

## Product Information

### BTN-DGK $\beta$ (DGKB)(235-803aa)

Product Number : 12-412-20N

#### Product description

Truncated human DGK $\beta$  [235-803(end) amino acids of accession number NP\_001337638.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (68 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

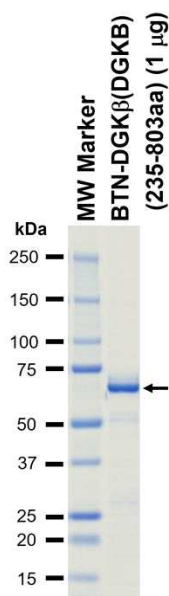
#### Storage buffer:

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.

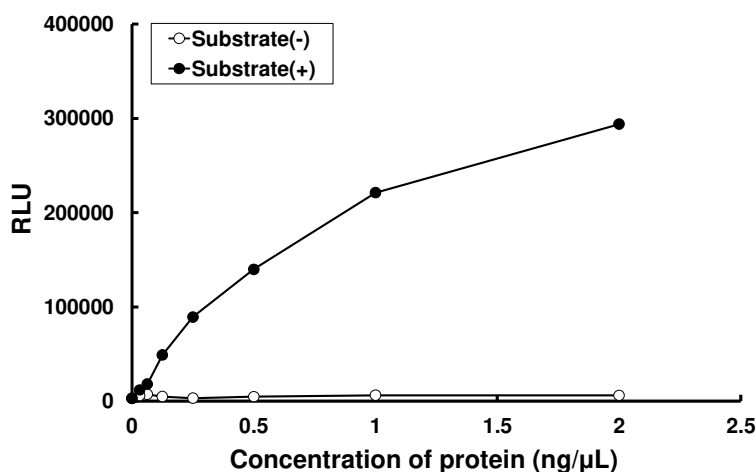
#### SDS-PAGE



Purity: 91 %

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  $\mu$ M