



Estrone ELISA Assay Kit

Catalog Number: ESN31-K01

For Research Use Only. Not for use in diagnostic procedures.

v. 13 (10 OCT 23)

EAGLE BIOSCIENCES, INC.
20A Northwest Blvd., Suite 112, Nashua, NH 03063
Phone: 617-419-2019 Fax: 617-419-1110
WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Estrone ELISA Assay Kit (enzyme-linked immunoassay kit) is intended for the quantitative measurement of Estrone in human serum. The Eagle Biosciences Estrone ELISA Assay Kit is for research use only and not to be used in diagnostic procedures.

For further information about this kit, its application, or the procedures in this insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.

LIMITATIONS RELATED TO INTENDED PURPOSE

1. This test is not intended to be used for screening purposes.
2. This test is not intended for home testing or self-testing
3. The kit is calibrated for the determination of estrone in human serum. The kit is not calibrated for the determination of pregnenolone in other specimens of human or animal origin.
4. The results obtained with this kit shall not be used for a clinical diagnosis and for therapeutic decisions.
5. Although common interfering substances have been evaluated with this test, other substances that have not been evaluated such as drugs and the occurrence of heterophilic antibodies in individuals regularly exposed to animals or animal products have the potential of causing interferences.

INTRODUCTION

Estrone is a steroid, a female sex hormone and, with estriol and estradiol, one of the three most important endogenous estrogens. Estrogens are involved in the development of female sex organs and secondary sex characteristics. Before the ovum is fertilized the main action of the estrogens is on the growth and function of the reproductive tract in order to prepare it for the fertilized ovum. During the follicular phase of the menstrual cycle the estrone level shows a slight increase. The production of estrone then increases markedly to peak at around day 13. The peak is of short duration and by day 16 of the cycle levels will be low. A second peak occurs at around day 21 of the cycle and if fertilization does not occur, then the production of estrone decreases.

PRINCIPLE OF THE ASSAY

The Estrone ELISA is a competitive immunoassay. Competition occurs between an estrone present in calibrators, controls and specimen samples and an enzyme-labelled antigen (HRP conjugate) for a limited number of antibody binding sites on the microplate. After a washing step that removes unbound materials, the TMB substrate (enzyme substrate) is added which reacts with HRP to form a blue-colored product that is inversely proportional to the amount of estrone. Following an incubation, the enzymatic reaction is terminated by addition of the stopping solution, converting the color from blue to yellow. The absorbance is measured on a microplate reader at 450 nm. A set of calibrators is used to plot a calibrator curve from which the amount of estrone in specimen samples and controls can be directly read.



PROCEDURAL CAUTIONS AND WARNINGS

1. This kit is for use by trained laboratory personnel (professional use only). For laboratory *in vitro* use only.
2. Practice good laboratory practices when handling kit reagents and specimens. This includes:
 - a. Do not pipette by mouth.
 - b. Do not smoke, drink, or eat in areas where specimens or kit reagents are handled.
 - c. Wear protective clothing and disposable gloves.
 - d. Wash hands thoroughly after performing test.
 - e. Avoid contact with eyes, use safety glasses; in case of contact with eyes, flush eyes with water immediately and contact a doctor.
3. Users should have a thorough understanding of this protocol for the successful use of this kit. Reliable performance will only be attained by strict and careful adherence to the instructions provided.
4. Do not use this kit beyond the expiry date stated on the label.
5. If the kit reagents are visibly damaged, do not use the test kit.
6. Do not use kit components from different kit lots within a test and do not use any component beyond the expiration date printed on the label.
7. All kit reagents and specimens must be brought to room temperature and mixed gently but thoroughly before use. Avoid repeated freezing and thawing of specimens.
8. When the use of water is specified for dilution or reconstitution, use deionized or distilled water.
9. Immediately after use, each individual component of the kit must be returned to the recommended storage temperature stated on the label.
10. A calibrator curve must be established for every run.
11. It is recommended to all customers to prepare their own control materials or serum pools which should be included in every run at a high and low level for assessing the reliability of results.
12. The controls (included in kit) must be included in every run and their results must fall within the ranges stated in the quality control certificate; a failed control result might indicate improper reagent storage.
13. When dispensing the substrate and stopping solutions, do not use pipettes in which these liquids will come into contact with any metal parts.
14. The TMB Substrate is sensitive to light and should remain colorless if properly stored. Instability or contamination may be indicated by the development of a blue color, in which case it should not be used.
15. Do not use grossly hemolyzed, grossly lipemic, icteric or improperly stored serum.
16. Samples or controls containing azide or thimerosal are not compatible with this kit, they may lead to false results.
17. Serum samples with a known low estrone concentration (< 60 pg/mL) may be used to dilute serum samples with values higher than the highest calibrator. Otherwise, results may be reported as "> 2000 pg/mL". The use of any other reagent will lead to false results.
18. Avoid microbial contamination of reagents.
19. To prevent the contamination of reagents, use a new disposable pipette tip for dispensing each reagent, sample, calibrator, and control.



20. To prevent contamination of reagents, do not pour reagents back into the original containers.
21. Kit reagents must be regarded as hazardous waste and disposed of according to local and/or national regulations.
22. Consumables used with the kit that are potentially biohazardous (e.g., pipette tips, bottles or containers containing human materials) must be handled according to biosafety practices to minimize the risk of infection and disposed of according to local and/or national regulations relating to biohazardous waste.
23. This kit contains 1 M sulfuric acid in the stopping solution component. Do not combine acid with waste material containing sodium azide or sodium hypochlorite.
24. The use of safety glasses, and disposable plastic, is strongly recommended when manipulating biohazardous or bio-contaminated solutions.
25. Proper calibration of the equipment used with the test, such as the pipettes and absorbance microplate reader, is required.
26. If a microplate shaker is required for the assay procedure, the type and speed of shaker required is stated in the REAGENTS AND EQUIPMENT NEEDED BUT NOT PROVIDED section. Both the type and speed of shaker used can influence the optical densities and test results. If a different type of shaker and/or speed is used, the user is responsible for validating the performance of the kit.
27. Do not reuse the microplate wells, they are for SINGLE USE only.
28. To avoid condensation within the microplate wells in humid environments, do not open the pouch containing the microplate until it has reached room temperature.

SAFETY CAUTIONS AND WARNINGS

Biohazards

The reagents should be considered a potential biohazard and handled with the same precautions applied to blood specimens. All human specimens should be considered a potential biohazard and handled as if capable of transmitting infections and in accordance with good laboratory practices.

The calibrators and controls provided with the kit contain processed human serum/plasma that has been tested by approved methods and found to be negative for the presence of HBsAg and antibodies to HCV, HIV 1/2 and HIV NAT. However, no test method can offer complete assurance that any viable pathogens are absent. Therefore, these components should be considered a potential biohazard and handled with the same precautions as applied to any blood specimen, following good laboratory practices.

Chemical Hazards

Avoid direct contact with any of the kit reagents. Specifically avoid contact with the TMB Substrate (contains tetramethylbenzidine) and Stopping Solution (contains sulfuric acid). If contacted with any of these reagents, wash with plenty of water and refer to SDS for additional information.

SPECIMEN COLLECTION, STORAGE AND PRE-TREATMENT

Specimen Collection and Storage

Approximately 0.15 mL of serum is required per duplicate determination. Collect 4–5 mL of venous blood into an appropriately labelled tube and allow it to clot. Centrifuge at room temperature and carefully transfer the serum into a new storage tube or container. Serum



samples may be stored at room temperature for up to 3 days, at 2-8°C for up to 7 days hours or at -20°C or lower for up to 1 month. Consider all human specimens as possible biohazardous materials and take appropriate precautions when handling.

Specimen Pre-Treatment

Specimen pre-treatment is not required

REAGENTS AND EQUIPMENT NEEDED BUT NOT SUPPLIED

1. Calibrated single-channel pipette to dispense 50 µL.
2. Calibrated multi-channel pipette to dispense 50 µL, 100 µL, and 150 µL.
3. Calibrated multi-channel pipettes to dispense 350 µL (if washing manually).
4. Automatic microplate washer (recommended).
5. Disposable pipette tips.
6. Distilled or deionized water.
7. Calibrated absorbance microplate reader with a 450 nm filter and an upper OD limit of 3.0 or greater.

MATERIALS PROVIDED

Microplate – Ready to use

Contents	One anti-estrone polyclonal antibody-coated 96-well (12x8) microplate in a resealable pouch with desiccant.
Storage	2-8°C
Stability	Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

HRP Conjugate – Ready to use

Contents	One bottle containing Estrone-Horse Radish Peroxidase (HRP) conjugate in a protein-based buffer with a non-mercury preservative.
Volume	15mL/bottle
Storage	2-8°C
Stability	Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

Calibrators (A – F) – Ready to use

Contents	Six bottles of calibrator containing specified estrone concentrations. Human serum-based buffer with a non-mercury preservative. Prepared by spiking buffer with define quantities of estrone. Listed below are approximate concentrations, please refer to vial labels for exact concentrations. Concentrations: 0, 20, 60, 200, 600, 2000 pg/mL
Volume	1mL/bottle
Storage	2-8°C
Stability	Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

**Controls (1 – 2) – Ready to use**

Contents	Two bottles of control containing different estrone concentrations. Human serum-based buffer with a non-mercury preservative. Prepared by spiking buffer with defined quantities of pregnenolone. Refer to the QC certificate for the target values and acceptable ranges.
Volume	1mL/bottle
Storage	2-8°C
Stability	Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

TMB Substrate – Ready to use

Contents	One bottle containing tetramethylbenzidine and hydrogen peroxide in a non-DMF or DMSO containing buffer.
Volume	16mL/bottle
Storage	2-8°C
Stability	Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

Stopping Solution – Ready to use

Contents	One bottle containing 1M sulfuric acid.
Volume	6mL/bottle
Storage	2-8°C
Stability	Unopened: Stable until the expiry date printed on the label. After Opening: Stable for four weeks.

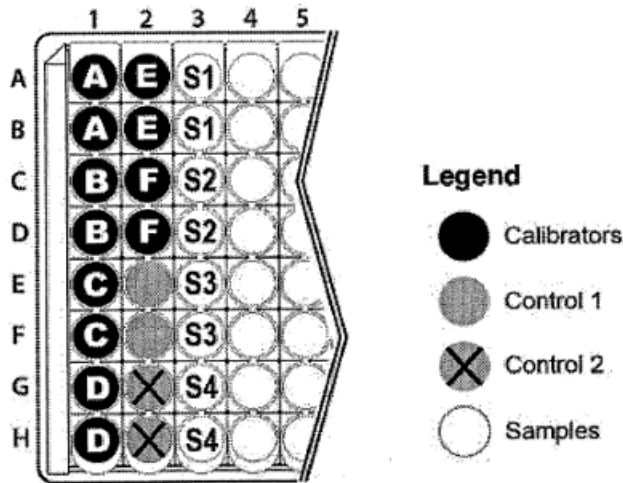
Wash Solution – Concentrated; Requires preparation

Contents	One bottle containing buffer with a non-ionic detergent and a non-mercury preservative.
Volume	50mL/bottle
Storage	2-8°C
Stability	Unopened: Stable until expiry date printed on the label. After Opening: Stable for four weeks. Following Preparation: The wash buffer working solution is stable for 2 weeks following preparation, assuming Good Laboratory Practices are adhered to. To prevent microbial growth, prepare the wash buffer working solution in a clean container and store under refrigerated conditions (2-8°C) when not in use.
Preparation	Dilute 1:10 in distilled or deionized water before use. If the whole microplate is to be used dilute 50 mL of the wash buffer concentrate in 450 mL of distilled or deionized water



RECOMMENDED ASSAY LAYOUT

RECOMMENDED ASSAY LAYOUT



ASSAY PROCEDURE

Specimen Pretreatment: None

All kit components, controls, and specimen samples must reach room temperature prior to use. Calibrators, controls, and specimen samples should be assayed in duplicate. Once the procedure has been started, all steps should be completed without interruption.

1. After all kit components have reached room temperature, mix gently by inversion.
2. Prepare the Wash Buffer Working Solution (See section Reagents Provided, HRP Conjugate Concentration, Wash Buffer Concentrate).
3. Plan the microplate wells to be used for calibrators, controls, and samples. See Recommended Assay Layout. Remove the strips from the microplate frame that will not be used and place them in the bag with desiccant. Reseal the bag with the unused strips and return it to the refrigerator.
4. Pipette 50 μ L of each calibrator, control, and pre-treated specimen sample into assigned wells.
5. Pipette 100 μ L of the HRP conjugate Working Solution into each well (the use of a multi-channel pipette is recommended).
6. Gently tap the microplate frame for 10 seconds to mix the contents of the wells and incubate the microplate at room temperature (no shaking) for 60 minutes.
7. Wash the microplate wells with an automatic microplate washer (preferred) or manually as state below.
 - a. Automatic: Using an automatic microplate washer, perform a 3-cycle wash using 350 μ L /well of Wash Buffer Working Solution (3 x 350 μ L). One cycle consists of aspirating all wells then filling each well with 350 μ L of Wash Buffer Working Solution. After the final wash cycle, aspirate all wells and then tap the microplate firmly against absorbent paper to remove any residual liquid.



- b. Manually: For manual washing, perform a 3-cycle wash using 350 μL /well of Wash Buffer Working Solution (3 x 350 μL). One cycle consists of aspirating all wells by briskly emptying the contents of the wells over a waster container, then pipetting 350 μL of Wash Buffer Working Solution into each well using a multi-channel pipette. After the final wash cycle, aspirate all wells by briskly emptying the contents over a waste container and then tap the microplate firmly against absorbent paper to remove any residual liquid.
8. Pipette 150 μL of TMB Substrate into each well (the use of a multi-channel pipette is recommended).
9. Incubate the microplate at room temperature (no shaking) for 20 minutes. Do not tap the microplate and avoid placing in intense light or air currents.
10. Pipette 50 μL of Stopping Solution into each well (the use of a multi-channel pipette is recommended) in the same order and speed as was used for the addition of TMB Substrate. Gently tap the microplate frame to mix the contents of the wells.
11. Measure the optical density (absorbance) in the microplate wells using an absorbance microplate reader set to 450 nm, within 20 minutes after addition of the Stopping Solution.

CALCULATIONS

1. Calculate the mean optical density of each calibrator, control, and specimen sample duplicate.
2. Use a 4-parameter or 5-parameter curve fit with immunoassay software to generate a calibrator curve.
3. The immunoassay software will calculate the concentrations of the controls and specimen samples using the mean optical density values and the calibrator curve.
4. If a sample reads more than 2000 pg/mL and needs to be diluted and retested, then dilute it with a known low estrone value (<60 pg/mL) serum sample not more than 1:10. The result obtained should be multiplied by the dilution factor.

QUALITY CONTROL

When assessing the validity of the test results, the following criteria should be evaluated:

1. The Calibrator A mean optical density meets the acceptable range as stated in the QC Certificate
2. The calibrator with the highest concentration meets the % binding acceptable range as stated in the QC Certificate. % Binding = (OD of calibrator / OD of calibrator A) X 100.
3. The values obtained for the kit controls are within the acceptable ranges as stated in the QC Certificate.
4. The results of any external controls that were used meet the acceptable ranges.

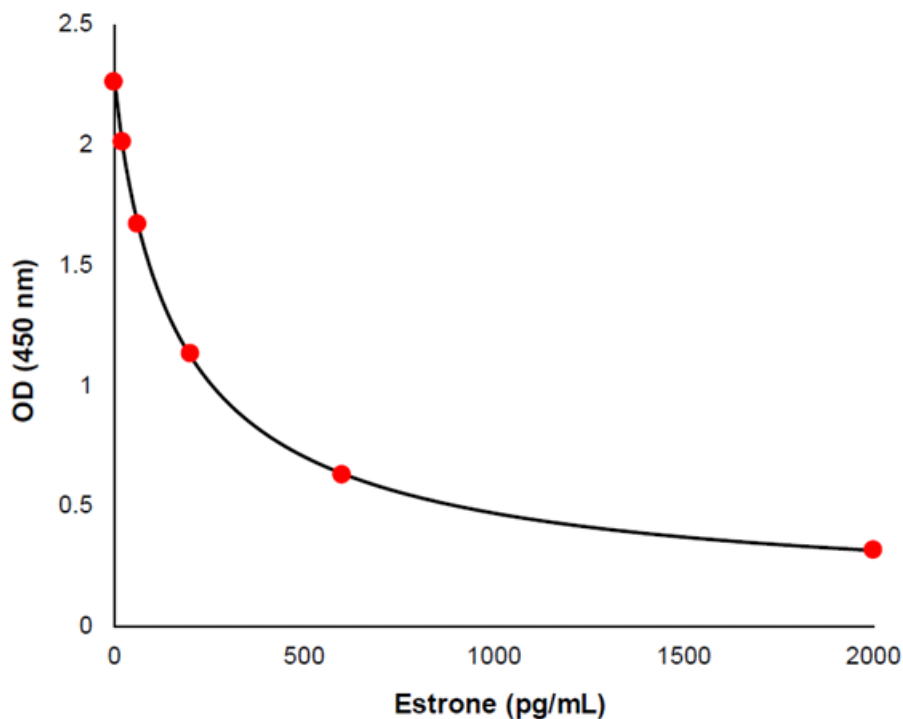


TYPICAL TABULATED DATA

Sample Data. Do not use to calculate results.

Calibrator	Mean OD	% Binding	Value (pg/mL)
A	2.261	100	0
B	2.012	89	20
C	1.671	75	60
D	1.132	50	200
E	0.632	28	600
F	0.319	14	2000
Unknown	1.481	-	96

TYPICAL CALIBRATOR CURVE



PERFORMANCE CHARACTERISTICS

Sensitivity

The analytical sensitivity study was performed according to the CLSI EP17-A2 guideline. The Limit of Background (LoB), limit of detection (LoD), limit of Quantitation (LoQ) are summarized in the table below.

Parameter	Estrone (pg/mL)
LoB	5.6
LoD	14.8
LoQ	17.7



SPECIFICITY (CROSS-REACTIVITY)

The following compounds were tested for cross-reactivity with the Direct Estrone ELISA kit with estrone cross-reacting at 100%:

Steroid	% Cross Reactivity
Estrone	100
11-Deoxycorticosterone	<0.1
17-Hydroxyprogesterone	<0.1
17 α -Estradiol	3.6
17 β -Estradiol	7.9
Aldosterone	<0.1
Androstenedione	<0.1
Androsterone	<0.1
Cholesterol	<0.01
Corticosterone	<0.1
Cortisol	0.2
Danazol	<0.1
DHEA	0.1
DHEAS	<0.1
DHT	<0.1
Equilin	19.1
Estradiol Sulfate	<2.9
Ethisterone	<0.1
Prednisone	<0.1
Pregnenolone Sulfate	<0.1
Progesterone	<0.1
Testosterone	<0.1
17 β -Estradiol-3-Glucuronide	0.14

INTERFERENCES

Potential interferents were spiked into human serum samples to determine the effect on the measured estrone values. Hemoglobin up to 10 g/L, Bilirubin conjugated up to 20 mg/dL, Bilirubin unconjugated up to 10mg/dL, Triglycerides up to 1500 mg/dL, Biotin up to 2.4 μ g/mL, HAMAS up to 1.2 μ g/mL, Rheumatoid Factor (RF) up to 1688 IU/mL, Fulvestrant up to 100 ng/mL, and Mifepristone up to 4.6 μ g/mL did not interfere with the assay. Interferences were observed for bilirubin unconjugated at levels of 20 mg/dL or higher.

PRECISION

The precision study was performed according to the CLSI EP05-A3 guideline



Repeatability

The experimental protocol used a nested components-of-variance design with 8 serum samples, 10 testing days, two lots and two scientists per day. Each scientist ran two tests with two lots per day and two replicate measurements per run (a 10 x 2 x 2 x 2 design) for each sample. The results were analyzed with a two-way nested ANOVA and are summarized in the table below.

Sample	Mean (pg/mL)	Within Run		Between Run		Total	
		SD (pg/mL)	CV %	SD (pg/mL)	CV %	SD (pg/mL)	CV %
1	91.5	8.5	9.2%	11.7	12.8%	14.4	15.8%
2	40.7	5.1	12.4%	6.2	15.1%	8.0	19.6%
3	144.8	11.9	8.2%	15.2	10.5%	20.1	13.9%
4	744.4	33.4	4.5%	31.7	4.3%	46.7	6.3%
5	632.8	26.3	4.2%	41.9	6.6%	56.2	8.9%
6	1027	55.1	5.4%	26.1	2.5%	73.5	7.2%
7	381.0	18.2	4.8%	25.5	6.7%	34.1	8.9%
8	1211.7	53.0	4.4%	71.2	5.9%	106.2	8.8%

Reproducibility

The reproducibility study evaluated the precision performance of the device following EP05-A3 experimental design model 3 x 5 x 5 (3 locations x five testing days x five replicates per day) across laboratories located in Italy, the USA, and Canada. The results were analyzed with a two-way nested ANOVA and are summarized in the table below.

Sample	Mean (pg/mL)	Repeatability		Within Location		Reproducibility	
		SD (pg/mL)	CV %	SD (pg/mL)	CV %	SD (pg/mL)	CV %
Control 1	88.3	6.7	7.5%	9.8	10.5%	10.8	12.2%
Control 2	515.5	25.9	5.0%	33.6	6.5%	46.1	8.9%
1	43.0	6.6	15.4%	7.0	16.3%	7.1	16.6%
2	75.1	6.8	15.4%	7.0	16.3%	7.1	16.6%
3	122.6	9.7	7.9%	8.8	11.8%	9.8	13%
4	129.4	8.7	6.7%	9.7	7.5%	11.9	9.2%
5	447.4	24.5	5.5%	31.3	7.0%	36.8	8.5%
6	912.3	51.0	5.6%	64.8	7.1%	66.9	7.3%



	Sample	Measured (pg/mL)	Expected	Recovery %
Low value: Sample A High Value: Sample B	100% Sample A	45.5	-	-
	100% Sample B	878.8	-	-
	90% Sample A / 10% Sample B	135.7	128.8	105.4
	70% Sample A/ 30% Sample B	285.7	295.3	96.8
	50% Sample A / 50% Sample B	397.3	461.9	86.0
	30% Sample A / 70% Sample B	616.9	628.4	98.2
	10% Sample A / 90% Sample B	838.3	795.0	105.5
Low Value: Sample C High Value: Sample D	100% Sample C	56.8	-	-
	100% Sample D	768.5	-	-
	90% Sample C / 10% Sample D	152.2	127.9	119.0
	70% Sample C/ 30% Sample D	318.4	270.3	117.8
	50% Sample C / 50% Sample D	482.8	412.6	117.0
	30% Sample C/ 70% Sample D	553.7	555.0	99.8
	10% Sample C / 90% Sample D	641.3	697.4	92.0
Low Value: Sample E High Value: Sample F	100% Sample E	45.2	-	-
	100% Sample F	1113.2	-	-
	90% Sample E / 10% Sample F	146.0	152.0	96.1
	70% Sample E/ 30% Sample F	373.4	365.6	102.1
	50% Sample E / 50% Sample F	676.3	579.2	116.8
	30% Sample E / 70% Sample F	928.4	792.8	117.1
	10% Sample E / 90% Sample F	994.7	1006.4	98.8

LINEARITY

The linearity study was performed according to the CLSI EP06-Ed2 guideline using six human serum samples covering the range of the assay. The samples were diluted in low



estrone value (<60 pg/mL) serum samples up to ten percent (1:10), tested in duplicate, and the regression equation of the results (y) compared to the concentration predicted (x) from the dilution factor was $y = 1.001x + 10.2$, $r=0.999$

The relative non-linearity ranged between -10.6% and 10.5% across all samples and measurement dilution points. The statistical analysis shows that the assay is sufficiently linear up to a 1:10 dilution when using low estrone value (<60 pg/mL) serum samples as the diluent.

COMPARITIVE STUDIES

The Estrone ELISA kit (y) was compared against a Liquid Chromatography-Mass Spectrometry (LC-MS/MS) method (x) and yielded the following linear regression results: $y = 0.80x + 25.82$, 105 samples, $r=0.92$, slope = 0.80

REFERENCE RANGES

Reference ranges (95%) were estimated using samples obtained from individuals of diverse races (all values reported in pg/mL). Each laboratory shall establish their own range of reference values.

Cohort	N	Mean	Median	95%	
				2.5%	97.5%
Adult Female Premenopausal*	140	93.9	83.3	19.5	231.9
<i>Adult Female, Menstrual Cycle</i>					
1 – 10 days	40	84.4	81.5	29.8	146.7
11 –20 days	40	87.7	79.6	20.9	232.0
21 – 30 days	40	82.2	73.2	27.2	173.8
Adult Female Postmenopausal*	205	31.9	42.5	ND	166.4
Adult Male	202	59.1	52.1	ND	187.2

*The menopausal status was classified according to age.
ND = Non-Detectable; results below the LoD (14.8 pg/mL)

REFERENCES

1. Fraser CG. Biological Variation: From Principles to Practice. AACCC Press, 2013.
2. Faslodex® Product Monograph. AstraZeneca Canada, 2017.
3. Robertson JFR and Harrison M. Fulvestrant Pharmacokinetics and pharmacology. Br J Cancer. 2004;90 Suppl 1(Suppl 1): S7–S10.
4. Siemens. Urgent Field Safety Notice. Advia Centaur Systems. Cross reactivity of Fulvestrant in Siemens Healthcare Diagnostics Estradiol Assays. 2016.
5. Berkane N, Liere P, Oudinet JP, et al. From Pregnancy to Preeclampsia: A Key Role for Estrogens. Endocr Rev. 2017; 38:123–144.
6. Fleck SC, Twaddle NC, Churchwell MI, Doerge DR, Pande P, Teeguarden JG. Comparative estrogenicity of endogenous, environmental, and dietary estrogens in pregnant women I: Serum levels, variability and the basis for urinary biomonitoring of serum estrogenicity. Food Chem Toxicol. 2018; 115:511–22.
7. Hauptmann H, et al. Concepts for the Synthesis of Biotinylated Steroids. Part II: 17β-Estradiol Derivatives as Immunochemical Probes. Bioconjug Chem. 2000;11(4):537–48.



8. Meyer HH, et al. Immunoaffinity Chromatography and a Biotin-Streptavidin Amplified Immunoassay for Sensitive and Specific Estimation of Estradiol-17 β . *J Steroid Biochem.* 1990; 35(2):263–69.
9. Folan J, et al. Solid-Phase Enzymoimmunoassay of Estrone in Serum. *Clin Chem.* 1988; 34(9):1843–6.
10. Folan J, et al. Solid-Phase Enzymoimmunoassay of Estrone in Saliva. *Clin Chem.* 1989; 35(4):569–72.
11. Speroff L, et al. Hormone Biosynthesis, Metabolism, and Mechanism of Action. In: *Clinical Gynecologic Endocrinology & Infertility*, 3rd ed. Baltimore: Willams & Wilkins; 1983:1–41.
12. Kim MH, et al. Plasma Levels of Estrogen, Androgens and Progesterone During Normal and Dexamethasone-treated Cycles. *J Clin Endocrinol Metab.* 1974; 39(4):706–12.
13. Speight AC, et al. Non-Protein Bound Oestrogens in Plasma and Urinary Excretion of Unconjugated Oestrogens in Non- Pregnant Women. *J Endocrinol.* 1979; 83(3):385–91.
14. Kricka, L.J. Human Anti-Animal Antibody Interferences in Immunological Assays. *Clin Chem.* 1999; 45(7):942–56.
15. Check JH, et al. Falsely Elevated Steroidal Assay Levels Related to Heterophile Antibodies Against Various Animal Species. *Gynecol Obstet Invest.* 1995; 40(2):139–40.
16. Bidlingmaier F, Wagner-Barnack M, Butenandt O, Knorr D: Plasma estrogens in childhood and puberty under physiologic and pathologic conditions. *Pediatr Res.* 1973;7(11):901-907.
17. Nelson RE, et al. Liquid Chromatography-Tandem Mass Spectrometry Assay for Simultaneous Measurement of Estradiol and Estrone in Human Plasma. *Clin Chem.* 2004;50(2):373–84.
18. Masi CM, et al. Serum Estrogen Metabolites and Systolic Blood Pressure Among Middle-Aged and Older Women and Men. *Am J Hypertens.* 2009;22(11): 1148–53.

WARRANTY INFORMATION

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.



For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.