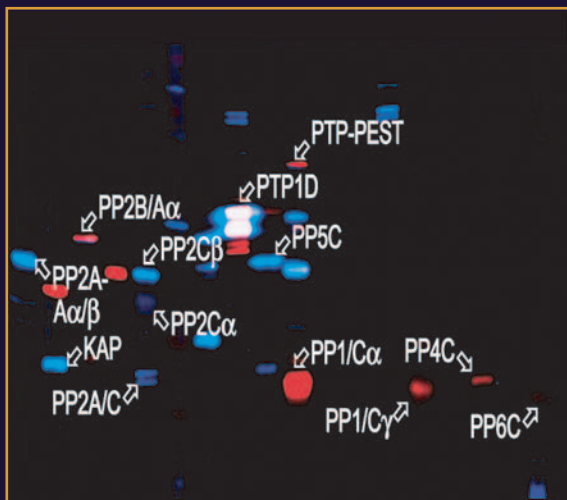
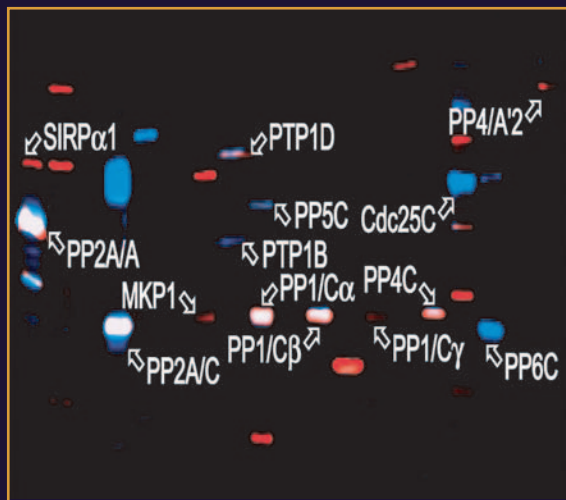


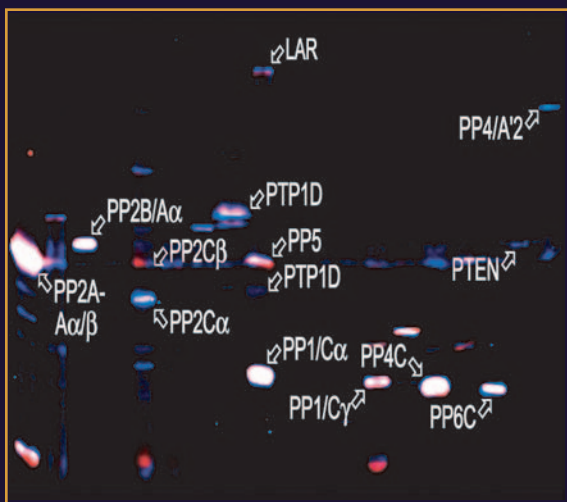
EXAMPLES OF KINETWORKS™ KPPS 1.2 PHOSPHATASE SCREENS



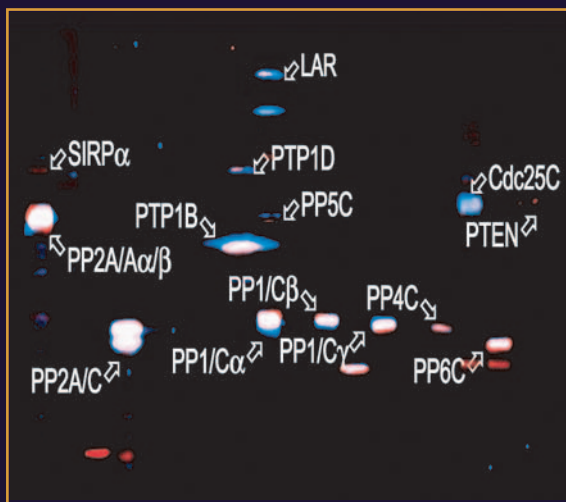
Human Spinal Cord (blue)
Mouse Spinal Cord (red)



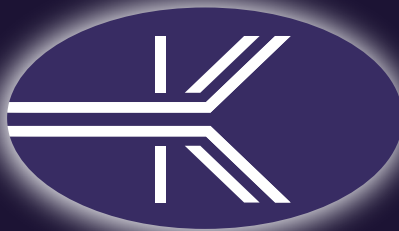
Human BG03 Embryo Stem Cells (blue)
Rat Corneal Endothelium Cells (red)



Mouse Skin Cells
+ Mechlorethamine



Human Embryonic Kidney HEK-293 Cells
+ Treatment



Effect of Treatment
on Phosphorylation

- Increased
- Unchanged
- Decreased