

Accelerate Discovery! Benefit from the assay expertise of Carna Biosciences...

QuickScout Screening Assist™ Kits

QuickScout Screening Assist™ Kits are designed to speed you through compound screening, particularly secondary- and counter- screening operations, by providing necessary reagents and detailed assay protocols, for more than 300 human kinases. Each kinase kit is made-to-order, with turnaround time of 2-3 weeks.



Advantages of Carna's Assay Kits...

- Prepared utilizing the extensive expertise of our profiling services
- Ready-To-Run products & protocols
- One kit allows you multiple assays

Designed for primary, in-house screening procedures
- Lead Generation through Lead Optimization!

Total
331
Kinases*

* As of April 14, 2015

Platform	Minimum Size	Available Target	Kit Components
<p>Mobility Shift Assay QSS Assist™ MSA</p> <p>This MSA kit works best using LabChip® technology from PerkinElmer, Inc.</p>	<p>400dp Equivalent to 1 x 384-well plate</p>	<p>287 Kinases</p>	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture (ATP, Metal included) ● Assay Buffer ● Termination Buffer ● Assay Protocol (Separation conditions included) <p>* Protocol sample is available online.</p> <p>After initial kit purchase, components may be purchased separately. Minimum order requirements apply.</p> <p>* The dilution ratio and MSA measuring parameters are different by kinase.</p>
<p>FP(IMAP™) QSS Assist™ FP</p>	<p>800dp Equivalent to 2 x 384-well plate</p>	<p>79 Kinases</p>	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture ● Assay Buffer ● Assay Protocol
<p>TR-FRET QSS Assist™ TR-FRET</p>	<p>800dp Equivalent to 2 x 384-well plate</p>	<p>28 Kinases</p>	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture ● Assay Buffer ● Assay Protocol
<p>ELISA QSS Assist™ ELISA</p>	<p>500dp Equivalent to 5 x 96-well plate</p>	<p>35 Kinases</p>	<ul style="list-style-type: none"> ● Protein Kinase ● Substrate Mixture ● Assay Buffer ● Antibody for ELISA (except for TTK & WEE1) ● Assay Protocol

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

Tyrosine Kinases	MSA	TR-FRET
ABL(ABL1)	●	
ABL(ABL1) [E255K]	●	
ABL(ABL1) [T315I]	●	
ACK(TNK2)	●	
ALK	●	●
ALK [C1156Y]	●	
ALK [F1174L]	●	
ALK [G1202R]	●	
ALK [G1269A]	●	
ALK [L1196M]	●	
ALK [R1275Q]	●	
ALK [T1151_L1152insT]	●	
EML4-ALK	●	
NPM1-ALK	●	
ALK2(ACVR1)	●	●
ALK4(ACVR1B)	●	●
ARG(ABL2)	●	●
AXL	●	●
BLK	●	
BMX	●	
BRK(PTK6)	●	
BTK	●	●
BTK [C481S]	●	
CSK	●	●
DDR1	●	
DDR2	●	
EGFR	●	
EGFR [d746-750]	●	
EGFR [d746-750/T790M]	●	
EGFR [L858R]	●	
EGFR [L861Q]	●	
EGFR [T790M/L858R]	●	
EGFR [T790M]	●	
EPHA1	●	
EPHA2	●	●
EPHA3	●	
EPHA4	●	
EPHA5	●	
EPHA6	●	
EPHA7	●	
EPHA8	●	
EPHB1	●	
EPHB2	●	
EPHB3	●	
EPHB4	●	
FAK(PTK2)	●	●
FER	●	
FES	●	●
FGFR1	●	
FGFR1 [V561M]	●	
FGFR2	●	
FGFR2 [N549H]	●	
FGFR3	●	●
FGFR3 [G697C]	●	
FGFR3 [K650E]	●	
FGFR3 [K650M]	●	
FGFR3 [V555M]	●	
FGFR4	●	
FGFR4 [N535K]	●	
FGFR4 [V550E]	●	
FGFR4 [V550L]	●	
FGR	●	
FLT1	●	
FLT3	●	●
FLT4	●	
FMS(CSF1R)	●	
FRK	●	
FYN[isoform a]	●	
FYN[isoform b]	●	
HCK	●	●
HER2(ERBB2)	●	
HER4(ERBB4)	●	
IGF1R	●	●
INSR	●	
IRR(INSSRR)	●	
ITK	●	
JAK1	●	●
JAK2	●	●
JAK2(JH1 JH2)	●	
JAK2(JH1 JH2) [V617F]	●	
JAK3	●	●
KDR	●	
KIT	●	
KIT [D816E]	●	
KIT [D816V]	●	
KIT [D816Y]	●	
KIT [T670I]	●	
KIT [V560G]	●	
KIT [V560G/D816V]	●	
KIT [V654A]	●	
LCK	●	●
LTK	●	
LYNa	●	●
LYNb	●	
MER(MERTK)	●	●
MET	●	●
MET [D1228H]	●	
MET [M1250T]	●	
MET [Y1235D]	●	
MJUSK	●	
PDGFRa(PDGFR)	●	
PDGFRa(PDGFR) [D842V]	●	
PDGFRa(PDGFR) [T674I]	●	
PDGFRa(PDGFR) [V561D]	●	
PDGFRb(PDGFRB)	●	●
PYK2(PTK2B)	●	
RET	●	
RET [G691S]	●	
RET [M918T]	●	
RET [S891A]	●	
RET [Y791F]	●	
RON(MST1R)	●	
ROS(ROS1)	●	
SRC	●	
SRM(SRMS)	●	
SYK	●	●
TEC	●	

Tyrosine Kinases	MSA	TR-FRET
TIE2(TEK)	●	●
TNK1	●	
TRKA(NTRK1)	●	●
TRKB(NTRK2)	●	
TRKC(NTRK3)	●	
TXK	●	
TYK2	●	●
TYRO3	●	
YES(YES1)	●	
YES(YES1) [T348I]	●	

Serine/Threonine Kinases	MSA	FP (IMAP™)	ELISA	TR-FRET
ACVR2A	●			●
AKT1	●	●		
AKT2	●	●		
AKT3	●	●		
AMPKα1/β1/γ1(PRKAA1/B1/G1)	●			
AMPKα2/β1/γ1(PRKAA2/B1/G1)	●			
AurA(AURKA)	●	●		
AurA(AURKA)/TPX2	●	●		
AurB(AURKB)/INCENP	●	●		
AurC(AURKC)	●	●		
BMPR1A	●		●	
BMPR1B	●			●
BRAF	●		●	
BRAF [V600E]	●		●	
BRSK1	●			
BRSK2	●	●		
CaMK1α(CAMK1)	●	●		
CaMK1δ(CAMK1D)	●	●		
CaMK2α(CAMK2A)	●	●		
CaMK2β(CAMK2B)	●	●		
CaMK2δ(CAMK2D)	●	●		
CaMK2γ(CAMK2G)	●	●		
CaMK4	●	●		
CDC2/CycB1	●	●		
CDC2L6/CycC	●		●	
CDC7/ASK	●		●	
CDK2/CycA2	●	●		
CDK2/CycE1	●	●		
CDK3/CycE1	●	●		
CDK4/CycD3	●	●		
CDK5/p25	●	●		
CDK6/CycD3	●	●		
CDK7/CycH/MAT1	●	●		
CDK8/CycC	●		●	
CDK9/CycT1	●	●		
CGK2(PRK2)	●	●		
CHK1(CHEK1)	●	●		
CHK2(CHEK2)	●	●		
CK1α(CSNK1A1)	●	●		
CK1δ(CSNK1D)	●	●		
CK1ε(CSNK1E)	●	●		
CK1γ1(CSNK1G1)	●	●		
CK1γ2(CSNK1G2)	●	●		
CK1γ3(CSNK1G3)	●	●		
CK2α1/β(CSNK2A1/B)	●	●		
CK2α2/β(CSNK2A2/B)	●	●		
CLK1	●	●		
CLK2	●	●		
CLK3	●	●		
COT(MAP3K8)	●		●	
CRIK(CIT)	●	●		
DAPK1	●	●		
DCAMK2	●	●		
DLK(MAP3K12)	●		●	
DYRK1A	●	●		
DYRK1B	●	●		
DYRK2	●	●		
DYRK3	●	●		
DYRK4	●	●		
EEF2K	●	●		
Erk1(MAPK3)	●	●		
Erk2(MAPK1)	●	●		
Erk5(MAPK7)	●	●		
GSK3α(GSK3A)	●	●		
GSK3β(GSK3B)	●	●		
Hspnin(GSG2)	●	●		
HGK(MAP4K4)	●	●		
HIPK1	●	●		
HIPK2	●	●		
HIPK3	●	●		
HIPK4	●	●		
IKKa(CHUK)	●	●		
IKKβ(IKBKB)	●	●		
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	●	●	●	

Serine/Threonine Kinases	MSA	FP (IMAP™)	ELISA	TR-FRET
MELK	●	●		
MGC42105	●	●		
MINK(MINK1)	●	●		
MLK1(MAP3K9)	●	●	●	
MLK2(MAP3K10)	●	●	●	
MLK3(MAP3K11)	●	●	●	
MNK1(MKMK1)	●	●	●	
MNK2(MKMK2)	●	●	●	
MOS	●	●	●	
MRCCKα(CDC42BPA)	●	●	●	
MRCCKβ(CDC42BPB)	●	●	●	
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)	●	●	●	
PAK1	●	●	●	
PAK2	●	●	●	
PAK4	●	●	●	
PAK5	●	●	●	
PAK6	●	●	●	
PASK	●	●	●	
PBK	●	●	●	
PDHK2(PDK2)	●	●	●	
PDHK4(PDK4)	●	●	●	
PEK	●	●	●	
PGK(PRK1)	●	●	●	
PHKG1	●	●	●	
PHKG2	●	●	●	
PIM1	●	●	●	
PIM2	●	●	●	
PIM3	●	●	●	
PKAcα(PRKACA)	●	●	●	
PKAcβ(PRKACB)	●	●	●	
PKAcγ(PRKACG)	●	●	●	
PKCa(PRKCA)	●	●	●	
PKCB1(PRKCB1)	●	●	●	
PKCB2(PRKCB2)	●	●	●	
PKCγ(PRKCG)	●	●	●	
PKCδ(PRKCD)	●	●	●	
PKCε(PRKCE)	●	●	●	
PKCζ(PRKCZ)	●	●	●	
PKCη(PRKH)	●	●	●	
PKCθ(PRKCO)	●	●	●	
PKCι(PRKCI)	●	●	●	
PKD1(PRKD1)	●	●	●	
PKD2(PRKD2)	●	●	●	
PKD3(PRKD3)	●	●	●	
PKR	●	●	●	
PLK1	●	●	●	
PLK2	●	●	●	
PLK3	●	●	●	
PRKX	●	●	●	
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	●	●	●	
TGFBR1(TGFBFR1)	●	●	●	●
TNIK	●	●	●	
TSSK1	●	●	●	
TSSK2	●	●	●	
TSSK3	●	●	●	
TTK	●	●	●	
WEE1	●	●	●	
WNK1	●	●	●	
WNK2	●	●	●	
WNK3	●	●	●	

Lipid Kinase	MSA	FP (IMAP™)	ELISA	TR-FRET
PIK3CA/PIK3R1	●	●		
SPHK1	●	●		
SPHK2	●	●		

MSA (Mobility Shift Assay)
FP (IMAP™)
TR-FRET
ELISA

287
79
28
35

Updated: 2015/4/14