

# MEK1 polyclonal antibody

MEK is a dual specificity kinase capable of phosphorylating both tyrosine and threonine residues. The MEK family, which is also known as MAP kinase kinase, phosphorylates MAP kinases on the conserved T-X-Y motif; phosphorylation of MAPK by MEK results in an increase in MAPK activity. MEK is involved in a diverse array of cellular processes such as stress-activated response, apoptosis, cytokine-induced cell proliferation, and DNA recombination during meiosis.

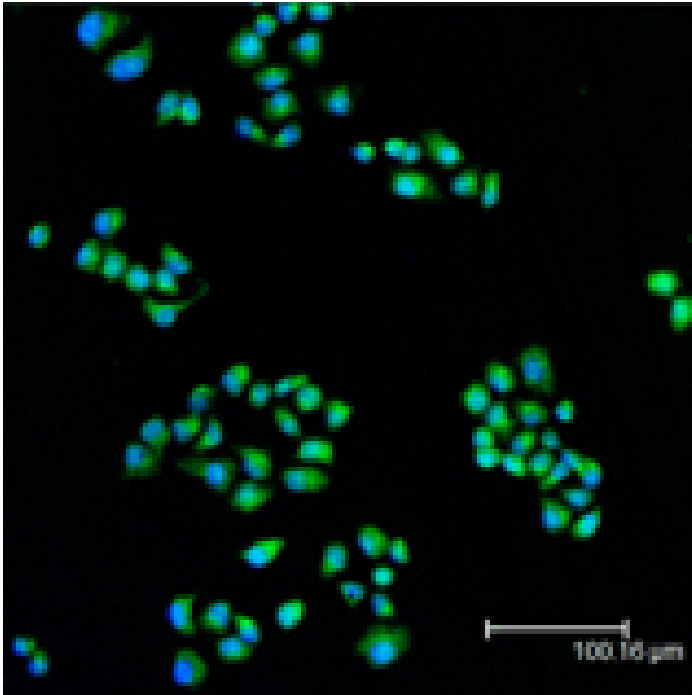
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

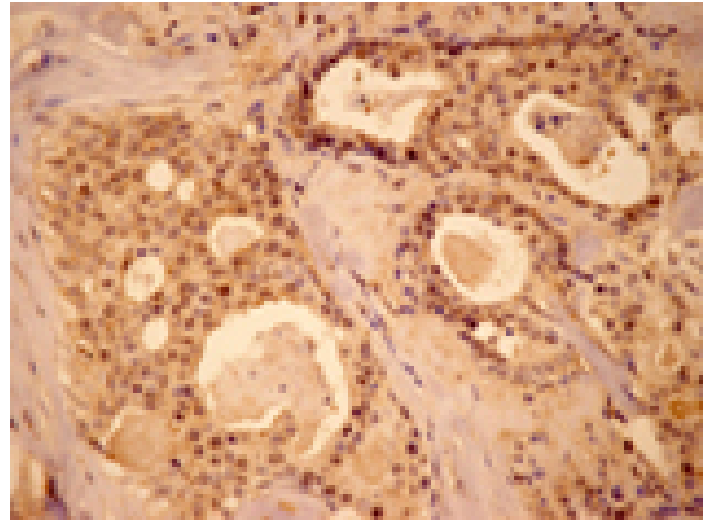
[View Online »](#)

Manuals, SDS & CofA

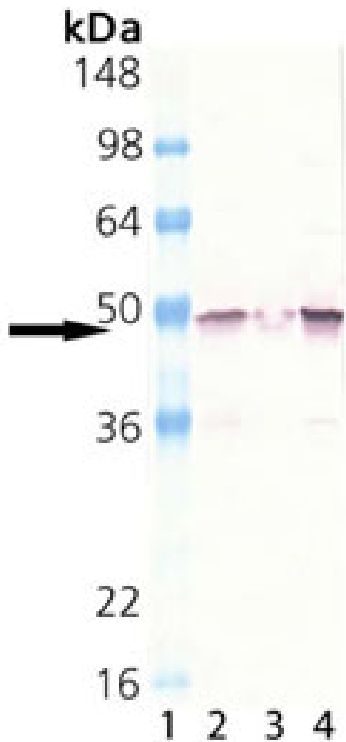
[View Online »](#)



Immunofluorescent analysis (confocal) staining of HeLa cells using MEK1 pAb (green); nuclei are stained in blue pseudocolor using DRAQ5.



Immunohistochemistry analysis of human breast cancer tissue with MEK1 pAb.



Western blot analysis of MEK1, pAb (Prod. No. ADI-KAP-MA010): Lane 1: MW marker; Lane 2: HeLa (heat shocked) (Prod. No. ADI-LYT-HL-101); Lane 3: Rat spleen extract; Lane 4: Mouse brain extract (Prod. No. ADI-LYT-MB100)

## Handling & Storage

**Handling** Avoid freeze/thaw cycles.

**Long Term Storage** -20°C

**Shipping** Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** MKK1, MAP kinase kinase 1

**Application** IF, IHC, WB

**Application Notes** Detects a band of ~48kDa by Western blot.

**Formulation** Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

**GenBank ID** L05624

**Host** Rabbit

**Immunogen** Synthetic peptide corresponding to the sequence near the N-terminus of human MEK1. The sequence is completely conserved in mouse, rat, hamster, rabbit and Xenopus.

**Purity Detail** Protein A affinity purified.

**Recommendation Dilutions/Conditions** Immunohistochemistry (1:50)Western Blot (1µg/ml, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

**Source** Purified from rabbit serum.

**Species Reactivity** Human, Mouse, Rat, Xenopus

**UniProt ID** Q02750

**Worry-free Guarantee** This antibody is covered by our [Worry-Free Guarantee](#)

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)