

Anti-phospho-MEK1/2 (Ser218 Ser222)

Rabbit Polyclonal Antibody

Catalog # M02-652R

Lot # J1274-20

Cited Applications

WB

Suggested Dilutions:
WB 1:1,000

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

Specificity

Recognizes the MEK1/2 protein phosphorylated at serine 218 and serine 222

Cross Reactivity

Human, Mouse, Rat, Bovine, Canine, Chicken, non-Human Primates, Xenopus and Zebrafish

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Ser218/222 conjugated to KLH

Formulation

100 µl in 10 mM HEPES (pH 7.5), 150 mM NaCl, 100 µg per ml BSA and 50% glycerol.

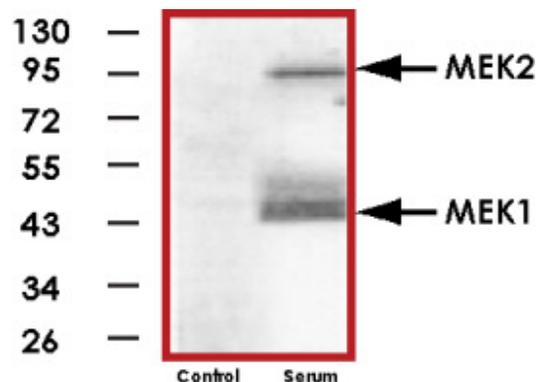
Scientific Background

MEK1 (MAP Kinase Kinase, also known as MKK) is a member of the dual specificity protein kinase family. MEK1 lies upstream of MAPK/ERK and stimulates the enzymatic activity of MAPK/ERK upon a wide variety of extra- and intracellular signals. As an essential component of the MAPK/ERK signal transduction pathway, MEK1 is involved in many cellular processes such as proliferation, differentiation, transcriptional regulation and development (1). This pathway plays a key role in synaptic plasticity in the brain (2). Constitutive activation of MEK1 results in cellular transformation. Thus, MEK1 represents a likely target for pharmacological intervention in proliferative diseases such as cancer (3). MEK1 can be phosphorylated by MAP kinase at Thr292 and Thr386.

References

1. Seger, R. et al: The MAPK signaling cascade. FASEB J. 1995 Jun;9(9):726-35.
2. Adams, J P. et al: Molecular psychology: Roles for the ERK MAP kinase cascade in memory. Annu Rev Pharmacol Toxicol 2002 42:135-163.
3. Sebolt-Leopold, J S. et al: Blockade of the MAP Kinase Pathway Suppresses growth of colon tumors in vivo. Nature Med. 1999 5: 810-816.

Sample Data



Western Blot of NIH 3T3 cell lysates showing specific immunolabeling of the ~45kDa MEK 1/2 protein phosphorylated at Ser218 and Ser222. The cells were either serum starved (Control) or incubated in the presence of serum (Serum). Immunolabeling of an additional band at ~95kDa was also observed.

Anti-phospho-MEK1/2 (Ser218 Ser222)

Rabbit Polyclonal Antibody

Catalog Number

M02-652R

Specific Lot Number

J1274-20

Purification
Stability
Storage & Shipping

Affinity chromatography
1yr at -20°C from date of shipment
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.