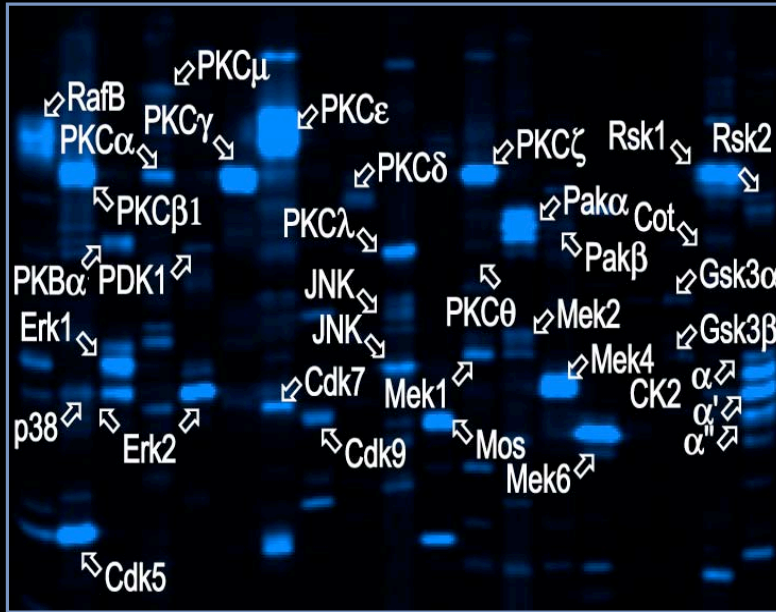


Kinetworks™ Multi-Immunoblotting



*Mouse Brain KPKS-1.2
Protein Kinase Screen Gel A*

- **Accurate** - based on immunoreactivity and M.W.
- **Quantitative** - ECL signals over 2000-fold range
- **Sensitive** - ~350 µg cell lysate protein needed
- **Reproducible** - ±10-20% S.D.
- **Broad utility** - works in human, rat, mouse, etc.