

# HSP70 High Sensitivity ELISA kit

**Only commercially available kit specifically validated and sensitive enough to quantify HSP70 (HSP72) in serum and plasma samples.**

The HSP70 high sensitivity EIA kit is a colorimetric immunometric enzyme immunoassay kit with results in 4.5 hours. Absorbance is read at 450 nm. This is the only commercially available kit specifically validated and sensitive enough to quantify HSP70 (HSP72) in serum and plasma samples, as referenced by the [Cell Stress Society International](#).

The Hsp70 family of heat shock proteins contains multiple homologs ranging in size from 66-78 kDa, and are the eukaryotic equivalents of the bacterial DnaK. The most studied Hsp70 members include the cytosolic stress-induced Hsp70 (Hsp72), the constitutive cytosolic Hsc70 (Hsp73), and the ER-localized BiP (Grp78). Hsp70 family members contain highly conserved N-terminal ATPase and C-terminal protein binding domains. Binding of peptide to Hsp70 is assisted by Hsp40, and stimulates the inherent ATPase activity of Hsp70, facilitating ATP hydrolysis and enhanced peptide binding. Hsp70 nucleotide exchange and substrate binding coordinates the folding of newly synthesized proteins, the re-folding of misfolded or denatured proteins, coordinates trafficking of proteins across cellular membranes, inhibits protein aggregation, and targets the degradation of proteins via the proteasomal pathway.

Citations: 70

[View Online »](#)

## Ordering Information

[Order Online »](#)

ADI-EKS-715

96 wells

## Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement, quantify as little as 90pg/ml Hsp70 (Hsp72)
- Fully validated for use with serum and plasma samples
- High specificity, measures stress-induced Hsp70 (Hsp72) with negligible reactivity from other Hsp70 family members (e.g. Hsc70/Hsp73, Grp78)
- Higher throughput format with results in just 4.5 hours for up to ~40 samples in duplicate

## 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 70, HSP72

**Application** Colorimetric detection, ELISA

**Application Notes** For the quantitative determination of HSP70 in serum and plasma of human, mouse, and rat origin. Cited sample(s) include cell lysates, urine, canine serum, and cerebrospinal fluid.

**Assay Time** 4.5 hours

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

**Contents** Microtiter plate, Assay buffer 28, Standard, Wash buffer concentrate, Antibody, Conjugate, TMB Substrate, Stop solution 2

**Crossreactivity** HSP72 (human) (100%), HSP70 (rat) (117.6%), HSP70 (salmon) (0.8%), HSP70B' (human) (5.4%), DnaK (E.coli) (0.5%), and <0.016%: Hsc70 (bovine), Grp78 (hamster), HSP71 (M. tuberculosis)

**Sensitivity** 90 pg/ml (range 0.20-12.5 ng/ml)

**Species Reactivity** Human, Mouse, Rat

**UniProt ID** P0DMV8 (HSPA1A), P0DMV9 (HSPA1B)

**Wavelength** 450 nm



Last modified: May 29, 2024

ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430

[info-us@enzolifesciences.com](mailto:info-us@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)