

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](#).

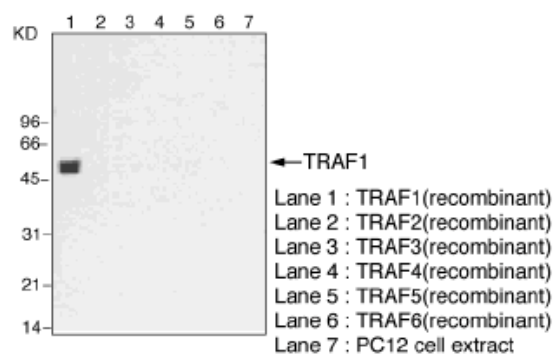
Ordering Information

[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 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com