

Product Information

TNIK

Product Number : **07-138**

Product description

Human TNIK, catalytic domain [1-314 amino acids of accession number NP_055843.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system. GST-TNIK was purified by using glutathione sepharose chromatography.

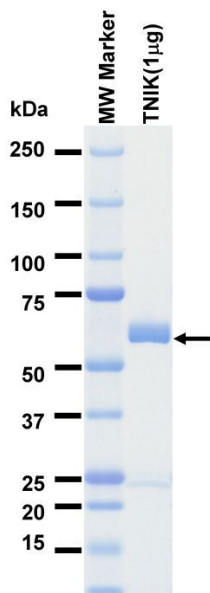
Storage buffer:

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

Storage and Handling:

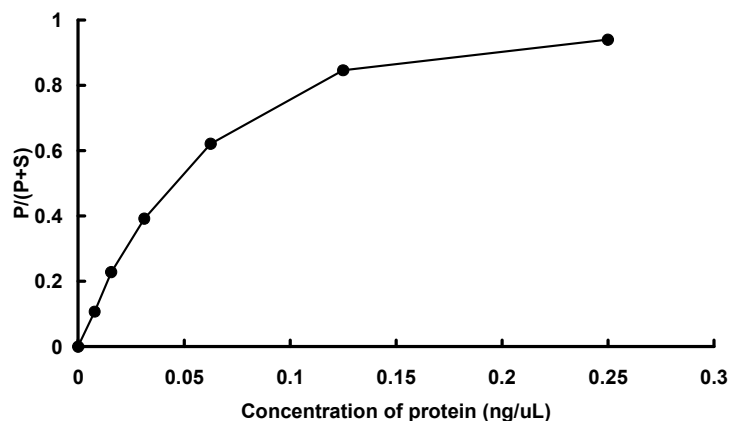
Store at -80C.
Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 90 %
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 : Moesin-derived peptide

ATP : 100 μ M