

Product datasheet

Human Heat Shock Protein 47 ELISA Kit (SERPINH1)
ab108861

2 References 1 Image

Overview

Product name	Human Heat Shock Protein 47 ELISA Kit (SERPINH1)			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Overall			6.1%
	Inter-assay			
	Sample	n	Mean	SD
	Overall			10.9%
Sample type	Serum, Plasma, Tissue, Cell Lysate, Cerebral Spinal Fluid			
Assay type	Sandwich (quantitative)			
Sensitivity	= 0.74 ng/ml			
Range	1.563 ng/ml - 25 ng/ml			
Recovery	98 %			
Assay time	5h 00m			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's Heat Shock Protein 47 (SERPINH1) Human <i>in vitro</i> 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.

Platform Microplate

Properties

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

Components	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

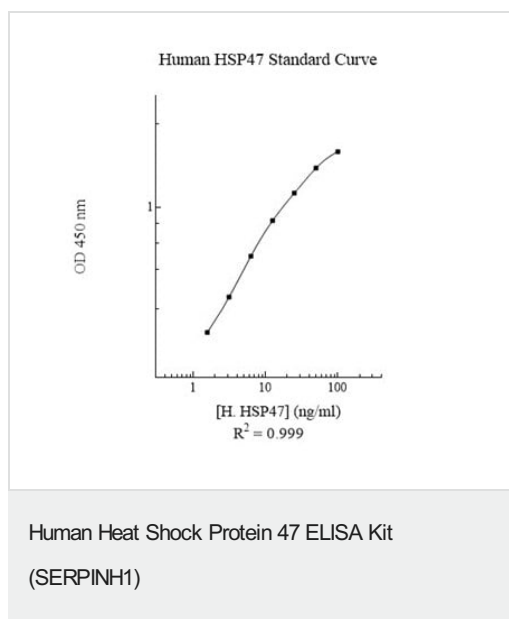
Function Binds specifically to collagen. Could be involved as a chaperone in the biosynthetic pathway of collagen.

Involvement in disease 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.

Sequence similarities Belongs to the serpin family.

Cellular localization Endoplasmic reticulum lumen.

Images



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.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