

Product Information

PIP5K1B

Product Number : 11-112

Product description

Full-length human PIP5K1B [1-540(end) amino acids of accession number NP_003549.1] was expressed as N-terminal GST-fusion protein (88 kDa) using baculovirus expression system. GST-PIP5K1B was purified by using glutathione sepharose chromatography.

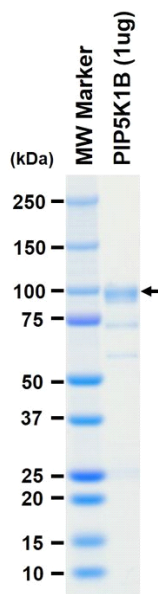
Storage buffer:

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.

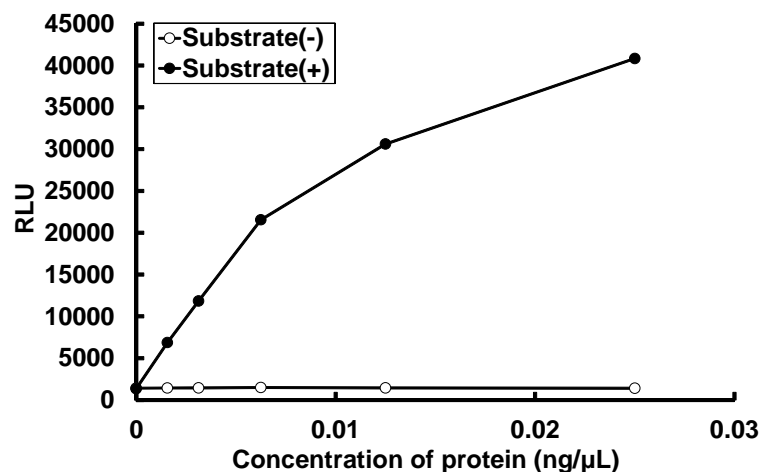
SDS-PAGE



Purity: 68 %

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(4)P and Phosphatidylserine

ATP : 100 μM