



The Cellular Landscape

Created for **Cell Signaling Technology** by Digizyme, and inspired by the stunning art of David Goodsell, this 3D rendering of a eukaryotic cell is modeled using X-ray, nuclear magnetic resonance (NMR), and cryo-electron microscopy datasets for all of its molecular actors. It is an attempt to recapitulate the myriad pathways involved in signal transduction, protein synthesis, endocytosis, vesicular transport, cell-cell adhesion, apoptosis, and other processes. Although dilute in its concentration relative to a real cell, this rendering is also an attempt to visualize the great complexity and beauty of the cell's molecular choreography. Explore: www.cst-science.com/cellularlandscapes

