



**EAGLE**  
BIOSCIENCES

# **Estrone LIA Assay Kit**

Catalog Number: ESN31-L01 (1 x 96 wells)

*For Research Use Only.*

*v. 1.1 (07.31.23)*

---

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



## INTENDED USE

The Eagle Biosciences Estrone LIA Assay Kit is for the direct quantitative determination of estrone in human serum by a chemiluminescence immunoassay (LIA). The Eagle Biosciences estrone LIA 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 kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023*

## INTRODUCTION

Estrone is a steroid like estriol and estradiol, belonging to the class of estrogens. The 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 principle of the following chemiluminescence immunoassay (LIA) test follows the typical competitive binding scenario. Competition occurs between an unlabeled antigen (present in standards, control and patient samples) and an enzyme-labelled antigen (conjugate) for a limited number of antibody binding sites on the microplate. The washing and decanting procedures remove unbound materials. After the washing step, the luminescence substrate solution is added. The relative luminescence units (RLUs) are measured on a microtiter plate luminometer. The RLU values are inversely proportional to the concentration of estrone in the sample. A set of calibrators are used to plot a standard curve from which the amount of estrone in patient samples and controls can be directly read.

## PROCEDURAL CAUTIONS AND WARNINGS

1. 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.
2. Control materials or serum pools should be included in every run at a high and low level for assessing the reliability of results.
3. When the use of water is specified for dilution or reconstitution, use deionized or distilled water.
4. In order to reduce exposure to potentially harmful substances, gloves should be worn when handling kit reagents and human specimens.
5. All kit reagents and specimens should be brought to room temperature and mixed gently but thoroughly before use. Avoid repeated freezing and thawing of reagents and specimens.
6. A calibrator curve must be established for every run.
7. The controls should be included in every run and fall within established confidence limits.
8. Improper procedural techniques, imprecise pipetting, incomplete washing as well as improper reagent storage may be indicated when assay values for the controls do not reflect established ranges.
9. The luminescence substrate solutions (A and B) are sensitive to light and should be stored in the original dark bottle away from direct sunlight.



10. When dispensing the substrate and stopping solution, do not use pipettes in which these liquids will come into contact with any metal parts.
11. To prevent contamination of reagents, use a new disposable pipette tip for dispensing each reagent, sample, standard and control.
12. Do not mix various lot numbers of kit components within a test and do not use any component beyond the expiration date printed on the label.
13. Kit reagents must be regarded as hazardous waste and disposed of according to national regulations.

## **LIMITATIONS**

1. All the reagents within the kit are calibrated for the determination of estrone in human serum. The kit is not calibrated for the determination of estrone in saliva, plasma or other specimens of human or animal origin.
2. Do not use grossly hemolyzed, grossly lipemic, icteric or improperly stored serum.
3. Any samples or control sera containing azide or thimerosal are not compatible with this kit, as they may lead to false results.
4. Only calibrator A may be used to dilute any high serum samples. The use of any other reagent may lead to false results.
5. This kit is intended for research use only and should not be used in diagnostic procedures.

## **SAFETY CAUTIONS AND WARNINGS**

### **POTENTIAL BIOHAZARDOUS MATERIAL**

Human serum that may be used in the preparation of the standards and control have been tested and found to be nonreactive for Hepatitis B surface antigen and have also been tested for the presence of antibodies to HCV and Human Immunodeficiency Virus (HIV) and found to be negative. No test method however, can offer complete assurance that HIV, HCV and Hepatitis B virus or any infectious agents are absent. The reagents should be considered a potential biohazard and handled with the same precautions as applied to any human specimen.

### **CHEMICAL HAZARDS**

Avoid contact with reagents containing TMB, hydrogen peroxide and sulfuric acid. If contacted with any of these reagents, wash with plenty of water. TMB is a suspected carcinogen.

### **SPECIMEN COLLECTION AND STORAGE**

Approximately 0.2 mL of serum is required per duplicate determination. Collect 4–5 mL of blood into an appropriately labelled tube and allow it to clot. Centrifuge and carefully remove the serum layer. Store at 4°C for up to 24 hours or at -10°C or lower if the analyses are to be done at a later date. Consider all human specimens as possible biohazardous materials and take appropriate precautions when handling.

### **SPECIMEN PRETREATMENT**

This assay is a direct system; no specimen pretreatment is necessary.

### **REAGENTS AND EQUIPMENT NEEDED BUT NOT PROVIDED**

1. Precision pipettes to dispense 20, 50, 100, 150 and 300  $\mu$ L
2. Disposable pipette tips
3. Distilled or deionized water



4. Plate shaker
5. Microplate luminometer.

## REAGENTS PROVIDED

### 1. Rabbit Anti-Estrone Antibody-Coated Break-Apart Well Microplate — Ready To Use

Contents: One 96-well (12x8) polyclonal antibody-coated microplate in a resealable pouch with desiccant.  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.

### 2. Estrone-Biotin Conjugate Concentrate — Requires Preparation X100

Contents: Estrone-biotin conjugate in a protein-based buffer with a non-mercury preservative.  
Volume: 0.2 mL/vial  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.  
Preparation: See preparation of conjugate working solution.

### 3. Avidin-Horse Radish Peroxidase (HRP) Conjugate Concentrate – Requires Preparation x100

Contents: Avidin-HRP conjugate in a protein-based buffer with a non-mercury preservative.  
Volume: 0.2 mL/vial  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.  
Preparation: See preparation of conjugate working solution.

#### Preparation of Conjugate Working Solution

Dilute both the estrone-biotin and avidin-HRP conjugate concentrates 1:100 into the same solution of assay buffer and mix thoroughly (example: To a tube containing 2 mL of assay buffer add 20 µL of estrone-biotin and 20 µL of avidin-HRP conjugate concentrates). If the whole plate is to be used add 120 µL of estrone-biotin and 120 µL of avidin-HRP conjugate concentrates to 12 mL of assay buffer. Discard any that is left over.

*It is essential that the conjugate working solution be prepared, mixed and allowed to stand for at least 15 minutes prior to use. Failure to do so may result in low optical densities and increased serum values.*

### 4. Estrone Calibrators — Ready To Use

Contents: Six vials containing estrone in a protein-based buffer with a non-mercury preservative. Prepared by spiking buffer with an exact quantity of estrone.

\* Listed below are approximate concentrations, please refer to bottle labels for exact concentrations.



Calibrator	Concentration	Volume
Calibrator A	0 pg/mL	2.0 mL
Calibrator B	15 pg/mL	0.5 mL
Calibrator C	50 pg/mL	0.5 mL
Calibrator D	200 pg/mL	0.5 mL
Calibrator E	800 pg/mL	0.5 mL
Calibrator F	2000 pg/mL	0.5 mL

Storage: Refrigerate at 2–8°C.  
Stability: 12 months in unopened vials or as indicated on label. Once opened, the standards should be used within 14 days or aliquoted and stored frozen. Avoid multiple freezing and thawing cycles.

#### 5. Controls — Ready To Use

Contents: Two vials containing estrone in a protein-based buffer with a non-mercury preservative. Prepared by spiking buffer with an exact quantity of estrone. Refer to vial labels for acceptable range.  
Volume: 0.5 mL/vial  
Storage: Refrigerate at 2–8°C  
Stability: 12 months in unopened vial or as indicated on label. Once opened, the controls should be used within 14 days or aliquoted and stored frozen. Avoid multiple freezing and thawing cycles.

#### 6. Wash Buffer Concentrate – Requires Preparation x10

Contents: One bottle containing buffer with a non-ionic detergent and a non-mercury preservative.  
Volume: 50 mL/bottle  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.  
Preparation: Dilute 1:10 in distilled or deionized water before use. If one whole plate is to be used dilute 50 mL of the wash buffer concentrate in 450 mL of water.

#### 7. Assay Buffer – Ready to Use

Contents: One bottle containing a protein-based buffer with a non-mercury preservative.  
Volume: 15 mL/bottle  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.

#### 8. LIA Substrate Reagent A – Requires Preparation

Contents: One vial containing luminol enhancer.  
Volume: 1.5 mL/vial  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.  
Preparation: See preparation of LIA working substrate solution.

#### 9. LIA Substrate Reagent B – Requires Preparation

Contents: One vial containing peroxide solution.



Volume: 1.5 mL/vial  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.  
Preparation: See preparation of LIA working substrate solution.

#### **10. LIA Substrate Reagent C – Requires Preparation**

Contents: One bottle containing buffer with a non-mercury preservative.  
Volume: 16 mL/bottle  
Storage: Refrigerate at 2–8°C  
Stability: 12 months or as indicated on label.  
Preparation: See preparation of LIA working substrate solution.

#### **PREPARATION OF LIA WORKING SUBSTRATE SOLUTION**

In a clean plastic container (glass is not suitable) mix 1 part of LIA substrate reagent A with 1 part of LIA substrate reagent B and 10 parts of LIA substrate reagent C. This gives the ready to use substrate solution. If the whole plate is to be used prepare working substrate solution as follows: Combine 1.4 mL of LIA substrate reagent A with 1.4 mL of LIA substrate reagent B and 14 mL of LIA substrate reagent C. It is suggested to wait at least 2 minutes prior to use after preparation of the working substrate solution. The working substrate solution is stable for up to 2 hours at room temperature. Discard the leftovers.

#### **ASSAY PROCEDURE**

Important Notes:

- All reagents must reach room temperature before use.
  - Once the procedure has been started, all steps should be completed without interruption to ensure equal elapsed time for each pipetting step.
  - The washing procedure influences the precision markedly; it is essential to ensure the washing is effective and thorough.
1. Prepare working solutions of the conjugate, wash buffer and LIA substrate (refer to reagents provided and preparation section).  
*After conjugate working solution is prepared, it is essential that it be mixed and allowed to stand for at least 15 minutes prior to use.*
  2. Remove the required number of well strips. Reseal the bag and return any unused strips to the refrigerator.
  3. Pipette 50 µL of each calibrator, control and specimen sample into correspondingly labelled wells in duplicate.
  4. Pipette 100 µL of the conjugate working solution into each well. (We recommend using a multichannel pipette.)
  5. Incubate on a plate shaker (approximately 200 rpm) for 1 hour at room temperature.
  6. Wash the wells 5 times with 300 µL of diluted wash buffer per well and tap the plate firmly against absorbent paper to ensure that it is dry. (The use of a washer is recommended).
  7. Pipette 150 µL of LIA working substrate solution into each well. (We recommend using a multichannel pipette.)
  8. Measure the RLU/second in each well on a microplate luminometer between 10–30 minutes after addition of the substrate.



## CALCULATIONS

1. Calculate the mean RLU of each calibrator duplicate.
2. Draw a calibrator curve on semi-log paper with the mean RLUs on the Y-axis and the calibrator concentrations on the X-axis. If immunoassay software is being used, a 4-parameter curve is recommended.
3. Calculate the mean RLU of each unknown duplicate.
4. Read the values of the unknowns directly off the calibrator curve.
5. If a sample reads more than 2000 pg/mL then dilute it with calibrator A at a dilution of no more than 1:8. The result obtained should be multiplied by the dilution factor.

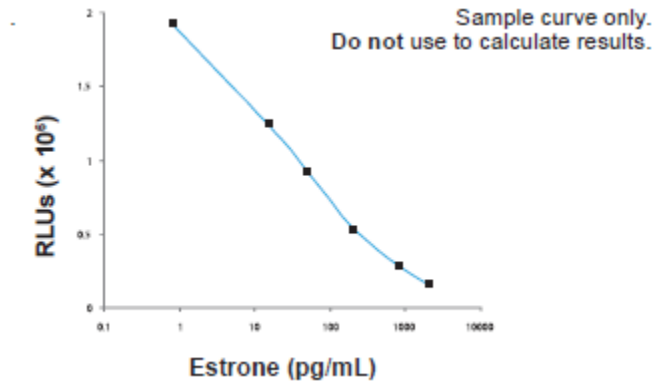
## TYPICAL TABULATED DATA

Sample data only. Do not use to calculate results.

Calibrator	RLU 1 x 10 <sup>3</sup>	RLU 2 x 10 <sup>3</sup>	Mean RLU x 10 <sup>3</sup>	RLU/RLU <sub>MAX</sub> (%)
A, 0 pg/mL	1980	1877	1928	100
B, 15 pg/mL	1256	1226	1241	64
C, 50 pg/mL	832	1004	918	48
D, 200 pg/mL	542	527	534	28
E, 800 pg/mL	311	276	294	15
F, 2000 pg/mL	169	161	165	9

\*\*It is recommended to use the RLU/RLU<sub>MAX</sub> values for comparative purposes since luminometers vary considerably between manufacturers. Results from different luminometers will show quite different RLU values, however, the RLU/RLU<sub>MAX</sub> values remain consistent.

## TYPICAL CALIBRATOR CURVE



## PERFORMANCE CHARACTERISTICS

### SENSITIVITY

The lower detection limit is calculated from the standard curve by determining the resulting concentration of the mean RLU of calibrator A (based on 10 replicate analyses) minus 2 SD. Therefore, the sensitivity of the estrone LIA kit is **8.8 pg/mL**.

### SPECIFICITY (CROSS-REACTIVITY)



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

<b>Steroid</b>	<b>% Cross Reactivity</b>
Estrone	100
Estrone-3-Sulfate	4.9
17 $\beta$ -Estradiol	2.2
Estrone-3-Glucuronide	1.2
17 $\beta$ -Estradiol-3-Glucuronide	0.14

The following steroids were tested but cross-reacted at less than 0.1%: Androstenedione, Cholesterol, Corticosterone, Cortisol, Cortisone, DHEAS, Diethylstilbesterol, Estriol, 17 $\beta$ -Estradiol-3-Glucuronide, Estradiol-Sulfate, Progesterone, 17-OH Progesterone and Testosterone.

### INTRA-ASSAY PRECISION

Three samples were assayed ten times each on the same calibrator curve. The results (in pg/mL) are tabulated below:

<b>Sample</b>	<b>Mean</b>	<b>SD</b>	<b>CV %</b>
1	70.4	4.12	6.6
2	278.7	16.16	5.8
3	787.5	76.9	9.8

### INTER-ASSAY PRECISION

Three samples were assayed ten times over a period of four weeks. The results (in pg/mL) are tabulated below:

<b>Sample</b>	<b>Mean</b>	<b>SD</b>	<b>CV %</b>
1	77.6	9.08	11.7
2	272.4	24.2	8.9
3	823.6	89.77	10.9

### RECOVERY

Spiked samples were prepared by adding defined amounts of estrone to three patient serum samples. The results (in pg/mL) are tabulated below:

<b>Sample</b>	<b>Obs. Result</b>	<b>Exp. Result</b>	<b>Recovery %</b>
1 Unspiked	52	-	-
+ 200	315	252	125
+ 400	557	452	120
+ 1000	1235	1052	117
2 Unspiked	75	-	-
+ 375	493	450	88.0
+ 750	505.23	559.81	90.3
+ 1500	712.44	794.88	89.6
3 Unspiked	720.11	-	-
+ 200	758.13	837.64	90.5
+ 400	856.46	955.17	89.7





+ 1000	1013.61	1190.24	85.1
--------	---------	---------	------

### LINEARITY

Three patient serum samples were diluted with calibrator A. The results (in pg/mL) are tabulated below:

Sample	Obs. Result	Exp. Result	Recovery %
1	340.67	-	-
1:2	165.35	170.34	97.1
1:4	95.39	85.17	112.0
1:8	48.47	42.58	113.8
2	1086.01	-	-
1:2	508.58	543.00	93.7
1:4	232.11	271.50	85.5
1:8	114.95	135.75	84.7
3	1313.21	-	-
1:2	612.98	656.61	93.4
1:4	318.63	328.30	97.1
1:8	134.98	164.15	82.2

### COMPARATIVE STUDIES

The estrone LIA kit (y) was compared with the estrone ELISA kit (x). The comparison of 50 serum samples yielded the following linear regression results:

$$y = 0.8872x - 2.382, \quad r^2 = 0.99$$

### EXPECTED NORMAL VALUES

As for all clinical assays each laboratory should collect data and establish their own range of expected normal values.

Group	Range (pg/mL)
Males	25-150
Females	25-350
Pregnancy	100-8000



## REFERENCES

1. Hauptmann H, et al. Concepts for the Synthesis of Biotinylated Steroids. Part 2:  $17\beta$ -estradiol Derivatives as Immunochemical Probes. *Bioconjug Chem.* 2000; 11(4):537–48.
2. Dressendörfer RA, et al. Synthesis of a Cortisol-biotin Conjugate and evaluation as a tracer in an immunoassay for salivary cortisol measurement. *J Steroid Biochem Mol Biol.* 1992; 43(7):683–92.
3. Meyer HH, et al. Immunoaffinity Chromatography and a Biotin-streptavidin Amplified Immunoassay for Sensitive and Specific Estimation of Estradiol-17. *J Steroid Biochem.* 1990; 35(2):263–9.
4. Folan J, et al. Solid-phase Enzymoimmunoassay of Estrone in Serum. *Clin Chem.* 1988; 34(9):1843–6.
5. Folan J, et al. Solid-phase Enzymoimmunoassay of Estrone in Saliva. *Clin Chem.* 1989; 35(4):569–72.
6. Speroff L, et al. Hormone biosynthesis, metabolism, and mechanism of action. In: *Clinical Gynecologic Endocrinology & Infertility*, 3rd ed. Philadelphia: Williams & Wilkins; 1983:1–41.
7. 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.
8. 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.
9. Kricka LJ. Human Anti-animal Antibody Interferences in Immunological Assays. *Clin Chem.* 1999; 45(7):942–56.
10. 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.

## 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](mailto:info@eaglebio.com) or at 866-411-8023.*