

## Product Information

### PKC $\gamma$ (PRKCG)

 Product Number : **01-137**

#### Product description

Full-length human PKC $\gamma$  [1-697(end) amino acids of accession number NP\_002730.1] was expressed as N-terminal GST-fusion protein (106 kDa) using baculovirus expression system. GST-PKC $\gamma$  was purified by using glutathione sepharose chromatography.

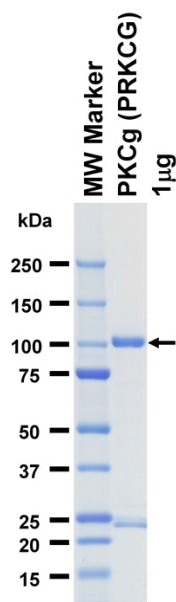
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
 1 mM DTT, 10% glycerol, pH 7.5

#### Storage and Handling:

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

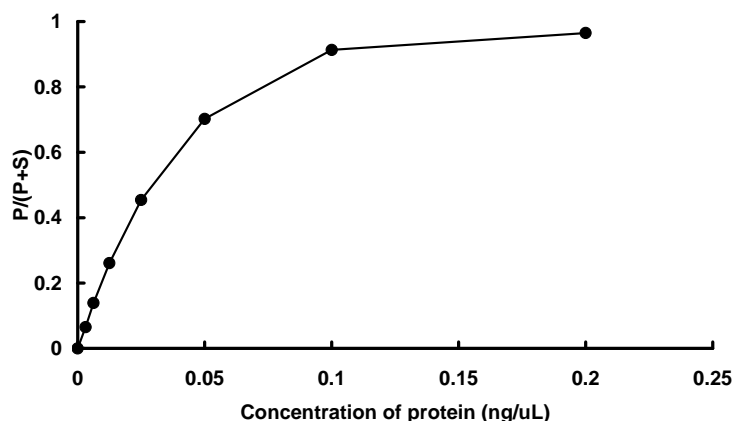
#### SDS-PAGE



Purity: 70 %

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  $\mu$ M