

BAFF monoclonal antibody (Buffy-2)

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

Citations: 19

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-131-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

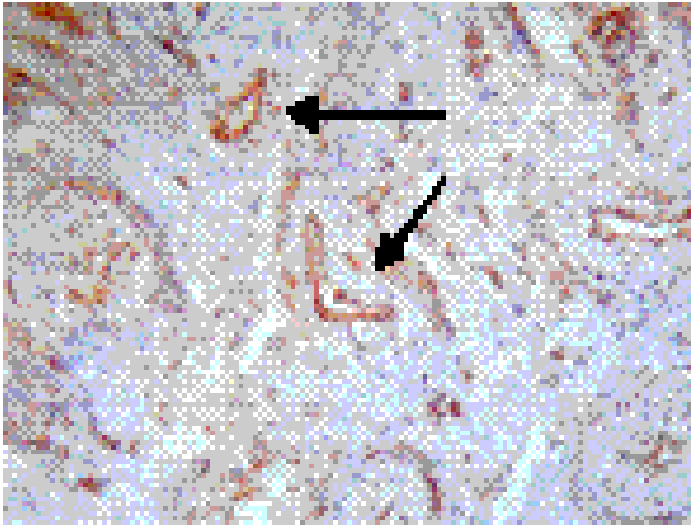
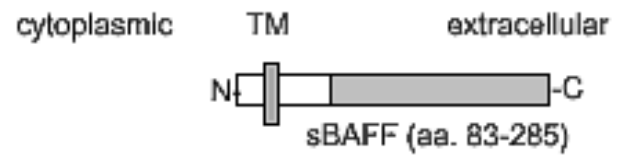


Figure: Staining of paraffin section of synovial tissue from a patient with Rheumatoid Arthritis. Arrows indicate vessels (20x) staining positive for BAFF using Buffy-2.

Method: Tissues are in paraffin sections. Sections are de-waxed and rehydrated, then cooked. Sections are stained in wet boxes, rehydrated and finally prepared as coverslip slides using Eukitt solution.



Schematic structure of rhsBAFF (aa. 83-285). MAb to BAFF (Buffy-2) recognizes the TNF-homology domain.

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles.
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, IHC (FS), IHC (PS), WB
Clone	Buffy-2
Crossreactivity	Weakly cross-reacts with mouse BAFF.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human soluble BAFF (aa 83-285) (Prod. No. ALX-522-025).
Isotype	IgM
Purity Detail	Purified by IgM purification HP column.
Source	Purified from ascites.
Species Reactivity	Human
Specificity	Recognizes membrane-bound and soluble BAFF.
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](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)