

## Product Information

### CK2 $\alpha$ 2/ $\beta$ (CSNK2A2/ $\beta$ )

 Product Number : **05-185**

#### Product description

Full-length human CK2 $\alpha$ 2 [1-350(end) amino acids of accession number NP\_001887.1] was co-expressed as N-terminal GST-fusion protein (68 kDa) with human His-tagged CK2 $\beta$  [1-215 amino acids of accession number NP\_001311.3] using baculovirus expression system. GST-CK2 $\alpha$ 2 was purified by using glutathione sepharose chromatography.

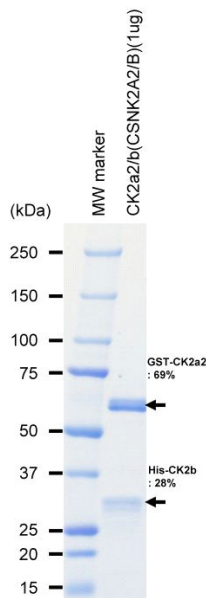
#### Storage buffer:

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.

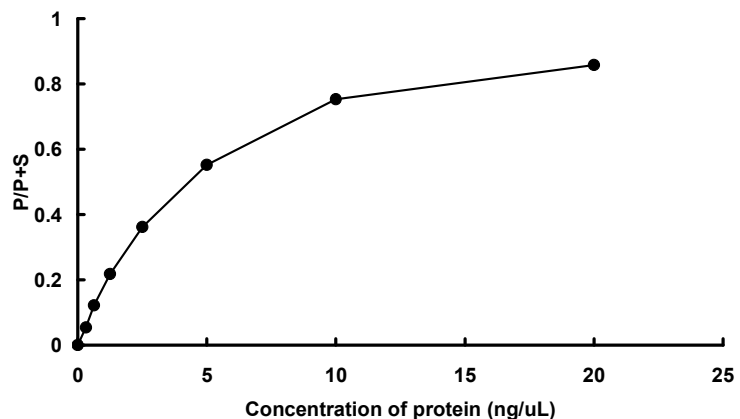
#### SDS-PAGE



Purity: 97 %

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 and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : CK2tide

 ATP : 100  $\mu$ M