

BAFF receptor (CT) BR3 polyclonal antibody

BAFF, or B-cell Activating Factor, is a member of the TNF superfamily also known as BLyS (B Lymphocyte Stimulator), TALL-1 (TNF- and ApoLrelated Leukocyte-expressed Ligand), and THANK (TNF Homologue that Activate Apoptosis, NF-κB and c-jun N-terminal Kinase). BAFF/BLyS was characterized as a B-cell activator since it induced B-cell proliferation and immunoglobulin secretion. Three receptors for BAFF were identified and designated TACI, BCMA, and BAFF-R. BAFF and its receptors are essential for B-cell development, survival, and humoral immune responses.

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

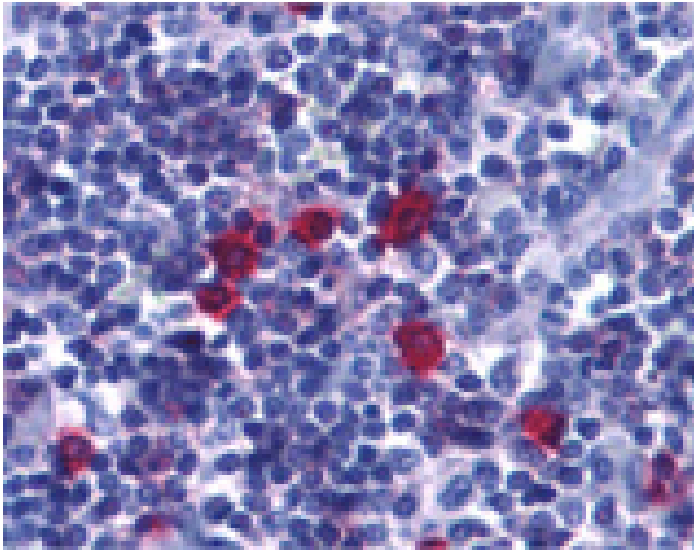
Ordering Information

[Order Online »](#)

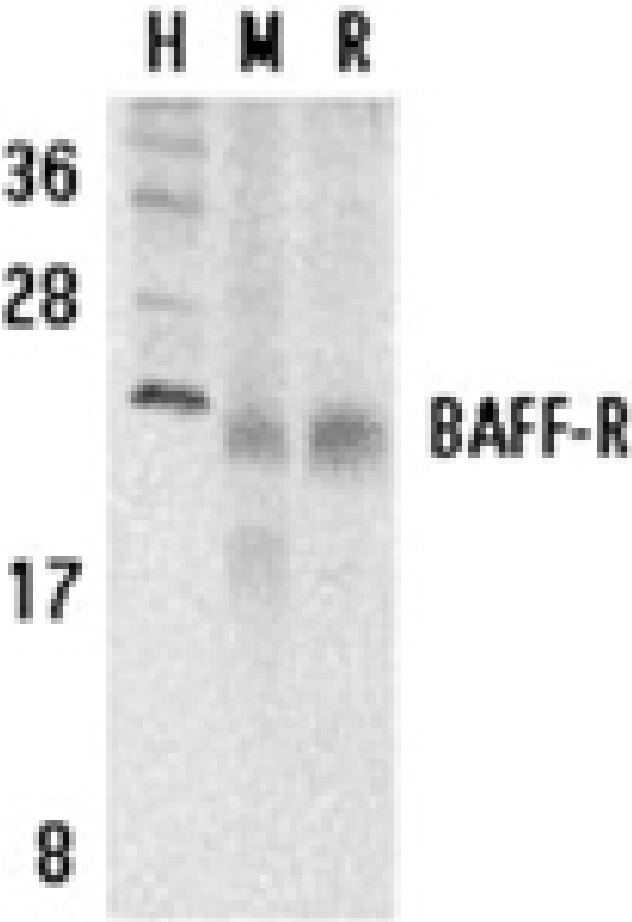
ADI-905-305-100	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of BAFF Receptor in human tonsil tissue with BAFF Receptor antibody at 5µg/ml.



Western blot analysis of BAFF Receptor in human (H), mouse (M), and rat (R) spleen tissue lysates with BAFF Receptor antibody at 5µg/ml.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	BlySR3, TNFRSF 13C, CD268
Application	IHC, WB
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rabbit
Immunogen	Synthetic peptide corresponding to the sequence near the C-terminus of human BAFF receptor.
Purity Detail	Peptide affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (5µg/ml)Western Blot (2-5µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human, Mouse, Rat
UniProt ID	Q96RJ3
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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@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)