

Product Information

CDC2L6/CycC

Product Number : **04-112**

Product description

Full-length human CDC2L6 [1-502(end) amino acids of accession number NP_055891.1] was co-expressed as N-terminal GST-fusion protein (85 kDa) with GST-CyclinC [1-283(end) amino acids of accession number NP_005181.2] using baculovirus expression system. GST-CDC2L6 was purified by using glutathione sepharose chromatography.

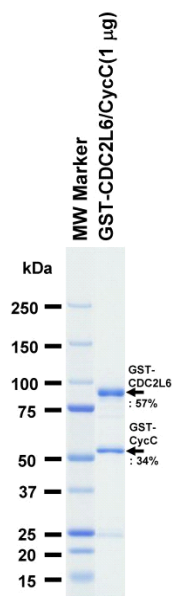
Storage buffer:

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.

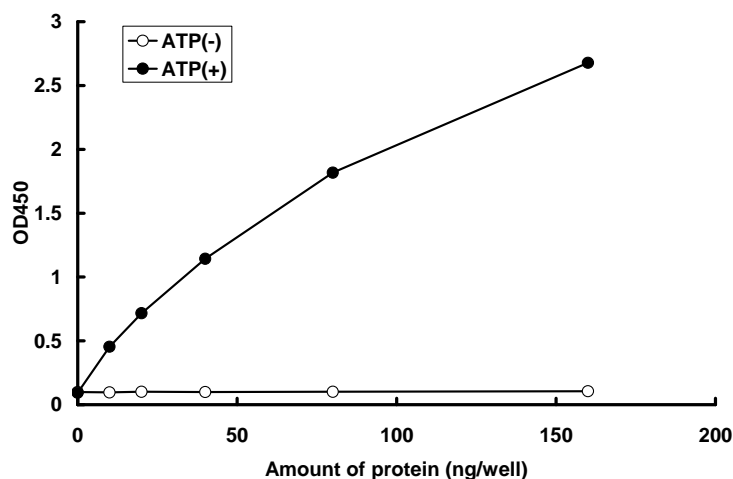
SDS-PAGE



Purity: 92 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was determined by ELISA. The enzyme was incubated with biotinylated peptide in streptavidin-coated ELISA plate. Phosphorylation was detected by anti-phospho antibody and HRP-labeled anti-rabbit IgG(or HRP-labeled anti-mouse IgG).

Substrate : RNA polymerase peptide

ATP : 150 μM