

Human Luteinizing hormone ELISA Kit (LH) ab178658

2 References 1 图像

概述

产品名称

检测方法

精确度

人Luteinizing hormone ELISA试剂盒(LH)

Colorimetric

批次内

样品	n	Mean	SD	CV%
Serum	20			<=9.21%

批次间

样品	n	Mean	SD	CV%
Serum	15			<=7.91%

样品类型

检测类型

灵敏度

范围

实验步骤

种属反应性

产品概述

Serum, Hep Plasma, Cit plasma

Sandwich (quantitative)

= 0.22 mU/ml

5 mU/ml - 200 mU/ml

Multiple steps standard assay

与反应: Human

Abcam's Luteinizing hormone (LH) Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of LH in serum.

A 96-well plate has been precoated with Streptavidin. Samples, standards and the LH-HRP and Bitoin conjugate are added to the wells. Biotinylated monoclonal and enzyme labeled antibodies are added and the reactants mixed: these antibodies have high affinity and specificity and are directed against distinct and different epitopes of LH. Reaction between the various LH antibodies and native LH occurs in the microwells without competition or steric hindrance forming a soluble sandwich 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 LH 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
50X Washing Solution	1 x 20ml
Cover Foil	1 unit
Luteinizing Hormone Coated Microplate (12 x 8 wells)	1 unit
Luteinizing Hormone Control	1 x 1ml
Luteinizing Hormone HRP and Biotin Conjugate	1 x 12ml
Luteinizing Hormone Standard 0 - 0 mU/mL	1 x 1ml
Luteinizing Hormone Standard 1 - 5 mU/mL	1 x 1ml
Luteinizing Hormone Standard 2 - 25 mU/mL	1 x 1ml
Luteinizing Hormone Standard 3 - 50 mU/mL	1 x 1ml
Luteinizing Hormone Standard 4 - 100 mU/mL	1 x 1ml
Luteinizing Hormone Standard 5 - 200 mU/mL	1 x 1ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

功能 Promotes spermatogenesis and ovulation by stimulating the testes and ovaries to synthesize steroids.

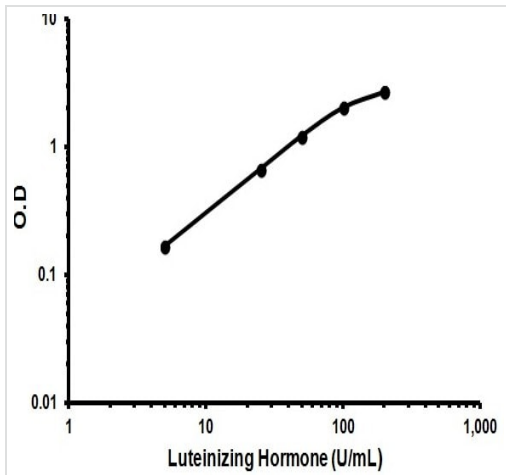
组织特异性 Pituitary gland.

疾病相关 Defects in LHB are a cause of hypogonadism LHB-related (HGON-LHB) [MIM:152780]. Hypogonadism is characterized by infertility and pseudohermaphroditism.

序列相似性 Belongs to the glycoprotein hormones subunit beta family.

细胞定位 Secreted.

图片



Representative standard curve using ab178658

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