

HSP60 (insect)

polyclonal antibody

The Hsp60 of *Heliothis virescens* belongs to a highly conserved family of molecular chaperones from several species, including plant Hsp60 (known as Rubisco binding protein), GroEL, the *E. coli* Hsp60, and 65 kDa major antigen of mycobacteria. In eukaryotes, Hsp60 is localized in the mitochondrial matrix, and in plants Hsp60 is localized in the chloroplast. Mitochondria, chloroplasts and bacteria share a common ancestry (>1 billion years), and this coupled with the high degree of homology between the divergent Hsp60s suggests that these proteins perform a primitive but vital function similar to all the different species. The common characteristics shared by the Hsp60s from the divergent species include high abundance; induction with environmental stress such as heat shock; homo-oligomeric structures of either 7 or 14 subunits which reversibly dissociate in the presence of Mg^{2+} and ATP; ATPase activity; and a role in folding and assembly of oligomeric protein structures. Studies support these similarities, showing expression of the single-ring human mitochondrial homolog Hsp60 with its co-chaperonin Hsp10, in a *E. coli* strain engineered so that the *groE* operon remained under strict regulatory control. The findings demonstrate that expression of Hsp60-Hsp10 enabled successful performance of all essential *in vivo* functions of GroEL and its co-chaperonin, GroES. Consistent with their functions as chaperones, Hsp60 and Hsp10 may act as docking molecules with a passive role in the maturation of caspase processing. Data indicates that recombinant Hsp60 and Hsp10 accelerate the activation of procaspase-3 by cytochrome c and dATP in an ATP-dependent manner. Hsps are intracellular proteins thought to serve protective functions against infection and cellular stress; however, several studies reveal a possible link between members of the Hsp60 and a number of autoimmune diseases, atherosclerosis, and chlamydial disease.

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

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

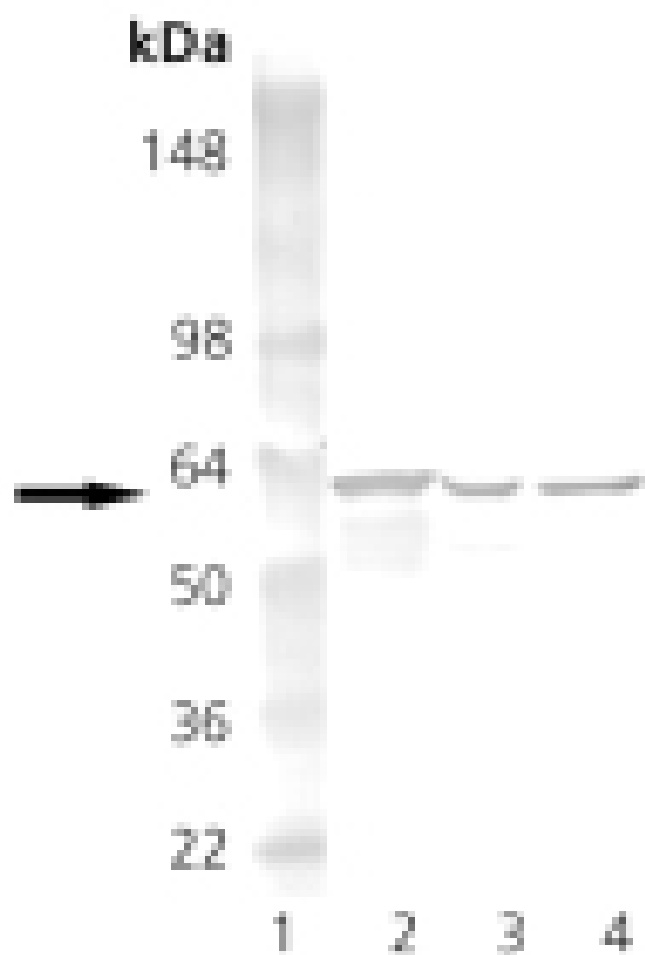
ADI-SPA-805-D

50µg

ADI-SPA-805-F	200µg
---------------	-------

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of HSP60: Lane 1: MW marker, Lane 2: HeLa (Heat Shocked), Lane 3: 3T3 (Heat Shocked), Lane 4: PC-12 (Heat Shocked).

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 IF, 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 X56034

Host Rabbit

Immunogen Native Hsp60 from Heliothis virescens (insect) sperm.

Purity Detail Protein A affinity purified.

Recommendation Dilutions/Conditions 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 rabbit serum.

Species Reactivity Beluga, Bovine, Chicken, Cockroach, Coral, Crab, Dog, Drosophila, E. coli, Ehrlichia, Fish, Grasshopper, Guinea pig, Hamster, Human, Lobster, Monkey, Moth, Mouse, Mussel, Porcine, Rabbit, Rat, Scallop, Sheep

UniProt ID P25420

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)