

QuickScout Screening Assist® Kits

カルナバイオサイエンスの **QuickScout Screening Assist® Kits** は希釈用の精製水を準備するだけで、受領後すぐに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る便利な製品です。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、通常で発注後2〜3週間でお届けが可能です。

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
- 1キットで最大4種 (1, 2, または4種) のキナーゼ選択が可能 (MSA 400dpのみ)



時間、手間、
コスト減に貢献!

4つの系で全312キナーゼ種に対応 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!

Mobility Shift Assay (MSA)

QSS Assist™ MSA

キャリパーライフサイエンス社のLabChip®テクノロジーを用いた機器でご利用可能なアッセイキットです



800dpセット

(384 well x 2 plate 相当)

272

キナーゼ

対象キナーゼから
1種をお選び下さい

400dpセット

(384 well x 1 plate 相当)

209

キナーゼ

対象キナーゼから
最大4種 (1, 2, または4種) を
お選び下さい

QSS Assist™ MSA キット構成

- 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でサンプルを公開しています

初回購入後は必要な試薬を単品で
追加で発注いただけます (MSAのみ)

詳しくはHPでご確認ください

FP (IMAP™)

QSS Assist™ FP

77

キナーゼ

キット構成

- 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

800dp/セット (384 well x 2 plate 相当)

TR-FRET

QSS Assist™ TR-FRET

22

キナーゼ

キット構成

- 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

800dp/セット (384 well x 2 plate 相当)

ELISA

QSS Assist™ ELISA

110

キナーゼ

キット構成

- 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

500dp/セット (96 well x 5 plate 相当)

2012年9月25日更新

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 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	●	●	●	
FGFR1 [V561M]	●	●	●	
FGFR2	●	●	●	
FGFR2 [N549H]	●	●	●	
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	●	●	●	●
JAK2(JH1 JH2)	●	●	●	
JAK2(JH1 JH2) [V617F]	●	●	●	
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)	●	●	●	
YES(YES1) [T348I]	●	●	●	
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)	●	●	●	
BMPRI1A				●
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)				●
CRK(CIT)	●			
DAPK1	●	●	●	
DCAMK12	●			
DLK(MAP3K12)				●
DYRK1A	●	●		
DYRK1B	●	●		
DYRK2	●	●		
DYRK3	●	●		
EEF2K	●			
Erf1(MAPK3)	●	●	●	
Erf2(MAPK1)	●	●	●	
Erf5(MAPK7)	●			●
GSK3α(GSK3A)	●	●		
GSK3β(GSK3B)	●	●		●
Haspin(GSG2)	●	●		
HGK(MAP4K4)	●	●	●	
HIPK1	●	●		
HIPK2	●	●		
HIPK3	●	●		
HIPK4	●	●		
IKKα(CHUK)	●		●	
IKKβ(IKKBK)	●	●		
IKKε(IKBE)	●		●	
IRAK1			●	
IRAK4	●		●	
JNK1(MAPK8)				●
JNK2(MAPK9)				●
JNK3(MAPK10)				●
LATS2	●	●		
LIHK1				●
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	●	●		
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				●
MRCkα(CDC42BPA)	●		●	
MRCkβ(CDC42BPB)	●			
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)	●	●	●	
PAK1	●	●		
PAK2	●	●		
PAK3	●	●		
PAK4	●	●		
PAK5	●	●		
PAK6	●	●		
PASK	●		●	
PBK	●	●	●	
PDHK2(PDK2)	●			
PDHK4(PDK4)	●			
PEK			●	
PGK(PRKG1)	●	●	●	
PHKG1	●	●	●	
PHKG2	●			
PIM1	●	●	●	
PIM2	●	●	●	
PIM3	●	●		
PKAα(PRKACA)	●		●	
PKAβ(PRKACB)	●	●		
PKAγ(PRKACG)	●			
PKCa(PRKCA)	●	●	●	
PKCB1(PRKCB1)	●	●	●	
PKCB2(PRKCB2)	●	●		
PKCγ(PRKCG)	●	●		
PKCδ(PRKCD)	●	●	●	
PKCe(PRKCE)	●	●	●	
PKCζ(PRKCZ)	●	●	●	
PKCη(PRKCH)	●	●	●	
PKCθ(PRKCQ)	●	●	●	
PKCι(PRKCI)	●	●	●	
PKD1(PRKD1)	●	●	●	
PKD2(PRKD2)	●	●	●	
PKD3(PRKD3)	●	●	●	
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 272
MSA (Mobility Shift Assay) - 400dp 209
FP (IMAP™) 77
ELISA 110
TR-FRET 22