

Anti-HSP60 IgG/A/M (human), ELISA kit

Most comprehensive portfolio of heat shock protein (chaperone) products commercially available.

This assay allows for reproducible, accurate and precise determination of IgG, IgA and IgM antibodies (total) to human Hsp60 in serum. The ELISA kit uses recombinant human Hsp60 bound to the wells of the plate to bind anti-human Hsp60 antibodies present in human serum. The captured anti-human Hsp60 antibodies are detected with a HRP conjugated goat polyclonal antibody specific for human IgG, IgA and IgM molecules. The assay is developed with tetramethylbenzidine substrate producing a blue color in proportion to the amount of captured anti-human Hsp60 antibodies. The color development is stopped with acid stop solution which converts the endpoint color to yellow. The intensity of the color is measured in a microplate reader at 450 nm.

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.

Citations: 9

[View Online »](#)

Ordering Information

[Order Online »](#)

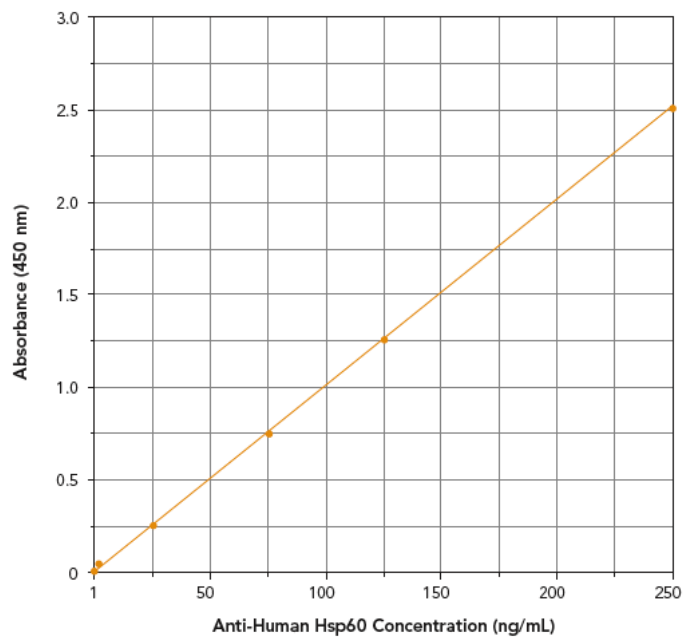
ADI-EKS-650

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Quantitative - fully quantitate levels of auto-antibodies to HSP60
- Higher throughput - obtain results from up to 41 samples in duplicate in less than 4 hours
- Easy-to-use - liquid color-coded reagents and pre-coated plates reduce errors and save time in the lab



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of anti-HSP60 IgG/A/M autoantibodies in human serum. Cited sample types include human plasma.

Assay Time <4 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents Microtiter plate, Standard, Sample diluent 1, Wash buffer concentrate, HRP Conjugate, TMB Substrate, Stop solution 2

Sensitivity 2.88 ng/ml (range 7.81 – 250 ng/ml)

Species Reactivity Human

UniProt ID P10809

Wavelength 450 nm



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-
fr@enzolifesciences.com](mailto:info-fr@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)