

Product Information

PIM1

Product Number : **02-054**

Product description

Full-length human PIM1 [1-313(end) amino acids of accession number NP_002639.1] was expressed as N-terminal His-tagged protein (39 kDa) using baculovirus expression system. His-tagged PIM1 was purified by using Ni-NTA affinity chromatography.

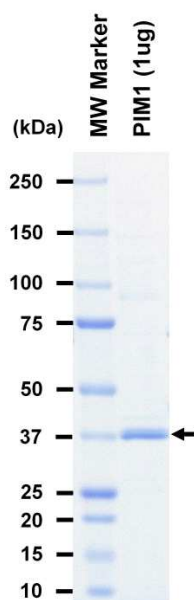
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

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

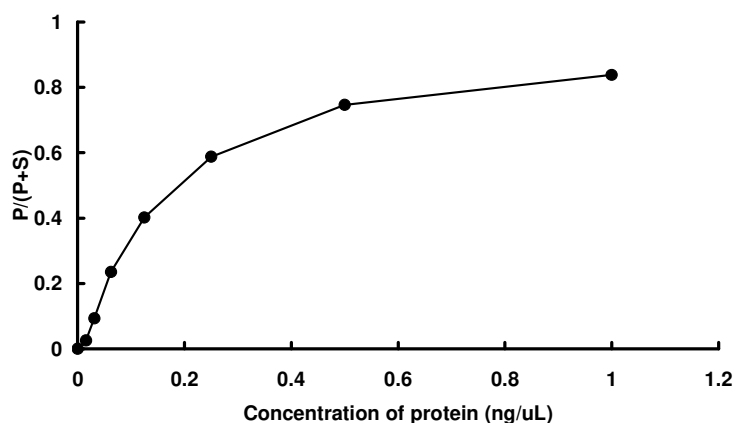
SDS-PAGE



Purity: 83 %

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 : S6K2 peptide

ATP : 100 μ M