

Product Information

CaMK1 β (PNCK)

 Product Number : **02-105**

Product description

Full-length human CaMK1 β [1-343(end) amino acid of accession number Q6P2M8] was expressed as N-terminal GST-fusion protein (66 kDa) using baculovirus expression system. GST-CaMK1 β was purified by using glutathione sepharose chromatography.

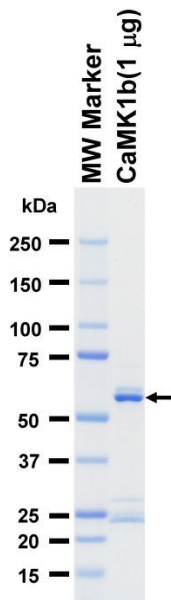
Storage buffer:

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

Storage and Handling:

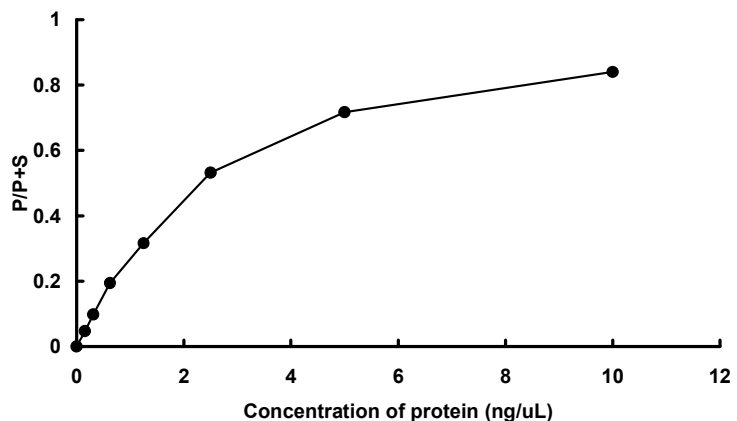
Store at -80C.
 Avoid repeating freeze-thaws.

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. 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 LabChip™ 3000.

Substrate : Synapsin peptide

ATP : 100 μ M