

DHEA ELISA kit

Highly sensitive ELISA, measuring DHEA in culture supernatants, plasma, serum, saliva, and urine from any species.

The DHEA ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 5 hours. Absorbance is read at 405 nm. This kit has low reactivity to structurally related molecules and offers sensitive measurement of DHEA in a variety of sample types.

Dehydroepiandrosterone (DHEA) is a 17-ketosteroid produced by side chain cleavage of 17-hydroxypregnenolone. DHEA is reversibly converted to its sulfate ester (DHEA-S) and Androstendiol. The human adrenal glands secrete both DHEA and DHEA-S which in turn serve as the precursors for both estrogenic and androgenic steroids. DHEA has been linked to a number of physiological conditions including depression, systemic lupus erythematosus, and rheumatoid arthritis.

Citations: 36

[View Online »](#)

Ordering Information

[Order Online »](#)

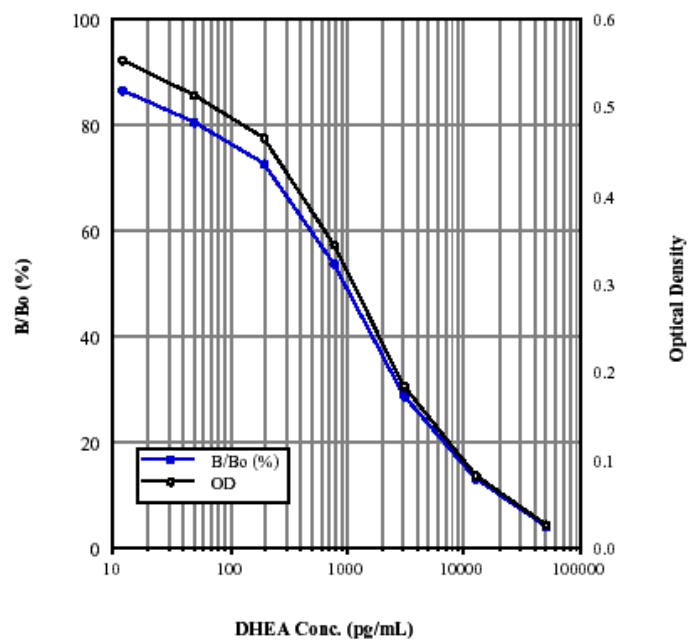
ADI-900-093

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Specific with low reactivity to structurally related molecules
- Highly reproducible with less than 9% inter- and intra-assay variation
- Highly sensitive, detecting low levels of DHEA
- Fast results for up to 37 samples in duplicate in 5 hours



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dehydroepiandrosterone

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of DHEA in culture supernatants, plasma, serum, saliva, and urine from any species.

Assay Time 5 hours

Compatibility This product is compatible with the [Absorbance 96 Plate Reader](#).

Contents GxR IgG Microtiter plate, Assay buffer, Conjugate, Antibody, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution

Crossreactivity DHEA (100%), DHEA sulfate (30.0%), Androstenedione (0.73%), Androsterone (0.29%), Pregnenolone (0.28%), Testosterone (0.10%), Progesterone (0.06%), Estradiol (0.04%) and <0.02%: 20-hydroxycholesterol, Androsterone sulphate, DHEA sodium sulphate, 17-hydroxyepipregnanolone, 19-hydroxyandrostendione, 17-hydroxyprogesterone, Cortisol

Sensitivity 2.9 pg/ml (range 12.21 – 50,000 pg/ml)

Species Reactivity Species independent

Technical Info / Product Notes **Cited samples:**
[Wildlife Endocrinology: Cited Sample Types](#)

Wavelength 405 nm



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)