

MFN2 polyclonal antibody

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

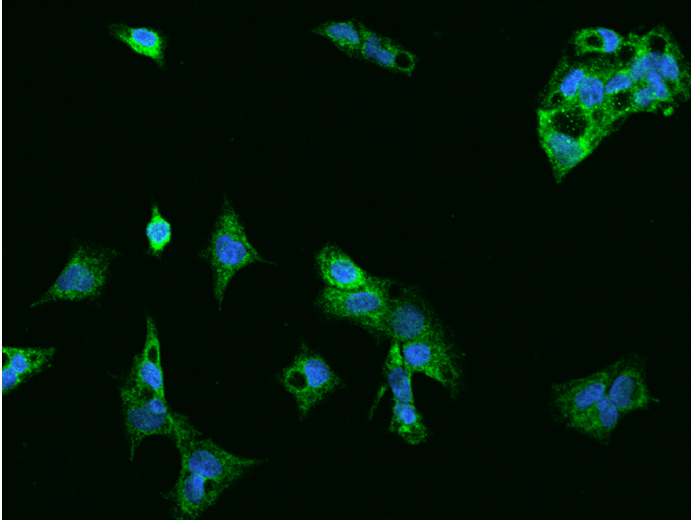
Ordering Information

[Order Online »](#)

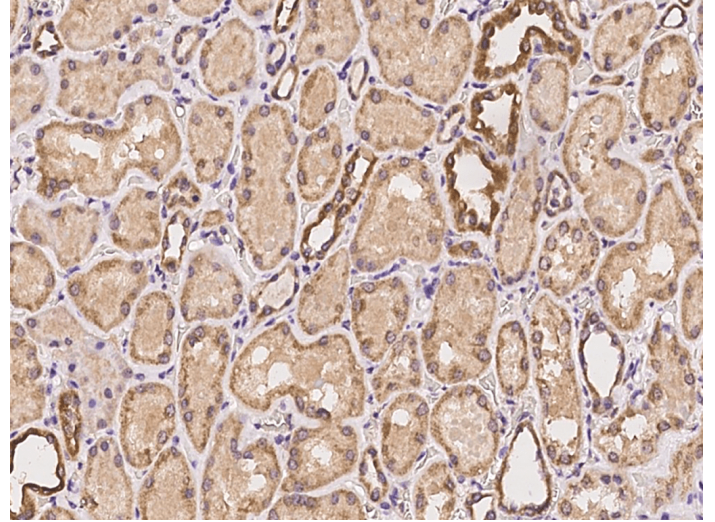
ENZ-ABS1002-0100	100µl
------------------	-------

Manuals, SDS & CofA

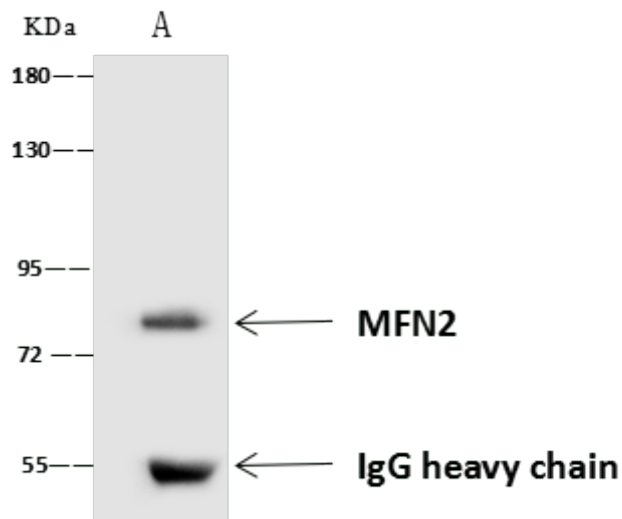
[View Online »](#)



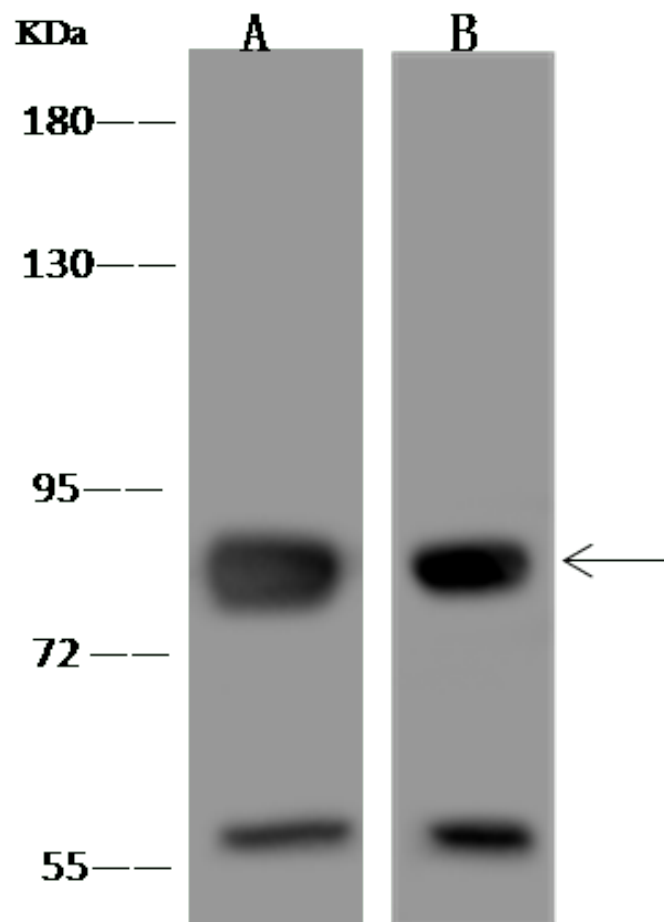
Immunofluorescence staining of MFN2 in HepG2 cells. Cells were fixed with 4% PFA, permeabilized with 0.1% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-Human MFN2 polyclonal antibody (dilution ratio 1:200) at 4 °C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to Cytoplasm.



Immunochemical staining of human MFN2 in human kidney with rabbit polyclonal antibody at 1:100 dilution, formalin-fixed paraffin embedded sections.



MFN2 was immunoprecipitated using: Lane A: 0.5 mg HeLa Jurkat Whole Cell Lysate, 4 μ L anti-MFN2 rabbit polyclonal antibody and 60 μ g of Immunomagnetic beads Protein A/G. Primary antibody: Anti-MFN2 rabbit polyclonal antibody, at 1:100 dilution. Secondary antibody: Goat Anti-Rabbit IgG (H+L)/HRP at 1/10000 dilution. Developed using the ECL technique. Performed under reducing conditions. Observed band size: 86 kDa.



Anti-MFN2 rabbit polyclonal antibody at 1:500. Lane A: Jurkat Whole Cell Lysate. Lysates/proteins at 30 μ g per lane. Secondary Goat Anti-Rabbit IgG (H+L)/HRP at 1/10000 dilution. Developed using the ECL technique. Performed under reducing conditions. Observed band size: 86 kDa.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Mitofusin 2

Application ICC, IF, IHC (PS), IP, WB

Formulation Liquid. In PBS, pH7.0 with 0.03% Proclin300. Preservative free.

Host Rabbit

Immunogen E. coli-derived human MFN2 fragment

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:1'000)Immunohistochemistry (paraffin sections) (1:50-1:200)Immunocytochemistry (1:100-1:500)Immunofluorescence (1:100-1:500)Immunoprecipitation (1-5 µl/ml of lysate)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Human

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



ENZO LIFE SCIENCES,
INC.
100 Morris Avenue
Rte. 100, Suite 300
Rutherford, NJ 07070
info-
usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(EU) AG
Phone: +41 61 926 8989
info-
eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-
be@enzolifesciences.com

France
Phone: +33 472 440 655
info-
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-
de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-
uk@enzolifesciences.com