

EXAMPLES OF KINETWORKS™

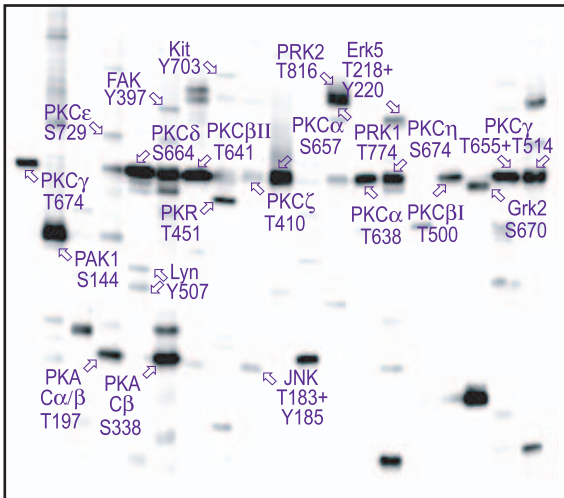
KPSS 11.0 PHOSPHO-SITE SCREENS



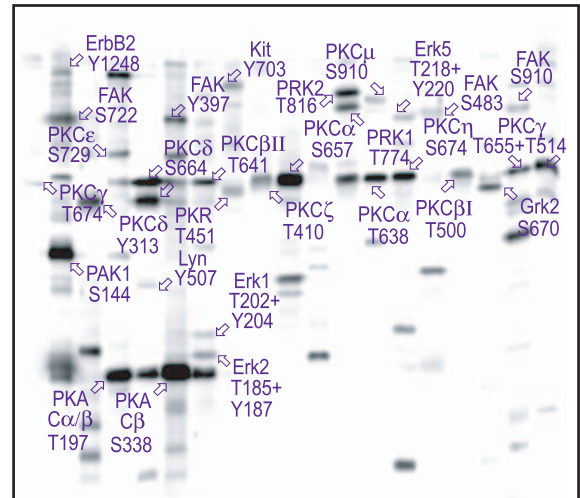
Human A431 Epidermoid Carcinoma Cells



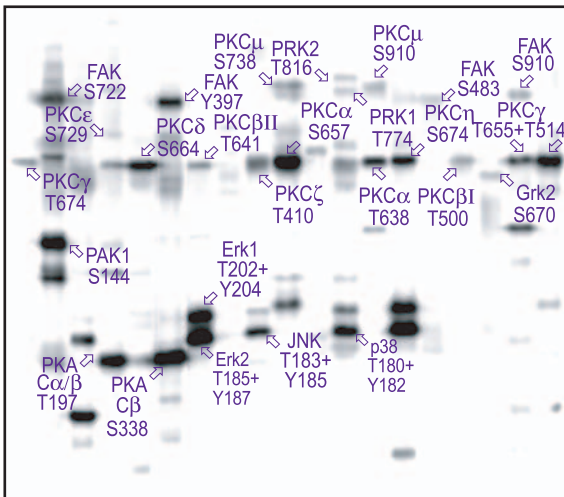
Human HeLa Cervical Carcinoma Cells



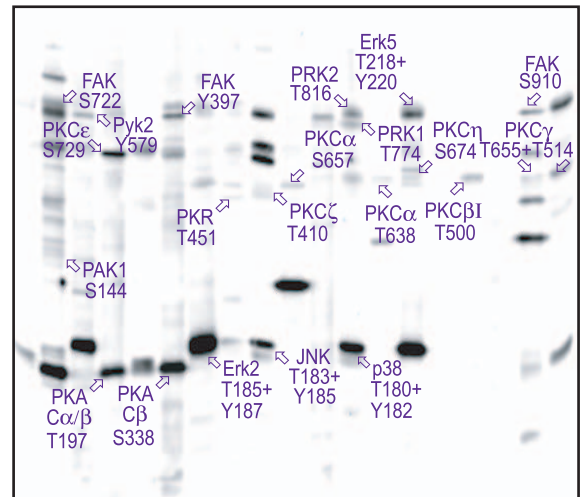
Human Jurkat T Cells



Mouse Pam212 Keratinocytes



Mouse NIH3T3 Fibroblasts



Xenopus laevis Maturing Oocytes

Suite 402, 6190 Agronomy Road, Vancouver, BC, Canada V6T 1Z3

Tel: 604 822 9963 • Fax: 604 822 9964 • www.kinexus.ca

