

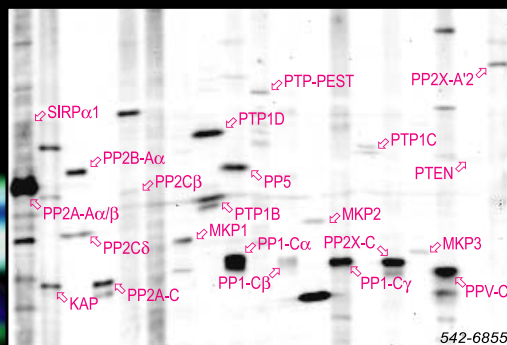
KINEXWORKS™

KPPS 1.2

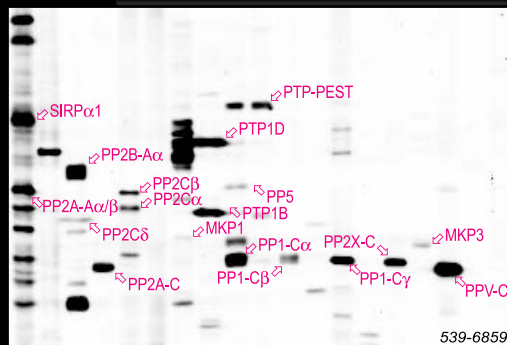
PROTEIN PHOSPHATASE PROFILING SERVICE

Using Highly Validated, Potent and
Specific Commercial Antibodies

CD45
Cdc25B
Cdc25C
KAP
LAR
MKP1
MKP2
PAC1
PP1/C α
PP1/C β
PP1/C γ
PP2A/A α /B'
PP2A/C α
PP2A/C β
PP2B/A α
PP2C α
PP2C β
PP2C δ
PP4/A'2
PP4C
PP5C
PP6C
PTEN
PTP1B
PTP1C
PTP1D
PTP-PEST
SIRP α 1
VHR



Human HC T116 Colon Carcinoma Cells



Rat Brain Cortex Tissue

Kinetworks™
is unmatched for:

- AFFORDABILITY
- ACCURACY
- SENSITIVITY
- QUANTITATION
- REPRODUCIBILITY
- TURN AROUND
- CONVENIENCE
- BREADTH

www.kinexus.ca

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