

## Product Information

### PI4KB

Product Number : 11-131

#### Product description

Full-length human PI4KB [1-828(end) amino acids of accession number NP\_002642.1] was expressed as N-terminal GST-fusion protein (121 kDa) using baculovirus expression system. GST-PI4KB was purified by using glutathione sepharose chromatography.

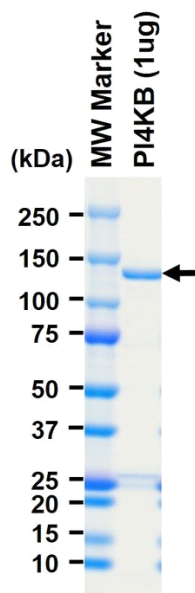
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.01% Brij35,  
 1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

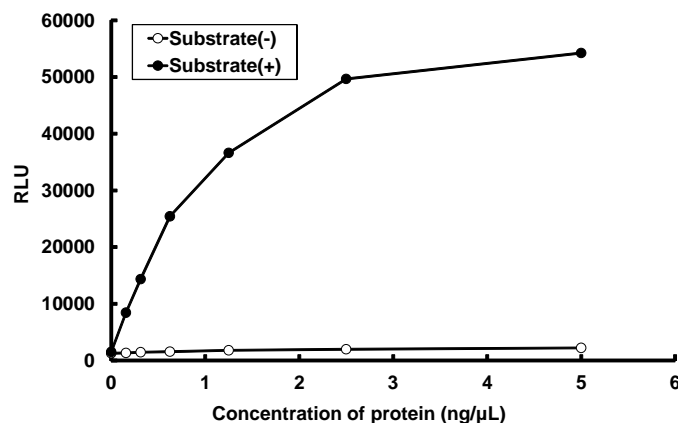
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 70 %  
 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 Mn/ATP. The phosphorylation was detected by the ADP-Glo™ Kinase Assay technology (luminescent ADP detection assay).

Substrate : PI and Phosphatidylserine

ATP : 100 μM