

Bim_{S/EL/L} monoclonal antibody (14A8)

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

Citations: 9

[View Online »](#)

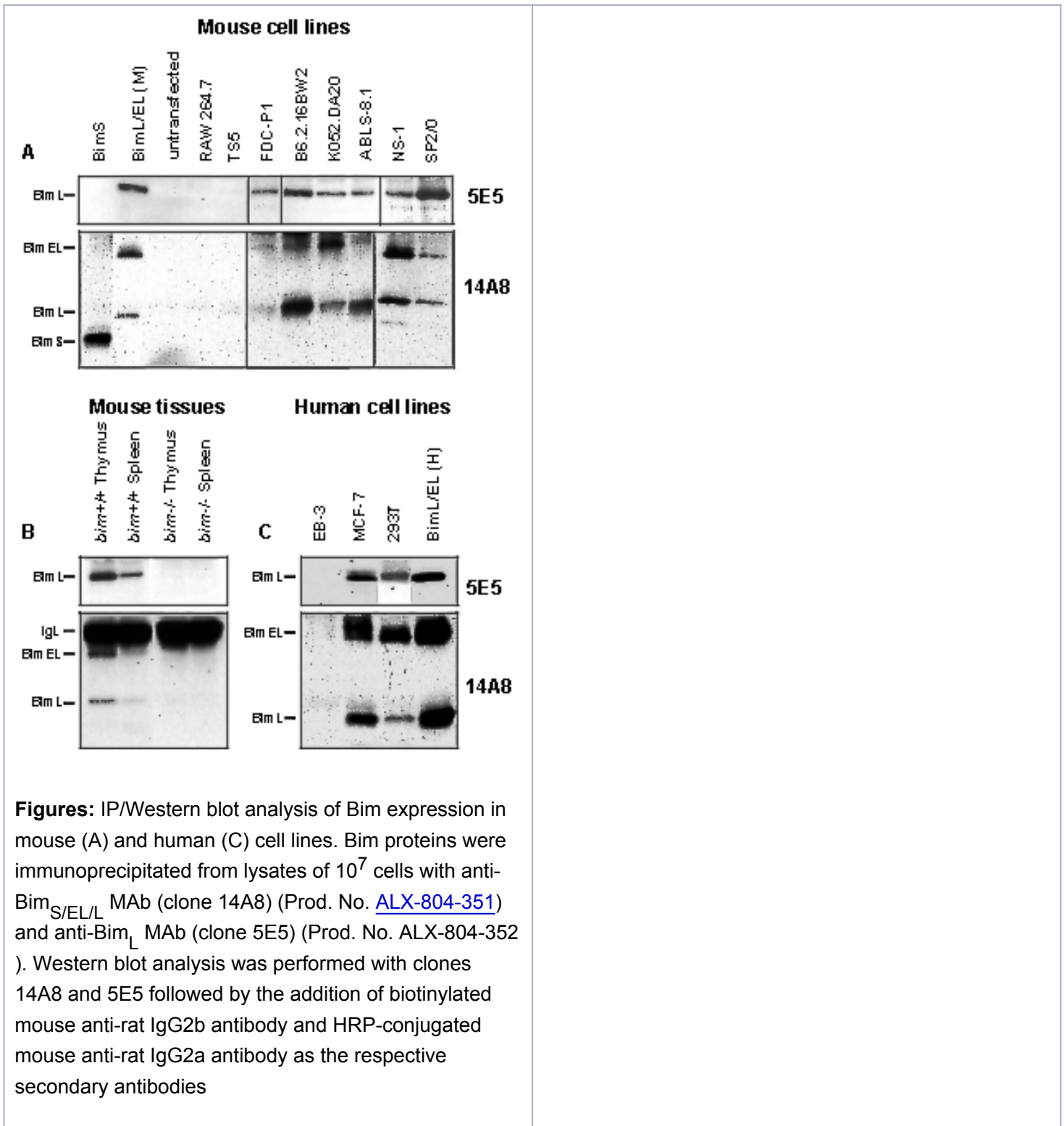
Ordering Information

[Order Online »](#)

ALX-804-351-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



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	Flow Cytometry, ICC, IHC (PS), IP, WB
Clone	14A8
Formulation	Liquid in PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse BimS.
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (0.5-1µg/ml; overexpressing cell lines)Immunocytochemistry (1-5µg/ml)Immunohistochemistry (paraffin sections, 1-5µg/ml)Immunoprecipitation (1-2µg/ml)Western Blot (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Monkey, Mouse
Specificity	Recognizes Bim _L , Bim _{EL} and Bim _S .

Technical Info / Product Notes

As Bim expression is generally low in non-transfected cell lines, please take care when choosing positive controls. Examples of cell lines which express detectable levels of Bim include: B6.2.16.BW2 or KO52.DA20 (T cell lymphomas) or ABL5.8.1 (B cell lymphoma) or MCF-7s. It is generally necessary to use a high amount of lysate (e.g. 1×10^7 cells). When using tissue samples, it is generally necessary to use 1-5mg of total protein from an appropriate tissue source.

UniProt ID

O54918

Worry-free Guarantee

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

Last modified: May 29, 2024



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

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-be@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