

Kinases Genes in Diverse Genomes

	Kinase Genes / Total Genes		
■ Human	515	/	30,000 (1.7%)
■ Fly	319	/	13,338 (2.4%)
■ Worm	437	/	18,266 (2.4%)
■ Yeast	122	/	6,144 (2.0%)
■ Mustard plant	1049	/	25,706 (4.1%)

*based on Intern. HGS Consortium (2001) Nature