

DHEA-S ELISA Kit ab178642

1 Image

Overview

Product name DHEA-S ELISA Kit

Detection method Colorimetric

Precision

Intra-assay

Sample	n	Mean	SD	CV%
Overall	14			<=7.8%

Inter-assay

Sample	n	Mean	SD	CV%
Overall	9			<=14.9%

Sample type Saliva

Assay type Competitive

Sensitivity 0.05 ng/ml

Range 0.2 ng/ml - 12 ng/ml

Recovery 108.86 %

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview

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

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

Platform Microplate

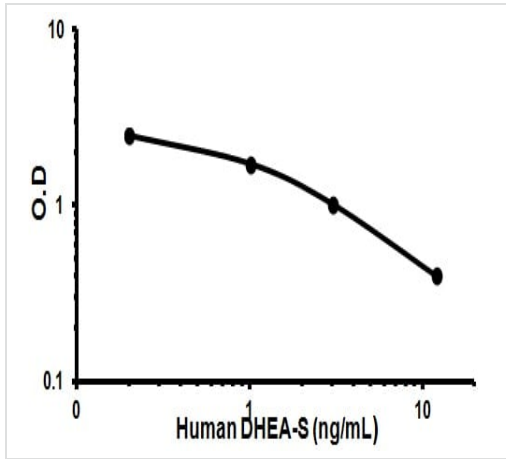
Properties

Storage instructions

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

Components	1 x 96 tests
50X Washing Solution	1 x 20ml
Anti-DHEA-S IgG Coated Microplate (12 x 8 wells)	1 unit
Cover Foil	1 unit
DHEA-S High Control	1 x 1ml
DHEA-S HRP Conjugate	1 x 1ml
DHEA-S Low Control	1 x 1ml
DHEA-S Standard 0 - 0 ng/mL	1 x 1ml
DHEA-S Standard 1 - 0.2 ng/mL	1 x 1ml
DHEA-S Standard 2 - 1.0 ng/mL	1 x 1ml
DHEA-S Standard 3 - 3.0 ng/mL	1 x 1ml
DHEA-S Standard 4 - 12.0 ng/mL	1 x 1ml
Incubation Buffer	1 x 30ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

Images



Representative standard curve using ab178642.

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