

17 beta Estradiol ELISA Kit ab108667

★★★★★ 2 Abreviews 45 References 2 图像

概述

产品名称17 beta Estradiol ELISA试剂盒

检测方法Colorimetric

批次内				
样品	n	Mean	SD	CV%
Overall	20			<=9%

批次间				
样品	n	Mean	SD	CV%
Overall	3			<=10%

样品类型Urine, Serum, Hep Plasma, EDTA Plasma, Cit plasma

检测类型Competitive

灵敏度8.68 pg/ml

范围20 pg/ml - 2000 pg/ml

回收率101.09 %

实验步骤Multiple steps standard assay

种属反应性与反应: Mammals

产品概述Abcam's 17 beta Estradiol *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of 17 beta Estradiol in serum and plasma (citrate).

A 96-well plate has been precoated with anti-Estradiol IgG. Samples and the Estradiol-HRP conjugate are added to the wells, where Estradiol in the sample competes with the added Estradiol-HRP for antibody binding. 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 inversely proportional to the amount of Estradiol in the sample and the intensity is measured at 450 nm.

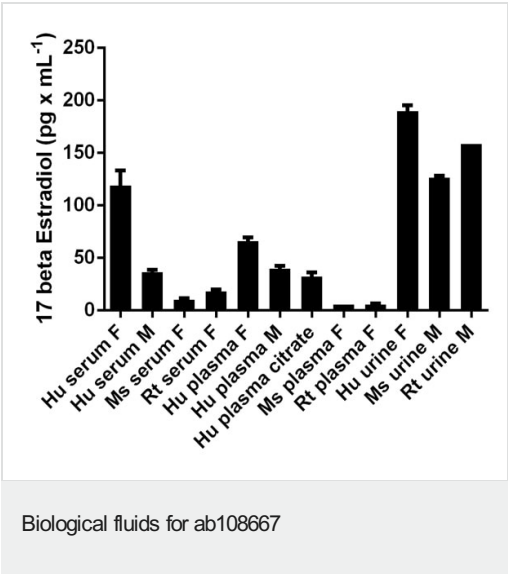
平台Microplate

性能

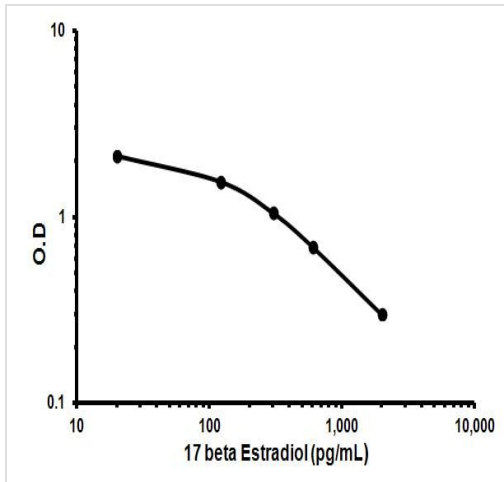
存放说明Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Washing Solution	1 x 50ml
17 beta Estradiol Control	1 x 0.5ml
17 beta Estradiol Standard 0	1 x 0.5ml
17 beta Estradiol Standard 1 (20 pg/mL)	1 x 0.5ml
17 beta Estradiol Standard 2 (120 pg/mL)	1 x 0.5ml
17 beta Estradiol Standard 3 (300 pg/mL)	1 x 0.5ml
17 beta Estradiol Standard 4 (600 pg/mL)	1 x 0.5ml
17 beta Estradiol Standard 5 (2,000 pg/mL)	1 x 0.5ml
17 beta Estradiol-HRP Conjugate	1 x 22ml
Anti-17 beta Estradiol IgG Coated Microplate (12 x 8 wells)	1 unit
Cover foils	1 unit
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

图片



17 beta Estradiol measured in biological fluids showing quantity (pg) per mL of tested sample. Samples were diluted 1-2 fold.



Representative Standard Curve using ab108667

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