

TRAF1 monoclonal antibody (3D4)

TRAF proteins are multifunctional signal adaptors that are involved in the activation of the kinase cascades which activate c-Jun N-terminal kinase, p38 mitogen activated protein kinase, and the transcription factor NF- κ B. Members of the TRAF family form homodimers as well as heterodimers, resulting in extensive diversity and specificity to their function in signal transduction pathways.

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

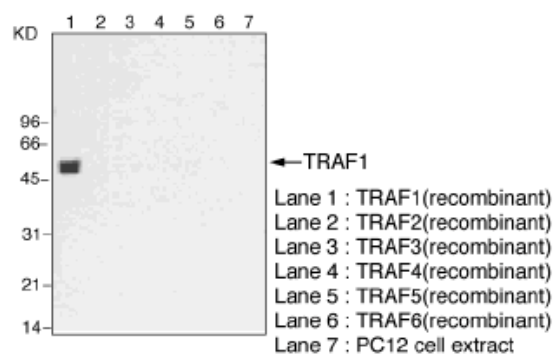
Ordering Information

[Order Online »](#)

ADI-AAM-421-E	100 μ g
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis probed with AAM-421.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TNF receptor-associated factor 1
Application	WB
Application Notes	Western Blot (1µg/ml, ECL) Optimal conditions must be determined individually for each application.
Clone	3D4
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol. No preservatives.
Host	Mouse
Immunogen	Recombinant mouse TRAF1.
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1µg/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
UniProt ID	P39428
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)