

## Product Information

### PIK3CD/PIK3R1

Product Number : 11-103

#### Product description

Full-length human PIK3CD[1-1044(end) amino acids of accession number NP\_005017.3] was co-expressed as N-terminal GST-fusion protein (147 kDa) with PIK3R1[1-724(end) amino acids of accession number NP\_852664.1] using baculovirus expression system. GST-PIK3CD/PIK3R1 was purified by using glutathione sepharose chromatography.

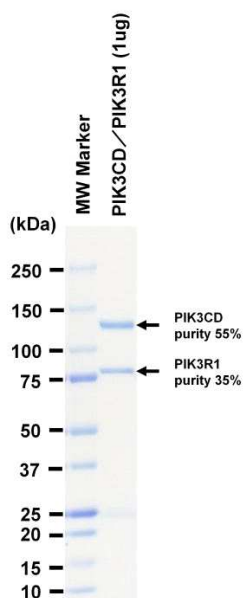
#### Storage buffer:

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

#### Storage and Handling:

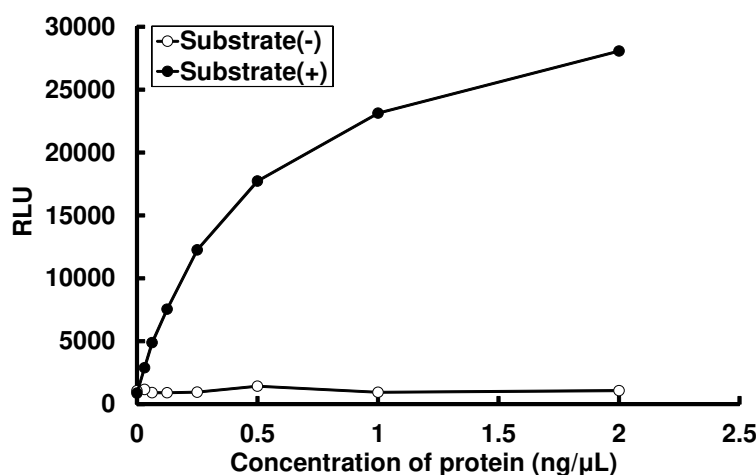
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 90 %  
 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(4,5)P2 and Phosphatidylserine

ATP : 100 μM