

# HSP70/HSP72 monoclonal antibody (C92F3A-5) (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 ATP-ase 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: 2

[View Online »](#)

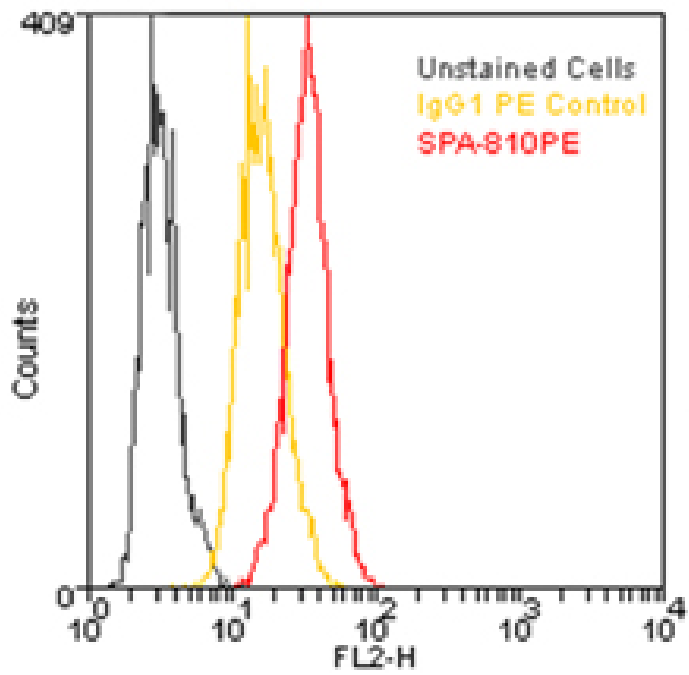
## Ordering Information

[Order Online »](#)

|                 |       |
|-----------------|-------|
| ADI-SPA-810PE-D | 50µg  |
| ADI-SPA-810PE-F | 200µg |

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometry analysis of  $10^6$  Jurkat cells stained using HSP70 (HSP72) mAb (C92F3A-5), R-Phycoerythrin Conjugate at a concentration of  $10\mu\text{g/ml}$ .

# 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                    | Hsp70, HspA1A, Heat shock protein 70, HspA1B, Hsp72                                                                                                                     |
| Application                         | Flow Cytometry                                                                                                                                                          |
| Application Notes                   | Detects a band of ~70kDa by Western blot.                                                                                                                               |
| Clone                               | C92F3A-5                                                                                                                                                                |
| Formulation                         | Liquid. In PBS containing 0.09% sodium azide.                                                                                                                           |
| GenBank ID                          | M11717                                                                                                                                                                  |
| Gene/Protein Identifier             | NP_005336.3 (RefSeq), NM_005345 (RefSeq), 3303 (Entrez GeneID), 140550 (OMIM)                                                                                           |
| Host                                | Mouse                                                                                                                                                                   |
| Immunogen                           | Native human Hsp70.                                                                                                                                                     |
| Isotype                             | IgG1                                                                                                                                                                    |
| Purity Detail                       | Protein G affinity purified.                                                                                                                                            |
| Recommendation Dilutions/Conditions | Flow Cytometry (1:100)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

Beluga, Bovine, Chicken, Dog, Drosophila, Fish, Guinea pig, Hamster, Human, Monkey, Mouse, Porcine, Rabbit, Rat, Sheep

## UniProt ID

P0DMV8 (HSPA1A), P0DMV9 (HSPA1B)

## 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)