

Product Information

CDK10(PISSLRE) / CycQ

Product Number : **04-171**

Product description

Full-length human CDK10 [1-360(end) amino acids of accession number NP_443714.3] was co-expressed as N-terminal GST-fusion protein (68 kDa) with His-tagged CyclinQ [1-248(end) amino acids of accession number NP_689487.2] using baculovirus expression system. GST-CDK10/CycQ was purified by using glutathione sepharose chromatography.

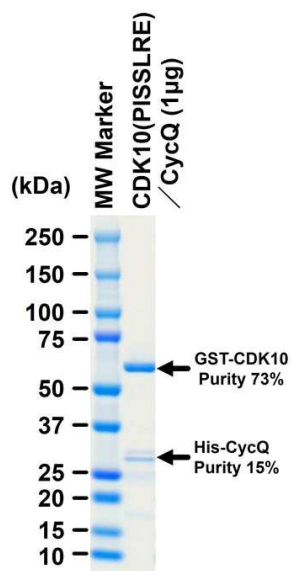
Storage buffer:

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.

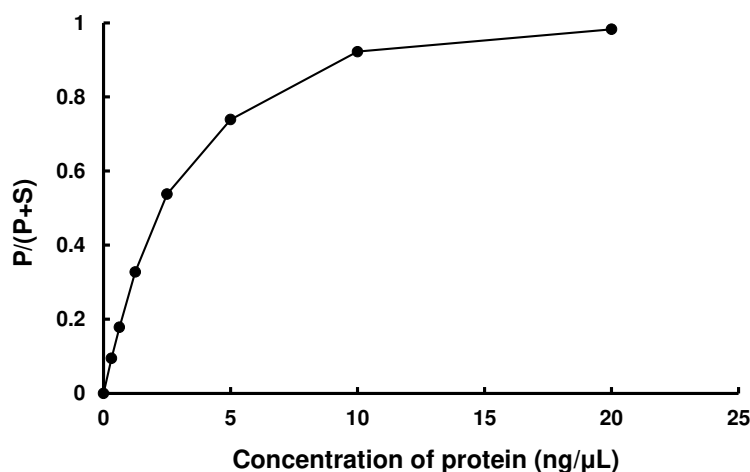
SDS-PAGE



Purity: 88 %

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