17β-Estradiol ELISA kit

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

The 17β -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, 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: 56

View Online »

Ordering Information

Order Online »

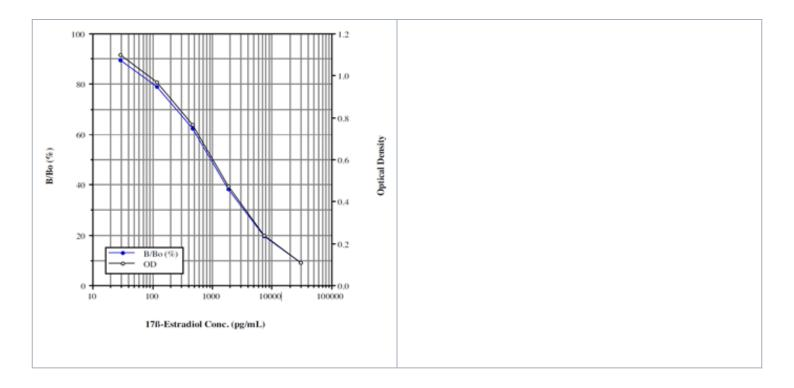
ADI-901-008

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 NotesFor the quantitative determination of 17β-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

