

Product datasheet

Estriol ELISA Kit ab178662

Overview

Product name Estriol ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Saliva	16			<=9.7%

Inter-assay

Sample	n	Mean	SD	CV%
Saliva	10			<=13.7%

Sample type Saliva

Assay type Competitive

Sensitivity 1 pg/ml

Recovery

Sample specific recovery

Sample type	Average %	Range
Spike	104	50pg/ml - 300pg/ml

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

Abcam's Estriol in vitro competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of Estriol in saliva. A 96-well plate has been precoated with anti-Estriol antibodies (solid-phase). Samples and the Estriol HRP Conjugate are added to the wells, where Estriol in the sample competes with the added Estriol 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 Estriol 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
anti-Estriol IgG Coated Microplate (12 x 8 wells)	1 unit
Cover Foil	1 unit
Estriol HRP Conjugate	1 x 1ml
Estriol Low Control	1 x 1ml
Estriol Control	1 x 1ml
Estriol Standard 0 - 0 pg/mL	1 x 1ml
Estriol Standard 1 - 2.5 pg/mL	1 x 1ml
Estriol Standard 2 - 15 pg/mL	1 x 1ml
Estriol Standard 3 - 100 pg/mL	1 x 1ml
Estriol Standard 4 - 600 pg/mL	1 x 1ml
Estriol Standard 5 - 4,000 pg/mL	1 x 1ml
Incubation Buffer	1 x 30ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

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