

Human Heat Shock Protein 27 ELISA Kit (HSP27) ab113334

1 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
サンプルの種類	Cell culture supernatant, Serum, Plasma
アッセイタイプ	Sandwich (quantitative)
検出感度	< 0.15 ng/ml
検出範囲	0.205 ng/ml - 50 ng/ml
添加回収試験	100 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Cell culture supernatant	133.1	123% - 151%
Serum	126.9	117% - 134%
Plasma	119.8	105% - 142%

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 ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> 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.

特記事項

Optimization may be required with urine samples.

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
20X Wash Buffer	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

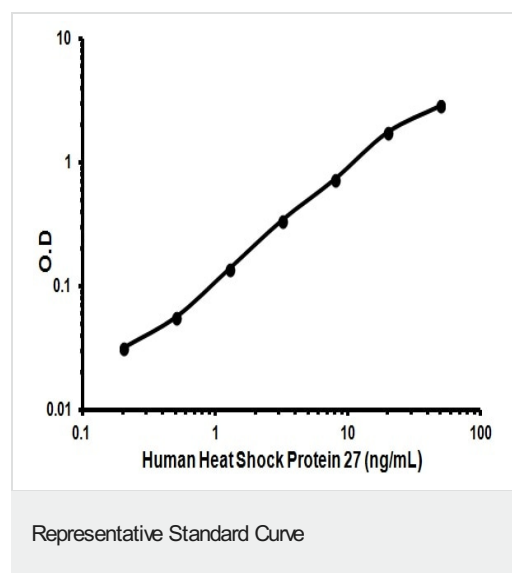
関連性

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 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.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