

---

---

## DGK $\delta$ (DGKD)

Product code: 12-404-20N

Catalytic domain of human DGK $\delta$  [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.

Assay platform	: ADP-Glo_Lipid
Substrate	: Diacylglycerol
ATP ( $\mu$ M) Km app / Bin	: 120 / 100
Metal	: Mg
Reference compound	: Non-disclosable
IC50 at ATP Bin (nM)	: Non-disclosable
IC50 at 1 mM ATP (nM)	: n.a.

---

---