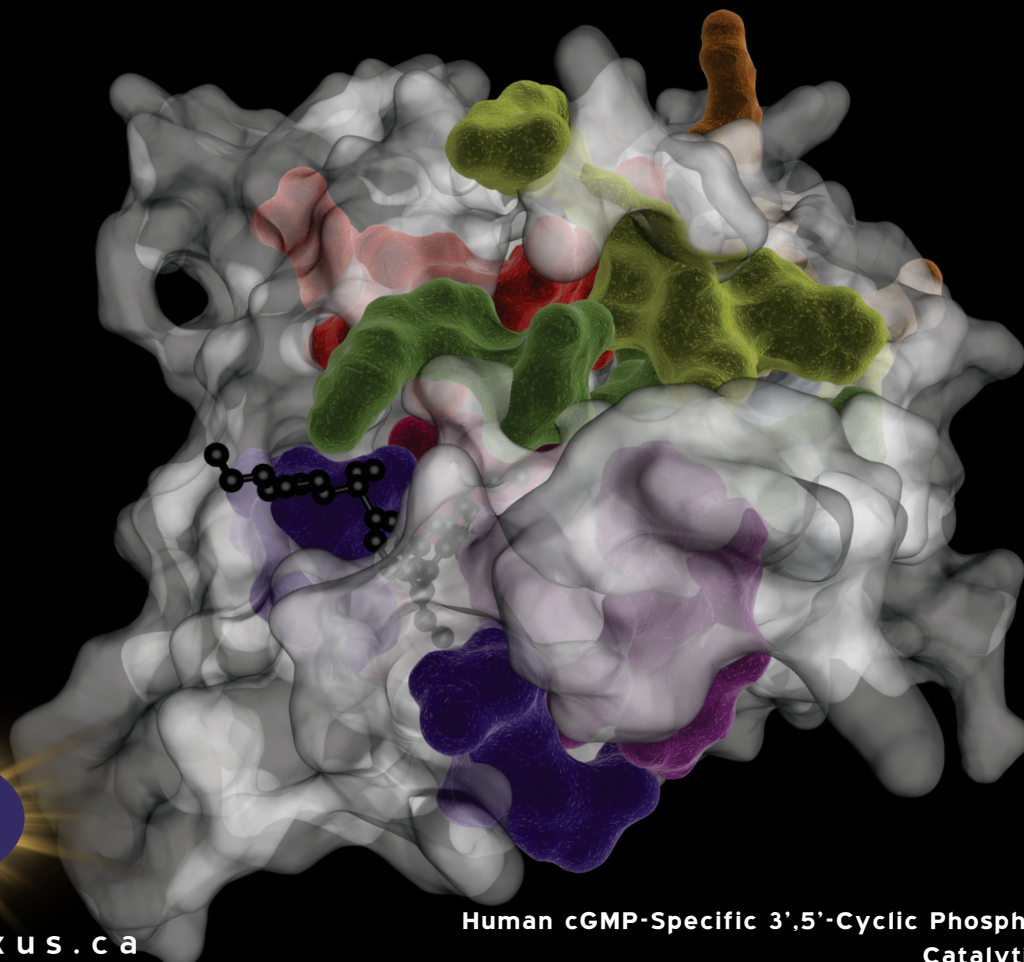




www.kinexus.ca

1-866-kinexus TOLL FREE PHONE



Human cGMP-Specific 3',5'-Cyclic Phosphodiesterase
Catalytic Fragment

Levitra (black) (Residues 537-860) PDB: 1UHO

Cyclic Nucleotide Phosphodiesterase Catalytic Domains

