

Product Information

BTN-DGK ζ (DGKZ)

Product Number : **12-410-20N**

Product description

Full-length human DGK ζ [1-929(end) amino acids of accession number NP_003637.2] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (109 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

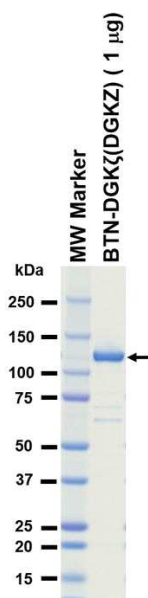
Storage buffer:

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.

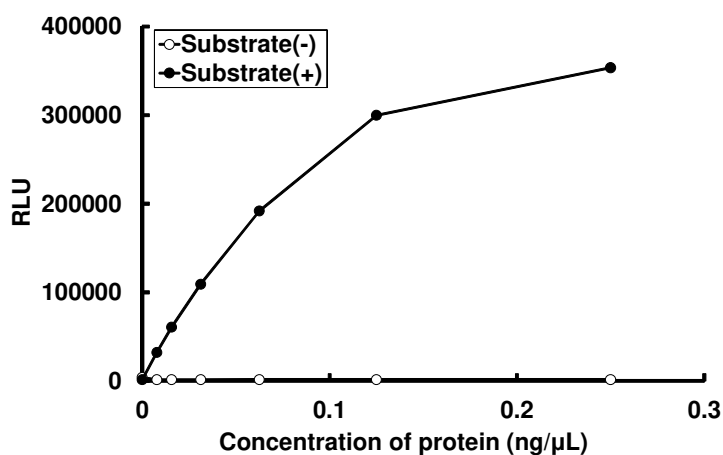
SDS-PAGE



Purity: 92 %

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 μ M