



www.kinexus.ca

1-866-kinexus TOLL FREE PHONE

Pig Phosphatidylinositol 3-Kinase p110

Catalytic Subunit

(Residues 141-1102) PDB: 1E7V

PI3K-Related Kinase Catalytic Domains

