



www.kinexus.ca



KINEXUS

Foreground [pdb 1CS4]: Adenylate cyclase C1A (yellow); & C2A (tan) domains & Gs-alpha subunit (teal).
Midground, right [pdb 1OMW]: G protein-coupled receptor kinase 2 (pink); G-beta subunit (green); & G-gamma subunit (ice blue); Cyclin A (blue); & p27 KIP1 CDK inhibitor (yellow). Background [pdb 1GOT]: Gs-alpha (teal); G-beta (green); & G-gamma (ice blue) subunits. ATP molecules (grey).