

Product Information

BTN-CDK4/CycD3

 Product Number : **04-405-20N**

Product description

Full-length human CDK4 [1-303(end) amino acids of accession number NP_000066.1] was co-expressed as N-terminal DYKDDDDK tagged, biotinylated protein (37 kDa) with human GST-CyclinD3 [1-292(end) amino acids and S259A of accession number NP_001751.1] using baculovirus expression system. The protein was purified by using DYKDDDDK tag antibody agarose.

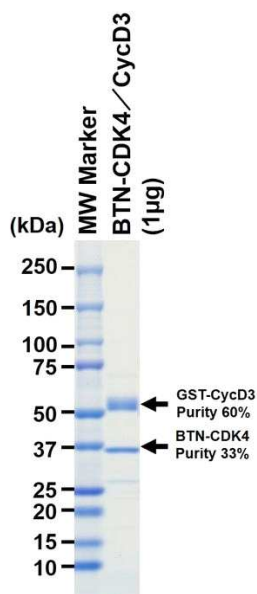
Storage buffer:

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.

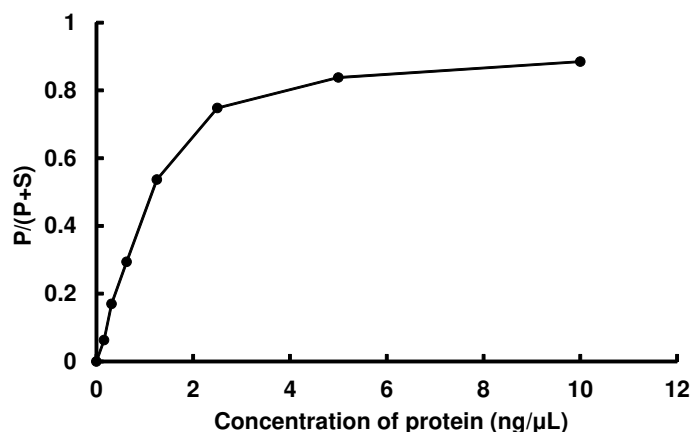
SDS-PAGE



Purity: 93 %

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 : DYRKtide-F

ATP : 200 μM