

Product datasheet

Human Chorionic Gonadotropin beta ELISA Kit ab108638

★★★★★ 2 Abreviews [画像数 2](#)

製品の概要

製品名	Human Chorionic Gonadotropin beta ELISA Kit
検出方法	Colorimetric
サンプルの種類	Serum
アッセイタイプ	Sandwich (quantitative)
検出感度	0.25 ng/ml
検出範囲	0.25 ng/ml - 50 ng/ml
ステップ	Multiple steps standard assay
種交差性	交差種: Human
製品の概要	<p>Abcam's Chorionic Gonadotropin beta <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Chorionic Gonadotropin beta in Human serum.</p> <p>A 96-well plate has been precoated with anti-Chorionic Gonadotropin beta antibodies. Samples and standards are added to the wells, where Chorionic Gonadotropin beta in the sample and standards binds to the precoated antibody. After incubation and washing, added anti-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.</p>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Microplate

製品の特性

保存方法 Store at +4°C. Please refer to protocols.

内容	1 x 96 tests
Anti-Chorionic Gonadotropin beta Coated Microplate (12 x 8 wells)	1 unit
Anti-Chorionic Gonadotropin beta HRP Conjugate	1 x 18ml

内容	1 x 96 tests
Chorionic Gonadotropin beta Standard 1 - 0 ng/mL (Lyophilized)	1 vial
Chorionic Gonadotropin beta Standard 2 - 2.5 ng/mL (Lyophilized)	1 vial
Chorionic Gonadotropin beta Standard 3 - 5 ng/mL (Lyophilized)	1 vial
Chorionic Gonadotropin beta Standard 4 - 10 ng/mL (Lyophilized)	1 vial
Chorionic Gonadotropin beta Standard 5 - 25 ng/mL (Lyophilized)	1 vial
Chorionic Gonadotropin beta Standard 6 - 50 ng/mL (Lyophilized)	1 vial
Stop Solution	1 x 11ml
TMB Substrate Solution	1 x 11ml
Zero Buffer	1 x 13ml

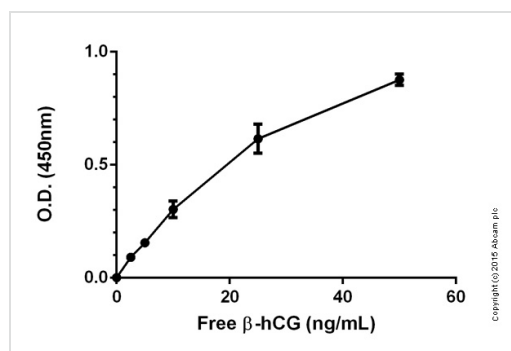
アプリケーション

Our [Abpromise guarantee](#) covers the use of **ab108638** in the following tested applications.

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

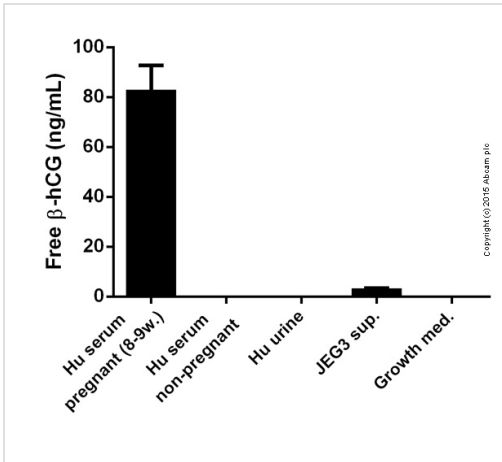
アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



Colorimetric standard curve: mean of duplicates (+/-) with background readings subtracted

Sandwich ELISA - Chorionic Gonadotropin beta
Human ELISA Kit ab108638



Free beta hCG measured in various samples showing concentration in ng/mL

Sandwich ELISA - Chorionic Gonadotropin beta
Human ELISA Kit ab108638

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.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