

Product Information

PKC α (PRKCA)

 Product Number : **01-133**

Product description

Full-length human PKC α [1-672(end) amino acids and V568I of accession number NP_002728.2] was expressed as N-terminal GST-fusion protein (104 kDa) using baculovirus expression system. GST-PKC α was purified by using glutathione sepharose chromatography.

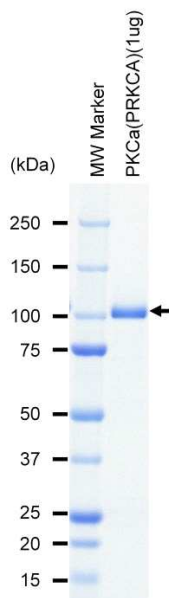
Storage buffer:

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

Storage and Handling:

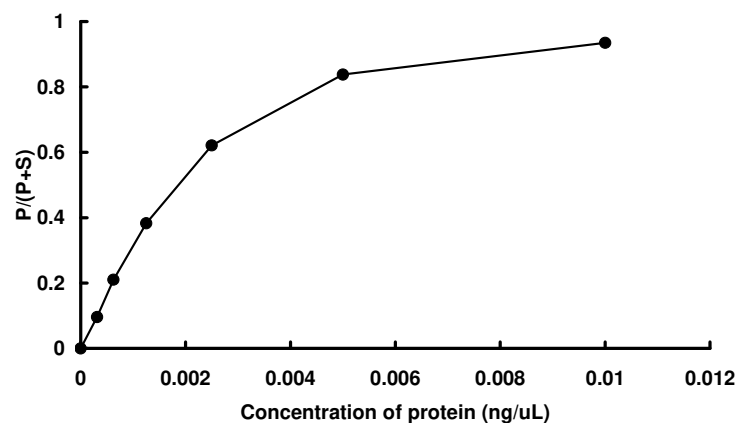
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 96 %
 The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). 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 MSA device.

Substrate : PKC peptide

ATP : 100 μ M