

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  
**327**  
Kinases\*

\* As of October 10, 2014

Platform	Minimum Size	Available Target	Kit Components
<p><b>Mobility Shift Assay</b> 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><b>284</b> Kinases</p>	<ul style="list-style-type: none"> <li>● Protein Kinase</li> <li>● Substrate Mixture (ATP, Metal included)</li> <li>● Assay Buffer</li> <li>● Termination Buffer</li> <li>● Assay Protocol (Separation conditions included)</li> </ul> <p>* Protocol sample is available online.</p> <p><b>After initial kit purchase, components may be purchased separately. Minimum order requirements apply.</b></p> <p>* The dilution ratio and MSA measuring parameters are different by kinase.</p>
<p><b>FP(IMAP™)</b> QSS Assist™ FP</p>	<p>800dp Equivalent to 2 x 384-well plate</p>	<p><b>79</b> Kinases</p>	<ul style="list-style-type: none"> <li>● Protein Kinase</li> <li>● Substrate Mixture</li> <li>● Assay Buffer</li> <li>● Assay Protocol</li> </ul>
<p><b>TR-FRET</b> QSS Assist™ TR-FRET</p>	<p>800dp Equivalent to 2 x 384-well plate</p>	<p><b>27</b> Kinases</p>	<ul style="list-style-type: none"> <li>● Protein Kinase</li> <li>● Substrate Mixture</li> <li>● Assay Buffer</li> <li>● Assay Protocol</li> </ul>
<p><b>ELISA</b> QSS Assist™ ELISA</p>	<p>500dp Equivalent to 5 x 96-well plate</p>	<p><b>35</b> Kinases</p>	<ul style="list-style-type: none"> <li>● Protein Kinase</li> <li>● Substrate Mixture</li> <li>● Assay Buffer</li> <li>● Antibody for ELISA (except for TTK &amp; WEE1)</li> <li>● Assay Protocol</li> </ul>

## 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 [L1196M]	●	
ALK [R1275Q]	●	
ALK [T1151_L1152insT]	●	
EML4-ALK	●	
NPM1-ALK	●	
ALK2(ACVR1)		●
ALK4(ACVR1B)		●
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/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(INSR)	●	
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]	●	
MUSK	●	
PDGFRa(PDGFR)	●	
PDGFRa(PDGFR) [D842V]	●	
PDGFRa(PDGFR) [T674I]	●	
PDGFRa(PDGFR) [V561D]	●	
PDGFRb(PDGFR)	●	●
PYK2(PTK2B)	●	
RET	●	
RET [G691S]	●	
RET [M918T]	●	
RET [S891A]	●	
RET [Y791F]	●	
RON(MST1R)	●	
ROS(ROS1)	●	
SRC	●	
SRM(SRMS)	●	
SYK	●	●
TEC	●	
TIE2(TEK)	●	●
TNK1	●	

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

Serine/Threonine Kinases	MSA	FP (IMAP™)	ELISA	TR-FRET
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]			●	●
BRK1	●			
BRK2	●			
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(PRKG2)	●	●		
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	●	●		
DCAMK12	●	●		
DLK(MAP3K12)	●		●	
DYRK1A	●			
DYRK1B	●			
DYRK2	●			
DYRK3	●			
DYRK4	●			
EEF2K	●			
Erk1(MAPK3)	●	●		
Erk2(MAPK1)	●	●		
Erk5(MAPK7)	●	●		
GSK3α(GSK3A)	●	●		
GSK3β(GSK3B)	●	●		
Haspin(GSG2)	●	●		
HGK(MAP4K4)	●	●		
HIPK1	●			
HIPK2	●			
HIPK3	●			
HIPK4	●			
IKKα(CHUK)	●	●		
IKKβ(IKKB)	●	●		
IKKε(IKBE)	●	●		
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(MKNK1)	●	●		
MNK2(MKNK2)	●	●		
MOS	●		●	
MRCRα(CDC42BPA)	●	●		
MRCRβ(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(PRKG1)	●	●		
PHKG1	●	●		
PHKG2	●	●		
PIM1	●	●		
PIM2	●	●		
PIM3	●	●		
PKAcα(PRKACA)	●	●		
PKAcβ(PRKACB)	●	●		
PKAcγ(PRKACG)	●	●		
PKCα(PRKCA)	●	●		
PKCβ1(PRKCB1)	●	●		
PKCβ2(PRKCB2)	●	●		
PKCγ(PRKCG)	●	●		
PKCδ(PRKCD)	●	●		
PKCε(PRKCE)	●	●		
PKCζ(PRKCZ)	●	●		
PKCη(PRKCH)	●	●		
PKCθ(PRKCQ)	●	●		
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(TGFB1)	●	●		●
TNIK	●	●		
TSSK1	●	●		
TSSK2	●	●		
TSSK3	●	●		
TTK	●			●
WEE1	●			●
WNK1	●	●		
WNK2	●	●		
WNK3	●	●		

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

MSA (Mobility Shift Assay) 284  
 FP (IMAP™) 79  
 TR-FRET 27  
 ELISA 35

Updated: 2014/10/10