

BCMA polyclonal antibody

Members of the TNF superfamily regulate immune responses and induce apoptosis. Two members of this family were identified and designated BAFF and April, respectively. Three receptors for BAFF were identified; these include: BCMA (B-cell maturation protein), TACI, and BAFF-R. BCMA, like TACI, binds BAFF and April. BAFF and its receptors induce B-Cell development and survival, activate NF- κ B and c-jun N-terminal kinase, and are involved in B-Cell associated autoimmune diseases.

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

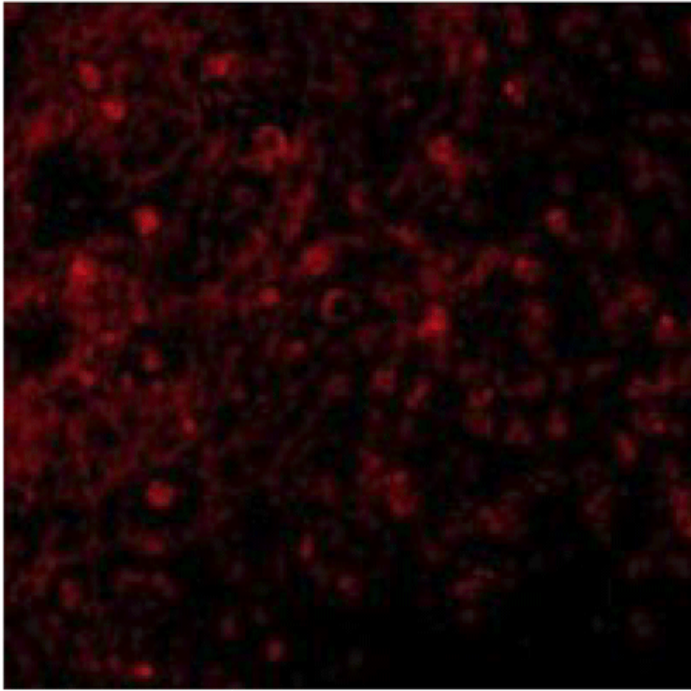
Ordering Information

[Order Online »](#)

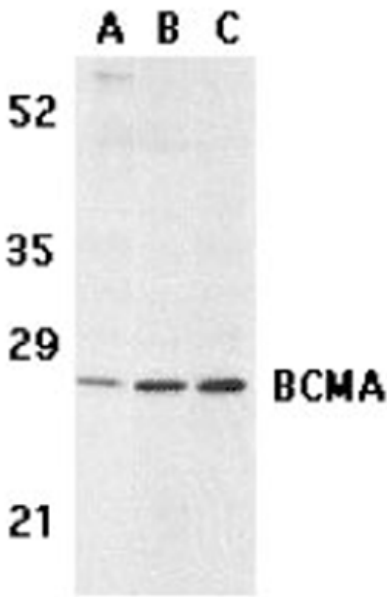
ADI-905-306-100	100 μ g
-----------------	-------------

Manuals, SDS & CofA

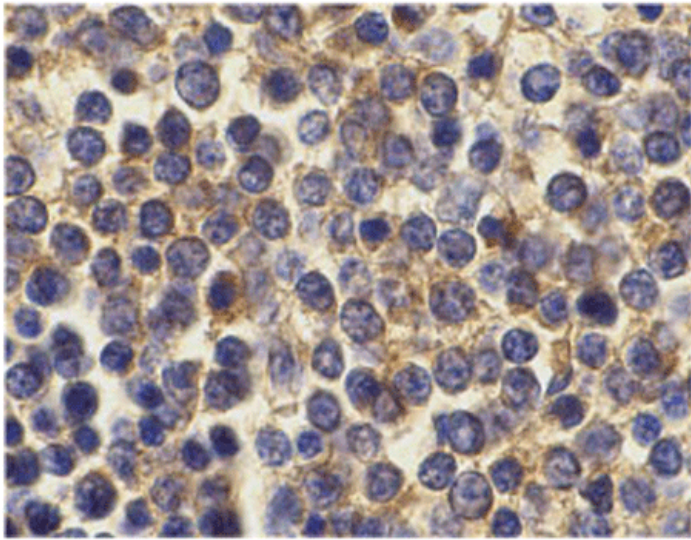
[View Online »](#)



Immunofluorescence of BCMA in human spleen cells with BCMA antibody at 10µg/ml.



Western blot analysis of BCMA in human spleen (A) tissue lysate, K562 (B), and U937 (C) cell lysates with BCMA antibody at 5µg/ml.



Immunohistochemistry analysis of BCMA in human spleen cells with BCMA antibody at 10µg/ml.

Handling & Storage

Use/Stability Stable at -20°C up to 1 year, at 4°C up to 3 months.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name B cell maturation protein, TNFRSF 17, CD269

Application ELISA, IF, IHC (PS), WB

Application Notes Detects a band of ~25kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to the sequence near the C-terminus of human BCMA.

Purity Detail Peptide affinity purified.

Recommendation Immunofluorescence (20µg/ml) Immunohistochemistry (paraffin sections,
Dilutions/Conditions 10µg/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

UniProt ID Q02223

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



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-
fr@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)