
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
Substrate	: Diacylglycerol
ATP (μ 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.
