

Product Information

CaMK2 γ (CAMK2G)

 Product Number : **02-112**

Product description

Full-length human CaMK2 γ [1-518(end) amino acids of accession number NP_751910.1] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-CaMK2 γ was purified by using glutathione sepharose chromatography.

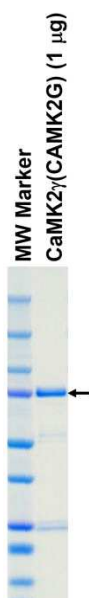
Storage buffer:

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.

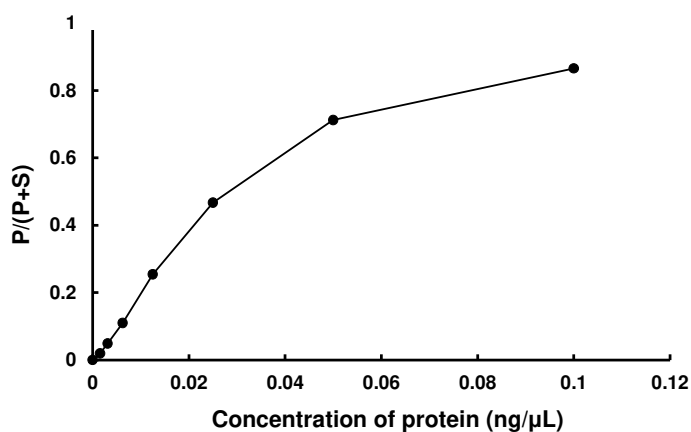
SDS-PAGE



Purity: 76 %

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