

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便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

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

* 2014年 10月 10日 現在

対象キナーゼ

327

種類*

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
Mobility Shift Assay QSS Assist™ MSA 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	400dp 1 x 384-well plate 相当	284 キナーゼ	● 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 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。 ※ 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。
FP(IMP™) QSS Assist™ FP	800dp 2 x 384-well plate 相当	79 キナーゼ	● 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
TR-FRET QSS Assist™ TR-FRET	800dp 2 x 384-well plate 相当	27 キナーゼ	● 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
ELISA QSS Assist™ ELISA	500dp 5 x 96-well plate 相当	35 キナーゼ	● 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 [L1196M]	●	
ALK [R1275Q]	●	
ALK [T1151_L1152insT]	●	
EML4-ALK	●	
NPM1-ALK	●	
ALK2(ACVR1)		●
ALK4(ACVR1B)		●
ARG(ABL2)	●	●
AXL	●	●
BLK	●	
BMX	●	
BRK(PTK6)	●	
BTX	●	●
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(INSRR)	●	
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(PDGFRB)	●	●
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	●	●		
AMPKa1/B1/y1(PRKAA1/B1/G1)	●			
AMPKa2/B1/y1(PRKAA2/B1/G1)	●			
AurA(AURKA)	●	●		
AurA(AURKA)/TPX2	●			
AurB(AURKB)/TNCENP	●	●		
AurC(AURKC)	●	●		
BMPR1A			●	
BMPR1B				●
BRAF			●	
BRAF [V600E]			●	
BRSK1	●			
BRSK2	●	●		
CaMK1a(CAMK1)	●			
CaMK1b(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)	●			
CK1b(CSNK1D)	●	●		
CK1e(CSNK1E)	●			
CK1y1(CSNK1G1)	●			
CK1y2(CSNK1G2)	●			
CK1y3(CSNK1G3)	●			
CK2a1/β(CSNK2A1/B)	●			
CK2a2/β(CSNK2A2/B)	●			
CLK1	●	●		
CLK2	●			
CLK3	●			
COT(MAP3K8)			●	
CRK(CIT)	●	●		
DAPK1	●	●		
DCAMK2	●			
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)		●		
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(MKNK1)	●	●		
MNK2(MKNK2)	●	●		
MOS			●	
MRCKa(CDC42BPA)	●	●		
MRCKb(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)	●	●		
p38b(MAPK13)	●	●		
p70S6K(RPS6KB1)	●	●		
PAK1	●			
PAK2	●			
PAK4	●			
PAK5	●			
PAK6	●			
PASK	●	●		
PBK	●	●		
PDHK2(PDK2)	●			
PDHK4(PDK4)	●			
PEK		●		
PGK(PRKG1)	●	●		
PHKG1	●	●		
PHKG2	●			
PIM1	●	●		
PIM2	●	●		
PIM3	●			
PKACa(PRKACA)	●	●		
PKACb(PRKACB)	●			
PKACy(PRKACG)	●			
PKCa(PRKCA)	●	●		
PKCb1(PRKCB1)	●	●		
PKCb2(PRKCB2)	●			
PKCy(PRKCG)	●			
PKCδ(PRKCD)	●	●		
PKCe(PRKCE)	●	●		
PKCζ(PRKCZ)	●	●		
PKCη(PRKCH)	●	●		
PKCθ(PRKCQ)	●	●		
PKCi(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	●			
TGFBF1(TGFBF1)				●
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