

Product Information

PKC β 1(PRKCB1)

 Product Number : **01-134**

Product description

Full-length human PKC β 1 [1-671(end) amino acids of accession number NP_997700.1] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-PKC β 1 was purified by using glutathione sepharose chromatography.

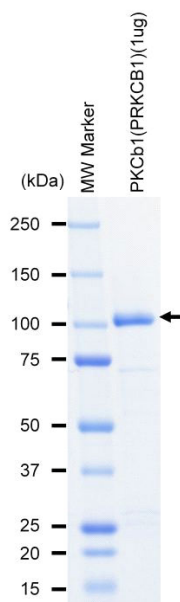
Storage buffer:

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.

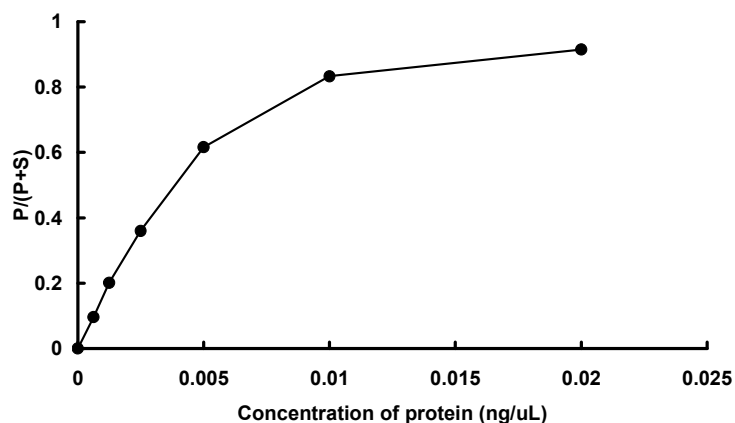
SDS-PAGE



Purity: 92 %

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 Ca/Lipid Activator. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : PKC peptide

 ATP : 100 μ M