

Product Information

PIK3C3

Product Number : 11-122

Product description

Full-length human PIK3C3 [1-887(end) amino acids of accession number NP_002638.2] was expressed as N-terminal GST-fusion protein (129 kDa) using baculovirus expression system. GST-PIK3C3 was purified by using glutathione sepharose chromatography.

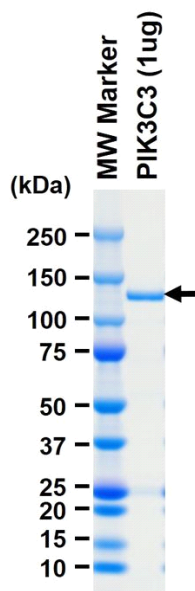
Storage buffer:

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

Storage and Handling:

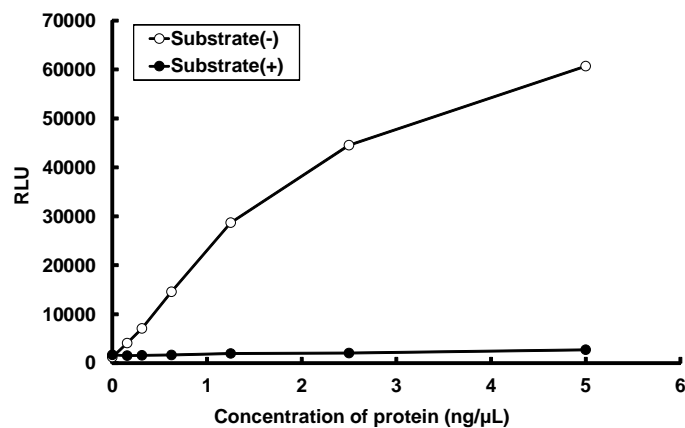
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 86 %
 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