

Product Information

NEK1

Product Number : **05-123**

Product description

Human NEK1, catalytic domain [1-505 amino acids of accession number NP_036356.1] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-NEK1 was purified by using glutathione sepharose chromatography.

Storage buffer:

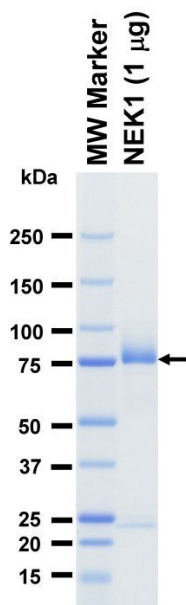
50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.

Avoid repeating freeze-thaws.

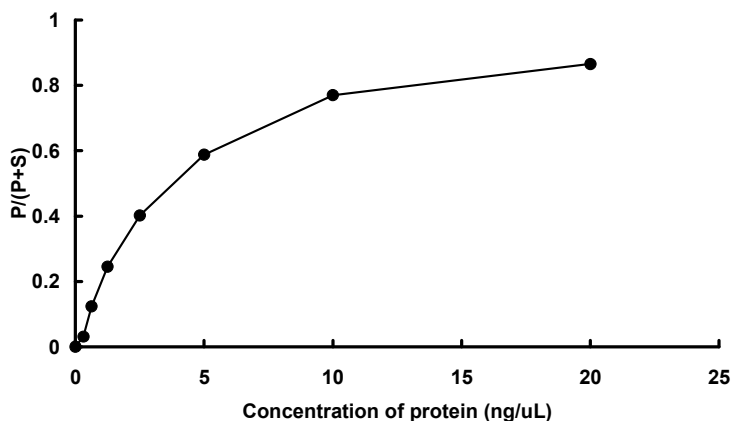
SDS-PAGE



Purity: 87 %

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 : CDK7 peptide

ATP : 100 µM