

HSC70/HSP73

monoclonal antibody

(1B5) (PE conjugate)

The Hsp70 family of heat shock proteins contains multiple homologs ranging in size from 66-78 kDa, and are the eukaryotic equivalents of the bacterial DnaK. The most studied Hsp70 members include the cytosolic stress-induced Hsp70 (Hsp72), the constitutive cytosolic Hsc70 (Hsp73), and the ER-localized BiP (Grp78). Hsp70 family members contain highly conserved N-terminal ATPase and C-terminal protein binding domains. Binding of peptide to Hsp70 is assisted by Hsp40, and stimulates the inherent ATPase activity of Hsp70, facilitating ATP hydrolysis and enhanced peptide binding. Hsp70 nucleotide exchange and substrate binding coordinates the folding of newly synthesized proteins, the re-folding of misfolded or denatured proteins, coordinates trafficking of proteins across cellular membranes, inhibits protein aggregation, and targets the degradation of proteins via the proteasomal pathway.

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

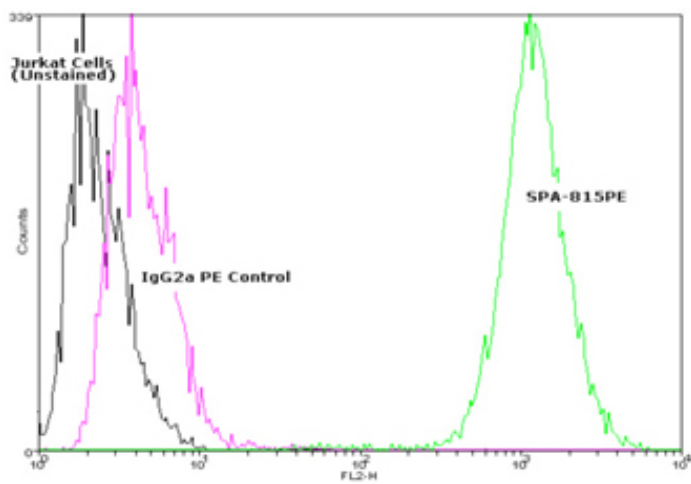
Ordering Information

[Order Online »](#)

ADI-SPA-815PE-E	100µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of 10⁶ Jurkat cells stained using HSC70 (HSP73) mAb (1B5), R-Phycoerythrin Conjugate at a concentration of 10µg/ml

Handling & Storage

Handling Avoid vigorous centrifugation and vortexing. Do not freeze. Protect from light.

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Heat shock protein 70, Hsc70, Hsp73

Application Flow Cytometry

Application Notes Detects a band of ~73kDa by Western blot.

Clone 1B5

Formulation Liquid. In PBS, pH 7.2, containing 0.09% sodium azide.

GenBank ID M34561

Host Rat

Immunogen Native hamster Hsc70 (Hsp73).

Isotype IgG2a

Purity Detail Protein G affinity purified.

Recommendation Flow Cytometry (10µg/ml) Suggested dilutions/conditions may not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from ascites.

Species Reactivity Bovine, Chicken, Dog, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep



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)

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee.

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.