Estriol ELISA kit

Certified for a large variety of sample types for steroid and hormone research.

The Estriol EIA kit is a colorimetric competitive enzyme immunoassay kit with results in 3 hours. Absorbance is read at 405 nm. This kit has a broad dynamic range suitable for a wide variety of sample matrices and the ability to < 60 pg/mL estriol.

Estriol is an important steroid hormone produced in large quantities during pregnancy. It is produced by the feto-placental unit and accounts for 90% of all estrogens produced during pregnancy.

Citations: 2

View Online »

Ordering Information

Order Online »

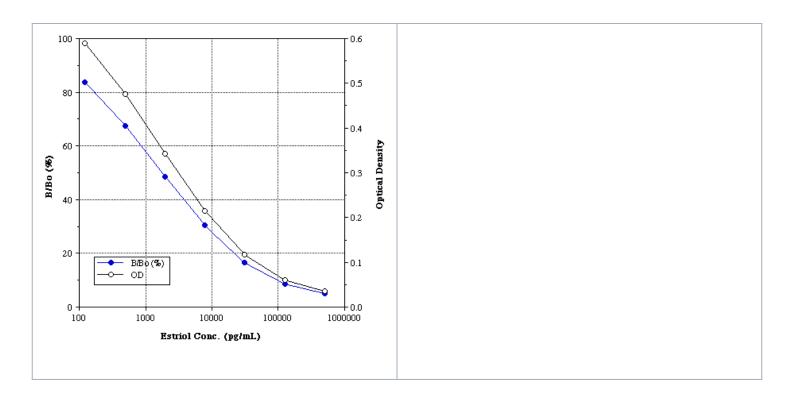
ADI-900-100

96 wells

Manuals, SDS & CofA

View Online »

- Fast results in from up to 37 samples in duplicate in just 3 hours
- Flexible broad dynamic range suitable for a wide variety of samples
- Sensitive measure < 60 pg/ml
 Estriol



Handling & Storage

Long Term Storage +4°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Application Colorimetric detection, ELISA

Application Notes For the quantitative determination of Estriol in culture

supernatants, plasma, serum, saliva, and urine from any

species.

Assay Time 3 hours

Compatibility This product is compatible with the Absorbance 96 Plate

Reader.

Contents GxR IgG Microtiter plate, Assay buffer concentrate,

Conjugate, Antibody, Wash buffer concentrate, Standard,

pNpp Substrate, Stop solution

Sensitivity 59.6 pg/ml (range 122 – 500,000 pg/ml)

Species Reactivity Species independent

Wavelength 405 nm

Last modified: May 29, 2024

