

Apaf-1 (human) monoclonal antibody (2E12)

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

Citations: 9

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

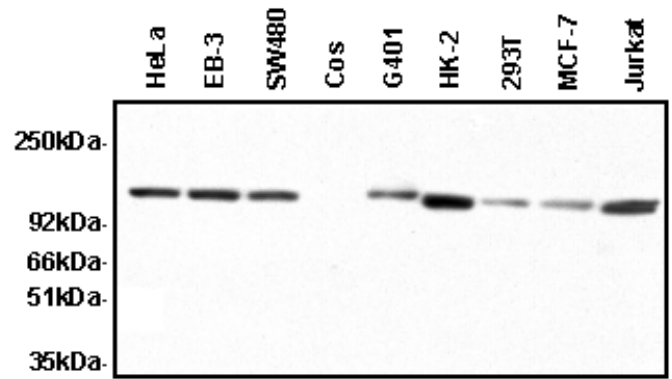
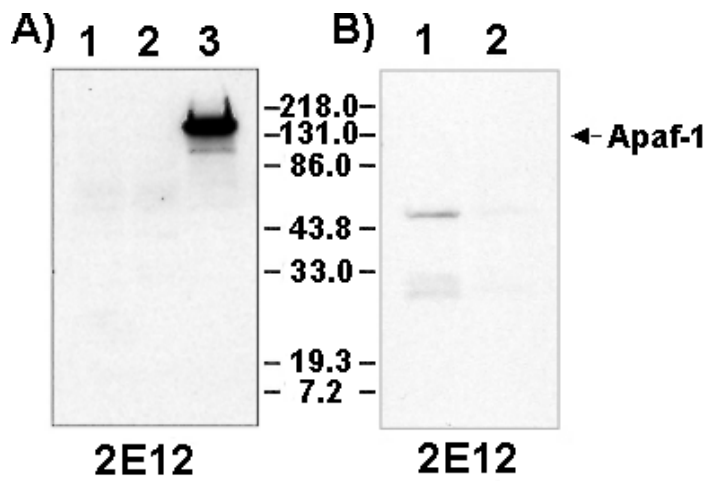


Figure: Western blot analysis with anti-Apaf-1 (human) MAb (clone 2E12) revealing endogenous Apaf-1. Lysates from 10^6 HeLa, MCF-7, 293T, COS, Jurkat, and activated primary human T cells were probed with clone 2E12. COS (monkey) cells were used as a negative control.

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, IHC (FS), IP, WB
Clone	2E12
Crossreactivity	Does not cross-react with mouse or rat Apaf-1.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Rat
Immunogen	Recombinant human Apaf-1 (apoptosis protease-activating factor 1) N-terminal CARD and CED-4 homologous domains (aa 1-464) .
Isotype	IgG2a
Purity Detail	Protein G-affinity purified.
Recommendation Dilutions/Conditions	Immunohistochemistry (EM on frozen sections; for details see G. Hausmann, et al., below.)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Human
Specificity	Recognizes the CARD domain of Apaf-1.

UniProt ID

O14727

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)