

Product Information

CK2 α 1/ β (CSNK2A1/ β)

 Product Number : **05-184**

Product description

Full-length human CK2 α 1 [1-391(end) amino acids of accession number NP_001886.1] was co-expressed as N-terminal GST-fusion protein (72 kDa) with human His-tagged CK2 β [1-215 amino acids of accession number NP_001311.3] using baculovirus expression system. GST-CK2 α 1 was purified by using glutathione sepharose chromatography.

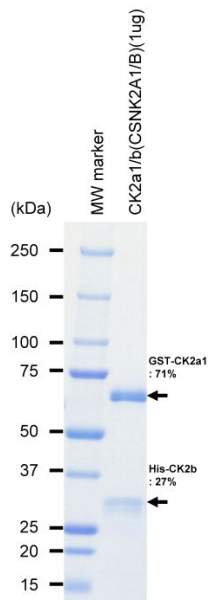
Storage buffer:

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

Storage and Handling:

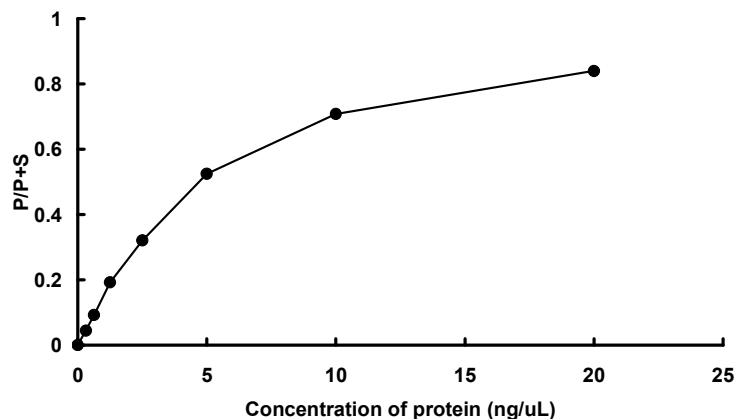
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 98 %
 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 μ M