

## Product Information

### DGKI(DGKI)

Product Number : **12-107**

#### Product description

Full-length human DGKI [1-1065(end) amino acids of accession number NP\_004708.1] was expressed as N-terminal GST-fusion protein (145 kDa) using baculovirus expression system. GST-DGKI was purified by using glutathione sepharose chromatography and gel filtration chromatography.

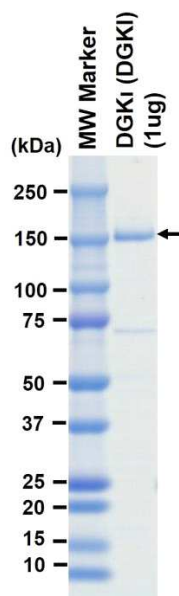
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

Store at -80C.  
Avoid repeating freeze-thaws.

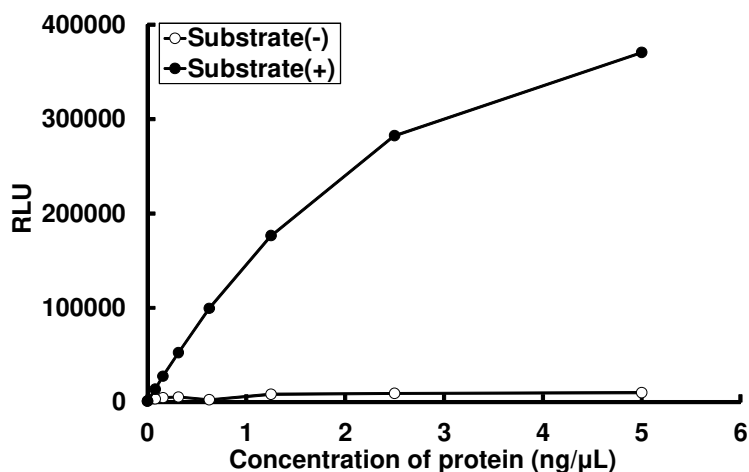
#### SDS-PAGE



Purity: 78 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was determined by ADP-Glo™ Assay. The enzyme was incubated with Lipid substrate and Mg(or Mn)/ATP. The phosphorylation was detected by the ADP-Glo™ Kinase Assay technology (luminescent ADP detection assay).

Substrate : Diacylglycerol and Phosphatidylserine

ATP : 100 μM