

## Product Information

### CDK6/CycD3

 Product Number : **04-107**

#### Product description

Full-length human CDK6 [1-326(end) amino acids of accession number NP\_001250.1] was co-expressed as N-terminal GST-fusion protein (64 kDa) with human GST-CyclinD3 [1-292(end) amino acids and S259A of accession number NP\_001751.1] using baculovirus expression system. GST-CDK6/CycD3 was purified by using glutathione sepharose chromatography and activated with His-CDK7/CycH/MAT1. Activated GST-CDK6/CycD3 was purified using glutathione sepharose chromatography.

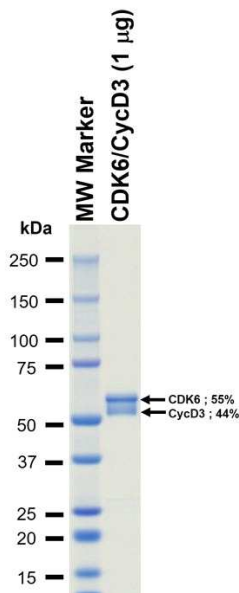
#### Storage buffer:

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.

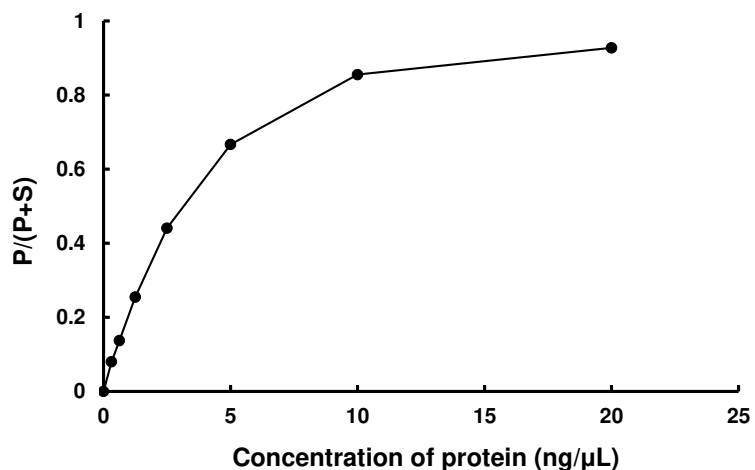
#### SDS-PAGE



Purity: 99 %

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 : 100 µM