

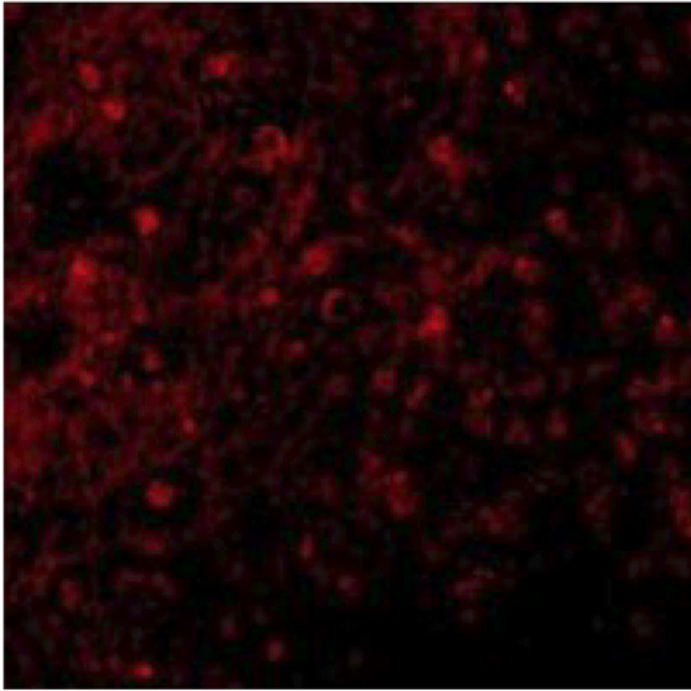
# BCMA polyclonal antibody

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- $\kappa$ B and c-jun N-terminal kinase, and are involved in B-Cell associated autoimmune diseases.

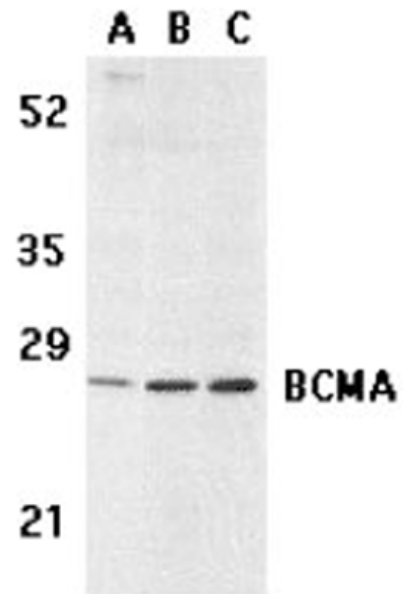
This antibody is covered by our [Worry-Free Guarantee](#).

**Manuals, SDS & CofA**

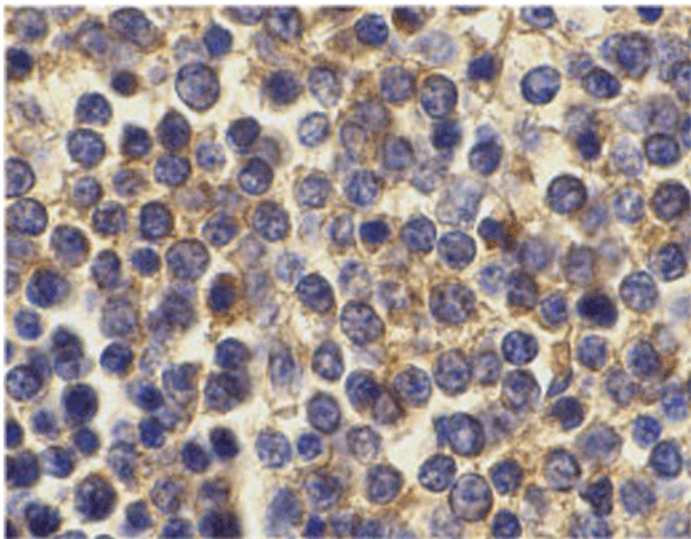
[View Online »](#)



Immunofluorescence of BCMA in human spleen cells with BCMA antibody at 10 $\mu$ g/ml.



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



Immunohistochemistry analysis of BCMA in human spleen cells with BCMA antibody at 10 $\mu$ g/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 Dilutions/Conditions** Immunofluorescence (20µg/ml) Immunohistochemistry (paraffin sections, 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: 800.948.0120  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)

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

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

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)