

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.

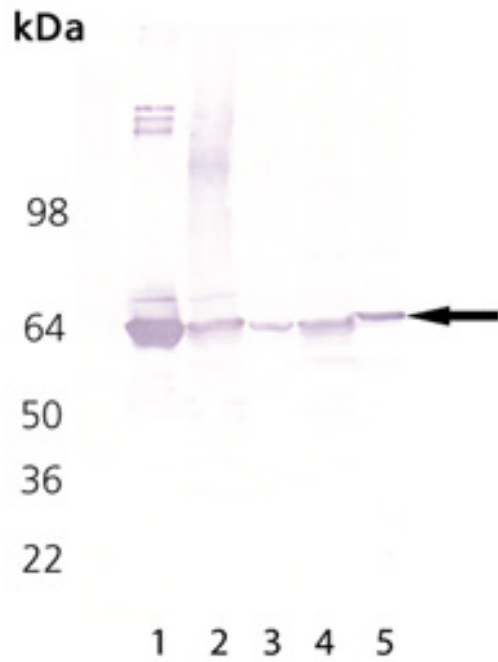
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 2

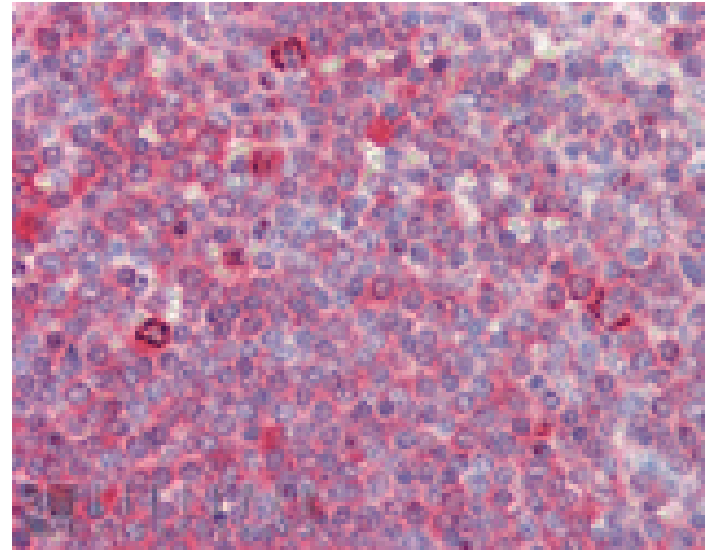
[View Online »](#)

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 µg/ml.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

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

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

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

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

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

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