

Product Information

BTN-CDK13(CHEd)/CycK

 Product Number : **04-483-20N**

Product description

Human CDK13, catalytic domain [694-1039 amino acids of accession number NP_003709.3] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (45 kDa) with His-CyclinK(CCNK) [1-580(end) amino acids of accession number NP_001092872.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

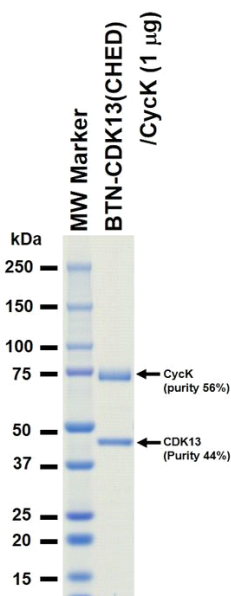
Storage buffer:

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.

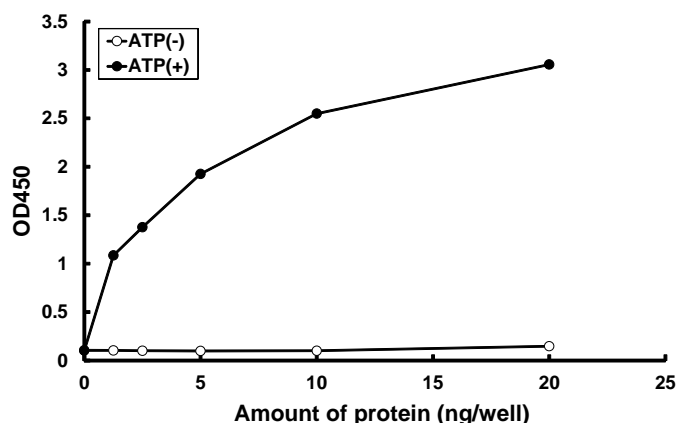
SDS-PAGE



Purity: 100 %

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 μ M