

## MAPKAPK5, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # **M42-14G**

Lot # U1739-6

### Product Description

Recombinant full-length human MAPKAPK5 was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_003668](#).

### Gene Aliases

PRAK

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°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.

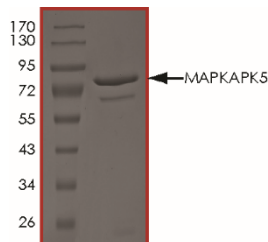
### Scientific Background

MAPKAPK5 is a member of the serine/threonine kinase family that responds to cellular stress and proinflammatory cytokines. MAPKAPK5 is activated through its phosphorylation by MAP kinases including MAPK1/ERK, MAPK14/p38-alpha, and MAPK11/p38-beta (1). MAPKAPK5 is activated in HeLa cells in response to cellular stress and proinflammatory cytokines. MAPKAPK5 activity is regulated by p38-alpha and p38-beta both *in vitro* and *in vivo*, and thr-182 is the regulatory phosphorylation site of MAPKAPK5 (2). *In vitro*, MAPKAPK5 kinase phosphorylates heat shock protein HSP27 at its physiologically relevant sites.

### References

1. New, L.; Jiang, Y. et al: PRAK, a novel protein kinase regulated by the p38 MAP kinase. EMBO J. 17: 3372-3384, 1998.
2. Ni, H.; Wang. et al: MAPKAPK5, a novel mitogen-activated protein kinase (MAPK)-activated protein kinase, is a substrate of the extracellular-regulated kinase (ERK) and p38 kinase. Biochem. Biophys. Res. Commun. 243: 492-496, 1998.

### Purity



The purity of MAPKAPK5 was determined to be **>85%** by densitometry, approx. MW **79kDa**.

## MAPKAPK5, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog # M42-14G

Lot # U1739-6

Purity >85%

Concentration 0.1 µg/µl

Stability 1yr at -70°C from date of shipment

Storage & Shipping Store product at -70°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.**