



Estrone ELISA Kit

Catalog Number KA1908

96 assays

Version: 09

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	5
Precautions for Use	5
Assay Protocol	7
Reagent Preparation	7
Sample Preparation	7
Assay Procedure	8
Data Analysis	9
Calculation of Results	9
Performance Characteristics	10
Resources	13
References	13
Plate Layout	14

Introduction

Intended Use

The Estrone ELISA Kit is an enzyme immunoassay for the quantitative measurement of Estrone in serum and EDTA plasma.

Principle of the Assay

The Estrone ELISA Kit is a solid phase enzyme-linked immunosorbent assay (ELISA), based on the principle of competitive binding.


The microtiter wells are coated with a polyclonal rabbit antibody directed towards an antigenic site of the Estrone molecule. Endogenous Estrone of a sample competes with an estrone-horseradish peroxidase conjugate for binding to the coated antibody. After incubation the unbound conjugate is washed off.

The amount of bound peroxidase conjugate is inversely proportional to the concentration of Estrone in the sample. After addition of the substrate solution, the intensity of colour developed is inversely proportional to the concentration of Estrone in the sample.

General Information

Materials Supplied

List of component

Component	Detail	Amount
Microtiterwells	12x8 (break apart) strips, 96 wells, wells coated with anti-Estrone antibody (polyclonal).	96 (8x12) wells
Enzyme Conjugate	Ready to use. Estrone conjugated to horseradish peroxidase contains non-mercury preservative.	14 mL
Substrate Solution	Ready to use. Tetramethylbenzidine (TMB)	14 mL
Stop Solution	Ready to use, contains 0.5 M H ₂ SO ₄ <div style="text-align: center;">  </div> Hazards identification: H290 May be corrosive to metals. H314 Causes severe skin burns and eye damage.	14 mL
Wash Solution	Concentrated 40x.	30 mL

Standards and Controls - Ready to use

Component	Concentration (pg/mL)	Volume/Vial
Standard A	0	1 mL
Standard B	15	1 mL
Standard C	30	1 mL
Standard D	90	1 mL
Standard E	270	1 mL
Standard F	810	1 mL
Standard G	2400	1 mL
Control 1	For control values and ranges please refer to vial label or	1 mL
Control 2	QC-Report	1 mL

Conversion: pg/mL x 3.69 = pmol/L

Contents: contain non mercury preservative.

The standards are calibrated against certified reference material code E-075 (Cerilliant).

Note: Additional Standard A for sample dilution is available upon request.

Storage Instruction

When stored at 2°C - 8°C unopened reagents will retain reactivity until expiration date. Do not use reagents beyond this date.

Opened reagents must be stored at 2°C - 8°C. Microtiter wells must be stored at 2°C - 8°C. Once the foil bag

has been opened, care should be taken to close it tightly again.

Opened kits retain activity for 8 weeks if stored as described above.

Materials Required but Not Supplied

- ✓ A microtiter plate calibrated reader (450 ± 10 nm)
- ✓ Calibrated variable precision micropipettes.
- ✓ Absorbent paper.
- ✓ Distilled or deionized water.
- ✓ Timer.
- ✓ Semi logarithmic graph paper or software for data reduction

Precautions for Use

1. This kit is for research use only.
2. All reagents of this test kit which contain human serum or plasma have been tested and confirmed negative for HIV I/II, HBsAg and HCV by FDA approved procedures. All reagents, however, should be treated as potential biohazards in use and for disposal.
3. Before starting the assay, read the instructions completely and carefully. Use the valid version of protocol provided with the kit. Be sure that everything is understood.
4. The microplate contains snap-off strips. Unused wells must be stored at 2-8°C in the sealed foil pouch and used in the frame provided.
5. Pipetting of samples and reagents must be done as quickly as possible and in the same sequence for each step.
6. Use reservoirs only for single reagents. This especially applies to the substrate reservoirs. Using a reservoir for dispensing a substrate solution that had previously been used for the conjugate solution may turn solution colored. Do not pour reagents back into vials as reagent contamination may occur.
7. Mix the contents of the microplate wells thoroughly to ensure good test results. Do not reuse microwells.
8. Do not let wells dry during assay; add reagents immediately after completing the rinsing steps.
9. Allow the reagents to reach room temperature (21-24°C) before starting the test. Temperature will affect the absorbance readings of the assay. However, values for the samples will not be affected.
10. Never pipet by mouth and avoid contact of reagents and specimens with skin and mucous membranes.
11. Do not smoke, eat, drink, or apply cosmetics in areas where specimens or kit reagents are handled.
12. Wear disposable latex gloves when handling specimens and reagents. Microbial contamination of reagents or specimens may give false results.
13. Handling should be done in accordance with the procedures defined by an appropriate national biohazard safety guideline or regulation.
14. Do not use reagents beyond expiry date as shown on the kit labels.
15. All indicated volumes have to be performed according to the protocol. Optimal test results are only

obtained when using calibrated pipettes and microtiterplate readers.

16. Do not mix or use components from kits with different lot numbers. It is advised not to exchange wells of different plates even of the same lot. The kits may have been shipped or stored under different conditions and the binding characteristic of the plates may result slightly different.
17. Avoid contact with Stop Solution containing 0.5 M H₂SO₄. It may cause skin irritation and burns.
18. Some reagents contain Proclin 300, BND and MIT as preservatives. In case of contact with eyes or skin, flush immediately with water.
19. TMB substrate has an irritant effect on skin and mucosa. In case of possible contact, wash eyes with an abundant volume of water and skin with soap and abundant water. Wash contaminated objects before reusing them. If inhaled, take the person to open air.
20. Chemicals and prepared or used reagents have to be treated as hazardous waste according to the national biohazard safety guideline or regulation.
21. For information on hazardous substances included in the kit please refer to Material Safety Data Sheet (MSDS). Material Safety Data Sheet for this product is available upon request directly from the manufacturer.

✓ Limitations of use

Reliable and reproducible results will be obtained when the assay procedure is performed with a complete understanding of the package insert instruction and with adherence to good laboratory practice. Any improper handling of samples or modification of this test might influence the results.

✓ Interfering Interferences

Hemoglobin (up to 4 mg/mL), Bilirubin (up to 0.5 mg/mL) and Triglyceride (up to 7.5 mg/mL) have no influence on the assay results.

✓ Drug interferences

Until today no substance (drugs) are known to us, which have an influence to the measurement of estrone in a sample.

✓ High-Dose-Hook effect

No hook effect was observed in this test.

✓ Disposal of the Kit

The disposal of the kit must be made according to the national regulations. Special Information for this product is given in the Material Safety Data Sheet.

✓ Damaged Test Kits

In case of any severe damage to the test kit or components, the manufacturer has to be informed in writing, at the latest, one week after receiving the kit. Severely damaged single components should not be used for a test run. They have to be stored until a final solution has been found. After this, they should be disposed according to the official regulations.

Assay Protocol

Reagent Preparation

Bring all reagents and required number of strips to room temperature (21°C - 24°C) prior to use.

✓ Wash Solution

Add deionized water to the 40X concentrated Wash Solution.

Dilute 30 mL of concentrated Wash Solution with 1170 mL deionized water to a final volume of 1200 mL.

The diluted Wash Solution is stable for 2 weeks at room temperature.

Sample Preparation

Serum or EDTA plasma can be used in this assay.

Do not use haemolytic, icteric or lipaemic specimens.

Please note: Samples containing sodium azide should not be used in the assay.

✓ Sample Collection

- Serum

Collect blood by venipuncture (e.g. Sarstedt Monovette for serum), allow to clot, and separate serum by centrifugation at room temperature. Do not centrifuge before complete clotting has occurred.

- Plasma

Whole blood should be collected into centrifuge tubes containing anti-coagulant (e.g. Sarstedt Monovette with the appropriate plasma preparation) and centrifuged immediately after collection.

✓ Specimen Storage and Preparation

Specimens should be capped and may be stored for up to 4 days at 2 °C - 8 °C prior to assaying.

Specimens held for a longer time (up to 18 months) should be frozen only once at -20 °C prior to assay.

Thawed samples should be inverted several times prior to testing.

✓ Specimen Dilution

If in an initial assay, a specimen is found to contain more than the highest standard, the specimens can be diluted with Standard A and re-assayed as described in Assay Procedure.

For the calculation of the concentrations this dilution factor has to be taken into account.

Example:

a) dilution 1:10: 10 µL sample + 90 µL Standard A (mix thoroughly)

b) dilution 1:100: 10 µL dilution a) 1:10 + 90 µL Standard A (mix thoroughly).

Assay Procedure

- ✓ All reagents and samples must be allowed to come to room temperature before use. All reagents must be mixed without foaming.
 - ✓ The test should not be performed above a temperature of 24°C.
 - ✓ Once the test has been started, all steps should be completed without interruption.
 - ✓ Use new disposal plastic pipette tips for each standard, control or sample in order to avoid cross contamination.
 - ✓ Absorbance is a function of the incubation time and temperature. Before starting the assay, it is recommended that all reagents are ready, caps removed, all needed wells secured in holder, etc. This will ensure equal elapsed time for each pipetting step without interruption.
 - ✓ As a general rule the enzymatic reaction is linearly proportional to time and temperature.
-
- ✓ Each run must include a standard curve.
 1. Secure the desired number of Microtiter wells in the frame holder.
 2. Dispense 25 µL of each Standard, Control and samples with new disposable tips into appropriate wells.
 3. Dispense 100 µL Enzyme Conjugate into each well.

Thoroughly mix for 10 seconds. It is important to have a complete mixing in this step.
 4. Incubate for 60 minutes at room temperature.
 5. Briskly shake out the contents of the wells.

Rinse the wells 5 times with diluted Wash Solution (400 µL per well). Strike the wells sharply on absorbent paper to remove residual droplets.

Important note: The sensitivity and precision of this assay is markedly influenced by the correct performance of the washing procedure!
 6. Add 100 µL of Substrate Solution to each well.
 7. Incubate for 15 minutes at room temperature.
 8. Stop the enzymatic reaction by adding 50 µL of Stop Solution to each well.
 9. Determine the absorbance (OD) of each well at 450 ± 10 nm with a microtiter plate reader.

It is recommended that the wells be read within 10 minutes after adding the Stop Solution.

Data Analysis

Calculation of Results

- ✓ Calculate the average absorbance values for each set of standards, controls and samples.
- ✓ Using semi-logarithmic graph paper, construct a standard curve by plotting the mean absorbance obtained from each standard against its concentration with absorbance value on the vertical (Y) axis and concentration on the horizontal (X) axis.
- ✓ Using the mean absorbance value for each sample determine the corresponding concentration from the standard curve.
- ✓ Automated method: The results in the Instructions for Use have been calculated automatically using a 4 Parameter curve fit. (4 Parameter Rodbard or 4 Parameter Marquardt are the preferred method.) Other data reduction functions may give slightly different results.
- ✓ The concentration of the samples can be read directly from this standard curve. Samples with concentrations higher than that of the highest standard have to be further diluted or reported as > 2400 pg/mL. For the calculation of the concentrations this dilution factor has to be taken into account.

- ✓ Example of typical standard curve

The following data is for demonstration only and cannot be used in place of data generations at the time of assay.

Standard	Optical Units (450 nm)
Standard A (0 pg/mL)	2.24
Standard B (15 pg/mL)	2.04
Standard C (30 pg/mL)	1.88
Standard D (90 pg/mL)	1.48
Standard E (270 pg/mL)	1.05
Standard F (810 pg/mL)	0.68
Standard G (2400 pg/mL)	0.43

- ✓ Expected Normal Values

It is strongly recommended that each laboratory should determine its own normal and abnormal values.

In a study conducted with apparently normal healthy adults, using the Estrone ELISA Kit the following values are observed:

Population	n	Range (min.-max.) (pg/mL)	Mean (pg/mL)	2.5 th – 97.5 th Percentile (pg/mL)	Median (pg/mL)
Males	60	9.0 – 79.1	47.6	15.6 – 77.0	47.9
Females					
premenopausal	110	11.2 – 338.3	69.4	15.5 – 220.2	52.8
postmenopausal	91	9.4 – 59.6	28.6	11.0 – 54.5	27.4

✓ Quality Control

Good laboratory practice requires that controls be run with each standard curve. A statistically significant number of controls should be assayed to establish mean values and acceptable ranges to assure proper performance.

It is recommended to use control samples according to state and federal regulations. The use of control samples is advised to assure the day to day validity of results. Use controls at both normal and abnormal levels.

The controls and the corresponding results of the QC-Laboratory are stated in the QC Report added to the kit. The values and ranges stated on the QC Report always refer to the current kit lot and should be used for direct comparison of the results.

Employ appropriate statistical methods for analysing control values and trends. If the results of the assay do not fit to the established acceptable ranges of control materials sample results should be considered invalid.

In this case, please check the following technical areas: Pipetting and timing devices; photometer, expiration dates of reagents, storage and incubation conditions, aspiration and washing methods.

After checking the above mentioned items without finding any error contact your distributor or the manufacturer directly.

Performance Characteristics

✓ Assay Dynamic Range

The range of the assay is between 8.1 - 2400 pg/mL.

✓ Specificity of Antibodies (Cross Reactivity)

The following substances were tested for cross reactivity of the assay.

Compound	Cross Reactivity (%)
17-OH-Progesterone	0.01
Androstenedione	0.04
Corticosterone	0.01
Cortisone	0.39
DHEA	0.01
Estradiol	1.19
Estriol	0.07
Estrone-3-β-D-glucuronide	0.35
Estrone-3-sulfate	0.44
Ethisterone	0.39
Hydrocortisone	0.30
Progesterone	< 0.01
Testosterone	0.03

✓ Sensitivity

The analytical sensitivity of the ELISA was calculated by subtracting 2 standard deviation from the mean of 20 replicate analyses of the Zero Standard (Standard A) and was found to be 8.1 pg/mL.

The limit of detection (LoD) of the assay is 14.6 pg/mL

The limit of quantification (LoQ) of the assay is 15.8 pg/mL.

✓ Reproducibility

• Intra-assay

The within assay variability is shown below:

Sample	n	Mean (pg/mL)	CV (%)
1	10	57.68	8.4
2	10	61.00	6.7
3	10	82.14	6.8
4	10	409.13	6.4

• Inter-assay

The between assay variability is shown below:

Sample	n	Mean (pg/mL)	CV (%)
1	30	32.98	11.9
2	30	61.91	9.7
3	30	81.96	7.4
4	30	451.6	8.7

• Inter-Lot

The inter-assay (between-lots) variation was determined by repeated measurements of 3 samples in 3 different kit lots.

Sample	n	Mean (pg/mL)	CV (%)
1	18	2005.49	4.8
2	18	1122.06	2.3
3	18	66.28	14.0
4	18	53.46	8.4

✓ Recovery

Samples have been spiked by adding Estrone solutions with known concentrations in a 1:1 ratio.

The % recovery has been calculated by multiplication of the ratio of the measurements and the expected values with 100 (expected value = (endogenous Estrone + added Estrone)/ 2; because of a 1:2 dilution of serum with spike material).

Sample	1	2	3	4	5	6	
Sample Type	Serum	EDTA	EDTA	Serum	Serum	Serum	
Concentration [pg/mL]	31.95	32.06	33.83	46.08	464.78	639.70	
Average Recovery [%]	102.0	103.4	93.3	108.0	105.6	92.4	
Range of Recovery [%]	from	96.0	96.1	85.7	101.7	95.9	89.2
	to	104.2	109.4	110.2	114.3	114.2	95.4

✓ Linearity

Sample	1	2	3	4	5	
Sample Type	EDTA	EDTA	Serum	Serum	Serum	
Concentration [pg/mL]	92.42	210.70	496.00	547.96	1649.26	
Average Recovery	110.3	101.4	106.0	96.6	101.6	
Range of Recovery [%]	from	108.4	95.7	88.9	85.0	95.7
	to	112.1	105.9	113.4	109.0	105.2

✓ Comparison Studies

For method comparison, 40 samples spanning the complete assay range were tested by the Estrone ELISA Kit and by the RIA.

The resulting regression equation was:

$$\text{ELISA} = 0.743 \text{ RIA} + 15.52; R^2 = 0.990$$

Resources

References

1. Resnik R, Killam AP, Battaglia FC et al: The stimulation of uterine blood flow by various estrogens. *Endocrinology*. 1974 Apr; 94(4):1192-6.
2. Fayman C, Winter JSD, Reyes FI. Patterns of gonadotropins and gonadal steroids throughout life. *Clin. Obstet. Gynecol*. 1976; 3:467-483.
3. Baird DT, Fraser IS. Blood production and ovarian secretion rates of estradiol-17beta and estrone in women throughout the menstrual cycle. *J Clin Endocrinol. Metab*. 1974; 38:1009-1017.
4. Lindbert BS, Johansson EDB, Nilsson BA: Plasma levels of non-conjugated oestrone, oestradiol-17b and oestriol during uncomplicated pregnancy. *Acta Obstet Gynecol Scand*. 1974; 32:21.
5. Drafta D, Schindler AE, Stroe EW, Neacsu E. Age-related changes of plasma steroids In normal adult males. *J Steroid Blochem*. 1982; 17:683-687.
6. Devane GW, Czekala NM, Judd HL, Yen SSC. Circulating gonadotropins, estrogens, and androgens in polycystic ovarian disease. *Am J Obstet Gynecol*. 1975; 121:496.
7. Kushnir MM et al. High-sensitivity tandem mass spectrometry assay for serum estrone and estradiol. *Am J Clin Pathol*. 2008; 129:530-9.
8. Jasuja GK et al. Age trends in estradiol and estrone levels measured using liquid chromatography tandem mass spectrometry in community-dwelling men of the Framingham Heart Study. *J Gerontol A Biol Sci Med Sci*. 2013; 68:733-40.
9. Kaaks R, Lukanova A, Kurzer MS. Obesity, endogenous hormones, and endometrial cancer risk: a synthetic review. *Cancer Epidemiol Biomarkers Prev*. 2002; 11(12):1531-43.
10. Kuiper GG, et al. Comparison of the ligand binding specificity and transcript tissue distribution of estrogen receptors alpha and beta. *Endocrinology*. 1997; 13B:863-B70.

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H