

# HSC70/HSP70 monoclonal antibody (N27F3-4) (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](#).

Citations: 1

[View Online »](#)

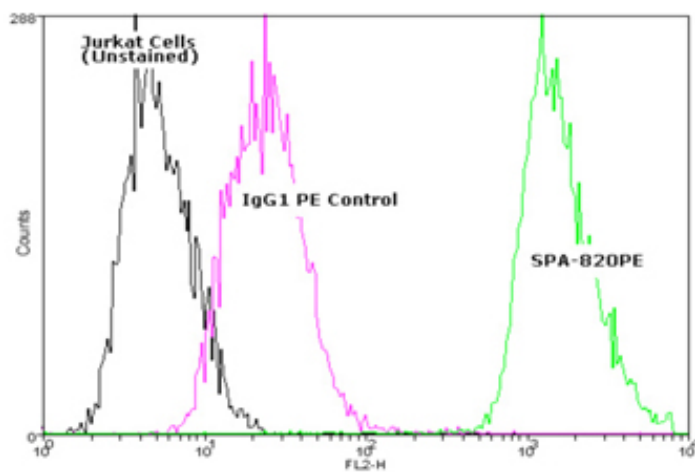
## Ordering Information

[Order Online »](#)

ADI-SPA-820PE-D	50µg
ADI-SPA-820PE-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of  $10^6$  Jurkat cells stained using HSP70/HSC70 mAb (N27F34), R-Phycoerythrin Conjugate at a concentration of  $10\mu\text{g/ml}$

## Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name Hsc70/Hsp73, Hsp70/Hsp72

Application Flow Cytometry

Clone N27F3-4

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

GenBank ID M11717 (HSP70), Y00371 (HSC70)

Host Mouse

Immunogen Purified Hsc70 (Hsp73) and Hsp70 (Hsp72) isolated from human HeLa cells.

Isotype IgG1

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 mouse ascites.

**Species Reactivity** Beluga, Bovine, Chicken, Dog, Fish, Guinea pig, Hamster, Human, Monkey, Mouse, Plant, Porcine, Rabbit, Rat, Sheep, Xenopus

**UniProt ID** P0DMV8 (HSP70/HSP72), P11142 (HSC70/HSP73)

**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)