

Product Information

BTN-CDK1(CDC2)/CycA2

 Product Number : **04-419-20N**

Product description

Full-length human CDK1 [1-297(end) amino acids of accession number NP_001777.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (38 kDa) with His-CyclinA2 [1-432(end) amino acids of accession number NP_001228.2] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

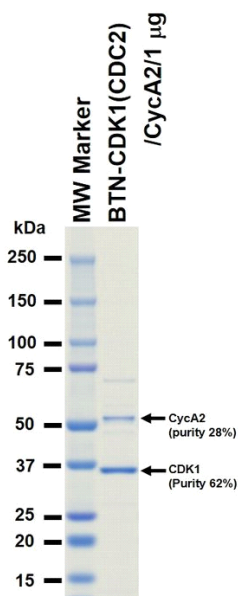
Storage buffer:

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.

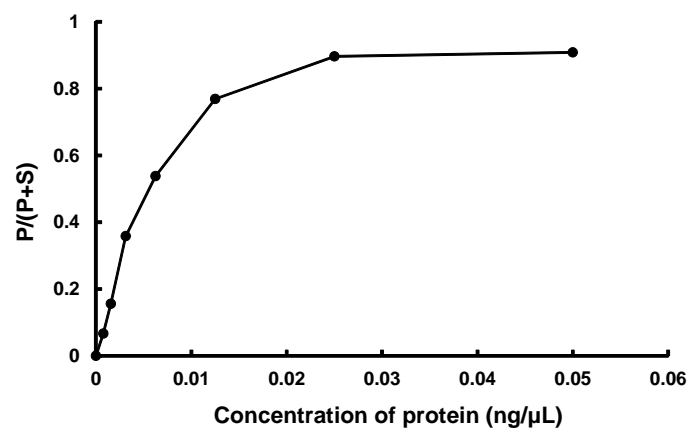
SDS-PAGE



Purity: 90 %

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

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : Modified Histone H1

ATP : 100 µM