

Product Information

PIP5KL1

Product Number : 11-114

Product description

Full-length human PIP5KL1 [1-394(end) amino acids of accession number NP_001128691.1] was expressed as N-terminal GST-fusion protein (72 kDa) using baculovirus expression system. GST-PIP5KL1 was purified by using glutathione sepharose chromatography.

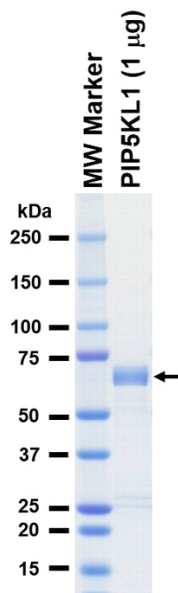
Storage buffer:

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.

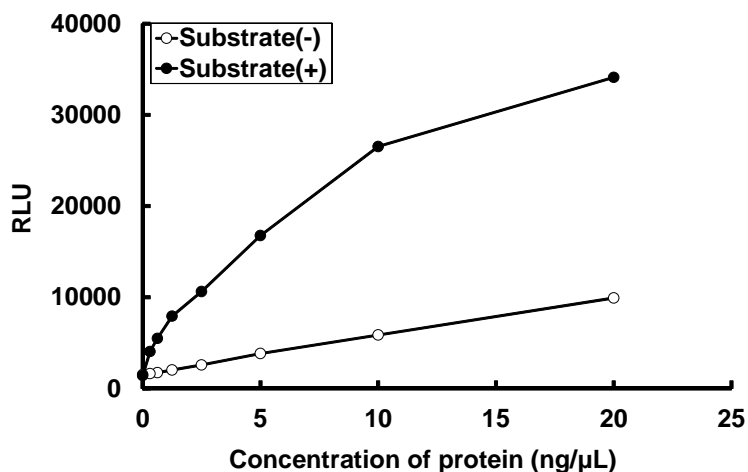
SDS-PAGE



Purity: 87 %

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