

## 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 $\beta$  [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.

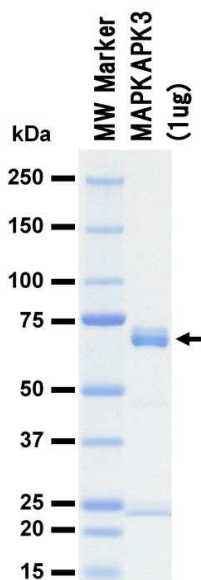
#### Storage buffer:

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.

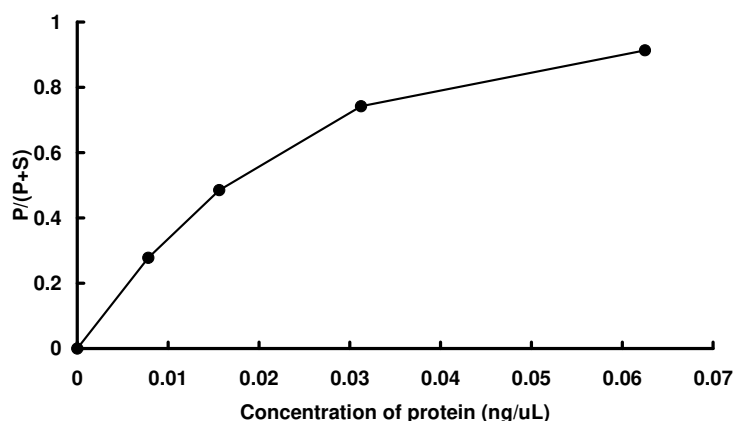
#### SDS-PAGE



Purity: 76 %

The purity was assessed by SDS-PAGE/CBB staining.

#### Activity data



The activity was measured by off-chip mobility shift assay(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 : GS peptide

ATP : 100  $\mu$ M