

Product Information

CDC7/ASK

Product Number : **05-109**

Product description

Full-length human CDC7 [1-574(end) amino acids of accession number NP_003494.1] was co-expressed as N-terminal GST-fusion protein (92 kDa) with Dbf4(ASK) [1-674(end) amino acids of accession number NP_006707.1] using baculovirus expression system. GST-CDC7 was purified by using glutathione sepharose chromatography.

Storage buffer:

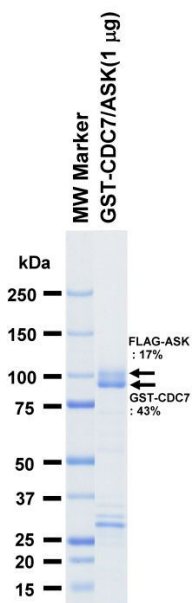
50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

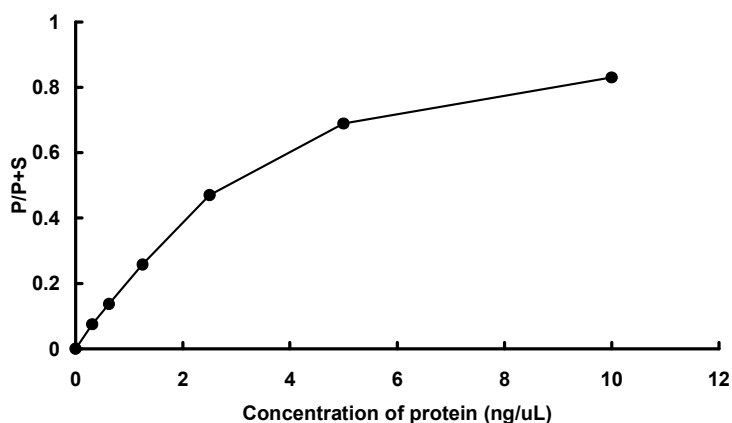
SDS-PAGE



Purity: 60 %

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 : MCM2 peptide

ATP : 100 µM