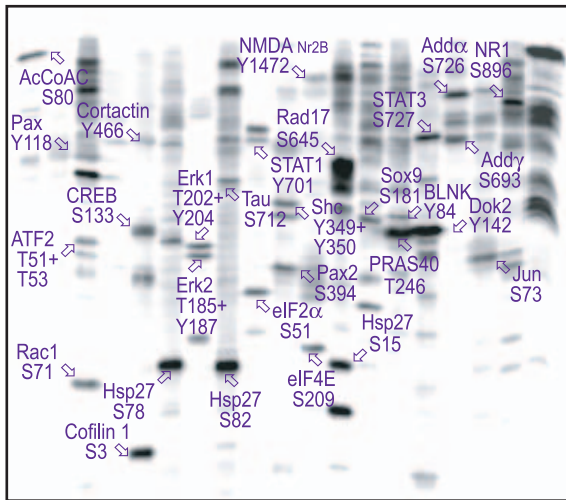


EXAMPLES OF KINETWORKS™

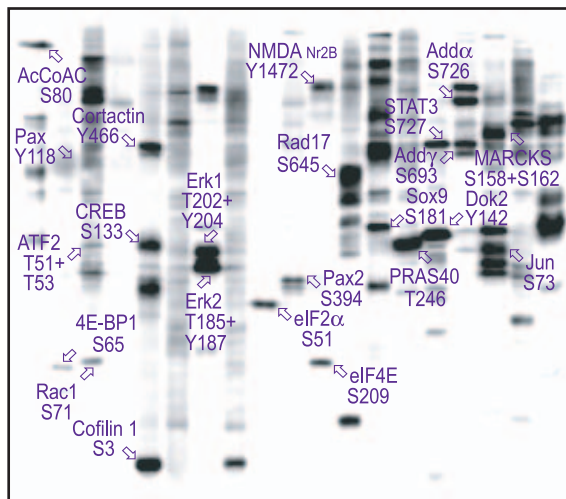
KPSS 12.0 PHOSPHO-SITE SCREENS



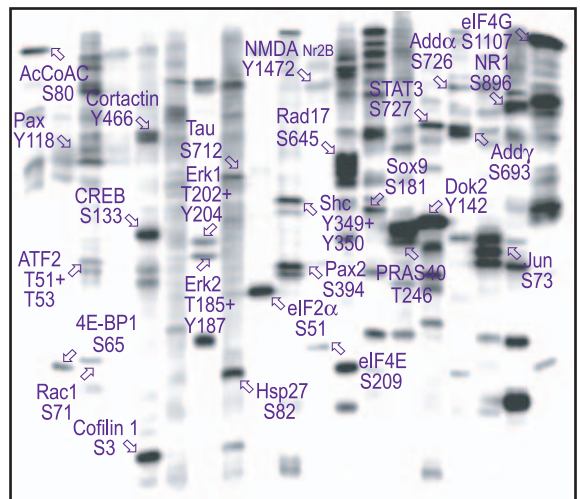
Human A431 Epidermoid Carcinoma Cells



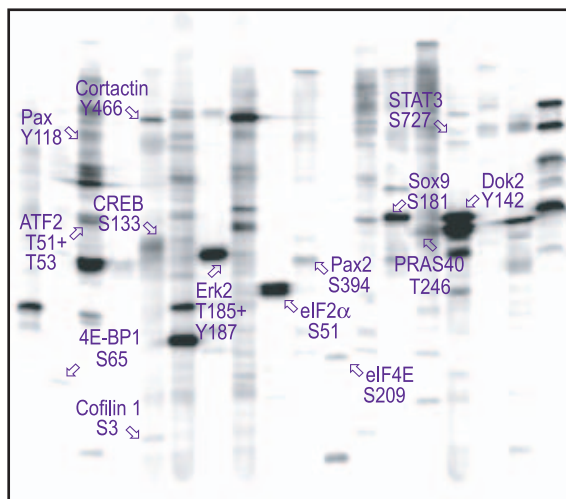
Human HeLa Cervical Carcinoma Cells



Mouse NIH-3T3 Fibroblasts



Mouse Pam212 Keratinocytes



Xenopus laevis Maturing Oocytes

S
 P
 X
 W
 Z
 T
 V