

Fis1 polyclonal antibody

The human Fis1, mitochondrial fission 1, is a novel molecule that was discovered as part of the mammalian fission machinery. Fis1 has been suggested to play a role in regulation of the fission processes that might be involved in apoptotic mechanisms.

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

Citations: 16

[View Online »](#)

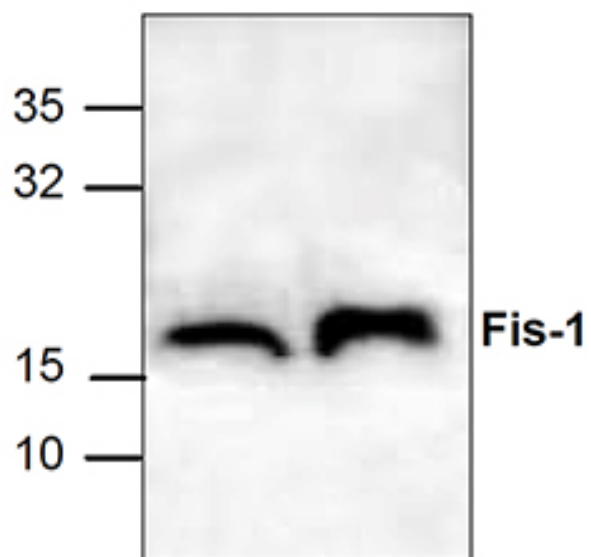
Ordering Information

[Order Online »](#)

ALX-210-1037-0100	100µg
-------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of Fis1 pAb (Prod. No. ALX-210-1037) with Jurkat cell lysates.

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Mitochondrial fission 1
Application	WB
Formulation	Liquid. In PBS containig 0.5% BSA, 0.01% thimerosal and 30% glycerol.
Host	Rabbit
Immunogen	Synthetic peptide surrounding aa 143 of human Fis1.
Isotype	IgG
Purity Detail	Affinity purified.
Recommendation Dilutions/Conditions	Western Blot (1-4 µg/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	Q9Y3D6
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-be@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-fr@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)