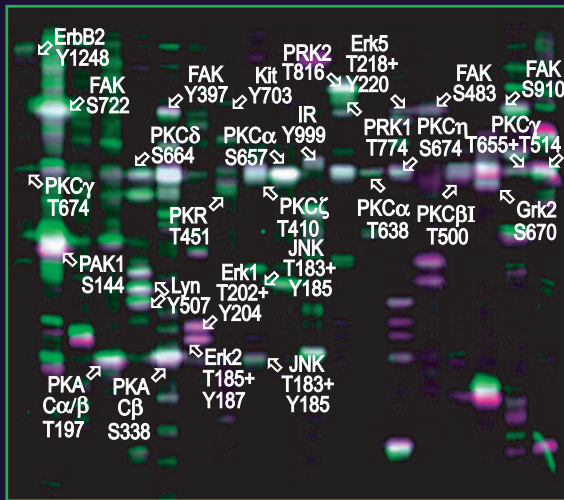


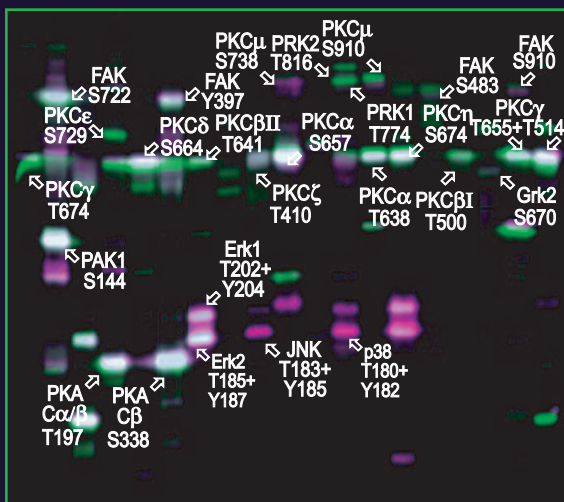
EXAMPLES OF KINETWORKS™ KPSS 11.0 PHOSPHO-SITE SCREENS



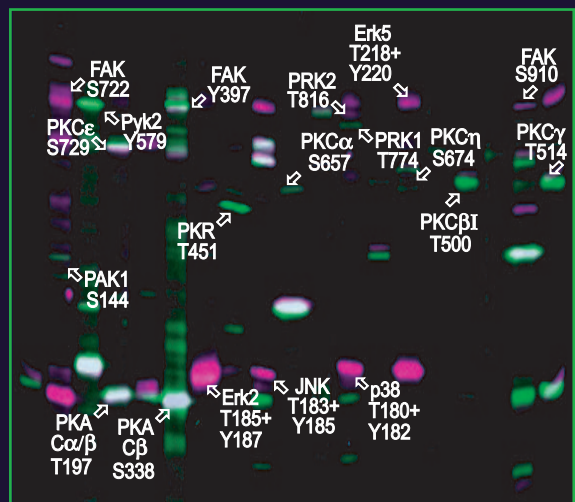
Human A431 Epidermoid Carcinoma Cells
+ 100 ng/ml Epidermal Growth Factor for 5 min



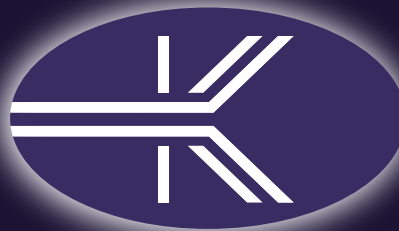
Human HeLa Cervical Carcinoma Cells
+ 20 ng/ml Tumour Necrosis Factor- α for 30 min



Mouse NIH3T3 Fibroblasts
+ 15 ng/ml Phorbol Myristate Acetate for 10 min



Xenopus laevis Maturing Oocytes
+ 1 μ M Progesterone for 6 h



Effect of Treatment on Phosphorylation

- Increased
- Unchanged
- Decreased