

Product Information

MAPKAPK3

 Product Number : **02-143**

Product description

Full-length human MAPKAPK3 [1-382(end) amino acids of accession number NP_004626.1] was co-expressed as N-terminal GST-fusion protein (70 kDa) with human His-tagged p38 β [1-364(end) amino acids of accession number NP_002742.3] and human His-tagged MAP2K6 [1-334(end) amino acids of accession number NP_002749.2] using baculovirus expression system. GST-MAPKAPK3 was purified by using glutathione sepharose chromatography and Ni-NTA chromatography.

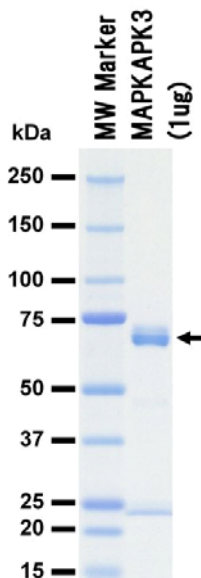
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
 1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

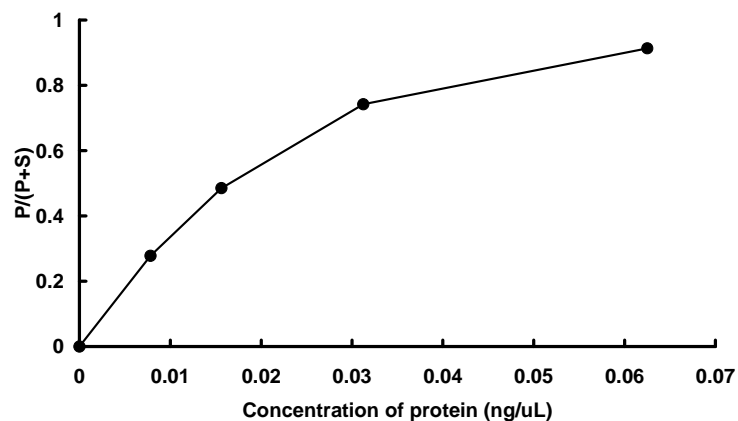
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 76 %
 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 : GS peptide

ATP : 100 μ M