

## Product Information

### PIP4K2A

Product Number : 11-115

#### Product description

Full-length human PIP4K2A [1-406(end) amino acids of accession number NP\_005019] was expressed as N-terminal GST-fusion protein (73 kDa) using baculovirus expression system. GST-PIP4K2A was purified by using glutathione sepharose chromatography.

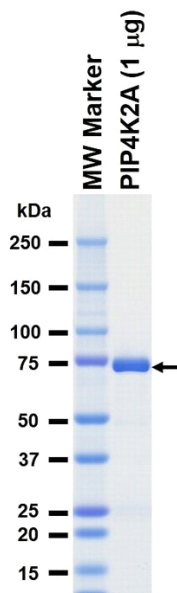
#### Storage buffer:

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.

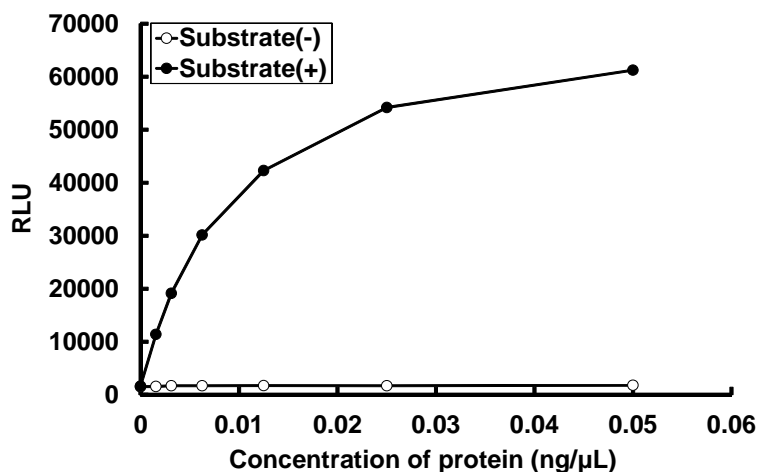
#### SDS-PAGE



Purity: 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 : PI(5)P and Phosphatidylserine

ATP : 100 µM