

Product Information

BTN-PKC η (PRKCH)

Product Number : **01-438-20N**

Product description

Full-length human PKC η [1-683(end) amino acids of accession number NP_006246] was expressed as N-terminal DYKDDDDK tagged, biotinylated protein (82 kDa) using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

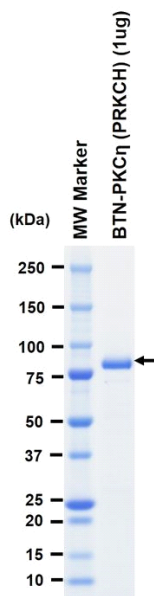
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

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

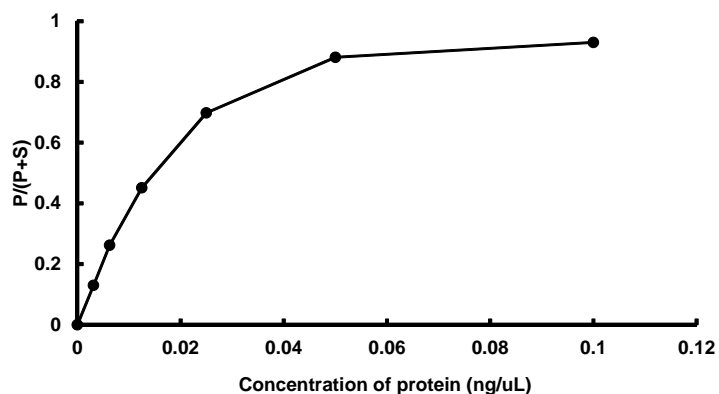
SDS-PAGE



Purity: 97 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate, Mg(or Mn)/ATP, and Lipid Activator. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : PKC peptide

ATP : 100 μ M