

# Apaf-1 monoclonal antibody (13F11)

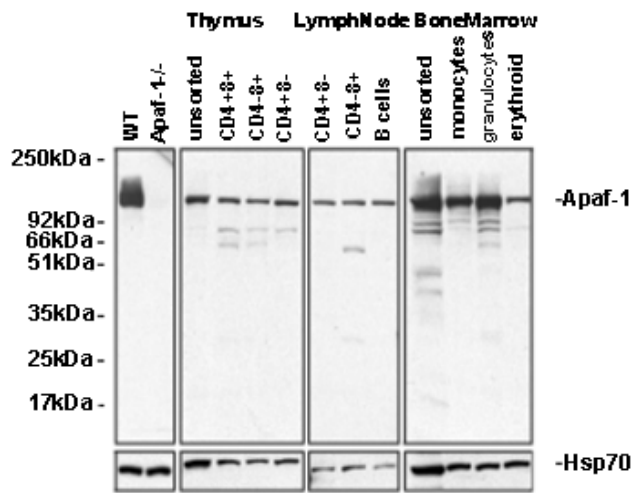
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 4

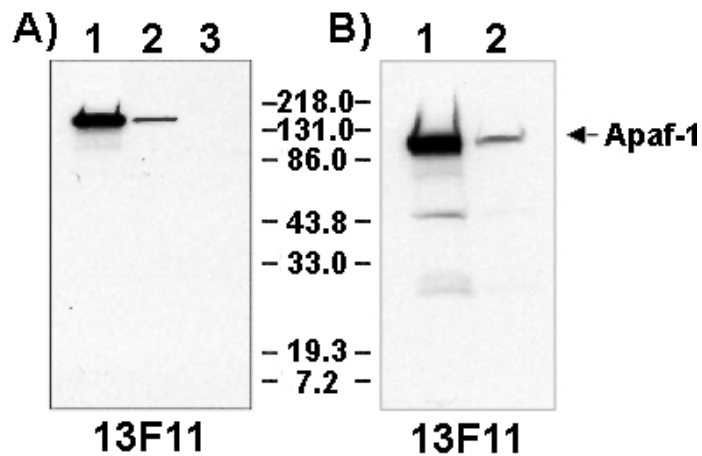
[View Online »](#)

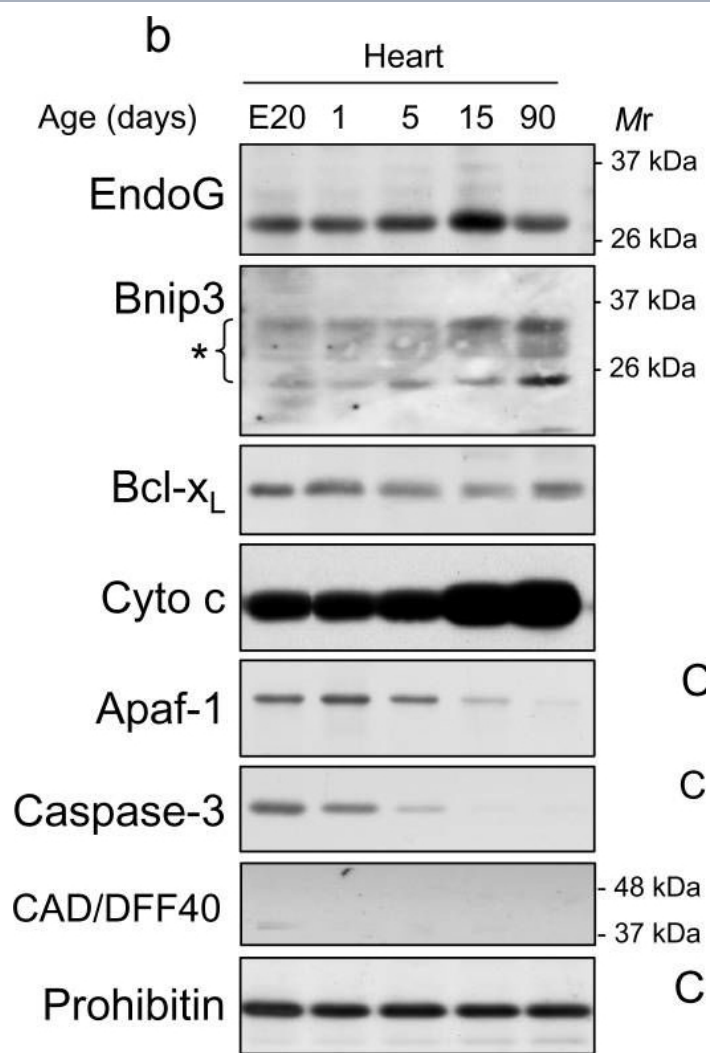
Manuals, SDS & CofA

[View Online »](#)



**Figure:** Western blot analysis with anti-Apaf-1 MAb clone 13F11 on purified cells from mouse tissues. The ~130kDa band corresponds to endogenous Apaf-1, which is absent in cells from Apaf-1- mice.





EndoG and Bnip3 are most highly expressed in rat heart, and are located to cardiac myocytes through development and during ischemia. a) Western blot analysis of EndoG, using a new anti-EndoG antibody (AntibodyBcn, BCN4778) and Bnip3 protein expression in total protein extracts from different adult rat tissues. EndoG is detected as a single band ( 27 kDa) and Bnip3 is detected as a group of bands around the 26 kDa mark (asterisk), in agreement with the information provided by the manufacturer. Both proteins are expressed mainly in the heart and skeletal muscle. Detection of EndoG and Bnip3 was performed in two different sets of independent samples with similar results. b) Western blot detection of EndoG, Bnip3 and proteins involved in caspase-dependent cell death in total protein extracts of ventricles from rats of different ages ranging from embryonic day 20 to adulthood. A representative image is shown from three independent experiments. c) Neonatal cardiomyocytes were treated with ischemia and the abundance of EndoG, Bnip3 and proteins involved in the caspase-dependent pathway was analyzed by Western blot in extracts obtained at different time points. Primary heart fibroblasts and neonatal liver samples were added for comparison. A representative image is shown from three independent experiments.



## 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	ICC, IP, WB
Clone	13F11
Crossreactivity	Does not cross-react with human Apaf-1.
Formulation	Liquid. 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
Purity Detail	Protein G-affinity purified.
Species Reactivity	Mouse, Rat
Specificity	Recognizes the CARD domain of Apaf-1.
UniProt ID	O88879
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a> .

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)