

BCMA (mouse) monoclonal antibody (Vicky-2)

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

Citations: 3

[View Online »](#)

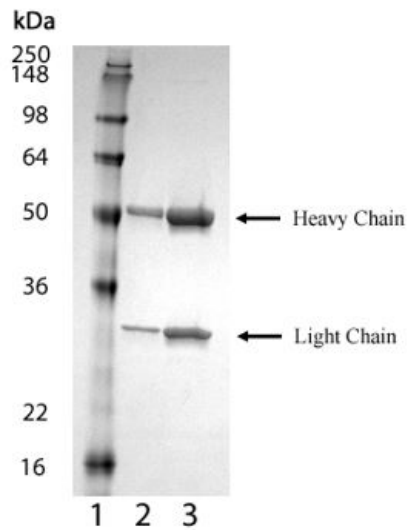
Ordering Information

[Order Online »](#)

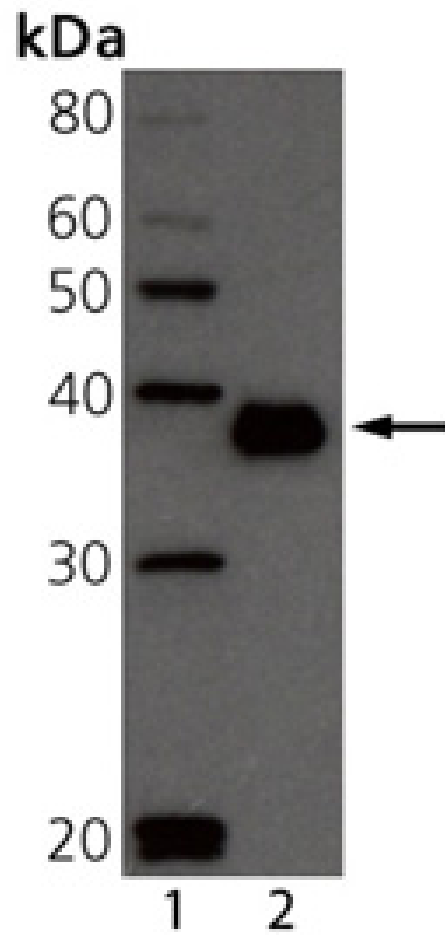
ALX-804-826-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis of Prod. No. ALX-804-826: Lane 1: MW marker, Lane 2: 1 μ g, Lane 3: 2 μ g BCMA (mouse) monoclonal antibody (Vicky-2), integrity of antibody is shown as both light chain and heavy chain are present.



Western blot analysis of BCMA: Lane 1: MW marker, Lane 2: Mouse BCMA Recombinant Protein.

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored at -20°C.
Handling	For long term storage, aliquot and freeze at -20°C. Avoid freeze/thaw cycles.
Short Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	B cell maturation protein, TNFRSF 17, CD269
Application	Flow Cytometry, IP, WB
Clone	Vicky-2
Crossreactivity	Does not cross-react with human BCMA.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse BCMA extracellular domain (aa 1-46).
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunoprecipitation (1:200)Flow Cytometry (1:1,000)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Mouse



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
088472 ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Worry-free Guarantee

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

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.