17β-Estradiol high sensitivity ELISA kit

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

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

Biologically active estrogen, 17β -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 »

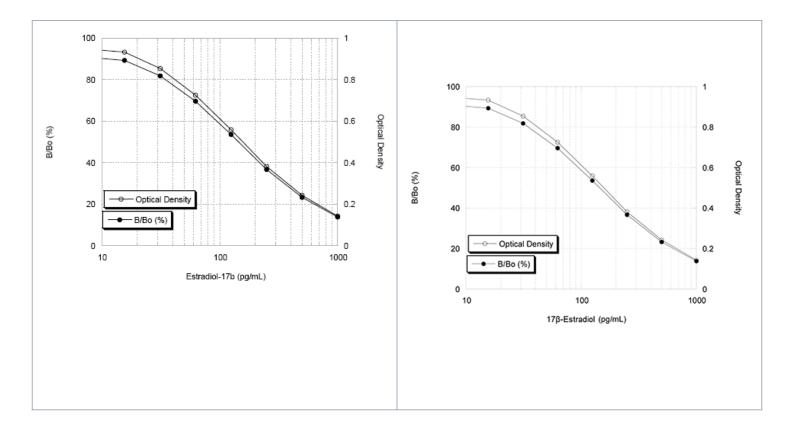
ADI-901-174

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 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 Microtiter plate, Conjugate, Antibody, Assay buffer

concentrate, Wash buffer concentrate, Standard, pNpp

Substrate, Stop solution

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

DHT (0.12%), Ethynylestradiol- 7α (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