

Product Information

CDK4/CycD3

 Product Number : **04-105**

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 human GST-CyclinD3 [1-292(end) amino acids of accession number AAA51927.1] using baculovirus expression system. GST-CDK4/CycD3 was purified by using glutathione sepharose chromatography and activated with His-tagged CDK7. Activated GST-CDK4/CycD3 was purified using glutathione sepharose chromatography.

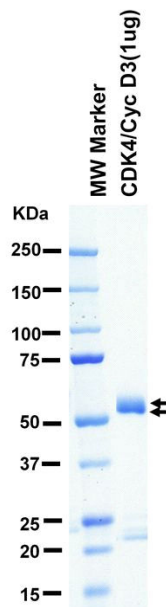
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

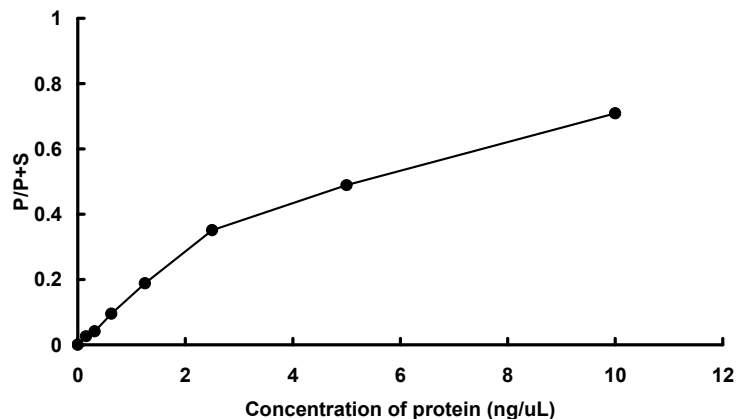
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 84 %
 The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay. The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by LabChip™3000.

Substrate : DYRKtide-F

ATP : 200 μM