

Product Information

BTN-CDK19(CDC2L6)/CycC

 Product Number : **04-412-20N**

Product description

Full-length human CDK19 [1-502(end) amino acids of accession number NP_055891.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (61 kDa) with GST-CyclinC [1-283(end) amino acids of accession number NP_005181.2] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

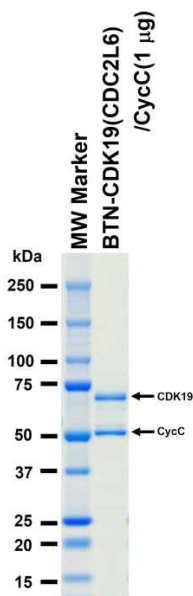
Storage buffer:

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.

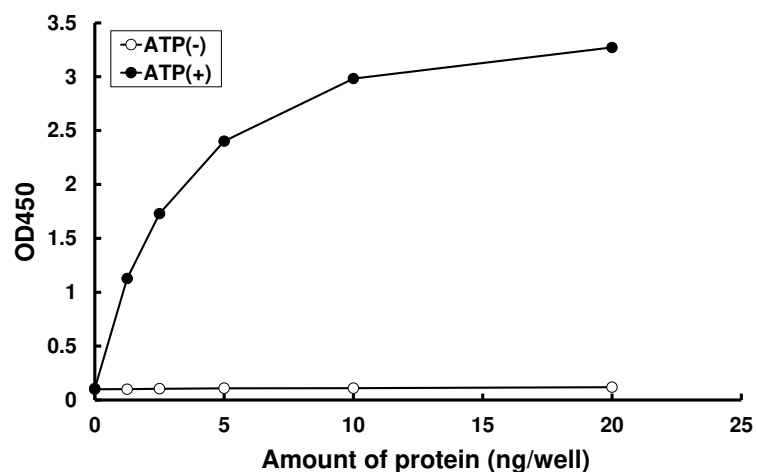
SDS-PAGE



Purity: 96 %

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