

Product Information

CaMK2 β (CAMK2B)

Product Number : 02-110

Product description

Full-length human CaMK2 β [1-503 amino acids of accession number NP_742078.1] was expressed as N-terminal GST-fusion protein (83 kDa) using baculovirus expression system. GST-CaMK2 β was purified by using glutathione sepharose chromatography.

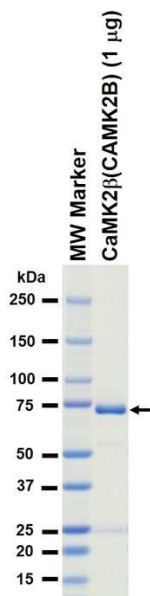
Storage buffer:

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.

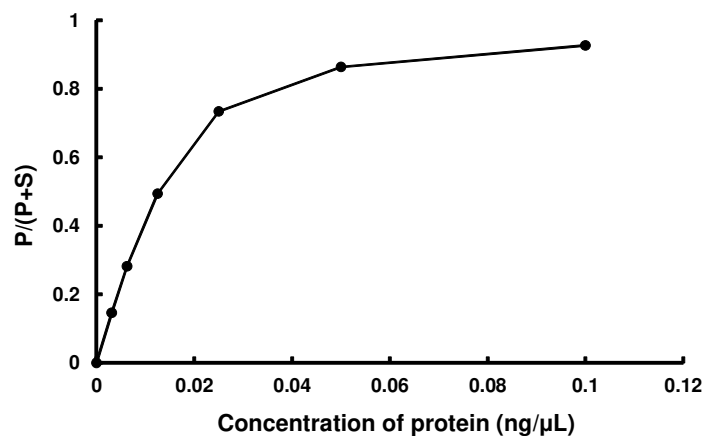
SDS-PAGE



Purity: 87 %

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 : GS peptide

ATP : 100 μ M