## 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 »

**Ordering Information** 

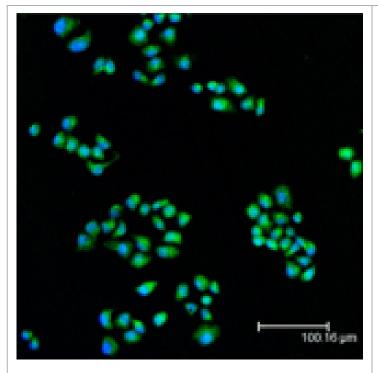
Order Online »

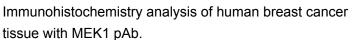
ADI-KAP-MA010-E

100µg

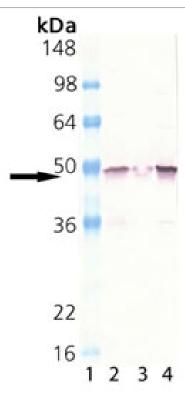
Manuals, SDS & CofA

**View Online** »





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



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



info-

eu@enzolifesciences.com