

# BAFF receptor monoclonal antibody (BAFFR/1557)

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

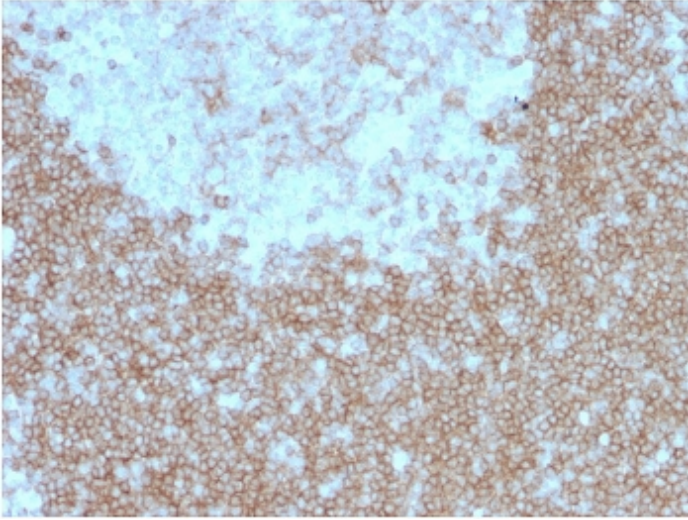
## Ordering Information

[Order Online »](#)

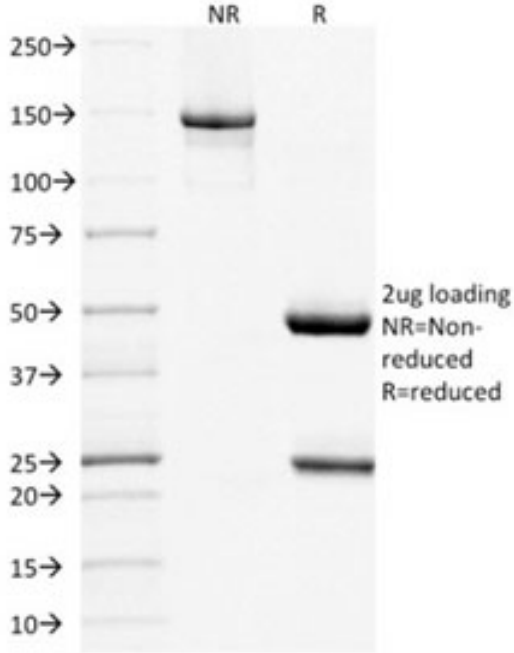
ENZ-ABS870-0100	100µg
-----------------	-------

## Manuals, SDS & CofA

[View Online »](#)



IHC staining of FFPE human tonsil with BAFF Receptor antibody (clone BAFFR/1557). HIER: boil tissue sections in pH 9 10mM Tris with 1mM EDTA for 10-20 min and allow to cool before testing.



SDS-PAGE analysis of purified, BSA-free BAFF Receptor antibody (clone BAFFR/1557) as confirmation of integrity and purity.

# Handling & Storage

Handling	Store the BAFF Receptor antibody at 2-8°C (with azide) or aliquot and store at -20°C or colder (without azide).
Long Term Storage	-20°C
Shipping	Blue Ice

## Regulatory Status

RUO - Research Use Only

# Product Details

Alternative Name	BAFF-R, B-cell-activating factor receptor, BLyS receptor 3, TNFRSF13C, Tumor necrosis factor receptor superfamily member 13C, CD268
Application	IHC (PS)
Clone	BAFFR/1557
Formulation	Liquid. 0.2 mg/ml in 1X PBS containing 0.1 mg/ml BSA (US sourced), 0.05% sodium azide.
Host	Mouse
Immunogen	Recombinant human full-length BAFF.
Isotype	IgG1κ
Purity Detail	Protein G affinity chromatography purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (FFPE) (1-2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
UniProt ID	Q96RJ3
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .



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)