

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

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

Ordering Information

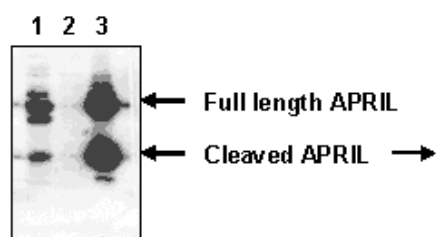
[Order Online »](#)

ALX-804-801B-C100	Biotin - 100µg
-------------------	----------------

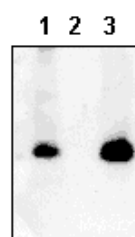
Manuals, SDS & CofA

[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 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



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)

This antibody is covered by our Worry-Free Guarantee.