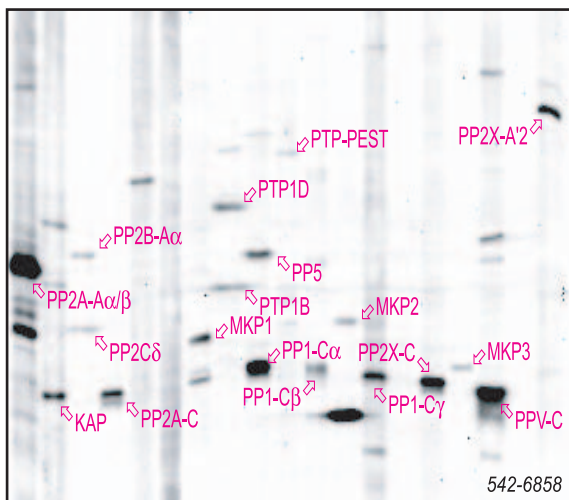


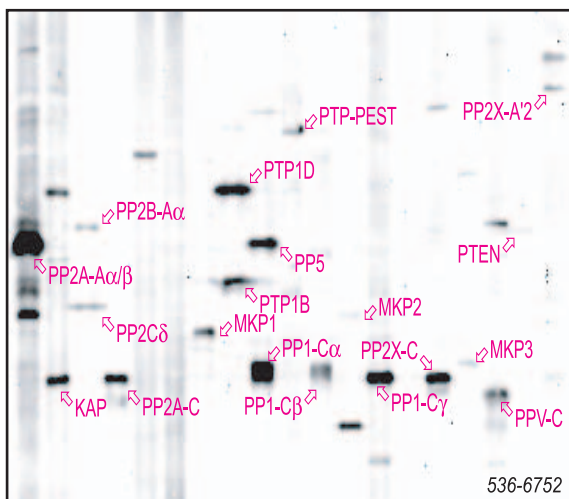
EXAMPLES OF KINETWORKS™ KPPS 1.1 PROTEIN PHOSPHATASE SCREENS



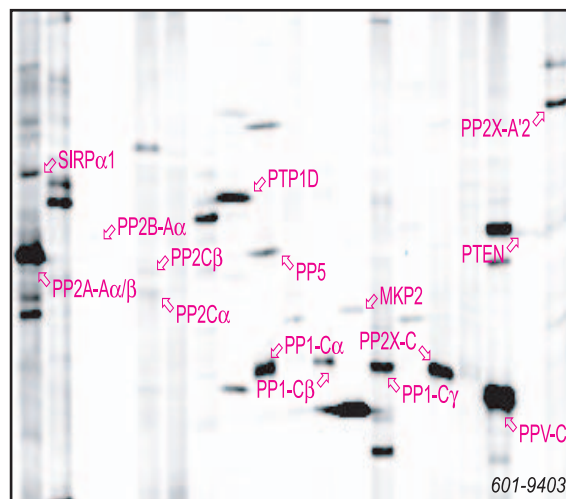
Human HeLa Cervical Carcinoma Cells



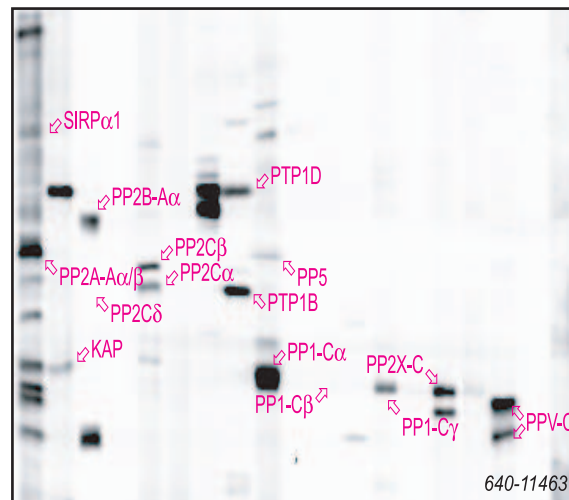
Human HCT116 Colon Tumour Cells



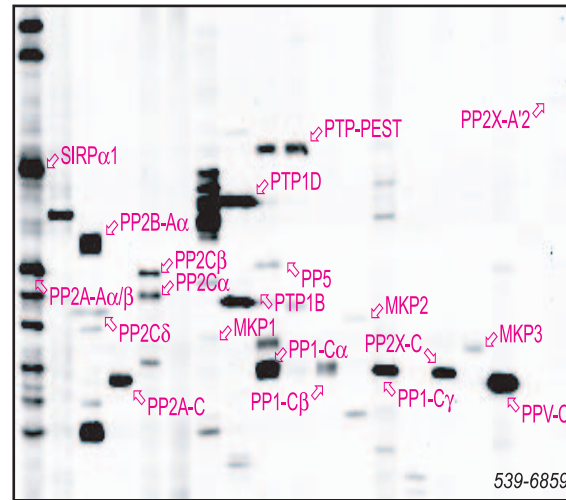
Human A375 Lung Tumour Cells



Chinese Hamster Ovary Cells

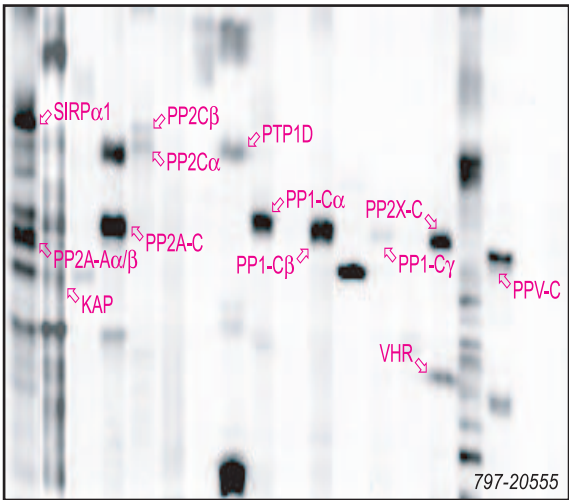


Mouse Whole Brain

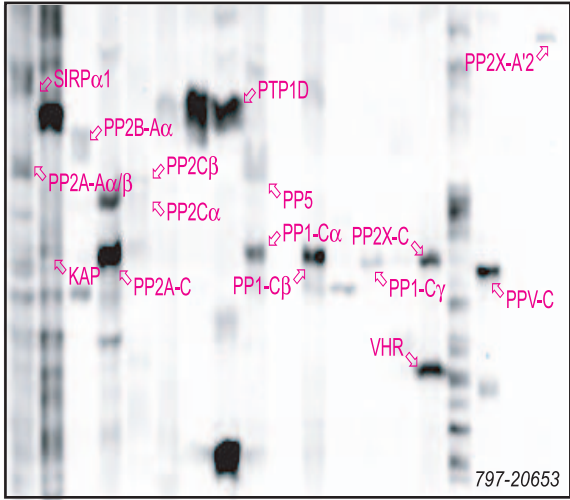


Rat Brain Cortex

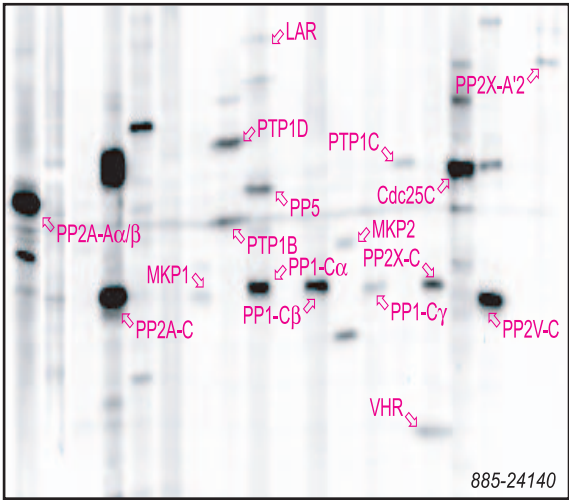
EXAMPLES OF KINETWORKS™ KPPS 1.1 PROTEIN PHOSPHATASE SCREENS



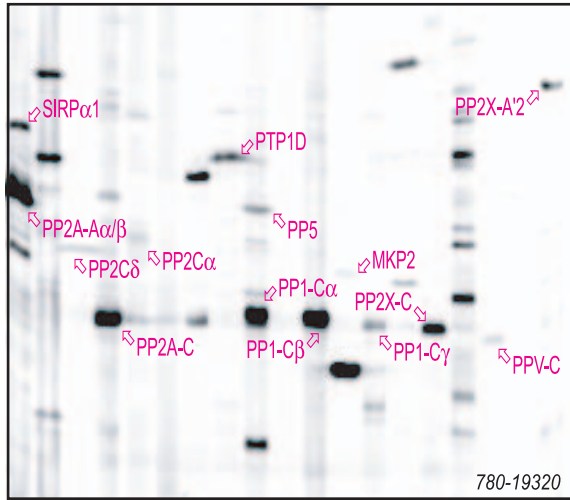
Human Glioblastoma Multiform Cells



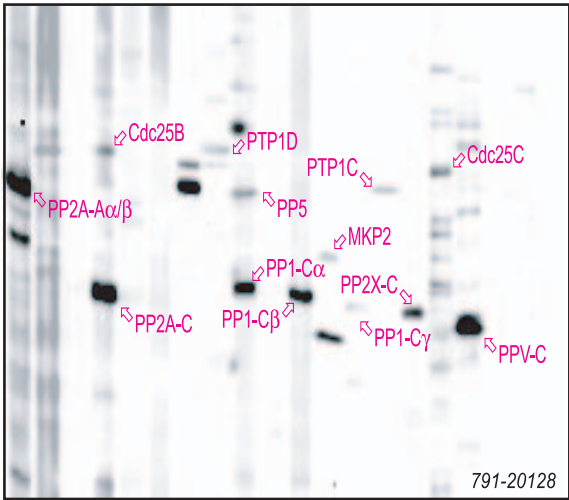
Rat Whole Brain



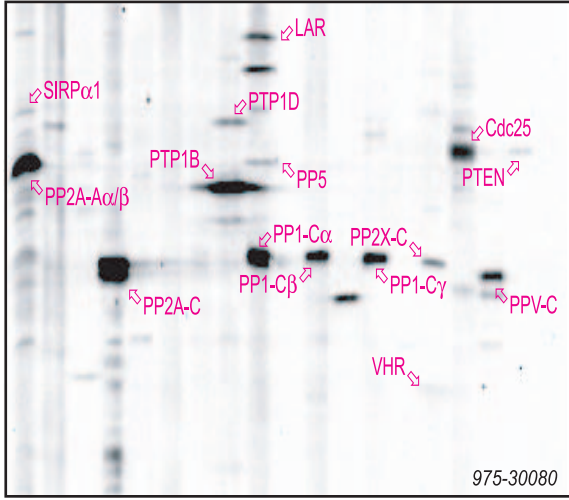
Human BG03 Embryo Stem Cells



Rat Corneal Endothelium Cells



Mouse Bone Marrow Cells



Human HEK 293 Embryonic Kidney Cells