

# Human Heat Shock Protein 47 ELISA Kit (SERPINH1) ab108861

[2 References](#) [1 图像](#)

### 概述

**产品名称** 人Heat Shock蛋白47 ELISA试剂盒(SERPINH1)

**检测方法** Colorimetric

**精确度**

批次内

样品	n	Mean	SD	CV%
Overall				6.1%

批次间

样品	n	Mean	SD	CV%
Overall				10.9%

**样品类型** Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid

**检测类型** Sandwich (quantitative)

**灵敏度** = 0.74 ng/ml

**范围** 1.563 ng/ml - 25 ng/ml

**回收率** 98 %

**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 47 (SERPINH1) Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of HSP47 levels in Human plasma, milk, serum, cell lysate, cerebrospinal fluid and tissue samples.

A HSP47 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently a HSP47 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 HSP47 captured in plate.

Get results in 90 minutes with Human Heat Shock Protein 47 ELISA Kit (SERPINH1) ([ab202407](#)) 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	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl	1 x 80µl
10X Diluent N Concentrate	1 x 30ml	1 x 30ml
1X Standard Diluent	1 x 2ml	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml	2 x 30ml
50X Biotinylated Human HSP47 Antibody	1 x 150µl	1 x 150µl
Chromogen Substrate	1 x 7ml	1 x 7ml
HSP47 Microplate (12 x 8 well strips)	1 unit	1 unit
HSP47 Standard	1 vial	1 vial
Sealing Tapes	3 units	3 units
Stop Solution	1 x 11ml	1 x 11ml

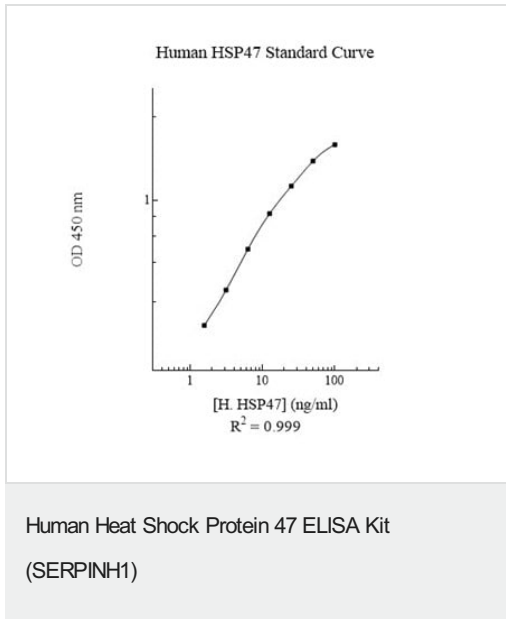
功能 Binds specifically to collagen. Could be involved as a chaperone in the biosynthetic pathway of collagen.

疾病相关 Note=Defects in SERPINH1 may cause severe autosomal recessive osteogenesis imperfecta (OI). Osteogenesis imperfecta defines a group of connective tissue disorders characterized by bone fragility and low bone mass.

序列相似性 Belongs to the serpin family.

细胞定位 Endoplasmic reticulum lumen.

## 图片



Representative Standard Curve using ab108861.

**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