

PRODUCT INFORMATION



MEK1 (Phospho-Thr³⁸⁶) Polyclonal Antibody

Item No. 10009517

Overview and Properties

Contents:	This vial contains 100 µl affinity-purified antibody in 10 mM HEPES, pH 7.5, containing 150 mM NaCl, 100 µg/ml BSA, and 50% glycerol.
Synonyms:	MAP Kinase Kinase 1, MAPKK1, Mitogen-activated Protein Kinase Kinase 1
Immunogen:	Phosphopeptide corresponding to amino acid residues surrounding Phospho-Thr ³⁸⁶ of human MEK1
Cross Reactivity:	(+) Rat MEK1; expected to react with bovine, canine, chicken, mouse, non-human primates, and <i>Xenopus</i> MEK1
Storage:	-20°C (as supplied)
Stability:	≥1 year
Host:	Rabbit
Applications:	Western blot (WB); the recommended starting dilution for WB is 1:1,000.

Description

MAP kinase kinase 1 (MEK1) is an integral component of the MAP kinase cascade that regulates cell growth and differentiation.^{1,2} This pathway also plays a key role in synaptic plasticity in the brain.³ Activated MEK1 acts as a dual specificity kinase phosphorylating both a threonine and a tyrosine residue on MAP kinase.⁴⁻⁶ Conversely, there also appears to be a feedback phosphorylation of MEK1 by MAP kinase. The sites on MEK1 that are phosphorylated by MAP kinase are Thr²⁹² and Thr³⁸⁶.⁷

References

1. Ahn, N.G. The MAP kinase cascade. Discovery of a new signal transduction pathway. *Mol. Cell Biochem.* **127-128**, 201-209 (1993).
2. Chong, H., Vikis, H.G., Guan, K.L. Mechanisms of regulating the Raf kinase family. *Cell. Signal.* **15(5)**, 463-469, (2003).
3. Adams, J.P., Sweatt, J.D. Molecular psychology: Roles for the ERK MAP kinase cascade in memory. *Annu. Rev. Pharmacol. Toxicol.* **42**, 135-163 (2002).
4. Kyriakis, J.M., Brautigan, D.L., Ingebritsen, T.S., *et al.* pp54 Microtubule-associated protein-2 kinase requires both tyrosine and serine/threonine phosphorylation for activity. *J. Biol. Chem.* **266(16)**, 10043-10046 (1991).
5. Seger, R., Ahn, N.G., Boulton, T.G., *et al.* Microtubule-associated protein 2 kinases, ERK1 and ERK2, undergo autophosphorylation on both tyrosine and threonine residues: Implications for their mechanism of activation. *Proc. Natl. Acad. Sci. U.S.A.* **88(14)**, 6142-6146 (1991).
6. Crews, C.M., Alessandrini, A. and Erikson, R.L. The primary structure of MEK a protein kinase that phosphorylates the ERK gene product (extracellular signal-regulated kinases). *Science* **258(5081)**, 478-480 (1992).
7. Mansour, S.J., Resing, K.A., Candi, J.M., *et al.* Mitogen-activated protein (MAP) kinase phosphorylation of MAP kinase kinase: Determination of phosphorylation sites by mass spectrometry and site-directed mutagenesis. *J. Biochem.* **116(2)**, 304-314 (1994).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 11/01/2017

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM