



DCM007-9  
Ed. 04/2013

# FREE ESTRIOL ELISA

for routine analysis

Direct immunoenzymatic determination of Free Estriol in human serum or plasma

IVD



LOT

See external label

2°C  8°C



Σ = 96 tests

REF DKO007

## INTENDED USE

The Eagle Biosciences [Free Estriol ELISA Assay Kit](#) is a competitive immunoenzymatic colorimetric method for quantitative determination of Free Estriol concentration in human serum or plasma.

The Free Estriol ELISA Assay Kit is intended for research use only and not to be used in diagnostic procedures.

## 1. CLINICAL SIGNIFICANCE

Estriol (also oestriol) is one of the three main estrogens produced by the human body. It is only produced in significant amounts during pregnancy as it is made by the fetus.

During pregnancy the production of estriol depends on an intact maternal-placental-fetal unit. Fetal-placental production of estriol leads to a progressive rise in maternal circulating levels reaching a late-gestational peak several orders of magnitude greater than non-pregnant levels. In the maternal circulation, estriol undergoes rapid conjugation in the liver followed by urinary excretion with a half-life of ~20 minutes. Since normal estriol production depends on an intact maternal-placental-fetal circulation and functional fetal metabolism, maternal estriol levels have been used to monitor fetal status during pregnancy, particularly during the third trimester.

DHEA is produced by the adrenal cortex of the fetus, this is converted to estriol by the placenta.

If levels are abnormally low in a pregnant woman, this may indicate a problem with the development in the child. Levels of estriol in non-pregnant women do not change much after menopause, and levels are not significantly different from levels in men.

## 2. PRINCIPLE

In the Free Estriol ELISA Assay Kit, the (antigen) in the sample competes with the antigenic estriol conjugated with horseradish peroxidase (HRP) for binding to the limited number of antibodies anti estriol coated on the microplate (solid phase).

After incubation, the bound/free separation is performed by a simple solid-phase washing.

Then, the enzyme HRP in the bound-fraction reacts with the Substrate (H<sub>2</sub>O<sub>2</sub>) and the TMB Substrate and develops a blue color that changes into yellow when the Stop Solution (H<sub>2</sub>SO<sub>4</sub>) is added. The color intensity is inversely proportional to the free estriol

concentration of in the sample. Free Estriol concentration in the sample is calculated through a calibration curve.

## 3. REAGENTS, MATERIALS AND INSTRUMENTATION

### 3.1. Reagents and materials supplied in the kit

#### 1. Calibrators (7 vials, 1 mL each)

CAL0	REF DCE002/0706-0
CAL1	REF DCE002/0707-0
CAL2	REF DCE002/0708-0
CAL3	REF DCE002/0709-0
CAL4	REF DCE002/0710-0
CAL5	REF DCE002/0711-0
CAL6	REF DCE002/0712-0

#### 2. Control (1 vial, 1 mL)

See concentration of Control on the Certificate of Analysis **REF DCE045/0703-0**

#### 3. Conjugate (1 vial, 15 mL, ready to use)

Estriol conjugated with Horseradish peroxidase (HRP) **REF DCE002/0702-0**

#### 4. Coated Microplate (1 breakable microplate)

Anti-Estriol antibody adsorbed on microplate **REF DCE002/0703-0**

#### 5. TMB Substrate (1 vial, 15 mL)

H<sub>2</sub>O<sub>2</sub>-TMB 0.26 g/L (avoid any skin contact) **REF DCE004-0**

#### 6. Stop Solution (1 vial, 15 mL)

Sulphuric acid 0.15 mol/L (avoid any skin contact) **REF DCE005-0**

#### 7. 10X Conc. Wash Solution (1 vial, 50 mL)

Phosphate buffer 0.2M, Proclin < 0,0015% **REF DCE054-0**

### 3.2. Reagents necessary not supplied

Distilled water.

### 3.3. Auxiliary materials and instrumentation

Automatic dispenser.

Microplates reader (450 nm, 620-630 nm)

## Note

Store all reagents between 2-8°C in the dark.

Open the bag of reagent 4 (Coated Microplate) only when it is at room temperature and close it immediately after use.

#### 4. WARNINGS

- This Free Estriol ELISA Assay Kit is intended for research use by professional persons only. Not for internal or external use in Humans or Animals.
- Use appropriate personal protective equipment while working with the reagents provided.
- Follow Good Laboratory Practice (GLP) for handling blood products.
- All human source material used in the preparation of the reagents has been tested and found negative for antibody to HIV 1&2, HbsAg, and HCV. No test method however can offer complete assurance that HIV, HBV, HCV or other infectious agents are absent. Therefore, the reagents should be handled in the same manner as potentially infectious material.
- Some reagents of the Free Estriol ELISA Assay Kit contain small amounts of Sodium Azide or Proclin 300<sup>R</sup> as preservative. Avoid the contact with skin or mucosa.
- Sodium Azide may be toxic if ingested or absorbed through the skin or eyes; moreover it may react with lead or copper plumbing to form potentially explosive metal azides. If you use a sink to remove the reagents, allow scroll through large amounts of water to prevent azide build-up.
- The TMB Substrate contains an irritant, which may be harmful if inhaled, ingested or absorbed through the skin. To prevent injury, avoid inhalation, ingestion or contact with skin and eyes.
- The Stop Solution consists of a diluted sulphuric acid solution. Sulphuric acid is poisonous and corrosive and can be toxic if ingested. To prevent chemical burns, avoid contact with skin and eyes.
- Avoid the exposure of reagent TMB/H<sub>2</sub>O<sub>2</sub> to directed sunlight, metals or oxidants. Do not freeze the solution.
- This Free Estriol ELISA Assay Kit allows the determination of Free Estriol from 0.15 ng/mL to 40 ng/mL.
- The clinical significance of Free Estriol determination can be invalidated if the sample was treated with natural or syntetic steroids.

#### 5. PRECAUTIONS

- Please adhere strictly to the sequence of pipetting steps provided in this protocol. The performance data represented here were obtained using specific reagents listed in this Instruction For Use.
- All reagents of the Free Estriol ELISA Assay Kit should be stored refrigerated at 2-8°C in their original container. Any exceptions are clearly indicated. The reagents are stable until the expiry date when stored and handled as indicated.
- Allow all Free Estriol ELISA Assay Kit components and specimens to reach room temperature (22-28°C) and mix well prior to use.
- Do not interchange Free Estriol ELISA Assay Kit components from different lots. The expiry date printed on box and vials labels must be observed.. Do not use any kit component beyond their expiry date.

- If you use automated equipment, the user has the responsibility to make sure that the kit has been appropriately tested.
- The incomplete or inaccurate liquid removal from the wells could influence the assay precision and/or increase the background.
- It is important that the time of reaction in each well is held constant for reproducible results. Pipetting of samples should not extend beyond ten minutes to avoid assay drift. If more than 10 minutes are needed, follow the same order of dispensation. If more than one plate is used, it is recommended to repeat the dose response curve in each plate
- Addition of the TMB Substrate solution initiates a kinetic reaction, which is terminated by the addition of the Stop Solution. Therefore, the TMB Substrate and the Stop Solution should be added in the same sequence to eliminate any time deviation during the reaction.
- Observe the guidelines for performing quality control in medical laboratories by assaying controls and/or pooled sera.
- Maximum precision is required for reconstitution and dispensation of the reagents.
- Samples microbiologically contaminated, highly lipemeic or haemolysed should not be used in the assay.
- Plate readers measure vertically. Do not touch the bottom of the wells.

#### 6. PROCEDURE

##### 6.1. Preparation of the Calibrators (C<sub>0</sub>...C<sub>6</sub>)

The Calibrators are ready to use and have the following concentration of Estriol:

	C <sub>0</sub>	C <sub>1</sub>	C <sub>2</sub>	C <sub>3</sub>	C <sub>4</sub>	C <sub>5</sub>	C <sub>6</sub>
ng/mL	0	0.15	0.75	2	5	15	40

The Calibrators are stable until the expiry date printed on the label. Once opened, the Calibrators are stable six months at 2÷8°C.

##### 6.2. Preparation of the Sample

The determination of Estriol can be performed in human plasma as well as in serum.

Store samples at -20°C if the determination is not performed on the same day of the sample connection. Avoid repetitive freezing and thawing of samples.

The Control is ready for use.

##### 6.3. Preparation of Wash Solution

Dilute the contents of each vial of "10X Conc. Wash Solution" with distilled water to a final volume of 500 mL prior to use. For smaller volumes respect the 1:10 dilution ratio. The diluted wash solution is stable for 30 days at 2-8°C. In concentrated wash solution is possible to observe the presence of crystals; in this case mix at room temperature until the complete dissolution of crystals; for greater accuracy, dilute the whole bottle of concentrated wash solution to 500 mL, taking care to transfer completely the crystals, then mix until crystals are completely dissolved.

#### 6.4. Procedure

- **Allow all reagents to reach room temperature (22-28°C) for at least 30 minutes.** At the end of the assay, store immediately the reagents at 2-8°C: avoid long exposure to room temperature.
- Unused coated microwell strips should be released securely in the foil pouch containing desiccant and stored at 2-8°C.
- To avoid potential microbial and/or chemical contamination, unused reagents should never be transferred into the original vials.
- As it is necessary to perform the determination in duplicate in order to improve accuracy of the test results, prepare two wells for each point of the calibration curve (C<sub>0</sub>-C<sub>6</sub>), two for each Control, two for each sample, one for Blank.

Reagent	Calibrator	Sample/Control	Blank
Calibrator C <sub>0</sub> -C <sub>6</sub>	25 µL		
Sample/Control		25 µL	
Conjugate	100 µL	100 µL	
Incubate at 37°C for 1 hour. Remove the contents from each well; wash the wells 6 times with 300 µL of diluted wash Solution. <b>Important note:</b> during each washing step, gently shake the plate for 5 seconds and remove excess solution by tapping the inverted plate on an absorbent paper towel.			
TMB Substrate	100 µL	100 µL	100 µL
Incubate at room temperature (22±28°C) for 15 minutes in the dark.			
Stop Solution	100 µL	100 µL	100 µL
Shake the microplate gently. Read the absorbance (E) at 450 nm against a reference wavelength of 620-630 nm or against Blank within 5 minutes.			

#### 7. QUALITY CONTROL

Each laboratory should assay controls at normal, high and low levels range of Estriol for monitoring assay performance. These controls should be treated as unknowns and values determined in every test procedure performed. Quality control charts should be maintained to follow the performance of the supplied reagents. Pertinent statistical methods should be employed to ascertain trends. The individual laboratory should set acceptable assay performance limits. Other parameters that should be monitored include the 80, 50 and 20% intercepts of the calibration curve for run-to-run reproducibility. In addition, maximum absorbance should be consistent with past experience. Significant deviation from established performance can indicate unnoticed change in experimental conditions or degradation of

kit reagents. Fresh reagents should be used to determine the reason for the variations.

#### 8. RESULTS

##### 8.1. Mean Absorbance

Calculate the mean of the absorbance (Em) for each point of the calibration curve (C<sub>0</sub>-C<sub>6</sub>) and of each sample.

##### 8.2. Calibration curve

Plot the values of absorbance (Em) of the Calibrators (C<sub>0</sub>-C<sub>6</sub>) against concentration. Draw the best-fit curve through the plotted points (es: Four Parameter Logistic).

##### 8.3. Calculation of Results

Interpolate the values of the samples on the calibration curve to obtain the corresponding values of the concentrations expressed in ng/mL.

#### 9. EXPECTED VALUES

47 samples from not pregnancy women were assayed with Diametra Free Estriol ELISA kit; the median of the values is:

N°	Median (Estriol)
47	0.08 ng/mL

Please pay attention to the fact that the determination of a range of expected values for a "normal" population in a given method is dependent on many factors, such as specificity and sensitivity of the method used and type of population under investigation. Therefore each laboratory should consider the range given by the Manufacturer as a general indication and produce their own range of expected values based on the indigenous population where the laboratory works.

#### 10. PERFORMANCE AND CHARACTERISTICS

##### 10.1. Precision

###### 10.1.1. Intra Assay Variation

Within run variation was determined by replicate the measurements (20x) of three different sera in one assay. The within assay variability is ≤ 7.4%.

###### 10.1.2. Inter Assay Variation

Between run variation was determined by replicate the measurements (10x) of three different sera in different lots of kit. The between assay variability is ≤ 7.7%.

##### 10.2. Accuracy

The recovery of 0.51 - 1.02 - 2.03 - 4.07 - 8.13 ng/mL of Estriol added to a sample gave an average value (±SD) of 96.96% ± 6.61% with reference to the original concentrations.

##### 10.3. Sensitivity

The lowest detectable concentration of free estriol that can be distinguished from the Calibrator 0 is 0.03 ng/mL at the 95% confidence limit.

#### 10.4. Specificity

The cross reaction of the antibody calculated at 50% according to Abraham are shown in the table:

Free Estriol	100 %
16 epi-estriol	10.5 %
15 $\alpha$ OH-estriol	7.0 %
Estriol 3 Sulphate	2.0 %
Estradiol	0.1 %
17 epi-estriol	$< 1 \times 10^{-2}$ %
Estriol 3 $\alpha$ -Glucuronate	$< 1 \times 10^{-2}$ %
Estriol 16 $\alpha$ -Glucuronate	$< 1 \times 10^{-2}$ %
Estrone	$< 1 \times 10^{-4}$ %

#### 10.5. Correlation

Diametra Free Estriol ELISA was compared to another commercially available Free Estriol assay. 80 serum samples were analysed.

The linear regression curve was calculated:

$$y = 0.90 x + 0.15$$

$$r^2 = 0.819$$

#### 11. WASTE MANAGEMENT

Reagents must be disposed off in accordance with local regulations.

---

#### BIBLIOGRAPHY

1. Fischer - Rasmussen, W., et al Acta Obstet Gynecol. Scand. 60-417 - 420 (1981)
2. Truran, P.L., et al Clin. Chem. 28/12, 2393 (1982)
3. Vining, R. F., et al J. Clin. Endoc. Metab. 56, 454 (1983)
4. Bagger, P. V., et al Acta Obstet Gynecol Scand 60, 187 (1981)
5. Osterman, T. M., et al Clin. Chem. 25 (5) 716 (1979)
6. Wisdom, G.B. Clin. Chem. 22 (8) 1243-1255 (1976)

Ed. 04/2013

DCM007-9

**DiaMetra S.r.l. Headquater:** Via Garibaldi, 18 –  
20090 SEGRATE (MI) Italy

Tel. 0039-02-2139184 – 02-26921595

Fax 0039-02-2133354.

**Manufactory:** Via Pozzuolo 14, 06034 SPELLO (PG)  
Italy

Tel. 0039-0742-24851

Fax 0039-0742-316197

E-mail: [info@diametra.com](mailto:info@diametra.com)

---

distributed in the US/Canada by:

**EAGLE BIOSCIENCES, INC.**

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

[www.EagleBio.com](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



## **ERROR POSSIBLE CAUSES / SUGGESTIONS**

### **No colorimetric reaction**

- no conjugate pipetted reaction after addition
- contamination of conjugates and/or of substrate
- errors in performing the assay procedure (e.g. accidental pipetting of reagents in a wrong sequence or from the wrong vial, etc.)

### **Too low reaction (too low ODs)**

- incorrect conjugate (e.g. not from original kit)
- incubation time too short, incubation temperature too low

### **Too high reaction (too high ODs)**

- incorrect conjugate (e.g. not from original kit)
- incubation time too long, incubation temperature too high
- water quality for wash buffer insufficient (low grade of deionization)
- insufficient washing (conjugates not properly removed)

### **Unexplainable outliers**

- contamination of pipettes, tips or containers
- insufficient washing (conjugates not properly removed) too high within-run
- reagents and/or strips not pre-warmed to CV% Room Temperature prior to use
- plate washer is not washing correctly (suggestion: clean washer head)
- too high between-run - incubation conditions not constant (time, CV % temperature)
- controls and samples not dispensed at the same time (with the same intervals) (check pipetting order)
- person-related variation