

Product Information

TLK1

Product Number : **05-166**

Product description

Full-length human TLK1 [1-766(end) amino acids of accession number NP_036422.3] was expressed as N-terminal GST-fusion protein (113 kDa) using baculovirus expression system. GST-TLK1 was purified by using glutathione sepharose chromatography.

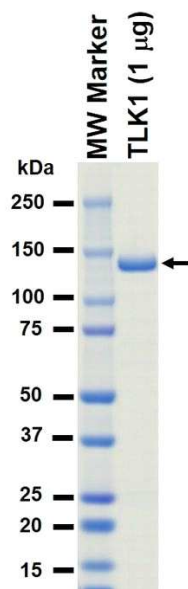
Storage buffer:

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.

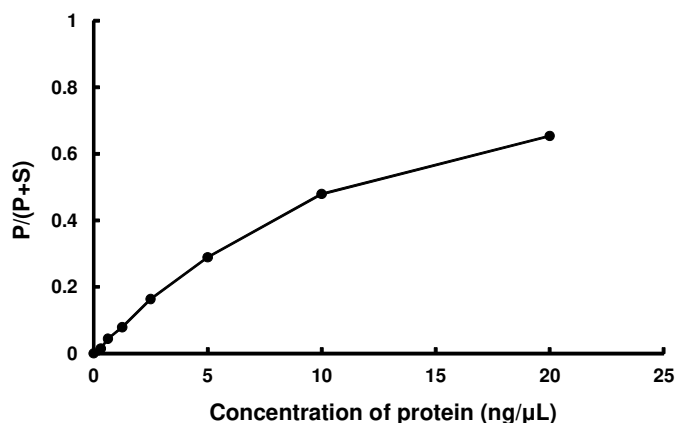
SDS-PAGE



Purity: 97 %

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 : Histone H3 peptide

ATP : 100 µM