

## Product Information

### DGK $\epsilon$ (DGKE)

Product Number : 12-115

#### Product description

Truncated human DGK $\epsilon$  [48-567(end) amino acids of accession number NP\_003638.1] was expressed as N-terminal GST-fusion protein (87 kDa) using baculovirus expression system. GST-DGK $\epsilon$  was purified by using glutathione sepharose chromatography.

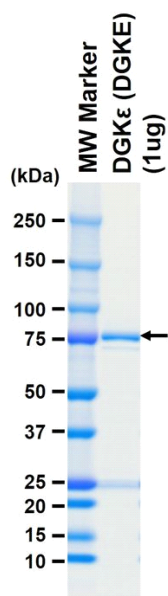
#### Storage buffer:

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.

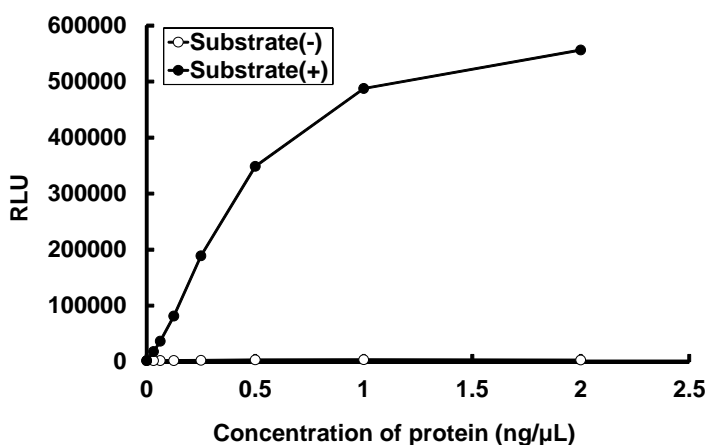
#### SDS-PAGE



Purity: 71 %

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