

## MAPKAPK2, Unactive

Recombinant protein expressed in E. coli cells

Catalog # M40-14G

Lot # K087-1

### Product Description

Recombinant human MAPKAPK2 (46-end) was expressed in E. coli cells using an N-terminal GST tag. The gene accession number is [NM\\_032960](#).

### Gene Aliases

(2 isoforms)

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at  $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

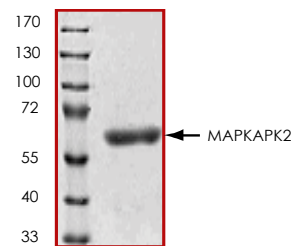
### Scientific Background

MAPKAPK2 (MAPKAP kinase2) is a Ser/Thr protein kinase which is regulated via direct phosphorylation by p38 MAP kinase (1). In conjunction with p38 MAP kinase, MAPKAPK2 is known to be involved in many cellular processes including stress and inflammatory responses, nuclear export, gene expression regulation and cell proliferation. Heat shock protein HSP27 has been shown to be one of the substrates of MAPKAPK2 *in vivo*.

### References

1. Stokoe, D. et al: The substrate specificity and structure of mitogen-activated protein (MAP) kinase-activated protein kinase-2. *Biochem. J.* 296: 843-849, 1993.

### Purity



The purity was determined to be **>90%** by densitometry. Approx. MW **68kDa**.

### Upstream Active Kinases

Active P38a	Cat # M39-10
Active P38 $\beta$	Cat # M36-10
Active P38 $\delta$	Cat # M38-10
Active P38 $\gamma$	Cat # M37-10

## MAPKAPK2, Unactive

Recombinant protein expressed in E. coli cells

Catalog Number	M40-14G
Specific Lot Number	K087-1
Purity	>90%
Concentration	0.2 $\mu\text{g}/\mu\text{l}$
Stability	1yr At $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**