

## Product datasheet

# Human Heat Shock Protein 27 ELISA Kit (HSP27) ab113334

[1 Image](#)

### Overview

<b>Product name</b>	Human Heat Shock Protein 27 ELISA Kit (HSP27)
<b>Detection method</b>	Colorimetric
<b>Sample type</b>	Cell culture supernatant, Serum, Plasma
<b>Assay type</b>	Sandwich (quantitative)
<b>Sensitivity</b>	< 0.15 ng/ml
<b>Range</b>	0.205 ng/ml - 50 ng/ml
<b>Recovery</b>	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	133.1	123% - 151%
Serum	126.9	117% - 134%
Plasma	119.8	105% - 142%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Abcam's Heat Shock Protein 27 (HSP27) Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human HSP27 in serum, plasma, cell culture supernatants.

This assay employs an antibody specific for Human Heat Shock Protein 27 coated on a 96-well plate. Standards and samples are pipetted into the wells and Heat Shock Protein 27 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human Heat Shock Protein 27 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of Heat Shock Protein 27 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human Heat Shock Protein 27 ELISA Kit (HSP27) ([ab193757](#)) from our SimpleStep ELISA® range.

**Notes** Optimization may be required with urine samples.

**Tested applications** **Suitable for:** Sandwich ELISA

**Platform** Microplate

### Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer Concentrate	1 x 25ml
300X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
Biotinylated anti-Human Heat Shock Protein 27	2 vials
Heat Shock Protein 27 Microplate (12 x 8 wells)	1 unit
Recombinant Human Heat Shock Protein 27 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

**Relevance** Involved in stress resistance and actin organization. Interacts with TGFB111. Associates with alpha- and beta-tubulin, microtubules and CRYAB. Interacts with HSPB8 and HSPBAP1.

**Cellular localization** Cytoplasm. Nucleus. Cytoplasm; cytoskeleton; spindle. Note: Cytoplasmic in interphase cells. Colocalizes with mitotic spindles in mitotic cells. Translocates to the nucleus during heat shock and resides in sub-nuclear structures known as SC35 speckles or nuclear splicing speckles.

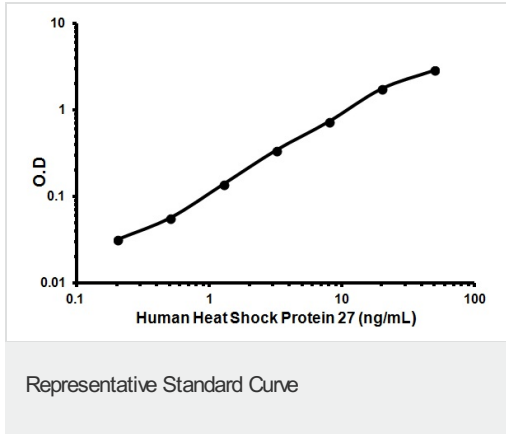
### Applications

Our [Abpromise guarantee](#) covers the use of **ab113334** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

### Images



Representative Standard Curve using ab113334

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

### Our Abpromise to you: Quality guaranteed and expert technical support

---

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
  
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

### Terms and conditions

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors