

Product Information

PKN3

 Product Number : **01-146**

Product description

Full-length human PKN3[1-889(end) amino acids of accession number NP_037487.2] was co-expressed as N-terminal GST-fusion protein (127 kDa) with CDC37 and HSP90a using baculovirus expression system. GST-PKN3 was purified by using glutathione sepharose chromatography and gel filtration chromatography.

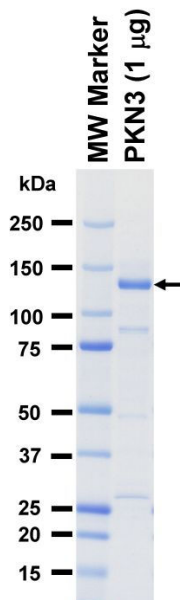
Storage buffer:

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

Storage and Handling:

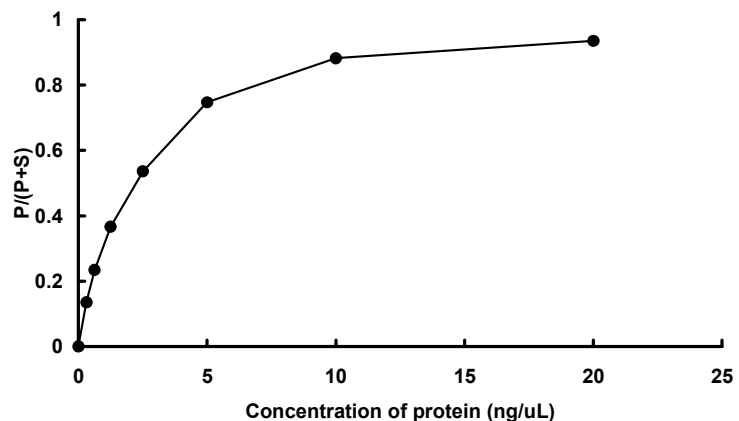
Store at -80C.
 Avoid repeating freeze-thaws.
 Use within 12 months from the date of shipment.

SDS-PAGE



Purity: 71 %
 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 : GS peptide

ATP : 100 µM