

Product Information

PCTAIRE1(CDK16)/CycY

Product Number : **04-116**

Product description

Human PCTAIRE1, catalytic domain [107-496(end) amino acids and S153A of accession number NP_006192.1] was co-expressed as N-terminal GST-fusion protein (71 kDa) with GST-CyclinY(CCN Y) [1-341(end) amino acids of accession number NP_659449.3] using baculovirus expression system. GST-PCTAIRE1 was purified by using glutathione sepharose chromatography.

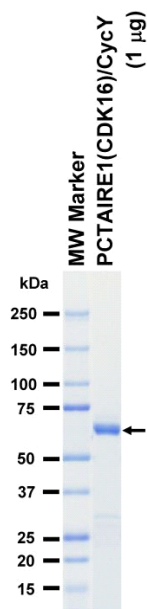
Storage buffer:

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.

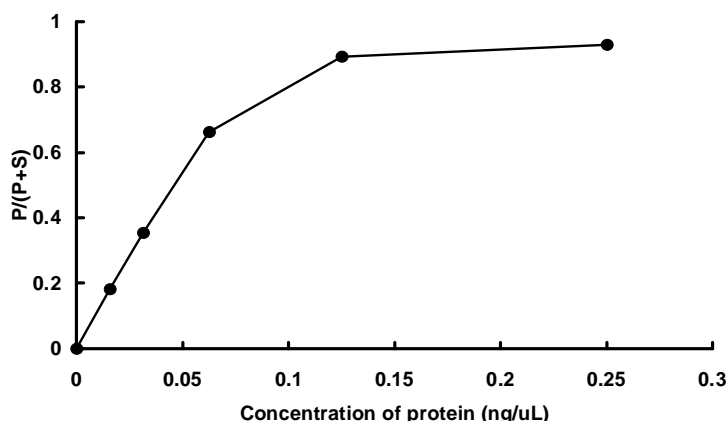
SDS-PAGE



Purity: 91 %

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 : PCTAIRE1 peptide

ATP : 100 µM