

## Product Information

### CaMK2δ(CAMK2D)

Product Number : **02-111**

#### Product description

Full-length human CaMK2δ [1-478 amino acids of accession number NP\_742113.1] was expressed as N-terminal GST-fusion protein (81 kDa) using baculovirus expression system. GST-CaMK2δ was purified by using glutathione sepharose chromatography.

#### Storage buffer:

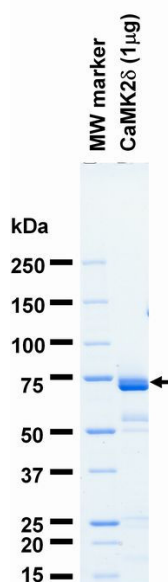
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,  
1 mM DTT, 10% glycerol, pH7.5

#### Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

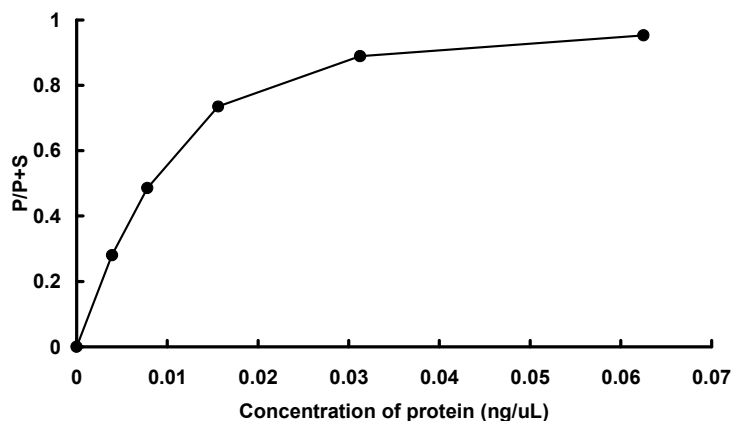
#### SDS-PAGE



Purity: 82 %

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 : GS peptide

ATP : 100 μM