

Kinase Products from Carna Biosciences, Inc.

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	●	●
LYN α	●	●
LYN β	●	●
PYK2	●	●
SRC	●	●
SRM	●	●
SYK	●	●
TEC	●	●
TNK1	●	●
TXK	●	●
TYK2	●	●
YES	●	●
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[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	●	●
FGFR2	●	●
FGFR3	●	●
FGFR3[K650E]	●	●
FGFR3[K650M]	●	●
FGFR4	●	●
FGFR4[N535K]	●	●
FGFR4[V550E]	●	●
FGFR4[V550L]	●	●

	Kinase	Profiling
FLT1	●	●
FLT3	●	●
FLT4	●	●
FMS	●	●
HER2	●	●
HER4	●	●
IGF1R	●	●
INSR	●	●
IRR	●	●
KDR	●	●
KIT	●	●
KIT[D816V]	●	●
KIT[T670I]	●	●
KIT[V560G]	●	●
KIT[V654A]	▲	●
LTK	●	●
MER	●	●
MET	●	●
MET[Y1235D]	●	●
MUSK	●	●
PDGFR α	●	●
PDGFR α [D842V]	●	▲
PDGFR α [T674I]	●	●
PDGFR α [V561D]	●	●
PDGFR β	●	●
RET	●	●
RET[G691S]	●	●
RET[M918T]	●	●
RET[S891A]	●	●
RET[Y791F]	●	●
RON	●	●
ROR1	●	▲
ROR2	●	▲
ROS	●	●
TIE2	●	●
TRKA	●	●
TRKB	●	●
TRKC	●	●
TYRO3	●	●

Lipid Kinases

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

Inactive Kinases

	Kinase	
Erk2	●	50ug/vial
Erk5	●	50ug/vial
JNK1	●	50ug/vial
JNK2 (MAPK9)	●	50ug/vial
MAP2K1	●	50ug/vial
MAP2K6	●	50ug/vial
MAP2K7	●	50ug/vial
p38 α	●	50ug/vial

Inactive Mutant Kinases

	Kinase	
Erk2	●	50ug/vial
Erk5	●	50ug/vial
JNK1	●	50ug/vial
MAP2K1	●	50ug/vial
MAP2K6	●	50ug/vial
MAP2K7	●	50ug/vial
p38 α	●	50ug/vial

Protein Substrates

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

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 - BTK	●	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 - 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 - IKK α (CHUK)	●	10, 100ug/vial
BTN - IKK α (CHUK) [inactive mutant]	●	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 - LYN α	●	10, 100ug/vial
BTN - MAP2K1	●	10, 100ug/vial
BTN - MET	●	10, 100ug/vial
BTN - PAK4	●	10, 100ug/vial
BTN - PIK3CA/PIK3R1	●	10, 100ug/vial
BTN - PLK1	●	10, 100ug/vial
BTN - PLK4	●	10, 100ug/vial
BTN - PYK2 (PTK2B)	●	10, 100ug/vial
BTN - RET	●	10, 100ug/vial
BTN - SPHK1	●	10, 100ug/vial
BTN - SRC	●	10, 100ug/vial
BTN - SYK	●	10, 100ug/vial
BTN - TGF β R1 (TGFB β R1)	●	10, 100ug/vial

	Kinase	
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

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

	Protein	Profiling (Km app.)	Profiling (1mM)
Cytoplasmic Tyrosine	33	34	34
Receptor Tyrosine	48	73	73
Serine/Threonine	259	201	59
Lipid	3	3	0
	343	311	166
Biotinylated	47		

To learn more, please contact:



Carna Biosciences, Inc.
BMA 3F, 1-5-5 Minatojima-Minami, Chuo, Kobe 650-0047 JAPAN
www.carnabio.com

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

To place an order :

Phone : +81-78-302-7091
FAX : +81-78-302-7086

To inquire :

E-mail : info@carnabio.com

Updated: 2012/2/1

Kinase Products from Carna Biosciences, Inc.

Serine/Threonine Kinases

	Kinase	Profiling		Kinase	Profiling		Kinase	Profiling	
ACTR2A	●	▲	HH498 (TNNI3K)	NEW	●	▲	PAK6	●	●
ACVR2B	●	▲	HIPK1	●	●	PASK	●	●	
AKT1	●	●	HIPK2	●	●	PBK	●	●	
AKT2	●	●	HIPK3	●	●	PCTAIRE1 (PCTK1)	●	▲	
AKT3	●	●	HIPK4	●	●	PKI1 (PDPK1)	▲	●	
ALK2 (ACVR1)	●	▲	HRI (EIF2AK1)	●	▲	PDHK1 (PDK1)	●	▲	
ALK4 (ACVR1B)	●	▲	ICK	●	▲	PDHK2 (PDK2)	●	●	
AMPKα1/β1/γ1(PRKAA1/B1/G1)	●	●	IKKα	●	●	PDHK3 (PDK3)	●	▲	
AMPKα2/β1/γ1(PRKAA2/B1/G1)	●	●	IKKβ	●	●	PDHK4 (PDK4)	●	●	
AurA	●	●	IKKε (IKBKE)	●	●	PEK (EIF2AK3)	●	●	
AurA (AURKA)/TPX2	▲	●	IRAK1	●	●	PGK (PRKG1)	●	●	
AurB (AURKB)/INCENP	●	●	IRAK4	●	●	PHKG1	●	●	
AurC (AURKC)	●	●	JNK1 (MAPK8)	●	●	PHKG2	●	●	
BARK1 (ADRBK1)	●	▲	JNK2 (MAPK9)	●	●	PIM1	●	●	
BARK2 (ADRBK2)	●	▲	JNK3 (MAPK10)	●	●	PIM2	●	●	
BMPR1A (ALK3)	●	●	LATS1	●	▲	PIM3	●	●	
BMPR1B	●	▲	LATS2	●	●	PKACα	●	●	
BMPR2	●	▲	LIMK1	●	●	PKACβ	●	●	
BRAF	●	●	LIMK2	●	▲	PKACγ	●	●	
BRAF[V600E]	●	●	LKB1	●	●	PKCα (PRKCA)	●	●	
BRAF[T599_V600insT]	●	▲	LOK	●	●	PKCβ1 (PRKCB1)	●	●	
BRSK1	●	●	LRRK2	●	▲	PKCβ2 (PRKCB2)	●	●	
BRSK2	●	●	LRRK2[G2019S]	●	▲	PKCγ (PRKCG)	●	●	
BUBR1 (BUB1B)	●	▲	LZK (MAP3K13)	●	▲	PKCδ (PRKCD)	●	●	
CAMKK1	●	▲	MAP2K1	●	●	PKCε (PRKCE)	●	●	
CAMKK2	●	▲	MAP2K1[P124L]	●	▲	PKCC (PRKCZ)	●	●	
CaMK1α	●	●	MAP2K2	●	●	PKCη (PRKCH)	●	●	
CaMK1β	●	▲	MAP2K3	●	●	PKCθ (PRKCO)	●	●	
CaMK1δ	●	●	MAP2K4	●	●	PKCι (PRKCI)	●	●	
CaMK2α	●	●	MAP2K5	●	●	PKD1 (PRKD1)	●	●	
CaMK2β	●	●	MAP2K6	●	●	PKD2 (PRKD2)	●	●	
CaMK2δ	●	●	MAP2K7	●	●	PKD3 (PRKD3)	●	●	
CaMK2γ	●	●	MAP3K1	●	●	PKN1	●	●	
CaMK4	●	●	MAP3K2	●	●	PKN2	●	▲	
CDC2/CycB1	●	●	MAP3K3	●	●	PKN3	●	▲	
CDC7/ASK	●	●	MAP3K4	●	●	PKR (EIF2AK2)	●	●	
CDK2/CycA2	●	●	MAP3K5	●	●	PLK1	●	●	
CDK2/CycE1	●	●	MAP3K6	●	▲	PLK2	●	●	
CDK3/CycE1	●	●	MAP3K14	●	▲	PLK3	●	●	
CDK4/CycD3	●	●	MAP4K2	●	▲	PLK4	●	●	
CDK5/p25	●	●	MAPKAPK2	●	●	PRKX	●	●	
CDK6/CycD3	●	●	MAPKAPK3	●	●	QIK (SNF1LK2)	●	●	
CDK7/CycH/MAT1	●	●	MAPKAPK5	●	●	RAF1	●	●	
CDK8/CycC	●	▲	MARK1	●	●	RIPK2	●	▲	
CDK9/CycT1	●	●	MARK2	●	●	ROCK1	●	●	
CDK11(CDC2L6)/CycC	●	▲	MARK3	●	●	ROCK2	●	●	
CGK2 (PRKG2)	●	●	MARK4	●	●	RSK1 (RPS6KA1)	●	●	
CHAK1 (TRPM7)	●	▲	MELK	●	●	RSK2 (RPS6KA3)	●	●	
CHK1	●	●	MGC42105	●	●	RSK3 (RPS6KA2)	●	●	
CHK2	●	●	MINK (MINK1)	●	●	RSK4 (RPS6KA6)	●	●	
CK1α (CSNK1A1)	●	●	MLK1 (MAP3K9)	●	●	SGK	●	●	
CK1δ	●	●	MLK2 (MAP3K10)	●	●	SGK2	●	●	
CK1ε (CSNK1E)	●	●	MLK3 (MAP3K11)	●	●	SGK3	●	●	
CK1γ1 (CSNK1G1)	●	●	MNK1 (MKNK1)	●	●	SIK (SNF1LK)	●	●	
CK1γ2 (CSNK1G2)	●	●	MNK2 (MKNK2)	●	●	skMLCK (MYLK2)	●	●	
CK1γ3 (CSNK1G3)	●	●	MOS	●	●	smMLCK (MYLK)	●	▲	
CK2α1/β (CSNK2A1/B)	●	●	MRCKα (CDC42BPA)	●	●	SLK	●	●	
CK2α2/β (CSNK2A2/B)	●	●	MRCKβ (CDC42BPB)	●	●	SRPK1	●	●	
CLK1	●	●	MSK1 (RPS6KA5)	▲	●	SRPK2	●	●	
CLK2	●	●	MSK2 (RPS6KA4)	●	●	SSTK (TSSK6)	●	▲	
CLK3	●	●	MSSK1 (STK23)	●	●	STK33	●	▲	
CLK4	●	▲	MST1 (STK4)	●	●	STLK3 (STK39)	●	▲	
COT (MAP3K8)	●	●	MST2 (STK3)	●	●	TAK1-TAB1 (MAP3K7)	●	●	
CRIK	●	●	MST3 (STK24)	●	●	TAOK2	●	●	
DAPK1	●	●	MST4	●	●	TAOK3	●	▲	
DAPK3	●	▲	MYT1 (PKMYT1)	●	▲	TBK1	●	●	
DCAMKL1	●	▲	NDR1 (STK38)	●	●	TESK1	●	▲	
DCAMKL2	●	●	NDR2 (STK38L)	●	●	TGFβR1 (TGFBR1)	●	▲	
DLK (MAP3K12)	●	●	NEK1	●	●	TGFβR2 (TGFBR2)	●	▲	
DMPK1 (DMPK)	●	▲	NEK2	●	●	TNIK	●	●	
DMPK2 (CDC42BPG)	●	▲	NEK3	●	▲	TSSK1	●	●	
DRAK1 (STK17A)	●	▲	NEK4	●	●	TSSK2	●	●	
DYRK1A	●	●	NEK6	●	●	TSSK3	●	●	
DYRK1B	●	●	NEK7	●	●	TTBK1	●	▲	
DYRK2	●	●	NEK9	●	●	TTBK2	●	▲	
DYRK3	●	●	NEK11	●	▲	TLK1	●	▲	
DYRK4	●	▲	NLK	●	●	TLK2	●	▲	
EEF2K	●	●	NuaK1	●	●	TTK	●	●	
Erk1 (MAPK3)	●	●	NuaK2	●	●	ULK3	●	▲	
Erk2 (MAPK1)	●	●	OSR1 (OXSR1)	●	▲	VRK1	●	▲	
Erk5 (MAPK7)	●	●	p38α (MAPK14)	●	●	VRK2	●	▲	
Erk7 (MAPK15)	●	▲	p38β (MAPK11)	●	●	WEE1	●	●	
GCN2 (EIF2AK4)	●	▲	p38γ (MAPK12)	●	●	WNK1	●	●	
GPRK4 (GRK4)	●	▲	p38δ (MAPK13)	●	●	WNK2	●	●	
GPRK6 (GRK6)	●	▲	p70S6K	●	●	WNK3	●	●	
GPRK7 (GRK7)	●	▲	p70S6Kβ	▲	●	WNK4	●	▲	
GSK3α	●	●	PAK1	●	●	YSK1	●	▲	
GSK3β	●	●	PAK2	●	●	ZAK	●	▲	
Haspin (GSG2)	●	●	PAK3	●	●				
HGK (MAP4K4)	●	●	PAK4	●	●				
			PAK5	●	●				

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