

# BAFF (human) monoclonal antibody (Buffy-1)

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

Citations: 5

[View Online »](#)

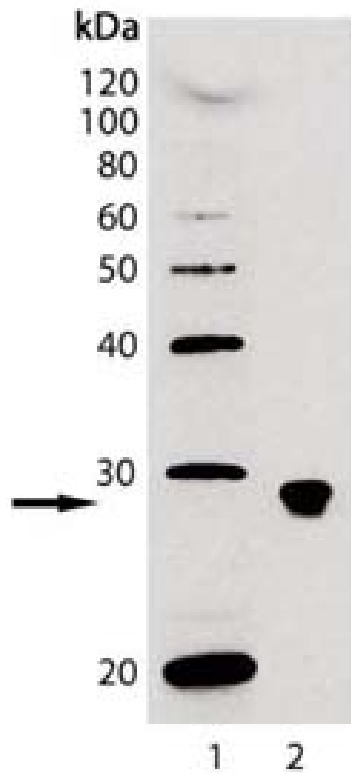
Ordering Information

[Order Online »](#)

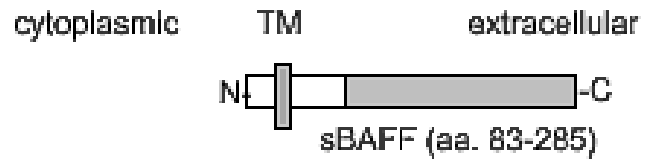
ALX-804-128-C100	100µg
------------------	-------

Manuals, SDS & CofA

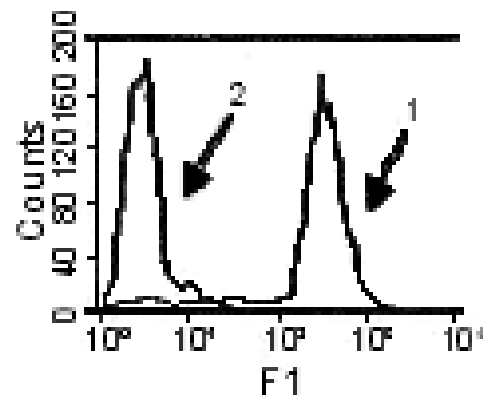
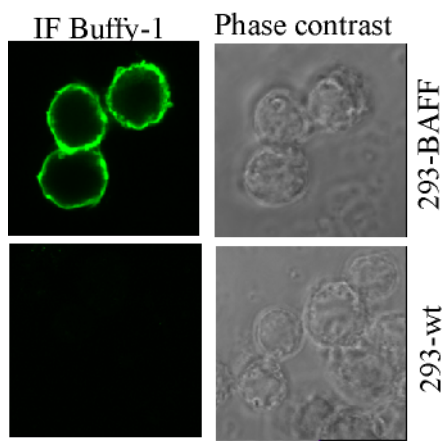
[View Online »](#)



Western blot analysis of BAFF (human), mAb (Buffy-1) (Prod. No. ALX-804-128): Lane 1: MW marker; Lane 2: BAFF (soluble) (human), (recombinant) (Prod. No. ALX-522-025).



Schematic structure of human soluble BAFF (rec.) (aa 83-285). Recognizes the stalk region present in the membrane-bound form of hBAFF.



Membrane-bound human BAFF detected by MAb to BAFF (human) (Buffy-1) (Prod. No. [ALX-804-128](#)). Left panel: ICC; Right panel: phase contrast. Lower panel: 293-wt control. BAFF (human) monoclonal antibody (Buffy-1) image, BAFF (human) monoclonal antibody (Buffy-1) image, yes

ALX-804-128, 804-128-

4.gif, <http://static.enzolifesciences.com/fileadmin/files/image7804-128-4.gif>, Flow cytometric profile of 293 cells stably transfected with a BAFF-expression plasmid (1). Control (2) are mock transfected 293 cells. Optimal concentration range is 0.5-5  $\mu\text{g/ml}$ . **Method:** 293T cells ( $5 \times 10^5$ ) expressing surface BAFF vector or Mock transfected

Flow cytometric profile of 293 cells stably transfected with a BAFF-expression plasmid (1). Control (2) are mock transfected 293 cells. Optimal concentration range is 0.5-5  $\mu\text{g/ml}$ . **Method:** 293T cells ( $5 \times 10^5$ ) expressing surface BAFF vector or Mock transfected, were incubated on ice for 30 min. in 50  $\mu\text{l}$  FACS buffer (PBS, 5% fetal calf serum, 0.02% azide) containing 4  $\mu\text{g/ml}$  of MAb to BAFF (human) (Buffy-1). After washing in FACS buffer, FITC-conjugated antibody to rat IgG (ALX-211-051) was added and cells were analyzed by flow cytometry.

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 6 months after receipt when stored as recommended.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	B cell-activating factor, BLyS, TALL-1, THANK, zTNF4, TNFSF 13B/20, CD257
<b>Application</b>	Flow Cytometry, ICC, WB
<b>Clone</b>	Buffy-1
<b>Formulation</b>	Liquid. In PBS containing 0.02% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human soluble BAFF (aa 83-285) (Prod. No. ALX-522-025).
<b>Isotype</b>	IgG2a
<b>Source</b>	Purified from mouse ascites.
<b>Species Reactivity</b>	Human
<b>Specificity</b>	Recognizes membrane-bound BAFF. Detects endogenous protein by FC.
<b>UniProt ID</b>	Q9Y275
<b>Worry-free Guarantee</b>	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



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)