

# 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: 139

[View Online »](#)

## Ordering Information

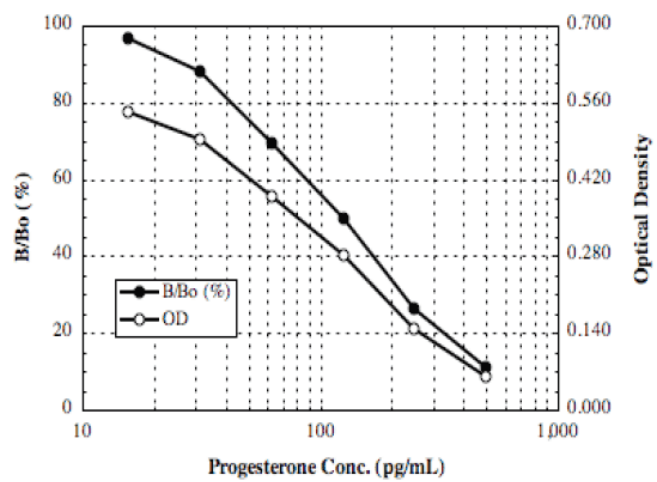
[Order Online »](#)

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

Last modified: May 29, 2024



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)