

## Product Information

### CaMK1 $\beta$ (PNCK)

Product Number : **02-105**

#### Product description

Full-length human CaMK1 $\beta$  [1-343(end) amino acid of accession number NP\_001353906.1] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system. GST-CaMK1 $\beta$  was purified by using glutathione sepharose chromatography.

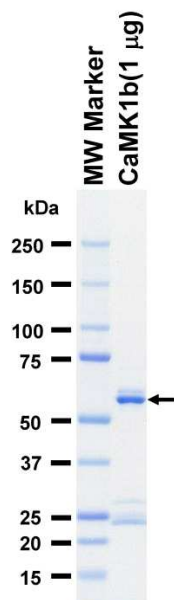
#### Storage buffer:

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.

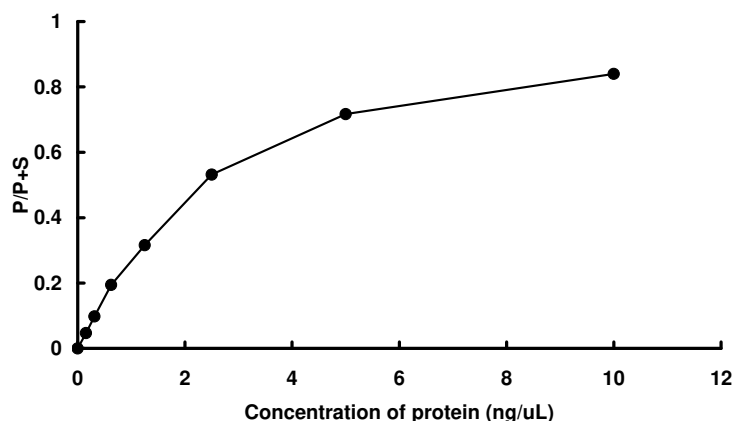
#### SDS-PAGE



Purity: 62 %

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/Calmodulin. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Synapsin peptide

ATP : 100  $\mu$ M