

Product Information

MAPKAPK2

Product Number : **02-142**

Product description

Full-length human MAPKAPK2 [1-400(end) amino acids of accession number NP_116584.2] was co-expressed as N-terminal GST-fusion protein (73 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-MAPKAPK2 was purified by using glutathione sepharose chromatography and Ni-NTA affinity chromatography.

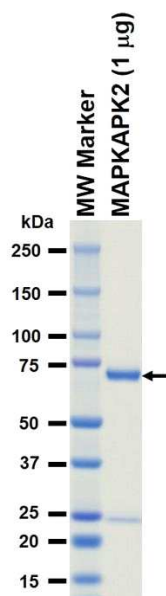
Storage buffer:

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.

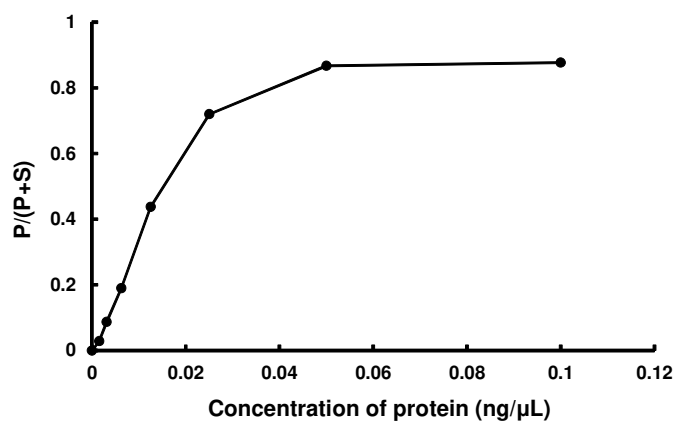
SDS-PAGE



Purity: 82 %

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 μ M