



Volume occupied by major intracellular compartments in a typical cell

Intracellular Compartment	Percent of Total Cell Volume	Approximate Number per Cell
Cytosol	54	1
Mitochondria	22	1700
Rough ER cisternae	9	3
Smooth ER cisternae + Golgi	6	6
Nucleus	6	1
Peroxisomes	1	400
Lysosomes	1	300
Endosomes	1	200