

## Product Information

### CDK4/CycD1

 Product Number : **04-120**

#### Product description

Full-length human CDK4 [1-303(end) amino acids of accession number NP\_000066.1] was co-expressed as N-terminal GST-fusion protein (61 kDa) with DYKDDDDK tagged CyclinD1(CCND1) [1-295(end) amino acids of accession number NP\_444284.1] using baculovirus expression system. GST-CDK4/CycD1 was purified by using glutathione sepharose chromatography.

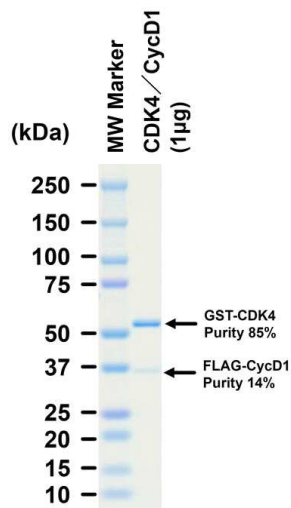
#### Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,  
 1 mM DTT, 10% glycerol, pH 7.5

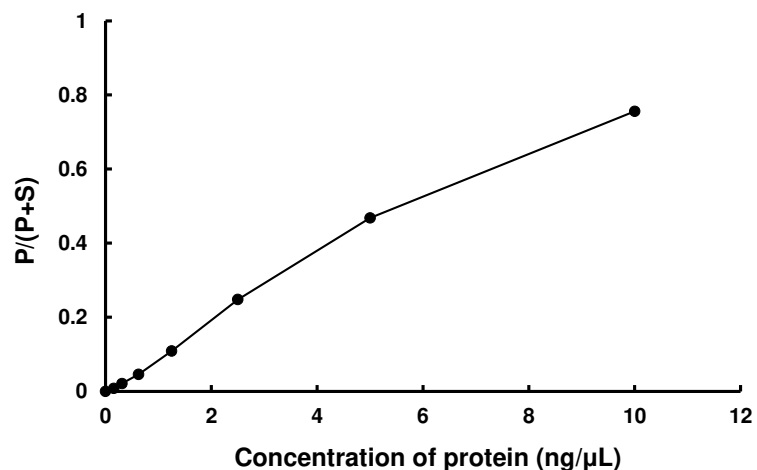
#### Storage and Handling:

Store at -80C.  
 Avoid repeating freeze-thaws.

#### SDS-PAGE



#### Activity data



Purity: 99 %  
 The purity was assessed by SDS-PAGE/CBB staining.

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