

DHEA sulfate (DHEA-S) ELISA Kit ab108669

4 References 1 Image

Overview

Product name	DHEA sulfate (DHEA-S) ELISA Kit			
Detection method	Colorimetric			
Precision	Intra-assay			
	Sample	n	Mean	SD
	Control	64		
				CV%
				< 7.9%
	Inter-assay			
	Sample	n	Mean	SD
	Control	48		
				CV%
				< 10.4%
Sample type	Serum, Hep Plasma, EDTA Plasma, Cit plasma			
Assay type	Competitive			
Sensitivity	= 0.04 ng/ml			
Range	0.1 µg/ml - 10 µg/ml			
Recovery	= 96.71 %			
Assay duration	Multiple steps standard assay			
Species reactivity	Reacts with: Human			
Product overview	Abcam's DHEA sulfate (DHEA-S) <i>in vitro</i> competitive ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate quantitative measurement of DHEA sulfate in plasma and serum.			

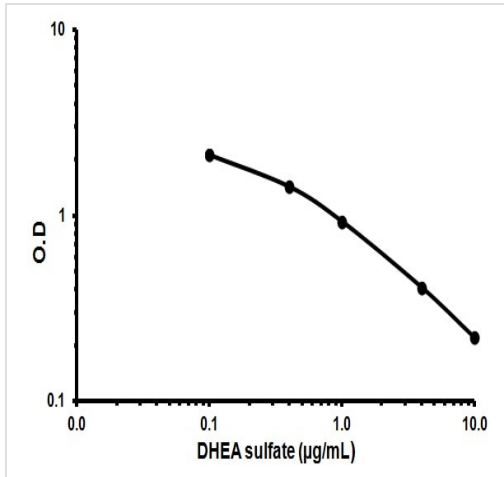
A 96-well plate has been precoated with anti-DHEA sulphate antibodies. Samples and the DHEA sulfate-HRP conjugate are added to the wells, where DHEA sulfate in the sample competes with the added DHEA sulfate-HRP for antibody binding. After incubation, the wells are washed to remove unbound material and 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 DHEA sulfate in the sample and the intensity is measured at 450 nm.

Properties**Storage instructions**

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

Components	1 x 96 tests
10X Wash Solution	1 x 50ml
Serum Diluent	1 x 100ml
Anti-DHEA sulfate antibody Coated Microplate (12 x 8 wells)	1 unit
Control A	1 x 1ml
Control B	1 x 1ml
Cover foils	1 unit
DHEA sulfate Standard 0 – 0 µg/mL	1 x 1ml
DHEA sulfate Standard 1 – 0.1 µg/mL	1 x 1ml
DHEA sulfate Standard 2 – 0.4 µg/mL	1 x 1ml
DHEA sulfate Standard 3 – 1.0 µg/mL	1 x 1ml
DHEA sulfate Standard 4 – 4.0 µg/mL	1 x 1ml
DHEA sulfate Standard 5 – 10.0 µg/mL	1 x 1ml
DHEA sulfate-HRP Conjugate	1 x 12ml
Stop Solution	1 x 15ml
Strip holder	1 unit
TMB Substrate Solution	1 x 15ml

Images



Representative Standard Curve using ab108669

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