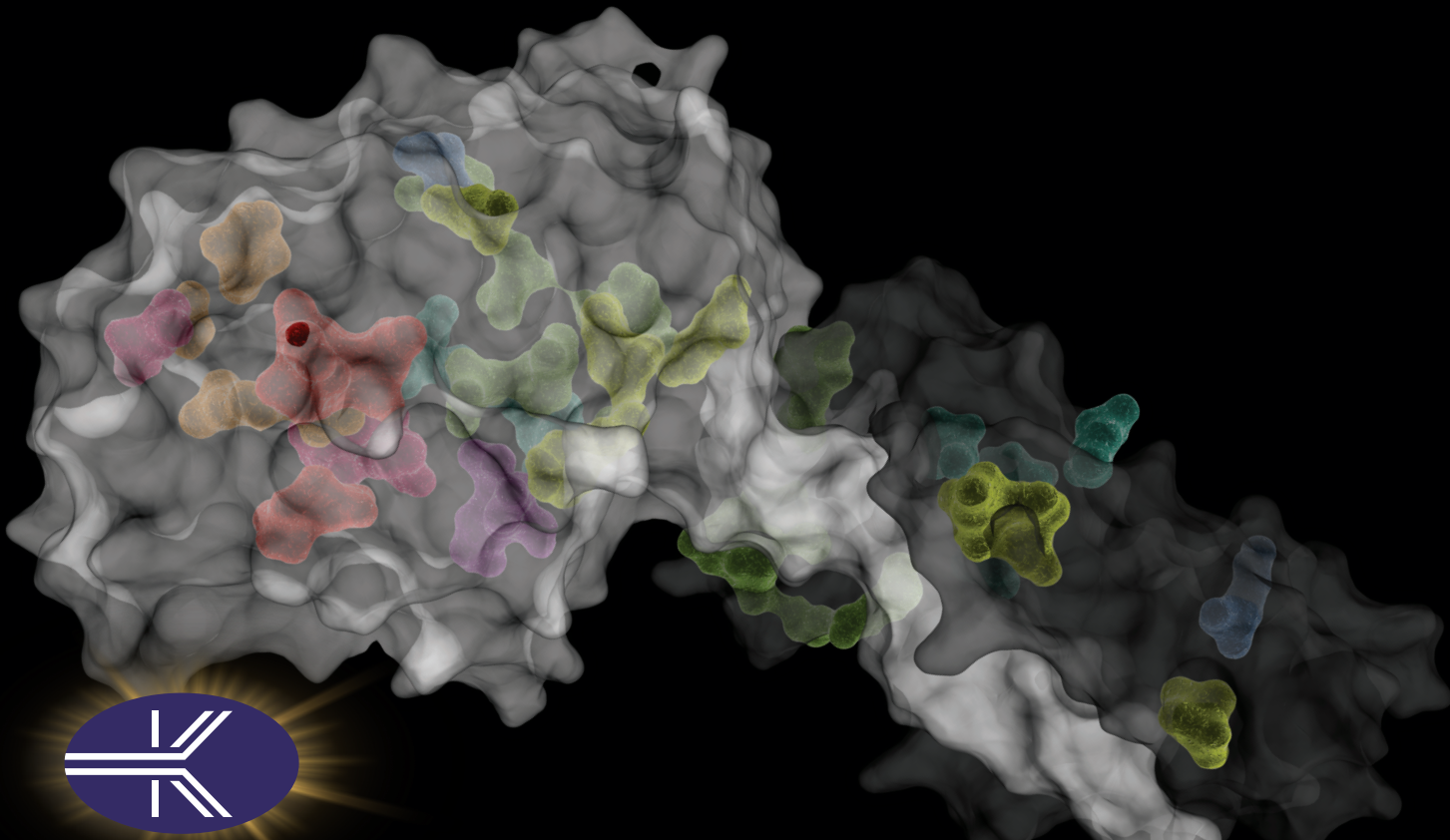




www.kinexus.ca

1-866-kinexus TOLL FREE PHONE

Human Protein-Serine Phosphatase 2B Calcineurin
complexed with Calmodulin (grey)
(Residues 2-185) PDB: 1AUI

Protein-Serine Phosphatase Catalytic Domains

