

Product Information

MNK2(MKNK2)

Product Number : **02-146**

Product description

Full-length human MNK2 [1-465(end) amino acids and T379D of accession number NP_951009.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system. GST-MNK2 was purified by using glutathione sepharose chromatography.

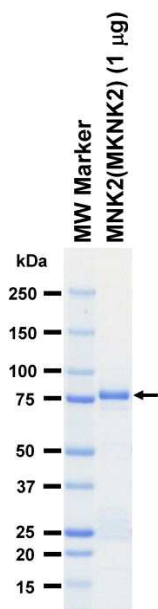
Storage buffer:

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.

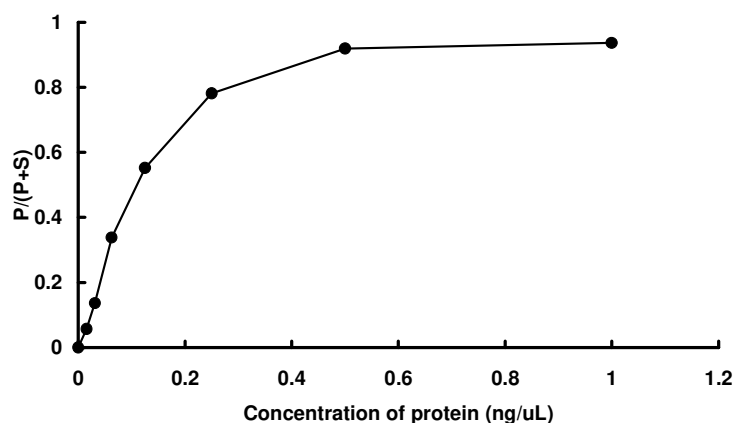
SDS-PAGE



Purity: 78 %

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 : RS peptide

ATP : 100 μM