

Product Information

DGK κ (DGKK)

 Product Number : **12-108**

Product description

Truncated human DGK κ [216-1271 amino acids and D1118N of accession number NP_001013764.1] was expressed as N-terminal GST-fusion protein (148 kDa) using baculovirus expression system. GST-DGK κ was purified by using glutathione sepharose chromatography.

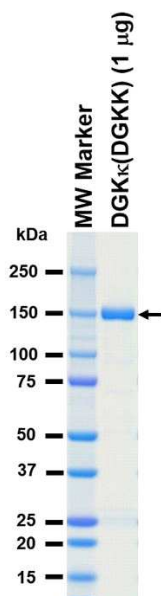
Storage buffer:

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

Storage and Handling:

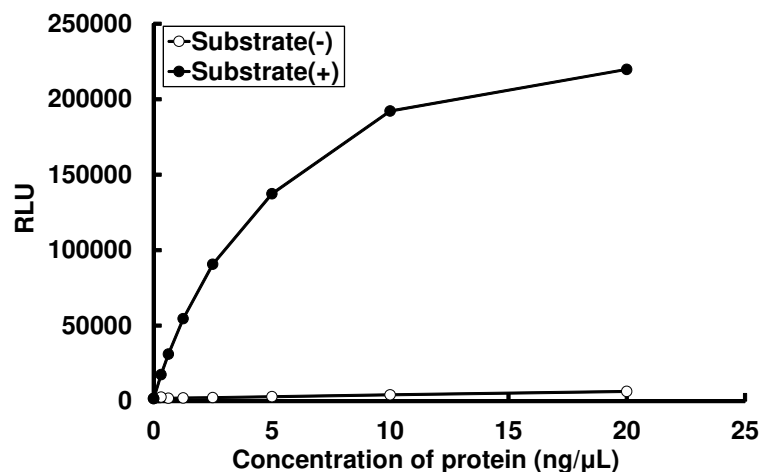
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 90 %
 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