

TRAF6 polyclonal antibody

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](#).

Citations: 3

[View Online »](#)

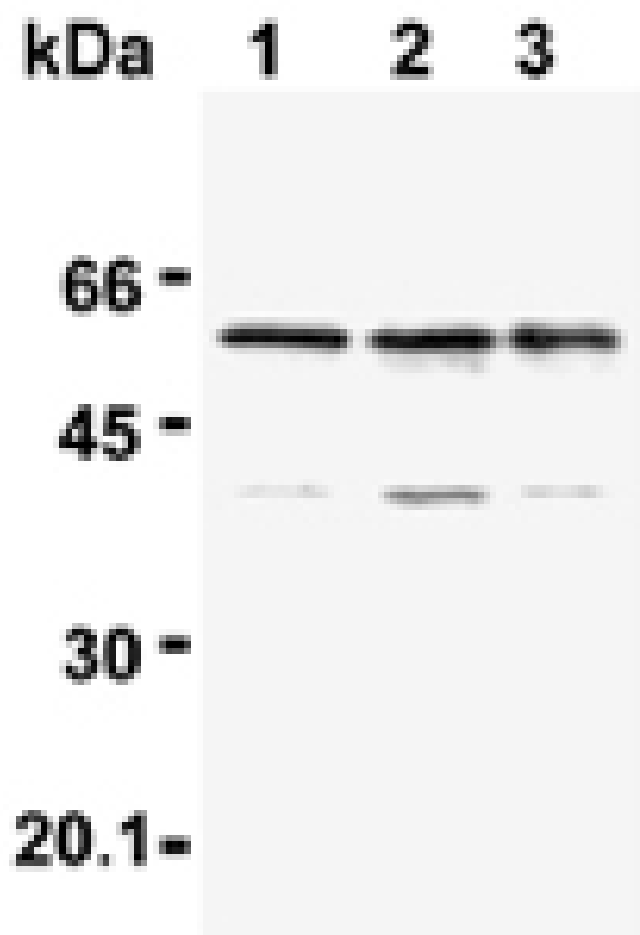
Ordering Information

[Order Online »](#)

ADI-AAP-426-E	100 μ l
---------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of L5178Y (1), WR19L (2) and NIH/3T3 (3) cells, probed with TRAF6 pAb.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name TNF receptor-associated factor 6

Application IP, WB

Application Notes Detects a band of ~58kDa by Western blot.

Formulation Liquid. In PBS containing 50% glycerol.

Host Rabbit

Immunogen Recombinant mouse TRAF6 (N-terminal fragment).

Purity Detail Immunoaffinity purified.

Recommendation Immunoprecipitation (2-4µL/100-500 µl cell extract from 5×10⁶ cells)Western Blot
Dilutions/Conditions (1:1000, 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 P70196

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)