

mtHSP70 monoclonal antibody (JG1)

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

Citations: 15

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)

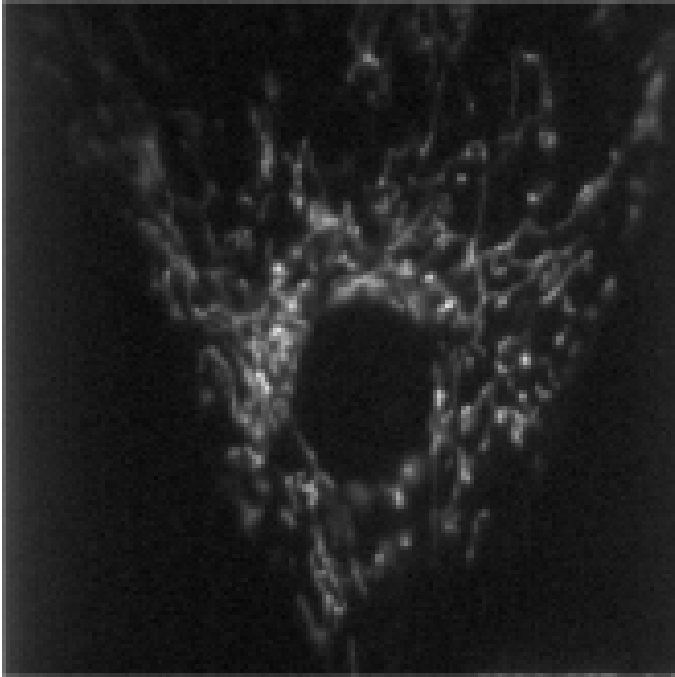


Figure: Immunolocalization of mtHSP70 in human fibroblasts using MAb to to mtHSP70 (JG1) (Prod. No. [ALX-804-077](#)).^{""},mtHSP70 monoclonal antibody (JG1) ICC,mtHSP70 monoclonal antibody (JG1)

Immunocytochemistry,yes

ALX-804-095,804-

095.gif,[http://static.enzolifesciences.com/fileadmin/files/image/804-](http://static.enzolifesciences.com/fileadmin/files/image/804-095.gif)

095.gif,**Figure:** Immunofluorescence staining of TGN38 (red) with MAb to trans-Golgi Network 38 (2F7.1) (Prod. No. [ALX-804-095](#)) and β -COP (green) with PAb to β -COP (Pro. No. ALX-210-127) in NRK cells.^{""""}

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Mortalin, Heat shock protein 70 (mitochondrial), GRP75

Application ICC, IP, WB

Application Notes Detects a band of ~75kDa by Western blot.

Clone JG1

Formulation Liquid. In PBS with 1 mg/mL BSA containing 0.05% sodium azide.

Host Mouse

Immunogen Synthetic peptide corresponding to C-terminal aa 615-633 (G615SGSSGTGEQKEDQKEEKQ633) of mouse mtHSP70 (mitochondrial heat shock protein 70).

Isotype IgG3

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunocytochemistry (1:50)Western Blot (1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Species Reactivity Dog, Human, Monkey, Mouse

UniProt ID P38647

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)