

Cortisol ELISA kit

Widely used sensitive cortisol ELISA kit with a large variety of sample types and highly reproducible results for stress, autoimmune disease and cancer research.

The Cortisol EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. This kit is highly cited in peer-reviewed publications. The ready-to-use liquid color-coded reagents provided are safe, non-radioactive, and reduce error in the lab. The broad dynamic range lets you accurately measure Cortisol levels in a variety of sample matrices.

Cortisol (hydrocortisone, compound F) is a steroid hormone synthesized from cholesterol. It is the primary glucocorticoid produced and secreted by the adrenal cortex. Cortisol is found in the blood either as free cortisol, or bound to corticosteroid-binding globulin (CBG). Serum levels are highest in the early morning and decrease throughout the day. Cortisol is known to regulate metabolism, promoting gluconeogenesis, liver glycogen deposition, and the reduction of glucose utilization. Cortisol production is positively correlated with stress, resulting in increased blood pressure, and exerts immunosuppressive effects on the immune system.

Citations: 168

[View Online »](#)

Ordering Information

[Order Online »](#)

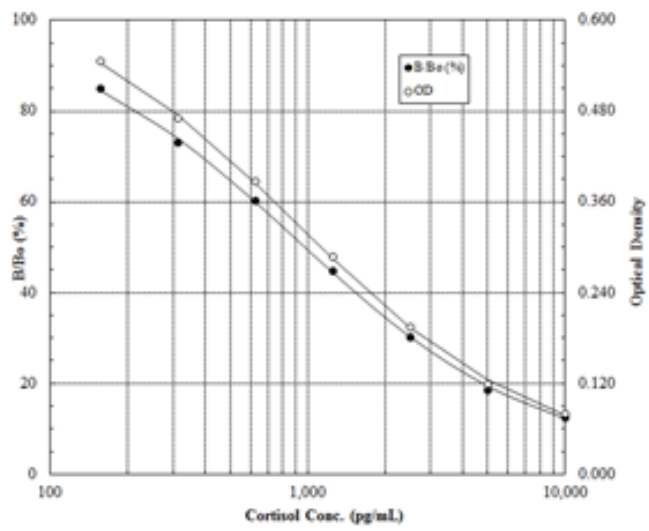
ADI-901-071

5x96 wells

Manuals, SDS & CofA

[View Online »](#)

- Sensitive measurement of cortisol, detecting as little as 56.72 pg/ml
- Highly cited in peer reviewed literature
- Reproducible results day-after-day and lot-after-lot
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Easy-to-use pre-coated plates and liquid color-coded reagents save time and reduce errors
- 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 Cortisol in culture supernatants, plasma, serum, saliva, and urine from any species. Cited sample type includes feces.

Assay Time 3 hours

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

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

Crossreactivity Cortisol (100%), Prednisolone (122.35%), Corticosterone (27.68%), 11-deoxycortisol (4.0%), Progesterone (3.64%), Prednisone (0.85%), Testosterone (0.12%) and <0.10%: Androstenedione, Cortisone, Estradiol

Sensitivity 56.72 pg/ml (range 156 – 10,000 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-
fr@enzolifesciences.com](mailto:info-fr@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)