
DGK γ (DGKG)

Product code: 12-403-20N

Full-length human DGK γ [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.

Assay platform : ADP-Glo
Substrate : Diacylglycerol

ATP (μ M) K_m app / B_{in} : 55 / 50
Metal : Mg
Reference compound : Non-disclosable
IC₅₀ at ATP B_{in} (nM) : Non-disclosable
IC₅₀ at 1 mM ATP (nM) : n.a.
