
PIP5K1C

Product code: 11-113

Full-length human PIP5K1C [1-668(end) amino acids of accession number NP_036530] was expressed as N-terminal GST-fusion protein (101 kDa) using baculovirus expression system.

Assay platform : ADP-Glo_Lipid

Substrate : PI(4)P

ATP (μ M) Km app / Bin : 33 / 50

Metal : Mg

Reference compound : AG-183

IC50 at ATP Bin (nM) : 1900

IC50 at 1 mM ATP (nM) : n.a.