
DGK ζ (DGKZ)

Product code: 12-410-20N

Full-length human DGK ζ [1-929(end) amino acids of accession number NP_003637.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (109 kDa) using baculovirus expression system.

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