

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

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

Citations: 6

[View Online »](#)

## Ordering Information

[Order Online »](#)

ALX-804-128F-T100	FITC - 100 tests
-------------------	------------------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as recommended.
Handling	Do not freeze. Protect from light.
Short Term Storage	+4°C
Long Term Storage	+4°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	Flow Cytometry, ICC, WB
Clone	Buffy-1
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human soluble BAFF (aa 83-285) (Prod. No. ALX-522-025).
Isotype	IgG2a
Source	Purified from mouse ascites.
Species Reactivity	Human
Specificity	Recognizes membrane-bound BAFF. Detects endogenous protein by FC.
Technical Info / Product Notes	1 test means: 1µl of a 1.0mg/ml stock solution was used to stain 200'000 cells in a sample volume of 50µl.

UniProt ID

Q9Y275



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**.