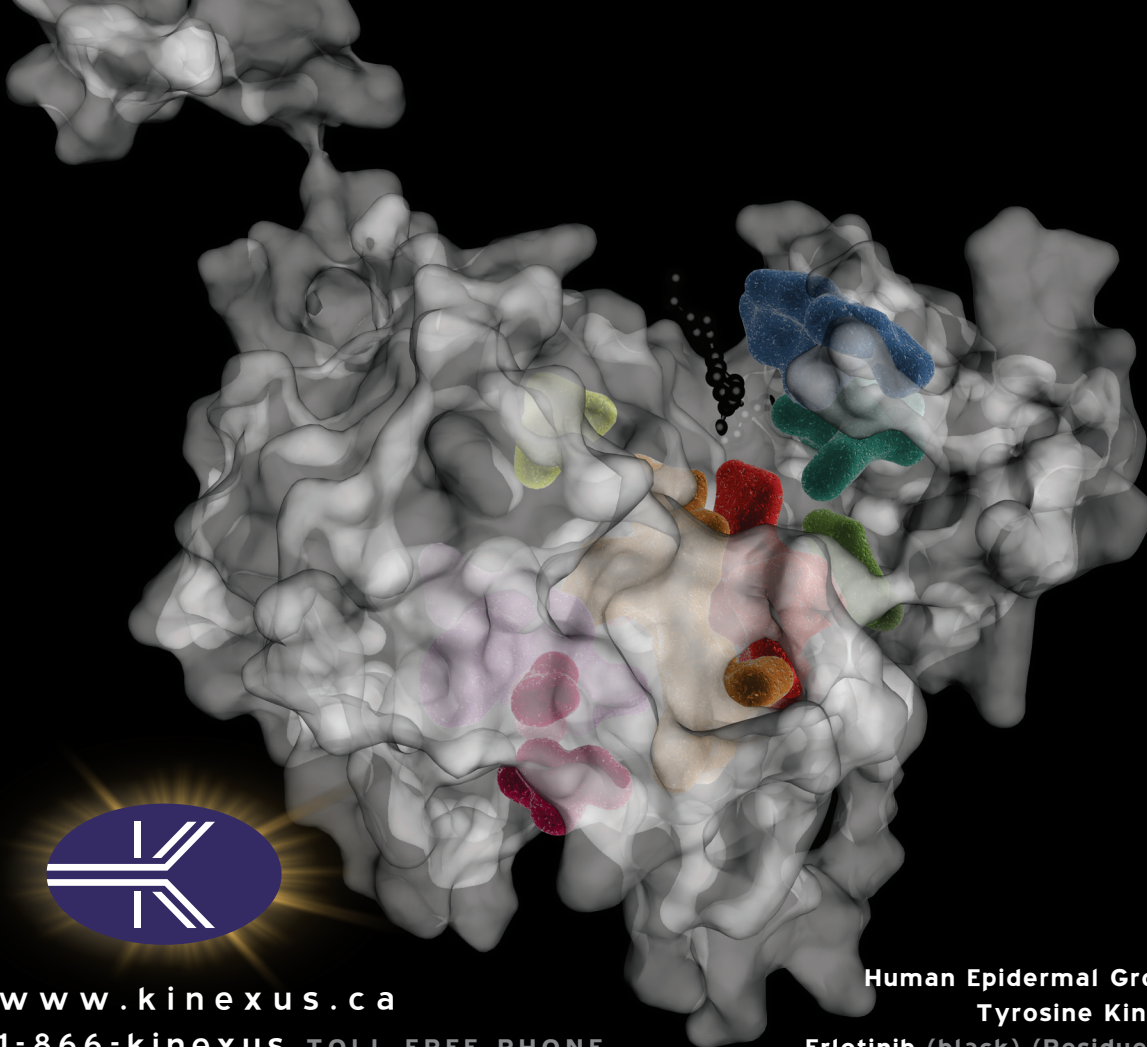




www.kinexus.ca
1-866-kinexus TOLL FREE PHONE

Human Epidermal Growth Factor Receptor
Tyrosine Kinase Catalytic Domain
Erlotinib (black) (Residues 671-998) PDB: 1M17

Protein Kinase Catalytic Domains

