
DGK β (DGKB)

Product code: 12-402-20N

Full-length human DGK β [1-803(end) amino acids of accession number NP_001337638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (95 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and activated with ATP. Activated protein was purified by using gel filtration.

Assay platform : ADP-Glo_Lipid

Substrate : Diacylglycerol

ATP (μ M) Km app / Bin : 61 / 50

Metal : Mg

Reference compound : Non-disclosable

IC50 at ATP Bin (nM) : Non-disclosable

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