
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. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform : ADP-Glo
Substrate : Diacylglycerol

ATP (μM) Km app / Bin : 37 / 50
Metal : Mg
Reference compound : Non-disclosable
IC50 at ATP Bin (nM) :
IC50 at 1 mM ATP (nM) :
