

KINEXUSTM

PHOSPHO-SITE PROFILING SERVICES

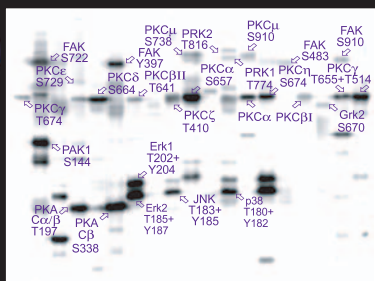
KPSS 10.0



Human A431 Cells

B23 (NPM) S4
B23 (NPM) T199
B23 (NPM) T234
BRCA1 S1497
CDK1/CDK2 T14+Y15
CDK1/CDK2 Y15
CDK1/CDK2 T161/T160
Erk1 T202+Y204
Erk2 T185+Y187
GSK3 α S21
GSK3 α Y279
GSK3 β S9
GSK3 β Y216
MEK1 S217+S221
MEK1 T291
MEK1 S297
MEK1 T385
MEK2 T394
mTOR S2448
p27 Kip1 T187
p53 S392
PDK1 S244
PKB (Akt) T308
PKB (Akt) S473
PTEN S380+S382+S385
Raf1 S259
Rb T356
Rb S612
Rb S780
Rb S807
Rb S811
Rb T821
Rb T826
Rsk1/2 S221/S227
Rsk1/2 S363/S369
Rsk1/2 S380/S386
S6 S235
S6K α T229
S6K α T421+S424
Src Y418
Src Y529

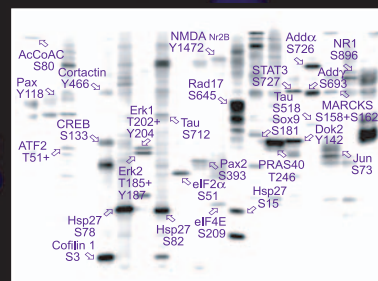
KPSS 11.0



Mouse NIH-3T3 Cells

EGFR Y1148
ErbB2 Y1248
Erk1 T202+Y204
Erk2 T185+Y187
Erk5 (BMK1) T218+Y220
FAK Y397
FAK S722
FAK S843
FAK S910
GRK2 S670
InsulinR Y999
IR/IGF1R T1189/Y1190
JNK T183+Y185
Kit Y703
Lyn Y507
p38 α T180+Y182
PAK1/2/3 S144/S141/S139
PKA C α / β T197
PKA C β S338
PKC α S657
PKC α / β 2 T638/T641
PKC β 1/ β 2 T500
PKC β 2 T641
PKC δ Y313
PKC δ S664
PKC ϵ S729
PKC γ T514
PKC γ T655
PKC γ T674
PKC η S674
PKC μ (PKD) S738+S742
PKC μ (PKD) S910
PKC ζ / λ T410/T403/T403
PKR1 T451
PRK1 T774
PRK2 T816
Pyk2 Y579

KPSS 12.0



Human HeLa Cells

4E-BP1 S65
AcCoA Carb S80
Adducin α S726
Adducin γ S693
ATF2 T51+T53
BLNK Y84
Cofilin S3
Cortactin Y470
CREB1 S133
Dok2 Y139
eIF2 α S51
eIF4E S209
eIF4G S1107
Erk1 T202+Y204
Erk2 T185+Y187
FKHRL1 T32
Hsp27 S15
Hsp27 S78
Hsp27 S82
IRS1 Y612
Jun S73
MARCKS S158+S162
NMDA Nr2B Y1474
NR1 S896
Paxillin Y31
Paxillin Y118
Pax2 S394
PRAS40 T246
Rac1/cdc42 S71
Rad17 S645
Shc Y349+Y350
SOX9 S181
STAT1 Y701
STAT3 S727
Synapsin 1a S9
Tau S515
Tau S518
Tau S712

Kinexus uses highly, validated commercial antibodies in our immunoblotting services.

www.kinexus.ca

Toll free (US): 1 866 KINEXUS Ph: 1 604 822 9963 Fax: 1 604 822 9964 info@kinexus.ca