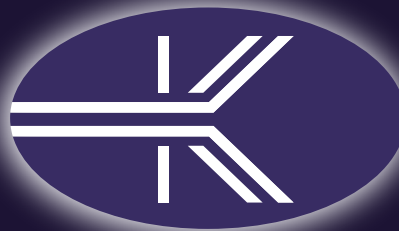
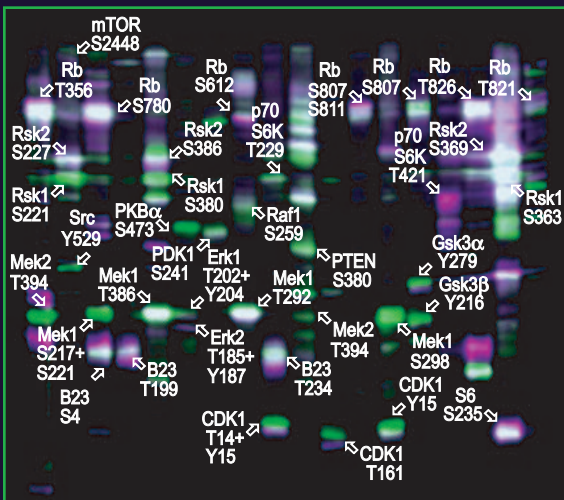
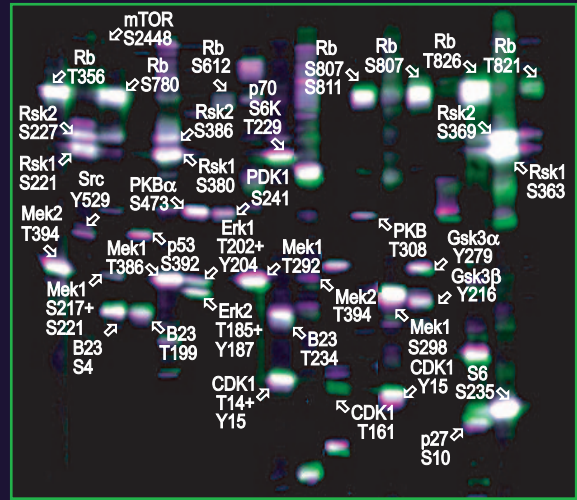


# EXAMPLES OF KINETWORKS™ KPSS 10.0 PHOSPHO-SITE SCREENS



**Effect of Treatment on Phosphorylation**

- Increased
- Unchanged
- Decreased