

Bcl-2 monoclonal antibody (83-8B)

The Bcl-2 family of proteins regulate the intrinsic apoptotic pathway and are characterized by the sharing of at least one of four conserved Bcl-2 Homology (BH) domains. Bcl-2 family members function as either pro- or anti-apoptotic factors, and are generally categorized into three groups based on structural homology and function. Bcl-2-like proteins (Bcl-2, Bcl-xL, Bcl-w, A1, and Mcl-1) are pro-survival. Pro-apoptotic members can be sub-divided into two groups: the multi-BH domain group, including Bax, Bak, and Bok; and the more numerous and diverse BH3-only group, which includes Bim, Bad, and Bid.

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

Citations: 4

[View Online »](#)

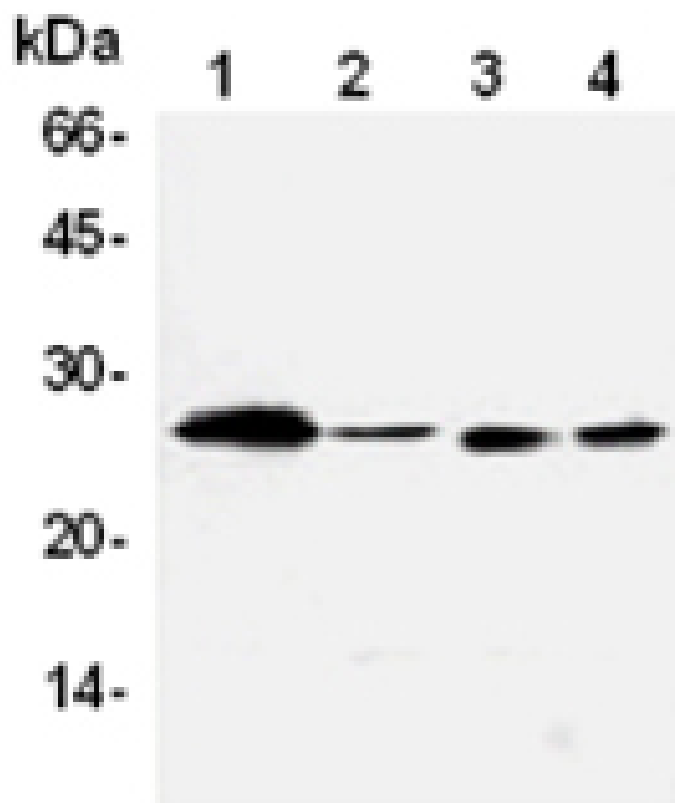
Ordering Information

[Order Online »](#)

ADI-AAM-072-E	100µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Jurkat (1), Raji (2), WR19L (3), and PC12 (4) cell lysates , probed with Bcl-2 mAb (83-8B).

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Flow Cytometry, WB

Application Notes Detects a band of ~26kDa by Western blot.

Clone 83-8B

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

Gene/Protein Identifier NP_058689 (RefSeq)

Host Mouse

Immunogen Recombinant rat Bcl-2 β .

Isotype IgG1

Purity Detail Protein A affinity purified.

Recommendation
Dilutions/Conditions Flow Cytometry (10 μ g/ml)Western Blot (1 μ g/ml, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Drosophila, Human, Mouse, Rat

UniProt ID P49950

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-
fr@enzolifesciences.com](mailto:info-fr@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)