
DGK γ (DGKG)

Product code: 12-403-20N

Full-length human DGK γ [1-791(end) amino acids and T142S and R316K of accession number NP_001337.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (94 kDa) using baculovirus expression system.

Assay platform : ADP-Glo_Lipid

Substrate : Diacylglycerol

ATP (μ M) Km app / Bin : 55 / 50

Metal : Mg

Reference compound : Non-disclosable

IC50 at ATP Bin (nM) : Non-disclosable

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