

# Bcl-w monoclonal antibody (16H12)

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

Citations: 3

[View Online »](#)

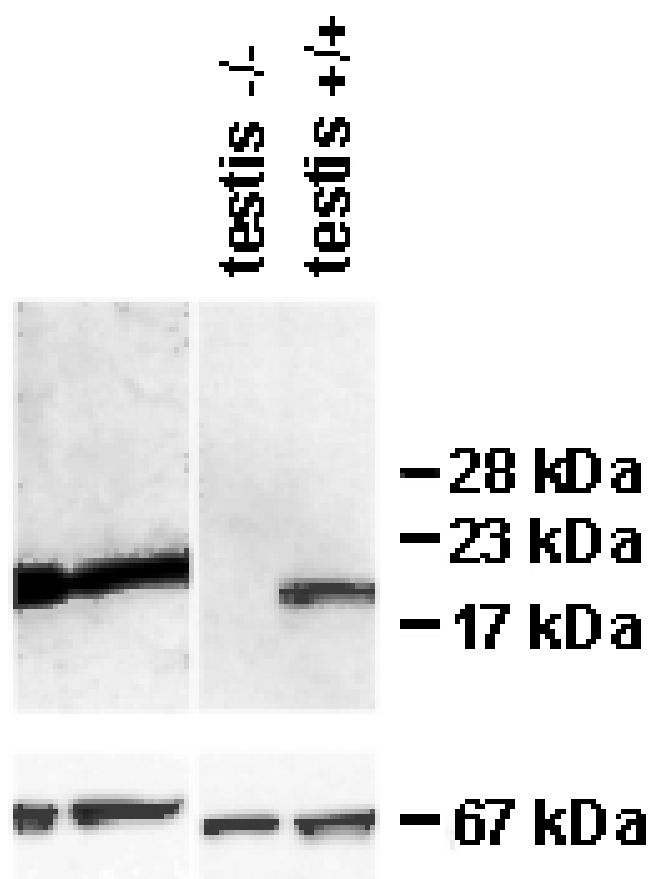
## Ordering Information

[Order Online »](#)

ALX-804-354-C100	100µg
------------------	-------

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 <a href="#">Worry-Free Guarantee</a> .



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)