

---

## DGKε(DGKE)

Product code 12-415-20N

Truncated human DGKε [48-567(end) amino acids of accession number NP\_003638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (63 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) :

---