

MEK5 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: 3

[View Online »](#)

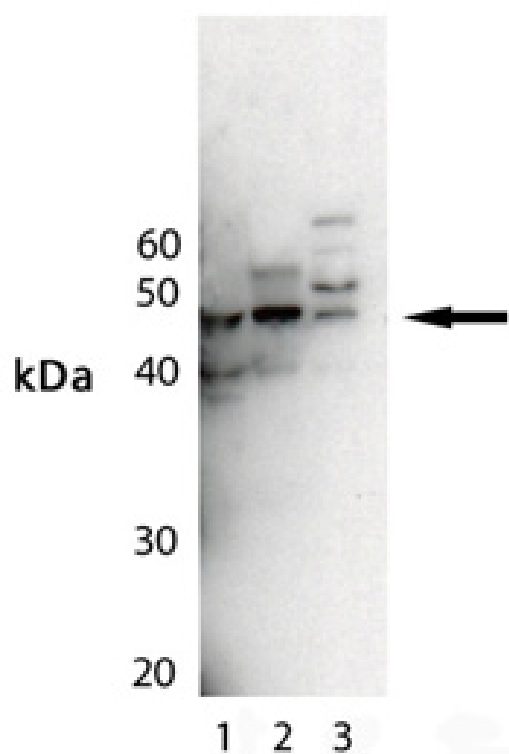
Ordering Information

[Order Online »](#)

ADI-KAP-MA003-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of MEK5, pAb (Prod. No. ADI-KAP-MA003): Lane 1: Jurkat, Lane 2: Mouse Testes, Lane 3: Rat Brain

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 MKK5, MAP kinase kinase 5

Application IHC, WB

Application Notes Detects a band of ~42kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

GenBank ID U71088

Host Rabbit

Immunogen Synthetic peptide corresponding to a portion of human MEK5. The sequence is completely conserved in rat.

Purity Detail Peptide affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:1,000, ECL)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 Dog, Human, Monkey, Mouse, Rabbit, Rat, Sheep, Xenopus

UniProt ID Q13163

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



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)