

Anti-MAPKAPK2/3

Rabbit Polyclonal Antibody

Catalog # M40-63R

Lot # H245-9

Cited Applications

- Western blot (1:1000)

Ideal working dilutions for each application should be empirically determined by the investigator.

Specificity

Recognizes the MAPKAPK2 and MAPKAPK3 protein

Cross Reactivity

- Western blot of human MAPKAPK2 and MAPKAPK3

MAPKAPK2 and MAPKAPK3 from other species may also be detectable

Host

Rabbit

Immunogen

Synthetic peptide corresponding to amino acids 332-336 of human MAPKAPK2

Formulation

PBS pH 7.4, 0.02% sodium azide, 50% glycerol

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles.

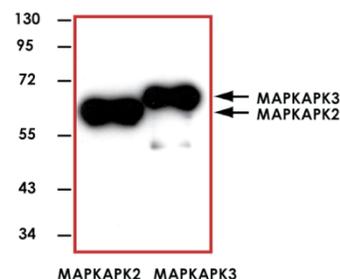
Scientific Background

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. MAPKAPK3 is closely related to MAPKAPK2 sharing 72% nucleotide and 75% amino acid identity (2). MAPKAPK3 is activated by growth inducers and stress stimulation of cells. In vitro studies demonstrated that ERK, p38 MAP kinase and Jun N-terminal kinase were all able to phosphorylate and activate this kinase, which suggested the role of MAPKAPK3 as an integrative element of signaling in both mitogen and stress responses (3).

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.
2. Sithanandam, G. et al: 3pK a new mitogen-activated protein kinase-activated protein kinase located in the small cell lung cancer tumor suppressor gene region. *Molec. Cell. Biol.* 16: 868-876, 1996.
3. Ludwig, S. et al: 3pK, a novel mitogen-activated protein (MAP) kinase-activated protein kinase, is targeted by three MAP kinase pathways. *Molec. Cell. Biol.* 16: 6687-6697, 1996.

Sample Data



Representative western blot with Anti-MAPKAPK2/3 (1:1000) using 20 ng of GST tagged human recombinant MAPKAPK2 and MAPKAPK3.

Anti-MAPKAPK2/3

Rabbit Polyclonal Antibody

Catalog Number

M40-63R

Specific Lot Number

H245-9

Purification

Affinity chromatography

Concentration

1.0 µg/µL

Stability

1 yr At -20°C from date of shipment

Storage & Shipping

Store product at -20°C. For optimal storage, aliquot antibody into smaller quantities after centrifugation and store at recommended temperature. For optimal performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on ice packs.

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

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