

## Product datasheet

# Androstenedione ELISA Kit ab108672

[1 References](#) [1 Image](#)

### Overview

**Product name** Androstenedione ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

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

Inter-assay

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

**Sample type** Serum, Heparin Plasma, Citrate Plasma

**Assay type** Competitive

**Sensitivity** = 0.01 ng/ml

**Range** 0.1 ng/ml - 10 ng/ml

**Recovery** 100.91 %

**Assay duration** Multiple steps standard assay

**Product overview** 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.

**Tested applications** **Suitable for:** Competitive ELISA

**Platform** Microplate

## Properties

### Storage instructions

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

Components	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

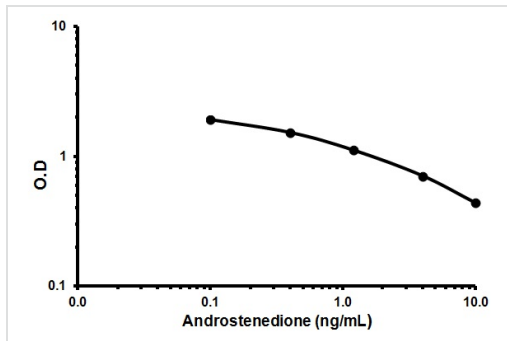
## Applications

Our [Abpromise guarantee](#) covers the use of **ab108672** in the following tested applications.

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

Application	Abreviews	Notes
Competitive ELISA		Use at an assay dependent dilution.

## Images



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.com/abpromise> or contact our technical team.

### Terms and conditions

---

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