

---

## DGKδ(DGKD)

Product code 12-404-20N

Catalytic domain of human DGKδ [1-1141 amino acids of accession number NP\_690618.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (131 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 : 120 / 100  
Metal : Mg  
Reference compound : Non-disclosable  
IC50 at ATP Bin (nM) :  
IC50 at 1 mM ATP (nM) :

---