
DGKθ(DGKQ)

Product code: 12-409-20N

Full-length human DGKθ [1-942(end) amino acids of accession number NP_001338.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (106 kDa) using baculovirus expression system.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP (μM) Km app / Bin	: 37 / 50
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	: Non-disclosable
IC50 at 1 mM ATP (nM)	: n.a.
