

BAFF (mouse) monoclonal antibody (5A8)

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

Citations: 3

[View Online »](#)

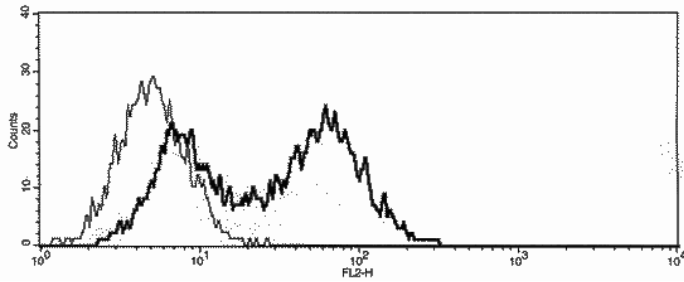
Ordering Information

[Order Online »](#)

ALX-804-158-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Method: HEK 293T cells were mock transfected (thin line) or transfected with an expression plasmid enabling surface expression of the extracellular domain of mouse BAFF (thick line). Cells (5×10^5) were incubated on ice for 30 min in 50 μ l FACS buffer (PBS, 5% fetal calf serum, 0.02% sodium azide) containing 1 μ g/ml of MAb to BAFF (mouse) (5A8). After washing in FACS buffer, R-PE-conjugated antibody to rat IgG (Prod. No. [ALX-211-053](#)) was added. Cells were incubated on ice for 30 min, washed and then analyzed by flow cytometry. """,BAFF (mouse) monoclonal antibody (5A8) Flow Cytometry,BAFF (mouse) monoclonal antibody (5A8) Flow Cytometry,yes ALX-804-192,804-192.gif,<http://static.enzolifesciences.com/fileadmin/files/image/804-192.gif>,**Figure:** Photomicrographs of thymus from a mouse treated with hydrocortisone (left) and of a control treated animal (right). Sections of formalin-fixed tissues were heated in formamide

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
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	B cell-activating factor, BLyS, TALL-1, THANK, zTNF4, TNFSF 13B/20, CD257
Application	ELISA, Flow Cytometry, WB
Application Notes	Excellent for Flow Cytometry.
Clone	5A8
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Rat RBL cells expressing mouse BAFF.
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (capture)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Mouse

UniProt ID

Q9WU72



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-
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)

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