



Pregnenolone ELISA Kit

Catalog Number KA1912

96 assays

Version: 04

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	5
Precautions for Use	5
Assay Protocol	7
Reagent Preparation	7
Sample Preparation	7
Assay Procedure	7
Data Analysis.....	9
Calculation of Results.....	9
Performance Characteristics	10
Resources	12
References	12
Plate Layout	13

Introduction

Intended Use

For the direct quantitative determination of Pregnenolone by enzyme immunoassay in human serum.

Background

Pregnenolone (3 β -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-450_{SCC} 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 17 α -hydroxylase or to progesterone via the enzymatic action of 3 β -hydroxysteroid dehydrogenase.

Elevated pregnenolone levels occur in forms of congenital adrenal hyperplasia (CAH), due to 3 β -hydroxysteroid dehydrogenase deficiency or 17 α -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 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 microwell plate. 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 amount of pregnenolone in human samples and controls can be directly read.

General Information

Materials Supplied

List of component

Component	State	Amount
Rabbit Anti-Pregnenolone Antibody Coated Microwell Plate-Break Apart Wells: Polyclonal antibody-coated microwell plate in a resealable pouch with desiccant.	Ready To Use	96 wells
Pregnenolone-Horseradish Peroxidase (HRP) Conjugate Concentrate (50x): Pregnenolone-HRP conjugate in a protein-based with a non-mercury preservative.	Concentrate	0.3 mL
Standards: Standard A: 0 ng/mL; Standard B: 0.1 ng/mL; Standard C: 0.4 ng/mL; Standard D: 1.6 ng/mL; Standard E: 6.4 ng/mL; Standard F: 25.6 ng/mL. Pregnenolone in a protein-based buffer with a non-mercury preservative. Prepared by spiking buffer with a defined quantity of pregnenolone.	Ready To Use	2.0 mL 0.5 mL 0.5 mL 0.5 mL 0.5 mL 0.5 mL
Controls (Control 1 and Control 2): Refer to vial labels for expected value and acceptable range! Prepared by spiking buffer with a defined quantity of pregnenolone.	Ready To Use	0.5 mL/vial
Wash Buffer Concentrate (10x): One bottle containing buffer with a non-ionic detergent and a non-mercury preservative.	Concentrate	50 mL
Assay Buffer: One vial containing a protein-based buffer with a non-mercury preservative.	Ready To Use	15 mL
TMB Substrate: One bottle containing tetramethylbenzidine and hydrogen peroxide in a non-DMF or DMSO containing buffer.	Ready To Use	16 mL
Stopping Solution: One vial containing 1 M sulfuric acid.	Ready To Use	6 mL

Storage Instruction

Store the complete kit at 2-8°C. Under these conditions, the kit is stable for 12 months or as indicated on label. Standards and Controls: once opened, the standards should be used within 14 days or aliquoted and stored frozen. Avoid multiple freezing and thawing cycles.

Materials Required but Not Supplied

- ✓ Precision pipettes to dispense 50, 100, 150 and 300 µL
- ✓ Disposable pipette tips
- ✓ Distilled or deionized water
- ✓ Plate shaker
- ✓ Microwell plate reader with a filter set at 450 nm and an upper OD limit of 3.0 or greater (see Assay Procedure step 10).

Precautions for Use

- ✓ Procedural Cautions and Warnings
- 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.
- Control materials or serum pools should be included in every run at a high and low level for assessing the reliability of results.
- When the use of water is specified for dilution or reconstitution, use deionized or distilled water.
- In order to reduce exposure to potentially harmful substances, gloves should be worn when handling kit reagents and human specimens.
- 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.
- A standard curve must be established for every run.
- The control should be included in every run and fall within established confidence limits.
- Improper procedural techniques, imprecise pipetting, incomplete washing as well as improper reagent storage may be indicated when assay values for the control do not reflect established ranges.
- When reading the microplate, the presence of bubbles in the microwells will affect the optical densities (ODs). Carefully remove any bubbles before performing the reading step.
- 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.
- When dispensing the substrate and stopping solution, do not use pipettes in which these liquids will come into contact with any metal parts.
- To prevent contamination of reagents, use a new disposable pipette tip for dispensing each reagent, sample, standard and control.
- 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.
- Kit reagents must be regarded as hazardous waste and disposed of according to national regulations.

✓ Limitations

- 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.
- Do not use grossly hemolyzed, grossly lipemic, icteric or improperly stored serum.
- Any samples or control sera containing azide or thimerosal are not compatible with this kit, as they may lead to false results.
- Only standard A may be used to dilute any high serum samples. The use of any other reagent may lead to false results.
- The results obtained with this kit should never be used as the sole basis for a clinical diagnosis. For example, the occurrence of heterophilic antibodies in patients regularly exposed to animals or animal products has the potential of causing interferences in immunological tests. Consequently, the clinical diagnosis should include all aspects of a patient's background including the frequency of exposure to animals/products if false results are suspected.

✓ Safety Cautions and Warnings

• Potential biohazards material

Human serum that may be used in the preparation of the standards and control has been tested and found to be non-reactive 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. However no test method 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.

Assay Protocol

Reagent Preparation

- ✓ Pregnenolone-Horseradish Peroxidase (HRP) Conjugate Concentrate - X50
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.
- ✓ Wash Buffer Concentrate (10x)
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.

Sample Preparation

- ✓ Specimen Collection and Storage
Approximately 0.2 mL of serum is required per duplicate determination. Collect 4-5 mL of blood into an appropriately labeled 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.

Assay Procedure

All reagents must reach room temperature before use. Standards, 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 microwell strips. Reseal the bag and return any unused strips to the refrigerator.
3. Pipette 50 μ L of each standard, control and specimen sample into the 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 prepared wash buffer (300 μ L/well for each wash) 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 at timed intervals.
8. Incubate the plate on a plate shaker for 10-15 minutes at room temperature (or until Standard 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 microwell plate 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 tested/control samples.

Data Analysis

Calculation of Results

- ✓ Calculate the mean optical density of each standard duplicate.
- ✓ Draw a standard curve on semi-log paper with the mean optical densities on the Y-axis and the standard concentrations on the X-axis. If immunoassay software is being used, a 4-parameter curve is recommended.
- ✓ Calculate the mean optical density of each unknown duplicate.
- ✓ Read the values of the unknowns directly off the standard curve.
- ✓ If a sample reads more than 25.6 ng/mL then dilute it with standard A at a dilution of no more than 1:8. The result obtained should be multiplied by the dilution factor.

- ✓ Typical tabulated data:

Standard	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 (Sample curve only. Do not use to calculate results):

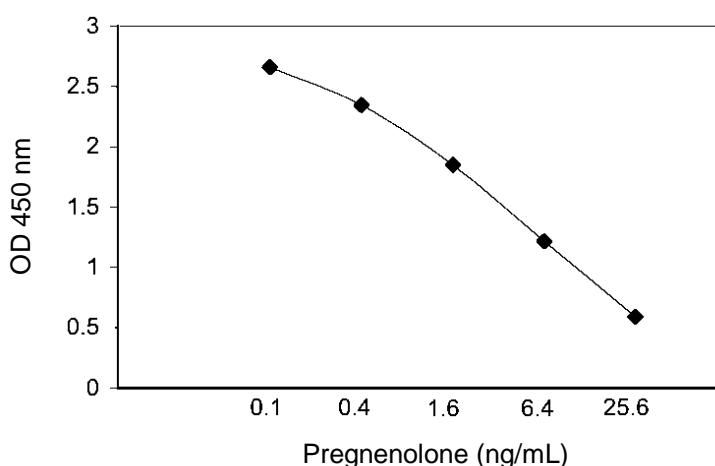


Figure 1: Typical Standard Curve for Pregnenolone ELISA Kit.

Performance Characteristics

✓ Sensitivity

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

✓ Specificity (cross reactivity)

The following compounds were tested for cross-reactivity with the 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

Spiked samples were prepared by adding defined amounts of pregnenolone to four human serum samples. The results (in ng/mL) are tabulated below:

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 human serum samples were diluted with standard A. The results (in ng/mL) are tabulated below:

Sample	Obs. Result	Exp. Result	Recovery%
1	5.31	-	-
1:2	2.89	2.66	108.6
1:4	1.26	1.33	94.7
1:8	0.71	0.66	107.6
2	6.51	-	-
1:2	2.75	3.26	84.4
1:4	1.54	1.63	94.5
1:8	0.80	0.81	98.8
3	8.34	-	-
1:2	3.78	4.17	90.6
1:4	2.15	2.09	102.9
1:8	1.05	1.04	101.0

✓ Expected normal values

As for all 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

Resources

References

1. Abraham, G.E., et al. Radioimmunoassay of Plasma Pregnenolone. J. Clin. Endocrinol. Metab. 37(1):40, 1973.
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 37(4):439, 1999 .
3. Robel, P., et al., Biosynthesis and Assay of Neurosteroids in Rats and Mice. J. Steroid Biochem. Molec. Biol. 53:335, 1995.
4. Weusten, J.J., et al., Early Time Sequence in Pregnenolone Metabolism to Testosterone in Homogenates of Human and Rat Testis. Endocrinology 120:1909, 1987.
5. Akwa, Y, et al., Neurosteroids: Biosynthesis, Metabolism and Function of Pregnenolone and Dehydroepiandrosterone in the brain. J. Steroid Biochem. Molec. Biol. 40(1-3):71, 1991.
6. McKenna, T.J. and Brown, R.D., Pregnenolone in Man: Plasma Levels in States of Normal and Abnormal Steroidogenesis. J. Clin. Endocrinol. Metab. 38(3):480, 1974.
7. McKenna, T.J., et al., Pregnenolone, 17-OH-Pregnenolone, and Testosterone in Plasma of Patients with Congenital Adrenal Hyperplasia. J. Clin. Endocrinol. Metab. 42(5):918, 1976.
8. McKenna, T.J., et al., Plasma Pregnenolone in Patients with Adrenal Tumors, ACTH Excess, or Idiopathic Hirsutism. J. Clin. Endocrinol. Metab. 44(2):231, 1977.
9. Wilson, J.D. and Foster, D.W. Williams Textbook of Endocrinology 8 th Edition. W.B. Saunders Company, London. pp 495-497, 1992.
10. Check, J.H., et al, Falsely elevated steroidal assay levels related to heterophile antibodies against various animal species. Gynecol Obstet Invest 40:139-140, 1995.

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H