

17 β -Estradiol high sensitivity ELISA kit

Highly sensitive ELISA, detecting as little as 14 pg/ml of 17 β -Estradiol

The 17 β -Estradiol (serum/plasma) EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours.

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: 81

[View Online »](#)

Ordering Information

[Order Online »](#)

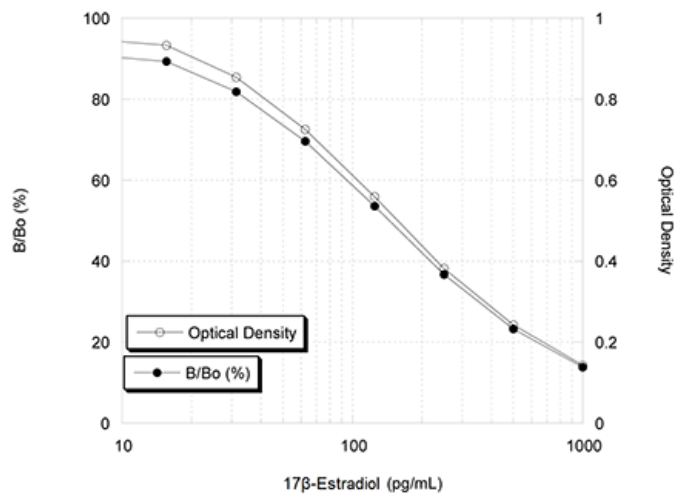
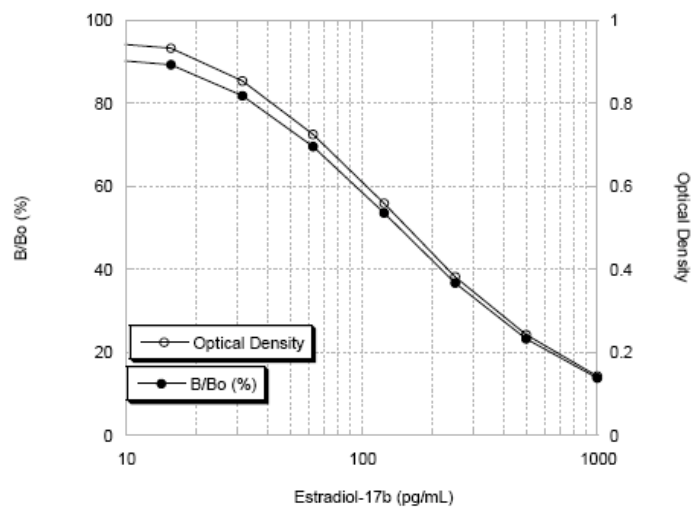
ADI-900-174

96 wells

Manuals, SDS & CofA

[View Online »](#)

- Highly sensitive measurement, as little as 14 pg/ml
- Fast and reliable results from up to 37 samples in duplicate in just 3 hours
- Fully quantitative results that surpass semi-quantitative Western blot analysis
- Convenient pre-coated plate and color-coded reagents reduces errors



Handling & Storage

Use/Stability	The 17 β -Estradiol Antibody and Conjugate are stable at -20°C. All other components of this kit are stable at +4°C until the kit's expiration date. The Antibody and Conjugate should be stored at -20°C upon receipt.
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 plasma and serum from any species.
Assay Time	3 hours
Compatibility	This product is compatible with the Absorbance 96 Plate Reader .
Contents	DxS IgG Microtiter plate, Conjugate, Antibody, Assay buffer, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution
Crossreactivity	17 β -Estradiol (100%), Estrone (17.8%), Estriol (0.9%), DHT (0.12%), Ethynylestradiol-7 α (0.043%), Dehydroisoandroesterone (0.38%), Estradiol-17 α (0.032%), Testosterone (0.01%), Progesterone (<0.001%), Danazol (<0.001%)
Sensitivity	14.0 pg/ml (range 15.6 – 1,000 pg/ml)
Species Reactivity	Species independent
Technical Info / Product Notes	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-
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-uk@enzolifesciences.com)