

Product Information

BTN-PIK3CA [R88Q]/PIK3R1

Product Number : 11-411-20N

Product description

Full-length human PIK3CA[1-1068(end) amino acids and R88Q of accession number NP_006209.2] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (128 kDa) with PIK3R1[1-724(end) amino acids of accession number NP_852664.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

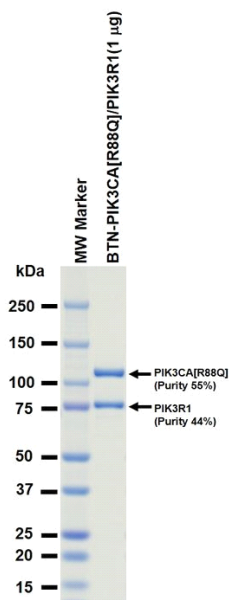
Storage buffer:

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

Storage and Handling:

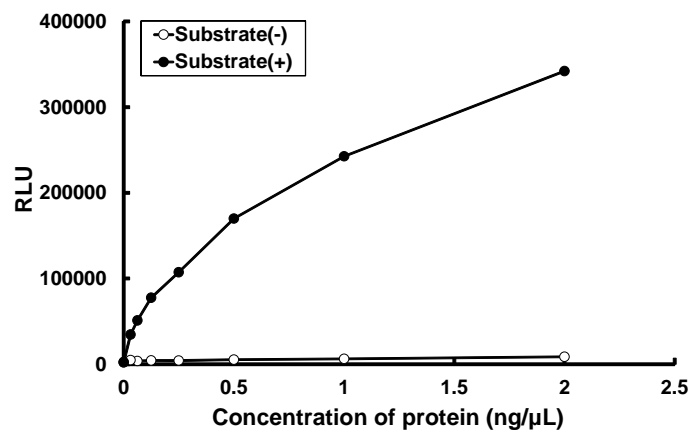
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 99 %
 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