

HSP60 (low endotoxin) (mouse), (recombinant)

Mouse Hsp60 belongs to a highly conserved family that includes molecular chaperones from a variety of species such as plant Hsp60 (known as Rubisco binding protein), GroEL, the *E.coli* Hsp60, and the 65kDa major antigen of mycobacteria. Hsp60 is localized in the mitochondrial matrix of eukaryotes, and in the chloroplast of plants. Hsp60s from divergent species share a number of common characteristics: 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. These similarities correspond with studies in which the single-ring human mitochondrial homolog Hsp60 and its co-chaperonin Hsp10 were expressed in an *E. coli* strain engineered to keep the *groE* operon 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 cochaperonin, GroES. Several studies reveal a possible link between members of the Hsp60 family and a number of autoimmune diseases, atherosclerosis, and chlamydial disease. Overexpression of self Hsp60 is seen in the synovial tissue of rheumatoid arthritic (RA) patients, and can accompany both cellular and humoral reactivity against Hsp60 in RA. Chlamydial heat shock protein Hsp60, a homolog of *E. coli* GroEL, appears capable of eliciting macrophage activation, and several studies reveal a correlation between Hsp60 responses and the immunopathologic manifestations of human chlamydial disease.

Citations: 2

[View Online »](#)

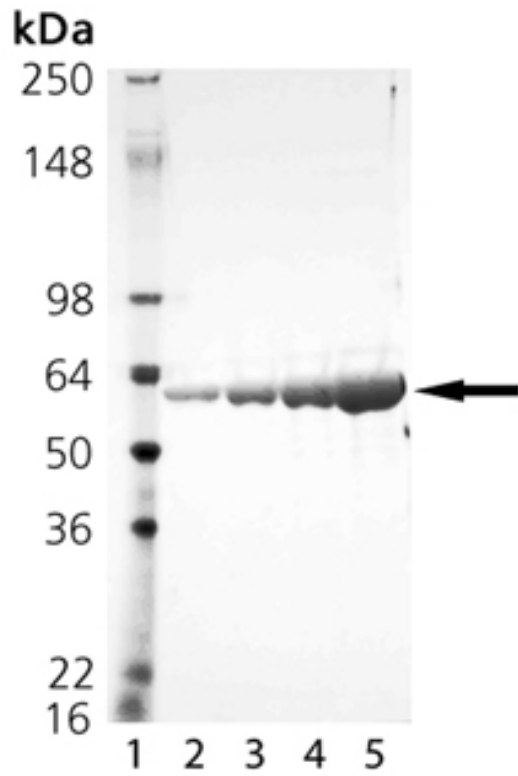
Ordering Information

[Order Online »](#)

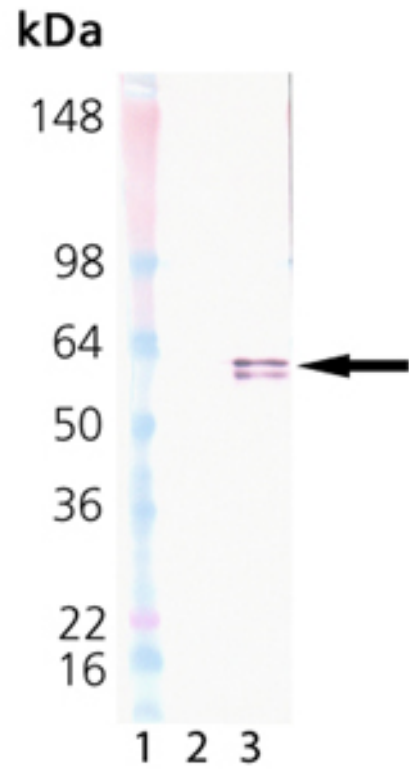
ADI-ESP-741-D	50µg
ADI-ESP-741-F	200µg

Manuals, SDS & CofA

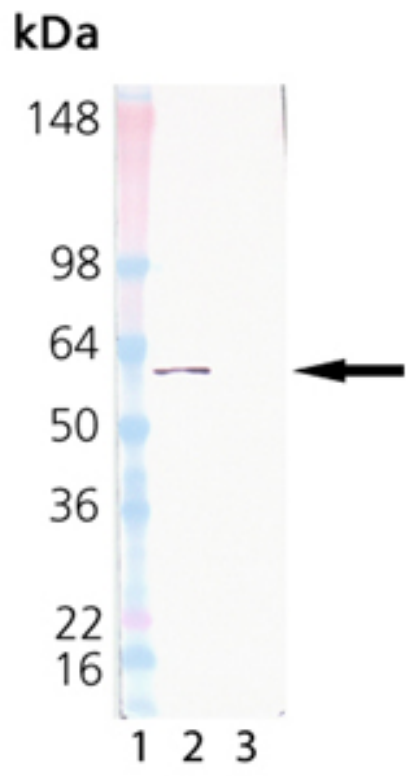
[View Online »](#)



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 0.5 μ g, Lane 3 : 1 μ g, Lane 4 : 2 μ g, Lane 5 : 5 μ g HSP60 (Prod. No. ADI-ESP-741).



Western Blot analysis: Lane 1: MWM, Lane 2: Hsp60 (Prod. No. ADI-ESP-741), Lane 3: GroEL (Prod. No. ADI-SPP-610) probed with anti-GroEL mAb (Prod. No. ADI-SPS-870).



Western Blot analysis: Lane 1: MWM, Lane 2: HSP60 (Prod. No. ADI-ESP-741), Lane 3: GroEL (Prod. No. ADI-SPP-610) probed with anti-HSP60 mAb (Prod. No. ADI-SPA-806).

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Chaperonin 60, CPN60, HspD1, Heat shock protein 60

Application Notes ATPase activity assay (positive). Western blot control.

Endotoxin Content <50EU/mg purified protein (LAL test)

Formulation Liquid. In 10mM TRIS, pH 7.5, containing 150mM sodium chloride, 1mM EDTA, and 1mM DTT.

MW ~60kDa

Purity ≥90% (SDS-PAGE; Western blot)

Purity Detail Purified by multi-step chromatography.

Source Produced in *E. coli*.

UniProt ID P63038



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