

Product Information

CDK9/CycT2

 Product Number : **04-121**

Product description

Full-length human CDK9 [1-372(end) amino acids of accession number NP_001252.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with His-CyclinT2(CCNT2) [1-730(end) amino acids of accession number NP_490595.1] using baculovirus expression system. GST-CDK9 was purified by using glutathione sepharose chromatography.

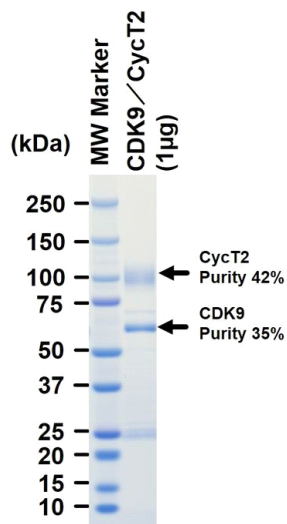
Storage buffer:

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.

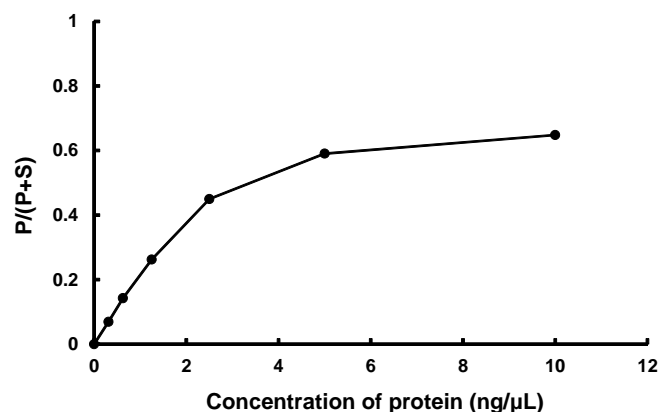
SDS-PAGE



Purity: 77 %

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 : CDK9 substrate

ATP : 100 μM