

17 β -Estradiol ELISA kit

Highly sensitive ELISA, measuring as little as 28.5 pg/ml of 17 β -Estradiol in < 3 hours

The 17 β -Estradiol ELISA kit is a colorimetric competitive enzyme immunoassay kit with results in < 3 hours. Absorbance is read at 405 nm. This kit offers specific and save non-radioactive quantification of estradiol with ready-to-use liquid color-coded reagents to reduce error.

Biologically active estrogen, 17 β -estradiol, is an important sex hormone, but also makes an important contribution to the high concentrations of estrogens, which are present in malignant breast tissues.

Citations: 55

[View Online »](#)

Ordering Information

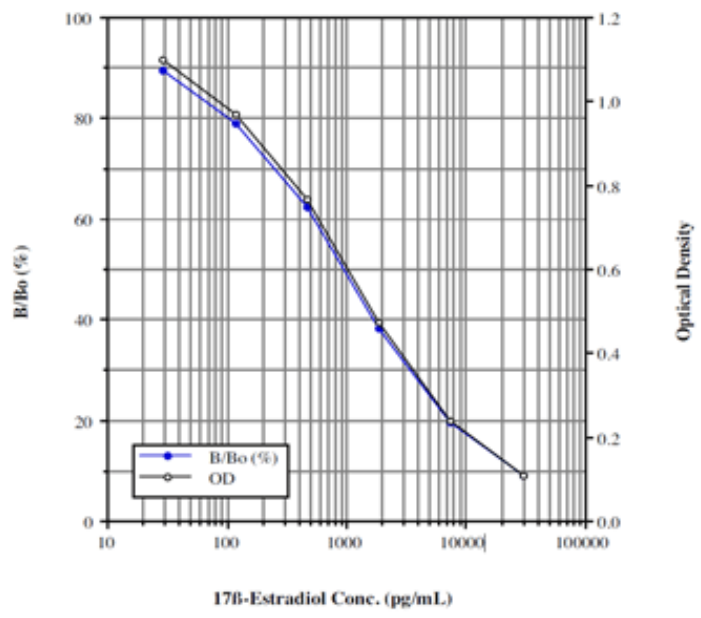
[Order Online »](#)

ADI-901-008	5x96 wells
-------------	------------

Manuals, SDS & CofA

[View Online »](#)

- Specific with low cross-reactivity to related steroids
- Safe and non-radioactive enzymatic assay
- Simple and color-coded reagents reduces error
- Reliable and reproducible results day-after-day and lot-after-lot
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Save with bulk kit package!



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Estrogen

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of 17 β -Estradiol in culture supernatants and saliva from any species. For serum and plasma use ADI-901-174.

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

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

Sensitivity 28.5 pg/ml (range 29.3 – 30,000 pg/ml)

Species Reactivity Species independent

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

Wavelength 405 nm

Last modified: August 19, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com