

# Bim<sub>S/EL/L</sub> monoclonal antibody (3C5)

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

Citations: 19

[View Online »](#)

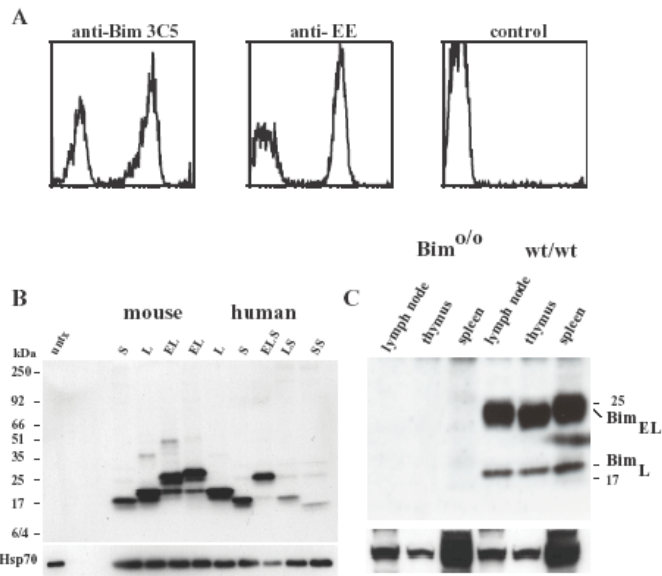
## Ordering Information

[Order Online »](#)

ALX-804-527-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



**Figure A:** Hybridomas producing MAbs to Bim (peptide) were identified using clones of the IL-3-dependent mouse myeloid cell line FDC-P1 stably expressing Bcl-2 alone, or Bcl-2 plus mouse BimL mixed at a 1:1 ratio. Transfected cells were fixed, permeabilized and stained with the rat anti-Bim MAb (clone 3C5) (Prod. No. ALX-804-527), the mouse anti-Glu-Glu (EE) MAb (positive control) or with secondary antibody alone (negative control). Staining was visualised by FITC-conjugated goat anti-rat or anti-mouse IgG antibodies. **Figure B:** Western blotting with anti-Bim MAb (clone 3C5) (Prod. No. ALX-804-527) recognizes all known isoforms of Bim. Western blotting of lysates of 293T cells transfected with EE-tagged forms of mouse or HA tagged forms of human Bim. Blots were probed with anti-Bim MAb (clone 3C5) (Prod. No. ALX-804-527), which detected EE mouse BimS, BimL, BimEL and EE human, BimSS, BimS, BimLS, BimL, BimELS, BimEL. Probing with an anti-Hsp-70 MAb was used as a loading control for all blots. **Figure C:** Anti-Bim MAb (clone 3C5) (Prod. No. ALX-804-527) detects endogenous Bim expression in mouse tissues but not those from Bim knockout mice. Bim protein was revealed in lysates from mouse tissues ( $5 \times 10^6$ /lane) by Western blotting with anti-Bim MAb (clone 3C5) (Prod. No. ALX-804-527) and detection by ECL. Probing with an anti-Hsp-70 MAb was used as a loading 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	BCL2L11
Application	ELISA, Flow Cytometry, ICC, IP, WB
Clone	3C5
Formulation	Liquid. In PBS containing 0.02% sodium azide and 25% glycerol.
Host	Rat
Immunogen	Synthetic peptide corresponding to aa 20-40 (Q20PAERPPQLRPGAPTSLQTEP40) of mouse BimL.
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (2µg/ml)Western Blot (1-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Dog, Human, Monkey, Mouse, Rat
Specificity	Recognizes an epitope within aa 20-40 of im <sub>L</sub> (~19kDa), Bim <sub>EL</sub> (~23kDa), and Bim <sub>S</sub> (~16kDa). Also recognizes the six novel isoforms of human Bim, generated by alternative splicing and lacking the C-terminal hydrophobic region recently identified.
UniProt ID	O54918

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)