## Bim<sub>S/EL/L</sub> 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** »

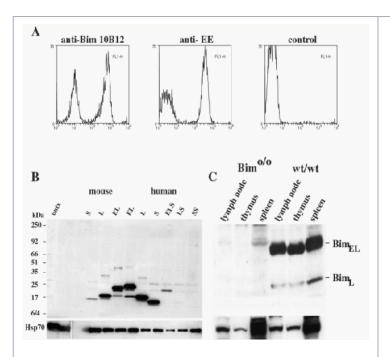


Figure A: Hybridomas producing MAbs to Bim (peptide) were identified using clones of the IL-3-dependent mouse myeloid cell line FDC-P1 stably expressing Bcl-2 alone, or Bcl-2 plus mouse BimL mixed at a 1:1 ratio. Transfected cells were fixed, permeabilized and stained with the rat anti-Bim MAb (clone 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 isforms of Bim. Western blotting of lysates of 293T cells 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 Dilutions/Conditions Immunohistochemistry (only EM)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 Bim<sub>L</sub>, Bim<sub>EL</sub>,

and Bim<sub>S</sub>. 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com