

Estrogen BPA Environmental ELISA Kit ab175820

3 References 1 Image

Overview

Product name Estrogen BPA Environmental ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	overall	9			0.08%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	overall	9			0.8%

Sample type Urine, Serum, Plasma, Tissue

Assay type Competitive

Sensitivity = 3 pg/ml

Range 0.01 ng/ml - 1000 ng/ml

Assay duration Multiple steps standard assay

Product overview Abcam's Estrogen BPA Environmental competitive *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Estrogen BPA Environmental in urine, serum, plasma, cells and tissues following proper isolation and purification.

This competitive ELISA kit is based on competition between the Estrogen BPA Environmental epitope and the Estrogen BPA Environmental-HRP conjugate for a limited number of binding sites available from the anti- Estrogen BPA Environmental antibody; which is coated on the bottom of the wells of the 96 well ELISA plate. The conjugate concentration is held constant in each well, while the concentration of the Estrogen BPA Environmental is variable, based on the concentration of the sample or standard. Thus the amount of the Estrogen BPA Environmental conjugate which is able to bind to each of the wells is inversely proportional to the concentration of Estrogen BPA Environmental in the standard or sample. The amount of the conjugate which is bound to each well is then determined by the amount of color obtained, when TMB is added. The TMB reacts with the HRP available in the well. With the addition of sulfuric acid, the blue colored

product is converted into a yellow colored product, which can be read on a plate reader at 450 nm.

Estrogen BPA Environmental is a phenolic environmental estrogen¹ which disrupts endocrine activity. In human, a Estrogen BPA Environmental glucuronide was a primary metabolite of BPA. In a recent study, the age group with highest Estrogen BPA Environmental exposure was 6-11 years old with a mean total (free + glucuronidated) Estrogen BPA Environmental level of 4.33 ng/g of creatinine. Urinary Estrogen BPA Environmental levels were correlated with cardiovascular diseases and diabetes. A recent study revealed that a 12-ounce serving of canned soup for 5 days increased urinary Estrogen BPA Environmental level 12-fold due to Estrogen BPA Environmental -containing epoxy resin lining of the cans.

The free Estrogen BPA Environmental level in urine or cell culture media can be measured using the Estrogen BPA Environmental ELISA, without ethyl acetate extraction after 4-fold dilution of the sample. The glucuronidated Estrogen BPA Environmental level can also be measured without extraction after beta-glucuronidase treatment using our specific protocol.

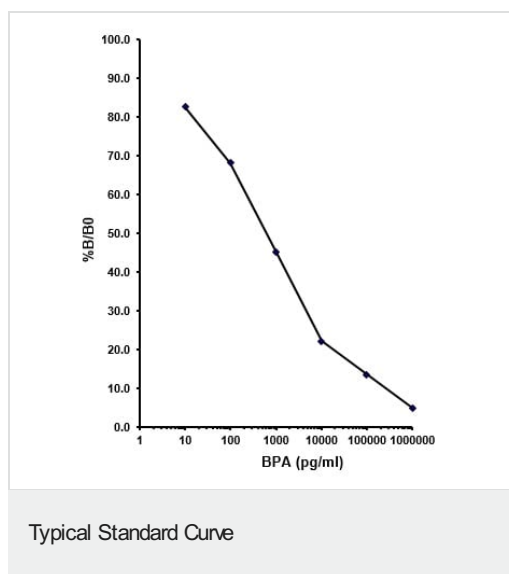
Platform Microplate

Properties

Storage instructions Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
BPA Standard	1 x 2µl
BPA-HRP Conjugates	1 x 12µl
HRP Buffer	1 x 15ml
Pre-coated 96-Well Microplate	1 unit
10X Sample Dilution Buffer	1 x 25ml
TMB Substrate	1 x 24ml
10X Wash Buffer Solution	1 x 25ml

Images



Example of standard curve obtained with ab175820.

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