

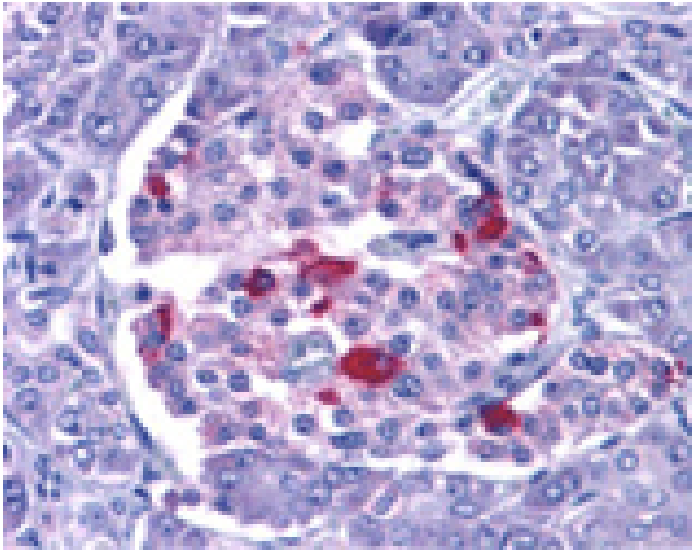
EndoG polyclonal antibody

The fragmentation of nuclear DNA is a hallmark of apoptotic cell death. EndoG is a mitochondrion specific nuclease that translocates to the nucleus and cleaves chromatin DNA during apoptosis in a caspase independent fashion.

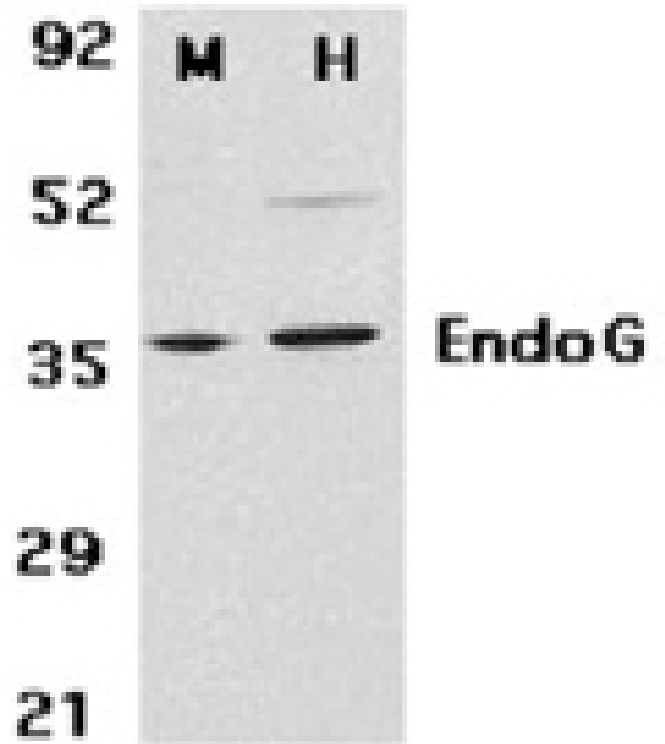
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Immunohistochemistry analysis of EndoG in human pancreas with EndoG antibody at 15 μ g/ml.



Western blot analysis of EndoG in mouse (M) 3T3 and human (H) HepG2 cell lysates with EndoG antibody at 2 μ g/ml.

Handling & Storage

Short Term Storage +4°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Endonuclease G

Application ELISA, IHC, IHC (PS), WB

Application Notes Detects a band of ~35kDa by Western blot.

Formulation Liquid. In PBS containing 0.02% sodium azide.

Host Rabbit

Immunogen Synthetic peptide corresponding to 15 amino acids near the N-terminus of human EndoG. The immunogen is located within amino acids 40 – 90 of EndoG.

Isotype IgG

Purity Detail Affinity chromatography purified via peptide column.

Recommendation Dilutions/Conditions Immunohistochemistry (2.5-15µg/ml)Western Blot (2µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human, Mouse, Rat

UniProt ID Q14249

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)