

Kinase Profiling Services

対応キナーゼ

311
種類

カルナバイオサイエンスの プロファイリングサービス

このような特長や利点があります

裏面リストを
ご覧下さい

■ 世界トップクラスのキナーゼ数

活性を指標にしたアッセイとして世界で最も多種のキナーゼでプロファイリングが可能です。
311種類からお客様ご自身で任意に、またはフルパネルを選択いただくことで、ご予算に応じて必要、十分なデータをご提供します。

■ 使用するキナーゼ蛋白質はすべて自社にて製造・品質管理

キナーゼの性質を十分把握できており、安定、且つご安心いただける材料を使用しています。同材料を5ugの少量包装からバルクまでご購入いただけます。

■ 独自のPre-セレクトパネル

網羅的なデータ取得には弊社が独自にキナーゼ種を選択したQuickScout®シリーズパネルがおすすめです。

■ ご研究のスピードアップに貢献

阻害率、IC₅₀測定共に結果を1~2週間*で報告します。
*MSA/IMAP™で測定するキナーゼ種が対象です。ELISAは最長で4週間かかりますこと予めご了承下さい。

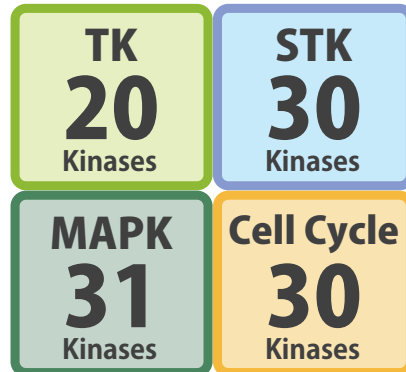
■ 高濃度設定(ATP=1mM)も可能

既存のKm値付近に加え、細胞内のATP濃度を想定した新しいサービスで、キナーゼに対する化合物の阻害作用、特に選択性を詳細に知る上で有効な手法の一つです。

■ メインのアッセイ系はMobility Shift Assay (MSA)

基質のリン酸化がダイレクトに見える Caliper社のLabChip®3000 を用いたアッセイプラットフォームで信頼性の高いデータをご提供します。

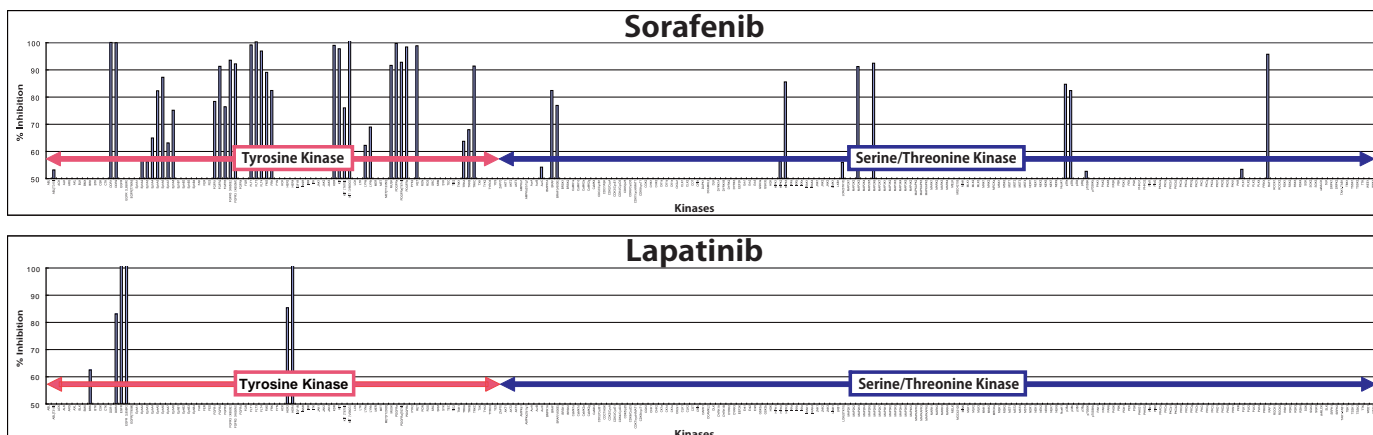
キナーゼを当社独自にセレクト
QuickScout®
Pre-Select パネルシリーズ



各パネルの詳細は別刷りのチラシ、
またはHPでご覧下さい

キナーゼ阻害剤のプロファイリング結果グラフ

自社製の256キナーゼ蛋白質に対する阻害率を濃度1uMで測定したものです(ATP=Km値付近)



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

QuickScout® Custom Profiling & Panel Profiling Series from Carna Biosciences, Inc.

Tyrosine Kinases	Custom		Panel
	Km bin.	1mM	TK
ABL(ABL1)	●	●	●
ABL(ABL1) [E255K]	●	●	●
ABL(ABL1) [T315I]	●	●	●
ACK(TNK2)	●	●	●
ALK	●	●	●
ALK [F1174L]	●	●	●
ALK [L1196M]	●	●	●
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]	●	●	●
EGFR [T790M/L858R]	●	●	●
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	●	●	●
NPM1-ALK	●	●	●
PDGFRα(PDGFR)	●	●	●
PDGFRα(PDGFR) [T674I]	●	●	●
PDGFRα(PDGFR) [V651D]	●	●	●
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	Custom		Panel		
	Km bin.	1mM	STK	MAPK	Cell
AKT1	●	●	●		
AKT2	●	●	●		
AKT3	●	●	●		
AMPKα1/β1/v1(PRKAA1/B1/G1)	●	●	●		
AMPKα2/β1/v1(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)	●	●	●		
CaMK2γ(CAMK2G)	●	●	●		
CaMK2δ(CAMK2D)	●	●	●		
CaMK4	●	●	●		
CDC2/CycB1	●	●	●		●
CDC7/ASK	●	●	●		●
CDK2/CycA2	●	●	●		●
CDK2/CycE1	●	●	●		●
CDK3/CycE1	●	●	●		●
CDK4/CycD3	●	●	●		●
CDK5/p25	●	●	●		●
CDK6/CycD3	●	●	●		●
CDK7/CycH/MAT1	●	●	●		●
CDK9/CycT1	●	●	●		●
CGK2(PRKG2)	●	●	●		●
CHK1(CHEK1)	●	●	●		●
CHK2(CHEK2)	●	●	●		●
CK1α(CSNK1A1)	●	●	●		●
CK1γ1(CSNK1G1)	●	●	●		●
CK1γ2(CSNK1G2)	●	●	●		●
CK1γ3(CSNK1G3)	●	●	●		●
CK1δ(CSNK1D)	●	●	●		●
CK1ε(CSNK1E)	●	●	●		●
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	●	●	●		
MAPKAPK2	●	●	●		
MAPKAPK3	●	●	●		
MAPKAPK5	●	●	●		
MARK1	●	●	●		
MARK2	●	●	●		
MARK3	●	●	●		
MARK4	●	●	●		
MELK	●	●	●		
MGC42105	●	●	●		
MINK(MINK1)	●	●	●		

Custom Profiling	Km bin.	311
	1mM	166
Panel Profiling	TK20	20
	STK30	30
	MAPK	31
	Cell Cycle	30

Serine/Threonine Kinases	Custom		Panel		
	Km bin.	1mM	STK	MAPK	Cell
MLK1(MAP3K9)	●	●		●	
MLK2(MAP3K10)	●	●		●	
MLK3(MAP3K11)	●	●		●	
MNK1(MKNK1)	●	●			
MNK2(MKNK2)	●	●			
MOS	●	●		●	
MRCRα(CDC42BPA)	●	●			
MRCRβ(CDC42BPB)	●	●			
MSK1(RPS6KA5)	●	●			
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)	●	●	●		●
p70S6KB(RPS6KB2)	●	●	●		●
PAK1	●	●	●		
PAK2	●	●	●		
PAK3	●	●	●		
PAK4	●	●	●		
PAK5(PAK7)	●	●	●		
PAK6	●	●	●		
PASK	●	●	●		
PBK	●	●	●		
PDHK2(PDK2)	●	●	●		
PDHK4(PDK4)	●	●	●		
PDK1(PDKP1)	●	●	●		
PEK(EIF2AK3)	●	●	●		
PGK(PRKG1)	●	●	●		
PHKG1	●	●	●		
PHKG2	●	●	●		
PIM1	●	●	●		●
PIM2	●	●	●		●
PIM3	●	●	●		●
PKAα(PRKACA)	●	●	●		
PKAβ(PRKACB)	●	●	●		
PKAγ(PRKACG)	●	●	●		
PKCα(PRKCA)	●	●	●		
PKCβ1(PRKCB1)	●	●	●		
PKCβ2(PRKCB2)	●	●	●		
PKCγ(PRKCG)	●	●	●		
PKCδ(PRKCD)	●	●	●		
PKCε(PRKCE)	●	●	●		
PKCζ(PRKCZ)	●	●	●		
PKCη(PRKCJ)	●	●	●		
PKCθ(PRCKO)	●	●	●		
PKCι(PRCKI)	●	●	●		
PKD1(PRKD1)	●	●	●		
PKD2(PRKD2)	●	●	●		
PKD3(PRKD3)	●	●	●		
PKN1	●	●	●		
PKR(EIF2AK2)	●	●	●		
PLK1	●	●	●		●
PLK2	●	●	●		●
PLK3	●	●	●		●
PLK4	●	●	●		●
PRXX	●	●	●		
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	●</				