

---

---

## DGK $\kappa$ (DGKK)

Product code: 12-408-20N

Full-length human DGK $\kappa$  [1-1271(end) amino acids and D1118N of accession number NP\_001013764.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (146 kDa) using baculovirus expression system.

Assay platform	: ADP-Glo
Substrate	: Diacylglycerol
ATP ( $\mu$ M) $K_m$ app / $B_{in}$	: 17 / 25
Metal	: Mg
Reference compound	: Non-disclosable
IC <sub>50</sub> at ATP $B_{in}$ (nM)	: Non-disclosable
IC <sub>50</sub> at 1 mM ATP (nM)	: n.a.

---

---