

Fas monoclonal antibody (ZB4)

Fas/APO-1/CD95 is a 36 kDa cell surface type I-membrane glycoprotein that is a member of the Tumor necrosis factor receptor (TNF-R) family, which includes TNF-R1, TNF-R2, CD27, CD30, and CD40. Binding of Fas-ligand (FasL) to Fas or crosslinking of Fas by anti-Fas monoclonal antibodies leads to apoptosis of Fas expressing cells. Apoptosis induced by Fas-Fas ligand interaction is inhibited by a number of cytosolic factors such as the ICE/CED3 family inhibitor CrmA, and the short and long form of FLIP.

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

Citations: 3

[View Online »](#)

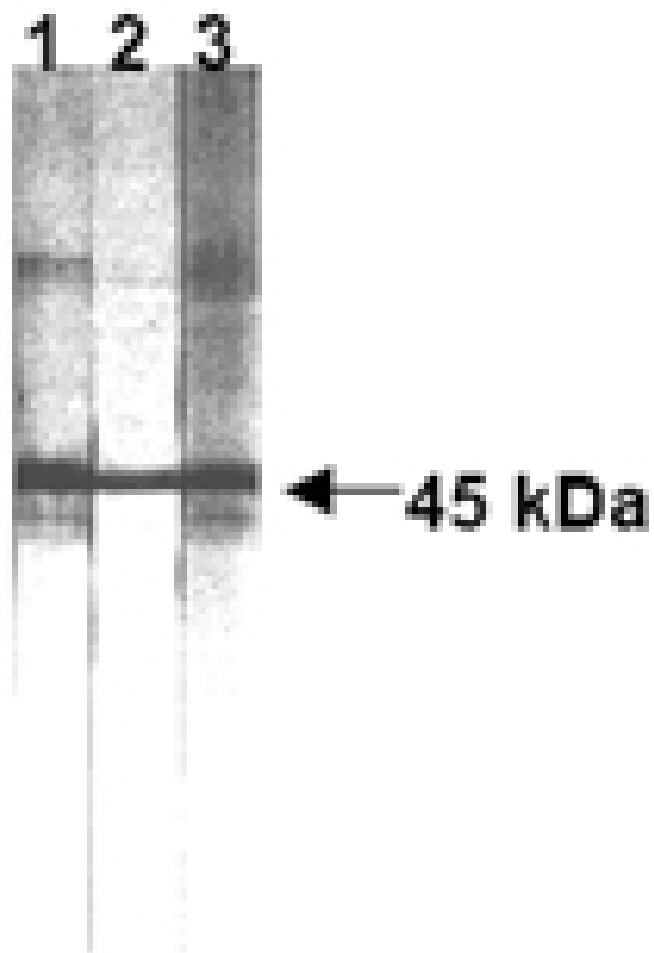
Ordering Information

[Order Online »](#)

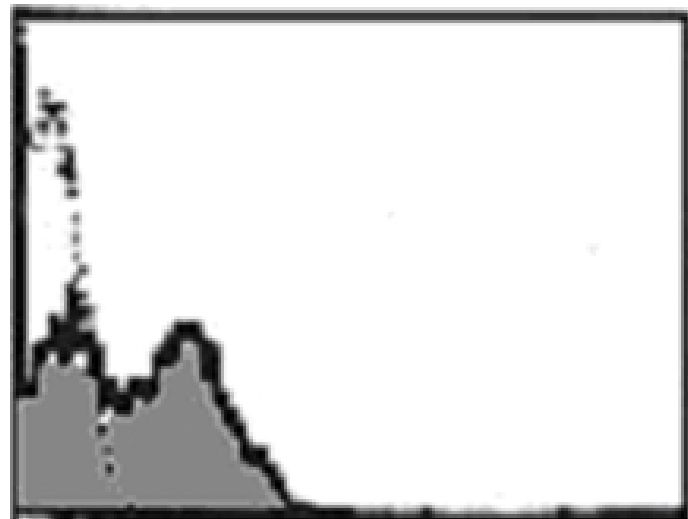
ADI-AAM-227-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of human Fas expression in human Fas transfected cells; Lane1: anti-human Fas (ZB4), Lane2: anti-human Fas (UB2) and Lane3: anti-human Fas (CH11).



Flow cytometric analysis of human Fas expression on normal human lymphocytes. Open histogram indicates the reaction of isotypic control to the cells. Shaded histogram indicates the reaction of Fas mAb (ZB4) to the cells.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name CD95, APO-1, TNFRSF 6

Application Flow Cytometry, FUNC, WB

Application Notes Detects a band of ~45kDa by Western blot.
Functional Application: neutralizing.

Clone ZB4

Formulation Liquid. In PBS, pH 7.2, with 50% glycerol.

Host Mouse

Immunogen Recombinant human Fas.

Isotype IgG1

Purity Detail Protein A affinity purified.

Recommendation Flow Cytometry (5-20µg/ml)Western Blot (5µg/ml, colorimetric)Suggested
Dilutions/Conditions dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

UniProt ID P25445

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)