

Product Information

PKCζ(PRKCZ)

 Product Number : **01-141**

Product description

Full-length human PKCζ [1-592(end) amino acids of accession number NP_002735.3] was expressed as N-terminal GST-fusion protein (94kDa) using baculovirus expression system. GST-PKCζ was purified by using glutathione sepharose chromatography.

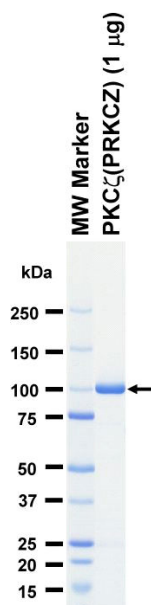
Storage buffer:

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.

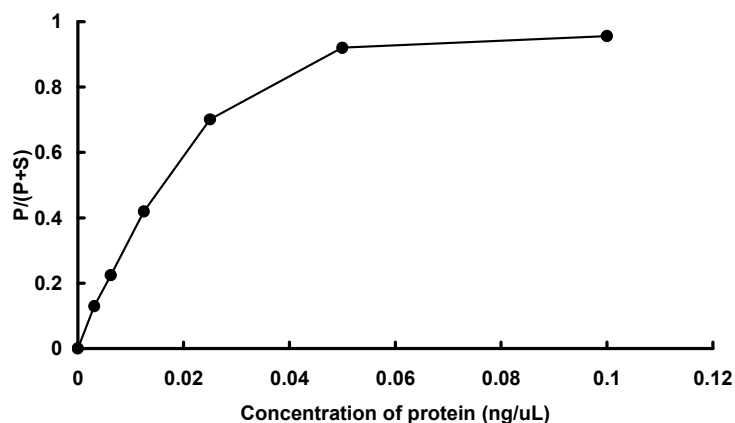
SDS-PAGE



Purity: 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 and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : PKC peptide

ATP : 100 μM