

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

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