

BCMA (human) monoclonal antibody (Vicky-1) (biotin conjugate)

BCMA (B cell maturation) is a nonglycosylated integral membrane type I protein that is preferentially expressed in mature B lymphocytes and belongs structurally to the TNF receptor family. BCMA associates with BAFF, APRIL, TNF receptor-associated factor (TRAF) 1, TRAF2, and TRAF3 and activates NF- κ B, Elk-1, c-Jun N-terminal kinase (JNK), and p38 mitogen-activated protein kinase. BCMA plays an important role in B cell development, function and regulation.

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

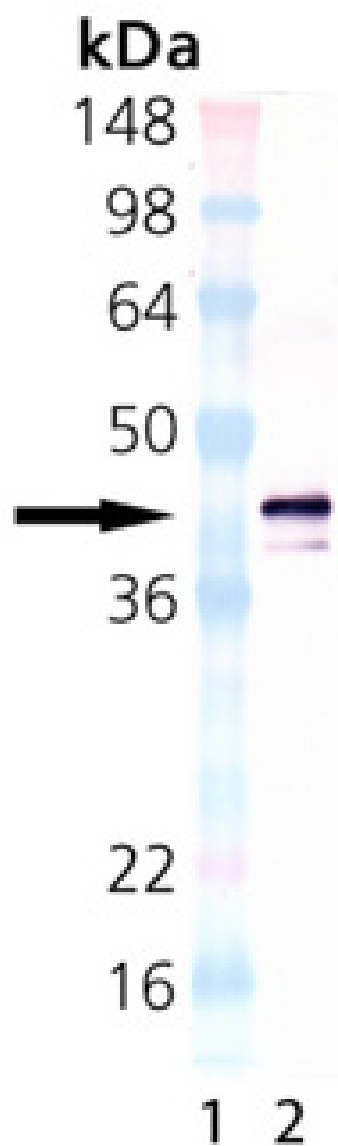
Ordering Information

[Order Online »](#)

ALX-804-151B-C100	100 μ g
-------------------	-------------

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of BCMA: Lane 1: MWM, Lane 2: BCMA (human):Fc (human), (recombinant) (Prod. No. ALX-522-026)

Handling & Storage

Handling	Avoid freeze/thaw cycles.
Short Term Storage	+4°C
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, WB
Clone	Vicky-1
Crossreactivity	Does not cross-react with mouse BCMA.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human BCMA extracellular domain (aa 2-54) (Prod. No. ALX-522-026).
Isotype	IgG1
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	ELISA (1:1,000)Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from ascites.
Species Reactivity	Human
UniProt ID	Q02223

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-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)