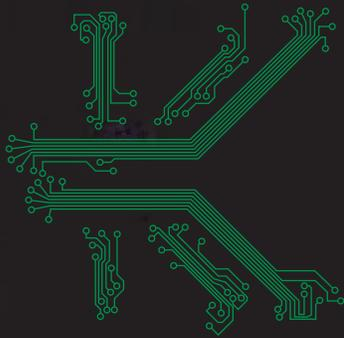


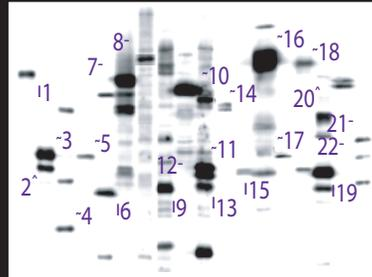


KINEXUS

CUSTOM IMMUNOBLOTTING PROFILING SERVICES

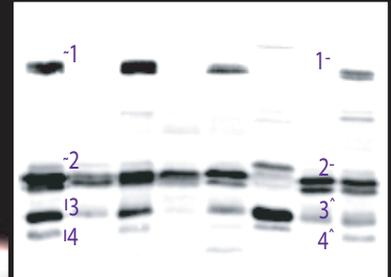


KCPS 1.0



Rat Brain Lysate Probed with 18 Protein Kinase Antibodies

KCSS 1.0



Monkey Tissue Lysates Probed with 3 Protein Kinase Antibodies

Whereas the Kinetworks™ Standard Signal Transduction Protein Profiling Services offer the highest value, the Kinetworks™ Custom Services provide the greatest flexibility. Clients can choose from combinations of at least 980 pan-specific and 1047 phosphosite-specific antibodies to have Kinexus test their tissue/cell lysates with two possible formats. The Kinetworks™ Custom Protein Screening Service (KCPS) permits probing of one sample with up to 18 different antibodies. The Kinetworks™ Custom Sample Screening Service (KCSS) allows probing of up to 8 samples with 3 different antibodies. These services can also accept antibodies that are provided by clients. The Custom Kinetworks™ screens are recommended to confirm results from initial standard Kinetworks™ or Kinex™ microarray analyses.

- PKCα 1
- ERK1 2
- ERK2 3
- CDK5 4
- MEK1 5
- MEK6 6
- PKCβ 7
- FAK 8
- Mos 9
- PKCζ 10
- CK2α 11
- CK2α' 12
- CK2α" 13
- PAKβ 14
- PKA 15
- PKCε 16
- MEK2 17
- RafB 18
- MEK4 19
- RSK1 20
- GSK3α 21
- GSK3β 22

- RSK1 1
- ERK1 2
- ERK2 3
- MEK6 4

- Brain - Lane 1
- Heart - Lane 2
- Kidney - Lane 3
- Liver - Lane 4
- Lung - Lane 5
- Sk. Muscle - Lane 6
- Spleen - Lane 7
- Testes - Lane 8

Kinetworks™ is unmatched for:

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

www.kinexus.ca