

APRIL (human) monoclonal antibody (Aprily-1)

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

Citations: 1

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-148-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

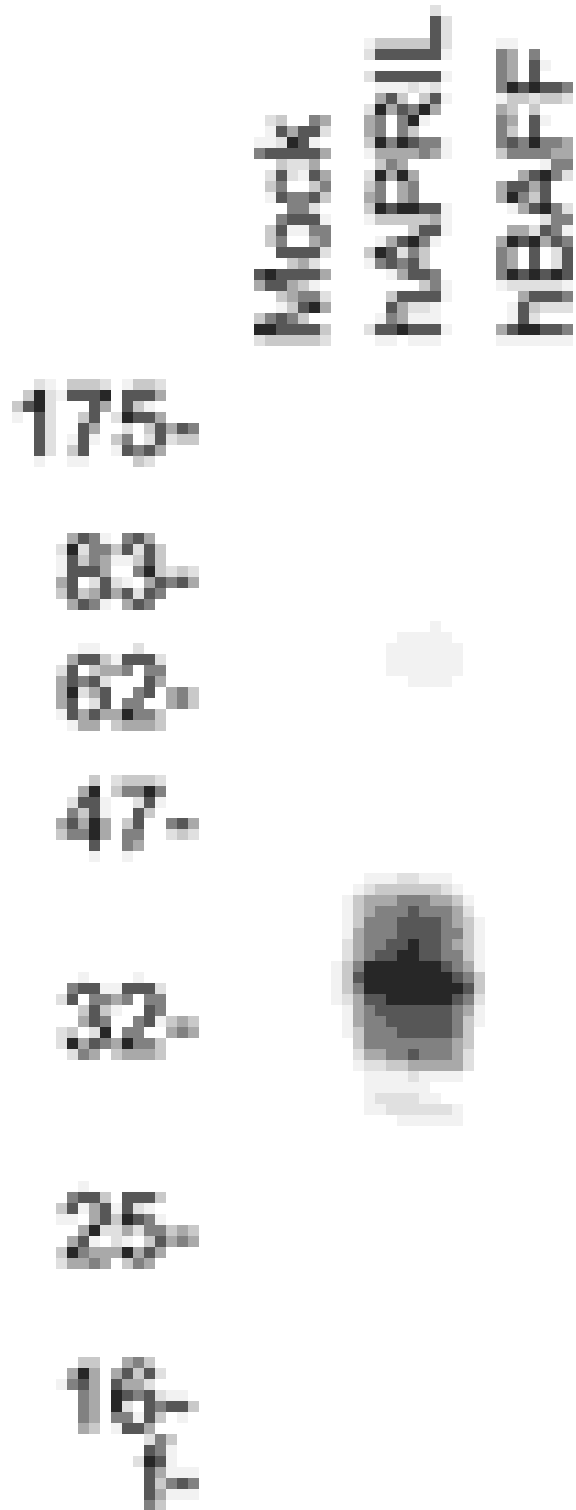


Figure 1: Western blot of total cell extracts from HEK 293 cells transfected with the indicated expression vector. Aprily-1 reacts specifically with human APRIL.
Method: 10µg of protein was applied to the gel. Revealed with Aprily-1 (1µg/ml) and HRP-coupled anti-mouse secondary antibody (1:4'000).
Note: The human APRIL construct used in this experiment is an uncleavable fusion protein between human BAFF (aa 1-132) and human APRIL (aa 93-233).

Handling & Storage

Use/Stability	Stable for at least 1 year after receipt when stored as recommended.
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	A-Proliferation-inducing ligand, TNFSF 13, CD256
Application	Flow Cytometry, IP, WB
Application Notes	Flow Cytometry: also works for intracellular FC staining. Excellent for Western Blot.
Clone	Aprily-1
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
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human

UniProt ID

075888



Worry-free Guarantee

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

European Sales Office
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

This antibody is covered by our **Worry-Free Guarantee**.