

---

---

## DGKθ(DGKQ)

Product code: 12-409-20N

Full-length human DGKθ [1-942(end) amino acids of accession number NP\_001338.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (106 kDa) using baculovirus expression system.

Assay platform : ADP-Glo\_Lipid

Substrate : Diacylglycerol

ATP (μM) Km app / Bin : 37 / 50

Metal : Mg

Reference compound : Non-disclosable

IC50 at ATP Bin (nM) : Non-disclosable

IC50 at 1 mM ATP (nM) : n.a.

---

---