

# Gene Families in Human Genome

■ Metabolic enzymes	3142	(9.2%)
■ Signalling proteins	6166	(21%)
■ Protein kinases	515	(1.7%)
■ Protein phosphatases	138	(0.5%)
■ Transcription factors	1850	(6.0%)
■ Receptors	1543	(5.0%)
■ Extracellular mediators	376	(1.2%)
■ Other signalling proteins	1722	(5.6%)

~30,000 genes (100%)

\*based on Venter et al. (2001) Science