

HSP60 (human), ELISA kit

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

The HSP60 (human), EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 450 nm. Save time, money, and precious sample with fully quantitative results and increased sensitivity compared to Western blot analysis.

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: 15

[View Online »](#)

Ordering Information

[Order Online »](#)

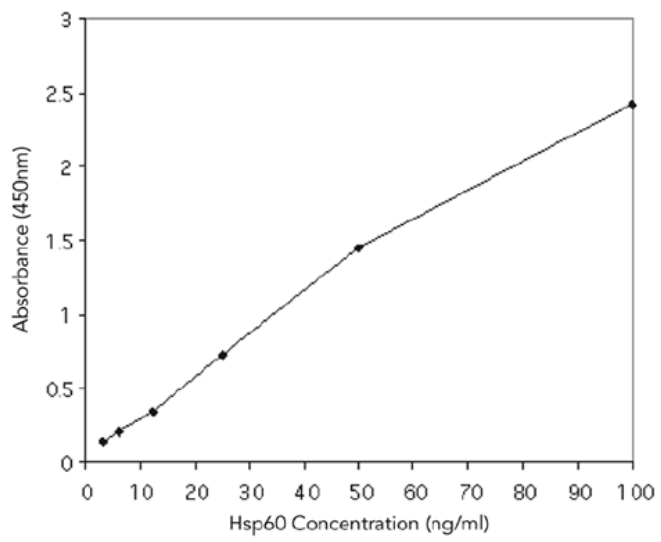
ADI-EKS-600

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive - use less precious sample per test compared to Western blot
- Higher throughput - test up to 40 samples in duplicate in just less than 3 hours
- Quantitative - obtain statistically significant results and compare data between multiple experiments



Handling & Storage

Use/Stability Store all components at 4°C, except standard at -20°C.

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Heat shock protein 60

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of HSP60 in cell lysates, serum, and tissue of human origin.

Assay Time <3 hours

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

Contents Microtiter plate, 5X Extraction reagent, Standard, Sample diluent, Wash buffer concentrate, Antibody, Antibody diluent, HRP Conjugate, Conjugate diluent, TMB Substrate, Stop solution 2

Crossreactivity HSP60 (100%). No reactivity with GroEL, HSP65, HSP27, HSP25, Cpn10, or GroES.

Sensitivity 3.125 ng/ml (range 3.125 – 100 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)