

Apaf-1 monoclonal antibody (18H2)

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

Citations: 4

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

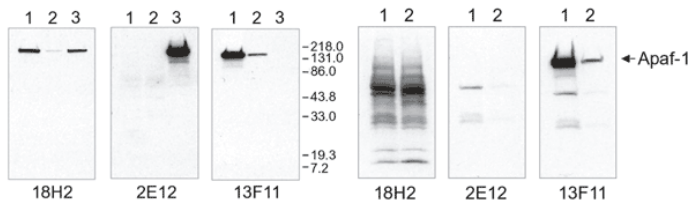
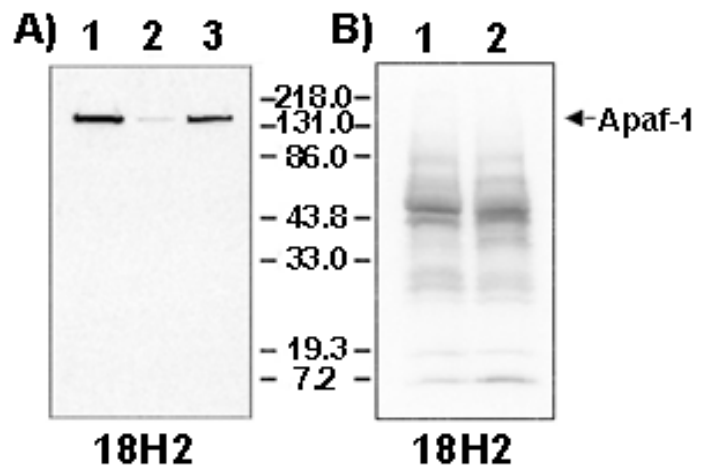


Figure: Species cross-reactivity of MABs to APAF-1 by WB. **left:** Lysates from 1-day-old mouse brain (lane 1), adult mouse brain (lane 2) and SH-SY5Y human neuroblastoma cell line (lane 3). **right:** Lysates from 2-days-old rat brain (lane 1) and adult rat brain (lane 2).



Figures: Species cross-reactivity of MAb to Apaf-1 (clone 18H2) (ALX-804-350) by Western blot. **A:** Lysates from 1-day-old **mouse** brain (lane 1), adult **mouse** brain (lane 2) and SH-SY5Y **human** neuroblastoma cell line (lane 3) were probed with clone 18H2. **B:** Lysates from 2-days-old **rat** brain (lane 1) and adult **rat** brain (lane 2) were probed with clone 18H2. **Results:** 18H2 is human and mouse specific but does not recognize rat Apaf-1.

Handling & Storage

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	Apoptosis protease-activating factor 1
Application	ELISA, ICC, IP, WB
Clone	18H2
Crossreactivity	Does not cross-react with rat Apaf-1.
Formulation	Liquid. Protein G purified antibody in PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant mouse Apaf-1 (apoptosis protease-activating factor-1) (aa 1-97) containing the N-terminal CARD domain.
Isotype	IgG2a
Species Reactivity	Human, Mouse
Specificity	Recognizes the CARD domain of Apaf-1.
UniProt ID	O88879
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)