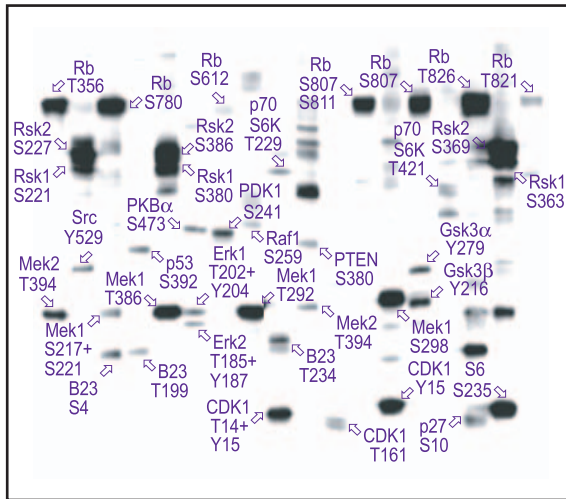
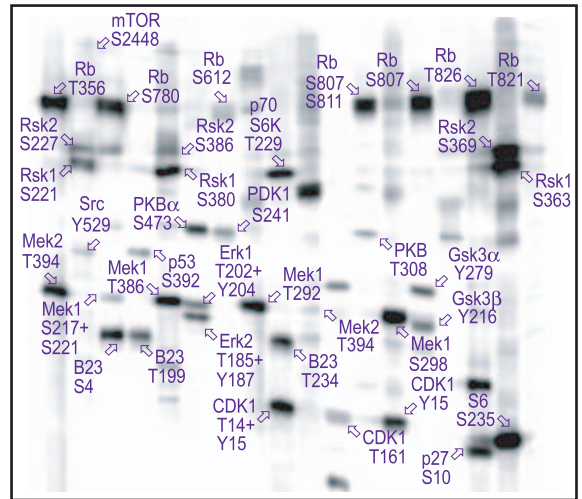


# EXAMPLES OF KINETWORKS™

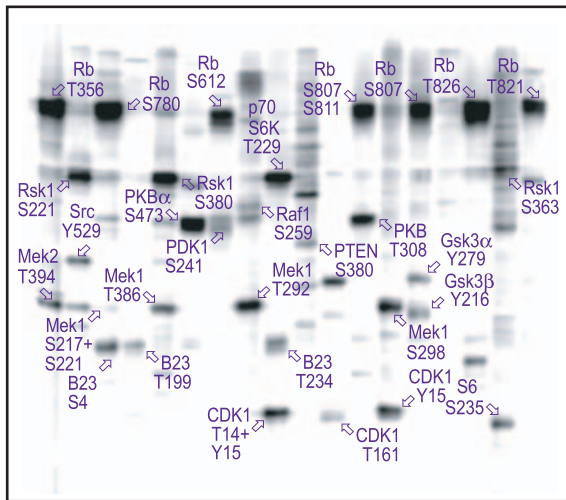
## KPSS 10.0 PHOSPHO-SITE SCREENS



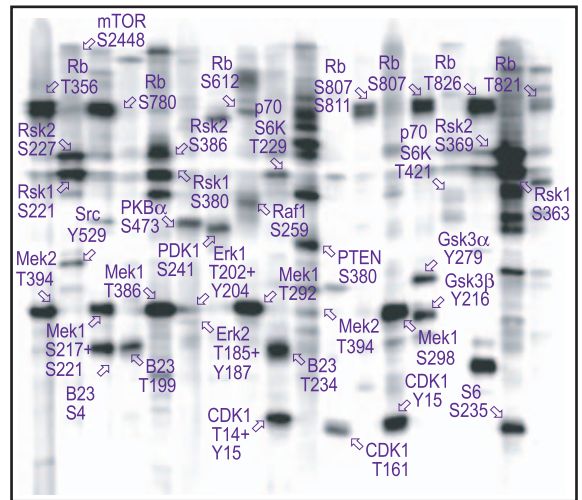
**Human A431 Epidermoid Carcinoma Cells**



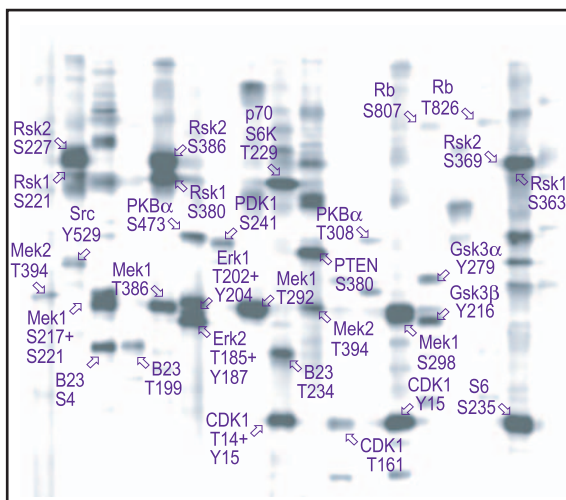
**Human HeLa Cervical Carcinoma Cells**



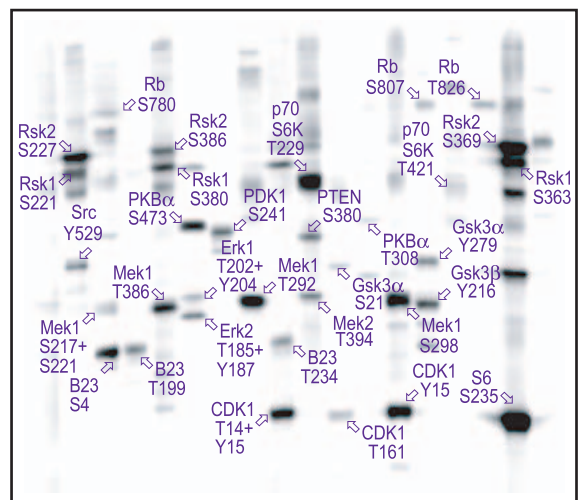
**Human Jurkat T Cells**



**Human HCT116 Colon Carcinoma Cells**



**Mouse NIH-3T3 Fibroblasts**



**Mouse Pam212 Keratinocytes**

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

Tel: 604 822 9963 • Fax: 604 822 9964 • [www.kinexus.ca](http://www.kinexus.ca)