

Product Information



MEK 1/2 (Phospho-Ser^{218,222}) Polyclonal Antibody

Catalog No. 10009178

Supplied as:	Affinity-purified IgG
Host:	Rabbit
Antigen:	Phosphopeptide corresponding to amino acid residues surrounding phospho-Ser ^{218,222} of human MEK 1/2
Cross Reactivity:	(+) NIH 3T3 cells
Stability:	≥1 year at -20°C
Application:	The recommended starting dilution for western blot is 1:1,000. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

MEK 1 (MAP kinase kinase, also known as MKK) 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 MEK 1 acts as a dual specificity kinase phosphorylating both a threonine and a tyrosine residue on MAP kinase.⁴⁻⁶

References

1. Ahn, N.G. Discovery of a new signal transduction pathway. *Mol. Cell. Biochem.* **127-128**, 201-209 (1993).
2. Chong, H., Vikis, H.G., and Guan, K.L. Mechanisms of regulating the Raf kinase family. *Cell. Signal.* **15**, 463-469 (2003).
3. Adams, J.P. and Sweatt, J.D. Molecular psychology: Roles for the ERK MAP kinase cascade in memory. *Annu. Rev. Pharmacol. Toxicol.* **42**, 135-163 (2002).
4. Crews, C.M., Alessandrini, A., and Erikson, R.L. The primary structure of MEK, a protein kinase that phosphorylates the ERK gene product. *Science* **258**, 478-480 (1992).
5. 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).
6. 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. USA* **88**, 6142-6146 (1991).

Related Products

5-Lipoxygenase (Phospho-Ser⁵²³) Polyclonal Antibody - Cat. No. 10007820 • p38 MAPK (Phospho-Thr¹⁸⁰/Tyr¹⁸²) Polyclonal Antibody - Cat. No. 10009177 • ERK/MAPK (Phospho-Thr²⁰²/Tyr²⁰⁴) Polyclonal Antibody - Cat. No. 10009179 • β-Catenin (Phospho-Ser^{33,37}) Polyclonal Antibody - Cat. No. 10009180 • CREB (Phospho-Ser¹³³) Polyclonal Antibody - Cat. No. 10009181

Cayman Chemical

Mailing address
1180 E. Ellsworth Road
Ann Arbor, MI
48108 USA

Phone
(800) 364-9897
(734) 971-3335

Fax
(734) 971-3640

E-Mail
custserv@caymanchem.com

Web
www.caymanchem.com

WARNING: THIS PRODUCT IS FOR LABORATORY RESEARCH ONLY: NOT FOR ADMINISTRATION TO HUMANS. NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman warrants only to the original customer that the material will meet our specifications at the time of delivery.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have any obligation or liability, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's exclusive remedy and Cayman's sole liability hereunder shall be limited to a refund of the purchase price, or at Cayman's option, the replacement, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our Warranty and Limitation of Remedy located on our website and in our catalog.

Copyright Cayman Chemical Company, 01/05/2010