

Product Information

NEK2

Product Number : **05-226**

Product description

Full-length human NEK2 [1-445(end) amino acids of accession number NP_002488.1] was expressed as N-terminal His-tagged protein (55 kDa) using baculovirus expression system. His-tagged NEK2 was purified by using Ni-NTA affinity chromatography. Purified His-NEK2 was digested by recombinant His-TEV protease, and His-tag free NEK2 (ca. 52 kDa) was collected as flow-through fraction from Ni-NTA affinity chromatography.

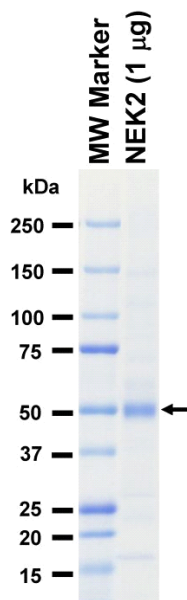
Storage buffer:

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.

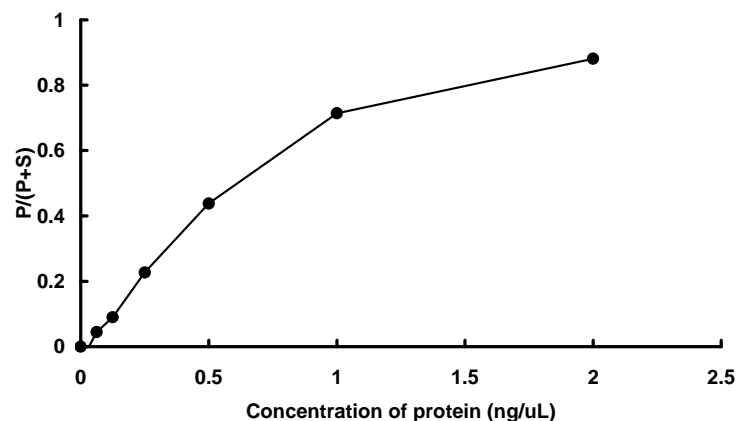
SDS-PAGE



Purity: 73 %

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 : CDK7 peptide

ATP : 100 μ M