

Estriol ELISA kit

Certified for a large variety of sample types for steroid and hormone research.

The Estriol EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. This kit has a broad dynamic range suitable for a wide variety of sample matrices and the ability to < 60 pg/mL estriol.

Estriol is an important steroid hormone produced in large quantities during pregnancy. It is produced by the feto-placental unit and accounts for 90% of all estrogens produced during pregnancy.

Citations: 2

[View Online »](#)

Ordering Information

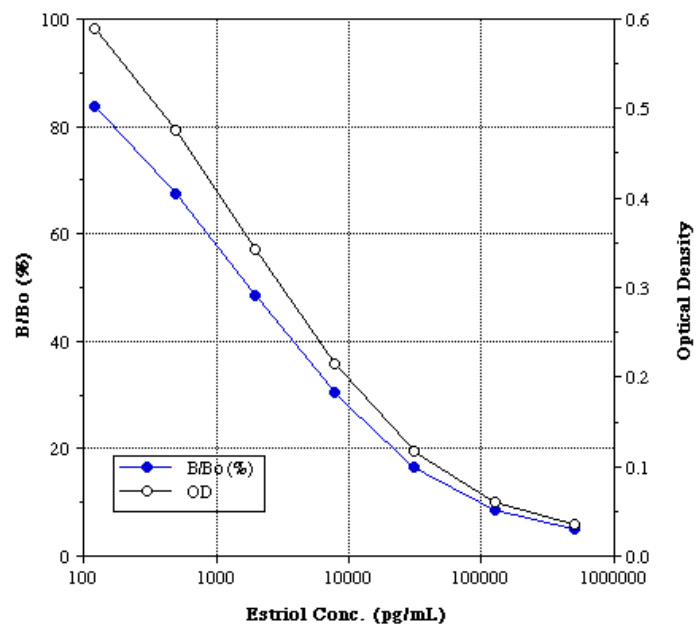
[Order Online »](#)

ADI-900-100	96 wells
-------------	----------

Manuals, SDS & CofA

[View Online »](#)

- Fast - results in from up to 37 samples in duplicate in just 3 hours
- Flexible - broad dynamic range suitable for a wide variety of samples
- Sensitive - measure < 60 pg/ml Estriol



Long Term Storage +4°C

Regulatory Status RUO - Research Use Only

Application	Colorimetric detection, ELISA
--------------------	-------------------------------

Assay Time 3 hours

Contents	GxR IgG Microtiter plate, Assay buffer concentrate, Conjugate, Antibody, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution
-----------------	---

Sensitivity 59.6 pg/ml (range 122 – 500,000 pg/ml)

Species Reactivity Species independent

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-
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@enzolifesciences.com)

For local distributors and detailed product information visit us online: enzo.com All product names, logos, and brands are the property of their respective owners.