

## Product Information

### BTN-DGK $\eta$ (DGKH)

 Product Number : **12-406-20N**

#### Product description

Catalytic domain of human DGK $\eta$  [1-1147 amino acids of accession number NP\_821077.1] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (131 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

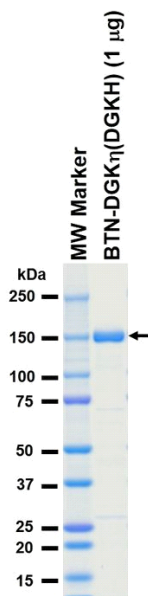
#### Storage buffer:

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.

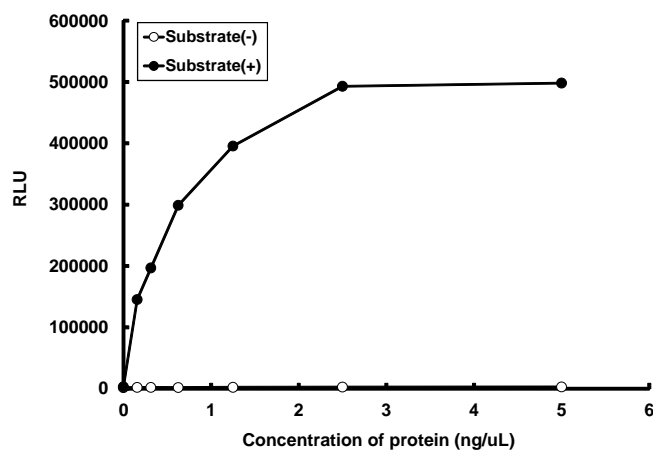
#### SDS-PAGE



Purity: 95 %

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