

## Product Information

### BTN-CDK5/p35

Product Number : **04-492-20N**

#### Product description

Full-length human CDK5 [1-292(end) amino acids of accession number NP\_004926.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (37 kDa) with GST-CDK5R1(p35) [1-307(end) amino acids of accession number NP\_003876.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

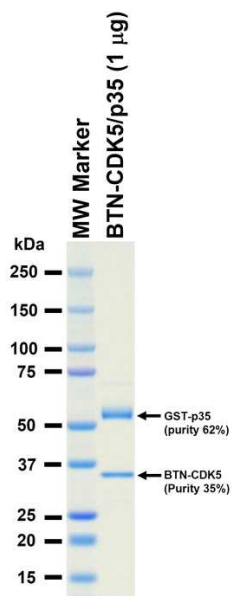
#### Storage buffer:

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.

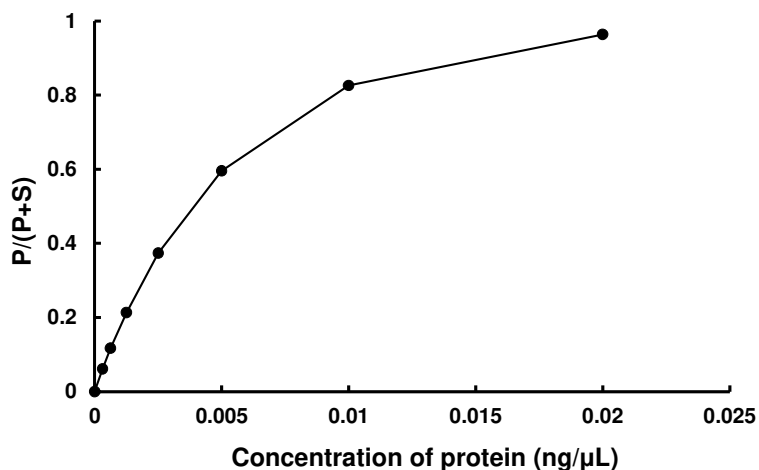
#### SDS-PAGE



Purity: 97 %

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