

---

## DGKζ(DGKZ)

Product code 12-410-20N

Full-length human DGKζ [1-929(end) amino acids of accession number NP\_003637.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (109 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 : 25 / 25  
Metal : Mg  
Reference compound : Non-disclosable  
IC50 at ATP Bin (nM) :  
IC50 at 1 mM ATP (nM) :

---