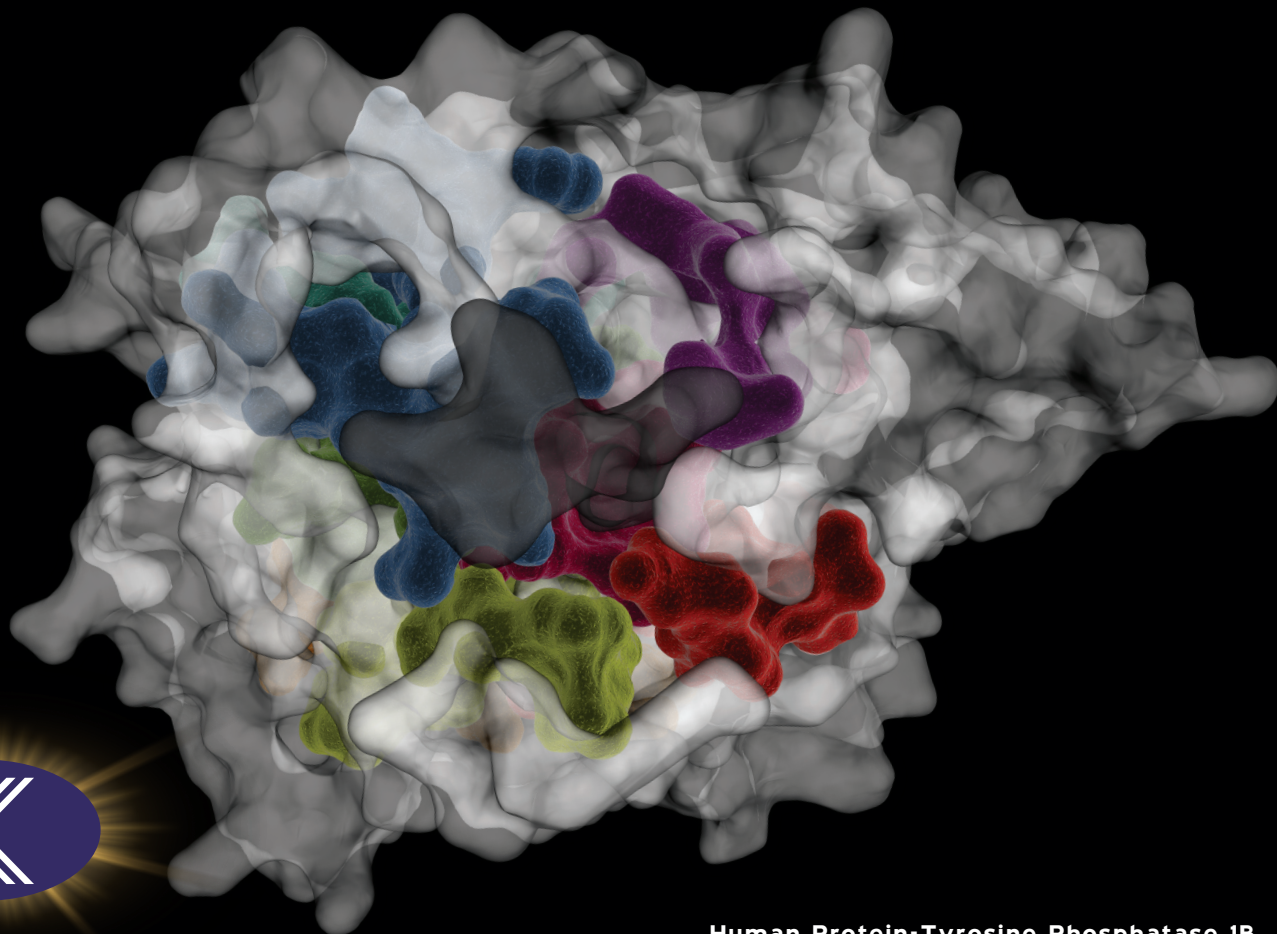




www.kinexus.ca

1-866-kinexus TOLL FREE PHONE



Human Protein-Tyrosine Phosphatase 1B
complexed with ETDPYRKGKGKLL peptide from
Insulin Receptor (dark grey) (Residues 1-299) PDB: 1G1G

Protein-Tyrosine Phosphatase Catalytic Domains

