

## Product Information

### CDK13(CHEd)/CycK

Product Number : **04-123**

#### Product description

Human CDK13, catalytic domain [694-1039 amino acids of accession number NP\_003709.3] was co-expressed as N-terminal GST-His-fusion protein (71 kDa) with His-CyclinK(CCNK) [1-580(end) amino acids of accession number NP\_001092872.1] using baculovirus expression system. GST-His-CDK13 was purified by using glutathione sepharose chromatography.

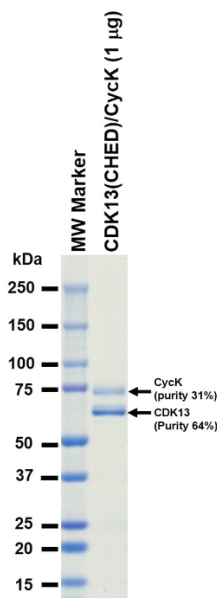
#### Storage buffer:

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.

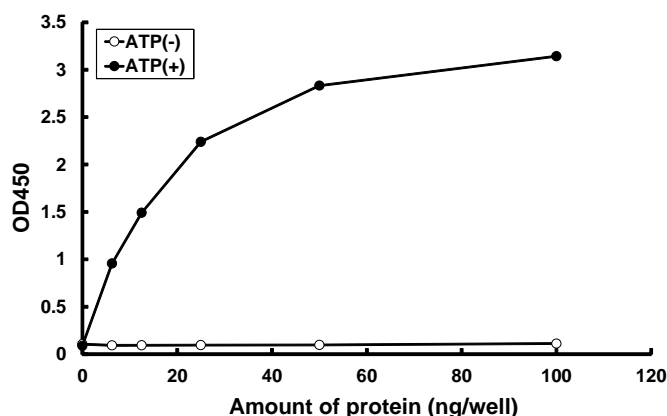
#### SDS-PAGE



Purity: 95 %

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 : 100 µM