

MEKK1 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 »](#)

ENZ-ABS293-0100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)

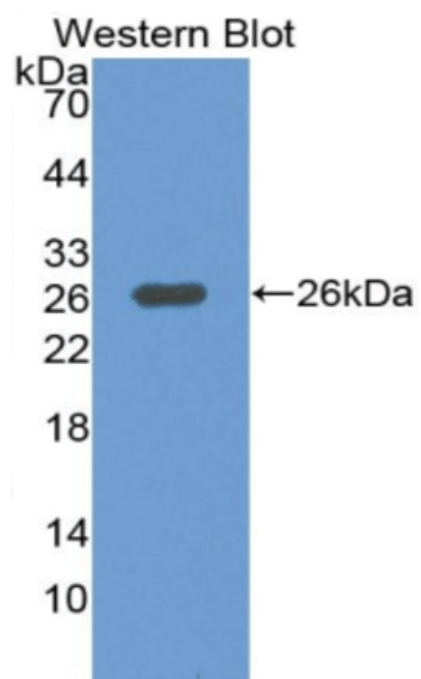


Figure 1. Western Blot analysis using MEKK1, pAb (Prod. No. ENZ-ABS293).

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Alternative Name	MAP3K1, MEK kinase 1, Mitogen-activated protein kinase kinase kinase 1, MAPK/ERK kinase kinase 1, MAPKKK1
Application	ELISA, ICC, IHC (FS), IHC (PS), WB
Application Notes	Detects a band of ~26KDa by Western blot.
Formulation	Liquid. In PBS, pH 7.4, containing 0.02% sodium azide and 50% glycerol.
Host	Rabbit
Immunogen	Recombinant MEKK1 (aa 1307-1509).
Positive Control	Included.
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	ELISA (1:100-200)Immunocytochemistry (formalin fixed cells) (1:100-500)Immunohistochemistry (frozen sections) (1:100-500)Immunohistochemistry (paraffin sections) (1:50-200)Western blot (1:100-400)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q13233
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)