



www.kinexus.ca



KINEXUS

14 Calmodulin-dependent protein kinase 2 subunits (pink) complexed with 14 Calmodulin subunits (blue) using pdb files 1HKX, 1COM & 1PHK (substituting for the catalytic domain with phosphorylase kinase gamma subunit).