)

Product code 03-104

Human CK1 ϵ , catalytic domain [1-348 amino acids of accession number NP_001885.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-CK1 ϵ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: 5-Iodotubercidin
IC50 at ATP Bin (nM)	: 300
IC50 at 1 mM ATP (nM)	: 5800

CK2 α 1/ β (CSNK2A1/B)

Product code 05-184

Full-length human CK2 α 1 [1-391(end) amino acids of accession number NP_001886.1] was co-expressed as N-terminal GST-fusion protein (72 kDa) with human His-tagged CK2 β [1-215 amino acids of accession number NP_001311.3] using baculovirus expression system. GST-CK2 α 1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CK2tide
ATP (μ M) Km app / Bin	: 2.9 / 5
Metal	: Mg
Reference compound	: TBB
IC50 at ATP Bin (nM)	: 60
IC50 at 1 mM ATP (nM)	: 4800

CK2 α 2/ β (CSNK2A2/B)

Product code 05-185

Full-length human CK2 α 2 [1-350(end) amino acids of accession number NP_001887.1] was co-expressed as N-terminal GST-fusion protein (68 kDa) with human His-tagged CK2 β [1-215 amino acids of accession number NP_001311.3] using baculovirus expression system. GST-CK2 α 2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CK2tide
ATP (μ M) Km app / Bin	: 2.1 / 5
Metal	: Mg
Reference compound	: TBB
IC50 at ATP Bin (nM)	: 50
IC50 at 1 mM ATP (nM)	: n.a.

CLK1

Product code 04-126

Human CLK1, catalytic domain [129-484(end) amino acids of accession number NP_004062.2] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-CLK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 9.6
IC50 at 1 mM ATP (nM)	: 60

CLK2

Product code 04-127

Full-length human CLK2 [1-499(end) amino acids of accession number AAH53603.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-CLK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 140 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.6
IC50 at 1 mM ATP (nM)	: 28

CLK3

Product code 04-128

Full-length human CLK3 [1-490(end) amino acids of accession number AAH02555.1] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system. GST-CLK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 75 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 820
IC50 at 1 mM ATP (nM)	: n.a.

CRK(CIT)

Product code 01-104

Human citron kinase (CRK), catalytic domain [1-449 amino acids of accession number NP_009105.1] was expressed as N-terminal GST fusion protein (77 kDa) using baculovirus expression system. GST-CRK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Histone H3 peptide
ATP (μ M) Km app / Bin	: 7.8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 31
IC50 at 1 mM ATP (nM)	: n.a.

CSK

Product code 08-111

Full-length human CSK [1-450(end) amino acids of accession number NP_004374.1] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system. GST-CSK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 4.8 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 53
IC50 at 1 mM ATP (nM)	: 1500

DAPK1

Product code 02-134

Human DAPK1, catalytic domain [1-289 amino acids of accession number NP_004929.1] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-DAPK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DAPK1 tide
ATP (μ M) Km app / Bin	: 1.1 / 1
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.9
IC50 at 1 mM ATP (nM)	: 490

DCAMKL2

Product code 02-140

Full-length human DCAMKL2 [1-695(end) amino acids of accession number NP_689832.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-DCAMKL2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 120 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 22
IC50 at 1 mM ATP (nM)	: n.a.

DDR1

Product code 08-113

Human DDR1, cytoplasmic domain [444-876(end) amino acids of accession number NP_001945.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-DDR1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μ M) Km app / Bin	: 94 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.0
IC50 at 1 mM ATP (nM)	: 3.1

DDR2

Product code 08-114

Human DDR2, cytoplasmic domain [422-855(end) amino acids of accession number NP_006173.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-DDR2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μ M) Km app / Bin	: 38 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 0.77

DGK α (DGKA)

Product code 12-401-20N

Full-length human DGK α [1-735(end) amino acids of accession number NP_958852.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (87 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μ M) Km app / Bin	: 130 / 100
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGK β (DGKB)

Product code 12-402-20N

Full-length human DGK β [1-803(end) amino acids of accession number NP_001337638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (95 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and activated with ATP. Activated protein was purified by using gel filtration.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μ M) Km app / Bin	: 61 / 50
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGK γ (DGKG)

Product code 12-403-20N

Full-length human DGK γ [1-791(end) amino acids and T142S and R316K of accession number NP_001337.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (94 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μ M) Km app / Bin	: 55 / 50
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGK δ (DGKD)

Product code 12-404-20N

Catalytic domain of human DGK δ [1-1141 amino acids of accession number NP_690618.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (131 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μ M) Km app / Bin	: 120 / 100
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGK ϵ (DGKE)

Product code 12-415-20N

Truncated human DGK ϵ [48-567(end) amino acids of accession number NP_003638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (63 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μ M) Km app / Bin	: 120 / 100
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGKζ(DGKZ)

Product code 12-410-20N

Full-length human DGKζ [1-929(end) amino acids of accession number NP_003637.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (109 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μM) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGKη(DGKH)

Product code 12-406-20N

Catalytic domain of human DGKη [1-1147 amino acids of accession number NP_821077.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (131 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μM) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGKθ(DGKQ)

Product code 12-409-20N

Full-length human DGKθ [1-942(end) amino acids of accession number NP_001338.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (106 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μM) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGKι(DGKI)

Product code 12-407-20N

Full-length human DGKι [1-1065(end) amino acids of accession number NP_004708.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (122 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μM) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DGKκ(DGKK)

Product code 12-408-20N

Full-length human DGKκ [1-1271(end) amino acids and D1118N of accession number NP_001013764.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (146 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μM) Km app / Bin	: 17 / 25
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	:
IC50 at 1 mM ATP (nM)	:

DYRK1A

Product code 04-130

Full-length human DYRK1A [1-763(end) amino acids of accession number NP_001387.2] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system. GST-DYRK1A was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 7.8
IC50 at 1 mM ATP (nM)	: 120

DYRK1B

Product code 04-131

Full-length human DYRK1B [1-629(end) amino acids of accession number NP_004705.1] was expressed as N-terminal GST-fusion protein (96 kDa) using baculovirus expression system. GST-DYRK1B was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 59 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 32

DYRK2

Product code 04-132

Full-length human DYRK2 [1-528(end) amino acids of accession number NP_003574.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-DYRK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 7.7 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 130
IC50 at 1 mM ATP (nM)	: 6300

DYRK3

Product code 04-133

Full-length human DYRK3 [1-588(end) amino acids of accession number NP_003573.2] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system. GST-DYRK3 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 6.8 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 17
IC50 at 1 mM ATP (nM)	: 1200

EEF2K

Product code 10-113

Full-length human EEF2K [1-725(end) amino acids of accession number NP_037434.1] was expressed as N-terminal GST-fusion protein (109 kDa) using E. coli expression system. GST-EEF2K was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: EEF2Ktide
ATP (μM) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: A-484954
IC50 at ATP Bin (nM)	: 330
IC50 at 1 mM ATP (nM)	: n.a.

EGFR

Product code 08-115

Human EGFR, cytoplasmic domain [669-1210(end) amino acids of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 2.7 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 53
IC50 at 1 mM ATP (nM)	: 7700

EGFR[C797S/L858R]

Product code 08-563

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and C797S/L858R of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[C797S/L858R] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 4.1 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 8.8
IC50 at 1 mM ATP (nM)	: 270

EGFR[d746-750]

Product code 08-527

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-EGFR[d746-750aa] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 19 / 25
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 13
IC50 at 1 mM ATP (nM)	: 93

EGFR[d746-750/C797S]

Product code 08-564

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and C797S of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-EGFR[d746-750aa/C797S] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 8.2 / 10
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 8.0
IC50 at 1 mM ATP (nM)	: 130

EGFR[d746-750/T790M]

Product code 08-528

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and T790M of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR [d746-750aa/T790M] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 5.4 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.52
IC50 at 1 mM ATP (nM)	: 9.7

EGFR[d746-750/T790M/C797S]

Product code 08-565

Human EGFR, cytoplasmic domain [669-745, 751-1210(end) amino acids and T790M/C797S of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-EGFR[d746-750aa/T790M/C797S] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 1.8 / 2
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.84
IC50 at 1 mM ATP (nM)	: 14

EGFR[D770_N771insNPG]

Product code 08-553

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and D770_N771insNPG of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[D770_N771insNPG] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 2.3 / 5
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 18
IC50 at 1 mM ATP (nM)	: 930

EGFR[L858R]

Product code 08-502

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and L858R of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[L858R] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 9.8 / 10
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 11
IC50 at 1 mM ATP (nM)	: 360

EGFR[L861Q]

Product code 08-513

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and L861Q of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[L861Q] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 7.5 / 10
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 68
IC50 at 1 mM ATP (nM)	: 2200

EGFR[T790M]

Product code 08-194

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and T790M of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[T790M] was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 0.9 / 1
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.8
IC50 at 1 mM ATP (nM)	: 190

EGFR[T790M/C797S/L858R]

Product code 08-559

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and T790M/C797S/L858R of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[T790M/C797S/L858R] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 0.85 / n.a.
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: n.a.
IC50 at 1 mM ATP (nM)	: 37

EGFR[T790M/L858R]

Product code 08-510

Human EGFR, cytoplasmic domain [669-1210(end) amino acids and T790M/L858R of accession number NP_005219.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-EGFR[T790M/L858R] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 1.9 / 2
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 56

EPHA1

Product code 08-119

Human EPHA1, cytoplasmic domain [586-976(end) amino acids of accession number NP_005223.3] was expressed as N-terminal GST-fusion protein (71 kDa) using baculovirus expression system. GST-EPHA1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyotide
ATP (μ M) Km app / Bin	: 22 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 20
IC50 at 1 mM ATP (nM)	: 340

EPHA2

Product code 08-121

Human EPHA2, cytoplasmic domain [572-976(end) amino acids of accession number NP_004422.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHA2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyotide
ATP (μ M) Km app / Bin	: 67 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 160
IC50 at 1 mM ATP (nM)	: 530

EPHA3

Product code 08-122

Human EPHA3, cytoplasmic domain [579-983(end) amino acids of accession number NP_005224.2] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-EPHA3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyotide
ATP (μ M) Km app / Bin	: 170 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 37
IC50 at 1 mM ATP (nM)	: 76

EPHA4

Product code 08-123

Human EPHA4, cytoplasmic domain [586-986(end) amino acids of accession number NP_004429.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-EPHA4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 52 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 50
IC50 at 1 mM ATP (nM)	: 330

EPHA5

Product code 08-124

Human EPHA5, catalytic domain [662-948 amino acids of accession number NP_004430.3] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system. GST-EPHA5 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 34
IC50 at 1 mM ATP (nM)	: 220

EPHA6

Product code 08-125

Human EPHA6, cytoplasmic domain [683-1130(end) amino acids of accession number NP_001073917.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-EPHA6 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 17
IC50 at 1 mM ATP (nM)	: 60

EPHA7

Product code 08-126

Human EPHA7, cytoplasmic domain [595-998(end) amino acids of accession number NP_004431.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHA7 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 58 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 48
IC50 at 1 mM ATP (nM)	: 480

EPHA8

Product code 08-127

Human EPHA8, catalytic domain [571-924 amino acids of accession number NP_065387.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-EPHA8 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 69 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 61
IC50 at 1 mM ATP (nM)	: 240

EPHB1

Product code 08-128

Human EPHB1, cytoplasmic domain [578-984(end) amino acids of accession number NP_004432.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHB1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 29 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 53
IC50 at 1 mM ATP (nM)	: 760

EPHB2

Product code 08-129

Human EPHB2, cytoplasmic domain [581-987(end) amino acids of accession number NP_004433.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHB2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 86 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 73
IC50 at 1 mM ATP (nM)	: 400

EPHB3

Product code 08-130

Human EPHB3, cytoplasmic domain [596-998(end) amino acids of accession number NP_004434.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-EPHB3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 49 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2000
IC50 at 1 mM ATP (nM)	: >10000

EPHB4

Product code 08-131

Human EPHB4, cytoplasmic domain [577-987(end) amino acids of accession number NP_004435.3] was expressed as N-terminal GST-protein (73 kDa) using baculovirus expression system. GST-EPHB4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 230
IC50 at 1 mM ATP (nM)	: 1500

Erk1(MAPK3)

Product code 04-142

Full-length human Erk1 [1-379(end) amino acids of accession number NP_002737.1] was expressed as N-terminal GST-fusion protein (70 kDa) using E.coli expression system. GST-Erk1 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K1. Activated GST-Erk1 was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 37
IC50 at 1 mM ATP (nM)	: 400

Erk2(MAPK1)

Product code 04-143

Full-length human Erk2 [1-360(end) amino acids of accession number NP_002736.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E.coli expression system. GST-Erk2 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K1. Activated GST-Erk2 was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 33 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 21
IC50 at 1 mM ATP (nM)	: 180

Erk5(MAPK7)

Product code 04-146

Human Erk5, catalytic domain [1-398 amino acids of accession number NP_002740.2] was expressed as N-terminal GST-fusion protein (72 kDa) using E. coli expression system. GST-Erk5 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K5. Activated GST-Erk5 was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: EGFR-derived peptide
ATP (μM) Km app / Bin	: 450 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 280
IC50 at 1 mM ATP (nM)	: 280

FAK(PTK2)

Product code 08-137

Truncated human FAK[376-1052(end) amino acids of accession number NP_722560.1] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-FAK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 47
IC50 at 1 mM ATP (nM)	: 230

FER

Product code 08-139

Full-length human FER [1-822(end) amino acids of accession number NP_005237.1] was expressed as N-terminal GST-fusion protein (122 kDa) using baculovirus expression system. GST-FER was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 12

FES

Product code 08-140

Full-length human FES [1-413, 416-822(end) amino acids of accession number NP_001996.1] was expressed as N-terminal GST-fusion protein (120 kDa) using baculovirus expression system. GST-FES was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 43 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.9
IC50 at 1 mM ATP (nM)	: 25

FGFR1

Product code 08-133

Human FGFR1, cytoplasmic domain [398-822(end) amino acids of accession number NP_075598.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 89 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.3
IC50 at 1 mM ATP (nM)	: 12

FGFR1[V561M]

Product code 08-536

Human FGFR1, cytoplasmic domain [398-822(end) amino acids and V561M of accession number NP_075598.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR1[V561M] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 33 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.14
IC50 at 1 mM ATP (nM)	: 1.3

FGFR2

Product code 08-134

Human FGFR2, cytoplasmic domain [399-821(end) amino acids of accession number NP_000132.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 66 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.4
IC50 at 1 mM ATP (nM)	: 5.4

FGFR2[V564I]

Product code 08-546

Human FGFR2, cytoplasmic domain [399-821(end) amino acids and V564I of accession number NP_000132.3] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-FGFR2[V564I] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.0
IC50 at 1 mM ATP (nM)	: 47

FGFR3

Product code 08-135

Human FGFR3, cytoplasmic domain [436-806(end) amino acids of accession number NP_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-FGFR3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 43 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.6
IC50 at 1 mM ATP (nM)	: 15

FGFR3[K650E]

Product code 08-501

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and K650E of accession number NP_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-FGFR3[K650E] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 41 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 14

FGFR3[K650M]

Product code 08-199

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and K650M of accession number NP_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-FGFR3[K650M] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 17 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.68
IC50 at 1 mM ATP (nM)	: 17

FGFR3[V555L]

Product code 08-548

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and V555L of accession number NP_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-FGFR3[V555L] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 29 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.49
IC50 at 1 mM ATP (nM)	: 9.4

FGFR3[V555M]

Product code 08-543

Human FGFR3, cytoplasmic domain [436-806(end) amino acids and V555M of accession number NP_000133.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-FGFR3[V555M] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.21
IC50 at 1 mM ATP (nM)	: 1.8

FGFR4

Product code 08-136

Human FGFR4, cytoplasmic domain [460-802(end) amino acids of accession number NP_002002.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-FGFR4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 230 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 43
IC50 at 1 mM ATP (nM)	: 120

FGFR4[N535K]

Product code 08-524

Human FGFR4, cytoplasmic domain [460-802(end) amino acids and N535K of accession number NP_002002.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-FGFR4[N535K] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 30 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 160
IC50 at 1 mM ATP (nM)	: 1200

FGFR4[V550E]

Product code 08-525

Human FGFR4, cytoplasmic domain [460-802(end) amino acids and V550E of accession number NP_002002.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-FGFR4[V550E] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 210 / 200
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 370
IC50 at 1 mM ATP (nM)	: 1300

FGFR4[V550L]

Product code 08-526

Human FGFR4, cytoplasmic domain [460-802(end) amino acids and V550L of accession number NP_002002.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-FGFR4[V550L] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 160 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 44

FGR

Product code 08-166

Full-length human FGR [1-529(end) amino acids of accession number NP_005239.1] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system. GST-FGR was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: 16

FLT1

Product code 08-189

Human FLT1, cytoplasmic domain [781-1338(end) amino acids of accession number NP_002010.1] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-FLT1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 140 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.2
IC50 at 1 mM ATP (nM)	: 6.8

FLT3

Product code 08-154

Human FLT3, cytoplasmic domain [564-993(end) amino acids of accession number NP_004110.2] was expressed as N-terminal GST-fusion protein (77 kDa) using baculovirus expression system. GST-FLT3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μM) Km app / Bin : 94 / 100
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 0.20
 IC50 at 1 mM ATP (nM) : 0.34

FLT4

Product code 08-190

Human FLT4, cytoplasmic domain [798-1298(end) amino acids of accession number NP_002011.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-FLT4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : CSK tide
 ATP (μM) Km app / Bin : 72 / 75
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 0.66
 IC50 at 1 mM ATP (nM) : 2.4

FMS(CSF1R)

Product code 08-155

Human FMS, cytoplasmic domain [538-972(end) amino acids of accession number NP_005202.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-FMS was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μM) Km app / Bin : 26 / 25
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 0.26
 IC50 at 1 mM ATP (nM) : 0.70

FRK

Product code 08-167

Human FRK, catalytic domain [223-505(end) amino acids of accession number NP_002022.1] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-FRK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μM) Km app / Bin : 62 / 75
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 3.4
 IC50 at 1 mM ATP (nM) : 40

FYN[isoform a]

Product code 08-168

Full-length human FYN [isoform a] [1-537(end) amino acids of accession number NP_002028.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-FYN [isoform a] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μM) Km app / Bin : 36 / 50
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 3.9
 IC50 at 1 mM ATP (nM) : 24

FYN[isoform b]

Product code 08-531

Full-length human FYN [isoform b] [1-534(end) amino acids of accession number NP_694592.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-FYN [isoform b] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.8
IC50 at 1 mM ATP (nM)	: 42

GSK3 α (GSK3A)

Product code 04-140

Full-length human GSK3 α [1-483(end) amino acids of accession number NP_063937.2] was expressed as N-terminal GST-fusion protein (78 kDa) using baculovirus expression system. GST-GSK3 α was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CREB tide-p
ATP (μ M) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 15
IC50 at 1 mM ATP (nM)	: 400

GSK3 β (GSK3B)

Product code 04-141

Full-length human GSK3 β [1-420(end) amino acids of accession number NP_001139628.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-GSK3 β was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CREB tide-p
ATP (μ M) Km app / Bin	: 9.1 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 9.2
IC50 at 1 mM ATP (nM)	: 240

Haspin(GSG2)

Product code 05-111

Full-length human Haspin [1-798(end) amino acids of accession number NP_114171.2] was expressed as N-terminal GST-fusion protein (116 kDa) using baculovirus expression system. GST-Haspin was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Histone H3 peptide
ATP (μ M) Km app / Bin	: 140 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 5.8
IC50 at 1 mM ATP (nM)	: n.a.

HCK

Product code 08-169

Truncated human HCK [25-526(end) amino acids of accession number NP_002101.2] was expressed as N-terminal GST-fusion protein (84 kDa) using baculovirus expression system. GST-HCK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 22

HER2(ERBB2)

Product code 08-016

Human HER2, cytoplasmic domain [676-1255(end) amino acids of accession number NP_004439.1] was expressed as N-terminal His-tagged protein (67 kDa) using baculovirus expression system. His-tagged HER2 was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 3.5 / 5
Metal	: Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 90
IC50 at 1 mM ATP (nM)	: >10000

HER4(ERBB4)

Product code 08-118

Human HER4, cytoplasmic domain [676-1308(end) amino acids of accession number NP_005226.1] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system. GST-HER4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 34
IC50 at 1 mM ATP (nM)	: 1000

HGK(MAP4K4)

Product code 07-137

Human HGK, catalytic domain [1-328 amino acids of accession number NP_004825.2] was expressed as N-terminal GST-fusion protein (64 kDa) using baculovirus expression system. GST-HGK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Moesin-derived peptide
ATP (μ M) Km app / Bin	: 9.4 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 23

HIPK1

Product code 04-135

Human HIPK1, catalytic domain [158-555 amino acids of accession number NP_689909.2] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-HIPK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRK tide-F
ATP (μ M) Km app / Bin	: 4.4 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 570
IC50 at 1 mM ATP (nM)	: n.a.

HIPK2

Product code 04-136

Full-length human HIPK2 [1-1198(end) amino acids of accession number NP_073577.3] was expressed as N-terminal GST-fusion protein (158 kDa) using baculovirus expression system. GST-HIPK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRK tide-F
ATP (μ M) Km app / Bin	: 5.9 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 170
IC50 at 1 mM ATP (nM)	: n.a.

HIPK3

Product code 04-137

Human HIPK3, catalytic domain [161-562 amino acids of accession number NP_005725.3] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-HIPK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 7.3 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 120
IC50 at 1 mM ATP (nM)	: >10000

HIPK4

Product code 04-138

Full-length human HIPK4 [1-616(end) amino acids of accession number NP_653286.2] was expressed as N-terminal GST-fusion protein (96 kDa) using baculovirus expression system. GST-HIPK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μM) Km app / Bin	: 7 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 71
IC50 at 1 mM ATP (nM)	: >10000

HPK1(MAP4K1)

Product code 07-410

Human HPK1, catalytic domain [1-346 amino acids of accession number NP_009112.1] was expressed as N-terminal DYKDDDDK tagged protein (41 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 22 / 25
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 6.9
IC50 at 1 mM ATP (nM)	: n.a.

IGF1R

Product code 08-141

Human IGF1R, cytoplasmic domain [959-1367(end) amino acids of accession number NP_000866.1] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-IGF1R was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μM) Km app / Bin	: 63 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 40
IC50 at 1 mM ATP (nM)	: 150

IKKα(CHUK)

Product code 05-112

Full-length human IKKα [1-745(end) amino acids of accession number NP_001269.3] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-IKKα was purified by using glutathione sepharose chromatography.

Assay platform	: IMAP
Substrate	: IκBα peptide
ATP (μM) Km app / Bin	: 41 / 40
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 310
IC50 at 1 mM ATP (nM)	: n.a.

IKK β (IKKBK)

Product code 05-084

Truncated human IKK β [1-662 amino acids of accession number NP_001547.1] was expressed as N-terminal His-tagged protein (77 kDa) using baculovirus expression system. His-tagged IKK β was purified by using Ni-NTA affinity chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified I κ B α -derived peptide
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 410
IC50 at 1 mM ATP (nM)	: >10000

IKK ϵ (IKBKE)

Product code 05-114

Full-length human IKK ϵ [1-716(end) amino acids of accession number NP_054721.1] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system. GST-IKK ϵ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: I κ B α peptide
ATP (μ M) Km app / Bin	: 9.5 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.79
IC50 at 1 mM ATP (nM)	: n.a.

INSR

Product code 08-142

Human INSR, catalytic domain [1005-1310 amino acids of accession number NP_000199.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-INSR was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μ M) Km app / Bin	: 58 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: 70

IRAK1

Product code 09-101

Truncated human IRAK1 [194-712(end) amino acids of accession number NP_001560.2] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-IRAK1 was purified by using glutathione sepharose chromatography.

Assay platform	: IMAP
Substrate	: SRPKtide
ATP (μ M) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 54
IC50 at 1 mM ATP (nM)	: n.a.

IRAK4

Product code 09-145

Full-length human IRAK4 [1-460(end) amino acids of accession number NP_057207.2] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-IRAK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRAK1 peptide
ATP (μ M) Km app / Bin	: 920 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 11
IC50 at 1 mM ATP (nM)	: 11

IRR(INSRR)

Product code 08-143

Human IRR, cytoplasmic domain [953-1297(end) amino acids of accession number NP_055030.1] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system. GST-IRR was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : IRS1
 ATP (μ M) Km app / Bin : 64 / 75
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 15
 IC50 at 1 mM ATP (nM) : 98

ITK

Product code 08-181

Full-length human ITK [2-620(end) amino acids of accession number NP_005537.3] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system. GST-ITK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μ M) Km app / Bin : 6.1 / 10
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 3.2
 IC50 at 1 mM ATP (nM) : 200

JAK1

Product code 08-144

Human JAK1, catalytic domain [850-1154(end) amino acids of accession number NP_002218.2] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-JAK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : JAK1 substrate peptide
 ATP (μ M) Km app / Bin : 68 / 75
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 0.71
 IC50 at 1 mM ATP (nM) : 10

JAK2

Product code 08-045

Human JAK2, catalytic domain [826-1132(end) amino acids of accession number NP_004963.1] was expressed as N-terminal His-tagged protein (39 kDa) using baculovirus expression system. His-tagged JAK2 was purified by using Ni-NTA affinity chromatography and gel filtration chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μ M) Km app / Bin : 13 / 10
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 0.34
 IC50 at 1 mM ATP (nM) : 6.0

JAK3

Product code 08-046

Human JAK3, catalytic domain [795-1124(end) amino acids of accession number NP_000206.2] was expressed as N-terminal His-tagged protein (41 kDa) using baculovirus expression system. His-tagged JAK3 was purified by using Ni-NTA affinity chromatography and gel filtration chromatography.

Assay platform : Mobility Shift Assay
 Substrate : Src tide
 ATP (μ M) Km app / Bin : 3.5 / 5
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at ATP Bin (nM) : 0.20
 IC50 at 1 mM ATP (nM) : 12

JNK1(MAPK8)

Product code 04-163

Human JNK1, catalytic domain [2-364 amino acids of accession number NP_620634.1] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-JNK1 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K4 and MAP2K7. Activated GST-JNK1 was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 29 / 100
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 99
IC50 at 1 mM ATP (nM)	: 770

JNK2(MAPK9)

Product code 04-164

Human JNK2, catalytic domain [1-364 amino acids of accession number NP_002743.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-JNK2 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K4 and MAP2K7. Activated GST-JNK2 was purified using Ni-NTA chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 21 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 110
IC50 at 1 mM ATP (nM)	: 1600

JNK3(MAPK10)

Product code 04-150

Full-length human JNK3 [1-426(end) amino acids of accession number NP_620446.1] was expressed as N-terminal GST-fusion protein (75 kDa) using E.coli expression system. GST-JNK3 was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K4 and MAP2K7. Activated GST-JNK3 was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 6 / 25
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 26
IC50 at 1 mM ATP (nM)	: 730

KDR

Product code 08-191

Human KDR, cytoplasmic domain [790-1356(end) amino acids of accession number NP_002244.1] was expressed as N-terminal GST-fusion protein (90 kDa) using baculovirus expression system. GST-KDR was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 74 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.1
IC50 at 1 mM ATP (nM)	: 13

KIT

Product code 08-156

Human KIT, cytoplasmic domain [544-976(end) amino acids of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 370 / 400
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 2.0

KIT[D816E]

Product code 08-541

Human KIT, cytoplasmic domain [544-976(end) amino acids and D816E of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[D816E] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 40 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.7
IC50 at 1 mM ATP (nM)	: 1.3

KIT[D816V]

Product code 08-505

Human KIT, cytoplasmic domain [544-976(end) amino acids and D816V of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[D816V] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.18
IC50 at 1 mM ATP (nM)	: 2.8

KIT[D816Y]

Product code 08-534

Human KIT, cytoplasmic domain [544-976(end) amino acids and D816Y of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[D816Y] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 22 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.27
IC50 at 1 mM ATP (nM)	: 2.1

KIT[T670I]

Product code 08-195

Human KIT, cytoplasmic domain [544-976(end) amino acids and T670I of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[T670I] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 100 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.80
IC50 at 1 mM ATP (nM)	: 3.4

KIT[V560G]

Product code 08-504

Human KIT, cytoplasmic domain [544-976(end) amino acids and V560G of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[V560G] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 110 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 1.6

KIT[V654A]

Product code 08-511

Human KIT, cytoplasmic domain [544-976(end) amino acids and V654A of accession number NP_000213.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-KIT[V654A] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 220 / 250
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP Bin (nM)	: 4.5
IC ₅₀ at 1 mM ATP (nM)	: 8.2

LATS1/MOBKL1A

Product code 01-523

Human LATS1, catalytic domain [589-1130(end) amino acids of accession number NP_004681.1] was co-expressed as N-terminal GST-fusion protein (90 kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP_775739.1] and human BTN-MST2 [1-491(end) amino acids of accession number NP_006272.2] using baculovirus expression system. GST-LATS1/MOBKL1A was purified by using glutathione sepharose chromatography and DYKDDDDK tag antibody agarose.

Assay platform	: Mobility Shift Assay
Substrate	: SGK tide
ATP (μ M) Km app / Bin	: 23 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP Bin (nM)	: 4.2
IC ₅₀ at 1 mM ATP (nM)	: 11

LATS2/MOBKL1A

Product code 01-524

Human LATS2, catalytic domain [553-1088(end) amino acids of accession number NP_055387.2] was co-expressed as N-terminal GST-fusion protein (89 kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP_775739.1] and human BTN-MST2 [1-491(end) amino acids of accession number NP_006272.2] using baculovirus expression system. GST-LATS2/MOBKL1A was purified by using glutathione sepharose chromatography and DYKDDDDK tag antibody agarose.

Assay platform	: Mobility Shift Assay
Substrate	: SGK tide
ATP (μ M) Km app / Bin	: 38 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP Bin (nM)	: 1.7
IC ₅₀ at 1 mM ATP (nM)	: 4.0

LCK

Product code 08-170

Full-length human LCK [1-509(end) amino acids of accession number NP_005347.2] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-LCK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP Bin (nM)	: 1.5
IC ₅₀ at 1 mM ATP (nM)	: 14

LOK(STK10)

Product code 07-315

Full-length human LOK [1-968(end) amino acids of accession number BAA35073.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-LOK was purified by using glutathione sepharose chromatography. GST-LOK was cleaved by PreScission protease and GST-free LOK (114 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Moesin-derived peptide
ATP (μ M) Km app / Bin	: 100 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC ₅₀ at ATP Bin (nM)	: 0.49
IC ₅₀ at 1 mM ATP (nM)	: n.a.

LTK

Product code 08-106

Human LTK, catalytic domain [498-796 amino acids of accession number NP_002335.2] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-LTK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 49 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.0
IC50 at 1 mM ATP (nM)	: 7.1

LYNa

Product code 08-171

Full-length human LYNa [1-512(end) amino acids of accession number NP_002341.1] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system. GST-LYNa was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.0
IC50 at 1 mM ATP (nM)	: 22

LYNb

Product code 08-172

Full-length human LYNb [1-491(end) amino acids of accession number NP_001104567.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-LYNb was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 18 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.0
IC50 at 1 mM ATP (nM)	: 21

MAP4K2

Product code 07-111

Full-length human MAP4K2 [1-820(end) amino acid of accession number NP_004570.2] was expressed as N-terminal GST-fusion protein (119 kDa) using baculovirus expression system. GST-MAP4K2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 93 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: n.a.

MAPKAPK2

Product code 02-142

Full-length human MAPKAPK2 [1-400(end) amino acids of accession number NP_116584.2] was co-expressed as N-terminal GST-fusion protein (73 kDa) with human His-tagged p38β [1-364(end) amino acids of accession number NP_002742.3] and human His-tagged MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] using baculovirus expression system. GST-MAPKAPK2 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 3.6 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 80
IC50 at 1 mM ATP (nM)	: 9300

MAPKAPK3

Product code 02-143

Full-length human MAPKAPK3 [1-382(end) amino acids of accession number NP_004626.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with human His-tagged p38 β [1-364(end) amino acids of accession number NP_002742.3] and human His-tagged MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] using baculovirus expression system. GST-MAPKAPK3 was purified by using glutathione sepharose chromatography and Ni-NTA chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 410
IC50 at 1 mM ATP (nM)	: n.a.

MAPKAPK5

Product code 02-144

Full-length human MAPKAPK5 [1-471(end) amino acids of accession number NP_003659.2] was co-expressed as N-terminal GST-fusion protein (81 kDa) with human His-tagged p38 β [1-364(end) amino acids of accession number NP_002742.3] and human His-tagged MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] using baculovirus expression system. GST-MAPKAPK5 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 320
IC50 at 1 mM ATP (nM)	: n.a.

MARK1

Product code 02-120

Full-length human MARK1 [1-795(end) amino acids of accession number AAF72103.1] was expressed as N-terminal GST-fusion protein (116 kDa) using baculovirus expression system. GST-MARK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.17
IC50 at 1 mM ATP (nM)	: n.a.

MARK2

Product code 02-121

Full-length human MARK2 [1-745(end) amino acids of accession number NP_059672.2] was expressed as N-terminal GST-fusion protein (110 kDa) using baculovirus expression system. GST-MARK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 8.8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.12
IC50 at 1 mM ATP (nM)	: n.a.

MARK3

Product code 02-122

Full-length human MARK3 [1-729(end) amino acids of accession number NP_002367.4] was expressed as N-terminal GST-fusion protein (108 kDa) using baculovirus expression system. GST-MARK3 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.20
IC50 at 1 mM ATP (nM)	: n.a.

MARK4

Product code 02-123

Full-length human MARK4 [1-688(end) amino acids of accession number NP_113605.2] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-MARK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μM) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.12
IC50 at 1 mM ATP (nM)	: 5.6

MELK

Product code 02-124

Truncated human MELK [1-493 amino acids of accession number NP_055606.1] was expressed as N-terminal GST-fusion protein (83 kDa) using E. coli expression system. GST-MELK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 38 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.81
IC50 at 1 mM ATP (nM)	: n.a.

MER(MERTK)

Product code 08-108

Human MER, cytoplasmic domain [528-999(end) amino acids of accession number NP_006334.2] was expressed as N-terminal GST-fusion protein (80 kDa) using baculovirus expression system. GST-MER was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 36 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.61
IC50 at 1 mM ATP (nM)	: 5.3

MET

Product code 08-151

Human MET, cytoplasmic domain [956-1390(end) amino acids of accession number NP_000236.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-MET was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 67
IC50 at 1 mM ATP (nM)	: 730

MET[D1228H]

Product code 08-540

Human MET, cytoplasmic domain [956-1390(end) amino acids and D1228H of accession number NP_000236.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-MET[D1228H] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 59
IC50 at 1 mM ATP (nM)	: 1200

MET[M1250T]

Product code 08-545

Human MET, cytoplasmic domain [956-1390(end) amino acids and M1250T of accession number NP_000236.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-MET[M1250T] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 17 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 86
IC50 at 1 mM ATP (nM)	: 1900

MET[Y1235D]

Product code 08-198

Human MET, cytoplasmic domain [956-1390(end) amino acids and Y1235D of accession number NP_000236.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-MET[Y1235D] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 71 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 79
IC50 at 1 mM ATP (nM)	: 390

MINK(MINK1)

Product code 07-139

Human MINK, catalytic domain [1-314 amino acids of accession number NP_056531.1] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-MINK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erk tide
ATP (μM) Km app / Bin	: 16 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 5.5
IC50 at 1 mM ATP (nM)	: 4.7

MNK1(MKNK1)

Product code 02-145

Full-length human MNK1 [1-424(end) amino acids and T344D of accession number NP_001129025.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-MNK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: RS peptide
ATP (μM) Km app / Bin	: 460 / 450
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 21
IC50 at 1 mM ATP (nM)	: n.a.

MNK2(MKNK2)

Product code 02-146

Full-length human MNK2 [1-465(end) amino acids and T379D of accession number NP_951009.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-MNK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: RS peptide
ATP (μM) Km app / Bin	: 110 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 7.5
IC50 at 1 mM ATP (nM)	: 44

MRCK α (CDC42BPA)

Product code 01-107

Truncated human MRCK α [1-574 amino acids of accession number NP_003598.2] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system. GST-MRCK α was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DAPK1tide
ATP (μ M) Km app / Bin	: 0.45 / 1
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: n.a.

MRCK β (CDC42BPB)

Product code 01-108

Truncated human MRCK β [1-473 amino acids of accession number NP_006026.3] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system. GST-MRCK β was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DAPK1tide
ATP (μ M) Km app / Bin	: 0.67 / 1
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.6
IC50 at 1 mM ATP (nM)	: n.a.

MSK1(RPS6KA5)

Product code 01-147

Full-length human MSK1 [1-802(end) amino acids of accession number NP_004746.2] was co-expressed as N-terminal GST-fusion protein (117 kDa) with human His-tagged Erk2 [1-360 amino acids of accession number NP_002736.3] using baculovirus expression system. GST-MSK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Crosstide
ATP (μ M) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.6
IC50 at 1 mM ATP (nM)	: 30

MSK2(RPS6KA4)

Product code 01-148

Full-length human MSK2 [1-772(end) amino acids of accession number NP_003933.1] was co-expressed as N-terminal GST-fusion protein (114 kDa) with human His-tagged Erk2 [1-360 amino acids of accession number NP_002736.3] using baculovirus expression system. GST-MSK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Crosstide
ATP (μ M) Km app / Bin	: 40 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.3
IC50 at 1 mM ATP (nM)	: 20

MSSK1(STK23)

Product code 04-159

Full-length human MSSK1 [1-567(end) amino acids of accession number NP_055185.2] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-MSSK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 220
IC50 at 1 mM ATP (nM)	: n.a.

MST1(STK4)

Product code 07-116

Full-length human MST1 [1-487(end) amino acids of accession number NP_006273.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-MST1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μM) Km app / Bin	: 50 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 0.55

MST2(STK3)

Product code 07-117

Full-length human MST2 [1-491(end) amino acids of accession number NP_006272.2] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-MST2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μM) Km app / Bin	: 69 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 6.7
IC50 at 1 mM ATP (nM)	: 3.1

MST3(STK24)

Product code 07-118

Full-length human MST3 [1-431(end) amino acids of accession number NP_001027467.2] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-MST3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Moesin-derived peptide
ATP (μM) Km app / Bin	: 66 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.9
IC50 at 1 mM ATP (nM)	: n.a.

MST4

Product code 07-119

Full-length human MST4 [1-416(end) amino acids of accession number NP_057626.2] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-MST4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Moesin-derived peptide
ATP (μM) Km app / Bin	: 76 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 6.3
IC50 at 1 mM ATP (nM)	: n.a.

MUSK

Product code 08-153

Human MUSK, catalytic domain [527-869(end) amino acids of accession number NP_005583.1] was expressed as N-terminal GST fusion protein (66 kDa) using baculovirus expression system. GST-MUSK was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 14 / 10
Metal	: Mg+Mn
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.1
IC50 at 1 mM ATP (nM)	: 2.6

NDR1(STK38)

Product code 01-125

Full-length human NDR1[1-465(end) amino acids of accession number NP_009202.1] was co-expressed as N-terminal GST-fusion protein (81kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP_775739.1] using baculovirus expression system. GST-NDR1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.8
IC50 at 1 mM ATP (nM)	: n.a.

NDR2(STK38L)

Product code 01-126

Full-length human NDR2 [1-464(end) amino acids of accession number NP_055815.1] was co-expressed as N-terminal GST-fusion protein (81 kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP_775739.1] using baculovirus expression system. GST-NDR2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 7.6 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.1
IC50 at 1 mM ATP (nM)	: n.a.

NEK1

Product code 05-123

Human NEK1, catalytic domain [1-505 amino acids of accession number NP_036356.1] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-NEK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDK7 peptide
ATP (μM) Km app / Bin	: 64 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 51
IC50 at 1 mM ATP (nM)	: 650

NEK2

Product code 05-226

Full-length human NEK2 [1-445(end) amino acids of accession number NP_002488.1] was expressed as N-terminal His-tagged protein (55 kDa) using baculovirus expression system. His-tagged NEK2 was purified by using Ni-NTA affinity chromatography. Purified His-NEK2 was digested by recombinant His-TEV protease, and His-tag free NEK2 (ca. 52 kDa) was collected as flow-through fraction from Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDK7 peptide
ATP (μM) Km app / Bin	: 65 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3700
IC50 at 1 mM ATP (nM)	: >10000

NEK4

Product code 05-128

Full-length human NEK4 [1-841(end) amino acids of accession number NP_003148.2] was expressed as N-terminal GST-fusion protein (122 kDa) using baculovirus expression system. GST-NEK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 51 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 120
IC50 at 1 mM ATP (nM)	: n.a.

NEK6

Product code 05-130

Full-length human NEK6 [1-313(end) amino acids of accession number NP_055212.2] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-NEK6 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDK7 peptide
ATP (μM) Km app / Bin	: 69 / 75
Metal	: Mg
Reference compound	: PKR Inhibitor
IC50 at ATP Bin (nM)	: 19000
IC50 at 1 mM ATP (nM)	: >10000

NEK7

Product code 05-131

Full-length human NEK7 [1-302(end) amino acids of accession number NP_598001.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-NEK7 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDK7 peptide
ATP (μM) Km app / Bin	: 40 / 50
Metal	: Mg
Reference compound	: PKR Inhibitor
IC50 at ATP Bin (nM)	: 8500
IC50 at 1 mM ATP (nM)	: >10000

NEK9

Product code 05-133

Truncated human NEK9 [1-346, 733-979(end) amino acids of accession number NP_149107.4] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system. GST-NEK9 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDK7 peptide
ATP (μM) Km app / Bin	: 190 / 200
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 150
IC50 at 1 mM ATP (nM)	: 400

NIM1K(MGC42105)

Product code 02-175

Full-length human NIM1K [1-436(end) amino acids of accession number NP_699192.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-NIM1K was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μM) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 300
IC50 at 1 mM ATP (nM)	: n.a.

Nuak1

Product code 02-126

Full-length human NuaK1 [1-661(end) amino acids of accession number NP_055655.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system. GST-NuaK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μM) Km app / Bin	: 59 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 3.4

NuaK2

Product code 02-127

Full-length human NuaK2 [1-628(end) amino acids of accession number NP_112214.1] was expressed as N-terminal GST-fusion protein (98kDa) using baculovirus expression system. GST-NuaK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CHKtide
ATP (μM) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 2.3

p38α(MAPK14)

Product code 04-152

Truncated human p38α [9-352 amino acids of accession number NP_620581.1] was expressed as N-terminal GST-fusion protein (66 kDa) using E. coli expression system. GST-p38α was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38α was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 150 / 150
Metal	: Mg
Reference compound	: SB202190
IC50 at ATP Bin (nM)	: 6.3
IC50 at 1 mM ATP (nM)	: 22

p38β(MAPK11)

Product code 04-153

Full-length human p38β [1-364(end) amino acids of accession number NP_002742.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38β was purified by using glutathione chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 63 / 75
Metal	: Mg
Reference compound	: SB202190
IC50 at ATP Bin (nM)	: 16
IC50 at 1 mM ATP (nM)	: 110

p38γ(MAPK12)

Product code 04-155

Full-length human p38γ [1-367(end) amino acids of accession number NP_002960.2] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38γ was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38γ was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 88
IC50 at 1 mM ATP (nM)	: 2800

p38δ(MAPK13)

Product code 04-154

Full-length human p38δ [1-365(end) amino acids of accession number NP_002745.1] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38δ was purified by using glutathione sepharose chromatography and activated with His-tagged MAP2K6. Activated GST-p38δ was purified using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Modified Erktide
ATP (μM) Km app / Bin	: 5.8 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 220
IC50 at 1 mM ATP (nM)	: >10000

p70S6K(RPS6KB1)

Product code 01-154

Human p70S6K, catalytic domain [1-421 amino acids and T412E of accession number NP_003152.1] was expressed as N-terminal GST-fusion protein (75 kDa) using baculovirus expression system. GST-p70S6K was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.1
IC50 at 1 mM ATP (nM)	: 9.8

p70S6Kβ(RPS6KB2)

Product code 01-155

Full-length human p70S6Kβ [1-482(end) amino acids of accession number NP_003943.2] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-p70S6Kβ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 3.3 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.7
IC50 at 1 mM ATP (nM)	: n.a.

PAK1

Product code 07-123

Full-length human PAK1 [1-545(end) amino acids of accession number NP_002567.3] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PAK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: LIMKtide
ATP (μM) Km app / Bin	: 300 / 300
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.0
IC50 at 1 mM ATP (nM)	: 11

PAK2

Product code 07-124

Full-length human PAK2 [1-524(end) amino acids of accession number NP_002568.2] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-PAK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DAPK1tide
ATP (μM) Km app / Bin	: 81 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.5
IC50 at 1 mM ATP (nM)	: 22

PAK4

Product code 07-126

Full-length human PAK4 [1-591(end) amino acids of accession number NP_005875.1] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system. GST-PAK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 2.5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: n.a.

PAK5(PAK7)

Product code 07-127

Human PAK5, catalytic domain [425-719(end) amino acids of accession number NP_065074.1] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-PAK5 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DAPK1tide
ATP (μM) Km app / Bin	: 1.9 / 1
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.5
IC50 at 1 mM ATP (nM)	: 290

PAK6

Product code 07-128

Full-length human PAK6 [1-681(end) amino acids of accession number NP_064553.1] was expressed as N-terminal GST-fusion protein (102 kDa) using baculovirus expression system. GST-PAK6 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 3.7 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: n.a.

PASK

Product code 02-128

Human PASK, catalytic domain [949-1323(end) amino acids of accession number NP_055963.2] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-PASK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 9.7 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 13
IC50 at 1 mM ATP (nM)	: 190

PBK

Product code 05-168

Full-length human PBK [1-322(end) amino acids of accession number NP_060962.2] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-PBK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Histone H3 peptide
ATP (μM) Km app / Bin	: 33 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 69
IC50 at 1 mM ATP (nM)	: 720

PDGFRα(PDGFRα)

Product code 08-157

Human PDGFRα, cytoplasmic domain [550-1089(end) amino acids of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-PDGFRα was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 28 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.30
IC50 at 1 mM ATP (nM)	: 1.4

PDGFR α (PDGFRA)[D842V]

Product code 08-506

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids and D842V of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-PDGFR α [D842V] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.25
IC50 at 1 mM ATP (nM)	: 1.9

PDGFR α (PDGFRA)[T674I]

Product code 08-503

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids and T674I of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-PDGFR α [T674I] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.12
IC50 at 1 mM ATP (nM)	: 1.1

PDGFR α (PDGFRA)[V561D]

Product code 08-507

Human PDGFR α , cytoplasmic domain [550-1089(end) amino acids and V561D of accession number NP_006197.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-PDGFR α [V561D] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 35 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.32
IC50 at 1 mM ATP (nM)	: 1.6

PDGFR β (PDGFRB)

Product code 08-158

Human PDGFR β , cytoplasmic domain [557-1106(end) amino acids of accession number NP_002600.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PDGFR β was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 23 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.27
IC50 at 1 mM ATP (nM)	: 0.62

PDHK2(PDK2)

Product code 10-140

Full-length human PDHK2 [1-407(end) amino acids of accession number NP_002602.2] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-PDHK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PDHKtide
ATP (μ M) Km app / Bin	: 28 / 25
Metal	: Mg+K
Reference compound	: VER-246608
IC50 at ATP Bin (nM)	: 33
IC50 at 1 mM ATP (nM)	: n.a.

PDHK4(PDK4)

Product code 10-125

Full-length human PDHK4 [1-411(end) amino acids of accession number NP_002603.1] was expressed as N-terminal GST-fusion protein (73 kDa) using E.coli expression system. GST-PDHK4 was purified by using glutathione affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PDHKtide
ATP (μM) Km app / Bin	: 19 / 25
Metal	: Mg+K
Reference compound	: VER-246608
IC50 at ATP Bin (nM)	: 120
IC50 at 1 mM ATP (nM)	: n.a.

PDK1(PDPK1)

Product code 01-132

Full-length human PDK1 [1-556(end) amino acids of accession number NP_002604.1] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system. GST-PDK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: T308tide
ATP (μM) Km app / Bin	: 9.6 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 9.2
IC50 at 1 mM ATP (nM)	: 12

PEK(EIF2AK3)

Product code 05-155

Human PEK, cytoplasmic domain [536-1116(end) amino acids of accession number NP_004827.3] was expressed as N-terminal GST-fusion protein (94 kDa) using E.coli expression system. GST-PEK was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: IMAP
Substrate	: SRPKtide
ATP (μM) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3600
IC50 at 1 mM ATP (nM)	: n.a.

PGK(PRKG1)

Product code 01-142

Full-length human PGK [1-686(end) amino acids of accession number NP_006249.1] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-PGK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μM) Km app / Bin	: 8.2 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.1
IC50 at 1 mM ATP (nM)	: n.a.

PHKG1

Product code 02-152

Full-length human PHKG1 [1-387(end) amino acids of accession number NP_006204.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-PHKG1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 71 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.22
IC50 at 1 mM ATP (nM)	: n.a.

PHKG2

Product code 02-153

Full-length human PHKG2 [1-406(end) amino acids of accession number NP_000285.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-PHKG2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 8.1 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.74
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CA/PIK3R1

Product code 11-401-20N

Full-length human PIK3CA[1-1068(end) amino acids of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 89 / 100
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 22
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CA[E542K]/PIK3R1

Product code 11-413-20N

Full-length human PIK3CA[1-1068(end) amino acids and E542K of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 42 / 50
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CA[E545K]/PIK3R1

Product code 11-414-20N

Full-length human PIK3CA[1-1068(end) amino acids and E545K of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 44 / 50
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 11
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CA[H1047R]/PIK3R1

Product code 11-415-20N

Full-length human PIK3CA[1-1068(end) amino acids and H1047R of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 78 / 75
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CA[P539R]/PIK3R1

Product code 11-412-20N

Full-length human PIK3CA[1-1068(end) amino acids and P539R of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 8.7
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CA[R88Q]/PIK3R1

Product code 11-411-20N

Full-length human PIK3CA[1-1068(end) amino acids and R88Q of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 42 / 50
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 16
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CB/PIK3R1

Product code 11-402-20N

Full-length human PIK3CB[1-1070(end) amino acids of accession number NP_006210.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (126 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] (84kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 88 / 100
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 22
IC50 at 1 mM ATP (nM)	: n.a.

PIK3CD/PIK3R1

Product code 11-403-20N

Full-length human PIK3CD[1-1044(end) amino acids of accession number NP_005017.3] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (123 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] (84kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform	: ADP-Glo
Substrate	: PI(4,5)P2
ATP (μM) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: PI-103
IC50 at ATP Bin (nM)	: 24
IC50 at 1 mM ATP (nM)	: n.a.

PIKFYVE(PIP5K3)

Product code 11-118

Full-length human PIKFYVE [1-2098(end) amino acids and S696N, L932S, Q995L, T998S, S1033A and Q1183K of accession number NP_055855.2] was expressed as N-terminal GST-fusion protein (265 kDa) using baculovirus expression system. GST-PIKFYVE was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(3)P
ATP (μM) Km app / Bin	: 36 / 50
Metal	: Mg
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 3900
IC50 at 1 mM ATP (nM)	: n.a.

PIM1

Product code 02-054

Full-length human PIM1 [1-313(end) amino acids of accession number NP_002639.1] was expressed as N-terminal His-tagged protein (39 kDa) using baculovirus expression system. His-tagged PIM1 was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 640 / 500
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 20

PIM2

Product code 02-155

Full-length human PIM2 [1-311(end) amino acids of accession number NP_006866.2] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system. GST-PIM2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 4 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 14
IC50 at 1 mM ATP (nM)	: 480

PIM3

Product code 02-156

Full-length human PIM3 [1-326(end) amino acids of accession number NP_001001852.1] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-PIM3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K2 peptide
ATP (μM) Km app / Bin	: 130 / 150
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.36
IC50 at 1 mM ATP (nM)	: 0.71

PIP4K2A

Product code 11-115

Full-length human PIP4K2A [1-406(end) amino acids of accession number NP_005019] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-PIP4K2A was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(5)P
ATP (μM) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 7600
IC50 at 1 mM ATP (nM)	: n.a.

PIP4K2B

Product code 11-116

Full-length human PIP4K2B [1-416(end) amino acids of accession number NP_003550] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system. GST-PIP4K2B was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(5)P
ATP (μM) Km app / Bin	: 18 / 25
Metal	: Mn
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 48000
IC50 at 1 mM ATP (nM)	: n.a.

PIP5K1A

Product code 11-111

Full-length human PIP5K1A [1-549(end) amino acids of accession number NP_003548.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PIP5K1A was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(4)P
ATP (μM) Km app / Bin	: 28 / 25
Metal	: Mg
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 10000
IC50 at 1 mM ATP (nM)	: n.a.

PIP5K1B

Product code 11-112

Full-length human PIP5K1B [1-540(end) amino acids of accession number NP_003549.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PIP5K1B was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(4)P
ATP (μM) Km app / Bin	: 95 / 100
Metal	: Mg
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 4300
IC50 at 1 mM ATP (nM)	: n.a.

PIP5K1C

Product code 11-113

Full-length human PIP5K1C [1-668(end) amino acids of accession number NP_036530] was expressed as N-terminal GST-fusion protein (101 kDa) using baculovirus expression system. GST-PIP5K1C was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(4)P
ATP (μM) Km app / Bin	: 33 / 50
Metal	: Mg
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 1900
IC50 at 1 mM ATP (nM)	: n.a.

PIP5KL1

Product code 11-114

Full-length human PIP5KL1 [1-394(end) amino acids of accession number NP_001128691.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-PIP5KL1 was purified by using glutathione sepharose chromatography.

Assay platform	: ADP-Glo
Substrate	: PI(4)P
ATP (μM) Km app / Bin	: 1 / 1
Metal	: Mg
Reference compound	: AG-183
IC50 at ATP Bin (nM)	: 2200
IC50 at 1 mM ATP (nM)	: n.a.

PKACα(PRKACA)

Product code 01-127

Full-length human PKACα [1-351(end) amino acids of accession number NP_002721.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-PKACα was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μM) Km app / Bin	: 2.6 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.80
IC50 at 1 mM ATP (nM)	: 86

PKAC β (PRKACB)

Product code 01-128

Full-length human PKAC β [1-351(end) amino acids of accession number NP_002722.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-PKAC β was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 4.7 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 37

PKAC γ (PRKACG)

Product code 01-129

Full-length human PKAC γ [1-351(end) amino acids of accession number NP_002723.2] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-PKAC γ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μ M) Km app / Bin	: 4.5 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3.1
IC50 at 1 mM ATP (nM)	: n.a.

PKC α (PRKCA)

Product code 01-133

Full-length human PKC α [1-672(end) amino acids of accession number NP_002728.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-PKC α was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 36 / 50
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.33
IC50 at 1 mM ATP (nM)	: 3.6

PKC β 1(PRKCB1)

Product code 01-134

Full-length human PKC β 1 [1-671(end) amino acids of accession number NP_997700.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-PKC β 1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 79 / 75
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.71
IC50 at 1 mM ATP (nM)	: n.a.

PKC β 2(PRKCB2)

Product code 01-165

Full-length human PKC β 2 [1-673(end) amino acids of accession number NP_002729.2] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-PKC β 2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 41 / 50
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.43
IC50 at 1 mM ATP (nM)	: n.a.

PKC γ (PRKCG)

Product code 01-137

Full-length human PKC γ [1-697(end) amino acids of accession number NP_002730.1] was expressed as N-terminal GST-fusion protein (106 kDa) using baculovirus expression system. GST-PKC γ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 74 / 75
Metal	: Mg+Ca
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 11

PKC δ (PRKCD)

Product code 01-135

Full-length human PKC δ [1-676(end) amino acids of accession number NP_006245.2] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-PKC δ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.22
IC50 at 1 mM ATP (nM)	: n.a.

PKC ϵ (PRKCE)

Product code 01-136

Full-length human PKC ϵ [1-737(end) amino acids of accession number NP_005391.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-PKC ϵ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.56
IC50 at 1 mM ATP (nM)	: 5.6

PKC ζ (PRKCZ)

Product code 01-141

Full-length human PKC ζ [1-592(end) amino acids of accession number NP_002735.3] was expressed as N-terminal GST-fusion protein (94kDa) using baculovirus expression system. GST-PKC ζ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 55
IC50 at 1 mM ATP (nM)	: n.a.

PKC η (PRKCH)

Product code 01-138

Full-length human PKC η [1-683(end) amino acids of accession number NP_006246.2] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-PKC η was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μ M) Km app / Bin	: 36 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.92
IC50 at 1 mM ATP (nM)	: n.a.

PKCθ(PRKCQ)

Product code 01-140

Full-length human PKCθ [1-706(end) amino acids of accession number NP_006248.1] was expressed as N-terminal GST-fusion protein (109 kDa) using baculovirus expression system. GST-PKCθ was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μM) Km app / Bin	: 18 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.72
IC50 at 1 mM ATP (nM)	: n.a.

PKCι(PRKCI)

Product code 01-139

Full-length human PKCι[1-587(end) amino acids of accession number NP_002731.3] was expressed as N-terminal GST-fusion protein (94 kDa) using baculovirus expression system. GST-PKCι was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: PKC peptide
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 13
IC50 at 1 mM ATP (nM)	: n.a.

PKD1(PRKD1)

Product code 02-157

Full-length human PKD1 [1-912(end) amino acids of accession number NP_002733.1] was expressed as N-terminal GST-fusion protein (129 kDa) using baculovirus expression system. GST-PKD1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 25 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: n.a.

PKD2(PRKD2)

Product code 02-158

Full-length human PKD2 [1-878(end) amino acids of accession number NP_057541.2] was expressed as N-terminal GST-fusion protein (124 kDa) using baculovirus expression system. GST-PKD2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 26 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 16

PKD3(PRKD3)

Product code 02-159

Full-length human PKD3 [1-890(end) amino acids of accession number NP_005804.1] was expressed as N-terminal GST-fusion protein (127 kDa) using baculovirus expression system. GST-PKD3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 34 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.80
IC50 at 1 mM ATP (nM)	: n.a.

PKN1

Product code 01-144

Full-length human PKN1 [1-942(end) amino acids of accession number NP_002732.3] was expressed as N-terminal GST-fusion protein (132 kDa) using baculovirus expression system. GST-PKN1 was purified by using glutathione sepharose chromatography.

Assay platform	: IMAP
Substrate	: S6K peptide
ATP (μ M) Km app / Bin	: 19 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.15
IC50 at 1 mM ATP (nM)	: n.a.

PKR(EIF2AK2)

Product code 05-156

Human PKR, catalytic domain [252-551(end) amino acids of accession number NP_002750.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-PKR was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: IMAP
Substrate	: SRPKtide
ATP (μ M) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 87
IC50 at 1 mM ATP (nM)	: n.a.

PLK1

Product code 05-157

Full-length human PLK1 [1-603(end) amino acids of accession number NP_005021.2] was expressed as N-terminal GST-fusion protein (95 kDa) using baculovirus expression system. GST-PLK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDC25ctide
ATP (μ M) Km app / Bin	: 5.6 / 5
Metal	: Mg
Reference compound	: GW843682X
IC50 at ATP Bin (nM)	: 3.6
IC50 at 1 mM ATP (nM)	: 47

PLK2

Product code 05-158

Full-length human PLK2 [1-685(end) amino acids of accession number NP_006613.2] was expressed as N-terminal GST-fusion protein (105 kDa) using baculovirus expression system. GST-PLK2 was purified by using glutathione sepharose chromatography.

Assay platform	: IMAP
Substrate	: CHK2 peptide
ATP (μ M) Km app / Bin	: 30 / 30
Metal	: Mg
Reference compound	: GW843682X
IC50 at ATP Bin (nM)	: 4.8
IC50 at 1 mM ATP (nM)	: n.a.

PLK3

Product code 05-159

Human PLK3, catalytic domain [58-340 amino acids of accession number NP_004064.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system. GST-PLK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CDC25ctide
ATP (μ M) Km app / Bin	: 6.8 / 5
Metal	: Mg
Reference compound	: GW843682X
IC50 at ATP Bin (nM)	: 33
IC50 at 1 mM ATP (nM)	: 450

PRKX

Product code 01-130

Full-length human PRKX [1-358(end) amino acids of accession number NP_005035.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-PRKX was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Kemptide
ATP (μM) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.59
IC50 at 1 mM ATP (nM)	: n.a.

PYK2(PTK2B)

Product code 08-138

Full-length human PYK2 [1-967(end) amino acids of accession number NP_775267.1] was expressed as N-terminal GST-fusion protein (138 kDa) using baculovirus expression system. GST-PYK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 56 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.2
IC50 at 1 mM ATP (nM)	: 4.9

QIK(SNF1LK2)

Product code 02-129

Full-length human QIK(SNF1LK2) [1-926(end) amino acids of accession number NP_056006.1] was expressed as N-terminal GST-fusion protein (132 kDa) using baculovirus expression system. GST-QIK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: AMARA peptide
ATP (μM) Km app / Bin	: 42 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.4
IC50 at 1 mM ATP (nM)	: 2.9

RET

Product code 08-159

Human RET, cytoplasmic domain [658-1114(end) amino acids of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 7.5 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: 20

RET[G691S]

Product code 08-522

Human RET, cytoplasmic domain [658-1114(end) amino acids and G691S of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[G691S] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.1
IC50 at 1 mM ATP (nM)	: 24

RET[M918T]

Product code 08-508

Human RET, cytoplasmic domain [658-1114(end) amino acids and M918T of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[M918T] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 4.2 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.4
IC50 at 1 mM ATP (nM)	: 81

RET[S891A]

Product code 08-523

Human RET, cytoplasmic domain [658-1114(end) amino acids and S891A of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[S891A] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.44
IC50 at 1 mM ATP (nM)	: 9.6

RET[Y791F]

Product code 08-521

Human RET, cytoplasmic domain [658-1114(end) amino acids and Y791F of accession number NP_066124.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-RET[Y791F] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μM) Km app / Bin	: 29 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.5
IC50 at 1 mM ATP (nM)	: 26

ROCK1

Product code 01-109

Human ROCK1, catalytic domain [1-477 amino acids of accession number NP_005397.1] was expressed as N-terminal GST-fusion protein (82 kDa) using baculovirus expression system. GST-ROCK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: LIMKtide
ATP (μM) Km app / Bin	: 3.1 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.6
IC50 at 1 mM ATP (nM)	: 73

ROCK2

Product code 01-110

Human ROCK2, catalytic domain [1-553 amino acids of accession number NP_004841.2] was expressed as N-terminal GST-fusion protein (91 kDa) using baculovirus expression system. GST-ROCK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: LIMKtide
ATP (μM) Km app / Bin	: 7.4 / 5
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.92
IC50 at 1 mM ATP (nM)	: 21

RON(MST1R)

Product code 08-152

Human RON, cytoplasmic domain [979-1400(end) amino acids of accession number NP_002438.1] was expressed as N-terminal GST-fusion protein (75kDa) using baculovirus expression system. GST-RON was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 27 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 37
IC50 at 1 mM ATP (nM)	: 550

ROS(ROS1)

Product code 08-163

Human ROS, cytoplasmic domain [1883-2347(end) amino acids of accession number NP_002935.2] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-ROS was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: IRS1
ATP (μM) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.4
IC50 at 1 mM ATP (nM)	: 1.0

RSK1(RPS6KA1)

Product code 01-149

Full-length human RSK1 [1-735(end) amino acids of accession number NP_002944.2] was expressed as N-terminal GST-fusion protein (110 kDa) using baculovirus expression system. GST-RSK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K peptide(N-FL)
ATP (μM) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.14
IC50 at 1 mM ATP (nM)	: 2.5

RSK2(RPS6KA3)

Product code 01-150

Full-length human RSK2 [1-740(end) amino acids of accession number NP_004577.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-RSK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K peptide(N-FL)
ATP (μM) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.10
IC50 at 1 mM ATP (nM)	: 2.1

RSK3(RPS6KA2)

Product code 01-151

Full-length human RSK3 [1-733(end) amino acids of accession number NP_066958.2] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-RSK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K peptide(N-FL)
ATP (μM) Km app / Bin	: 9.9 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.11
IC50 at 1 mM ATP (nM)	: 1.7

RSK4(RPS6KA6)

Product code 01-152

Full-length human RSK4 [1-745(end) amino acids of accession number NP_055311.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-RSK4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: S6K peptide(N-FL)
ATP (μM) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.051
IC50 at 1 mM ATP (nM)	: 0.56

SGK

Product code 01-158

Truncated human SGK [61-431(end) amino acids and S422D of accession number NP_005618.2] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-SGK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 52 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 10
IC50 at 1 mM ATP (nM)	: 99

SGK2

Product code 01-159

Full-length human SGK2 [1-367(end) amino acids and S356D of accession number NP_733794.1] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-SGK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 58 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 30
IC50 at 1 mM ATP (nM)	: n.a.

SGK3(SGKL)

Product code 01-160

Truncated human SGK3 [119-496(end) amino acids and S486D of accession number NP_037389.4] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged PDK1 [1-556(end) amino acids of accession number NP_002604.1] using baculovirus expression system. GST-SGK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SGKtide
ATP (μM) Km app / Bin	: 17 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 42
IC50 at 1 mM ATP (nM)	: n.a.

SIK(SNF1LK)

Product code 02-131

Full-length human SIK [1-783(end) amino acids of accession number NP_775490.2] was expressed as N-terminal GST-fusion protein (112 kDa) using baculovirus expression system. GST-SIK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: AMARA peptide
ATP (μM) Km app / Bin	: 47 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.8
IC50 at 1 mM ATP (nM)	: 1.0

skMLCK(MYLK2)

Product code 02-150

Full-length human skMLCK [1-596(end) amino acids of accession number NP_149109.1] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system. GST-skMLCK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MLCtide
ATP (μM) Km app / Bin	: 820 / 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 51
IC50 at 1 mM ATP (nM)	: 51

SLK

Product code 07-129

Full-length human SLK [1-1152(end) amino acids and S5N of accession number NP_055535.1] was expressed as N-terminal GST-fusion protein (160 kDa) using baculovirus expression system. GST-SLK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Moesin-derived peptide
ATP (μM) Km app / Bin	: 36 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.32
IC50 at 1 mM ATP (nM)	: n.a.

SPHK1

Product code 11-105

Full-length human SPHK1 [1-384(end) amino acids of accession number NP_001136074.1] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-SPHK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Sphingosine
ATP (μM) Km app / Bin	: 20 / 25
Metal	: Mg
Reference compound	: PF-543
IC50 at ATP Bin (nM)	: 3.9
IC50 at 1 mM ATP (nM)	: n.a.

SPHK2

Product code 11-106

Full-length human SPHK2 [1-618(end) amino acids of accession number NP_001191089.1] was expressed as N-terminal GST-fusion protein (92 kDa) using baculovirus expression system. GST-SPHK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Sphingosine
ATP (μM) Km app / Bin	: 620 / 600
Metal	: Mg
Reference compound	: PF-543
IC50 at ATP Bin (nM)	: 400
IC50 at 1 mM ATP (nM)	: n.a.

SRC

Product code 08-173

Full-length human SRC [1-536(end) amino acids of accession number NP_005408.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-SRC was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Srctide
ATP (μM) Km app / Bin	: 31 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 5.3
IC50 at 1 mM ATP (nM)	: 33

SRM(SRMS)

Product code 08-174

Human SRM, catalytic domain [215-488(end) amino acids of accession number NP_543013.1] was expressed as N-terminal GST-fusion protein (58kDa) using baculovirus expression system. GST-SRM was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 38 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 290
IC50 at 1 mM ATP (nM)	: 5000

SRPK1

Product code 04-160

Full-length human SRPK1 [1-655(end) amino acids and V504 deletion of accession number NP_003128.3] was expressed as N-terminal GST-fusion protein (101 kDa) using E. coli expression system. GST-SRPK1 was purified by using glutathione sepharose chromatography.

Assay platform	: IMAP
Substrate	: SRPKtide
ATP (μ M) Km app / Bin	: 200 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 85
IC50 at 1 mM ATP (nM)	: n.a.

SRPK2

Product code 04-161

Full-length human SRPK2 [1-688(end) amino acids of accession number NP_872633.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-SRPK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: DYRKtide-F
ATP (μ M) Km app / Bin	: 14 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 600
IC50 at 1 mM ATP (nM)	: n.a.

SYK

Product code 08-176

Full-length human SYK [1-635(end) amino acids of accession number NP_003168.2] was expressed as N-terminal GST-fusion protein (99 kDa) using baculovirus expression system. GST-SYK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μ M) Km app / Bin	: 59 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.29
IC50 at 1 mM ATP (nM)	: 0.63

TAK1-TAB1(MAP3K7)

Product code 09-019

Fused gene of human TAK1 [1-303 amino acids of accession number NP_663304.1] and human TAB1 [437-504 amino acids of accession number NP_006107.1] was expressed as N-terminal His-tagged protein (45kDa) using baculovirus expression system. His-tagged TAK1-TAB1 was purified by using Ni-NTA affinity chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: LRRKtide
ATP (μ M) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: 27

TAOK2

Product code 07-133

Human TAOK2, catalytic domain [1-319 amino acid of accession number NP_004774.1] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-TAOK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: TAOKtide
ATP (μM) Km app / Bin	: 39 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 24
IC50 at 1 mM ATP (nM)	: 100

TBK1

Product code 05-115

Full-length human TBK1 [1-729(end) amino acids of accession number NP_037386.1] was expressed as N-terminal GST-fusion protein (111 kDa) using baculovirus expression system. GST-TBK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CKtide
ATP (μM) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.2
IC50 at 1 mM ATP (nM)	: 6.2

TEC

Product code 08-182

Human TEC, catalytic domain [359-631(end) amino acids of accession number NP_003206.2] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system. GST-TEC was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 55 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 20
IC50 at 1 mM ATP (nM)	: 220

TIE2(TEK)

Product code 08-185

Human TIE2, cytoplasmic domain [771-1124(end) amino acids of accession number NP_000450.1] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-TIE2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Blk/Lyntide
ATP (μM) Km app / Bin	: 94 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 100
IC50 at 1 mM ATP (nM)	: 190

TNIK

Product code 07-138

Human TNIK, catalytic domain [1-314 amino acids of accession number NP_055843.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-TNIK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Moesin-derived peptide
ATP (μM) Km app / Bin	: 16 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 11

TNK1

Product code 08-104

Human TNK1, catalytic domain [106-390 amino acids of accession number Q13470-2] was expressed as N-terminal GST-fusion protein (58 kDa) using baculovirus expression system. GST-TNK1 was purified by using glutathione sepharose chromatography and gel filtration chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 71 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.55
IC50 at 1 mM ATP (nM)	: 1.7

TRKA(NTRK1)

Product code 08-186

Human TRKA, cytoplasmic domain [436-790(end) amino acids of accession number NP_001012331.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-TRKA was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSKtide
ATP (μ M) Km app / Bin	: 65 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.34
IC50 at 1 mM ATP (nM)	: 0.64

TRKB(NTRK2)

Product code 08-187

Human TRKB, cytoplasmic domain [456-822(end) amino acids of accession number NP_001018074.1] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-TRKB was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 80 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.29
IC50 at 1 mM ATP (nM)	: 0.55

TRKC(NTRK3)

Product code 08-197

Human TRKC, cytoplasmic domain [456-825(end) amino acids of accession number NP_002521.2] was expressed as N-terminal GST-fusion protein (69 kDa) using baculovirus expression system. GST-TRKC was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μ M) Km app / Bin	: 47 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.32
IC50 at 1 mM ATP (nM)	: 1.0

TSSK1

Product code 02-364

Full-length human TSSK1 [1-367(end) amino acids of accession number NP_114417.1] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-TSSK1 was purified by using glutathione sepharose chromatography. GST-TSSK1 was cleaved by PreScission protease and GST-free TSSK1 (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μ M) Km app / Bin	: 11 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 0.19
IC50 at 1 mM ATP (nM)	: 0.95

TSSK2

Product code 02-165

Full-length human TSSK2 [1-358(end) amino acids of accession number NP_443732.3] was expressed as N-terminal GST-fusion protein (68 kDa) using baculovirus expression system. GST-TSSK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 8.8 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 4.7
IC50 at 1 mM ATP (nM)	: n.a.

TSSK3

Product code 02-166

Full-length human TSSK3 [2-268(end) amino acids of accession number NP_443073.1] was expressed as N-terminal GST-fusion protein (57 kDa) using baculovirus expression system. GST-TSSK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: GS peptide
ATP (μM) Km app / Bin	: 45 / 50
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 12
IC50 at 1 mM ATP (nM)	: n.a.

TXK

Product code 08-183

Human TXK, catalytic domain [260-527(end) amino acids of accession number NP_003319.2] was expressed as N-terminal GST-fusion protein (58 kDa) using baculovirus expression system. GST-TXK was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 110 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 45
IC50 at 1 mM ATP (nM)	: 220

TYK2

Product code 08-147

Human TYK2, catalytic domain [871-1187(end) amino acids of accession number NP_003322.3] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-TYK2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 18 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.0
IC50 at 1 mM ATP (nM)	: 7.0

TYRO3

Product code 08-109

Human TYRO3, cytoplasmic domain of [453-890(end) amino acids of accession number NP_006284.2] was expressed as N-terminal GST fusion protein (76 kDa) using baculovirus expression system. GST-TYRO3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: CSK tide
ATP (μM) Km app / Bin	: 80 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.3
IC50 at 1 mM ATP (nM)	: 2.9

WNK1

Product code 05-179

Human WNK1, catalytic domain [1-491 amino acids of accession number NP_061852.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-WNK1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SPAktide
ATP (μM) Km app / Bin	: 140 / 150
Metal	: Mg+Mn
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 4500
IC50 at 1 mM ATP (nM)	: n.a.

WNK2

Product code 05-180

Human WNK2, catalytic domain [166-489 amino acids of accession number NP_006639.3] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-WNK2 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SPAktide
ATP (μM) Km app / Bin	: 48 / 50
Metal	: Mg+Mn
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 2300
IC50 at 1 mM ATP (nM)	: n.a.

WNK3

Product code 05-181

Human WNK3, catalytic domain [1-434 amino acids of accession number NP_065973.2] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-WNK3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: SPAktide
ATP (μM) Km app / Bin	: 48 / 50
Metal	: Mg+Mn
Reference compound	: K252a
IC50 at ATP Bin (nM)	: 1300
IC50 at 1 mM ATP (nM)	: n.a.

YES(YES1)

Product code 08-175

Full-length human YES [1-543(end) amino acids of accession number NP_005424.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-YES was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 13 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.4
IC50 at 1 mM ATP (nM)	: 23

YES(YES1)[T348I]

Product code 08-533

Full-length human YES [1-543(end) amino acids and T348I of accession number NP_005424.1] was expressed as N-terminal GST-fusion protein (89 kDa) using baculovirus expression system. GST-YES[T348I] was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: Src tide
ATP (μM) Km app / Bin	: 8.5 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.4
IC50 at 1 mM ATP (nM)	: 45

ZAP70

Product code 08-177

Full-length human ZAP70 [1-619(end) amino acids of accession number NP_001070] was expressed as N-terminal GST-fusion protein (97 kDa) using baculovirus expression system. GST-ZAP70 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Blk/Lyntide

ATP (μ M) Km app / Bin : 3.3 / 5
Metal : Mg+Mn
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 0.76
IC50 at 1 mM ATP (nM) : 34

BRAF

Product code 09-122

Human BRAF, catalytic domain [433-726 amino acid of accession number NP_004324.2] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-BRAF was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : MAP2K1
Cascade Assay*
Metal : Mg
Reference compound : ZM336372
IC50 at 1 mM ATP (nM) : >10000

**MAP2K1/Erk2/Modified Erktide*

BRAF[V600E]

Product code 09-144

Human BRAF, catalytic domain [433-726 amino acids and V600E of accession number NP_004324.2] was expressed as N-terminal GST-fusion protein (60 kDa) using baculovirus expression system. GST-BRAF[V600E] was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : MAP2K1
Cascade Assay*
Metal : Mg
Reference compound : ZM336372
IC50 at 1 mM ATP (nM) : 662

**MAP2K1/Erk2/Modified Erktide*

COT(MAP3K8)

Product code 07-301

Human COT, catalytic domain [30-397 amino acids of accession number NP_005195.2] was expressed as N-terminal GST-fusion protein using baculovirus expression system. GST-COT was purified by using glutathione sepharose chromatography. GST-COT was cleaved by PreScission protease and GST-free COT (42 kDa) was collected as flow-through fraction from glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : MAP2K1
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 120

**MAP2K1/Erk2/Modified Erktide*

DLK(MAP3K12)

Product code 09-111

Human DLK, catalytic domain [1-520 amino acid of accession number NP_006292.3] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system. GST-DLK was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : MAP2K4/MAP2K7
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 460

**(MAP2K4/MAP2K7)/JNK2/Modified Erktide*

MAP2K1

Product code 07-141

Full-length human MAP2K1 [1-393(end) amino acids of accession number NP_002746.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged RAF1 [306-648(end) amino acids and Y340D and Y341D of accession number NP_002871.1] using baculovirus expression system. GST-MAP2K1 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

Assay platform : Mobility Shift Assay
Substrate : Erk2
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 58

**Erk2/Modified Erktide*

MAP2K2

Product code 07-142

Full-length human MAP2K2 [1-400(end) amino acids of accession number NP_109587.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged RAF1 [306-648(end) amino acids and Y340D and Y341D of accession number NP_002871.1] using baculovirus expression system. GST-MAP2K2 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

Assay platform : Mobility Shift Assay
Substrate : Erk2
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 54

**Erk2/Modified Erktide*

MAP2K3

Product code 07-143

Full-length human MAP2K3 [1-347(end) amino acids of accession number NP_659731.1] was co-expressed as N-terminal GST-fusion protein (67 kDa) with human His-tagged MLK3 [99-398 amino acids of accession number NP_002410.1] using baculovirus expression system. GST-MAP2K3 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : p38α
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 790

**p38α/Modified Erktide*

MAP2K4

Product code 07-144

Full-length human MAP2K4 [1-399(end) amino acids of accession number NP_003001.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] using baculovirus expression system. GST-MAP2K4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : JNK2
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 4600

**JNK2/Modified Erktide*

MAP2K5

Product code 07-145

Full-length human MAP2K5 [1-448(end) amino acids of accession number NP_660143.1] was co-expressed as N-terminal GST-fusion protein (77 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP_002392.2], CDC37 and HSP90 using baculovirus expression system. GST-MAP2K5 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : Erk5
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 62

**Erk5/EGFR-derived peptide*

MAP2K6

Product code 07-146

Full-length human MAP2K6 [2-334(end) amino acids of accession number NP_002749.2] was co-expressed as N-terminal GST-fusion protein (64 kDa) with human His-tagged MLK3 [99-398 amino acids of accession number NP_002410.1] using baculovirus expression system. GST-MAP2K6 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : p38α
Cascade Assay*
Metal : Mg
Reference compound : Staurosporine
IC50 at 1 mM ATP (nM) : 140

**p38α/Modified Erktide*

MAP2K7

Product code 07-148

Full-length human MAP2K7 [1-419(end) amino acids of accession number NP_660186.1] was co-expressed as N-terminal GST-fusion protein (75 kDa) with human His-tagged MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] using baculovirus expression system. GST-MAP2K7 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: JNK2 Cascade Assay*
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 1100

**JNK2/Modified Erktide*

MAP3K1

Product code 07-103

Human MAP3K1, catalytic domain [1327-1646(end) amino acids of accession number XP_042066.8] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-MAP3K1 was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MAP2K1 Cascade Assay*
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 160

**MAP2K1/Erk2/Modified Erktide*

MAP3K2

Product code 07-104

Human MAP3K2, catalytic domain [336-619(end) amino acids of accession number NP_006600.3] was expressed as N-terminal GST-fusion protein (59 kDa) using baculovirus expression system. GST-MAP3K2 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MAP2K4/MAP2K7 Cascade Assay*
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 45

**(MAP2K4/MAP2K7)/JNK2/Modified Erktide*

MAP3K3

Product code 07-105

Full-length human MAP3K3 [1-626(end) amino acids of accession number NP_002392.2] was expressed as N-terminal GST-fusion protein (98 kDa) using baculovirus expression system. GST-fusion MAP3K3 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MAP2K6 Cascade Assay*
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 72

**MAP2K6/p38α/Modified Erktide*

MAP3K4

Product code 07-106

Human MAP3K4, catalytic domain [1312-1608(end) amino acids of accession number NP_005913.2] was expressed as N-terminal GST-fusion protein (61 kDa) using baculovirus expression system. GST-MAP3K4 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MAP2K6 Cascade Assay*
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 100

**MAP2K6/p38α/Modified Erktide*

MAP3K5

Product code 07-107

Human MAP3K5, catalytic domain [654-971 amino acids of accession number NP_005914.1] was expressed as N-terminal GST-tagged protein (62 kDa) using baculovirus expression system. GST-MAP3K5 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : MAP2K6
 Cascade Assay*
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at 1 mM ATP (nM) : 14

**MAP2K6/p38α/Modified Erktide*

MLK1(MAP3K9)

Product code 09-115

Human MLK1, catalytic domain [110-422 amino acids of accession number NP_149132.2] was expressed as N-terminal GST-fusion protein (62kDa) using baculovirus expression system. GST-MLK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
 Substrate : MAP2K1
 Cascade Assay*
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at 1 mM ATP (nM) : 11

**MAP2K1/Erk2/Modified Erktide*

MLK2(MAP3K10)

Product code 09-116

Human MLK2, catalytic domain and leucine-zipper domain [75-462 amino acids of accession number NP_002437.2] was expressed as N-terminal GST-fusion protein (71kDa) using baculovirus expression system. GST-MLK2 was purified by using glutathione sepharose chromatography and gel filtration chromatography.

Assay platform : Mobility Shift Assay
 Substrate : MAP2K1
 Cascade Assay*
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at 1 mM ATP (nM) : 45

**MAP2K1/Erk2/Modified Erktide*

MLK3(MAP3K11)

Product code 09-017

Human MLK3, catalytic domain [99-398 amino acids of accession number NP_002410.1] was expressed as N-terminal His-tagged protein (37kDa) using baculovirus expression system. His-tagged MLK3 was purified by using Ni-NTA affinity chromatography.

Assay platform : Mobility Shift Assay
 Substrate : MAP2K1
 Cascade Assay*
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at 1 mM ATP (nM) : 4.8

**MAP2K1/Erk2/Modified Erktide*

MOS

Product code 05-118

Full-length, human MOS [1-346(end) amino acids of accession number NP_005363.1] was expressed as N-terminal GST-fusion protein (65 kDa) using baculovirus expression system. GST-MOS was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform : Mobility Shift Assay
 Substrate : MAP2K1
 Cascade Assay*
 Metal : Mg
 Reference compound : Staurosporine
 IC50 at 1 mM ATP (nM) : 32

**MAP2K1/Erk2/Modified Erktide*

RAF1

Product code 09-125

Human RAF1, catalytic domain [306-648(end) amino acids and Y340D and Y341D of accession number NP_002871.1] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system. GST-RAF1 was purified by using glutathione sepharose chromatography.

Assay platform	: Mobility Shift Assay
Substrate	: MAP2K1 Cascade Assay*
Metal	: Mg
Reference compound	: ZM336372
IC50 at 1 mM ATP (nM)	: 2800

**MAP2K1/Erk2/Modified Erktide*

Assay conditions

Test compounds

The test compound is dissolved in and diluted with dimethylsulfoxide (DMSO) to achieve 100-fold higher concentration which is specified by the sponsor. Then the solution is further 25-fold diluted with assay buffer to make the final test compound solution. Reference compounds for the assay control are prepared similarly.

Assay reagents and procedures

IMAP™ assay

- 1) 4x Substrate/ATP/Metal solution and 2x kinase solution are prepared with assay buffer (20 mM HEPES, 0.01% Tween-20, 2 mM DTT, pH7.4).
- 2) 5 µL of 4x compound solution, 5 µL of 4x Substrate/ATP/Metal solution, and 10 µL of 2x kinase solution are mixed and incubated in a well of polystyrene 384 well black microplate for 1 hour at room temperature.
- 3) 60 µL of IMAP binding reagent (IMAP™ Screening Express kit; Molecular Devices) is added to the well, and incubated for over 30 minutes.
- 4) The kinase reaction is evaluated by the fluorescence polarization at 485 nm for excitation and 530 nm for emission of the well.

Off-chip Mobility Shift Assay (MSA)

- 1) 4x Substrate/ATP/Metal solution is prepared with kit buffer (20 mM HEPES, 0.01% Triton X-100, 5 mM DTT, pH7.5), and 2x kinase solution is prepared with assay buffer (20 mM HEPES, 0.01% Triton X-100, 1 mM DTT, pH7.5).
- 2) 5 µL of 4x compound solution, 5 µL of 4x Substrate/ATP/Metal solution, and 10 µL of 2x kinase solution are mixed and incubated in a well of polypropylene 384 well microplate for 1 or 5 hour(s)* at room temperature. (*; depend on kinase)
- 3) 70 µL of Termination Buffer (127 mM HEPES, 0.01% Triton X-100, 26.7 mM EDTA-2Na, 1% DMSO, pH7.5) is added to the well.
- 4) The reaction mixture is applied to LabChip™ system (PerkinElmer), and the product and substrate peptide peaks are separated and quantitated.
- 5) The kinase reaction is evaluated by the product ratio calculated from peak heights of product(P) and substrate(S) peptides ($P/(P+S)$).

Off-chip Mobility Shift Assay (MSA) with pre-incubation

- 1) 4x Substrate/ATP/Metal solution is prepared with kit buffer (20 mM HEPES, 0.01% Triton X-100, 5 mM DTT, pH7.5), and 2x kinase solution is prepared with assay buffer (20 mM HEPES, 0.01% Triton X-100, 1 mM DTT, pH7.5).
- 2) 5 µL of 4x compound solution and 10 µL of 2x kinase solution are mixed and incubated in a well of polypropylene 384 well microplate for 30 minutes at room temperature.
- 3) 5 µL of 4x Substrate/ATP/Metal solution is added to the well, and incubated for 1 hour at room temperature.
- 4) 70 µL of Termination Buffer (127 mM HEPES, 0.01% Triton X-100, 26.7 mM EDTA-2Na, 1% DMSO, pH7.5) is added to the well.
- 5) The reaction mixture is applied to LabChip™ system (PerkinElmer), and the product and substrate peptide peaks are separated and quantitated.
- 6) The kinase reaction is evaluated by the product ratio calculated from peak heights of product(P) and substrate(S) peptides ($P/(P+S)$).

ADP-Glo™ Kinase Assay

- 1) 4x compound solution and 4x ATP solution are prepared with assay buffer (50 mM MOPS, 1 mM DTT, pH7.2). 4x Substrate solution and 4x kinase/Metal solution are prepared with MOPS based buffer containing individual kinase specific additives.
- 2) 5 μ L of 4x compound solution, 5 μ L of 4x Substrate solution, 5 μ L of 4x ATP solution, and 5 μ L of 4x kinase/Metal solution are mixed and incubated in a well of polystyrene 384 well black microplate for 1 hour at room temperature.
- 3) 20 μ L of ADP-Glo™ Reagent (Promega) is added to the well, and incubated for over 40 minutes.
- 4) 40 μ L of Kinase Detection Reagent (Promega) is added to the well, and incubated for over 40 minutes.
- 5) The kinase reaction is evaluated by the endpoint luminescence of the well.

Reaction conditions

ATP Km Bin

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
ABL	MSA	ABLTide	1000	16	25	Mg	5	Staurosporine
ABL[E255K]	MSA	ABLTide	1000	17	25	Mg	5	Staurosporine
ABL[T315I]	MSA	ABLTide	1000	4.0	5	Mg	5	Staurosporine
ACK ¹⁾	MSA	WASP peptide	1000	97	100	Mg	5	Staurosporine
AKT1	MSA	Crosstide	1000	31	50	Mg	5	Staurosporine
AKT2	MSA	Crosstide	1000	110	100	Mg	5	Staurosporine
AKT3	MSA	Crosstide	1000	54	50	Mg	5	Staurosporine
ALK	MSA	Srctide	1000	57	50	Mg	5	Staurosporine
ALK[C1156Y]	MSA	Srctide	1000	64	75	Mg	5	Staurosporine
ALK[F1174L]	MSA	Srctide	1000	49	50	Mg	5	Staurosporine
ALK[G1202R]	MSA	Srctide	1000	31	50	Mg	5	Staurosporine
ALK[G1269A]	MSA	Srctide	1000	27	25	Mg	5	Staurosporine
ALK[L1196M]	MSA	Srctide	1000	57	75	Mg	5	Staurosporine
ALK[R1275Q]	MSA	Srctide	1000	84	100	Mg	5	Staurosporine
ALK[T1151_L1152insT]	MSA	Srctide	1000	110	100	Mg	5	Staurosporine
EML4-ALK ¹⁾	MSA	Srctide	1000	43	50	Mg	5	Staurosporine
NPM1-ALK	MSA	Srctide	1000	57	50	Mg	5	Staurosporine
AMPKα1/β1/γ1	MSA	SAMS peptide	1000	130	150	Mg	5	Staurosporine
AMPKα2/β1/γ1	MSA	SAMS peptide	1000	100	100	Mg	5	Staurosporine
ARG	MSA	ABLTide	1000	24	25	Mg	5	Staurosporine
AurA	MSA	Kemptide	1000	27	25	Mg	5	Staurosporine
AurA/TPX2 ⁹⁾	MSA	Kemptide	1000	1.7	2	Mg	5	Staurosporine
AurB/INCENP	MSA	Kemptide	1000	16	25	Mg	5	Staurosporine
AurC	MSA	Kemptide	1000	24	25	Mg	5	Staurosporine
AXL	MSA	CSKTide	1000	32	50	Mg	5	Staurosporine
BLK	MSA	Srctide	1000	62	75	Mg	5	Staurosporine
BMX	MSA	Srctide	1000	75	75	Mg	5	Staurosporine
BRK ¹⁾	MSA	Blk/Lyntide	1000	250	250	Mg	5	Staurosporine
BRSK1	MSA	CHKTide	1000	30	25	Mg	5	Staurosporine
BRSK2	MSA	CHKTide	1000	31	50	Mg	5	Staurosporine
BTK	MSA	Srctide	1000	22	25	Mg	5	Staurosporine
BTK[C481S]	MSA	Srctide	1000	27	25	Mg	5	Staurosporine
BUB1/BUB3	MSA	H2A peptide	1000	2.9	5	Mg	5	Staurosporine
CaMK1α ¹⁾³⁾	MSA	GS peptide	1000	750	1000	Mg	5	Staurosporine
CaMK1δ ¹⁾³⁾	MSA	Synapsin peptide	1000	11	10	Mg	5	Staurosporine
CaMK2α ³⁾	MSA	GS peptide	1000	33	50	Mg	5	Staurosporine
CaMK2β ³⁾	MSA	GS peptide	1000	19	25	Mg	5	Staurosporine
CaMK2γ ³⁾	MSA	GS peptide	1000	23	25	Mg	5	Staurosporine
CaMK2δ ³⁾	MSA	GS peptide	1000	6.3	5	Mg	5	Staurosporine
CaMK4 ³⁾	MSA	GS peptide	1000	20	25	Mg	5	Staurosporine
CDC7/ASK ¹⁾	MSA	MCM2 peptide	1000	2.8	5	Mg	10	Staurosporine
CDK1(CDC2)/CycB1	MSA	Modified Histone H1	1000	34	50	Mg	5	Staurosporine
CDK2/CycA2	MSA	Modified Histone H1	1000	27	25	Mg	5	Staurosporine
CDK2/CycE1	MSA	Modified Histone H1	1000	130	150	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
CDK3/CycE1	MSA	Modified Histone H1	1000	1000	1000	Mg	5	Staurosporine
CDK4/CycD3 ¹⁾	MSA	DYRKtide-F	1000	200	200	Mg	5	Staurosporine
CDK5/p25	MSA	Modified Histone H1	1000	10	10	Mg	5	Staurosporine
CDK6/CycD3 ¹⁾	MSA	DYRKtide-F	1000	330	300	Mg	5	Staurosporine
CDK7/CycH/MAT1 ¹⁾	MSA	CTD3 peptide	1000	32	50	Mg	5	Staurosporine
CDK9/CycT1 ¹⁾	MSA	CDK9 substrate	1000	9.4	10	Mg	5	Staurosporine
CGK2 ⁴⁾	MSA	Kemptide	1000	24	25	Mg	5	Staurosporine
CHK1	MSA	CHKtide	1000	50	50	Mg	5	Staurosporine
CHK2	MSA	CHKtide	1000	51	50	Mg	5	Staurosporine
CK1α ¹⁾	MSA	CKtide	1000	4.1	5	Mg	5	5-Iodotubercidin
CK1γ1	MSA	CKtide	1000	6.3	5	Mg	5	5-Iodotubercidin
CK1γ2	MSA	CKtide	1000	10	10	Mg	5	5-Iodotubercidin
CK1γ3	MSA	CKtide	1000	3.2	5	Mg	5	5-Iodotubercidin
CK1δ	MSA	CKtide	1000	7.7	10	Mg	5	5-Iodotubercidin
CK1ε ¹⁾	MSA	CKtide	1000	16	25	Mg	5	5-Iodotubercidin
CK2α1/β	MSA	CK2tide	1000	2.9	5	Mg	5	TBB
CK2α2/β	MSA	CK2tide	1000	2.1	5	Mg	5	TBB
CLK1	MSA	DYRKtide-F	1000	11	10	Mg	5	Staurosporine
CLK2	MSA	DYRKtide-F	1000	140	150	Mg	5	Staurosporine
CLK3	MSA	DYRKtide-F	1000	75	75	Mg	5	Staurosporine
CRIK ¹⁾	MSA	Histone H3 peptide	1000	7.8	10	Mg	5	Staurosporine
CSK ¹⁾	MSA	Srctide	1000	4.8	5	Mg, Mn	5, 1	Staurosporine
DAPK1	MSA	DAPK1tide	1000	1.1	1	Mg	5	Staurosporine
DCAMKL2 ¹⁾	MSA	GS peptide	1000	120	150	Mg	5	Staurosporine
DDR1 ¹⁾	MSA	IRS1	1000	94	100	Mg	5	Staurosporine
DDR2 ¹⁾	MSA	IRS1	1000	38	50	Mg	5	Staurosporine
DGKα	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	130	100	Mg	5	-
DGKβ	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	61	50	Mg	5	-
DGKγ	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	55	50	Mg	5	-
DGKδ	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	120	100	Mg	5	-
DGKε	ADP-Glo	Diacylglycerol, POPS	775000, 1600000	120	100	Mg	5	-
DGKζ	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	25	25	Mg	5	-
DGKη	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	24	25	Mg	5	-
DGKθ	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	37	50	Mg	5	-
DGKι	ADP-Glo	Diacylglycerol, POPS	850000, 6750000	34	50	Mg	5	-
DGKκ	ADP-Glo	Diacylglycerol, POPS	67500, 800000	17	25	Mg	1	-
DYRK1A	MSA	DYRKtide-F	1000	16	25	Mg	5	Staurosporine
DYRK1B	MSA	DYRKtide-F	1000	59	50	Mg	5	Staurosporine
DYRK2	MSA	DYRKtide-F	1000	7.7	10	Mg	5	Staurosporine
DYRK3	MSA	DYRKtide-F	1000	6.8	5	Mg	5	Staurosporine
EEF2K ¹⁾³⁾	MSA	EEF2Ktide	1000	12	10	Mg	5	A-484954
EGFR	MSA	Srctide	1000	2.7	5	Mg, Mn	5, 1	Staurosporine
EGFR[C797S/L858R]	MSA	Srctide	1000	4.1	5	Mg, Mn	5, 1	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
EGFR[d746-750]	MSA	Srctide	1000	19	25	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750/C797S]	MSA	Srctide	1000	8.2	10	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750/T790M]	MSA	Srctide	1000	5.4	5	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750/T790M/C797S]	MSA	Srctide	1000	1.8	2	Mg, Mn	5, 1	Staurosporine
EGFR[D770_N771insNPG]	MSA	Srctide	1000	2.3	5	Mg, Mn	5, 1	Staurosporine
EGFR[L858R]	MSA	Srctide	1000	9.8	10	Mg, Mn	5, 1	Staurosporine
EGFR[L861Q]	MSA	Srctide	1000	7.5	10	Mg, Mn	5, 1	Staurosporine
EGFR[T790M]	MSA	Srctide	1000	0.90	1	Mg, Mn	5, 1	Staurosporine
EGFR[T790M/L858R]	MSA	Srctide	1000	1.9	2	Mg, Mn	5, 1	Staurosporine
EPHA1	MSA	Blk/Lyntide	1000	22	25	Mg	5	Staurosporine
EPHA2	MSA	Blk/Lyntide	1000	67	75	Mg	5	Staurosporine
EPHA3	MSA	Blk/Lyntide	1000	170	150	Mg	5	Staurosporine
EPHA4	MSA	Blk/Lyntide	1000	52	50	Mg	5	Staurosporine
EPHA5	MSA	Blk/Lyntide	1000	56	50	Mg	5	Staurosporine
EPHA6	MSA	Blk/Lyntide	1000	27	25	Mg	5	Staurosporine
EPHA7	MSA	Blk/Lyntide	1000	58	50	Mg	5	Staurosporine
EPHA8	MSA	Blk/Lyntide	1000	69	75	Mg	5	Staurosporine
EPHB1	MSA	Blk/Lyntide	1000	29	25	Mg	5	Staurosporine
EPHB2	MSA	Blk/Lyntide	1000	86	100	Mg	5	Staurosporine
EPHB3	MSA	Blk/Lyntide	1000	49	50	Mg	5	Staurosporine
EPHB4	MSA	Blk/Lyntide	1000	56	50	Mg	5	Staurosporine
Erk1	MSA	Modified Erktide	1000	34	50	Mg	5	K252a
Erk2	MSA	Modified Erktide	1000	33	50	Mg	5	K252a
Erk5 ¹⁾	MSA	EGFR-derived peptide	1000	450	1000	Mg	5	Staurosporine
FAK ¹⁾	MSA	Blk/Lyntide	1000	25	25	Mg	5	Staurosporine
FER	MSA	Srctide	1000	26	25	Mg	5	Staurosporine
FES	MSA	Srctide	1000	43	50	Mg	5	Staurosporine
FGFR1	MSA	CSKtide	1000	89	100	Mg	5	Staurosporine
FGFR1[V561M]	MSA	CSKtide	1000	33	50	Mg	5	Staurosporine
FGFR2	MSA	CSKtide	1000	66	75	Mg	5	Staurosporine
FGFR2[V564I]	MSA	CSKtide	1000	21	25	Mg	5	Staurosporine
FGFR3	MSA	CSKtide	1000	43	50	Mg	5	Staurosporine
FGFR3[K650E]	MSA	CSKtide	1000	41	50	Mg	5	Staurosporine
FGFR3[K650M]	MSA	CSKtide	1000	17	25	Mg	5	Staurosporine
FGFR3[V555L]	MSA	CSKtide	1000	29	25	Mg	5	Staurosporine
FGFR3[V555M]	MSA	CSKtide	1000	37	50	Mg	5	Staurosporine
FGFR4	MSA	CSKtide	1000	230	250	Mg	5	Staurosporine
FGFR4[N535K]	MSA	CSKtide	1000	30	25	Mg	5	Staurosporine
FGFR4[V550E]	MSA	CSKtide	1000	210	200	Mg	5	Staurosporine
FGFR4[V550L]	MSA	CSKtide	1000	160	150	Mg	5	Staurosporine
FGR	MSA	Srctide	1000	34	50	Mg	5	Staurosporine
FLT1	MSA	CSKtide	1000	140	150	Mg	5	Staurosporine
FLT3	MSA	Srctide	1000	94	100	Mg	5	Staurosporine
FLT4	MSA	CSKtide	1000	72	75	Mg	5	Staurosporine
FMS	MSA	Srctide	1000	26	25	Mg	5	Staurosporine
FRK	MSA	Srctide	1000	62	75	Mg	5	Staurosporine
FYN[isoform a]	MSA	Srctide	1000	36	50	Mg	5	Staurosporine
FYN[isoform b]	MSA	Srctide	1000	20	25	Mg	5	Staurosporine
GSK3α	MSA	CREBtide-p	1000	12	10	Mg	5	Staurosporine
GSK3β	MSA	CREBtide-p	1000	9.1	10	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
Haspin	MSA	Histone H3 peptide	1000	140	150	Mg	5	Staurosporine
HCK	MSA	Srctide	1000	11	10	Mg	5	Staurosporine
HER2	MSA	Srctide	1000	3.5	5	Mn	5	Staurosporine
HER4	MSA	Srctide	1000	27	25	Mg	5	Staurosporine
HGK	MSA	Moesin-derived peptide	1000	9.4	10	Mg	5	Staurosporine
HIPK1	MSA	DYRKtide-F	1000	4.4	5	Mg	5	Staurosporine
HIPK2	MSA	DYRKtide-F	1000	5.9	5	Mg	5	Staurosporine
HIPK3	MSA	DYRKtide-F	1000	7.3	5	Mg	5	Staurosporine
HIPK4	MSA	DYRKtide-F	1000	7.0	5	Mg	5	Staurosporine
HPK1	MSA	S6K2 peptide	1000	22	25	Mg	1.25	K252a
IGF1R	MSA	IRS1	1000	63	75	Mg	5	Staurosporine
IKKα	IMAP	IκBα peptide	100	41	40	Mg	10	Staurosporine
IKKβ	MSA	Modified IκBα-derived peptide	1000	16	25	Mg	5	Staurosporine
IKKε ¹⁾	MSA	IκBα peptide	1000	9.5	10	Mg	5	Staurosporine
INSR	MSA	IRS1	1000	58	50	Mg	5	Staurosporine
IRAK1	IMAP	SRPKtide	100	27	25	Mg	2.5	Staurosporine
IRAK4 ¹⁾	MSA	IRAK1 peptide	1000	920	1000	Mg	5	Staurosporine
IRR	MSA	IRS1	1000	64	75	Mg	5	Staurosporine
ITK	MSA	Srctide	1000	6.1	10	Mg	5	Staurosporine
JAK1 ¹⁽⁶⁾	MSA	JAK1 substrate peptide	1000	68	75	Mg	5	Staurosporine
JAK2	MSA	Srctide	1000	13	10	Mg	5	Staurosporine
JAK3	MSA	Srctide	1000	3.5	5	Mg	5	Staurosporine
JNK1	MSA	Modified Erktide	1000	29	100	Mg	5	K252a
JNK2	MSA	Modified Erktide	1000	21	50	Mg	5	K252a
JNK3	MSA	Modified Erktide	1000	6.0	25	Mg	5	K252a
KDR	MSA	CSKtide	1000	74	75	Mg	5	Staurosporine
KIT ⁽⁶⁾	MSA	Srctide	1000	370	400	Mg	5	Staurosporine
KIT[D816E] ⁽⁶⁾	MSA	Srctide	1000	40	50	Mg	5	Staurosporine
KIT[D816V] ⁽⁶⁾	MSA	Srctide	1000	14	10	Mg	5	Staurosporine
KIT[D816Y] ⁽⁶⁾	MSA	Srctide	1000	22	25	Mg	5	Staurosporine
KIT[T670I] ⁽⁶⁾	MSA	Srctide	1000	100	100	Mg	5	Staurosporine
KIT[V560G] ⁽⁶⁾	MSA	Srctide	1000	110	250	Mg	5	Staurosporine
KIT[V654A] ⁽⁶⁾	MSA	Srctide	1000	220	250	Mg	5	Staurosporine
LATS1/MOBKL1A ¹⁾	MSA	SGKtide	1000	23	25	Mg	5	Staurosporine
LATS2/MOBKL1A ¹⁾	MSA	SGKtide	1000	38	50	Mg	5	Staurosporine
LCK	MSA	Srctide	1000	14	10	Mg	5	Staurosporine
LOK ¹⁾	MSA	Moesin-derived peptide	1000	100	100	Mg	5	Staurosporine
LTK	MSA	Srctide	1000	49	50	Mg	5	Staurosporine
LYNa	MSA	Srctide	1000	14	10	Mg	5	Staurosporine
LYNb	MSA	Srctide	1000	18	25	Mg	5	Staurosporine
MAP4K2	MSA	S6K2 peptide	1000	93	100	Mg	5	Staurosporine
MAPKAPK2	MSA	GS peptide	1000	3.6	5	Mg	5	Staurosporine
MAPKAPK3	MSA	GS peptide	1000	13	10	Mg	5	K252a
MAPKAPK5	MSA	GS peptide	1000	12	10	Mg	5	Staurosporine
MARK1	MSA	CHKtide	1000	8.0	10	Mg	5	Staurosporine
MARK2	MSA	CHKtide	1000	8.8	10	Mg	5	Staurosporine
MARK3	MSA	CHKtide	1000	5.0	5	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
MARK4	MSA	CHKtide	1000	12	10	Mg	5	Staurosporine
MELK ¹⁾	MSA	GS peptide	1000	38	50	Mg	5	Staurosporine
MER	MSA	CSKtide	1000	36	50	Mg	5	Staurosporine
MET	MSA	Srctide	1000	27	25	Mg	5	Staurosporine
MET[D1228H]	MSA	Srctide	1000	25	25	Mg	5	Staurosporine
MET[M1250T]	MSA	Srctide	1000	17	25	Mg	5	Staurosporine
MET[Y1235D]	MSA	Srctide	1000	71	75	Mg	5	Staurosporine
MINK ¹⁾	MSA	Modified Erktide	1000	16	50	Mg	5	K252a
MNK1	MSA	RS peptide	1000	460	450	Mg	5	Staurosporine
MNK2	MSA	RS peptide	1000	110	100	Mg	5	Staurosporine
MRCKα ¹⁾	MSA	DAPK1tide	1000	0.45	1	Mg	5	Staurosporine
MRCKβ	MSA	DAPK1tide	1000	0.67	1	Mg	5	Staurosporine
MSK1	MSA	Crosstide	1000	13	10	Mg	5	Staurosporine
MSK2 ¹⁾	MSA	Crosstide	1000	40	50	Mg	5	Staurosporine
MSSK1 ¹⁾	MSA	DYRKtide-F	1000	56	50	Mg	5	K252a
MST1 ¹⁾²⁾	MSA	IRS1	1000	50	50	Mg	5	Staurosporine
MST2 ¹⁾⁷⁾	MSA	IRS1	1000	69	75	Mg	5	Staurosporine
MST3 ¹⁾	MSA	Moesin-derived peptide	1000	66	75	Mg	5	Staurosporine
MST4 ¹⁾	MSA	Moesin-derived peptide	1000	76	75	Mg	5	Staurosporine
MUSK ¹⁾	MSA	CSKtide	1000	14	10	Mg, Mn	5, 1	Staurosporine
NDR1 ¹⁾	MSA	SGKtide	1000	12	10	Mg	5	Staurosporine
NDR2 ¹⁾	MSA	SGKtide	1000	7.6	10	Mg	5	Staurosporine
NEK1 ¹⁾	MSA	CDK7 peptide	1000	64	75	Mg	5	Staurosporine
NEK2	MSA	CDK7 peptide	1000	65	75	Mg	5	Staurosporine
NEK4	MSA	GS peptide	1000	51	50	Mg	5	Staurosporine
NEK6 ¹⁾	MSA	CDK7 peptide	1000	69	75	Mg	5	PKR Inhibitor
NEK7 ¹⁾	MSA	CDK7 peptide	1000	40	50	Mg	5	PKR Inhibitor
NEK9 ¹⁾	MSA	CDK7 peptide	1000	190	200	Mg	5	Staurosporine
NIM1K	MSA	CHKtide	1000	21	25	Mg	5	Staurosporine
NuaK1	MSA	CHKtide	1000	59	50	Mg	5	Staurosporine
NuaK2	MSA	CHKtide	1000	26	25	Mg	5	Staurosporine
p38α	MSA	Modified Erktide	1000	150	150	Mg	5	SB202190
p38β	MSA	Modified Erktide	1000	63	75	Mg	5	SB202190
p38γ	MSA	Modified Erktide	1000	13	10	Mg	5	Staurosporine
p38δ	MSA	Modified Erktide	1000	5.8	5	Mg	5	Staurosporine
p70S6K	MSA	S6K2 peptide	1000	14	10	Mg	5	Staurosporine
p70S6Kβ	MSA	S6K2 peptide	1000	3.3	5	Mg	5	Staurosporine
PAK1	MSA	LIMKtide	1000	300	300	Mg	5	Staurosporine
PAK2	MSA	DAPK1tide	1000	81	100	Mg	5	Staurosporine
PAK4 ¹⁾	MSA	SGKtide	1000	2.5	5	Mg	5	Staurosporine
PAK5	MSA	DAPK1tide	1000	1.9	1	Mg	5	Staurosporine
PAK6 ¹⁾	MSA	SGKtide	1000	3.7	5	Mg	5	Staurosporine
PASK ¹⁾	MSA	GS peptide	1000	9.7	10	Mg	5	Staurosporine
PBK ¹⁾	MSA	Histone H3 peptide	1000	33	50	Mg	5	Staurosporine
PDGFRα	MSA	CSKtide	1000	28	25	Mg	5	Staurosporine
PDGFRα[D842V]	MSA	CSKtide	1000	21	25	Mg	5	Staurosporine
PDGFRα[T674I] ¹⁾	MSA	CSKtide	1000	11	10	Mg	5	Staurosporine
PDGFRα[V561D]	MSA	CSKtide	1000	35	50	Mg	5	Staurosporine
PDGFRβ	MSA	CSKtide	1000	23	25	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
PDHK2 ¹⁾	MSA	PDHKtide	1000	28	25	Mg, K	5, 3	VER-246608
PDHK4 ¹⁾	MSA	PDHKtide	1000	19	25	Mg, K	5, 25	VER-246608
PDK1 ¹⁾⁸⁾	MSA	T308tide	1000	9.6	10	Mg	5	Staurosporine
PEK	IMAP	SRPKtide	100	13	10	Mg	5	Staurosporine
PGK ¹⁾⁴⁾	MSA	Kemptide	1000	8.2	10	Mg	5	Staurosporine
PHKG1 ¹⁾	MSA	GS peptide	1000	71	75	Mg	5	Staurosporine
PHKG2	MSA	GS peptide	1000	8.1	10	Mg	5	Staurosporine
PIK3CA/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	89	100	Mg	5	PI-103
PIK3CA[E542K]/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	42	50	Mg	5	PI-103
PIK3CA[E545K]/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	44	50	Mg	5	PI-103
PIK3CA[H1047R]/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	78	75	Mg	5	PI-103
PIK3CA[P539R]/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	37	50	Mg	5	PI-103
PIK3CA[R88Q]/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	42	50	Mg	5	PI-103
PIK3CB/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	88	100	Mg	5	PI-103
PIK3CD/PIK3R1	ADP-Glo	PI(4,5)P2, POPS	10000, 5000	37	50	Mg	5	PI-103
PIKFYVE	ADP-Glo	PI(3)P, POPS	10000, 20000	36	50	Mg	5	AG-183
PIM1	MSA	S6K2 peptide	1000	640	500	Mg	5	Staurosporine
PIM2 ¹⁾	MSA	S6K2 peptide	1000	4.0	5	Mg	5	Staurosporine
PIM3	MSA	S6K2 peptide	1000	130	150	Mg	5	Staurosporine
PIP4K2A	ADP-Glo	PI(5)P, POPS	10000, 20000	20	25	Mg	5	AG-183
PIP4K2B	ADP-Glo	PI(5)P, POPS	10000, 20000	18	25	Mn	0.25	AG-183
PIP5K1A	ADP-Glo	PI(4)P, POPS	10000, 20000	28	25	Mg	5	AG-183
PIP5K1B	ADP-Glo	PI(4)P, POPS	10000, 20000	95	100	Mg	5	AG-183
PIP5K1C	ADP-Glo	PI(4)P, POPS	10000, 20000	33	50	Mg	5	AG-183
PIP5KL1	ADP-Glo	PI(4)P, POPS	10000, 20000	1.0	1	Mg	5	AG-183
PKAα	MSA	Kemptide	1000	2.6	5	Mg	5	Staurosporine
PKAβ	MSA	Kemptide	1000	4.7	5	Mg	5	Staurosporine
PKAγ ¹⁾	MSA	Kemptide	1000	4.5	5	Mg	5	Staurosporine
PKCα ⁵⁾	MSA	PKC peptide	1000	36	50	Mg, Ca	5, 0.05	Staurosporine
PKCβ1 ⁵⁾	MSA	PKC peptide	1000	79	75	Mg, Ca	5, 0.05	Staurosporine
PKCβ2 ⁵⁾	MSA	PKC peptide	1000	41	50	Mg, Ca	5, 0.05	Staurosporine
PKCγ ⁵⁾	MSA	PKC peptide	1000	74	75	Mg, Ca	5, 0.05	Staurosporine
PKCδ ⁵⁾	MSA	PKC peptide	1000	26	25	Mg	5	Staurosporine
PKCε ⁵⁾	MSA	PKC peptide	1000	16	25	Mg	5	Staurosporine
PKCζ	MSA	PKC peptide	1000	11	10	Mg	5	Staurosporine
PKCη ⁵⁾	MSA	PKC peptide	1000	36	50	Mg	5	Staurosporine
PKCθ ⁵⁾	MSA	PKC peptide	1000	18	25	Mg	5	Staurosporine
PKCι	MSA	PKC peptide	1000	27	25	Mg	5	Staurosporine
PKD1	MSA	GS peptide	1000	25	25	Mg	5	Staurosporine
PKD2	MSA	GS peptide	1000	26	25	Mg	5	Staurosporine
PKD3	MSA	GS peptide	1000	34	50	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
PKN1	IMAP	S6K peptide	100	19	25	Mg	1	Staurosporine
PKR	IMAP	SRPKtide	100	13	10	Mg	5	Staurosporine
PLK1 ¹⁾	MSA	CDC25ctide	1000	5.6	5	Mg	5	GW843682X
PLK2	IMAP	CHK2 peptide	50	30	30	Mg	10	GW843682X
PLK3	MSA	CDC25ctide	1000	6.8	5	Mg	5	GW843682X
PRKX ¹⁾	MSA	Kemptide	1000	20	25	Mg	5	Staurosporine
PYK2	MSA	Blk/Lyntide	1000	56	50	Mg	5	Staurosporine
QIK	MSA	AMARA peptide	1000	42	50	Mg	5	Staurosporine
RET	MSA	CSKtide	1000	7.5	10	Mg	5	Staurosporine
RET[G691S]	MSA	CSKtide	1000	13	10	Mg	5	Staurosporine
RET[M918T]	MSA	CSKtide	1000	4.2	5	Mg	5	Staurosporine
RET[S891A]	MSA	CSKtide	1000	11	10	Mg	5	Staurosporine
RET[Y791F]	MSA	CSKtide	1000	29	25	Mg	5	Staurosporine
ROCK1	MSA	LIMKtide	1000	3.1	5	Mg	5	Staurosporine
ROCK2	MSA	LIMKtide	1000	7.4	5	Mg	5	Staurosporine
RON	MSA	Srctide	1000	27	25	Mg	5	Staurosporine
ROS	MSA	IRS1	1000	37	50	Mg	5	Staurosporine
RSK1	MSA	S6K peptide (N-FL)	1000	21	25	Mg	5	Staurosporine
RSK2	MSA	S6K peptide (N-FL)	1000	14	10	Mg	5	Staurosporine
RSK3	MSA	S6K peptide (N-FL)	1000	9.9	10	Mg	5	Staurosporine
RSK4	MSA	S6K peptide (N-FL)	1000	20	25	Mg	5	Staurosporine
SGK	MSA	SGKtide	1000	52	50	Mg	5	Staurosporine
SGK2	MSA	SGKtide	1000	58	50	Mg	5	Staurosporine
SGK3	MSA	SGKtide	1000	17	25	Mg	5	Staurosporine
SIK ¹⁾	MSA	AMARA peptide	1000	47	50	Mg	5	Staurosporine
skMLCK ³⁾	MSA	MLCtide	1000	820	1000	Mg	5	Staurosporine
SLK ¹⁾	MSA	Moesin-derived peptide	1000	36	50	Mg	5	Staurosporine
SPHK1	MSA	Sphingosine	1000	20	25	Mg	5	PF-543
SPHK2	MSA	Sphingosine	1000	620	600	Mg	5	PF-543
SRC	MSA	Srctide	1000	31	50	Mg	5	Staurosporine
SRM	MSA	Blk/Lyntide	1000	38	50	Mg	5	Staurosporine
SRPK1	IMAP	SRPKtide	100	200	100	Mg	10	Staurosporine
SRPK2 ¹⁾	MSA	DYRKtide-F	1000	14	10	Mg	5	Staurosporine
SYK	MSA	Blk/Lyntide	1000	59	50	Mg	5	Staurosporine
TAK1-TAB1	MSA	LRRKtide	1000	37	50	Mg	1.25	Staurosporine
TAOK2 ¹⁾⁷⁾	MSA	TAOKtide	1000	39	50	Mg	5	Staurosporine
TBK1	MSA	CKtide	1000	21	25	Mg	5	Staurosporine
TEC	MSA	Srctide	1000	55	50	Mg	5	Staurosporine
TIE2	MSA	Blk/Lyntide	1000	94	100	Mg	5	Staurosporine
TNIK	MSA	Moesin-derived peptide	1000	16	25	Mg	5	Staurosporine
TNK1 ¹⁾	MSA	CSKtide	1000	71	75	Mg	5	Staurosporine
TRKA	MSA	CSKtide	1000	65	75	Mg	5	Staurosporine
TRKB	MSA	Srctide	1000	80	75	Mg	5	Staurosporine
TRKC	MSA	Srctide	1000	47	50	Mg	5	Staurosporine
TSSK1	MSA	GS peptide	1000	11	10	Mg	5	Staurosporine
TSSK2 ¹⁾	MSA	GS peptide	1000	8.8	10	Mg	5	Staurosporine
TSSK3 ¹⁾	MSA	GS peptide	1000	45	50	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
TXK ¹⁾	MSA	Srctide	1000	110	100	Mg	5	Staurosporine
TYK2 ¹⁾	MSA	Srctide	1000	18	25	Mg	5	Staurosporine
TYRO3	MSA	CSKtide	1000	80	75	Mg	5	Staurosporine
WNK1 ¹⁾	MSA	SPAKtide	1000	140	150	Mg, Mn	5, 3	K252a
WNK2 ¹⁾	MSA	SPAKtide	1000	48	50	Mg, Mn	5, 3	K252a
WNK3 ¹⁾	MSA	SPAKtide	1000	48	50	Mg, Mn	5, 3	K252a
YES	MSA	Srctide	1000	13	10	Mg	5	Staurosporine
YES[T348I]	MSA	Srctide	1000	8.5	10	Mg	5	Staurosporine
ZAP70	MSA	Blk/Lyntide	1000	3.3	5	Mg, Mn	5, 1	Staurosporine

ATP 1mM

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
ABL	MSA	ABLTide	1000	16	1000	Mg	5	Staurosporine
ABL[E255K]	MSA	ABLTide	1000	17	1000	Mg	5	Staurosporine
ABL[T315I]	MSA	ABLTide	1000	4.0	1000	Mg	5	Staurosporine
ACK ¹⁾	MSA	WASP peptide	1000	97	1000	Mg	5	Staurosporine
AKT1	MSA	Crosstide	1000	31	1000	Mg	5	Staurosporine
ALK	MSA	Srctide	1000	57	1000	Mg	5	Staurosporine
ALK[C1156Y]	MSA	Srctide	1000	64	1000	Mg	5	Staurosporine
ALK[F1174L]	MSA	Srctide	1000	49	1000	Mg	5	Staurosporine
ALK[G1202R]	MSA	Srctide	1000	31	1000	Mg	5	Staurosporine
ALK[G1269A]	MSA	Srctide	1000	27	1000	Mg	5	Staurosporine
ALK[L1196M]	MSA	Srctide	1000	57	1000	Mg	5	Staurosporine
ALK[R1275Q]	MSA	Srctide	1000	84	1000	Mg	5	Staurosporine
ALK[T1151_L1152insT]	MSA	Srctide	1000	110	1000	Mg	5	Staurosporine
EML4-ALK ¹⁾	MSA	Srctide	1000	43	1000	Mg	5	Staurosporine
NPM1-ALK	MSA	Srctide	1000	57	1000	Mg	5	Staurosporine
AMPKα1/β1/γ1	MSA	SAMS peptide	1000	130	1000	Mg	5	Staurosporine
ARG	MSA	ABLTide	1000	24	1000	Mg	5	Staurosporine
AurA	MSA	Kemptide	1000	27	1000	Mg	5	Staurosporine
AurB/INCENP	MSA	Kemptide	1000	16	1000	Mg	5	Staurosporine
AurC	MSA	Kemptide	1000	24	1000	Mg	5	Staurosporine
AXL	MSA	CSKtide	1000	32	1000	Mg	5	Staurosporine
BLK	MSA	Srctide	1000	62	1000	Mg	5	Staurosporine
BMX	MSA	Srctide	1000	75	1000	Mg	5	Staurosporine
BRK ¹⁾	MSA	Blk/Lyntide	1000	250	1000	Mg	5	Staurosporine
BRSK1	MSA	CHKtide	1000	30	1000	Mg	5	Staurosporine
BTK	MSA	Srctide	1000	22	1000	Mg	5	Staurosporine
BTK[C481S]	MSA	Srctide	1000	27	1000	Mg	5	Staurosporine
CaMK2α ³⁾	MSA	GS peptide	1000	33	1000	Mg	5	Staurosporine
CaMK2β ³⁾	MSA	GS peptide	1000	19	1000	Mg	5	Staurosporine
CaMK2γ ³⁾	MSA	GS peptide	1000	23	1000	Mg	5	Staurosporine
CaMK2δ ³⁾	MSA	GS peptide	1000	6.3	1000	Mg	5	Staurosporine
CaMK4 ³⁾	MSA	GS peptide	1000	20	1000	Mg	5	Staurosporine
CDC7/ASK ¹⁾	MSA	MCM2 peptide	1000	2.8	1000	Mg	10	Staurosporine
CDK1(CDC2)/CycB1	MSA	Modified Histone H1	1000	34	1000	Mg	5	Staurosporine
CDK2/CycA2	MSA	Modified Histone H1	1000	27	1000	Mg	5	Staurosporine
CDK2/CycE1	MSA	Modified Histone H1	1000	130	1000	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
CDK4/CycD3 ¹⁾	MSA	DYRKtide-F	1000	200	1000	Mg	5	Staurosporine
CDK5/p25	MSA	Modified Histone H1	1000	10	1000	Mg	5	Staurosporine
CDK6/CycD3 ¹⁾	MSA	DYRKtide-F	1000	330	1000	Mg	5	Staurosporine
CDK7/CycH/MAT1 ¹⁾	MSA	CTD3 peptide	1000	32	1000	Mg	5	Staurosporine
CDK9/CycT1 ¹⁾	MSA	CDK9 substrate	1000	9.4	1000	Mg	5	Staurosporine
CHK1	MSA	CHKtide	1000	50	1000	Mg	5	Staurosporine
CHK2	MSA	CHKtide	1000	51	1000	Mg	5	Staurosporine
CK1α ¹⁾	MSA	CKtide	1000	4.1	1000	Mg	5	5-Iodotubercidin
CK1δ	MSA	CKtide	1000	7.7	1000	Mg	5	5-Iodotubercidin
CK1ε ¹⁾	MSA	CKtide	1000	16	1000	Mg	5	5-Iodotubercidin
CK2α1/β	MSA	CK2tide	1000	2.9	1000	Mg	5	TBB
CLK1	MSA	DYRKtide-F	1000	11	1000	Mg	5	Staurosporine
CLK2	MSA	DYRKtide-F	1000	140	1000	Mg	5	Staurosporine
CSK ¹⁾	MSA	Srctide	1000	4.8	1000	Mg, Mn	5, 1	Staurosporine
DAPK1	MSA	DAPK1tide	1000	1.1	1000	Mg	5	Staurosporine
DDR1 ¹⁾	MSA	IRS1	1000	94	1000	Mg	5	Staurosporine
DDR2 ¹⁾	MSA	IRS1	1000	38	1000	Mg	5	Staurosporine
DYRK1A	MSA	DYRKtide-F	1000	16	1000	Mg	5	Staurosporine
DYRK1B	MSA	DYRKtide-F	1000	59	1000	Mg	5	Staurosporine
DYRK2	MSA	DYRKtide-F	1000	7.7	1000	Mg	5	Staurosporine
DYRK3	MSA	DYRKtide-F	1000	6.8	1000	Mg	5	Staurosporine
EGFR	MSA	Srctide	1000	2.7	1000	Mg, Mn	5, 1	Staurosporine
EGFR[C797S/L858R]	MSA	Srctide	1000	4.1	1000	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750]	MSA	Srctide	1000	19	1000	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750/C797S]	MSA	Srctide	1000	8.2	1000	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750/T790M]	MSA	Srctide	1000	5.4	1000	Mg, Mn	5, 1	Staurosporine
EGFR[d746-750/T790M/C797S]	MSA	Srctide	1000	1.8	1000	Mg, Mn	5, 1	Staurosporine
EGFR[D770_N771insNPG]	MSA	Srctide	1000	2.3	1000	Mg, Mn	5, 1	Staurosporine
EGFR[L858R]	MSA	Srctide	1000	9.8	1000	Mg, Mn	5, 1	Staurosporine
EGFR[L861Q]	MSA	Srctide	1000	7.5	1000	Mg, Mn	5, 1	Staurosporine
EGFR[T790M]	MSA	Srctide	1000	0.90	1000	Mg, Mn	5, 1	Staurosporine
EGFR[T790M/C797S/L858R]	MSA	Srctide	1000	0.85	1000	Mg, Mn	5, 1	Staurosporine
EGFR[T790M/L858R]	MSA	Srctide	1000	1.9	1000	Mg, Mn	5, 1	Staurosporine
EPHA1	MSA	Blk/Lyntide	1000	22	1000	Mg	5	Staurosporine
EPHA2	MSA	Blk/Lyntide	1000	67	1000	Mg	5	Staurosporine
EPHA3	MSA	Blk/Lyntide	1000	170	1000	Mg	5	Staurosporine
EPHA4	MSA	Blk/Lyntide	1000	52	1000	Mg	5	Staurosporine
EPHA5	MSA	Blk/Lyntide	1000	56	1000	Mg	5	Staurosporine
EPHA6	MSA	Blk/Lyntide	1000	27	1000	Mg	5	Staurosporine
EPHA7	MSA	Blk/Lyntide	1000	58	1000	Mg	5	Staurosporine
EPHA8	MSA	Blk/Lyntide	1000	69	1000	Mg	5	Staurosporine
EPHB1	MSA	Blk/Lyntide	1000	29	1000	Mg	5	Staurosporine
EPHB2	MSA	Blk/Lyntide	1000	86	1000	Mg	5	Staurosporine
EPHB3	MSA	Blk/Lyntide	1000	49	1000	Mg	5	Staurosporine
EPHB4	MSA	Blk/Lyntide	1000	56	1000	Mg	5	Staurosporine
Erk1	MSA	Modified Erktide	1000	34	1000	Mg	5	K252a
Erk2	MSA	Modified Erktide	1000	33	1000	Mg	5	K252a
FAK ¹⁾	MSA	Blk/Lyntide	1000	25	1000	Mg	5	Staurosporine
FER	MSA	Srctide	1000	26	1000	Mg	5	Staurosporine
FES	MSA	Srctide	1000	43	1000	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μ M)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
FGFR1	MSA	CSKtide	1000	89	1000	Mg	5	Staurosporine
FGFR1[V561M]	MSA	CSKtide	1000	33	1000	Mg	5	Staurosporine
FGFR2	MSA	CSKtide	1000	66	1000	Mg	5	Staurosporine
FGFR2[V564I]	MSA	CSKtide	1000	21	1000	Mg	5	Staurosporine
FGFR3	MSA	CSKtide	1000	43	1000	Mg	5	Staurosporine
FGFR3[K650E]	MSA	CSKtide	1000	41	1000	Mg	5	Staurosporine
FGFR3[K650M]	MSA	CSKtide	1000	17	1000	Mg	5	Staurosporine
FGFR3[V555L]	MSA	CSKtide	1000	29	1000	Mg	5	Staurosporine
FGFR3[V555M]	MSA	CSKtide	1000	37	1000	Mg	5	Staurosporine
FGFR4	MSA	CSKtide	1000	230	1000	Mg	5	Staurosporine
FGFR4[N535K]	MSA	CSKtide	1000	30	1000	Mg	5	Staurosporine
FGFR4[V550E]	MSA	CSKtide	1000	210	1000	Mg	5	Staurosporine
FGFR4[V550L]	MSA	CSKtide	1000	160	1000	Mg	5	Staurosporine
FGR	MSA	Srctide	1000	34	1000	Mg	5	Staurosporine
FLT1	MSA	CSKtide	1000	140	1000	Mg	5	Staurosporine
FLT3	MSA	Srctide	1000	94	1000	Mg	5	Staurosporine
FLT4	MSA	CSKtide	1000	72	1000	Mg	5	Staurosporine
FMS	MSA	Srctide	1000	26	1000	Mg	5	Staurosporine
FRK	MSA	Srctide	1000	62	1000	Mg	5	Staurosporine
FYN[isoform a]	MSA	Srctide	1000	36	1000	Mg	5	Staurosporine
FYN[isoform b]	MSA	Srctide	1000	20	1000	Mg	5	Staurosporine
GSK3 α	MSA	CREBtide-p	1000	12	1000	Mg	5	Staurosporine
GSK3 β	MSA	CREBtide-p	1000	9.1	1000	Mg	5	Staurosporine
HCK	MSA	Srctide	1000	11	1000	Mg	5	Staurosporine
HER2	MSA	Srctide	1000	3.5	1000	Mn	5	Staurosporine
HER4	MSA	Srctide	1000	27	1000	Mg	5	Staurosporine
HGK	MSA	Moesin-derived peptide	1000	9.4	1000	Mg	5	Staurosporine
HIPK3	MSA	DYRKtide-F	1000	7.3	1000	Mg	5	Staurosporine
HIPK4	MSA	DYRKtide-F	1000	7.0	1000	Mg	5	Staurosporine
IGF1R	MSA	IRS1	1000	63	1000	Mg	5	Staurosporine
IKK β	MSA	Modified I κ B α -derived peptide	1000	16	1000	Mg	5	Staurosporine
INSR	MSA	IRS1	1000	58	1000	Mg	5	Staurosporine
IRR	MSA	IRS1	1000	64	1000	Mg	5	Staurosporine
ITK	MSA	Srctide	1000	6.1	1000	Mg	5	Staurosporine
JAK1 ^{1/6)}	MSA	JAK1 substrate peptide	1000	68	1000	Mg	5	Staurosporine
JAK2	MSA	Srctide	1000	13	1000	Mg	5	Staurosporine
JAK3	MSA	Srctide	1000	3.5	1000	Mg	5	Staurosporine
JNK1	MSA	Modified Erktide	1000	29	1000	Mg	5	K252a
JNK2	MSA	Modified Erktide	1000	21	1000	Mg	5	K252a
JNK3	MSA	Modified Erktide	1000	6.0	1000	Mg	5	K252a
KDR	MSA	CSKtide	1000	74	1000	Mg	5	Staurosporine
KIT ⁶⁾	MSA	Srctide	1000	370	1000	Mg	5	Staurosporine
KIT[D816E] ⁶⁾	MSA	Srctide	1000	40	1000	Mg	5	Staurosporine
KIT[D816V] ⁶⁾	MSA	Srctide	1000	14	1000	Mg	5	Staurosporine
KIT[D816Y] ⁶⁾	MSA	Srctide	1000	22	1000	Mg	5	Staurosporine
KIT[T670I] ⁶⁾	MSA	Srctide	1000	100	1000	Mg	5	Staurosporine
KIT[V560G] ⁶⁾	MSA	Srctide	1000	110	1000	Mg	5	Staurosporine
KIT[V654A] ⁶⁾	MSA	Srctide	1000	220	1000	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
LATS1/MOBKL1A ¹⁾	MSA	SGKtide	1000	23	1000	Mg	5	Staurosporine
LATS2/MOBKL1A ¹⁾	MSA	SGKtide	1000	38	1000	Mg	5	Staurosporine
LCK	MSA	Srctide	1000	14	1000	Mg	5	Staurosporine
LTK	MSA	Srctide	1000	49	1000	Mg	5	Staurosporine
LYNa	MSA	Srctide	1000	14	1000	Mg	5	Staurosporine
LYNb	MSA	Srctide	1000	18	1000	Mg	5	Staurosporine
MAPKAPK2	MSA	GS peptide	1000	3.6	1000	Mg	5	Staurosporine
MARK4	MSA	CHKtide	1000	12	1000	Mg	5	Staurosporine
MER	MSA	CSKtide	1000	36	1000	Mg	5	Staurosporine
MET	MSA	Srctide	1000	27	1000	Mg	5	Staurosporine
MET[D1228H]	MSA	Srctide	1000	25	1000	Mg	5	Staurosporine
MET[M1250T]	MSA	Srctide	1000	17	1000	Mg	5	Staurosporine
MET[Y1235D]	MSA	Srctide	1000	71	1000	Mg	5	Staurosporine
MINK ¹⁾	MSA	Modified Erktide	1000	16	1000	Mg	5	K252a
MNK2	MSA	RS peptide	1000	110	1000	Mg	5	Staurosporine
MSK1	MSA	Crosstide	1000	13	1000	Mg	5	Staurosporine
MSK2 ¹⁾	MSA	Crosstide	1000	40	1000	Mg	5	Staurosporine
MST1 ¹⁾²⁾	MSA	IRS1	1000	50	1000	Mg	5	Staurosporine
MST2 ¹⁾⁷⁾	MSA	IRS1	1000	69	1000	Mg	5	Staurosporine
MUSK ¹⁾	MSA	CSKtide	1000	14	1000	Mg, Mn	5, 1	Staurosporine
NEK1 ¹⁾	MSA	CDK7 peptide	1000	64	1000	Mg	5	Staurosporine
NEK2	MSA	CDK7 peptide	1000	65	1000	Mg	5	Staurosporine
NEK6 ¹⁾	MSA	CDK7 peptide	1000	69	1000	Mg	5	PKR Inhibitor
NEK7 ¹⁾	MSA	CDK7 peptide	1000	40	1000	Mg	5	PKR Inhibitor
NEK9 ¹⁾	MSA	CDK7 peptide	1000	190	1000	Mg	5	Staurosporine
NuaK1	MSA	CHKtide	1000	59	1000	Mg	5	Staurosporine
NuaK2	MSA	CHKtide	1000	26	1000	Mg	5	Staurosporine
p38α	MSA	Modified Erktide	1000	150	1000	Mg	5	SB202190
p38β	MSA	Modified Erktide	1000	63	1000	Mg	5	SB202190
p38γ	MSA	Modified Erktide	1000	13	1000	Mg	5	Staurosporine
p38δ	MSA	Modified Erktide	1000	5.8	1000	Mg	5	Staurosporine
p70S6K	MSA	S6K2 peptide	1000	14	1000	Mg	5	Staurosporine
PAK1	MSA	LIMKtide	1000	300	1000	Mg	5	Staurosporine
PAK2	MSA	DAPK1tide	1000	81	1000	Mg	5	Staurosporine
PAK5	MSA	DAPK1tide	1000	1.9	1000	Mg	5	Staurosporine
PASK ¹⁾	MSA	GS peptide	1000	9.7	1000	Mg	5	Staurosporine
PBK ¹⁾	MSA	Histone H3 peptide	1000	33	1000	Mg	5	Staurosporine
PDGFRα	MSA	CSKtide	1000	28	1000	Mg	5	Staurosporine
PDGFRα[D842V]	MSA	CSKtide	1000	21	1000	Mg	5	Staurosporine
PDGFRα[T674I] ¹⁾	MSA	CSKtide	1000	11	1000	Mg	5	Staurosporine
PDGFRα[V561D]	MSA	CSKtide	1000	35	1000	Mg	5	Staurosporine
PDGFRβ	MSA	CSKtide	1000	23	1000	Mg	5	Staurosporine
PDK1 ¹⁾⁸⁾	MSA	T308tide	1000	9.6	1000	Mg	5	Staurosporine
PIM1	MSA	S6K2 peptide	1000	640	1000	Mg	5	Staurosporine
PIM2 ¹⁾	MSA	S6K2 peptide	1000	4.0	1000	Mg	5	Staurosporine
PIM3	MSA	S6K2 peptide	1000	130	1000	Mg	5	Staurosporine
PKACα	MSA	Kemptide	1000	2.6	1000	Mg	5	Staurosporine
PKACβ	MSA	Kemptide	1000	4.7	1000	Mg	5	Staurosporine
PKCα ⁵⁾	MSA	PKC peptide	1000	36	1000	Mg, Ca	5, 0.05	Staurosporine
PKCγ ⁵⁾	MSA	PKC peptide	1000	74	1000	Mg, Ca	5, 0.05	Staurosporine
PKCε ⁵⁾	MSA	PKC peptide	1000	16	1000	Mg	5	Staurosporine

Kinase	Platform	Substrate		ATP (μ M)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
PKD2	MSA	GS peptide	1000	26	1000	Mg	5	Staurosporine
PLK1 ¹⁾	MSA	CDC25ctide	1000	5.6	1000	Mg	5	GW843682X
PLK3	MSA	CDC25ctide	1000	6.8	1000	Mg	5	GW843682X
PYK2	MSA	Blk/Lyntide	1000	56	1000	Mg	5	Staurosporine
QIK	MSA	AMARA peptide	1000	42	1000	Mg	5	Staurosporine
RET	MSA	CSKtide	1000	7.5	1000	Mg	5	Staurosporine
RET[G691S]	MSA	CSKtide	1000	13	1000	Mg	5	Staurosporine
RET[M918T]	MSA	CSKtide	1000	4.2	1000	Mg	5	Staurosporine
RET[S891A]	MSA	CSKtide	1000	11	1000	Mg	5	Staurosporine
RET[Y791F]	MSA	CSKtide	1000	29	1000	Mg	5	Staurosporine
ROCK1	MSA	LIMKtide	1000	3.1	1000	Mg	5	Staurosporine
ROCK2	MSA	LIMKtide	1000	7.4	1000	Mg	5	Staurosporine
RON	MSA	Srctide	1000	27	1000	Mg	5	Staurosporine
ROS	MSA	IRS1	1000	37	1000	Mg	5	Staurosporine
RSK1	MSA	S6K peptide (N-FL)	1000	21	1000	Mg	5	Staurosporine
RSK2	MSA	S6K peptide (N-FL)	1000	14	1000	Mg	5	Staurosporine
RSK3	MSA	S6K peptide (N-FL)	1000	9.9	1000	Mg	5	Staurosporine
RSK4	MSA	S6K peptide (N-FL)	1000	20	1000	Mg	5	Staurosporine
SGK	MSA	SGKtide	1000	52	1000	Mg	5	Staurosporine
SIK ¹⁾	MSA	AMARA peptide	1000	47	1000	Mg	5	Staurosporine
SRC	MSA	Srctide	1000	31	1000	Mg	5	Staurosporine
SRM	MSA	Blk/Lyntide	1000	38	1000	Mg	5	Staurosporine
SYK	MSA	Blk/Lyntide	1000	59	1000	Mg	5	Staurosporine
TAK1-TAB1	MSA	LRRKtide	1000	37	1000	Mg	1.25	Staurosporine
TAOK2 ¹⁷⁾	MSA	TAOKtide	1000	39	1000	Mg	5	Staurosporine
TBK1	MSA	CKtide	1000	21	1000	Mg	5	Staurosporine
TEC	MSA	Srctide	1000	55	1000	Mg	5	Staurosporine
TIE2	MSA	Blk/Lyntide	1000	94	1000	Mg	5	Staurosporine
TNIK	MSA	Moesin-derived peptide	1000	16	1000	Mg	5	Staurosporine
TNK1 ¹⁾	MSA	CSKtide	1000	71	1000	Mg	5	Staurosporine
TRKA	MSA	CSKtide	1000	65	1000	Mg	5	Staurosporine
TRKB	MSA	Srctide	1000	80	1000	Mg	5	Staurosporine
TRKC	MSA	Srctide	1000	47	1000	Mg	5	Staurosporine
TSSK1	MSA	GS peptide	1000	11	1000	Mg	5	Staurosporine
TXK ¹⁾	MSA	Srctide	1000	110	1000	Mg	5	Staurosporine
TYK2 ¹⁾	MSA	Srctide	1000	18	1000	Mg	5	Staurosporine
TYRO3	MSA	CSKtide	1000	80	1000	Mg	5	Staurosporine
YES	MSA	Srctide	1000	13	1000	Mg	5	Staurosporine
YES[T348I]	MSA	Srctide	1000	8.5	1000	Mg	5	Staurosporine
ZAP70	MSA	Blk/Lyntide	1000	3.3	1000	Mg, Mn	5, 1	Staurosporine

- 1) Reaction time is 5 hours.
- 2) Cantharidin mixed in Substrate/ATP/Metal solution is added at the final concentration of 20 μ M.
- 3) CaCl_2 and Calmodulin mixed in Substrate/ATP/Metal solution are added at the final concentration of 1 mM and 10 μ g/ml, respectively.
- 4) cGMP mixed in Substrate/ATP/Metal solution is added at the final concentration of 5 μ M.
- 5) Phosphatidylserine and Diacyl Glycerol mixed in Substrate/ATP/Metal solution are added at the final concentration of 50 μ g/mL and 5 μ g/mL, respectively.
- 6) Sodium orthovanadate mixed in Substrate/ATP/Metal solution is added at the final concentration of 25 μ M.
- 7) Cantharidin mixed in Substrate/ATP/Metal solution is added at the final concentration of 10 μ M.
- 8) PIFtide and Cantharidin mixed in Substrate/ATP/Metal solution are added at the final concentration of 2 μ M and 20 μ M, respectively.
- 9) TPX2 peptide mixed in Substrate/ATP/Metal solution is added at the final concentration of 200 nM.

Cascade assay

Kinase	Platform	Substrate		ATP (μM)		Metal		Positive control
		Name	(nM)	Km	Assay	Name	(mM)	
BRAF	MSA	MAP2K1	1	-	1000	Mg	5	ZM336372
BRAF[V600E]	MSA	MAP2K1	1	-	1000	Mg	5	ZM336372
COT	MSA	MAP2K1	1	-	1000	Mg	5	Staurosporine
DLK ¹⁾	MSA	MAP2K4, MAP2K7	0.5, 0.5	-	1000	Mg	5	Staurosporine
MAP2K1	MSA	Erk2	2.5	-	1000	Mg	5	Staurosporine
MAP2K2	MSA	Erk2	2.5	-	1000	Mg	5	Staurosporine
MAP2K3	MSA	p38a	10	-	1000	Mg	5	Staurosporine
MAP2K4 ¹⁾	MSA	JNK2	50	-	1000	Mg	5	Staurosporine
MAP2K5 ¹⁾	MSA	Erk5	50	-	1000	Mg	5	Staurosporine
MAP2K6	MSA	p38a	10	-	1000	Mg	5	Staurosporine
MAP2K7 ¹⁾	MSA	JNK2	50	-	1000	Mg	5	Staurosporine
MAP3K1	MSA	MAP2K1	1	-	1000	Mg	5	Staurosporine
MAP3K2 ¹⁾	MSA	MAP2K4, MAP2K7	0.5, 0.5	-	1000	Mg	5	Staurosporine
MAP3K3	MSA	MAP2K6	1	-	1000	Mg	5	Staurosporine
MAP3K4	MSA	MAP2K6	1	-	1000	Mg	5	Staurosporine
MAP3K5	MSA	MAP2K6	1	-	1000	Mg	5	Staurosporine
MLK1	MSA	MAP2K1	1	-	1000	Mg	5	Staurosporine
MLK2	MSA	MAP2K1	1	-	1000	Mg	5	Staurosporine
MLK3	MSA	MAP2K1	1	-	1000	Mg	5	Staurosporine
MOS	MSA	MAP2K1	1	-	1000	Mg	5	Staurosporine
RAF1	MSA	MAP2K1	1	-	1000	Mg	5	ZM336372

1) Reaction time is 5 hours.

Substrate information of cascade assay

Kinase	Substrate					
	MAP2K	(nM)	MAPK	(nM)	peptide	(nM)
BRAF	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
BRAF[V600E]	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
COT	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
DLK	MAP2K4, MAP2K7	0.5, 0.5	JNK2	50	Modified Erktide	1000
MAP2K1	-	-	Erk2	2.5	Modified Erktide	1000
MAP2K2	-	-	Erk2	2.5	Modified Erktide	1000
MAP2K3	-	-	p38 α	10	Modified Erktide	1000
MAP2K4	-	-	JNK2	50	Modified Erktide	1000
MAP2K5	-	-	Erk5	50	EGFR-derived peptide	1000
MAP2K6	-	-	p38 α	10	Modified Erktide	1000
MAP2K7	-	-	JNK2	50	Modified Erktide	1000
MAP3K1	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
MAP3K2	MAP2K4, MAP2K7	0.5, 0.5	JNK2	50	Modified Erktide	1000
MAP3K3	MAP2K6	1	p38 α	10	Modified Erktide	1000
MAP3K4	MAP2K6	1	p38 α	10	Modified Erktide	1000
MAP3K5	MAP2K6	1	p38 α	10	Modified Erktide	1000
MLK1	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
MLK2	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
MLK3	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
MOS	MAP2K1	1	Erk2	2.5	Modified Erktide	1000
RAF1	MAP2K1	1	Erk2	2.5	Modified Erktide	1000

Data analysis

The readout value of reaction control (complete reaction mixture) is set as a 0% inhibition, and the readout value of background (Enzyme(-)) is set as a 100% inhibition, then the percent inhibition of each test solution is calculated.

IC₅₀ value is calculated from concentration vs. %Inhibition curves by fitting to a four parameter logistic curve.

<http://www.carnabio.com>

Carna Biosciences, Inc.

BMA 3F, 1-5-5 Minatojima-Minamimachi, Chuo-ku, Kobe 650-0047 JAPAN

TEL: 078-302-7091 / FAX: 078-302-7086

E-mail: info@carnabio.com

Carna's wholly-owned subsidiary

CarnaBio USA, Inc.

209 West Central Street, Suite 307, Natick, MA 01760 USA

TEL: +1-508-650-1244 / Toll-Free: +1-888-645-1233

E-mail: orders@carnabio.com / FAX: +1-508-650-1722