

Human hCG ELISA Kit ab100533

14 References 2 图像

概述

产品名称	人hCG ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 50 pg/ml
范围	54.87 pg/ml - 40000 pg/ml
回收率	> 90 %

特定样本回收率

样品类型	平均 %	范围
Cell culture supernatant	130	120% - 135%
Serum	113.7	103% - 124%
Plasma	90.57	83% - 103%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's hCG Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an <i>in vitro</i> enzyme linked immunosorbent assay for the quantitative measurement of Human hCG in serum, plasma, and cell culture supernatants.</p> <p>This assay employs an antibody specific for Human hCG coated on a 96-well plate. Standards and samples are pipetted into the wells and hCG present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human hCG 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 hCG bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.</p>
说明	Optimization may be required with urine samples
平台	Microplate

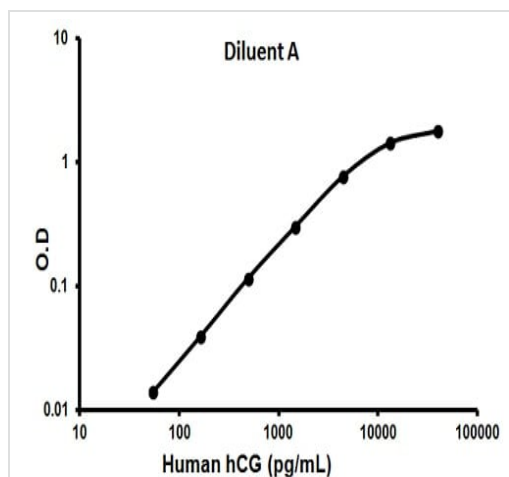
性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
700X HRP-Streptavidin Concentrate	1 x 200µl
Assay Diluent A	1 x 30ml
Biotinylated anti-Human hCG	2 vials
hCG Microplate (12 x 8 wells)	1 unit
Human hCG Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

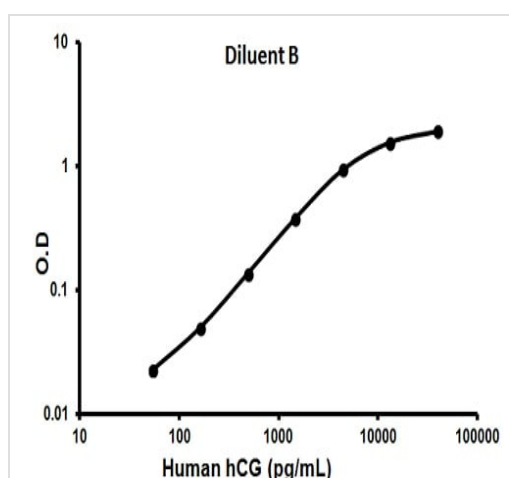
相关性 Human chorionic gonadotropin (hCG) is a glycoprotein hormone produced by trophoblastic cells of the placenta beginning 10 to 12 days after conception. Maintenance of the fetus in the first trimester of pregnancy requires the production of hCG, which binds to the corpus luteum of the ovary which is stimulated to produce progesterone which in turn maintains the secretory endometrium. hCG is present only in trace amounts in non pregnant urine and sera. It rises sharply during pregnancy. HCG is composed of two non identical, non covalently linked polypeptide chains designated as the a and b subunits. The a subunit of HCG is nearly identical to that of thyroid stimulating hormone (TSH), follicle stimulating hormone (FSH) and luteinizing hormone (LH). A germ cell tumor which is positive for cytokeratin, placental alkaline phosphatase (PLAP) and HCG but negative for EMA and AFP is probably a choriocarcinoma.

图片



Representative standard curve using ab100533

Typical standard curve



Representative standard curve using ab100533

Typical standard curve

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.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors