

Human Chorionic Gonadotropin beta ELISA Kit ab178633

2 References

Overview

Product name Human Chorionic Gonadotropin beta ELISA Kit

Detection method Colorimetric

Precision Intra-assay

Sample	n	Mean	SD	CV%
Overall	48			< 7.6%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	3			< 8.8%

Sample type Serum, Hep Plasma, Cit plasma

Assay type Sandwich (quantitative)

Sensitivity > 0.09 mU/ml

Range 1 mU/ml - 400 mU/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's Chorionic Gonadotropin beta *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Chorionic Gonadotropin beta in Human serum and plasma.

A 96-well plate has been precoated with anti-Chorionic Gonadotropin beta IgG antibodies. Samples, standards and the Chorionic Gonadotropin beta-HRP conjugate are added to the wells, where Chorionic Gonadotropin beta in the sample and standards binds to the precoated antibody and added Chorionic Gonadotropin beta-HRP conjugate binds to this antibody-Chorionic Gonadotropin beta 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 beta in the sample and the intensity is measured at 450 nm.

Platform

Microplate

Properties**Storage instructions**

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

Components	1 x 96 tests
10X Anti-Chorionic Gonadotropin beta HRP Conjugate	1 x 1ml
50X Washing Solution	1 x 20ml
Anti-Chorionic Gonadotropin beta Coated Microplate (12 x 8 wells)	1 unit
Chorionic Gonadotrophin beta Control	1 x 1ml
Chorionic Gonadotropin beta Standard 0 - 0 mU/mL	1 x 1ml
Chorionic Gonadotropin beta Standard 1 - 1 mU/mL	1 x 1ml
Chorionic Gonadotropin beta Standard 2 - 5 mU/mL	1 x 1ml
Chorionic Gonadotropin beta Standard 3 - 20 mU/mL	1 x 1ml
Chorionic Gonadotropin beta Standard 4 - 100 mU/mL	1 x 1ml
Chorionic Gonadotropin beta Standard 5 - 400 mU/mL	1 x 1ml
Cover Foil	1 unit
Incubation Buffer	1 x 50ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

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