

Estrogen ELISA Kit ab285313

1 Image

Overview

Product name	Estrogen ELISA Kit
Detection method	Colorimetric
Sample type	Serum, Plasma, Other biological fluids, Tissue Homogenate
Assay type	Quantitative
Sensitivity	< 9.375 pg/ml
Range	15.625 pg/ml - 1000 pg/ml
Assay duration	Multiple steps standard assay
Product overview	<p>Estrogen ELISA Kit (ab285313, E4616) is a competitive ELISA assay for the quantitative measurement of estrogen in Human serum, plasma and cell culture supernatants. The density of color is proportional to the amount of Estrogen captured from the samples.</p> <p>Estrogen is the primary female sex hormone. It is responsible for the development and regulation of the female reproductive system and secondary sex characteristics. Estrogen may also refer to any substance, natural or synthetic, that mimics the effects of the natural hormone. The steroid 17-beta-estradiol is the most potent and prevalent endogenous estrogen, although several metabolites of estradiol also have estrogenic hormonal activity. Synthetic estrogens are used as part of some oral contraceptives, in estrogen replacement therapy for postmenopausal women, and in hormone replacement therapy for trans women.</p>
Platform	Pre-coated microplate (12 x 8 well strips)

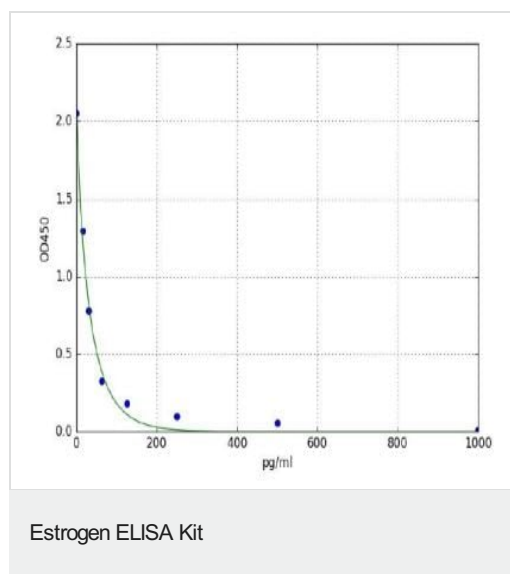
Properties

Storage instructionsStore at +4°C. Please refer to protocols.

Components	96 tests
25X Wash buffer	1 x 30ml
Antibody dilution buffer	1 x 10ml
Biotin/Detection antibody (Lyophilized)	1 vial

Components	96 tests
HRP-Streptavidin Conjugate (SABC)	1 x 120µl
Lyophilized Standard	2 vials
Micro ELISA Plate	1 unit
Plate Sealer	5 units
Purified water	1 x 200µl
SABC dilution buffer	1 x 10ml
Sample / Standard Dilution Buffer	1 x 20ml
Stop Solution	1 x 10ml
TMB Substrate	1 x 10ml

## Images



Typical Standard Curve: These standard curves are for demonstration only. A standard curve must be run with each assay

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

## Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you

- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

#### **Terms and conditions**

---

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors