

Human Heat Shock Protein 27 ELISA Kit (HSP27) ab108862

5 References 1 图像

概述

产品名称	人Heat Shock蛋白27 ELISA试剂盒(HSP27)				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Overall				4.8%
	批次间				
	样品	n	Mean	SD	CV%
	Overall				9.9%
样品类型	Milk, Serum, Plasma, Tissue, Cell Lysate				
检测类型	Sandwich (quantitative)				
灵敏度	> 0.16 ng/ml				
范围	1.25 ng/ml - 20 ng/ml				
回收率	97 %				
检测时间	5h 00m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's Heat Shock Protein 27 (HSP27) Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of HSP27 levels in cell culture samples, milk, serum, plasma and tissue extract.				

A HSP27 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a HSP27 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of HSP27 captured in plate.

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

**The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.**

平台 Microplate

## 性能

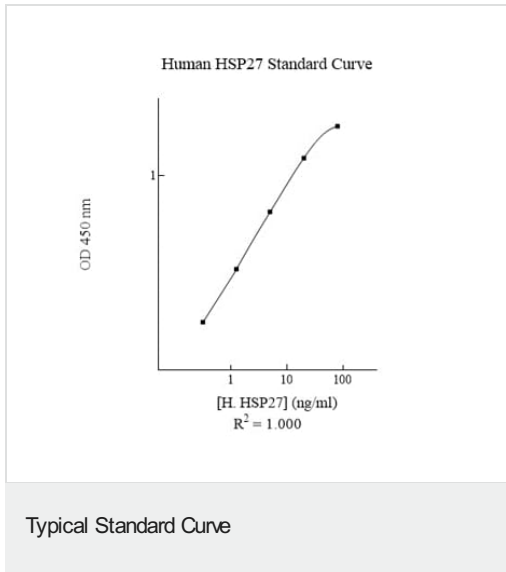
存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent M Concentrate	1 x 20ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
50X Biotinylated Human HSP27 Antibody	1 x 120µl
Chromogen Substrate	1 x 7ml
HSP27 Microplate (12 x 8 well strips)	1 unit
HSP27 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

相关性 Involved in stress resistance and actin organization. Interacts with TGFB1I1. Associates with alpha- and beta-tubulin, microtubules and CRYAB. Interacts with HSPB8 and HSPBAP1.

细胞定位 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.

## 图片



Representative Standard Curve using ab108862.

**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.cn/abpromise> or contact our technical team.

### Terms and conditions

---

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