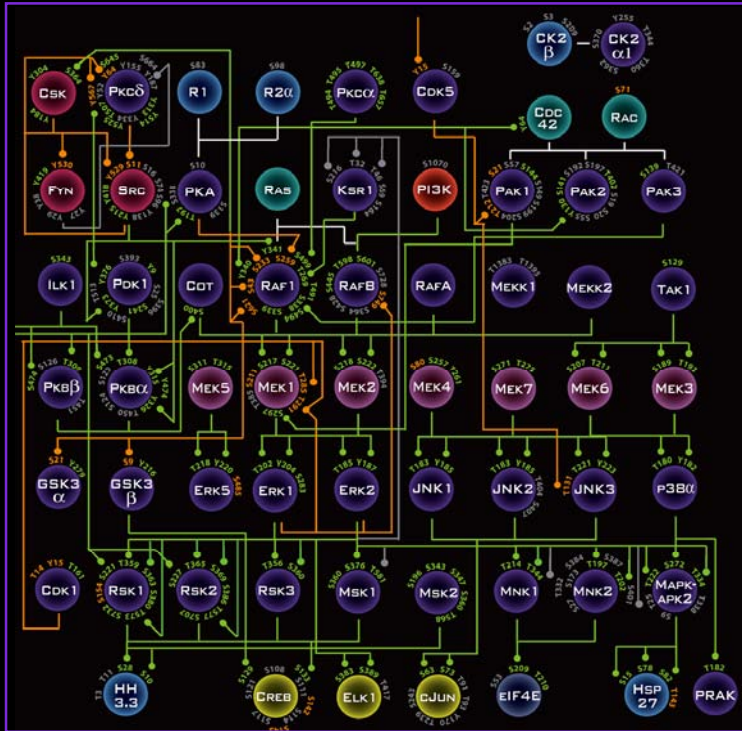


# Protein Kinases are Major Drug Targets



- More than 30% of pharmaceutical drug discovery is focused on protein kinases
- Over 400 human diseases linked to defects in kinase signalling
- Only ~ 24 of 515 protein kinases targeted for drug screening in humans