

Androstenedione ELISA Kit ab178609

1 Image

Overview

Product name Androstenedione ELISA Kit

Detection method Colorimetric

Precision

Inter-assay

Sample	n	Mean	SD	CV%
Saliva	10			<=11%

Sample type Saliva

Assay type Competitive

Sensitivity 5 pg/ml

Range 20 pg/ml - 1000 pg/ml

Recovery 102.6 %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's Androstenedione *in vitro* competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Androstenedione in saliva.

A 96-well plate has been precoated with anti-Androstenedione antibodies. Samples and the Androstenedione-HRP conjugate are added to the wells, where Androstenedione in the sample competes with the added Androstenedione-HRP conjugate 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.

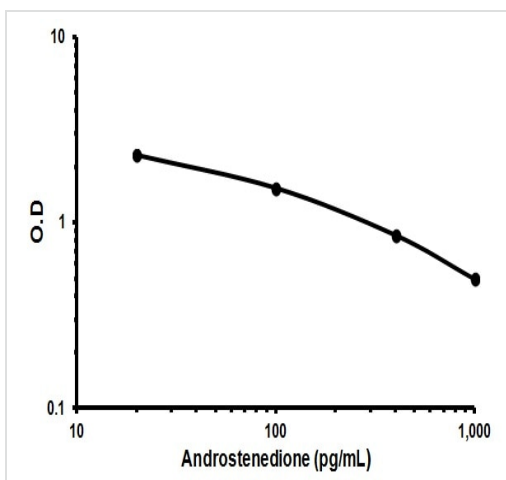
Platform Microplate

Properties

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

Components	1 x 96 tests
50X Washing Solution	1 x 20ml
Androstenedione HRP Conjugate	1 x 1ml
Androstenedione Standard 0 - 0 pg/mL	1 x 1ml
Androstenedione Standard 1 - 20 pg/mL	1 x 1ml
Androstenedione Standard 2 - 100 pg/mL	1 x 1ml
Androstenedione Standard 3 - 400 pg/mL	1 x 1ml
Androstenedione Standard 4 - 1,000 pg/mL	1 x 1ml
Anti-Androstenedione IgG Coated Microplate (12 x 8 wells)	1 unit
Control A	1 vial
Control B	1 vial
Cover Foil	1 unit
Incubation Buffer	1 x 30ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

Images



Typical Standard Curve

Representative standard curve using ab178609.

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