

## Product Information

### BTN-CDK9/CycT2

 Product Number : **04-421-20N**

#### Product description

Full-length human CDK9 [1-372(end) amino acids of accession number NP\_001252.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (47 kDa) with His-CyclinT2(CCNT2) [1-730(end) amino acids of accession number NP\_490595.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

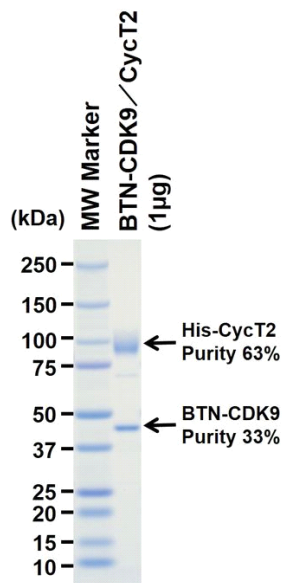
#### Storage buffer:

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.

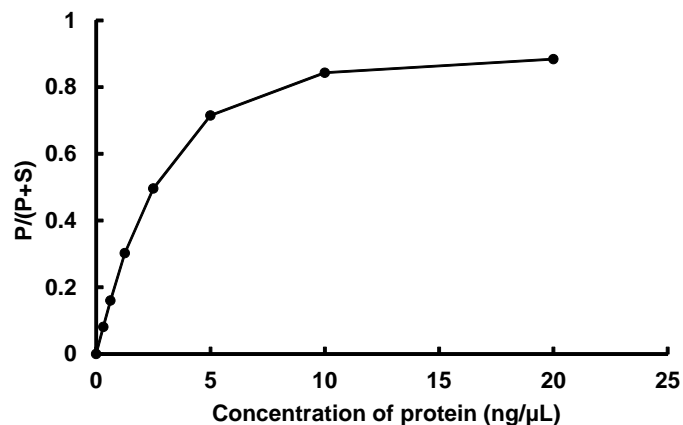
#### SDS-PAGE



Purity: 96 %

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 : CDK9 substrate

ATP : 100 µM