
DGKy(DGKG)

Product code 12-403-20N

Full-length human DGKy [1-791(end) amino acids and T142S and R316K of accession number NP_001337.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (94 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

Assay platform : ADP-Glo
Substrate : Diacylglycerol

ATP (μ M) Km app / Bin : 55 / 50
Metal : Mg
Reference compound : Non-disclosable
IC50 at ATP Bin (nM) :
IC50 at 1 mM ATP (nM) :
