

## Product Information

### CaMK1δ(CAMK1D)

 Product Number : **02-106**

#### Product description

Full-length human CaMK1δ[1-357(end) amino acid of accession number NP\_065130.1] was expressed as N-terminal GST-fusion protein (67 kDa) using baculovirus expression system. GST-CaMK1δ was purified by using glutathione sepharose chromatography.

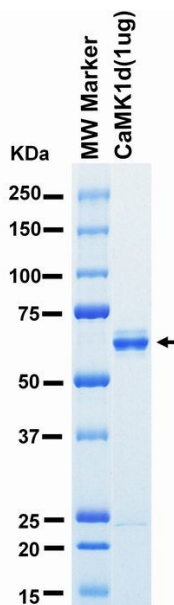
#### Storage buffer:

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

#### Storage and Handling:

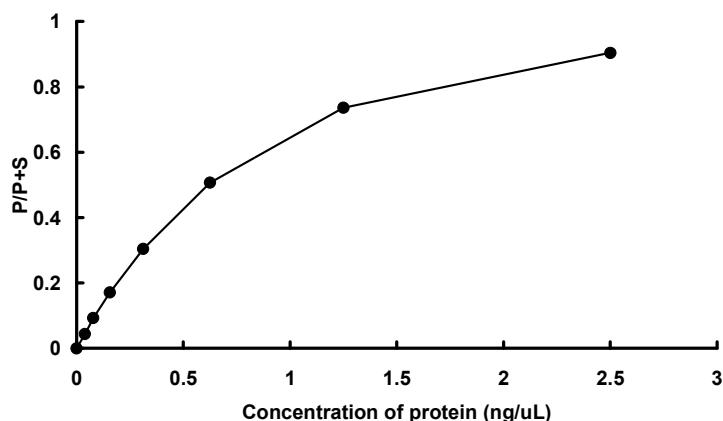
Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



Purity: 89 %  
 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 μM