

APRIL (human) monoclonal antibody (Aprily-8)

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

Citations: 4

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-804-149-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)

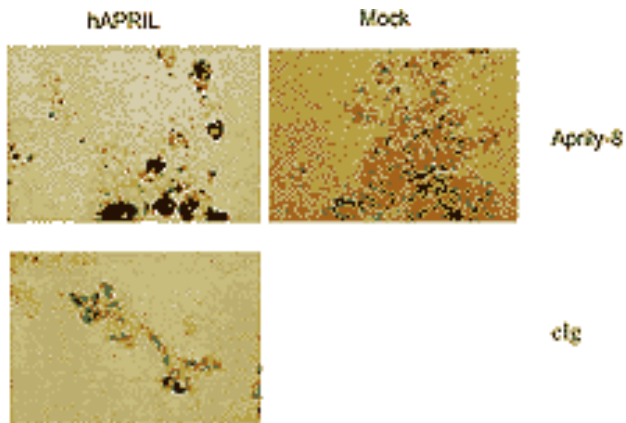


Figure 3: Immunostaining of HEK 293 cells transfected with a human APRIL expression plasmid (left panels), or mock transfected (right panel) by Aprily-8. **Method:** 3 days after transfection of cells with the indicated constructs, cells were fixed with 4% formaldehyde 5 min. at RT. After a wash in PBS, samples were dehydrated by washes in 60%, 80%, 90%, 100% EtOH and xylol. Cells were then dried and embedded in paraffin. Sections were cut, mounted on slides and dried overnight at 50°C. Slides were then successively washed 2x 10 min. in xylol, 2x 10 min. in 100% ethanol, and then treated 10 min. in 100% methanol/0.6% H₂O₂ to inhibit endogenous peroxidase. Samples were rehydrated by washes in 90%, 80%, 60% ethanol and PBS. After micro-wave treatment, slides were washed 3x in PBS, blocked with IgG, and incubated for 1 hour with 5µg/ml Aprily-8 or control mouse IgG (isotype control) in 1%BSA / 1x PBS for 1 hour. After PBS washes, samples were incubated with the secondary Ab for 1 hour, washed in PBS and revealed with StreptABComplex/HRP (Vector) and AEC.

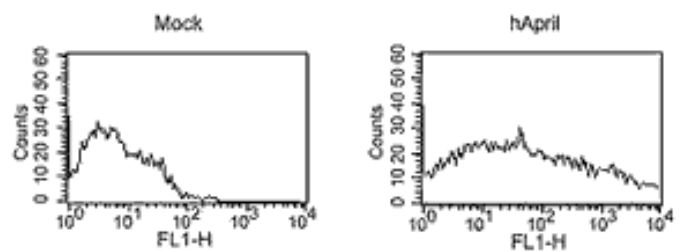


Figure 2: FACS analysis of cells with MAbs to APRIL (Aprily-8). **Method:** HEK 293 cells were mock transfected or transfected with an expression plasmid coding for a non-cleavable human APRIL. Cells (5×10^5) were incubated on ice for 30 min. in 50µl FACS buffer (PBS, 5% fetal calf serum, 0.02% azide) containing 10µg/ml of Aprily-8 antibody. After washing in FACS buffer, FITC-conjugated antibody to mouse IgG was added. Cells were incubated on ice for 30 min., washed and analyzed by flow cytometry.

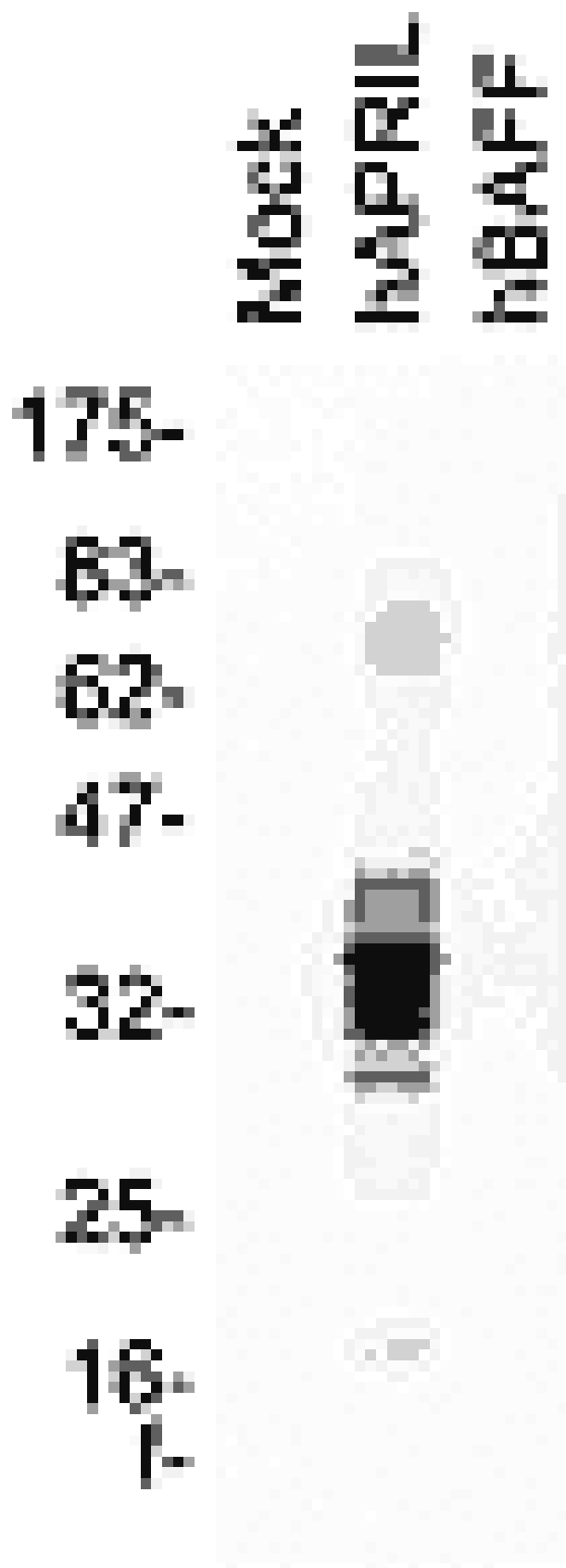


Figure 1: Western blot of total cell extracts from HEK 293 cells transfected with the indicated expression vector. Aprily-8 reacts specifically with human APRIL.
Method: 10µg of protein was applied to the gel. Revealed with Aprily-8 (1µg/ml) and HRP-coupled anti-mouse secondary antibody (1:4'000).
Note: 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 one year after receipt when stored at -20°C.
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, ICC, IHC (PS), IP, WB
Application Notes	Excellent for Immunocytochemistry and Western Blot.
Clone	Aprily-8
Crossreactivity	Does not cross-react with mouse APRIL or TWE-PRIL.
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 supernatant.
Species Reactivity	Human
Specificity	Recognizes APRIL and TWE-PRIL.
UniProt ID	O75888

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)