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: 37

View Online »

Ordering Information

Order Online »

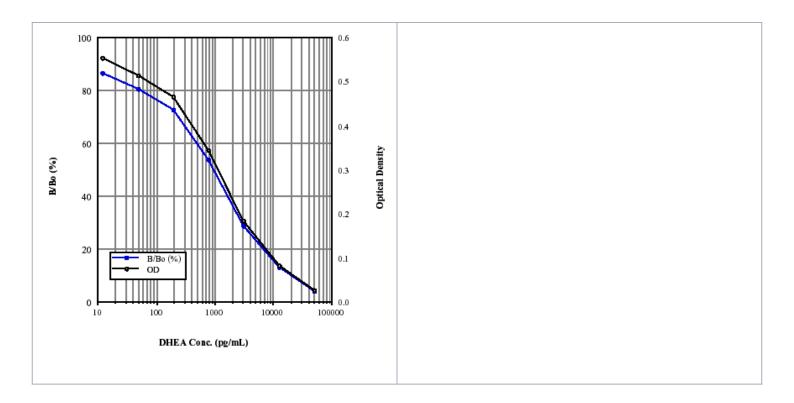
ADI-900-093

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 NotesFor 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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com