

QuickScout Screening Assist[®] Kits from Carna Biosciences, Inc.

Tyrosine Kinases	MSA 800	MSA 400	TK-ELISA	TR-FRET
ABL(ABL1)	●	●	●	
ABL(ABL1) [E255K]	●	●	●	
ABL(ABL1) [T315I]	●	●	●	
ACK(TNK2)	●	●	●	
ALK	●	●	●	●
ALK [F1174L]	●	●	●	
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/L858R]	●	●	●	
EGFR [T790M]	●	●	●	
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	●	●	●	
PDGFRα(PDGFR)	●	●	●	
PDGFRα(PDGFR) [T674I]	●	●	●	
PDGFRα(PDGFR) [V561D]	●	●	●	
PDGFRβ(PDGFRB)	●	●	●	●
PYK2(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)	●	●	●	
TXK	●	●	●	
TYK2	●	●	●	●
TYRO3	●	●	●	
YES(YES1)	●	●	●	
ZAP70	●	●	●	

Serine/Threonine Kinases	MSA 800	MSA 400	FP (IMAP™)	STK-ELISA
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	●	●	●	●
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(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	●	●	●	
DCAMKL2	●	●	●	
DLK(MAP3K12)	●	●	●	●
DYRK1A	●	●	●	
DYRK1B	●	●	●	
DYRK2	●	●	●	
DYRK3	●	●	●	
EEF2K	●	●	●	
Erk1(MAPK3)	●	●	●	
Erk2(MAPK1)	●	●	●	
Erk5(MAPK7)	●	●	●	●
GSK3α(GSK3A)	●	●	●	
GSK3β(GSK3B)	●	●	●	●
Haspin(GSG2)	●	●	●	
HGK(MAP4K4)	●	●	●	
HIPK1	●	●	●	
HIPK2	●	●	●	
HIPK3	●	●	●	
HIPK4	●	●	●	
IKKα(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	●	●	●	
MAPKAP2	●	●	●	
MAPKAP3	●	●	●	
MAPKAP5	●	●	●	
MARK1	●	●	●	
MARK2	●	●	●	
MARK3	●	●	●	
MARK4	●	●	●	
MELK	●	●	●	
MGC42105	●	●	●	
MINK(MINK1)	●	●	●	
MLK1(MAP3K9)	●	●	●	●
MLK2(MAP3K10)	●	●	●	●
MLK3(MAP3K11)	●	●	●	●

Serine/Threonine Kinases	MSA 800	MSA 400	FP (IMAP™)	STK-ELISA
MNK1(MKNK1)	●	●	●	
MNK2(MKNK2)	●	●	●	
MOS	●	●	●	●
MRCα(CDC42BPA)	●	●	●	
MRCβ(CDC42BFB)	●	●	●	
MSK2(RPS6KA4)	●	●	●	
MSSK1(STK23)	●	●	●	
MST1(STK4)	●	●	●	
MST2(STK3)	●	●	●	
MST3(STK24)	●	●	●	
MST4	●	●	●	
NDR1(STK38)	●	●	●	
NDR2(STK38L)	●	●	●	
NEK1	●	●	●	
NEK2	●	●	●	
NEK4	●	●	●	
NEK6	●	●	●	
NEK7	●	●	●	
NEK9	●	●	●	
NPM1-ALK	●	●	●	
NuaK1	●	●	●	
NuaK2	●	●	●	
p38α(MAPK14)	●	●	●	
p38β(MAPK11)	●	●	●	
p38γ(MAPK12)	●	●	●	
p38δ(MAPK13)	●	●	●	
p70S6K(RPS6KB1)	●	●	●	
p70S6Kβ(RPS6KB2)	●	●	●	
PAK1	●	●	●	
PAK2	●	●	●	
PAK3	●	●	●	
PAK4	●	●	●	
PAK5	●	●	●	
PAK6	●	●	●	
PASK	●	●	●	
PBK	●	●	●	
PDHK2(PDK2)	●	●	●	
PDHK4(PDK4)	●	●	●	
PEK	●	●	●	
PGK(PRKG1)	●	●	●	
PHK1	●	●	●	
PHK2	●	●	●	
PIM1	●	●	●	
PIM2	●	●	●	
PIM3	●	●	●	
PKAα(PRKACA)	●	●	●	
PKAβ(PRKACB)	●	●	●	
PKAγ(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)	●	●	●	
PKN1	●	●	●	
PKR	●	●	●	
PLK1	●	●	●	
PLK2	●	●	●	
PLK3	●	●	●	
PLK4	●	●	●	●
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	●	●	●	
TNIK	●	●	●	
TSSK1	●	●	●	
TSSK2	●	●	●	
TSSK3	●	●	●	
TTK(TK-ELISA)	●	●	●	●
WEE1(TK-ELISA)	●	●	●	●
WNK1	●	●	●	
WNK2	●	●	●	
WNK3	●	●	●	

Lipid Kinase	MSA 800	MSA 400	FP (IMAP™)	STK-ELISA
PIK3CA/PIK3R1	●	●	●	
SPHK1	●	●	●	

MSA (Mobility Shift Assay) - 800dp 268
MSA (Mobility Shift Assay) - 400dp 205
FP (IMAP™) 78
ELISA 110
TR-FRET 22