



Estradiol Saliva ELISA Kit

Catalog Number: EST32-K01

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

V. 9 (12 APR 24)

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 Estradiol Saliva ELISA kit is designed and validated for the quantitative measurement of 17 β -Estradiol in human saliva samples. This assay is intended for research use only and should not be used for diagnostic procedures.

Warning: Estradiol immunoassays have been reported to demonstrate significant cross-reactivity with the drug Fulvestrant (FaslodexR). This cross-reactivity can cause falsely elevated estradiol levels in those being treated with Fulvestrant. Due to the possibility of this cross-reactivity, this Estradiol Immunoassay should not be used for those being treated with the drug Fulvestrant. Fulvestrant (FaslodexR) is used to treat a certain type of breast cancer in postmenopausal women.

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.

ASSAY BACKGROUND

17 β -Estradiol (E2) is a steroid hormone produced mainly by the Graafian follicle of the ovary in females and in small amounts by the testes in male subjects. E2 is biologically the most active and naturally produced human estrogens. The majority of E2 (98%) is bound to the sex hormone binding globulin (SHBG) and to the lesser extent to the other serum proteins such as albumin. Only a small fraction circulates in the free form is conjugated to sulfates and glucuronides. In non-pregnant women there is a cyclic variation in the concentration of E2, the highest values being measured usually the day before ovulation. Positive feedback influence of this peak value is considered essential for occurrence of the midcycle luteinizing hormone (hLH) peak and consequently, for ovulation. During pregnancy the E2 concentration increases considerably and remains high throughout pregnancy.

This assay of E2 is a valuable tool for assessing and researching the etiology of amenorrhea and/or infertility in female samples. It is also a useful aid in monitoring ovulation induction treatment with clomiphene citrate, LH-RH (LH-releasing hormone) or exogenous gonadotropins (8,9). In male samples, serum E2 measurements are used for researching feminizing syndromes.

ASSAY PRINCIPLE

The Eagle Biosciences, Inc. Estradiol Saliva ELISA Kit, is based on the competition principle and microplate separation. An unknown amount of estradiol present in a saliva sample and a fixed amount of estradiol conjugated to horse radish peroxidase (E2-HRP) compete for binding sites with a rabbit monoclonal estradiol antiserum bound to GARGG (goat anti-rabbit gamma globulin) coated wells of a microplate. After incubation, unbound components are washed away. Enzyme substrate solution is then added, and a blue color formed. This reaction is stopped with an acid solution to produce a yellow color. The optical density is then read at 450 nm. The amount of E2-HRP detected is inversely proportional to the amount of estradiol in a sample.

MATERIALS PROVIDED

1. **GARGG Plate:** One 96 well microplate (12x8 breakable strip wells) coated with goat anti-rabbit gamma globulin (GARGG) placed in a resealable foil bag with desiccant. One (1) 96 well kit is sufficient for 78 singlicate or 39 duplicate patient measurements.



2. **Stock Estradiol (17- β E2) in BSA buffer. Concentration:** 3200 pg/mL, 1 bottle, 0.250 mL

Preparation: From stock **3200 pg/mL**, dilute **1:100** with **assay buffer** to obtain the **32 pg/mL** calibrator. Make serial dilutions (in assay buffer) starting with the **32 pg/mL** calibrator to obtain the following calibrator concentrations: **16 pg/mL, 8 pg/mL, 4 pg/mL, 2 pg/mL, 1 pg/mL**. "**0**" calibrator is assay buffer.

3. **Assay buffer.** 1 bottle, 20 mL.
4. Salivary **Estradiol EIA Control #1.** 1 bottle, 1.0 mL. Concentration is on the label and is traceable to Estradiol (17 β -Estradiol) from USP.
5. Salivary **Estradiol EIA Control #2:** 1 bottle, 1.0 mL. Concentration is on the label and is traceable to Estradiol (17 β -Estradiol) from USP.
6. Salivary **Estradiol EIA Antibody.** 1 bottle, 3 mL. The solution is blue.
7. Salivary **Estradiol Horse Radish Peroxidase Conjugate concentrate (E2-HRP concentrate)** 1 amber bottle, 0.200 mL. The solution is yellow and light sensitive.
8. **E2-HRP Conjugate Buffer.** 4 mL. To be **used for E2-HRP working reagent preparation ONLY.**

E2-HRP working reagent preparation: Determine the amount of working E2HRP working reagent needed and dilute 1:20 with conjugate buffer. For example, mix 125 μ L (0.125 mL) of E2-HRP concentrate (number 7) to a total volume of 2.5 mL (0.125 mL + 2.375 mL) with conjugate buffer (number 8). This is sufficient for 100 EIA wells. Immediately after use, store the unused portion of the Estradiol-HRP working reagent at 2-8°C. Discard if not used within 3 weeks of mixing.

9. **Wash solution (10X concentrated) EIA #1:** 1 bottle, 50 mL. Prior to use dilute 1:10 with deionized water.
10. **Color Development Reagent EIA #1:** 1 amber plastic bottle, 15 mL of Tetramethyl-Benzidine (TMB) plus hydrogen peroxide. Light sensitive.
11. **Stopping Solution EIA #1:** 1 bottle 15 mL diluted acid solution.

STORAGE AND STABILITY

- When stored at 2° – 8° C, unopened reagents will retain activity until the expiration date. Do not use the reagents beyond this date.
- Use only reagents supplied with this kit. Do not interchange reagents with different lot numbers.
- Opened reagents must be stored at 2° – 8° C.
- Microtiter wells must be stored at 2° – 8° C. Once the foil bag has been opened, care should be taken to reseal tightly.
- Opened kits retain activity for 3 weeks if stored as described above.
- Expiration dates and lot numbers are printed on the labels.



MATERIALS NEEDED BUT NOT PROVIDED

- Device to dispense very accurately 100 μ L of saliva.
- Multichannel pipettors.
- Microplate or orbital shake
- Vortex mixer
- Microplate washer (not required, plates can be washed manually).
- Microplate reader capable of reading 450 nm with 4 parameter data reduction or comparable software.
- Plate sealers.
- Collection device

INTERFERENCES

Commonly consumed products such as alcohol, caffeine, nicotine and food and gum extracts may interfere with Estradiol measurements. As a precautionary measure, follow the sample collection instructions as stated in section VIII. 1, under Collection.

SAMPLE COLLECTION AND PROCESSING

Collection: This sample collection and processing procedure must be followed.

- Avoid food consumption, drinking coffee or alcohol, smoking or chewing gum one (1) hour prior to sample collection.
- Rinse mouth thoroughly with water 15 minutes prior to collection.
- Collect whole saliva by unstimulated passive drool by allowing saliva to drip off the lower lip into a graduated plastic test or by allowing saliva to accumulate in the floor of the mouth and spitting it into the recommended collection device.
- Time and date specimen. Refrigerate then freeze (-20°C or below) samples until day of assay. On day of assay, Thaw samples to facilitate precipitation of mucins. Centrifuge at 1500g for ten minutes. Bring samples to room temperature and assay.

SAMPLE STABILITY

Storage	Room Temperature 20 - 30°C	37°C	2 - 8°C	$\leq -15^{\circ}\text{C}$ (7 Freeze / thaw cycles)	$\leq -15^{\circ}\text{C}$ (Long term)
Stability	Up to 7 days	Up to 7 days	Up to 7 days	Up to 7 days	Up to 9 months



PROCEDURE SUMMARY FLOW SHEET

Calibrator Control Sample (I.D.)	Calibrator control (µL)	Cortisol-HRP Working Reagent (µL)	Anti-Cortisol (µL)		Diluted 10X Wash Solution (µL)		Color Developer (µL)		Stopping Solution (µL)	
0	100	25	25	Mix. Incubate 30 min. at room temperature. with shaking	300	Wash 3x	125	Mix. Incubate 30 min. At room temperature	125	Mix. Read at 450nm
1	100	25	25		300		125			
2	100	25	25		300		125			
4	100	25	25		300	125				
8	100	25	25		300	125				
16	100	25	25		300	125				
32	100	25	25		300	125				
Control #1	100	25	25		300	125				
Control #2	100	25	25		300	125				
Sample #1	100	25	25		300	125				
Sample #2	100	25	25		300	125				
Sample #3	100	25	25		300	125				

ASSAY PROCEDURE

1. It is recommended that the **calibrators, controls** and **samples** should be tested in duplicate and the mean value should be used to report results.
2. To the GARGG microplate **dispense 100 µL** of the previously prepared working **Estradiol Calibrators** (0, 1, 2, 4, 8, 16 and 32 pg/mL), **controls**, and **saliva** samples. Mix the microplate by shaking gently (manual) for a few seconds.
3. Add **25 µL** of **Estradiol HRP Working Reagent**.
4. Add **25 µL** of **Anti-Estradiol EIA Antibody**.
5. Cover the microplate with plastic sealer. **Incubate by shaking** on a microplate orbital shaker set at **500 – 900 rpm** for **2 hours** at room temperature.
6. After incubation decant content of the wells. Wash 3 times with 300 µL of diluted Wash Solution (**10 mL** of 10X Wash solution EIA #1 diluted with **90 mL** of D.I. water) after the 3rd wash invert GARGG microplate on absorbent paper and tap dry.
7. Dispense **125 µL** of Color Development reagent EIA #1 into each well. Shake briefly (manual). Cover microplate with plastic sealer. Incubate for 30 minutes at room temperature.



8. Dispense **125µL** of Stopping Solution EIA #1 into each microtiter well of the GARGG plate. Shake briefly (manual). Color changes from blue to yellow.
9. **Read at 450 nm** on a microplate reader **within 10 minutes**.

TYPICAL RESULTS

Calibrators (pg/mL)	Mean Absorbance	%B/Bo	Value (pg/mL)
0	2.16	100.0	0
1	1.82	84.3	1
2	1.65	76.4	2
4	1.33	61.6	4
8	0.86	39.8	8
16	0.50	23.1	16
32	0.32	14.8	32
Control #1	1.53	70.8	2.5
Control #2	0.73	33.8	10.7
Sample #1	1.65	76.4	1.9
Sample #2	0.33	15.3	28.2
Sample #3	0.97	44.9	6.9

CALULATION OF RESULTS

1. Compute the average optical density (OD) for the zero (Bo) calibrator.
2. Calculate the percent bound (B/Bo) for each calibrator, control and unknown by dividing the average OD (B) by the average OD for the zero (Bo) x 100.
3. Plot percent bound (B/Bo) versus the calibrator concentrations and draw the best fit for the curve.
4. Plot percent bound (B/Bo) of the controls and unknowns to determine saliva estradiol concentrations.
5. Alternately determine the concentrations of the controls and unknowns by interpolation using software capable of logistics using a 4-parameter sigmoid minus curve fit.

Analytical Measuring range (AMR)	1 pg/mL – 32 pg/mL
---	--------------------

Samples with estradiol values greater than 32 pg/mL should be diluted 1:10 with zero (0) calibrator and re-run for accuracy. Obtain the final estradiol concentration by multiplying the diluted sample by the dilution factor.

QUALITY CONTROL

The expected values for the controls are stated on the label of each control which are included in the kit. The results can only be accepted if the expected values are met.

EXPECTED VALUES

The following AM values were obtained from apparently healthy subjects with the Salivary Estradiol Elisa Kit

Estradoil Saliva ELISA Kit
Catalog Number: EST32-K01



Estradiol (pg/mL)		Median	Range
Females Premenopausal (not using contraceptives) N=90, Age 21 – 50 yrs	Follicular Phase	3.16	2.13 – 5.6
	Luteal Phase	3.15	1.42 – 14.43
Females Postmenopausal n=51, – 75 Yrs		2.53	0.80 – 4.07
Males n=43, Age 20-70 Yrs		2.68	1.07 – 5.00

It is recommended that each laboratory establish its own normal ranges

PERFORMANCE CHARACTERISTICS

C-18 Steroids	Spiked Concentration	% Cross Reactivity
Estradiol-17 β	10,000 pg/mL	100.00
Estradiol-17a	10,000 pg/mL	0.0112
Estriol	10,000 pg/mL	0.0615
Estrone	10,000 pg/mL	0.9225
Estrone-3-SO4	10,000 pg/mL	0.3398
Estriol-3-SO4	10,000 pg/mL	0.0066
C-19 Steroids		
Androstenedione	10,000 pg/mL	0.0008
Testosterone	10,000 pg/mL	0.0082
5 α -Dihydrotestosterone	10,000 pg/mL	0.0050
Dehydroepiandrosterone SO4	10,000 pg/mL	0.0030
Androstenedione	10,000 pg/mL	0.0148
C-21 Steroids		
Progesterone	10,000 pg/mL	<0.001
17-OH Progesterone	10,000 pg/mL	0.0048
Pregnenolone	10,000 pg/mL	0.0044
17-OH-Pregnenolone	10,000 pg/mL	0.0096
Deoxycorticosterone	10,000 pg/mL	0.0063
11-Desoxycortisol	10,000 pg/mL	0.0085
Corticosterone	10,000 pg/mL	0.0171
Aldosterone	10,000 pg/mL	0.0072
Cortisol	10,000 pg/mL	0.0104

DETECTION LIMITS

The LOB (limit of the blank) and the LOD (limit of detection) were determined by generating one hundred twenty (120) measurements each of “estradiol free saliva” and low level (<1 pg/mL) estradiol samples (Reference, CLSI EP 17-A, protocols for Determination of Limits of Detection and Limits of Quantitation).

Limit of the Blank (LoB) pg/mL	Limit of Detection (LoD) pg/mL
0.3680	0.5239



INTRA-ASSAY PRECISION

The intra-assay precision was determined from the mean of 20 replicates of low, medium and high samples.

Sample	N	Mean	SD	CV %
Low	20	3.29	0.352	10.7
Medium	20	27.24	0.999	3.7
High	20	52.56	3.103	5.9

INTER-ASSAY PRECISION

The inter-assay precision was determined from the mean of the average duplicates of 12 separate assays with low, medium and high samples

Sample	N	Mean	SD	CV %
Low	12	3.01	0.266	8.8
Medium	12	26.13	0.939	3.6
High	12	49.19	4.552	9.3

REPEATABILITY

This study was conducted during 5 days of a familiarization period and 20 days of testing. Two assays were performed daily. Three (3) different reagent lots and three (3) saliva pools were used for the study (low, medium and high). The pools were aliquoted and frozen until day of assay.

PRECISION LOW CONCENTRATION POOL

	SD	CV %
Within Run	0.047	1.7
Between Run	0.052	1.9
Repeatability	0.105	3.9
Total Device Precision	0.126	4.6

PRECISION MEDIUM CONCENTRATION POOL

	SD	CV %
Within Run	0.323	1.2
Between Run	0.096	0.4
Repeatability	0.552	2.1
Total Device Precision	0.647	2.5

PRECISION HIGH CONCENTRATION POOL

	SD	CV %
Within Run	0.321	0.7
Between Run	0.964	2.1
Repeatability	0.859	1.8
Total Device Precision	1.331	2.8



LINEARITY

S=10 samples (dilutions)

Concentration=(C1*V1+C10*V10)/(V1+V10)

	C1 (pg/mL)	V1 (mL)	C10 (pg/mL)	V10 (mL)	Calculated Concentration (pg/mL)	Obtained Concentration (pg/mL)	Recovery (%)
1				*	0.40	0.42	105
2	0.42	0.889	81.57	0.111	9.43	8.44	90
3	0.42	0.778	81.57	0.222	18.44	16.61	90
4	0.42	0.667	81.57	0.333	27.44	24.88	91
5	0.42	0.556	81.57	0.444	36.45	36.32	100
6	0.42	0.444	81.57	0.556	45.54	45.50	100
7	0.42	0.333	81.57	0.667	54.55	56.08	103
8	0.42	0.222	81.57	0.778	63.55	65.69	103
9	0.42	0.111	81.57	0.889	72.56	76.66	106
10				*	80.00	81.57	102

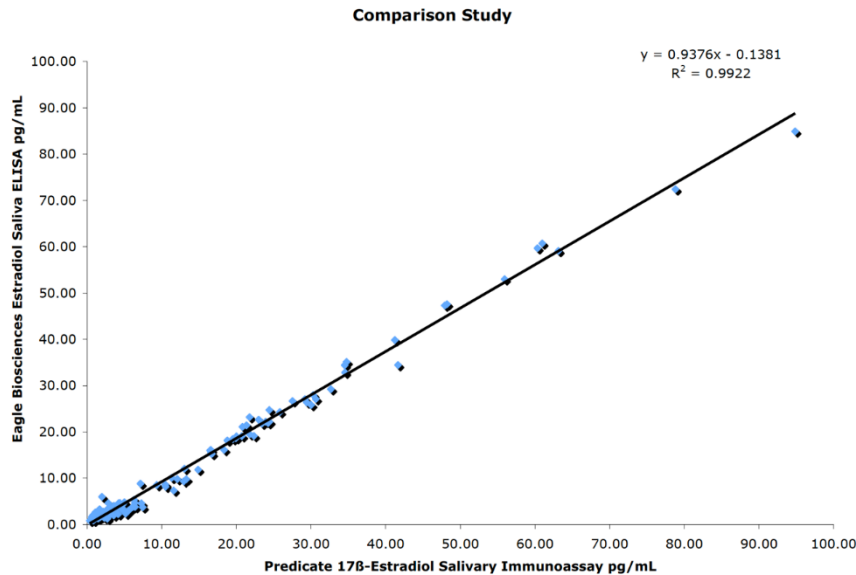
RECOVERY

Seven (7) saliva samples containing different levels of endogenous estradiol were spiked with known quantities of estradiol, assayed and the percent recovery determined.

Sample	Endogenous (pg/mL)	Estradiol Added (pg/mL)	Expected (pg/mL)	Observed (pg/mL)	Recovery (%)
1	1.00	64.0	65.00	60.37	92.9
2	3.20	32.0	35.20	34.53	98.1
3	4.42	16.0	20.42	19.55	95.7
4	4.24	16.0	12.24	11.53	94.2
5	24.61	4.0	28.61	26.82	93.7
6	40.90	2.0	42.90	42.02	97.9
7	4.86	1.0	5.86	5.30	90.4

COMPARISON STUDY

One hundred twenty-five (125) saliva samples spanning the analytical measuring range (AMR) and beyond were analyzed by comparing the Eagle Biosciences Direct Estradiol Saliva ELISA Kit, Cat #EST31-K01 with a commercially available predicate Estradiol salivary Immunoassay. See correlation results below.



REFERENCES

1. Baird, D.T. (1976): Ovarian steroid secretion and metabolism in women. In the Endocrine Function of the Human Ovary, Eds. V.H.T. James, M., Serio, G. Giusti. Academic Press, London, New York, San Francisco, pp. 125-133.
2. Tsang, B.K., et al. (1980) Steroid biosyntheses by isolated human ovarian follicular cells in vitro, *J.Clin. Endocrinol. Metab.* 51, 1407-11.
3. Gore-Langton, R.E. et.al (1988): Follicular steroidogenesis and its control. In: The physiology of Reproduction. Ed.: Knobil et al. p.331-385, Raven press, New York.
4. Abraham, G.E., Odell, W.D., Swerdloff, R.S. and Hopper, K (1972): Simultaneous radioimmunoassay of plasma FSH, LH, progesterone, 17-hydroxyprogesterone, and estradiol-17 during the menstrual cycle. *J. Clin. Endocrinol. Metab.* 34, 312-318.
5. Punnonen, R., Nummi, S., Ylikorkala, O., Alapiessa, U., Karvonen, P. and Viinikka, L. (1974): A composite picture of the normal menstrual cycle. *Acta Obstet. Gynecol. Scand. Suppl.* 51, 65-70.
6. Jaffe, R. (1974): Regulation of the human menstrual cycle. In *Physiology and Genetics of Reproduction, Part A*. Eds. E.M. Countinho and F. Fuchs. Plenum Press, New York and London, pp. 371-383.
7. Klopper, A., Jandial, V. and Wilson, G. (1975): Plasma steroid assay in the assessment of foetoplacental function. *J. Steroid Biochem.* 6, 651-656.
8. Kerin, J.F., Warnes, G.M., Quinn, P., Kirby, C., Godefrey, B. and Cox, L.W.(1984): Endocrinology of ovarian stimulation for in vitro fertilization. *Aust. NZ. J. Obstet. Gynaec.* 24, 121-124.
9. Ronnberg, L., Martikainen, H. and Kirkinen, P. (1988): The effect of the mode of clomiphene citrate administration on ovarian response in an in vitro fertilization program. *Int. J. Fertil.* 33, 334-337.
10. Odell, W.D., and Swerdloff, R.S. (1978): Abnormalities of gonadal function in men. *Clin. Endocrinol.* 8, 149-180.



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 insert, please contact the Technical Service Team at Eagle Biosciences, Inc at www.EagleBio.com or at 866-411-8023.