

---

# DGK $\beta$ (DGKB)

Product code 12-402-20N

Full-length human DGK $\beta$  [1-803(end) amino acids of accession number NP\_001337638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (95 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose and activated with ATP. Activated protein was purified by using gel filtration.

Assay platform : ADP-Glo  
Substrate : Diacylglycerol

ATP ( $\mu$ M) Km app / Bin : 61 / 50  
Metal : Mg  
Reference compound : Non-disclosable  
IC50 at ATP Bin (nM) :  
IC50 at 1 mM ATP (nM) :

---