

Human Heat Shock Protein 27 ELISA Kit (HSP27) ab108862

5 References 画像数 1

製品の概要

製品名	Human Heat Shock Protein 27 ELISA Kit (HSP27)			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				4.8%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				9.9%
サンプルの種類	Milk, Serum, Plasma, Tissue, Cell Lysate			
アッセイタイプ	Sandwich (quantitative)			
検出感度	> 0.16 ng/ml			
検出範囲	1.25 ng/ml - 20 ng/ml			
添加回収試験	97 %			
全工程の試験時間	5h 00m			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	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.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	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

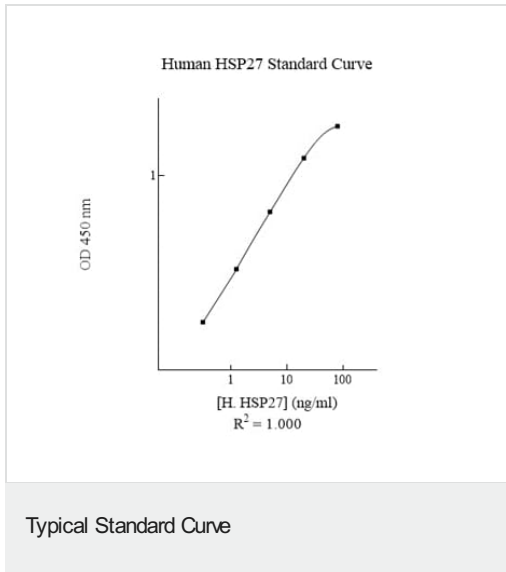
関連性

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

細胞内局在

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.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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