HSP70/HSP72 monoclonal antibody (C92F3A-5) (DyLight™ 488 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: 1

View Online »

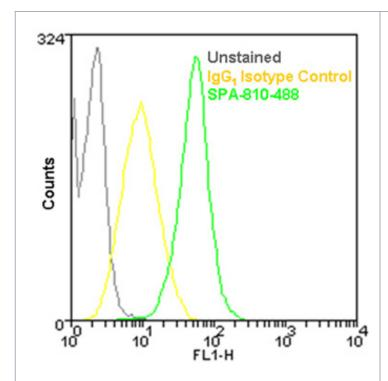
Ordering Information

Order Online »

ADI-SPA-810-488-D	50µg
ADI-SPA-810-488-F	200μg

Manuals, SDS & CofA

View Online »



Flow cytometry analysis of 10⁶ Jurkat cells stained using HSP70 (HSP72) mAb (C92F3A -5), DyLight™ 488 Conjugate at a concentration of 50µg/ml

Handling & Storage

Handling Avoid freeze/thaw cycles. 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 NotesDetects 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.

lsotype lgG1

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:20)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, C. elegans, Chicken, Dog, Drosophila, Fish,

Human, Mouse, Rat

Worry-free Guarantee

This antibody is covered by our Worry-Free Guarantee

Last modified: May 29, 2024



info-

European Sales Office ENZO LIFE SCIENCES Phone: 800.942.0430 (ELS) AG Phone: +41 61 926 8989 usa@enzolifesciences.com

eu@enzolifesciences.com

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

Phone: +33 472 440 655 fr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 de@enzolifesciences.com

UK & Ireland Phone (UK customers): 0845 601 1488 Phone: +44 1392 825900 info-uk@enzolifesciences.com