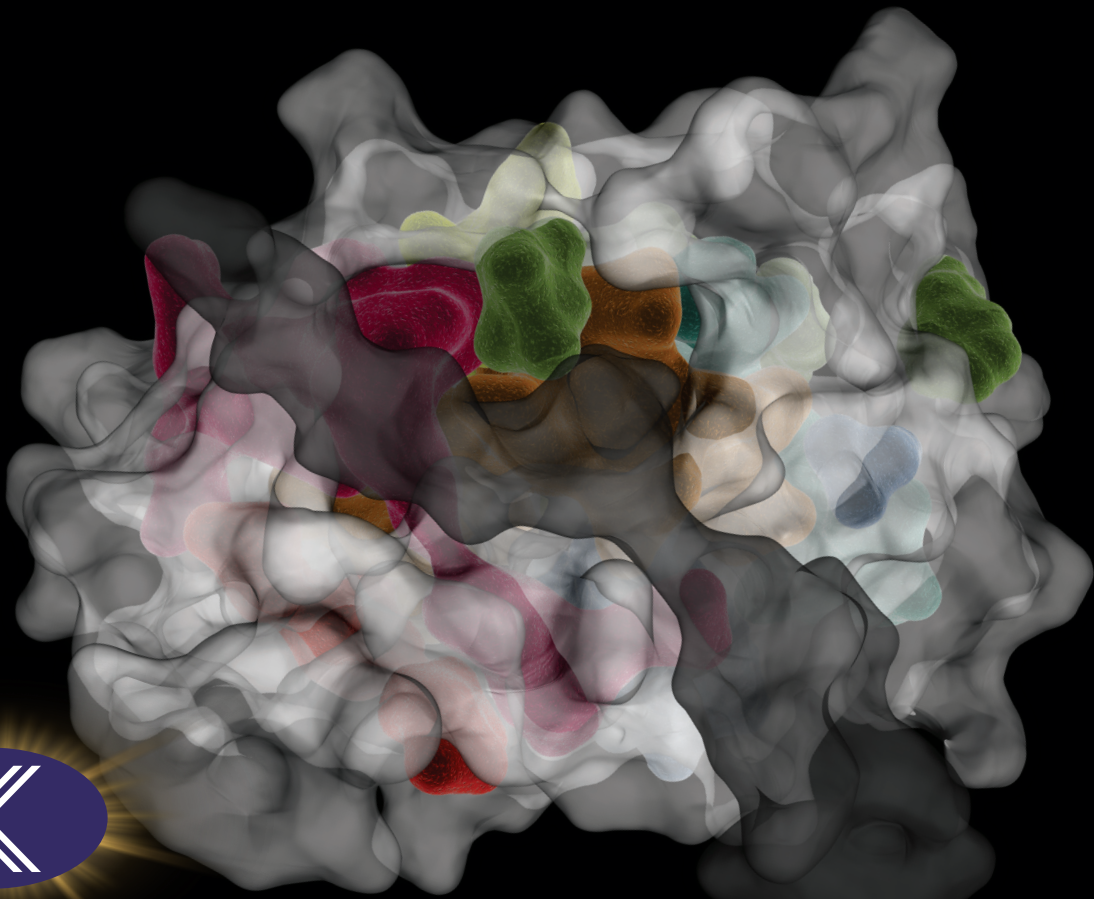




www.kinexus.ca

1-866-kinexus TOLL FREE PHONE

Human VHR-Related Dual-Specificity Phosphatase VHR
complexed with DDEXGpYVATR substrate peptide (grey)
(Residues 2-185) PDB: 1J4X

Protein-Dual Phosphatase Catalytic Domains

