

# Bcl-w monoclonal antibody (13F9)

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

Citations: 4

[View Online »](#)

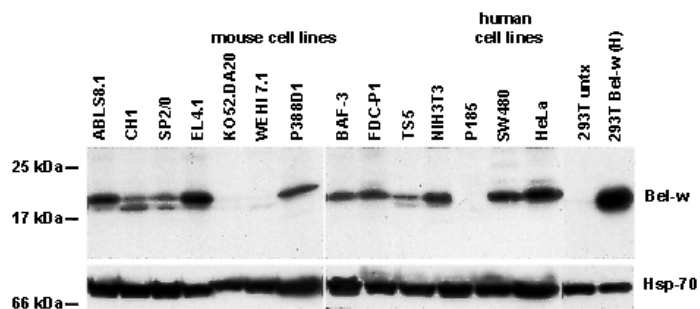
## Ordering Information

[Order Online »](#)

ALX-804-353-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



**Figure:** Immunoprecipitation/Western blot showing the detection of Bcl-w in human and mouse cell lines. Lysates from  $10^7$  cells were immunoprecipitated by clone 13F9 and the fractionated proteins were then blotted with the same antibody. Hsp-70 expression was used as an equal loading control, and 293T cells transiently transfected with a human *bcl-w* expression construct were used as a positive control.

## Handling & Storage

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
Application Notes	Excellent for Immunoprecipitation.
Clone	13F9
Formulation	Liquid. In PBS containing 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
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-  
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)