

Anti-ERK1

Mouse Monoclonal Antibody

Catalog # M29-60M

Lot # V058-4

Cited Applications

- Immunoprecipitation (0.1 µg)
- Western blot (1:1000 dilution)

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

Specificity

Recognizes only the ERK1 protein

Cross Reactivity

- Immunoprecipitation of ERK1 from human cells
- Western blot ERK1 from human and mouse cells

ERK1 from other species may also be detectable

Host/Isotype/Clone#

Mouse, IgG1, E32

Immunogen

Recombinant full length human ERK1 protein

Formulation

10mM PBS pH 7.2

Stability

Store at 4°C (add 0.1% NaN₃) for several months, and at -20°C for longer periods. 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.

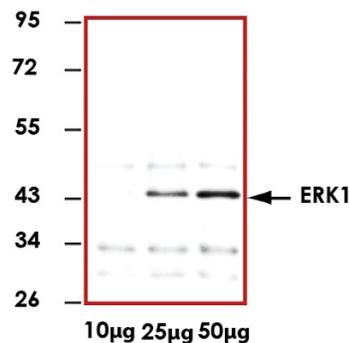
Scientific Background

ERK1 is a protein serine/threonine kinase that is a member of the extracellular signal-regulated kinases (ERKs) which are activated in response to numerous growth factors and cytokines (1). Activation of ERK1 requires both tyrosine and threonine phosphorylation that is mediated by MEK. ERK1 is ubiquitously distributed in tissues with the highest expression in heart, brain and spinal cord. Activated ERK1 translocates into the nucleus where it phosphorylates various transcription factors (e.g., Elk-1, c-Myc, c-Jun, c-Fos, and C/EBP beta).

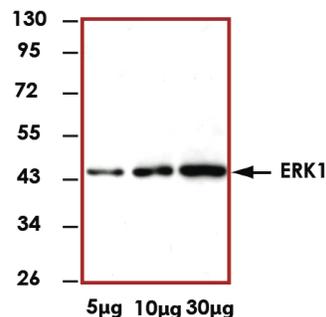
References

1. Boulton, TG. et al: Purification and properties of extracellular signal-regulated kinase 1, an insulin-stimulated microtubule-associated protein 2 kinase. *Biochemistry*. 1991 Jan 8;30(1):278-86.

Sample Data



Representative immunoprecipitation with 0.1 µg of Anti-ERK1 using 10 µg, 25 µg, and 50 µg of HeLa cell lysate.



Representative western blot with Anti-ERK1 (1:1000) using 5 µg, 10 µg, and 30 µg HEK-293 cell lysate.

Anti-ERK1

Mouse Monoclonal Antibody

Catalog Number	M29-60M
Specific Lot Number	V058-4
Purification	Affinity purified using Protein G
Concentration	0.1 µg/µL
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

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.