
DGK ϵ (DGKE)

Product code: 12-415-20N

Truncated human DGK ϵ [48-567(end) amino acids of accession number NP_003638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (63 kDa) using baculovirus expression system.

Assay platform	:	ADP-Glo_Lipid
Substrate	:	Diacylglycerol
ATP (μ M) Km app / Bin	:	120 / 100
Metal	:	Mg
Reference compound	:	Non-disclosable
IC50 at ATP Bin (nM)	:	Non-disclosable
IC50 at 1 mM ATP (nM)	:	n.a.
