

## Product Information

### TNK1

Product Number : **08-104**

#### Product description

Human TNK1, catalytic domain [106-390 amino acids of accession number NP\_003976.2] was expressed as N-terminal GST-fusion protein (58 kDa) using baculovirus expression system. GST-TNK1 was purified by using glutathione sepharose chromatography and gel filtration chromatography.

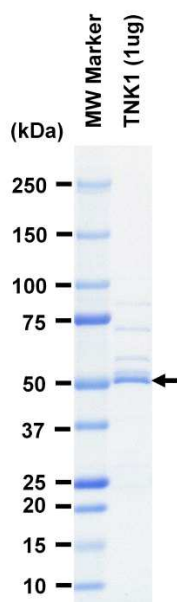
#### Storage buffer:

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.

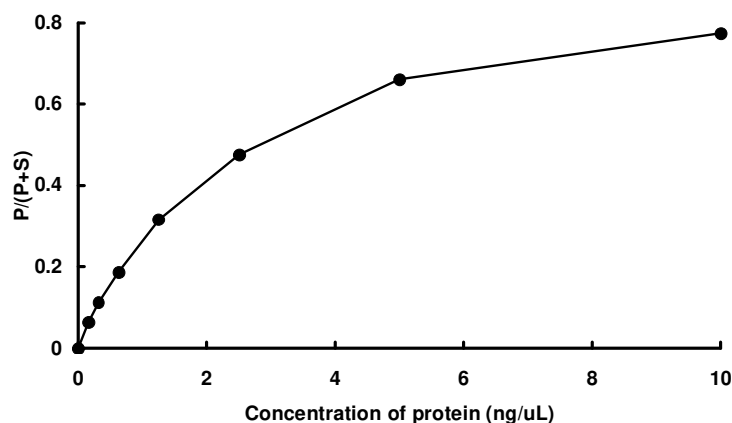
#### SDS-PAGE



Purity: 72 %

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 : CSKtide

ATP : 100  $\mu$ M