

Kinase Profiling Services

311
validated targets!

Customizable and Pre-Selected Profiling Services to fit Your Budget and Needs!

See reverse side for available kinases!

The Carna Advantage...

- **Massive Kinome Coverage,** Activity Based Profiling & Screening.
- Confirm your results at **higher ATP concentrations** to identify potential Type II/III inhibitors.
- Build your own panel or consider our **unique Pre-Selected Panels.** BOTH offering **Time and Cost Savings!**
- Opt for Modern Technology... **Mobility Shift Assay** measures kinase activity efficiently and effectively.

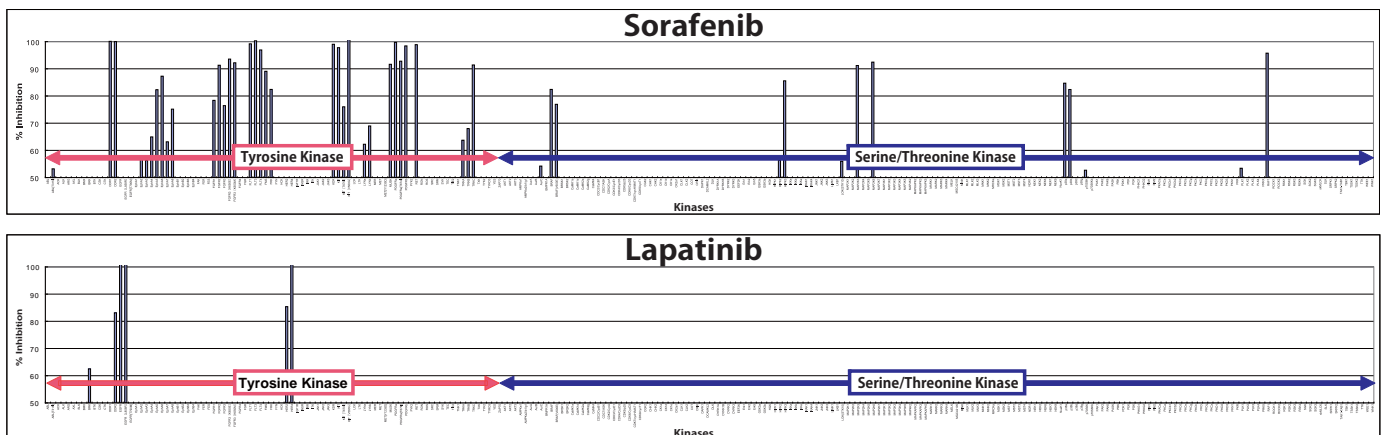
Unique Pre-Selected QuickScout™ Panel Profiling Series

TK 20 Kinases	STK 30 Kinases
MAPK 31 Kinases	Cell Cycle 30 Kinases

For more details, please request our printed brochure or visit our website.

Therapeutic Kinase Inhibitors Profiled against 256 Kinases

% Inhibition measured at 1uM against 256 kinases made in-house at Carna Biosciences (ATP conc. = Km bin.)



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

QuickScout™ Custom Profiling & Panel Profiling Series from Carna Biosciences, Inc.

Tyrosine Kinases	Custom		Panel
	Km bin.	1mM	TK
ABL(ABL1)	●	●	●
ABL(ABL1) [E255K]	●	●	●
ABL(ABL1) [T315I]	●	●	●
ACK(TNK2)	●	●	●
ALK	●	●	●
ALK [F1174L]	●	●	●
ALK [L1196M]	●	●	●
ALK [R1275Q]	●	●	●
ARG(ABL2)	●	●	●
AXL	●	●	●
BLK	●	●	●
BMX	●	●	●
BRK(PTK6)	●	●	●
BTK	●	●	●
CSK	●	●	●
DDR1	●	●	●
DDR2	●	●	●
EGFR	●	●	●
EGFR [d746-750]	●	●	●
EGFR [d746-750/T790M]	●	●	●
EGFR [L858R]	●	●	●
EGFR [L861Q]	●	●	●
EGFR [T790M]	●	●	●
EGFR [T790M/L858R]	●	●	●
EML4-ALK	●	●	●
EPHA1	●	●	●
EPHA2	●	●	●
EPHA3	●	●	●
EPHA4	●	●	●
EPHA5	●	●	●
EPHA6	●	●	●
EPHA7	●	●	●
EPHA8	●	●	●
EPHB1	●	●	●
EPHB2	●	●	●
EPHB3	●	●	●
EPHB4	●	●	●
FAK(PTK2)	●	●	●
FER	●	●	●
FES	●	●	●
FGFR1	●	●	●
FGFR2	●	●	●
FGFR3	●	●	●
FGFR3 [K650E]	●	●	●
FGFR3 [K650M]	●	●	●
FGFR4	●	●	●
FGFR4 [N535K]	●	●	●
FGFR4 [V550E]	●	●	●
FGFR4 [V550L]	●	●	●
FGR	●	●	●
FLT1	●	●	●
FLT3	●	●	●
FLT4	●	●	●
FMS(CSF1R)	●	●	●
FRK	●	●	●
FYN	●	●	●
HCK	●	●	●
HER2(ERBB2)	●	●	●
HER4(ERBB4)	●	●	●
IGF1R	●	●	●
INSR	●	●	●
IRR(INSRR)	●	●	●
ITK	●	●	●
JAK1	●	●	●
JAK2	●	●	●
JAK3	●	●	●
KDR	●	●	●
KIT	●	●	●
KIT [D816V]	●	●	●
KIT [T670I]	●	●	●
KIT [V560G]	●	●	●
KIT [V654A]	●	●	●
LCK	●	●	●
LTK	●	●	●
LYNa	●	●	●
LYNb	●	●	●
MER(MERTK)	●	●	●
MET	●	●	●
MET [Y1235D]	●	●	●
MUSK	●	●	●
NPM1-ALK	●	●	●
PDGFRa(PDGFR)	●	●	●
PDGFRa(PDGFR) [T674I]	●	●	●
PDGFRa(PDGFR) [V651D]	●	●	●
PDGFRb(PDGFRb)	●	●	●
PK2(PTK2B)	●	●	●
RET	●	●	●
RET [G691S]	●	●	●
RET [M918T]	●	●	●
RET [S891A]	●	●	●
RET [Y791F]	●	●	●
RON(MST1R)	●	●	●
ROS(ROS1)	●	●	●
SRC	●	●	●
SRM(SRMS)	●	●	●
SYK	●	●	●
TEC	●	●	●
TIE2(TEK)	●	●	●
TNK1	●	●	●
TRKA(NTRK1)	●	●	●
TRKB(NTRK2)	●	●	●
TRKC(NTRK3)	●	●	●
TYK	●	●	●
TYK2	●	●	●
TYRO3	●	●	●
YES(YES1)	●	●	●
ZAP70	●	●	●

Serine/Threonine Kinases	Custom		Panel		
	Km bin.	1mM	STK	MAPK	Cell
AKT1	●	●	●		
AKT2	●	●	●		
AKT3	●	●	●		
AMPKa1/β1/γ1(PRKAA1/B1/G1)	●	●	●		
AMPKa2/β1/γ1(PRKAA2/B1/G1)	●	●	●		
AurA(AURKA)	●	●	●		●
AurA(AURKA)/TPX2	●	●	●		●
AurB(AURKB)/INCENP	●	●	●		●
AurC(AURKC)	●	●	●		●
BMPRI1A	●	●	●		
BRAF	●	●	●	●	
BRAF [V600E]	●	●	●	●	
BRSK1	●	●	●		
BRSK2	●	●	●		●
CaMK1a(CAMK1)	●	●	●		
CaMK1δ(CAMK1D)	●	●	●		
CaMK2a(CAMK2A)	●	●	●		
CaMK2β(CAMK2B)	●	●	●		
CaMK2γ(CAMK2G)	●	●	●		
CaMK2δ(CAMK2D)	●	●	●		
CaMK4	●	●	●		
CDC2/CycB1	●	●	●		●
CDC7/ASK	●	●	●		●
CDK2/CycA2	●	●	●		●
CDK2/CycE1	●	●	●		●
CDK3/CycE1	●	●	●		●
CDK4/CycD3	●	●	●		●
CDK5/p25	●	●	●		●
CDK6/CycD3	●	●	●		●
CDK7/CycH/MAT1	●	●	●		●
CDK9/CycT1	●	●	●		●
CGK2(PRKG2)	●	●	●		●
CHK1(CHEK1)	●	●	●		●
CHK2(CHEK2)	●	●	●		●
CK1a(CSNK1A1)	●	●	●		●
CK1γ1(CSNK1G1)	●	●	●		●
CK1γ2(CSNK1G2)	●	●	●		●
CK1γ3(CSNK1G3)	●	●	●		●
CK1δ(CSNK1D)	●	●	●		●
CK1ε(CSNK1E)	●	●	●		●
CK2a1/β(CSNK2A1/B)	●	●	●		●
CK2a2/β(CSNK2A2/B)	●	●	●		●
CLK1	●	●	●		
CLK2	●	●	●		
CLK3	●	●	●		
COT(MAP3K8)	●	●	●	●	
CRIK(CIT)	●	●	●	●	
DAPK1	●	●	●		
DCAMK12	●	●	●		
DLK(MAP3K12)	●	●	●	●	
DYRK1A	●	●	●		
DYRK1B	●	●	●		
DYRK2	●	●	●		
DYRK3	●	●	●		
EEF2K	●	●	●	●	
Erk1(MAPK3)	●	●	●	●	
Erk2(MAPK1)	●	●	●	●	
Erk5(MAPK7)	●	●	●	●	
GSK3a(GSK3A)	●	●	●		●
GSK3β(GSK3B)	●	●	●		●
Haspin(GSG2)	●	●	●		
HGK(MAP4K4)	●	●	●		
HIPK1	●	●	●		
HIPK2	●	●	●		
HIPK3	●	●	●		
HIPK4	●	●	●		
IKKα(CHUK)	●	●	●		
IKKβ(IKBK)	●	●	●		
IKKε(IKBKE)	●	●	●		
IRAK1	●	●	●		
IRAK4	●	●	●		
JNK1(MAPK8)	●	●	●	●	
JNK2(MAPK9)	●	●	●	●	
JNK3(MAPK10)	●	●	●	●	
LATS2	●	●	●		
LIMK1	●	●	●		
LKB1(STK11)/MO25a/STRADA	●	●	●		
LOK(STK10)	●	●	●		
MAP2K1	●	●	●	●	
MAP2K2	●	●	●	●	
MAP2K3	●	●	●	●	
MAP2K4	●	●	●	●	
MAP2K5	●	●	●	●	
MAP2K6	●	●	●	●	
MAP2K7	●	●	●	●	
MAP3K1	●	●	●	●	
MAP3K2	●	●	●	●	
MAP3K3	●	●	●	●	
MAP3K4	●	●	●	●	
MAP3K5	●	●	●	●	
MAP4K2	●	●	●	●	
MAPKAPK2	●	●	●	●	
MAPKAPK3	●	●	●	●	
MAPKAPK5	●	●	●	●	
MARK1	●	●	●		
MARK2	●	●	●		
MARK3	●	●	●		
MARK4	●	●	●		
MELK	●	●	●		
MGC42105	●	●	●		
MINK(MINK1)	●	●	●		

Custom Profiling	Km bin.	311
	1mM	166
Panel Profiling	TK20	20
	STK30	30
	MAPK	31
	Cell Cycle	30

Serine/Threonine Kinases	Custom		Panel		
	Km bin.	1mM	STK	MAPK	Cell
MLK1(MAP3K9)	●	●	●	●	
MLK2(MAP3K10)	●	●	●	●	
MLK3(MAP3K11)	●	●	●	●	
MNK1(MKNK1)	●	●	●		
MNK2(MKNK2)	●	●	●		
MOS	●	●	●	●	
MRCkα(CDC42BPA)	●	●	●		
MRCkβ(CDC42BPB)	●	●	●		
MSK1(RPS6KA5)	●	●	●		
MSK2(RPS6KA4)	●	●	●		
MSSK1(STK23)	●	●	●		
MST1(STK4)	●	●	●		
MST2(STK3)	●	●	●		
MST3(STK24)	●	●	●		
MST4	●	●	●		
NDR1(STK38)	●	●	●		
NDR2(STK38L)	●	●	●		
NEK1	●	●	●		●
NEK2	●	●	●		●
NEK4	●	●	●		
NEK6	●	●	●		●
NEK7	●	●	●		●
NEK9	●	●	●		●
Nuak1	●	●	●		
Nuak2	●	●	●		
p38α(MAPK14)	●	●	●	●	●
p38β(MAPK11)	●	●	●	●	
p38γ(MAPK12)	●	●	●	●	
p38δ(MAPK13)	●	●	●	●	
p70S6K(RPS6KB1)	●	●	●		●
p70S6Kβ(RPS6KB2)	●	●	●		
PAK1	●	●	●		
PAK2	●	●	●		
PAK3	●	●	●		
PAK4	●	●	●		
PAK5(PAK7)	●	●	●		
PAK6	●	●	●		
PASK	●	●	●		
PBK	●	●	●		
PDHK2(PDK2)	●	●	●		
PDHK4(PDK4)	●	●	●		
PDK1(PDPK1)	●	●	●		
PEK(EIF2AK3)	●	●	●		
PGK(PRKG1)	●	●	●		
PHKG1	●	●	●		
PHKG2	●	●	●		
PIM1	●	●	●		●
PIM2	●	●	●		●
PIM3	●	●	●		
PKAcα(PRKACA)	●	●	●		
PKAcβ(PRKACB)	●	●	●		
PKAcγ(PRKACG)	●	●	●		
PKCα(PRKCA)	●	●	●		
PKCβ1(PRKCβ1)	●	●	●		
PKCβ2(PRKCβ2)	●	●	●		
PKCγ(PRKCγ)	●	●	●		
PKCδ(PRKCD)	●	●	●		
PKCε(PRKCE)	●	●	●		
PKCζ(PRK CZ)	●	●	●		
PKCη(PRKCH)	●	●	●		
PKCθ(PRK Cθ)	●	●	●		
PKCi(PRKCI)	●	●	●		
PKD1(PRKD1)	●	●	●		
PKD2(PRKD2)	●	●	●		
PKD3(PRKD3)	●	●	●		
PKN1	●	●	●		
PKR(EIF2AK2)	●	●	●		
PLK1	●	●	●		●
PLK2	●	●	●		●
PLK3	●	●	●		
PLK4	●	●	●		
PRXK	●	●	●		
QIK(SNF1LK2)	●	●	●		
RAF1	●	●	●	●	
ROCK1	●	●	●		
ROCK2	●	●	●		
RSK1(RPS6KA1)	●	●	●		●
RSK2(RPS6KA3)	●	●	●		
RSK3(RPS6KA2)	●	●	●		
RSK4(RPS6KA6)	●	●	●		
SGK	●	●	●		
SGK2	●	●	●		
SGK3(SGKL)	●	●	●		
SIK(SNF1LK)	●	●	●		
skMLCK(MYLK2)	●	●	●		
SLK	●	●	●		
SRPK1	●	●	●		
SRPK2	●	●	●		
TAK1-TAB1(MAP3K7)	●	●	●	●	
TAOK2	●	●	●		
TBK1	●	●	●		
TNIK	●	●	●		
TSSK1	●	●	●		
TSSK2	●	●	●	</	