

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	●	●
PIK3CB/PIK3R1	●	▲
PIK3CD/PIK3R1	NEW	▲
SPHK1	●	●
SPHK2	▲	●

バイオチン化キナーゼタンパク質

2012年10月1日 **大幅に価格改定**を実施！
よりお求めやすくなりました！

詳しくは営業部までお気軽にお問合せください

	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
	342	311	166
Biotinylated	55		

Updated: 2012/10/1

Biotinylated Kinases

	Kinase
BTN - ABL (ABL1)	●
BTN - ACTR2B(ACVR2B)	●
BTN - AKT1	●
BTN - AKT2	●
BTN - ALK	●
BTN - ALK[L1196M]	●
BTN - AurA (AURKA)	●
BTN - AXL	●
BTN - BMX	●
BTN - BTK	●
BTN - CHK1 (CHEK1)	●
BTN - DDR2	NEW
BTN - DLK (MAP3K12)	●
BTN - EGFR	●
BTN - EGFR[T790M/L858R]	●
BTN - FGFR1	●
BTN - FGFR3	●
BTN - FAK (PTK2)	●
BTN - FLT3	●
BTN - FMS (CSF1R)	●
BTN - FYN	●
BTN - HER2 (ERBB2)	●
BTN - IGF1R	●
BTN - IKKa (CHUK)	●
BTN - IKKb	●
BTN - IRAK4	●
BTN - ITK	●
BTN - JAK1	●
BTN - JAK2	●
BTN - KDR (VEGFR2)	●
BTN - LCK	●
BTN - LYNa	●
BTN - MAP2K1	●
BTN - MAP3K5	●
BTN - MET	●
BTN - PAK4	●
BTN - PIK3CA/PIK3R1	●
BTN - PKN1	●

	Kinase
BTN - PLK1	●
BTN - PLK4	●
BTN - PYK2 (PTK2B)	●
BTN - RET	●
BTN - ROCK1	●
BTN - SPHK1	●
BTN - SRC	●
BTN - SYK	●
BTN - TGFbR1 (TGFB1R1)	●
BTN - TIE2	●
BTN - TNK1	●
BTN - WNK1	●
BTN - WNK4	●
BTN - TRKA (NTRK1)	●
BTN - TRKB (NTRK2)	●
BTN - TRKC (NTRK3)	●
BTN - TYK2	●

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

Serine/Threonine Kinases

	Kinase	Profiling
ACTR2A	●	▲
ACVR2B	●	▲
AKT1	●	●
AKT2	●	●
AKT3	●	●
ALK2 (ACVR1)	●	▲
ALK4 (ACVR1B)	●	▲
AMPKa1/b1/g1 (PRKAA1/B1/G1)	●	●
AMPKa2/b1/g1 (PRKAA2/B1/G1)	●	●
AurA	●	●
AurA (AURKA)/TPX2	▲	●
AurB (AURKB)/INCENP	●	●
AurC (AURKC)	●	●
BARK1 (ADRBK1)	●	▲
BARK2 (ADRBK2)	●	▲
BMPR1A (ALK3)	●	●
BMPR1B	●	▲
BMPR2	●	▲
BRAF	●	●
BRAF[V600E]	●	●
BRAF[T599_V600insT]	●	▲
BRSK1	●	●
BRSK2	●	●
BUBR1 (BUB1B)	●	▲
CAMKK1	●	▲
CAMKK2	●	▲
CaMK1a	●	●
CaMK1b	●	▲
CaMK1d	●	●
CaMK2a	●	●
CaMK2b	●	●
CaMK2d	●	●
CaMK2g	●	●
CaMK4	●	●
CDC2/CycB1	●	●
CDC7/ASK	●	●
CDK2/CycA2	●	●
CDK2/CycE1	●	●
CDK3/CycE1	●	●
CDK4/CycD3	●	●
CDK5/p25	●	●
CDK6/CycD3	●	●
CDK7/CycH/MAT1	●	●
CDK8/CycC	●	▲
CDK9/CycT1	●	●
CDK11(CDC2L6)/CycC	●	▲
CGK2 (PRKG2)	●	●
CHAK1 (TRPM7)	●	▲
CHK1	●	●
CHK2	●	●
CK1a (CSNK1A1)	●	●
CK1d	●	●
CK1e (CSNK1E)	●	●
CK1g1 (CSNK1G1)	●	●
CK1g2 (CSNK1G2)	●	●
CK1g3 (CSNK1G3)	●	●
CK2a1/b (CSNK2A1/B)	●	●
CK2a2/b (CSNK2A2/B)	●	●
CLK1	●	●
CLK2	●	●
CLK3	●	●
CLK4	●	▲
COT (MAP3K8)	●	●
CRIK	●	●
DAPK1	●	●
DAPK3	●	▲
DCAMKL1	●	▲
DCAMKL2	●	●

	Kinase	Profiling
DLK (MAP3K12)	●	●
DMPK1 (DMPK)	●	▲
DMPK2 (CDC42BPB)	●	▲
DRAK1 (STK17A)	●	▲
DYRK1A	●	●
DYRK1B	●	●
DYRK2	●	●
DYRK3	●	●
DYRK4	●	▲
EEF2K	●	●
Erk1 (MAPK3)	●	●
Erk2 (MAPK1)	●	●
Erk5 (MAPK7)	●	●
Erk7 (MAPK15)	●	▲
GCN2 (EIF2AK4)	●	▲
GPRK4 (GRK4)	●	▲
GPRK6 (GRK6)	●	▲
GPRK7 (GRK7)	●	▲
GSK3a	●	●
GSK3b	●	●
Haspin (GSG2)	●	●
HGK (MAP4K4)	●	●
HH498 (TNNI3K)	●	▲
HIPK1	●	●
HIPK2	●	●
HIPK3	●	●
HIPK4	●	●
HRI (EIF2AK1)	●	▲
ICK	●	▲
IKKa	●	●
IKKb	●	●
IKKe (IKBKE)	●	●
IRAK1	●	●
IRAK4	●	●
JNK1 (MAPK8)	●	●
JNK2 (MAPK9)	●	●
JNK3 (MAPK10)	●	●
LATS1	●	▲
LATS2	●	●
LIMK1	●	●
LIMK2	●	▲
LKB1	●	●
LOK	●	●
LRRK2	●	▲
LRRK2[G2019S]	●	▲
LZK (MAP3K13)	●	▲
MAP2K1	●	●
MAP2K1[F129L]	●	▲
MAP2K1[P124L]	●	▲
MAP2K2	●	●
MAP2K3	●	●
MAP2K4	●	●
MAP2K5	●	●
MAP2K6	●	●
MAP2K7	●	●
MAP3K1	●	●
MAP3K2	●	●
MAP3K3	●	●
MAP3K4	●	●
MAP3K5	●	●
MAP3K6	●	▲
MAP3K14	●	▲
MAP4K2	●	●
MAPKAPK2	●	●
MAPKAPK3	●	●
MAPKAPK5	●	●
MARK1	●	●
MARK2	●	●

	Kinase	Profiling
MARK3	●	●
MARK4	●	●
MELK	●	●
MGC42105	●	●
MINK (MINK1)	●	●
MLK1 (MAP3K9)	●	●
MLK2 (MAP3K10)	●	●
MLK3 (MAP3K11)	●	●
MLK4	●	▲
MNK1 (MKNK1)	●	●
MNK2 (MKNK2)	●	●
MOS	●	●
MRCKa (CDC42BPA)	●	●
MRCKb (CDC42BPB)	●	●
MSK1 (RPS6KA5)	▲	●
MSK2 (RPS6KA4)	●	●
MSSK1 (STK23)	●	●
MST1 (STK4)	●	●
MST2 (STK3)	●	●
MST3 (STK24)	●	●
MST4	●	●
MYT1 (PKMYT1)	●	▲
NDR1 (STK38)	●	●
NDR2 (STK38L)	●	●
NEK1	●	●
NEK2	●	●
NEK3	●	▲
NEK4	●	●
NEK6	●	●
NEK7	●	●
NEK9	●	●
NEK11	●	▲
NLK	●	▲
NuaK1	●	●
NuaK2	●	●
OSR1 (OXSR1)	●	▲
p38a (MAPK14)	●	●
p38b (MAPK11)	●	●
p38g (MAPK12)	●	●
p38d (MAPK13)	●	●
p70S6K	●	●
p70S6Kb	▲	●
PAK1	●	●
PAK2	●	●
PAK3	●	●
PAK4	●	●
PAK5	●	●
PAK6	●	●
PASK	●	●
PBK	●	●
PCTAIRE1 (PCTK1)	●	▲
PDK1 (PDPK1)	▲	●
PDHK1 (PDK1)	●	▲
PDHK2 (PDK2)	●	●
PDHK3 (PDK3)	●	▲
PDHK4 (PDK4)	●	●
PEK (EIF2AK3)	●	●
PGK (PRKG1)	●	●
PHKG1	●	●
PHKG2	●	●
PIM1	●	●
PIM2	●	●
PIM3	●	●
PKACa	●	●
PKACb	●	●
PKACg	●	●
PKCa (PRKCA)	●	●
PKCb1 (PRKCB1)	●	●

	Kinase	Profiling
PKCb2 (PRKCB2)	●	●
PKCg (PRKCG)	●	●
PKCd (PRKCD)	●	●
PKCe (PRKCE)	●	●
PKCz (PRKCZ)	●	●
PKCh (PRKCH)	●	●
PKCq (PRKCQ)	●	●
PKCi (PRKCI)	●	●
PKD1 (PRKD1)	●	●
PKD2 (PRKD2)	●	●
PKD3 (PRKD3)	●	●
PKN1	●	●
PKN2	●	▲
PKN3	●	▲
PKR (EIF2AK2)	●	●
PLK1	●	●
PLK2	●	●
PLK3	●	●
PLK4	●	●
PRKX	●	●
QIK (SNF1LK2)	●	●
RAF1	●	●
RIPK2	●	▲
ROCK1	●	●
ROCK2	●	●
RSK1 (RPS6KA1)	●	●
RSK2 (RPS6KA3)	●	●
RSK3 (RPS6KA2)	●	●
RSK4 (RPS6KA6)	●	●
SGK	●	●
SGK2	●	●
SGK3	●	●
SIK (SNF1LK)	●	●
skMLCK (MYLK2)	●	●
smMLCK (MYLK)	●	▲
SLK	●	●
SRPK1	●	●
SRPK2	●	●
SSTK (TSSK6)	●	▲
STK33	●	▲
STLK3 (STK39)	●	▲
TAK1-TAB1 (MAP3K7)	●	●
TAOK2	●	●
TAOK3	●	▲
TBK1	●	●
TESK1	●	▲
TGFbR1 (TGFB1)	●	▲
TGFbR2 (TGFB2)	●	▲
TNIK	●	●
TSSK1	●	●
TSSK2	●	●
TSSK3	●	●
TTBK1	●	▲
TTBK2	●	▲
TLK1	●	▲
TLK2	●	▲
TTK	●	●
ULK3	●	▲
VRK1	●	▲
VRK2	●	▲
WEE1	●	●
WNK1	●	●
WNK2	●	●
WNK3	●	●
WNK4	●	▲
YSK1	●	▲
ZAK	●	▲

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

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
p38a	●	50ug/vial

Inactive Mutant Kinases

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

Protein Substrates

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

To learn more, please contact:

カルナバイオサイエンス株式会社

〒650-0047 神戸市中央区港島南町1-5-5 BMA3F

TEL: 078-302-7091 (営業部直通) / FAX: 078-302-7086

E-mail: info@carnabio.com