

Product Information

p38 α (MAPK14)[inactive mutant]

Product Number : **04-152-11**

Product description

Truncated human p38 α [9-352 amino acids and K53M of accession number NP_620581.1] was expressed as N-terminal GST-fusion protein (66 kDa) using E. coli expression system. GST-p38 α was purified by using glutathione sepharose chromatography.

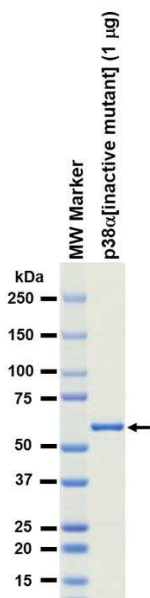
Storage buffer:

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.

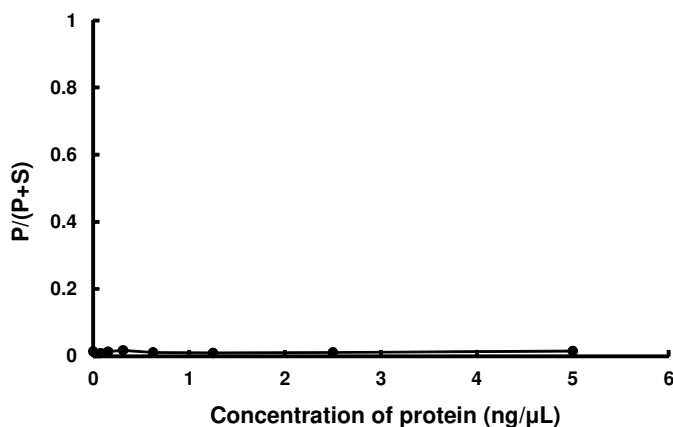
SDS-PAGE



Purity: 99 %

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 Erktide

ATP : 100 μ M