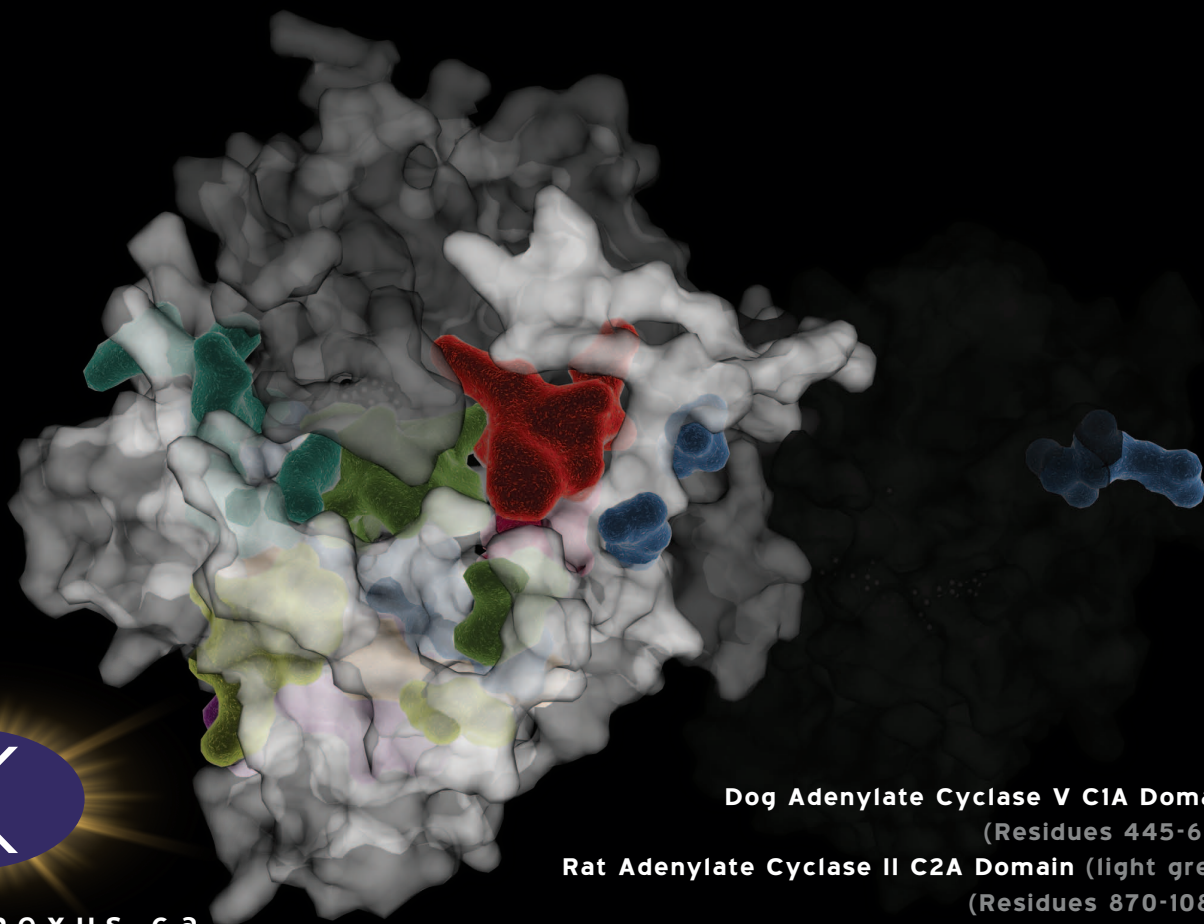




www.kinexus.ca

1-866-kinexus TOLL FREE PHONE



Dog Adenylate Cyclase V C1A Domain
(Residues 445-661)

Rat Adenylate Cyclase II C2A Domain (light grey)
(Residues 870-1082)

Cow G Protein Gsα Subunit (dark grey)
(Residues 1-394) PDB: 1CJK

Adenylate and Guanylate Cyclase Catalytic Domains

