
DGK α (DGKA)

Product code: 12-401-20N

Full-length human DGK α [1-735(end) amino acids of accession number NP_958852.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (87 kDa) using baculovirus expression system.

Assay platform : ADP-Glo_Lipid

Substrate : Diacylglycerol

ATP (μ M) Km app / Bin : 130 / 100

Metal : Mg

Reference compound : Non-disclosable

IC50 at ATP Bin (nM) : Non-disclosable

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