

Cytoplasmic Tyrosine Kinases

	Kinase	Profiling
ABL	●	●
ABL[E255K]	●	●
ABL[T315I]	●	●
ACK	●	●
ARG	●	●
BLK	●	●
BMX	●	●
BRK	●	●
BTK	●	●
CSK	●	●
FAK	●	●
FER	●	●
FES	●	●
FGR	●	●
FRK	●	●
FYN	●	●
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.

Receptor Tyrosine Kinases

	Kinase	Profiling
ALK	●	●
ALK[C1156Y]	●	▲
ALK[F1174L]	●	●
ALK[G1269A]	NEW	▲
ALK[L1196M]	●	●
ALK[R1275Q]	●	●
EML4-ALK	●	●
NPM1-ALK	●	●
AXL	●	●
DDR1	●	●
DDR2	●	●
EGFR	●	●
EGFR[d746-750]	●	●
EGFR[d746-750/T790M]	●	●
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]	●	▲
FGFR3	●	●
FGFR3[G697C]	●	▲
FGFR3[K650E]	●	●
FGFR3[K650M]	●	●
FGFR3[V555M]	●	▲
FGFR4	●	●
FGFR4[N535K]	●	●
FGFR4[V550E]	●	●
FGFR4[V550L]	●	●

	Kinase	Profiling
FLT1	●	●
FLT3	●	●
FLT4	●	●
FMS	●	●
HER2	●	●
HER4	●	●
IGF1R	●	●
INSR	●	●
IRR	●	●
KDR	●	●
KIT	●	●
KIT[D816V]	●	●
KIT[D816Y]	●	▲
KIT[T670I]	●	●
KIT[V560G]	●	●
KIT[V560G/D816V]	●	▲
KIT[V654A]	▲	●
LTK	●	●
MER	●	●
MET	●	●
MET[D1228H]	NEW	▲
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	●	●

Lipid Kinases

	Kinase	Profiling	
PIK3CA/PIK3R1	●	●	20ug/vial
PIK3CB/PIK3R1	●	▲	20ug/vial
PIK3CD/PIK3R1	NEW	▲	20ug/vial
SPHK1	●	●	5ug/vial
SPHK2	▲	●	---

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 - AXL	●	10, 100ug/vial
BTN - BMX	●	10, 100ug/vial
BTN - BTK	●	10, 100ug/vial
BTN - CHK1 (CHEK1)	●	10, 100ug/vial
BTN - DDR2	NEW	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 - FAK (PTK2)	●	10, 100ug/vial
BTN - FLT3	●	10, 100ug/vial
BTN - FMS (CSF1R)	●	10, 100ug/vial
BTN - FYN	●	10, 100ug/vial
BTN - HER2 (ERBB2)	●	10, 100ug/vial
BTN - IGF1R	●	10, 100ug/vial
BTN - IKKa (CHUK)	●	10, 100ug/vial
BTN - IKKb	●	10, 100ug/vial
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 - LCK	●	10, 100ug/vial
BTN - LYNa	●	10, 100ug/vial
BTN - MAP2K1	●	10, 100ug/vial
BTN - MAP3K5	●	10, 100ug/vial
BTN - MET	●	10, 100ug/vial
BTN - PAK4	●	10, 100ug/vial
BTN - PIK3CA/PIK3R1	●	10, 100ug/vial
BTN - PKN1	●	10, 100ug/vial

	Kinase	
BTN - PLK1	●	10, 100ug/vial
BTN - PLK4	●	10, 100ug/vial
BTN - PYK2 (PTK2B)	●	10, 100ug/vial
BTN - RET	●	10, 100ug/vial
BTN - ROCK1	●	10, 100ug/vial
BTN - SPHK1	●	10, 100ug/vial
BTN - SRC	●	10, 100ug/vial
BTN - SYK	●	10, 100ug/vial
BTN - TGFbR1 (TGFB1)	●	10, 100ug/vial
BTN - TIE2	●	10, 100ug/vial
BTN - TNIK	●	10, 100ug/vial
BTN - WNK1	●	10, 100ug/vial
BTN - WNK4	●	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

Biotinylated Kinases
PRICE REDUCED
as of October 1, 2012!

	Protein	Profiling (Km app.)	Profiling (1mM)
Cytoplasmic Tyrosine	32	34	34
Receptor Tyrosine	46	73	73
Serine/Threonine	260	201	59
Lipid	4	3	0
Biotinylated	342	311	166

Updated: 2012/10/1

Kinase

- Available
- Available as 5ug only
- ▲ Under development
- Specification change

Profiling

- Available at [ATP] Km app.
- Available at both [ATP] Km app. and 1mM
- ▲ Under development

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
ACTR2A	● ▲	DLK (MAP3K12)	● ●	MARK3	● ●	PKCb2 (PRKCB2)	● ●
ACVR2B	● ▲	DMPK1 (DMPK)	● ▲	MARK4	● ●	PKCg (PRKCG)	● ●
AKT1	● ●	DMPK2 (CDC42BPG)	● ▲	MELK	● ●	PKCd (PRKCD)	● ●
AKT2	● ●	DRAK1 (STK17A)	● ▲	MGC42105	● ●	PKCe (PRKCE)	● ●
AKT3	● ●	DYRK1A	● ●	MINK (MINK1)	● ●	PKCz (PRKCZ)	● ●
ALK2 (ACVR1)	● ▲	DYRK1B	● ●	MLK1 (MAP3K9)	● ●	PKCh (PRKCH)	● ●
ALK4 (ACVR1B)	● ▲	DYRK2	● ●	MLK2 (MAP3K10)	● ●	PKCq (PRKCO)	● ●
AMPKa1/b1/g1 (PRKAA1/B1/G1)	● ●	DYRK3	● ●	MLK3 (MAP3K11)	● ●	PKCi (PRKCI)	● ●
AMPKa2/b1/g1 (PRKAA2/B1/G1)	● ●	DYRK4	● ▲	MLK4	● ▲	PKD1 (PRKD1)	● ●
AurA	● ●	EEF2K	● ●	MNK1 (MKNK1)	● ●	PKD2 (PRKD2)	● ●
AurA (AURKA)/TPX2	● ●	Erk1 (MAPK3)	● ●	MNK2 (MKNK2)	● ●	PKD3 (PRKD3)	● ●
AurB (AURKB)/INCENP	● ●	Erk2 (MAPK1)	● ●	MOS	● ●	PKN1	● ●
AurC (AURKC)	● ●	Erk5 (MAPK7)	● ●	MRCKa (CDC42BPA)	● ●	PKN2	● ▲
BARK1 (ADRBK1)	● ▲	Erk7 (MAPK15)	● ▲	MRCKb (CDC42BPB)	● ●	PKN3	● ▲
BARK2 (ADRBK2)	● ▲	GCN2 (EIF2AK4)	● ▲	MSK1 (RPS6KA5)	● ●	PKR (EIF2AK2)	● ●
BMPR1A (ALK3)	● ●	GPRK4 (GRK4)	● ▲	MSK2 (RPS6KA4)	● ●	PLK1	● ●
BMPR1B	● ▲	GPRK6 (GRK6)	● ▲	MSSK1 (STK23)	● ●	PLK2	● ●
BMPR2	● ▲	GPRK7 (GRK7)	● ▲	MST1 (STK4)	● ●	PLK3	● ●
BRAF	● ●	GSK3a	● ●	MST2 (STK3)	● ●	PLK4	● ●
BRAF[V600E]	● ●	GSK3b	● ●	MST3 (STK24)	● ●	PRKX	● ●
BRAF[T599_V600InsT]	● ▲	Haspin (GSG2)	● ●	MST4	● ●	QIK (SNF1LK2)	● ●
BRSK1	● ●	HGK (MAP4K4)	● ●	MYT1 (PKMYT1)	● ▲	RAF1	● ●
BRSK2	● ●	HH498 (TNNI3K)	● ▲	NDR1 (STK38)	● ●	RIPK2	● ▲
BUBR1 (BUB1B)	● ▲	HIPK1	● ●	NDR2 (STK38L)	● ●	ROCK1	● ●
CAMKK1	● ▲	HIPK2	● ●	NEK1	● ●	ROCK2	● ●
CAMKK2	● ▲	HIPK3	● ●	NEK2	● ●	RSK1 (RPS6KA1)	● ●
CaMK1a	● ●	HIPK4	● ●	NEK3	● ▲	RSK2 (RPS6KA2)	● ●
CaMK1b	● ▲	HRI (EIF2AK1)	● ▲	NEK4	● ●	RSK3 (RPS6KA2)	● ●
CaMK1d	● ●	ICK	● ▲	NEK6	● ●	RSK4 (RPS6KA6)	● ●
CaMK2a	● ●	IKKa	● ●	NEK7	● ●	SGK	● ●
CaMK2b	● ●	IKKb	● ●	NEK9	● ●	SGK2	● ●
CaMK2d	● ●	IKKe (IKBKE)	● ●	NEK11	● ▲	SGK3	● ●
CaMK2g	● ●	IRAK1	● ●	NLK	● ▲	SIK (SNF1LK)	● ●
CaMK4	● ●	IRAK4	● ●	NuaK1	● ●	skMLCK (MYLK2)	● ●
CDC2/CycB1	● ●	JNK1 (MAPK8)	● ●	NuaK2	● ●	smMLCK (MYLK)	● ▲
CDC7/ASK	● ●	JNK2 (MAPK9)	● ●	OSR1 (OXSR1)	● ▲	SLK	● ●
CDK2/CycA2	● ●	JNK3 (MAPK10)	● ●	p38a (MAPK14)	● ●	SRPK1	● ●
CDK2/CycE1	● ●	LATS1	● ▲	p38b (MAPK11)	● ●	SRPK2	● ●
CDK3/CycE1	● ●	LATS2	● ●	p38g (MAPK12)	● ●	SSTK (TSSK6)	● ▲
CDK4/CycD3	● ●	LIMK1	● ●	p38d (MAPK13)	● ●	STK33	● ▲
CDK5/p25	● ●	LIMK2	● ▲	p70S6K	● ●	STLK3 (STK39)	● ▲
CDK6/CycD3	● ●	LKB1	● ●	p70S6Kb	● ●	TAK1-TAB1 (MAP3K7)	● ●
CDK7/CycH/MAT1	● ●	LOK	● ●	PAK1	● ●	TAOK2	● ●
CDK8/CycC	● ▲	LRRK2	● ▲	PAK2	● ●	TAOK3	● ▲
CDK9/CycT1	● ●	LRRK2[G2019S]	● ▲	PAK3	● ●	TBK1	● ●
CDK11(CDC2L6)/CycC	● ▲	LZK (MAP3K13)	● ▲	PAK4	● ●	TESK1	● ▲
CGK2 (PRKG2)	● ●	MAP2K1	● ●	PAK5	● ●	TGFbR1 (TGFB1)	● ▲
CHAK1 (TRPM7)	● ▲	MAP2K1[F129L]	● ▲	PAK6	● ●	TGFbR2 (TGFB2)	● ▲
CHK1	● ●	MAP2K1[P124L]	● ▲	PASK	● ●	TNIK	● ●
CHK2	● ●	MAP2K2	● ●	PBK	● ●	TSSK1	● ●
CK1a (CSNK1A1)	● ●	MAP2K3	● ●	PCTAIRE1 (PCTK1)	● ▲	TSSK2	● ●
CK1d	● ●	MAP2K4	● ●	PDK1 (PDK1)	● ▲	TSSK3	● ●
CK1e (CSNK1E)	● ●	MAP2K5	● ●	PDHK1 (PDK1)	● ▲	TTBK1	● ▲
CK1g1 (CSNK1G1)	● ●	MAP2K6	● ●	PDHK2 (PDK2)	● ●	TTBK2	● ▲
CK1g2 (CSNK1G2)	● ●	MAP2K7	● ●	PDHK3 (PDK3)	● ▲	TLK1	● ▲
CK1g3 (CSNK1G3)	● ●	MAP3K1	● ●	PDHK4 (PDK4)	● ●	TLK2	● ▲
CK2a1/b (CSNK2A1/B)	● ●	MAP3K2	● ●	PEK (EIF2AK3)	● ●	TTK	● ●
CK2a2/b (CSNK2A2/B)	● ●	MAP3K3	● ●	PGK (PRKG1)	● ●	ULK3	● ●
CLK1	● ●	MAP3K4	● ●	PHKG1	● ●	VRK1	● ▲
CLK2	● ●	MAP3K5	● ●	PHKG2	● ●	VRK2	● ▲
CLK3	● ●	MAP3K6	● ▲	PIM1	● ●	WEE1	● ●
CLK4	● ▲	MAP3K14	● ▲	PIM2	● ●	WNK1	● ●
COT (MAP3K8)	● ●	MAP4K2	● ●	PIM3	● ●	WNK2	● ●
CRIK	● ●	MAPKAPK2	● ●	PKACa	● ●	WNK3	● ●
DAPK1	● ●	MAPKAPK3	● ●	PKACb	● ●	WNK4	● ▲
DAPK3	● ▲	MAPKAPK5	● ●	PKACg	● ●	YSK1	● ▲
DCAMKL1	● ▲	MARK1	● ●	PKCa (PRKCA)	● ●	ZAK	● ▲
DCAMKL2	● ●	MARK2	● ●	PKCb1 (PRKCB1)	● ●		

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

Inactive Kinases

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

Inactive Mutant Kinases

Kinase	50ug/vial
Erk2	●
Erk5	●
JNK1	●
MAP2K1	●
MAP2K6	●
MAP2K7	●
p38a	●

Protein Substrates

Protein Substrate	50ug/vial
Cofilin2	●
EIF2S1	●
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