

Product Information

MAPKAPK5

 Product Number : **02-144**

Product description

Full-length human MAPKAPK5 [1-471(end) amino acids of accession number NP_003659.2] was co-expressed as N-terminal GST-fusion protein (81 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] using baculovirus expression system. GST-MAPKAPK5 was purified by using glutathione sepharose chromatography.

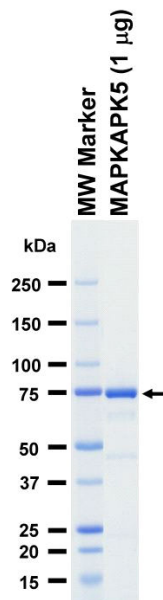
Storage buffer:

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

Storage and Handling:

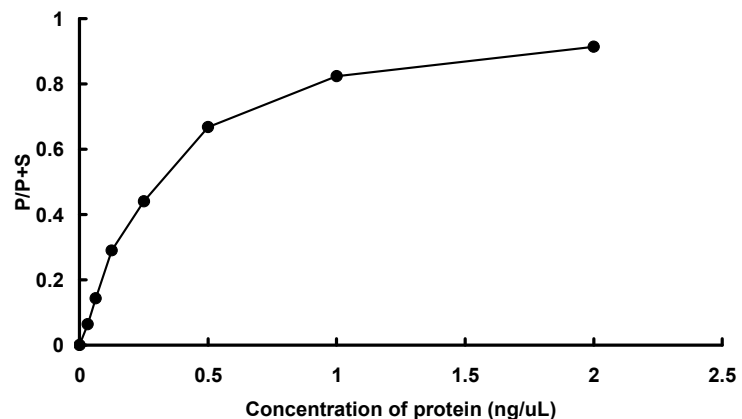
Store at -80C.
 Avoid repeating freeze-thaws.

SDS-PAGE



Purity: 90 %
 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