

# 17 $\beta$ -Estradiol high sensitivity ELISA kit

Highly sensitive ELISA, detecting as little as 14 pg/ml of 17 $\beta$ -Estradiol

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

Biologically active estrogen, 17 $\beta$ -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 »](#)

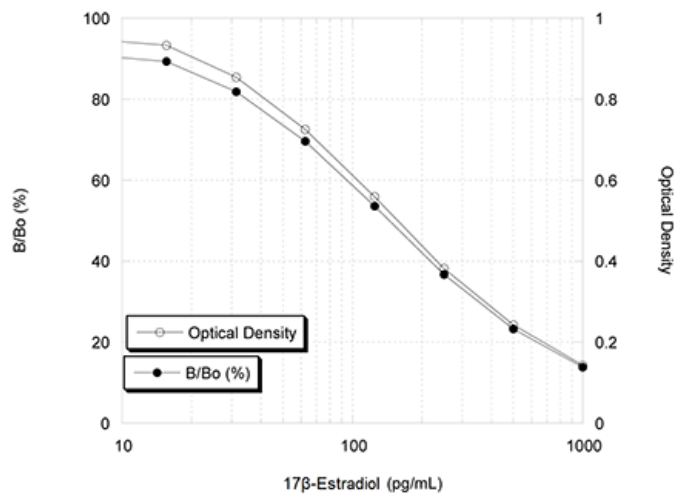
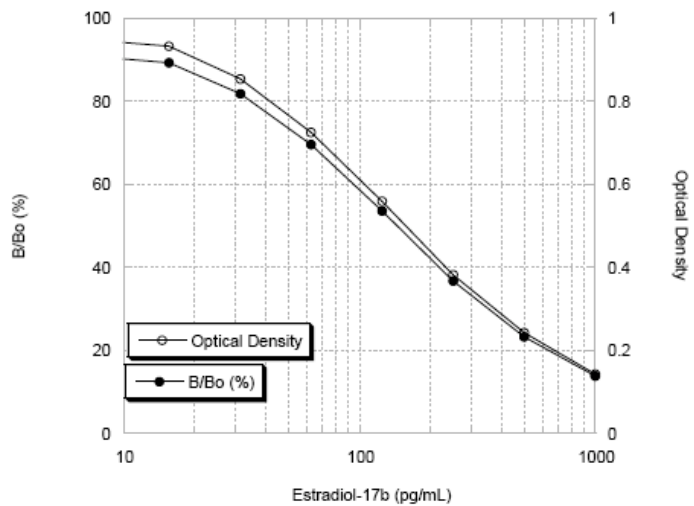
ADI-901-174

5x96 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
- Save with bulk kit package!



## Handling & Storage

Use/Stability	The 17 $\beta$ -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 $\beta$ -Estradiol in plasma and serum from any species.
Assay Time	3 hours
Compatibility	This product is compatible with the <a href="#">Absorbance 96 Plate Reader</a> .
Contents	DxS Microtiter plate, Conjugate, Antibody, Assay buffer concentrate, Wash buffer concentrate, Standard, pNpp Substrate, Stop solution
Crossreactivity	17 $\beta$ -Estradiol (100%), Estrone (17.8)%, Estriol (0.9%), DHT (0.12%), Ethynylestradiol-7 $\alpha$ (0.043%), Dehydroisoandroesterone (0.38%), Estradiol-17 $\alpha$ (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	<b>Cited samples:</b> <a href="#">Wildlife Endocrinology: Cited Sample Types</a>
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)