

Anti-phospho-MEK1 (Thr386)

Rabbit Polyclonal Antibody

Catalog # M02-65CR

Lot # J1274-22

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 protein phosphorylated at threonine 386

Cross Reactivity

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

Host/Isotype/Clone#

Rabbit, IgG

Immunogen

Synthetic phospho-peptide corresponding to amino acid residues surrounding Thr386 conjugated to KLH

Formulation

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

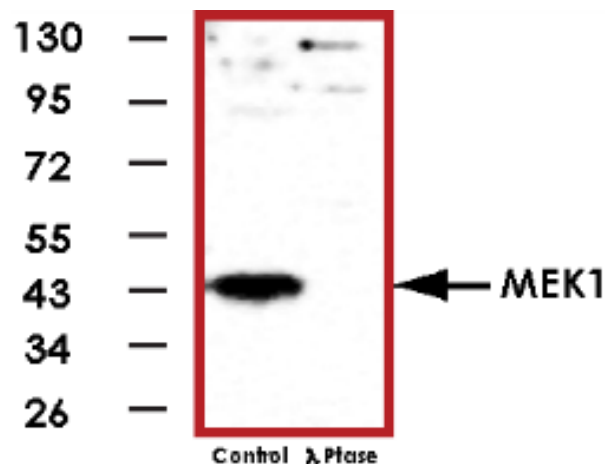
Scientific Background

MEK1 is a member of the dual specificity protein kinase family and is also referred to as MAP Kinase Kinase or MKK. 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 including proliferation, differentiation, transcriptional regulation and development (1). MEK1 plays a key role in synaptic plasticity in the brain (2). Constitutive activation of MEK1 results in cellular transformation. Thus, MEK 1 represents a likely target for pharmacological intervention in proliferative diseases such as cancer (3). MAP kinase can phosphorylate MEK1 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 PharmacolToxicol 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 human T47D cells showing specific immunolabeling of the ~45kDa MEK1 (Control). The phosphospecificity of this labeling is shown in the second lane (lambda phosphatase: lambdaPtase). The blot is identical to the control except that it was incubated in lambda Ptase (1200 units for 30 min) before being exposed to the Anti-phospho-MEK1 (Thr386) antibody. The immunolabeling of MEK1 is completely eliminated by lambda Ptase.

Anti-phospho-MEK1 (Thr386)

Rabbit Polyclonal Antibody

Catalog Number

M02-65CR

Specific Lot Number

J1274-22

Purification

Affinity chromatography

Stability

1yr 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.