

Product datasheet

Human HTRA2 ELISA Kit ab155434

1 Image

Overview

Product name Human HTRA2 ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| overall | | | | < 10% |

Inter-assay

| Sample | n | Mean | SD | CV% |
|---------|---|------|----|-------|
| overall | | | | < 12% |

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 0.2 ng/ml

Range 0.205 ng/ml - 50 ng/ml

Recovery

Sample specific recovery

| Sample type | Average % | Range |
|--------------------|-----------|------------|
| Serum | = 100.1 | 93% - 107% |
| Plasma | = 103.4 | 91% - 117% |
| Cell culture media | = 97.26 | 91% - 103% |

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's HTRA2 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay designed for the quantitative measurement of Human HTRA2 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for Human HTRA2 coated on a 96-well plate. Standards

and samples are pipetted into the wells and HTRA2 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human HTRA2 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 HTRA2 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Notes Optimization may be required with urine samples.

Tested applications **Suitable for:** Sandwich ELISA

Platform Microplate

Properties

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

| Components | 1 x 96 tests |
|--|--------------|
| 20X Wash Buffer Concentrate | 1 x 25ml |
| 400X HRP-Streptavidin Concentrate | 1 x 200µl |
| 5X Assay Diluent B | 1 x 15ml |
| 5X Assay Diluent D | 1 x 15ml |
| Biotinylated anti-Human HTRA2 | 2 vials |
| HTRA2 Microplate (12 x 8 wells) | 1 unit |
| Recombinant Human HTRA2 Standard (lyophilized) | 2 vials |
| Stop Solution | 1 x 8ml |
| TMB One-Step Substrate Reagent | 1 x 12ml |

Relevance This gene encodes a serine protease. The protein has been localized in the endoplasmic reticulum and interacts with an alternatively spliced form of mitogen-activated protein kinase 14. The protein has also been localized to the mitochondria with release to the cytosol following apoptotic stimulus. The protein is thought to induce apoptosis by binding the apoptosis inhibitory protein baculoviral IAP repeat-containing 4. Nuclear localization of this protein has also been observed. Alternate splicing of this gene results in two transcript variants encoding different isoforms. Additional transcript variants have been described, but their full-length sequences have not been determined.

Cellular localization Mitochondrial, Endoplasmic reticulum, Cytoplasmic.

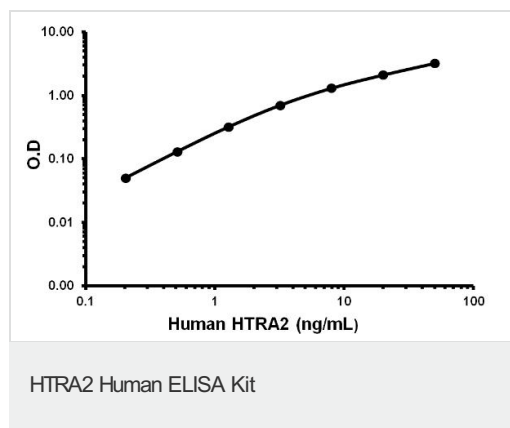
Applications

Our [Abpromise guarantee](#) covers the use of **ab155434** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

| Application | Abreviews | Notes |
|----------------|-----------|--|
| Sandwich ELISA | | Use at an assay dependent concentration. |

Images



Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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