

HSP70 recombinant monoclonal antibody (001)

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

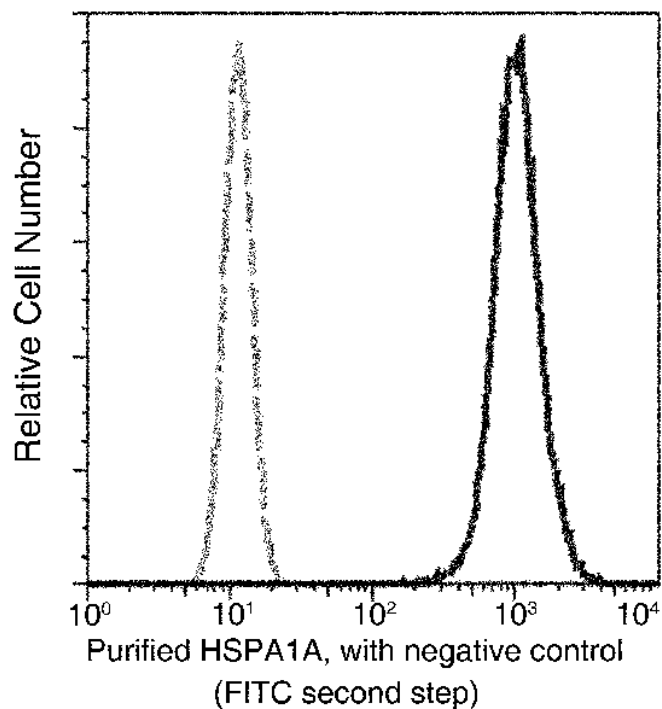
Ordering Information

[Order Online »](#)

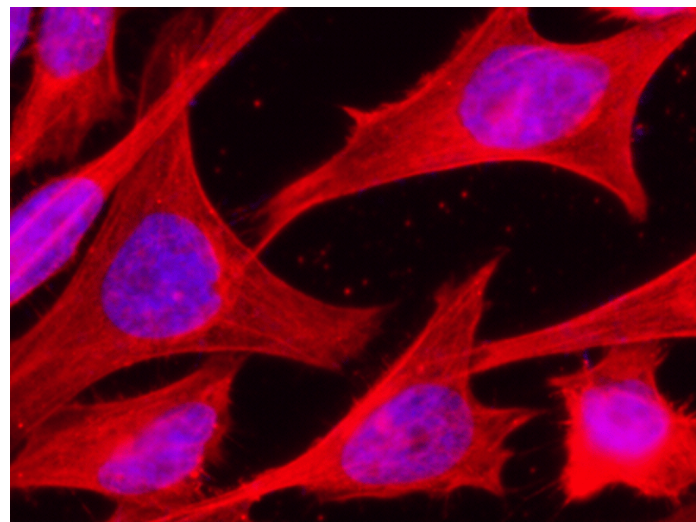
ENZ-ABS982-0100	100µl
-----------------	-------

Manuals, SDS & CofA

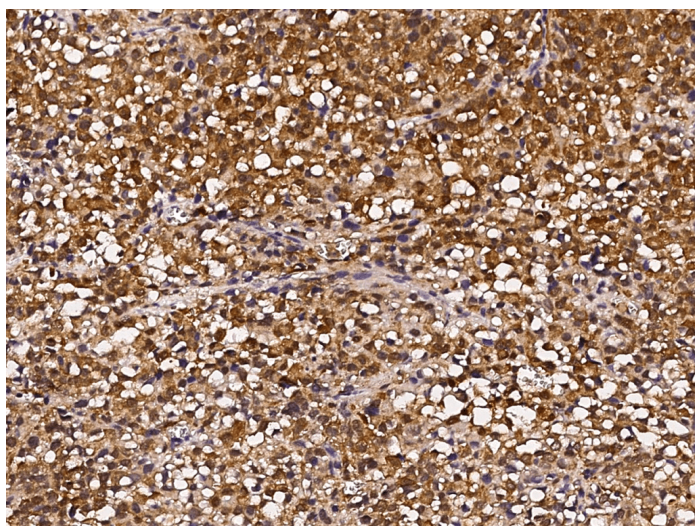
[View Online »](#)



Flow cytometric analysis of Human HSPA1A expression on HeLa cells. The cells were treated and stained with purified anti-Human HSPA1A, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.

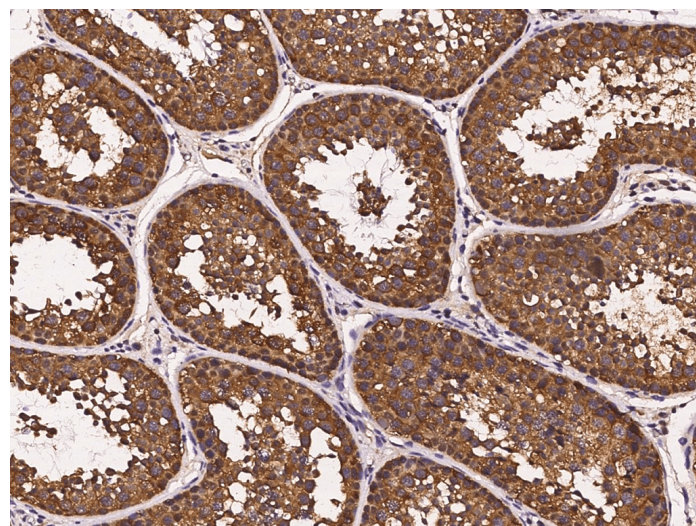


Immunofluorescence staining of Human HSPA1A in HeLa cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with rabbit anti-Human HSPA1A monoclonal antibody (1:60) at 37 °C for 1 hour. Then cells were stained with the Alexa Fluor® 594-conjugated goat Anti-rabbit IgG secondary antibody (red) and counterstained with DAPI (blue). Positive staining was localized to cytoplasm.

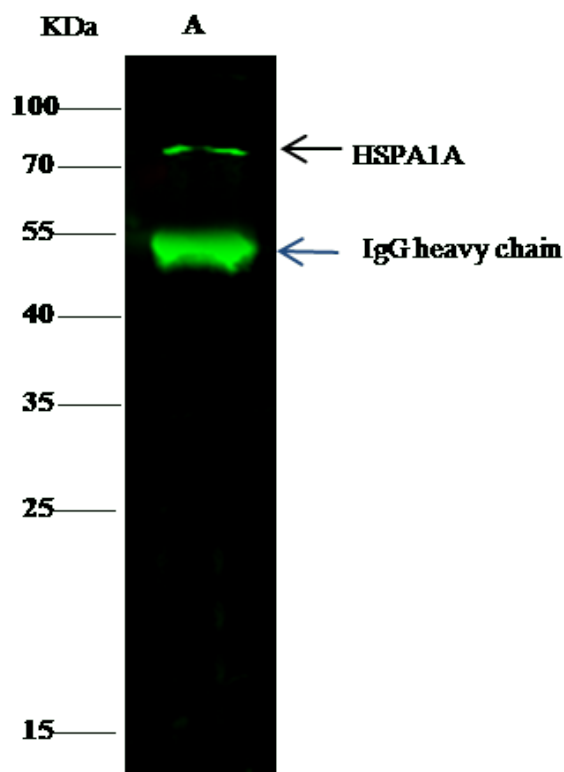


© Sino Biological Inc. All rights reserved

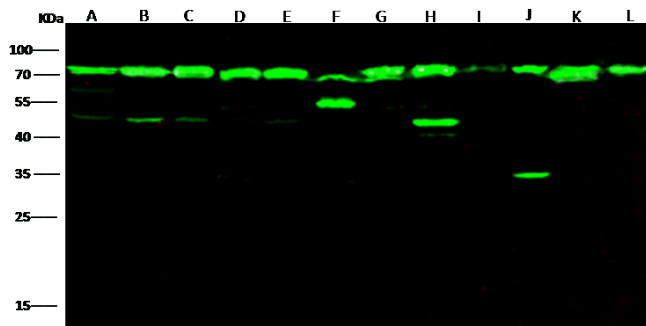
Immunochemical staining of HSPA1A in human hepatoma with rabbit monoclonal antibody (1:1000, formalin-fixed paraffin embedded sections).



Immunochemical staining of HSPA1A in cynomolgus macaque testis with rabbit monoclonal antibody (1:1000, formalin-fixed paraffin embedded sections).



HSPA1A was immunoprecipitated using: Lane A: 0.5 mg HeLa Whole Cell Lysate, 2 μ L anti-HSPA1A rabbit monoclonal antibody and 15 μ L of 50 % Protein G agarose. Primary antibody: Anti-HSPA1A rabbit monoclonal antibody, at 1:200 dilution. Secondary antibody: Dylight 800-labeled antibody to rabbit IgG (H+L), at 1:5000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size: 70 kDa.



Anti-HSPA1A rabbit monoclonal antibody at 1:500 dilution. Lane A: A549 whole cell lysate. Lane B: HCT116 whole cell lysate. Lane C: HeLa whole cell lysate. Lane D: HepG2 whole cell lysate. Lane E: HL60 whole cell lysate. Lane F: A431 whole cell lysate. Lane G: MCF7 whole cell lysate. Lane H: K562 whole cell lysate. Lane I: C6 whole cell lysate. Lane J: 293T whole cell lysate. Lane K: MOLT4 whole cell lysate. Lane L: Jurkat whole cell lysate. Lysates/proteins at 30 μ g per lane. Secondary Goat Anti-Rabbit IgG H&L (Dylight800) at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size: 70 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 HSPA1A, Heat shock protein 70

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

Clone 001

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Host Rabbit

Immunogen Recombinant human HSP70 / HSPA1A protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:1'000)ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:500-1:2'000)Immunocytochemistry (1:20-1:100)Immunofluorescence (1:20-1:100)Immunoprecipitation (1-4µl/mg 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.
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-
fr@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)

Last modified: September 17, 2024

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.