

Product datasheet

Human Chorionic Gonadotropin ELISA Kit ab108637

Overview

| | |
|----------------------------|---|
| Product name | Human Chorionic Gonadotropin ELISA Kit |
| Detection method | Colorimetric |
| Sample type | Urine, Serum |
| Assay type | Sandwich (qualitative) |
| Sensitivity | = 20 mU/ml |
| Assay duration | Multiple steps standard assay |
| Species reactivity | Reacts with: Human |
| Product overview | <p>Abcam's Chorionic Gonadotropin <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of Chorionic Gonadotropin in Human urine or serum.</p> <p>A 96-well plate has been precoated with anti-Chorionic Gonadotropin antibodies. Samples and standards are added to the wells, where Chorionic Gonadotropin in the sample and standards binds to the precoated antibody. After incubation and washing, added Anti-Chorionic Gonadotropin HRP conjugate binds to the antibody-Chorionic Gonadotropin complex. After incubation, the wells are washed to remove unbound material and TMB substrate is then added which is catalyzed by HRP to produce blue coloration. The reaction is terminated by addition of Stop Solution which stops the color development and produces a color change from blue to yellow. The intensity of signal is directly proportional to the amount of Chorionic Gonadotropin in the sample and the intensity is measured at 450 nm.</p> |
| Tested applications | Suitable for: Sandwich ELISA |
| Platform | Microplate |

Properties

Storage instructions Store at +4°C. Please refer to protocols.

| Components | Identifier | 1 x 96 tests |
|--|------------|--------------|
| Anti-Chorionic Gonadotropin Coated Microplate (12 x 8 wells) | | 1 unit |
| Anti-Chorionic Gonadotropin HRP Conjugate | | 1 x 7ml |
| Chorionic Gonadotropin Standard 0 - 0 mU/mL | White cap | 1 x 1ml |

| Components | Identifier | 1 x 96 tests |
|---|------------|--------------|
| Chorionic Gonadotropin Standard 1 - 20 mU/mL | Yellow cap | 1 x 1ml |
| Chorionic Gonadotropin Standard 2 - 150 mU/mL | Black cap | 1 x 1ml |
| Stop Solution | | 1 x 7ml |
| TMB Reagent | | 1 x 7ml |

Applications

Our [Abpromise guarantee](#) covers the use of **ab108637** 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 dilution. |

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