

Corticosterone ELISA kit

Highly sensitive corticosterone ELISA kit, species independent, used for animal stress research.

The Corticosterone ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. The Corticosterone ELISA kit has low cross-reactivity with related steroids. The broad dynamic range makes this kit ideal for a wide variety of sample matrices from any species.

Corticosterone is a glucocorticoid secreted by the cortex of the adrenal gland. Corticosterone is produced in response to the stimulation of the adrenal cortex by adrenocorticotrophic hormone (ACTH) and is the precursor of aldosterone. Corticosterone production has been shown to increase with stress, with elevated corticosterone levels being associated with impairment of long term memory retrieval and following burn injury. In addition to stress levels, corticosterone is believed to play a decisive role in sleep-wake patterns.

Citations: 500

[View Online »](#)

Ordering Information

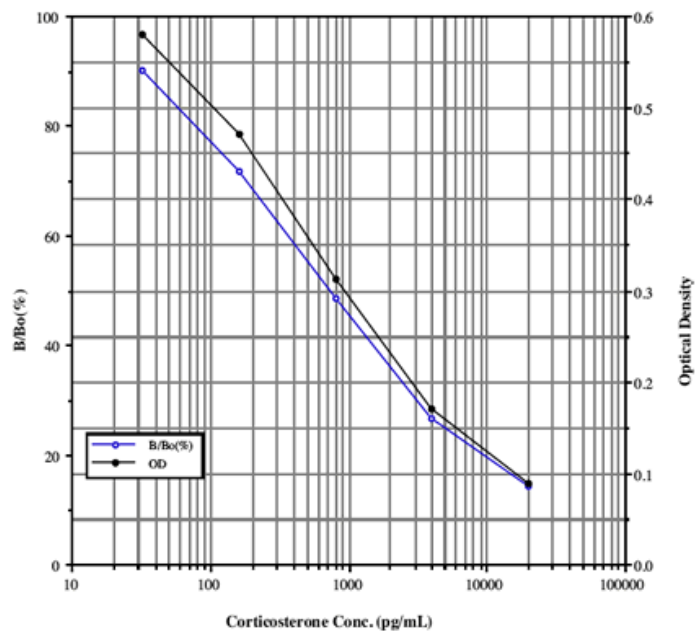
[Order Online »](#)

ADI-901-097	5x96 wells
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of Corticosterone, detecting as little as 27 pg/ml
- High throughput format with results in 3 hours for up to 39 samples
- Broad dynamic range
- Species independent
- Routinely used to monitor stress levels in zoo animals
- Save with bulk kit package!



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of Corticosterone in culture supernatants, feces, plasma, serum, and saliva from any species. Cited sample types include cell lysate, urine and whole blood.

Assay Time 3 hours

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

Contents DxS Microtiter plate, Conjugate, Antibody, Assay buffer 15 concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution, Steroid displacement reagent

Crossreactivity Corticosterone (100%), Deoxycorticosterone (21.3%), Desoxycorticosterone (21.0%), Progesterone (0.46%), Testosterone (0.31%), Tetrahydrocorticosterone (0.28%), Aldosterone (0.18%), Cortisol (0.046%) and <0.03%: Pregnenolone, Estradiol, Cortisone, 11-dehydrocorticosterone acetate

Sensitivity 27.0 pg/ml (range 32 – 20,000 pg/ml)

Species Reactivity Species independent

Technical Info / Product Notes **Cited samples:**
[Hormone ELISAs: 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)