

Human Genome Sequencing

- Human genome is 2.9 billion base pairs long
- 50 terabytes of data = 80,000 compact disks worth of data
- Only about 3-4% of the genome encodes genes

