

Bcl-w monoclonal antibody (16H12)

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

Citations: 3

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

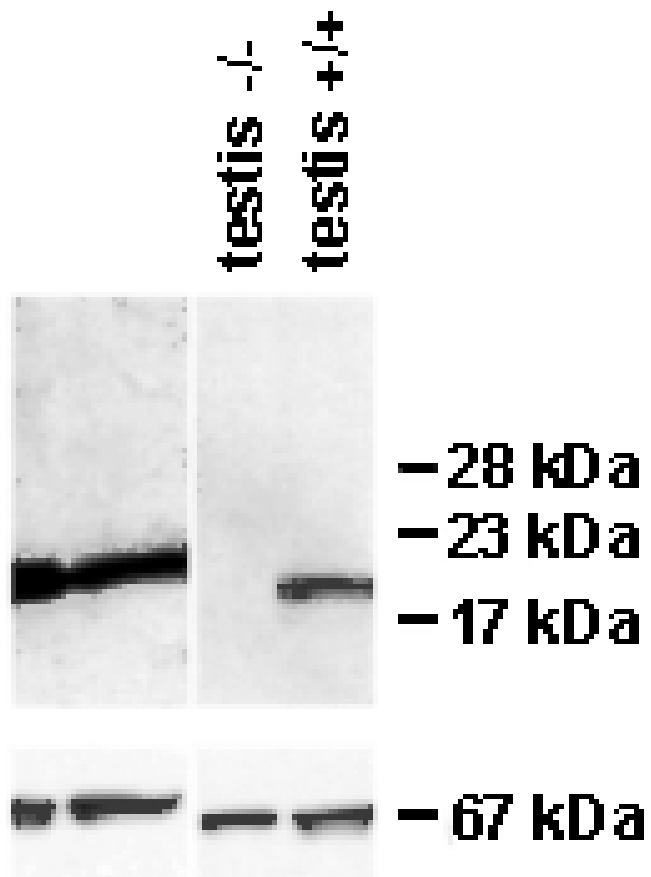


Figure: Western blot showing the detection of Bcl-w in fractionated lysates of mouse tissues with clone 16H12. Antibody specificity is indicated by the absence of the expected 21 kDa protein in *Bcl-w*^{-/-} testis control tissues. The 67 kDa band is Hsp-70 used as an equal loading control.

Handling & Storage

Use/Stability	Stable for up to 3 months when stored at +4°C.
Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	BCL2L2
Application	IP, WB
Clone	16H12
Formulation	Liquid. In PBS containing 25% glycerol and 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human Bcl-w.
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Species Reactivity	Dog, Human, Monkey, Mouse, Rat
Technical Info / Product Notes	Western Blot works for transfected and Bcl-w overproducing cells. For detection of endogenous Bcl-w, immunoprecipitation is required.
UniProt ID	Q92843
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



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)