

---

---

## DGKI(DGKI)

Product code: 12-407-20N

Full-length human DGKI [1-1065(end) amino acids of accession number NP\_004708.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (122 kDa) using baculovirus expression system.

Assay platform	:	ADP-Glo_Lipid
Substrate	:	Diacylglycerol
ATP ( $\mu$ M) Km app / Bin	:	34 / 50
Metal	:	Mg
Reference compound	:	Non-disclosable
IC50 at ATP Bin (nM)	:	Non-disclosable
IC50 at 1 mM ATP (nM)	:	n.a.

