## 17β-Estradiol high sensitivity ELISA kit

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

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: 79

View Online »

**Ordering Information** 

Order Online »

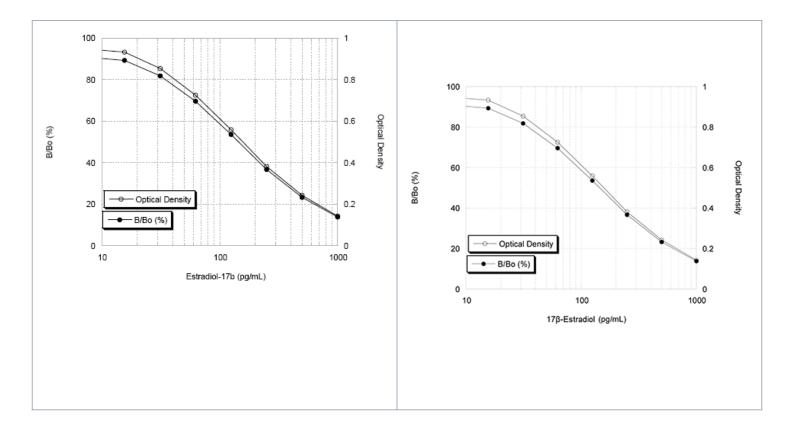
ADI-900-174

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 17β-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 17β-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** 17β-Estradiol (100%), Estrone (17.8%), Estriol (0.9%),

DHT (0.12%), Ethynylestradiol- $7\alpha$  (0.043%),

Dehydroisoandroesterone (0.38%), Estradiol-17α

(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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com