

Product Information

CDK2/CycE1

 Product Number : **04-165**

Product description

Full-length human CDK2 [1-298(end) amino acids of accession number NP_001789.2] was co-expressed as N-terminal GST-tagged protein (61 kDa) with CyclinE1 [1-410(end) amino acids of accession number NP_001229.1] using baculovirus expression system. GST-CDK2 was purified by using glutathione sepharose chromatography.

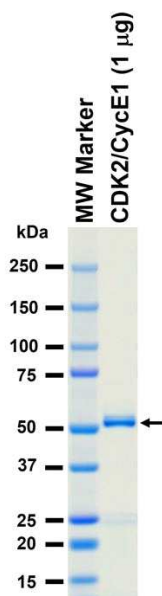
Storage buffer:

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.

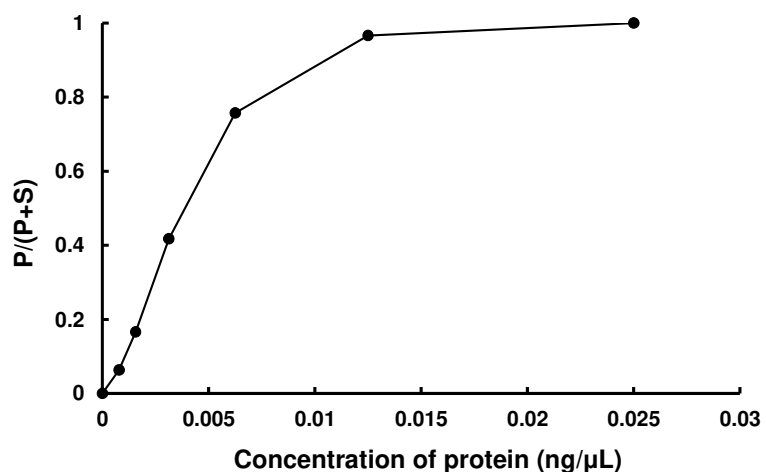
SDS-PAGE



Purity: 90 %

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 : Modified Histone H1

ATP : 100 µM