

# BAFF (mouse) monoclonal antibody (1C9)

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

Citations: 2

[View Online »](#)

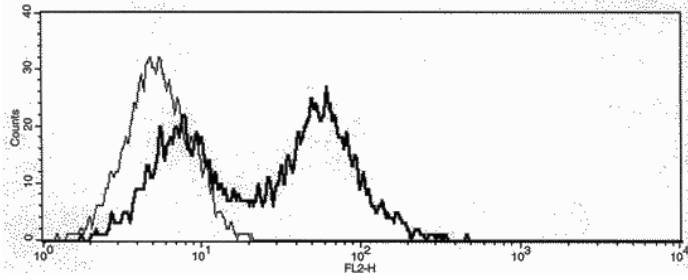
## Ordering Information

[Order Online »](#)

ALX-804-157-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 ( $5 \times 10^5$ ) were incubated on ice for 30 min in 50  $\mu$ l FACS buffer (PBS, 5% fetal calf serum, 0.02% sodium azide) containing 1  $\mu$ g/ml of MAb to BAFF (mouse) (1C9). 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.

## 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
Clone	1C9
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Rat RBL cells expressing mouse BAFF.
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (detection)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
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-  
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)

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