

Bim_{S/EL/L} monoclonal antibody (10B12)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-528-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

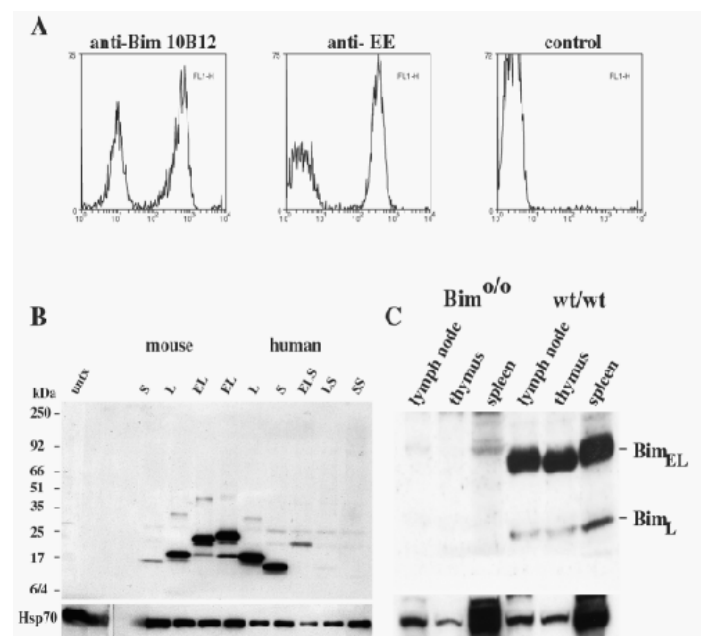


Figure A: Hybridomas producing MAb to Bim (peptide) were identified using clones of the IL-3-dependent mouse myeloid cell line FDC-P1 stably expressing Bcl-2 alone, or Bcl-2 plus mouse BimL mixed at a 1:1 ratio. Transfected cells were fixed, permeabilized and stained with the rat anti-Bim MAb (clone 10B12) ([Prod. No. ALX-804-528](#)), 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 10B12) ([Prod. No. ALX-804-528](#)) 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 10B12) ([Prod. No. ALX-804-528](#))

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, IHC, IP, WB
Clone	10B12
Formulation	Liquid. Protein G affinity purified antibody in PBS containing 25% glycerol and 0.02% sodium azide.
Host	Rat
Immunogen	Synthetic peptide corresponding to aa 20-40 (Q20PAERPPQLRPGAPTSLQTEP40) of mouse BimL.
Isotype	IgG2a
Recommendation	Immunohistochemistry (only EM)Immunoprecipitation (2µg/ml)Western Blot (1-
Dilutions/Conditions	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 Bim _L , Bim _{EL} , and Bim _S . Also recognizes the six novel isoforms of human Bim, generated by alternative splicing and lacking the C-terminal hydrophobic region recently identified.



ENZO LIFE SCIENCES
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

054918

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com

Worry-free Guarantee

This antibody is covered by our **Worry-Free Guarantee.**