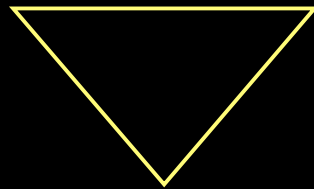
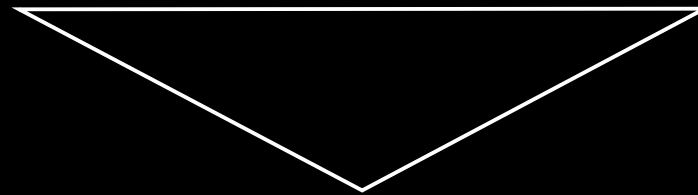


Human Protein Phosphorylation By Kinases

PROTEOME



~30,000 total protein genes



KINEOME

~515 protein kinase genes



>1000 catalytically active
protein kinase isoforms

**PHOSPHO-
PROTEOME**

>1,000,000 phospho-sites