

### Androstenedione ELISA Kit ab108672

[3 References](#) [1 图像](#)

#### 概述

**产品名称** Androstenedione ELISA试剂盒

**检测方法** Colorimetric

**精确度** 批次内

样品	n	Mean	SD	CV%
Control Sera	60			<=10%

批次间

样品	n	Mean	SD	CV%
Control Sera	30			<=9.5%

**样品类型** Serum, Hep Plasma, Cit plasma

**检测类型** Competitive

**灵敏度** = 0.01 ng/ml

**范围** 0.1 ng/ml - 10 ng/ml

**回收率** 100.91 %

**实验步骤** Multiple steps standard assay

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

A 96-well plate has been precoated with anti-Androstenedione IgG. Samples and the Androstenedione-HRP conjugate are added to the wells, where Androstenedione in the sample competes with the added Androstenedione-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 Androstenedione in the sample and the intensity is measured at 450 nm.

**平台** Microplate

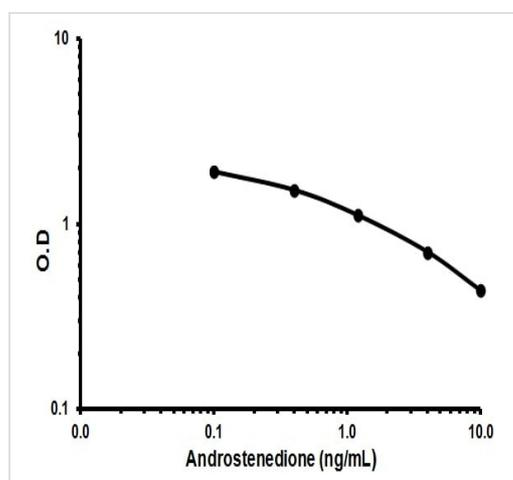
## 性能

### 存放说明

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

组件	1 x 96 tests
10X Washing Solution	1 x 50ml
Androstenedione Control	1 x 1ml
Androstenedione Standard 0 – 0 ng/mL	1 x 1ml
Androstenedione Standard 1 – 0.1 ng/mL	1 x 1ml
Androstenedione Standard 2 – 0.4 ng/mL	1 x 1ml
Androstenedione Standard 3 – 1.2 ng/mL	1 x 1ml
Androstenedione Standard 4 – 4.0 ng/mL	1 x 1ml
Androstenedione Standard 5 – 10.0 ng/mL	1 x 1ml
Androstenedione-HRP Conjugate	1 x 21ml
Anti-Androstenedione 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



Representative Standard Curve using ab108672

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