

Progesterone ELISA kit

Highly specific and sensitive ELISA kit with results in < 3h

The Progesterone EIA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. This non-radioactive enzyme immunoassay is safe and reliable with ready-to-use liquid color-coded reagents to reduce errors.

Progesterone is the major female sex hormone, and is responsible for reproductive-related activities such as breast glandular development, the endometrial aspects of the menstrual cycle, and the establishment and maintenance of pregnancy. In addition, progesterone also directs pregnancy support physiology including changes in carbohydrate, protein and lipid metabolism, thermoregulation, sodium reabsorption in renal tubules and the reduction of alveolar and atrial carbon dioxide partial pressures (pCO_2). Progesterone is secreted in large amounts by the corpus luteum and by the extracellular conversion of cholesterol, cholesteryl esters, adrenal steroids, pregnenolone and pregnenolone sulfate. Small quantities are also secreted directly from the adrenal glands.

Citations: 135

[View Online »](#)

Ordering Information

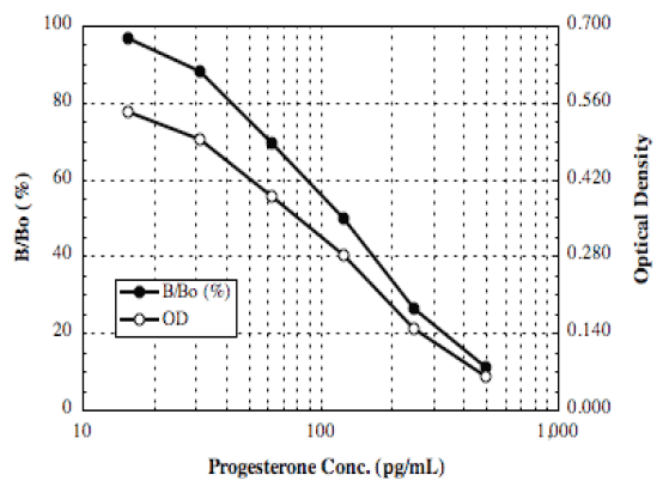
[Order Online »](#)

ADI-900-011	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- High sensitivity, measuring as little as 8.57 pg/ml of progesterone
- Highly specific with negligible cross-reactivity to other steroids
- High throughput format with results in < 3 hours for up to 38 samples in duplicate
- Easy-to-use liquid color-coded reagents reduce errors



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 progesterone in culture supernatants, serum, and saliva from any species. Cited sample type includes plasma.

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

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

Sensitivity 8.57 pg/ml (range 15.62 – 500 pg/ml)

Species Reactivity Species independent

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

Application Note:
[Validation of Two Steroid Hormone ELISA Kits with Ovine Blood Samples.](#)

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)