

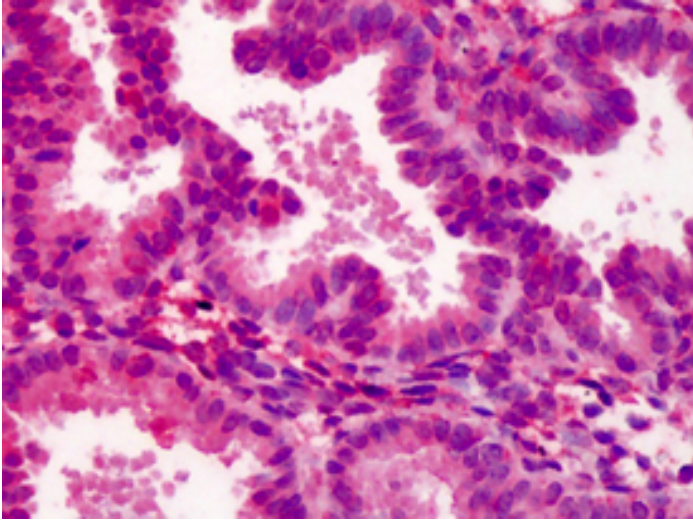
MEKK3 polyclonal antibody

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](#).

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