

Cytoplasmic Tyrosine Kinases

	Kinase	Profiling
ABL	●	●
ABL[E255K]	●	●
ABL[T315I]	●	●
ACK	●	●
ARG	●	●
BLK	●	●
BMX	●	●
BRK	●	●
BTK	●	●
BTK [C481S]	●	●
CSK	●	●
FAK	●	●
FER	●	●
FES	●	●
FGR	●	●
FRK	●	●
FYN [isoform a]	●	●
FYN [isoform b]	●	●
HCK	●	●
ITK	●	●
JAK1	●	●
JAK2	●	●
JAK2 (JH1 JH2)	●	▲
JAK2 (JH1 JH2)[V617F]	●	▲
JAK3	●	●
LCK	●	●
LYNa	●	●
LYNb	●	●
PYK2	●	●
SRC	●	●
SRM	●	●
SYK	●	●
TEC	●	●
TNK1	●	●
TXK	●	●
TYK2	●	●
YES	●	●
YES (YES1)[T348I]	●	●
ZAP70	●	●

Pack sizes are 5, 100, 200, 500 and 1000ug.

Bulk production is available upon receipt of purchase order.

Please see Serine/Threonine Kinases on reverse side.

Lipid Kinases

	Kinase	Profiling	
DGKa (DGKA)	●	●*	5ug/vial
DGKβ (DGKB)	●	●*	5ug/vial
DGKy (DGKG)	●	●*	5ug/vial
DGKδ (DGKD)	●	●*	5ug/vial
DGKe (DGKE)	●	●*	5ug/vial
DGKη (DGKH)	●	●*	5ug/vial
DGKi (DGKI)	●	●*	5ug/vial
DGKk (DGKK)	●	●*	5ug/vial
DGKθ (DGKQ)	●	●*	5ug/vial
DGKζ (DGKZ)	●	●*	5ug/vial
PIK3CA/PIK3R1	●	▲	20ug/vial
PIK3CB/PIK3R1	●	▲	20ug/vial
PIK3CD/PIK3R1	●	▲	20ug/vial
PIK3C3	●	▲	20ug/vial
PIKFYVE(PIP5K3)	●	▲	5ug/vial
PI4KB	●	▲	20ug/vial
PIP4K2A	●	▲	5ug/vial
PIP4K2B	●	▲	5ug/vial
PIP4K2C	●	▲	5ug/vial
PIP5K1A	●	▲	5ug/vial
PIP5K1B	●	▲	5ug/vial
PIP5K1C	●	▲	5ug/vial
SPHK1	●	●	5ug/vial
SPHK2	●	●	5ug/vial

\* Provided by biotinylated protein

	Protein	Profiling (Km app.)	Profiling (1mM)
Cytoplasmic Tyrosine	32	37	37
Receptor Tyrosine	46	86	86
Serine/Threonine	266	200	69
Lipid	24	13	0
<b>Total</b>	<b>368</b>	<b>336</b>	<b>192</b>
Biotinylated	<b>84</b>		

Updated: 2018/07/10

Receptor Tyrosine Kinases

	Kinase	Profiling
ALK	●	●
ALK[C1156Y]	●	●
ALK[F1174L]	●	●
ALK[G1202R]	●	●
ALK[G1269A]	●	●
ALK[L1196M]	●	●
ALK[R1275Q]	●	●
ALK[T1151_L1152insT]	●	●
EML4-ALK	●	●
NPM1-ALK	●	●
AXL	●	●
DDR1	●	●
DDR2	●	●
EGFR	●	●
EGFR[d746-750]	●	●
EGFR[d746-750/T790M]	●	●
EGFR[D770_N771insNPG]	●	▲
EGFR[L858R]	●	●
EGFR[L861Q]	●	●
EGFR[T790M]	●	●
EGFR[T790M/L858R]	●	●
EphA1	●	●
EphA2	●	●
EphA3	●	●
EphA4	●	●
EphA5	●	●
EphA6	●	●
EphA7	●	●
EphA8	●	●
EphB1	●	●
EphB2	●	●
EphB3	●	●
EphB4	●	●
FGFR1	●	●
FGFR1[V561M]	●	●
FGFR2	●	●
FGFR2[N549H]	●	▲
FGFR2[V564I]	●	●
FGFR3	●	●
FGFR3[G697C]	●	▲
FGFR3[K650E]	●	●
FGFR3[K650M]	●	●
FGFR3[V555L]	●	●
FGFR3[V555M]	●	●
FGFR4	●	●
FGFR4[N535K]	●	●
FGFR4[V550E]	●	●

	Kinase	Profiling
FGFR4[V550L]	●	●
FLT1	●	●
FLT3	●	●
FLT4	●	●
FMS	●	●
HER2	●	●
HER4	●	●
IGF1R	●	●
INSR	●	●
IRR	●	●
KDR	●	●
KIT	●	●
KIT[D816E]	●	●
KIT[D816V]	●	●
KIT[D816V]	●	●
KIT[T670I]	●	●
KIT[T670I/ΔD816V]	●	▲
KIT[V560G]	●	●
KIT[V560G/D816V]	●	▲
KIT[V654A]	●	▲
KIT[V654A/ΔD816V]	●	▲
LTK	●	●
MER	●	●
MET	●	●
MET[D1228H]	●	●
MET[M1250T]	●	●
MET[Y1235D]	●	●
MUSK	●	●
PDGFRa	●	●
PDGFRa[D842V]	●	●
PDGFRa[T674I]	●	●
PDGFRa[V561D]	●	●
PDGFRb	●	●
RET	●	●
RET[G691S]	●	●
RET[M918T]	●	●
RET[S891A]	●	●
RET[Y791F]	●	●
RON	●	●
ROS	●	●
TIE2	●	●
TRKA	●	●
TRKB	●	●
TRKC	●	●
TYRO3	●	●

Biotinylated Kinases

	Kinase	
BTN - ABL (ABL1)	●	10, 100ug/vial
BTN - ACTR2B (ACVR2B)	●	10, 100ug/vial
BTN - AKT1	●	10, 100ug/vial
BTN - AKT2	●	10, 100ug/vial
BTN - ALK	●	10, 100ug/vial
BTN - ALK[L1196M]	●	10, 100ug/vial
BTN - AurA (AURKA)	●	10, 100ug/vial
BTN - AurB (AURKB)/INCENP	●	10, 100ug/vial
BTN - AXL	●	10, 100ug/vial
BTN - BMX	●	10, 100ug/vial
BTN - BRAF	●	10, 100ug/vial
BTN - BTK	●	10, 100ug/vial
BTN - BTK[C481S]	●	10, 100ug/vial
BTN - BTK[T316A]	●	10, 100ug/vial
BTN - BTK[T474I]	●	10, 100ug/vial
BTN - BTK[T474S]	●	10, 100ug/vial
BTN - CHK1 (CHEK1)	●	10, 100ug/vial
BTN - DDR2	●	10, 100ug/vial
BTN - DGKa (DGKA)	●	10, 100ug/vial
BTN - DGKβ (DGKB)	●	10, 100ug/vial
BTN - DGKy (DGKG)	●	10, 100ug/vial
BTN - DGKδ (DGKD)	●	10, 100ug/vial
BTN - DGKe (DGKE)	●	10, 100ug/vial
BTN - DGKη (DGKH)	●	10, 100ug/vial
BTN - DGKi (DGKI)	●	10, 100ug/vial
BTN - DGKk (DGKK)	●	10, 100ug/vial
BTN - DGKθ (DGKQ)	●	10, 100ug/vial
BTN - DGKζ (DGKZ)	●	10, 100ug/vial
BTN - DLK (MAP3K12)	●	10, 100ug/vial
BTN - EGFR	●	10, 100ug/vial
BTN - EGFR[T790M/L858R]	●	10, 100ug/vial
BTN - FGFR1	●	10, 100ug/vial
BTN - FGFR3	●	10, 100ug/vial
BTN - FGFR4	●	10, 100ug/vial
BTN - FAK (PTK2)	●	10, 100ug/vial
BTN - FLT3	●	10, 100ug/vial
BTN - FMS (CSF1R)	●	10, 100ug/vial
BTN - FYN[isoform a]	●	10, 100ug/vial
BTN - IGF1R	●	10, 100ug/vial
BTN - IKKa (CHUK)	●	10, 100ug/vial
BTN - IKKa (CHUK) [inactive mutant]	●	10, 100ug/vial
BTN - IKKb	●	10, 100ug/vial

	Kinase	
BTN - IRAK4	●	10, 100ug/vial
BTN - ITK	●	10, 100ug/vial
BTN - JAK1	●	10, 100ug/vial
BTN - JAK2	●	10, 100ug/vial
BTN - KDR (VEGFR2)	●	10, 100ug/vial
BTN - KIT	●	10, 100ug/vial
BTN - LCK	●	10, 100ug/vial
BTN - LYNa	●	10, 100ug/vial
BTN - MAP2K1 [inactive]	●	10, 100ug/vial
BTN - MAP3K5	●	10, 100ug/vial
BTN - MET	●	10, 100ug/vial
BTN - MNK1 (MKNK1)	●	10, 100ug/vial
BTN - NEK9	●	10, 100ug/vial
BTN - p38α (MAPK14)	●	10, 100ug/vial
BTN - p38α (MAPK14)[inactive]	●	10, 100ug/vial
BTN - PAK4	●	10, 100ug/vial
BTN - PIK3CA/PIK3R1	●	10, 100ug/vial
BTN - PIK3CB/PIK3R1	●	10, 100ug/vial
BTN - PIK3CD/PIK3R1	●	10, 100ug/vial
BTN - PIK3CG/PIK3R1	●	10, 100ug/vial
BTN - PKCη (PRKCH)	●	10, 100ug/vial
BTN - PKN1	●	10, 100ug/vial
BTN - PLK1	●	10, 100ug/vial
BTN - PYK2 (PTK2B)	●	10, 100ug/vial
BTN - RAF1	●	10, 100ug/vial
BTN - RET	●	10, 100ug/vial
BTN - ROCK1	●	10, 100ug/vial
BTN - SPHK1	●	10, 100ug/vial
BTN - SPHK2	●	10, 100ug/vial
BTN - SRC	●	10, 100ug/vial
BTN - SRPK1	●	10, 100ug/vial
BTN - SYK	●	10, 100ug/vial
BTN - TGFbR1 (TGFBFR1)	●	10, 100ug/vial
BTN - TIE2	●	10, 100ug/vial
BTN - TNK1	●	10, 100ug/vial
BTN - WNK1	●	10, 100ug/vial
BTN - WNK4	●	10, 100ug/vial
BTN - TAK1-TAB1(MAP3K7)	●	10, 100ug/vial
BTN - TRKA (NTRK1)	●	10, 100ug/vial
BTN - TRKB (NTRK2)	●	10, 100ug/vial
BTN - TRKC (NTRK3)	●	10, 100ug/vial
BTN - TYK2	●	10, 100ug/vial

Note: Please contact us for availability or further information. Information may be changed without notice.

Serine/Threonine Kinases

Kinase		Profiling	Kinase		Profiling	Kinase		Profiling	Kinase		Profiling
ACTR2B	●	▲	DAPK1	●	●	MAPKAPK5	●	●	PKCb1 (PRKCB1)	●	●
ACVR2A	●	▲	DAPK3	●	▲	MARK1	●	●	PKCb2 (PRKCB2)	●	●
ACVRL1(ALK1)	●	▲	DCAMKL1	●	▲	MARK2	●	●	PKCg (PRKCG)	●	●
AKT1	●	●	DCAMKL2	●	●	MARK3	●	●	PKCd (PRKCD)	●	●
AKT2	●	●	DLK (MAP3K12)	●	●	MARK4	●	●	PKCe (PRKCE)	●	●
AKT3	●	●	DMPK1 (DMPK)	●	▲	MELK	●	●	PKCz (PRKCZ)	●	●
ALK2 (ACVR1)	●	▲	DMPK2 (CDC42BPG)	●	▲	MINK (MINK1)	●	●	PKCh (PRKCH)	●	●
ALK2 (ACVR1)[Q207D]	●	▲	DRAK1 (STK17A)	●	▲	MLK1 (MAP3K9)	●	●	PKCq (PRKCO)	●	●
ALK2 (ACVR1)[R206H]	●	▲	DYRK1A	●	●	MLK2 (MAP3K10)	●	●	PKCi (PRKCI)	●	●
ALK4 (ACVR1B)	●	▲	DYRK1B	●	●	MLK3 (MAP3K11)	●	●	PKD1 (PRKD1)	●	●
AMPKa1/b1/g1(PRKAA1/B1/G1)	●	●	DYRK2	●	●	MLK4	●	▲	PKD2 (PRKD2)	●	●
AMPKa2/b1/g1(PRKAA2/B1/G1)	●	●	DYRK3	●	●	MNK1 (MKNK1)	●	●	PKD3 (PRKD3)	●	●
AMPKa1/b2/g1(PRKAA1/B2/G1)	●	▲	DYRK4	●	▲	MNK2 (MKNK2)	●	●	PKN1	●	●
AMPKa2/b2/g1(PRKAA2/B2/G1)	●	▲	EEF2K	●	●	MOS	●	●	PKN2	●	▲
AurA	●	●	Erk1 (MAPK3)	●	●	MRCka (CDC42BPA)	●	●	PKN3	●	▲
AurA (AURKA)/TPX2	▲	●	Erk2 (MAPK1)	●	●	MRCkb (CDC42BPB)	●	●	PKR (EIF2AK2)	●	●
AurB (AURKB)/INCENP	●	●	Erk5 (MAPK7)	●	●	MSK1 (RPS6KA5)	●	●	PLK1	●	●
AurC (AURKC)	●	●	Erk7 (MAPK15)	●	▲	MSK2 (RPS6KA4)	●	●	PLK2	●	●
BARK1 (ADRBK1)	●	▲	GCN2 (EIF2AK4)	●	▲	MSSK1 (STK23)	●	●	PLK3	●	●
BARK2 (ADRBK2)	●	▲	GPRK4 (GRK4)	●	▲	MST1 (STK4)	●	●	PLKX	●	●
BMPR1A (ALK3)	●	●	GPRK6 (GRK6)	●	▲	MST2 (STK3)	●	●	QIK (SNF1LK2)	●	●
BMPR1B	●	▲	GPRK7 (GRK7)	●	▲	MST3 (STK24)	●	●	RAF1	●	●
BMPR2	●	▲	GSK3a	●	●	MST4	●	●	RIPK2	●	▲
BRAF	●	●	GSK3b	●	●	MYT1 (PKMYT1)	●	▲	ROCK1	●	●
BRAF[V600E]	●	●	Haspin (GSG2)	●	●	NDR1 (STK38)	●	●	ROCK2	●	●
BRAF[T599_V600insT]	●	▲	HGK (MAP4K4)	●	●	NDR2 (STK38L)	●	●	RSK1 (RPS6KA1)	●	●
BRSK1	●	●	HH498 (TNNI3K)	●	▲	NEK1	●	●	RSK2 (RPS6KA3)	●	●
BRSK2	●	●	HIPK1	●	●	NEK2	●	●	RSK3 (RPS6KA2)	●	●
BUB1/BUB3	●	●	HIPK2	●	●	NEK3	●	▲	RSK4 (RPS6KA6)	●	●
BUBR1 (BUB1B)	●	▲	HIPK3	●	●	NEK4	●	●	SGK	●	●
CAMKK1	●	▲	HIPK4	●	●	NEK6	●	●	SGK2	●	●
CAMKK2	●	▲	HPK1 (MAP4K1)	●	▲	NEK7	●	●	SGK3	●	●
CaMK1a	●	●	HRI (EIF2AK1)	●	▲	NEK9	●	●	SIK (SNF1LK)	●	●
CaMK1b	●	▲	ICK	●	▲	NEK11	●	▲	skMLCK (MYLK2)	●	●
CaMK1d	●	●	IKKa	●	●	NIM1K(MGC42105)	●	●	smMLCK (MYLK)	●	▲
CaMK2a	●	●	IKKb	●	●	NLK	●	▲	SLK	●	●
CaMK2b	●	●	IKKe (IKBKE)	●	●	NuaK1	●	●	SRPK1	●	●
CaMK2d	●	●	IRAK1	●	●	NuaK2	●	●	SRPK2	●	●
CaMK2g	●	●	IRAK4	●	●	OSR1 (OXSR1)	●	▲	SSTK (TSSK6)	●	▲
CaMK4	●	●	JNK1 (MAPK8)	●	●	p38a (MAPK14)	●	●	STK33	●	▲
CDC2/CycB1	●	●	JNK2 (MAPK9)	●	●	p38b (MAPK11)	●	●	STK39	●	▲
CDC7/ASK	●	●	JNK3 (MAPK10)	●	●	p38g (MAPK12)	●	●	TAK1-TAB1 (MAP3K7)	●	●
CDK2/CycA2	●	●	LATS1	●	▲	p38d (MAPK13)	●	●	TAOK2	●	●
CDK2/CycE1	●	●	LATS2	●	●	p70S6K	●	●	TAOK3	●	▲
CDK3/CycE1	●	●	LIMK1	●	●	p70S6Kb	●	●	TBK1	●	●
CDK4/CycD3	●	●	LIMK2	●	▲	PAK1	●	●	TESK1	●	▲
CDK5/p25	●	●	LKB1	●	●	PAK2	●	●	TGFbR1 (TGFBFR1)	●	▲
CDK6/CycD1	●	▲	LOK	●	●	PAK3	●	▲	TGFbR2 (TGFBFR2)	●	▲
CDK6/CycD3	●	●	LRRK2	●	▲	PAK4	●	●	TNIK	●	●
CDK7/CycH/MAT1	●	●	LRRK2[G2019S]	●	▲	PAK5	●	●	TSSK1	●	●
CDK8/CycC	●	▲	LZK (MAP3K13)	●	▲	PAK6	●	●	TSSK2	●	●
CDK9/CycK	●	▲	MAP2K1	●	●	PASK	●	●	TSSK3	●	●
CDK9/CycT1	●	●	MAP2K1[F129L]	●	▲	PBK	●	●	TTBK1	●	▲
CDK11(CDC2L6)/CycC	●	▲	MAP2K1[P124L]	●	▲	PCTAIRE1 (CDK16)/CycY	●	▲	TTBK2	●	▲
CGK2 (PRKG2)	●	●	MAP2K2	●	●	PDK1 (PDPK1)	●	▲	TLK1	●	▲
CHAK1 (TRPM7)	●	▲	MAP2K3	●	●	PDHK1 (PDK1)	●	▲	TLK2	●	▲
CHK1	●	●	MAP2K4	●	●	PDHK2 (PDK2)	●	●	TTK	●	●
CHK2	●	●	MAP2K5	●	●	PDHK3 (PDK3)	●	▲	ULK3	●	▲
CK1a (CSNK1A1)	●	●	MAP2K6	●	●	PDHK4 (PDK4)	●	●	VRK1	●	▲
CK1d	●	●	MAP2K7	●	●	PEK (EIF2AK3)	●	●	VRK2	●	▲
CK1e (CSNK1E)	●	●	MAP3K1	●	●	PGK (PRKG1)	●	●	WEE1	●	●
CK1g1 (CSNK1G1)	●	●	MAP3K2	●	●	PHKG1	●	●	WNK1	●	●
CK1g2 (CSNK1G2)	●	●	MAP3K3	●	●	PHKG2	●	●	WNK2	●	●
CK1g3 (CSNK1G3)	●	●	MAP3K4	●	●	PIM1	●	●	WNK3	●	●
CK2a1/b (CSNK2A1/B)	●	●	MAP3K5	●	●	PIM2	●	●	WNK4	●	▲
CK2a2/b (CSNK2A2/B)	●	●	MAP3K6	●	▲	PIM3	●	●	YSK1	●	▲
CLK1	●	●	MAP3K14	●	▲	PKACa	●	●	ZAK	●	▲
CLK2	●	●	MAP4K2	●	●	PKACb	●	●			
CLK3	●	●	MAPKAPK2	●	●	PKACg	●	●			
CLK4	●	▲	MAPKAPK3	●	●	PKCa (PRKCA)	●	●			
COT (MAP3K8)	●	●									
CRIK	●	●									

Pack sizes are 5, 100, 200, 500 and 1000uq. Bulk production is available upon receipt of purchase order.

Inactive Kinases

Kinase	50ug/vial
Erk2 (MAPK1)	●
Erk5 (MAPK7)	●
JNK1 (MAPK8)	●
JNK2 (MAPK9)	●
MAP2K1	●
MAP2K6	●
MAP2K7	●
p38a (MAPK14)	●

Inactive Mutant Kinases

Kinase	50ug/vial
Erk2 (MAPK1)	●
Erk5 (MAPK7)	●
JNK1 (MAPK8)	●
MAP2K1	●
MAP2K6	●
MAP2K7	●
p38a (MAPK14)	●

Protein Substrates

Protein Substrate	50ug/vial
Cofilin2	●
EIF2S1	●
EIF4EBP1	●
Smad1	●
Smad3	●

To learn more, please contact:

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