

APRIL (human) monoclonal antibody (Aprily-5) (biotin conjugate)

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

Ordering Information

[Order Online »](#)

ALX-804-801B-C100

Biotin - 100µg

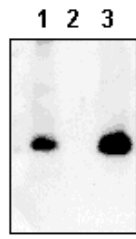
Manuals, SDS & CofA

[View Online »](#)

A.



B.



Handling & Storage

Short Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	A-Proliferation-inducing ligand, TNFSF 13, CD256
Application	ELISA, IP, WB
Application Notes	Excellent for Western Blot.
Clone	Aprily-5
Crossreactivity	Does not cross-react with mouse APRIL.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human APRIL (aa 93-233).
Isotype	IgG1
Recommendation Dilutions/Conditions	ELISA (detection)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	Human
Specificity	Recognizes full-length and cleaved APRIL.
UniProt ID	O75888
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)