

Product Information



ERK/MAPK (Phospho-Thr²⁰²/Tyr²⁰⁴) Polyclonal Antibody

Catalog No. 10009179

Supplied as:	Affinity-purified IgG
Host:	Rabbit
Antigen:	Phosphopeptide corresponding to amino acid residues surrounding the phospho-Thr ²⁰² and Tyr ²⁰⁴ of rat ERK/MAPK. The amino acid residues surrounding the Thr/Tyr phosphorylation sites of ERK/MAPK of Rat, primate(s), mouse, dog, cow, frog, and zebrafish (for example, G F L T E Y V A T) are identical.
Cross Reactivity:	(+) Human and rat ERK/MAPK, others not tested
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.

Extracellular-signal regulated kinase/mitogen-activated protein kinase (ERK/MAPK) is an integral component of cellular signaling during mitogenesis and differentiation of mitotic cells and also is thought to play a key role in learning and memory.¹⁻⁴ The activity of this kinase is regulated by dual phosphorylation at Thr²⁰² and Tyr²⁰⁴.³

References

1. Johnson, G.L. and Lapadat, R. Mitogen-activated protein kinase pathways mediated by ERK, JNK, and p38 protein kinases. *Science* **298**, 1911-1912 (2002).
2. 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).
3. Ahn, N.G. Discovery of a new signal transduction pathway. *Mol. Cell. Biochem.* **127-128**, 201-209 (1993).
4. Tanoue, T.J. and Nishida, E. Molecular recognitions in the MAP kinase cascades. *Cell. Signal.* **15**, 455-462 (2003).

Related Products

5-Lipoxygenase (Phospho-Ser⁵²³) Polyclonal Antibody - Cat. No. 10007820 • p38 MAPK (Phospho-Thr¹⁸⁰/Tyr¹⁸²) Polyclonal Antibody - Cat. No. 10009177 • MEK 1/2 (Phospho-Ser^{218,222}) Polyclonal Antibody - Cat. No. 10009178 • β-Catenin (Phospho-Ser^{33,37}) Polyclonal Antibody - Cat. No. 10009180 • CREB (Phospho-Ser¹³³) Polyclonal Antibody - Cat. No. 10009181

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, 03/23/2010

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