

Product Information

p38 β (MAPK11)

Product Number : **04-153**

Product description

Full-length human p38 β [1-364(end) amino acids of accession number NP_002742.3] was expressed as N-terminal GST-fusion protein (69 kDa) using E. coli expression system. GST-p38 β was purified by using glutathione chromatography.

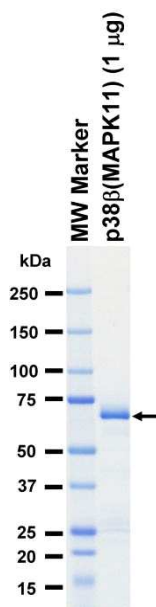
Storage buffer:

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.

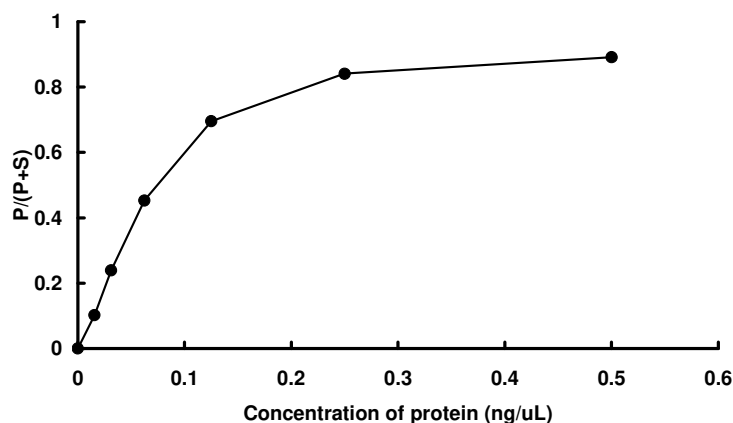
SDS-PAGE



Purity: 88 %

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 : Modified Erkide

ATP : 100 μ M