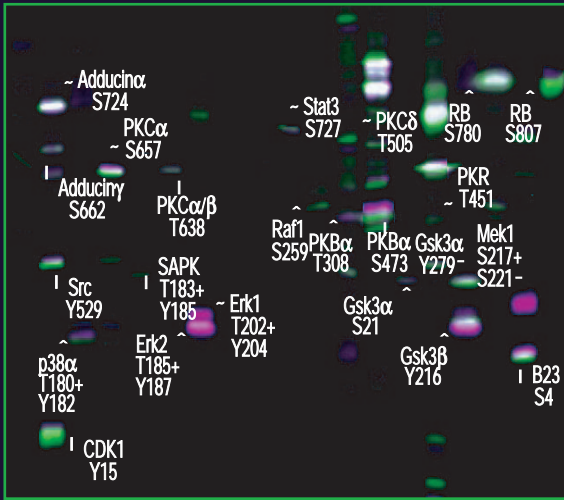
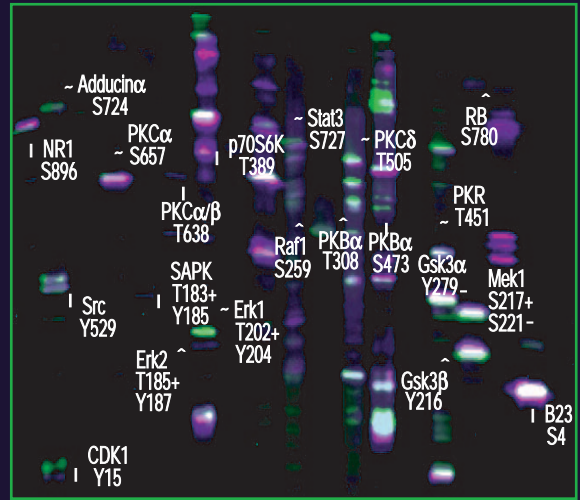


EXAMPLES OF KINETWORKS^a

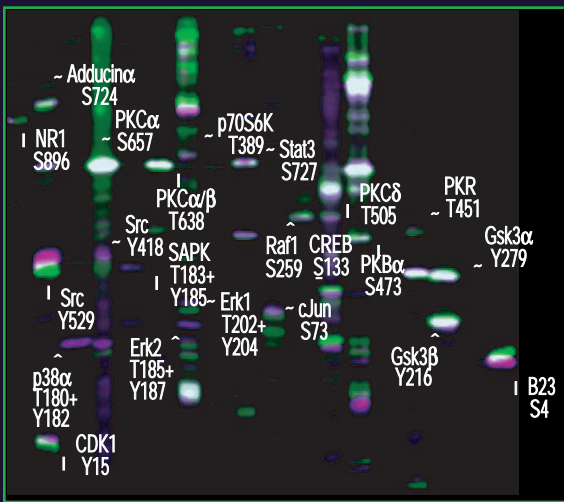
KPSS 1.1 PHOSPHO-SITE SCREENS



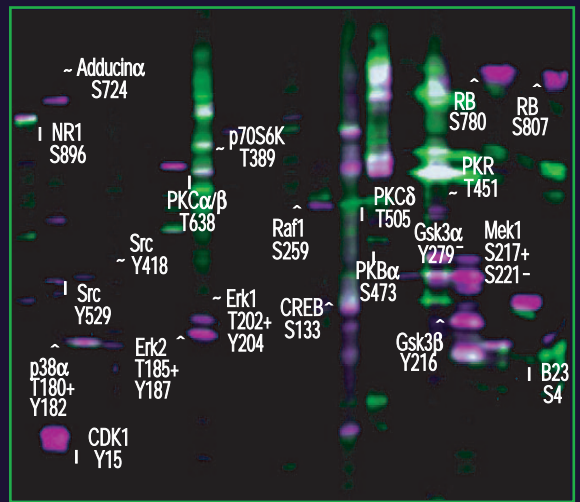
Human CaCo Colon Carcinoma Cells
+ 20 ng/ml Epidermal Growth Factor for 5 min



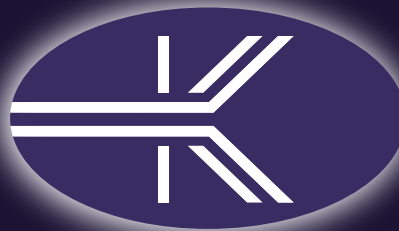
Human HCT-116 Colon Carcinoma Cells
+ 200 ng/ml Nocadazole for 72 h



Mouse Pam-212 Keratinocytic Cells
+ 1 µg/ml Anisomycin for 6 h



Human THP1 Thymocyte Cells
+ 10 nM Phorbol Ester for 1 h



Effect of Treatment
on Phosphorylation

- Increased
- Unchanged
- Decreased