

MEKK3 polyclonal antibody

MEKKs (Mitogen activated protein kinase kinase kinases) are serine-threonine kinases that act as the first tier of cellular MAP kinase pathways by activation of MAP/ERK kinases, or MEKs. Many enzymes with MEKK activity have been identified, including MEKK1-4, Raf, MLK3, TAK, and DLK. MEKKs generally display little similarity outside of their catalytic kinase domains. MEKK1-4 are nearly 50% identical within their catalytic domains, and are known to regulate Erk, Jnk, and p38 MAP kinase pathways. MEKK2 and MEKK3 bind MEK5 via conserved PB1 domains, leading to downstream activation of Erk5.

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

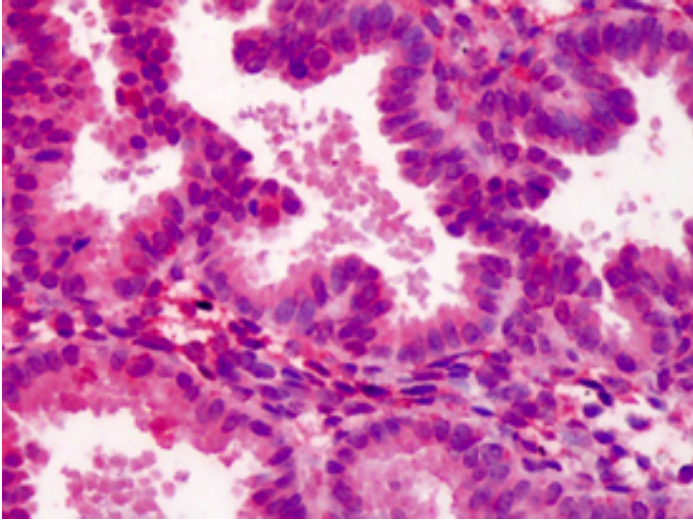
Ordering Information

[Order Online »](#)

ADI-KAS-MA013-E	100µl
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



IHC Analysis: Human uterus tissue stained with MEKK3, pAb at 1:100.

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 MAP3K3, MEK kinase 3

Application IHC (PS), WB

Formulation Liquid. Whole rabbit serum.

GenBank ID U43187

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the N-terminus of mouse MEKK3.

**Recommendation
Dilutions/Conditions** Western Blot (1:5,000) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Whole rabbit serum.

Species Reactivity Dog, Drosophila, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep, Xenopus

UniProt ID Q61084

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)