

QuickScout Screening Assist® Kits

カルナバイオサイエンスの QuickScout Screening Assist® Kits はインハウスでキナーゼ阻害剤の特異性を調べるスクリーニング、プロファイリングにご利用いただけます。ATP濃度があらかじめKm値付近にされており、短期間で測定が可能です。本製品はカスタムメイドで、通常、ご発注後2~3週間でお届けいたします。



キナーゼの活性測定が簡単にできるアッセイキットには
このような特長や利点があります

- 当社プロファイリングサービスのノウハウを盛り込んだ信頼性の高い製品
- 必要な試薬とプロトコルがセットになった Ready-To-Run タイプ
- 1キットを複数回に分けてアッセイ実施が可能

対象キナーゼ
331
種類*

リピート購入率の高い **簡便さを実感** いただける製品です

*2015年 4月 14日 現在

プラットフォーム	最少販売単位	対象キナーゼ数	キット構成
<p>Mobility Shift Assay QSS Assist™ MSA</p> <p>パーキンエルマー社のLabChip® テクノロジーを用いた機器でご利用いただけます</p>	<p>400dp 1 x 384-well plate 相当</p>	<p>287 キナーゼ</p>	<ul style="list-style-type: none"> ● キナーゼ蛋白質 ● 基質ミクスチャー (ATP, Metal 含む) ● アッセイバッファー ● ターミネーションバッファー ● プロトコル (分離条件込み) HPでサンプル*公開 <p>MSAキット400dpを初回ご購入後は必要な試薬を単品でお求めいただけます。ATP, Metalを含まない基質単体も用意しておりますので、お客様で最適化を行ないアッセイいただくことも可能です。</p> <p>※ キナーゼの希釈倍率とMSAの測定パラメータはキナーゼ毎に異なります。</p>
<p>FP (IMAP™) QSS Assist™ FP</p>	<p>800dp 2 x 384-well plate 相当</p>	<p>79 キナーゼ</p>	<ul style="list-style-type: none"> ● キナーゼ蛋白質 ● 基質ミクスチャー ● アッセイバッファー ● プロトコル
<p>TR-FRET QSS Assist™ TR-FRET</p>	<p>800dp 2 x 384-well plate 相当</p>	<p>28 キナーゼ</p>	<ul style="list-style-type: none"> ● キナーゼ蛋白質 ● 基質ミクスチャー ● アッセイバッファー ● プロトコル
<p>ELISA QSS Assist™ ELISA</p>	<p>500dp 5 x 96-well plate 相当</p>	<p>35 キナーゼ</p>	<ul style="list-style-type: none"> ● キナーゼ蛋白質 ● 基質ミクスチャー ● アッセイバッファー ● ELISA用抗体 (TTK, WEE1は除く) ● プロトコル

[製造元]

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

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(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]	●	
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	●	●		
AMPKa1/B1/g1(PRKAA1/B1/G1)	●	●		
AMPKa2/B1/g1(PRKAA2/B1/G1)	●	●		
AurA(AURKA)	●	●		
AurA(AURKA)/TPX2	●	●		
AurB(AURKB)/INCENP	●	●		
AurC(AURKC)	●	●		
BMPR1A			●	
BMPR1B				●
BRAF			●	
BRAF [V600E]			●	
BRSK1	●			
BRSK2	●	●		
CaMK1a(CAMK1)	●	●		
CaMK1d(CAMK1D)	●	●		
CaMK2a(CAMK2A)	●	●		
CaMK2b(CAMK2B)	●	●		
CaMK2d(CAMK2D)	●	●		
CaMK2y(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)	●	●		
CK1a(CSNK1A1)	●	●		
CK1d(CSNK1D)	●	●		
CK1e(CSNK1E)	●	●		
CK1y1(CSNK1G1)	●	●		
CK1y2(CSNK1G2)	●	●		
CK1y3(CSNK1G3)	●	●		
CK2a1/B(CSNK2A1/B)	●	●		
CK2a2/B(CSNK2A2/B)	●	●		
CLK1	●	●		
CLK2	●	●		
CLK3	●	●		
COT(MAP3K8)			●	
CR1(CIT)	●	●		
DAPK1	●	●		
DCAMKL2	●	●		
DLK(MAP3K12)	●	●		
DYRK1A	●	●		
DYRK1B	●	●		
DYRK2	●	●		
DYRK3	●	●		
DYRK4	●	●		
EEF2K	●	●		
Erk1(MAPK3)	●	●		
Erk2(MAPK1)	●	●		
Erk5(MAPK7)	●	●		
GSK3a(GSK3A)	●	●		
GSK3b(GSK3B)	●	●		
Haspin(GSG2)	●	●		
HGK(MAP4K4)	●	●		
HIPK1	●	●		
HIPK2	●	●		
HIPK3	●	●		
HIPK4	●	●		
IKKa(CHUK)	●	●		
IKKb(IKKB)	●	●		
IKKe(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			●	
MRCCKa(CDC42BPA)	●	●		
MRCCKb(CDC42BPB)	●	●		
MSK2(RPS6KA4)	●	●		
MSSK1(STK23)	●	●		
MST1(STK4)	●	●		
MST2(STK3)	●	●		
MST3(STK24)	●	●		
MST4	●	●		
NDR1(STK38)	●	●		
NDR2(STK38L)	●	●		
NEK1	●	●		
NEK2	●	●		
NEK4	●	●		
NEK6	●	●		
NEK7	●	●		
NEK9	●	●		
NuaK1	●	●		
NuaK2	●	●		
p38a(MAPK14)	●	●		
p38b(MAPK11)	●	●		
p38y(MAPK12)	●	●		
p38d(MAPK13)	●	●		
p70S6K(RPS6KB1)	●	●		
PAK1	●	●		
PAK2	●	●		
PAK4	●	●		
PAK5	●	●		
PAK6	●	●		
PASK	●	●		
PBK	●	●		
PDHK2(PDK2)	●	●		
PDHK4(PDK4)	●	●		
PEK	●	●		
PGK(PRKG1)	●	●		
PHKG1	●	●		
PHKG2	●	●		
PIM1	●	●		
PIM2	●	●		
PIM3	●	●		
PKAcg(PRKACA)	●	●		
PKAcB(PRKACB)	●	●		
PKAcY(PRKACG)	●	●		
PKCa(PRKCA)	●	●		
PKCB1(PRKCB1)	●	●		
PKCB2(PRKCB2)	●	●		
PKCy(PRKCG)	●	●		
PKCδ(PRKCD)	●	●		
PKCe(PRKCE)	●	●		
PKCζ(PRKCF)	●	●		
PKCη(PRKCH)	●	●		
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) 287
 FP (IMAP™) 79
 TR-FRET 28
 ELISA 35

Updated: 2015/4/14