

# HSP60 polyclonal antibody

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: 6

[View Online »](#)

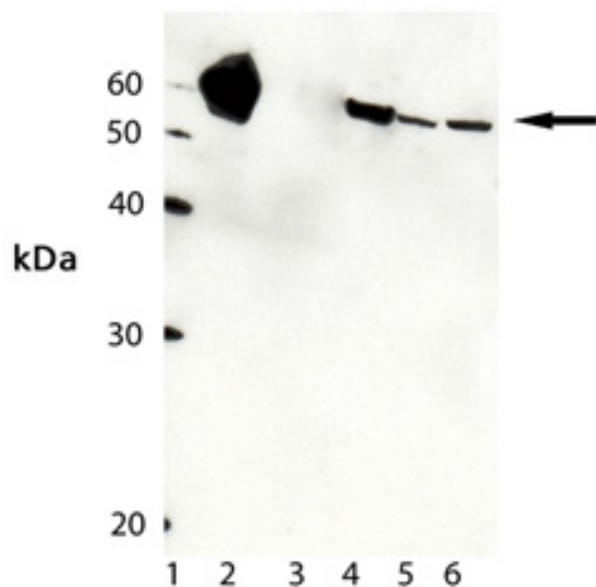
## Ordering Information

[Order Online »](#)

ADI-SPA-828-J	1mg
ADI-SPA-828-D	50µg
ADI-SPA-828-F	200µg

## Manuals, SDS & CofA

[View Online »](#)



Western Blot analysis of HSP60 pAb (Prod. No. ADI-SPA-828) Lane1: MW Marker; Lane 2:HSP60 protein (human) (recombinant) (Prod. No. ADI-NSP-540); Lane 3: HSP65 protein (M. bovis) (recombinant); Lane 4: HeLa Cell Lysate (Prod. No. ADI-LYC-HL100); Lane 5: 3T3 Cell Lysate; and Lane 6: RK-13 Cell Lysate.

# 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 IHC, IP, WB

Application Notes Detects a band of ~60kDa by Western blot.

Formulation Liquid. In PBS, pH 7.2, containing 50% glycerol and 0.09% sodium azide.

GenBank ID M34664

Host Goat

Immunogen Recombinant human Hsp60.

Purity Detail Protein G affinity purified.

Recommendation Dilutions/Conditions Immunoprecipitation (1:100)Western Blot (1:1,000, ECL)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.

Source Purified from goat serum.

Species Reactivity Beluga, Bovine, Chicken, Dog, Drosophila, Fish, Guinea pig, Hamster, Human, Monkey, Mouse, Mycobacterium bovis, Porcine, Rabbit, Rat, Sheep, Xenopus

UniProt ID P10809

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-  
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)