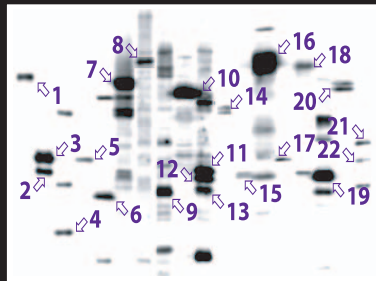
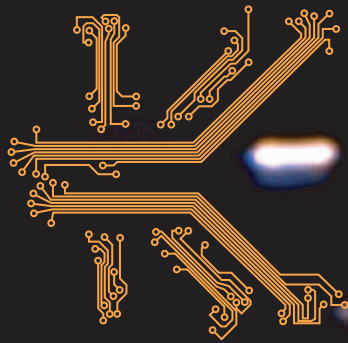


KINETWORKS™

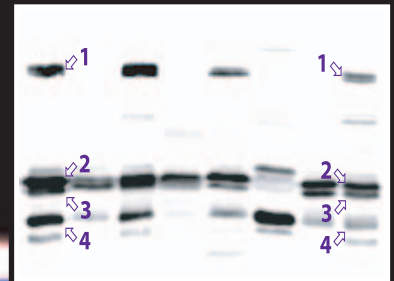
CUSTOM PROTEIN PROFILING SERVICES

KCPS 1.0

KCSS 1.0



Rat Brain Lysate Probed with 18 Protein Kinase Antibodies



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 over 300 pan-specific and 200 phospho-site-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.

PKC α	1	RSK1	1
ERK1	2	ERK1	2
ERK2	3	ERK2	3
CDK5	4	MEK6	4
MEK1	5		
MEK6	6	Brain - Lane 1	
PKC β	7	Heart - Lane 2	
FAK	8	Kidney - Lane 3	
Mos	9	Liver - Lane 4	
PKC ζ	10	Lung - Lane 5	
CK2 α	11	Sk. Muscle - Lane 6	
CK2 α'	12	Spleen - Lane 7	
CK2 α''	13	Testes - Lane 8	
PAK β	14		
PKA	15		
PKC ϵ	16		
MEK2	17		
RafB	18		
MEK4	19		
RSK1	20		
GSK3 α	21		
GSK3 β	22		

Kinetworks™ is unmatched for:

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

www.kinexus.ca