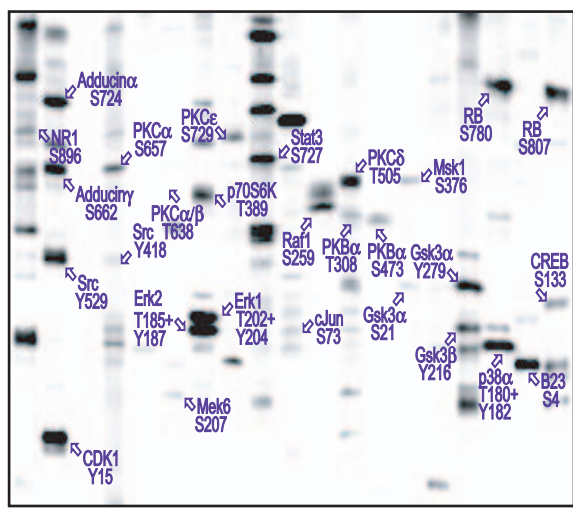


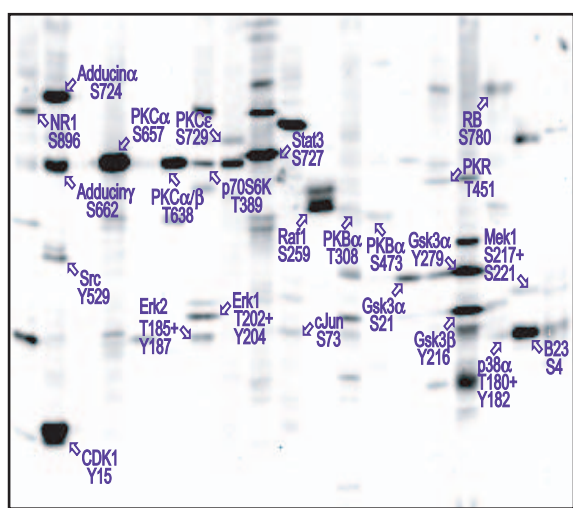
EXAMPLES OF KINETWORKS™ KPSS 1.3 PHOSPHO-SITE SCREENS



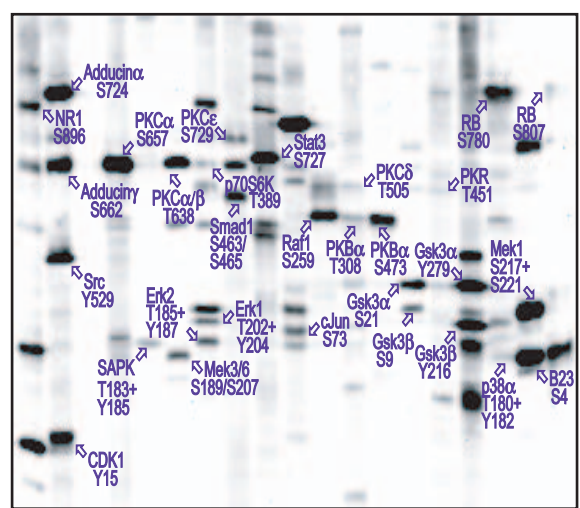
Human MCF7 Breast Tumour Cells



Human T47D Breast Tumour Cells



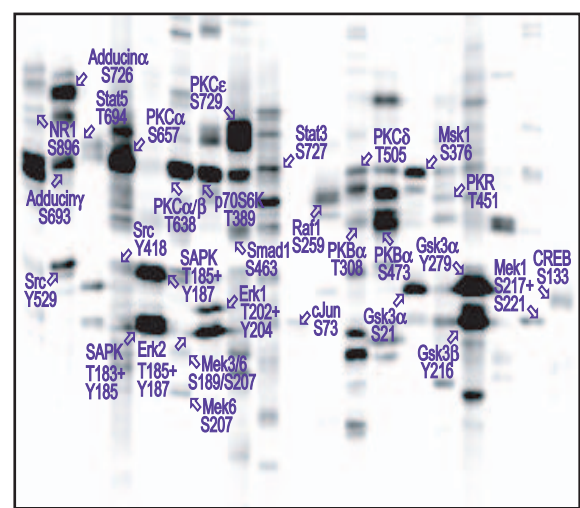
Human HeLa Cervical Carcinoma Cells



Human A375 Melanoma Cells



Human NCI-H69 Lung Carcinoma Cells



Rat Brain Cortex