

Product Information

JNK2(MAPK9)[inactive]

 Product Number : **04-164-10**

Product description

Human JNK2, catalytic domain [1-364 amino acids of accession number NP_002743.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-JNK2 was purified by using glutathione sepharose chromatography.

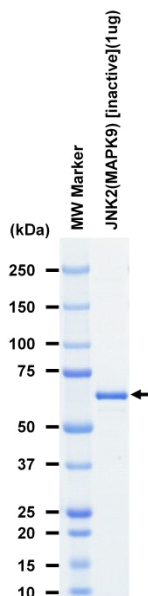
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.1% CHAPS,
 1 mM DTT, 10% glycerol, pH 7.5

Storage and Handling:

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

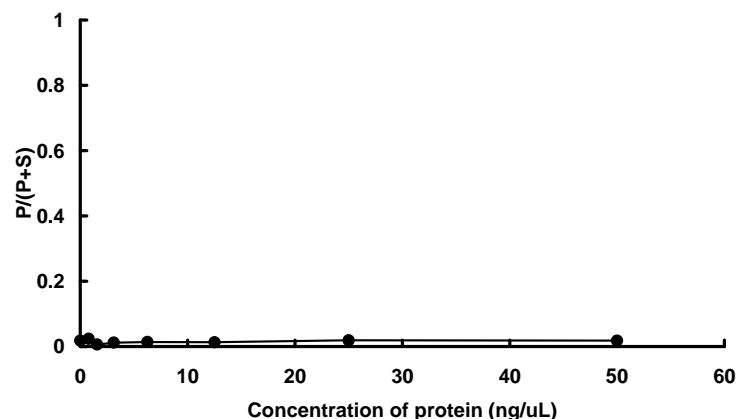
SDS-PAGE



Purity: 96 %

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 : Modified Erktide

 ATP : 1000 μ M