

HSP60 monoclonal antibody (Mab11-13)

Hsp60 is a member of the chaperonin family of heat shock proteins, with homologs functioning in the cytosol and mitochondria to fold nascent and aggregated proteins. Hsp60 is the eukaryotic homolog of the *E. coli* GroEL protein, and forms a multimeric complex in the mitochondria with Hsp10 (Cpn10) to form a large central cavity in which ATP-dependent protein folding takes place. TRiC/CCT, a eukaryotic relative of Hsp60, is expressed in the cytosol and participates in the folding of actin and tubulin substrates, but lacks any association with an Hsp10-like co-factor.

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

Citations: 1

[View Online »](#)

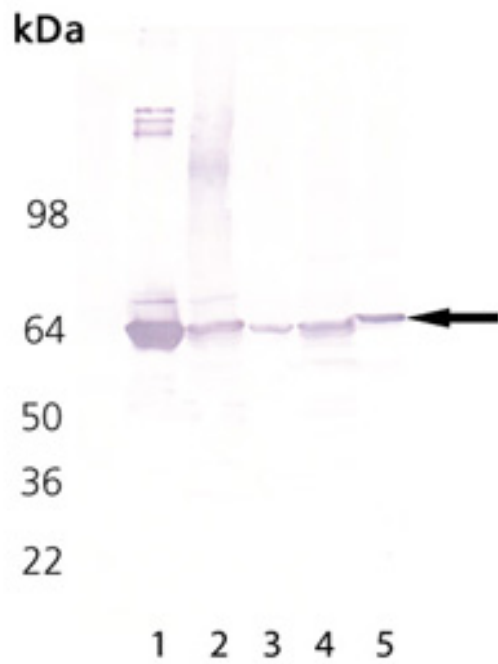
Ordering Information

[Order Online »](#)

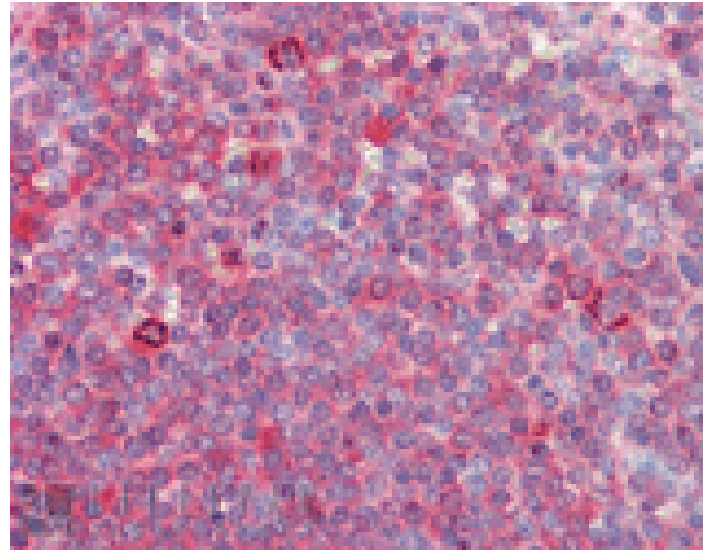
ADI-SPA-829-D	50µg
ADI-SPA-829-F	200µg

Manuals, SDS & CofA

[View Online »](#)



Western Blot Analysis of ADI-SPA-829: Lane 1: HSP60 protein(Human)(recombinant) (ADI-NSP-540), Lane 2: HeLa (Heat Shocked) Cell Lysate (ADI-LYC-HL101), Lane 3: PC-12 (Heat Shocked) Cell Lysate (ADI-LYC-PC101), Lane 4: 3T3 (Heat Shocked) Cell Lysate (ADI-LYC-3T101), Lane 5: GPC-16 (Heat Shocked) Cell Lysate



Immunohistochemistry analysis of human spleen tissue stained with HSP60, mAb (Mab11-13) at 5 μ g/ml.

Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Product Details

Alternative Name	Chaperonin 60, CPN60, HspD1, heat shock protein 60
Application	Flow Cytometry, IF, IHC (PS), WB
Application Notes	Detects a band of ~60kDa by Western blot.
Clone	Mab-11-13
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.
GenBank ID	M34664
Host	Mouse
Immunogen	Recombinant human Hsp60.
Isotype	IgG2a
Purity Detail	Protein G affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (10µg/ml)Western Blot (1:1,000, colorimetric)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, Dog, Dolphin, Drosophila, Fish, Guinea pig, Hamster, Human, Monkey, Mosquito, Mouse, Porcine, Rabbit, Rat, Snake

UniProt ID

P10809

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)