

Product Information

TNK1

Product Number : **08-104**

Product description

Human TNK1, catalytic domain [106-390 amino acids of accession number Q13470-2] 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.

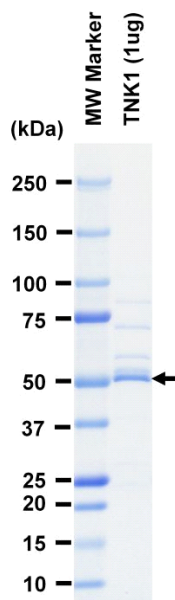
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

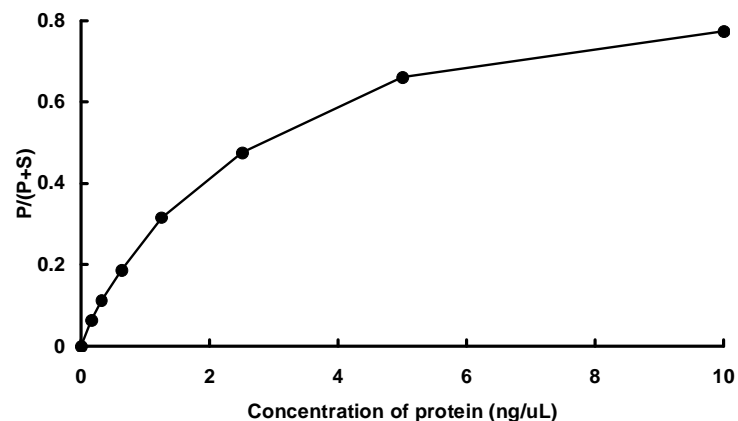
Store at -80C.
 Avoid repeating freeze-thaws.

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. 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 : CSKtide

ATP : 100 μ M