## 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-кB 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.

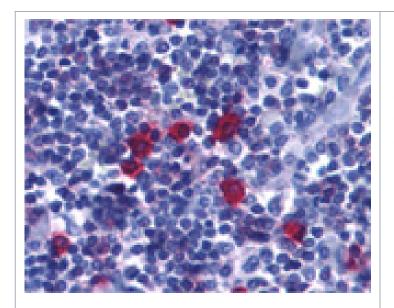
## **Ordering Information**

Order Online »

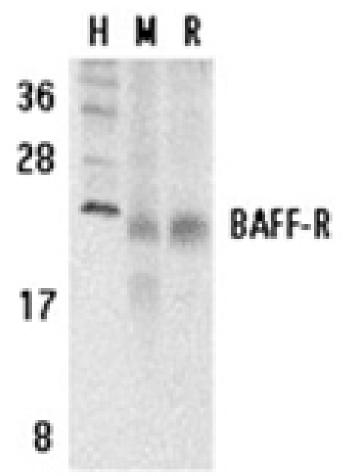
**ADI-905-305-100** 100μg

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  $5\mu$ g/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

