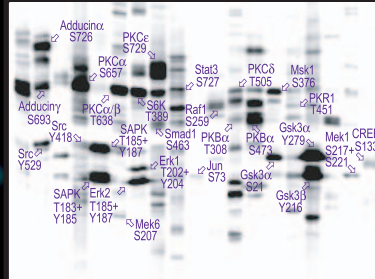




KINEXUS™

PHOSPHO-SITE PROFILING SERVICES

KPSS 1.3



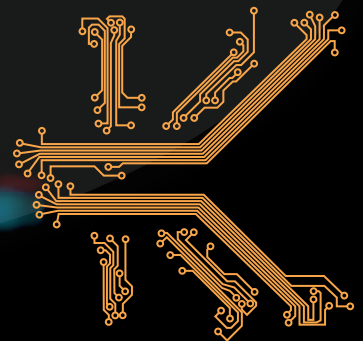
Rat Brain Cortex

Reversible phosphorylation is the major form of post-translational modification and regulation of proteins in eukaryotic cells. Kinexus has tested over 800 commercial phospho-site antibodies for incorporation of the best of these into our Kinetworks™ KPSS services. The KPSS screens utilize from 34 to 42 phospho-site antibodies, and work in human, mouse, rat and many other species due to the high conservation of phospho-sites. In addition to tracking over 200 known phospho-sites, hundreds of other phosphorylation sites can also be visualized by cross-reactivity. With antibody-based phosphoprotein discovery, it is often possible to quickly follow up with Western blotting, immunoprecipitation, immunohistochemistry and flow cytometry studies. Kinetworks™ is unmatched for breadth, accuracy, reproducibility, convenience, turnaround time and affordability.

- Adducin-α S726
- Adducin-γ S693
- B23 (NPM) S4
- CDK1/2 T14+Y15
- CREB1 S133
- Erk1 T202+Y204
- Erk2 T185+Y187
- GSK3α S21
- GSK3α Y279
- GSK3β S9
- GSK3β Y216
- JNK T183+Y185
- Jun S73
- MEK1 S217+S221
- MEK3 S189
- MEK6 S207
- Msk1 S376
- NR1 S896
- p38α MAPK T180+Y182
- PKBα (Akt1) T308
- PKBα (Akt1) S473
- PKCα S657
- PKCα/β2 T638/T641
- PKCδ T507
- PKCε S729
- PKR1 T451
- Raf1 S259
- Rb S780
- Rb S807+S811
- RSK1/2 T359/T365
- S6Kα T389
- Smad1/5/9 S463+S465
- Src Y418
- Src Y529
- STAT1 Y701
- STAT3 S727
- STAT5 Y694

Kinetworks™
is unmatched for:

- CONVENIENCE
- ACCURACY
- HIGH SENSITIVITY
- FAST TURN AROUND
- REPRODUCIBILITY
- AFFORDABILITY
- QUANTIFICATION
- BREADTH



www.kinexus.ca

KINEXUS™

PHOSPHO-SITE PROFILING SERVICES

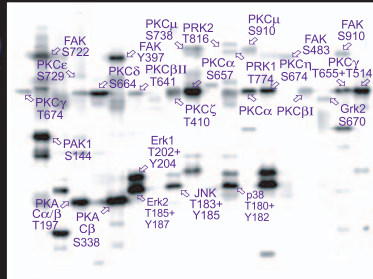
KPSS 10.0



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

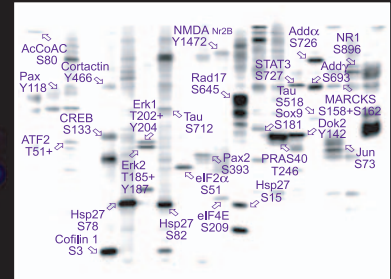
KPSS 11.0



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
 IIR/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 / η / λ T410/T403/T403
 PKR1 T451
 PRK1 T774
 PRK2 T816
 Pyk2 Y579

KPSS 12.0



Human HeLa Cells

4E-BP1 S65
 AcCoACarb 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