



EAGLE
BIOSCIENCES

Pregnenolone ELISA Assay Kit

Catalog Number:
PGN31-K01 (1 x 96 wells)
For Research Use Only
v. 2.0 (11.17.22)

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

INTRODUCTION

Pregnenolone (3β -hydroxypregn-5-en-20-one) is the first steroid to be derived from cholesterol in the pathway of steroidogenesis, and it is the common precursor for all of the adrenal and gonadal steroids. Its production occurs in the mitochondrion by cleavage of the C-20 side chain of cholesterol by the P-450SCC enzyme. Once produced, pregnenolone may be utilized by two pathways of steroidogenesis. Pregnenolone may either be converted to 17-OH pregnenolone via the enzymatic action of 17α -hydroxylase or to progesterone via the enzymatic action of 3β hydroxysteroid dehydrogenase. Elevated pregnenolone levels occur in forms of congenital adrenal hyperplasia (CAH), due to 3β -hydroxysteroid dehydrogenase deficiency or 17α hydroxylase deficiencies. Higher levels have also been reported in women with idiopathic hirsutism. Studies on pregnenolone levels in regard to sex and age differences indicate that maximum levels occur at approximately 17 and 16 years of age for women and men, while minimum levels occur at approximately 37 and 38 years of age for women and men, respectively. In general, women were found to have slightly higher values when compared to men. Many areas of pregnenolone physiology remain to be investigated. Current research indicates that the determination of pregnenolone in serum may be useful for studying its metabolite, pregnenolone sulfate, which has been reported to have various effects in the mammalian brain and central nervous system.

PRINCIPLE OF THE ASSAY

The principle of the following enzyme immunoassay test follows the typical competitive binding scenario. Competition occurs between an unlabelled antigen (present in standards, controls 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 enzyme substrate is added. The enzymatic reaction is terminated by addition of the stopping solution. The absorbance is measured on a microtiter plate reader. The intensity of the colour formed is inversely proportional to the concentration of pregnenolone in the sample. A set of standards is used to plot a standard curve from which the concentration of pregnenolone 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. It is recommended to all customers to prepare their own control materials or serum pools that 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 provided in the kit 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. When reading the microplate, the presence of bubbles in the wells will affect the optical densities (ODs). Carefully remove any bubbles before performing the reading step.
10. The substrate solution (TMB) is sensitive to light and should remain colourless if properly stored. Instability or contamination may be indicated by the development of a blue colour, in which case it should not be used.
11. When dispensing the substrate and stopping solution, do not use pipettes in which these liquids will come into contact with any metal parts.
12. To prevent contamination of reagents, use a new disposable pipette tip for dispensing each reagent, sample, standard and control.
13. 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.
14. 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 direct determination of pregnenolone in human serum. The kit is not calibrated for the determination of pregnenolone 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 for research use only and should not be used as a diagnostic tool.

SAFETY CAUTIONS AND WARNINGS

POTENTIAL BIOHAZARDOUS MATERIAL

Human serum that may be used in the preparation of the standards and controls has been tested and found to be nonreactive for Hepatitis B surface antigen and has 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 blood 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 50, 100, 150 and 300 μ L
2. Disposable pipette tips
3. Distilled or deionized water
4. Plate shaker
5. Microplate reader with a filter set at 450 nm and an upper OD limit of 3.0 or greater* (see assay procedure step 10)

REAGENTS PROVIDED 1. Rabbit Anti-Pregnenolone 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. Pregnenolone-Horseradish Peroxidase (HRP) Conjugate Concentrate — Requires Preparation X51

Contents: Pregnenolone-HRP conjugate in a protein-based buffer with a non-mercury preservative. Volume: 300 μ L/vial Storage: Refrigerate at 2–8°C
Stability: 12 months or as indicated on label.
Preparation: Dilute 1:50 in assay buffer before use (eg. 40 μ L of HRP in 2 mL of assay buffer). If the whole plate is to be used dilute 240 μ L of HRP in 12 mL of assay buffer. Discard any that is left over.

3. Pregnenolone Calibrators — Ready To Use

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

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

| Calibrator | Concentration | Volume/Vial |
|-------------------|----------------------|--------------------|
| Calibrator A | 0 ng/mL | 2.0 mL |
| Calibrator B | 0.1 ng/mL | 0.5 mL |
| Calibrator C | 0.4 ng/mL | 0.5 mL |
| Calibrator D | 1.6 ng/mL | 0.5 mL |
| Calibrator E | 6.4 ng/mL | 0.5 mL |
| Calibrator F | 25.6 ng/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.

4. **Controls** — Ready To Use

Contents: Two vials containing pregnenolone in a protein-based buffer with a non-mercury preservative. Prepared by spiking buffer with defined quantities of pregnenolone. Refer to vial labels for the 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.

5. **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 the whole plate is to be used dilute 50 mL of the wash buffer concentrate in 450 mL of water.

6. **Assay Buffer** – Ready to Use

Contents: One bottle containing 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.

7. **TMB Substrate** – Ready to Use

Contents: One bottle containing tetramethylbenzidine and hydrogen peroxide in a non-DMF or DMSO containing buffer.

Volume: 16 mL/vial

Storage: Refrigerate at 2–8°C

Stability: 12 months or as indicated on label.

8. **Stopping Solution** – Ready to Use

Contents: One bottle containing 1M sulfuric acid. Volume: 6 mL/bottle

Storage: Refrigerate at 2–8°C

Stability: 12 months or as indicated on label.



ASSAY PROCEDURE

Specimen Pretreatment: None

All reagents must reach room temperature before 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. Prepare working solutions of the pregnenolone-HRP conjugate and wash buffer.
2. Remove the required number of well strips. Reseal the bag and return any unused strips to the refrigerator.
3. Pipette 50 μ L of each calibrator, control and specimen sample into correspondingly labelled wells in duplicate.
4. Pipette 100 μ L 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 3 times with 300 μ L 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 μ L of TMB substrate into each well.
8. Incubate on a plate shaker for 10–15 minutes at room temperature (or until calibrator A attains dark blue colour for desired OD).
9. Pipette 50 μ L of stopping solution into each well at the same timed intervals as in step 7.
10. Read the plate on a microplate reader at 450 nm within 20 minutes after addition of the stopping solution.

* If the OD exceeds the upper limit of detection or if a 450 nm filter is unavailable, a 405 or 415 nm filter may be substituted. The optical densities will be lower, however, this will not affect the results of patient/control samples.

CALCULATIONS

1. Calculate the mean optical density of each calibrator duplicate.
2. Draw a calibrator curve on semi-log paper with the mean optical densities on the Y-axis and the calibrator concentrations on the X-axis. If immunoassay software is being used, a 4-parameter or 5-parameter curve is recommended.
3. Calculate the mean optical density of each unknown duplicate.
4. Read the values of the unknowns directly off the calibrator curve.
5. If a sample reads more than 25.6 ng/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

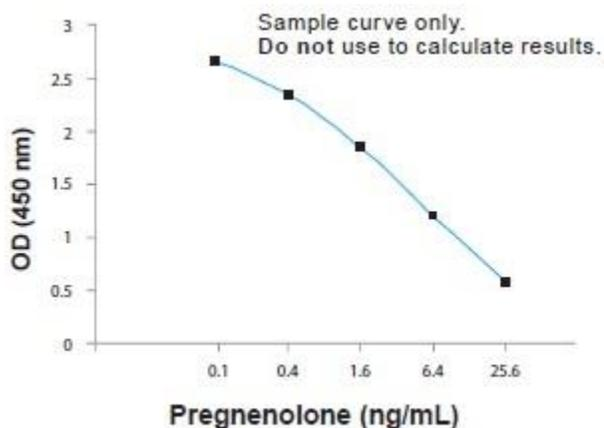
Sample data only. Do not use to calculate results.

| Calibrator | OD 1 | OD 2 | Mean OD | Value (ng/mL) |
|------------|-------|-------|---------|---------------|
| A | 2.891 | 2.808 | 2.850 | 0 |
| B | 2.613 | 2.651 | 2.632 | 0.1 |



| | | | | |
|---------|-------|-------|-------|------|
| C | 2.350 | 2.343 | 2.347 | 0.4 |
| D | 1.823 | 1.879 | 1.851 | 1.6 |
| E | 1.237 | 1.197 | 1.217 | 6.4 |
| F | 0.589 | 0.594 | 0.591 | 25.6 |
| Unknown | 1.431 | 1.451 | 1.441 | 4.0 |

TYPICAL CALIBRATOR CURVE



PERFORMANCE CHARACTERISTICS SENSITIVITY

The lower detection limit is calculated from the standard curve by determining the resulting concentration of the mean OD of Calibrator A (based on 10 replicate analyses) minus 2 SD. Therefore, the sensitivity of the Direct Pregnenolone ELISA kit is 0.05 ng/mL.

SPECIFICITY (CROSS-REACTIVITY)

The following compounds were tested for cross-reactivity with the Direct Pregnenolone ELISA kit with pregnenolone cross-reacting at 100%.

| Steroid | % Cross Reactivity |
|---------------------------|--------------------|
| Pregnenolone | 100 |
| Progesterone | 6.0 |
| Dehydroisoandrosterone | 5.2 |
| 5 α -Androstandiol | 4.7 |
| Epiandrosterone | 1.0 |
| Pregnenolone Sulfate | 0.4 |
| Androstandione | 0.3 |
| 5 α -Androsterone | 0.3 |
| DHEAS | 0.2 |
| Etiocholanolone | 0.1 |

The following steroids were tested but cross-reacted at less than 0.1%: Adrenosterone, Aldosterone, Androstenedione, Cholesterol, Corticosterone, 5 α -DHT, 17 β -Estradiol, Estriol and Testosterone.



INTRA-ASSAY PRECISION

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

| Sample | Mean | SD | CV % |
|--------|------|------|------|
| 1 | 0.19 | 0.02 | 10.6 |
| 2 | 1.04 | 0.85 | 8.2 |
| 3 | 4.77 | 0.37 | 7.8 |

INTER-ASSAY PRECISION

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

| Sample | Mean | SD | CV % |
|--------|------|------|------|
| 1 | 0.22 | 0.03 | 14.5 |
| 2 | 1.14 | 0.14 | 12.3 |
| 3 | 4.56 | 0.44 | 9.6 |

RECOVERY

| Sample | Obs. Result | Exp. Result | Recovery % |
|------------|-------------|-------------|------------|
| 1 Unspiked | 0.37 | - | - |
| +4.14 | 5.31 | 4.51 | 117.7 |
| 2 Unspiked | 0.77 | - | - |
| +4.01 | 5.69 | 4.78 | 119.0 |
| 3 Unspiked | 0.85 | - | - |
| +3.98 | 5.18 | 4.83 | 107.2 |
| 4 Unspiked | 1.47 | - | - |
| +3.78 | 6.31 | 5.25 | 120.2 |

LINEARITY

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

| Sample | Obs. Result | Exp. Result | Recovery % |
|---------|-------------|-------------|------------|
| 1 | 5.31 | - 2.66 | - |
| 1:2 | 2.89 | 1.33 | 108.6 |
| 1:4 1:8 | 1.26 | 0.66 | 94.7 |
| | 0.71 | | 107.6 |
| 2 | 6.51 | - 3.26 | - |
| 1:2 | 2.75 | 1.63 | 84.4 |
| 1:4 1:8 | 1.54 | 0.81 | 94.5 |
| | 0.80 | | 98.8 |



| | | | |
|---------|-----------|--------|-------------|
| 3 | 8.34 3.78 | - 4.17 | - 90.6 |
| 1:2 | 2.15 | 2.09 | 102.9 101.0 |
| 1:4 1:8 | 1.05 | 1.04 | |

EXPECTED NORMAL VALUES

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

| Group | N | Mean (ng/mL) | Abs. Range (ng/mL) |
|---------|----|--------------|--------------------|
| Males | 30 | 0.50 | 0.1-3.4 |
| Females | 50 | 0.55 | 0.1-3.8 |

REFERENCES

1. Abraham GE, et al. Radioimmunoassay of Plasma Pregnenolone, 17-hydroxypregnenolone and Dehydroepiandrosterone Under Various Physiological Conditions. *J Clin Endocrinol Metab.* 1973; 37(1):140–4.
2. Hill M, et al. Age Relationships and Sex Differences in Serum Levels of Pregnenolone and 17-hydroxypregnenolone in Healthy Subjects; *Clin Chem Lab Med.* 1999; 37(4):439–47.
3. Robel P, et al. Biosynthesis and Assay of Neurosteroids in Rats and Mice. *J Steroid Biochem Mol Biol.* 1995; 53:355–60.
4. Weusten JJ, et al. Early Time Sequence in Pregnenolone Metabolism to Testosterone in Homogenates of Human and Rat Testis. *Endocrinology.* 1987; 120(5):1909–13.
5. Akwa Y, et al. Neurosteroids: Biosynthesis, Metabolism and Function of Pregnenolone and Dehydroepiandrosterone in the brain. *J Steroid Biochem Mol Biol.* 1991; 40(1-3):7181.
6. McKenna TJ, Brown RD. Pregnenolone in Man: Plasma Levels in States of Normal and Abnormal Steroidogenesis. *J Clin Endocrinol Metab.* 1974; 38(3):480–5.
7. McKenna TJ, et al. Pregnenolone, 17-OH-Pregnenolone, and Testosterone in Plasma of Patients with Congenital Adrenal Hyperplasia. *J Clin Endocrinol Metab.* 1976; 42(5):918–25.
8. McKenna TJ, et al. Plasma Pregnenolone in Patients with Adrenal Tumors, ACTH Excess, or Idiopathic Hirsutism. *J Clin Endocrinol Metab.* 1977; 44(2):231–6.
9. Wilson J D, Foster, DW, eds. *Williams Textbook of Endocrinology* 8th Edition. London: W.B. Saunders Company; 1992:495-7.
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 or at 866-411-8023.