

# BAFF monoclonal antibody (Buffy-2)

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

Citations: 19

[View Online »](#)

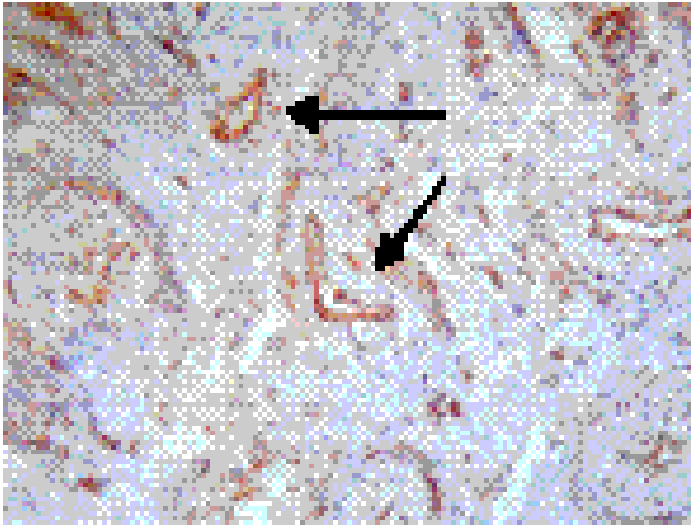
## Ordering Information

[Order Online »](#)

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

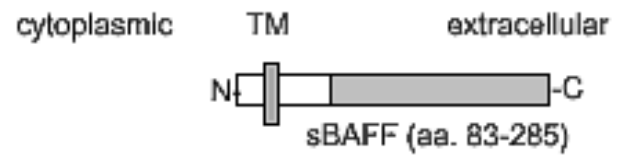
Manuals, SDS & CofA

[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



ENZO LIFE SCIENCES,  
INC.  
Phone: +1 940.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(UK) LTD  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
ENZO LIFE SCIENCES  
Phone: +32 3 966 0426  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[de@enzolifesciences.com](mailto: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**