
DGK κ (DGKK)

Product code: 12-408-20N

Full-length human DGK κ [1-1271(end) amino acids and D1118N of accession number NP_001013764.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (146 kDa) using baculovirus expression system.

Assay platform : ADP-Glo_Lipid

Substrate : Diacylglycerol

ATP (μ M) Km app / Bin : 17 / 25

Metal : Mg

Reference compound : Non-disclosable

IC50 at ATP Bin (nM) : Non-disclosable

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