
DGK η (DGKH)

Product code: 12-406-20N

Catalytic domain of human DGK η [1-1147 amino acids of accession number NP_821077.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (131 kDa) using baculovirus expression system.

Assay platform	: ADP-Glo_Lipid
Substrate	: Diacylglycerol
ATP (μ M) Km app / Bin	: 24 / 25
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	: Non-disclosable
IC50 at 1 mM ATP (nM)	: n.a.
