

Product Information

PIKFYVE(PIP5K3)

Product Number : 11-118

Product description

Full-length human PIKFYVE [1-2098(end) amino acids and S696N, L932S, Q995L, T998S, S1033A and Q1183K of accession number NP_055855.2] was expressed as N-terminal GST-fusion protein (265 kDa) using baculovirus expression system. GST-PIKFYVE was purified by using glutathione sepharose chromatography.

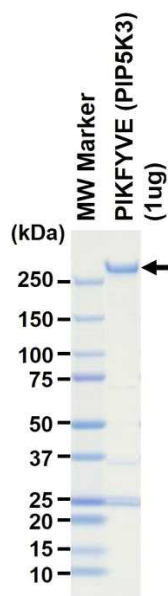
Storage buffer:

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.

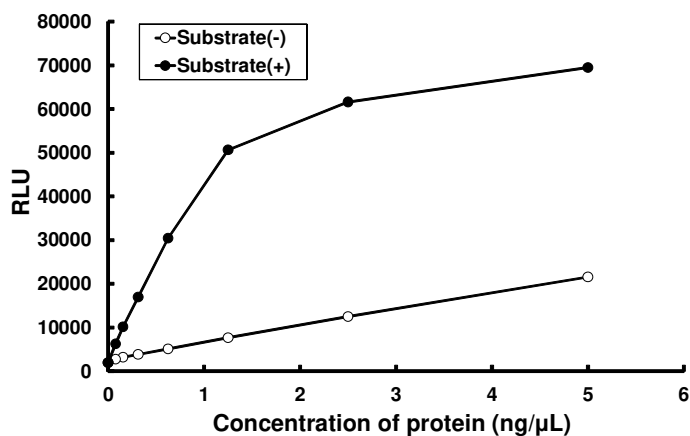
SDS-PAGE



Purity: 57 %

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(3)P and Phosphatidylserine

ATP : 100 μM