



Aldosterone ELISA Kit

Catalog Number KA1883

96 assays

Version: 09

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	9
Data Analysis.....	10
Calculation of Results.....	10
Performance Characteristics	12
Resources.....	15
References	15
Plate Layout	16

Introduction

Intended Use

The Aldosterone ELISA Kit is an enzyme immunoassay for the measurement of aldosterone in serum, plasma (EDTA-, heparin- or citrate plasma) and urine.

Background

The steroid hormone aldosterone is a potent mineral corticoid that is produced by the zona glomerulosa of the adrenal cortex in the adrenal gland. The synthesis and release are controlled by the renin-angiotensin-aldosterone system (RAAS)¹, as well as by plasma potassium concentration², the pituitary peptide ACTH, and by the blood pressure via pressure sensitive baroreceptors in the vessel walls of nearly all large arteries of the body³. Aldosterone binds to mineralocorticoid receptors (MR) and triggers the transcription of hormone-responsive genes.

Principle of the Assay


The Aldosterone ELISA Kit is a solid phase enzyme-linked immunosorbent assay (ELISA), based on the principle of competitive binding.

The microtiter wells are coated with a polyclonal rabbit antibody directed towards an antigenic site of the aldosterone molecule. Endogenous aldosterone of a sample competes with an aldosterone-horseradish peroxidase conjugate for binding to the coated antibody. After incubation the unbound conjugate is washed off. After addition of the substrate solution, the intensity of colour developed is inversely proportional to the concentration of aldosterone in the sample.

General Information

Materials Supplied

List of component

Component	Detail	Amount
Microtiterwells	12 x 8 (break apart) strips. Wells coated with anti-aldosterone antibody (polyclonal rabbit).	96 wells
Enzyme Conjugate	Ready to use. Aldosterone conjugated to horseradish peroxidase, contain non-mercury preservative.	14 mL
Substrate Solution	Ready to use. Tetramethylbenzidine (TMB).	14 mL
Stop Solution	Ready to use. Contains 0.5 M H ₂ SO ₄ . Avoid contact with the stop solution. It may cause skin irritations and burns.  Hazards identification: H290 May be corrosive to metals. H314 Causes severe skin burns and eye damage.	14 mL
Wash Solution	40x concentrated	30 mL

Standards and Controls - lyophilized

Component	Concentration (pg/mL)	Amount
Standard A	0	1.0 mL
Standard B	20	1.0 mL
Standard C	80	1.0 mL
Standard D	200	1.0 mL
Standard E	500	1.0 mL
Standard F	1000	1.0 mL
Control 1	For control values and ranges please refer to vial label or QC-Report.	1.0 mL
Control 2		1.0 mL

Contents: contain non-mercury preservative.

Conversion: 1 pg/mL corresponds to 2.77 pmol/L

Note: Additional Standard A for sample dilution is available upon request.

Storage Instruction

When stored at 2 °C to 8 °C unopened reagents will retain reactivity until expiration date. Do not use reagents beyond this date.

Opened reagents must be stored at 2 °C to 8 °C. Microtiter wells must be stored at 2 °C to 8 °C. Once the foil bag has been opened, care should be taken to close it tightly again.

Opened kits retain activity for two months if stored as described above.

Materials Required but Not Supplied

- ✓ A microtiter plate calibrated reader (450 ± 10 nm)
- ✓ Calibrated variable precision micropipettes.
- ✓ Absorbent paper.
- ✓ Distilled or deionized water
- ✓ Timer
- ✓ Scale paper or semi-logarithmic graph paper or software for data reduction

Precautions for Use

1. This kit is for research use only. For professional use only.
2. All reagents of this test kit which contain human serum or plasma have been tested and confirmed negative for HIV I/II, HBsAg and HCV by FDA approved procedures. All reagents, however, should be treated as potential biohazards in use and for disposal.
3. Before starting the assay, read the instructions completely and carefully. Use the valid version of protocol provided with the kit. Be sure that everything is understood.
4. The microplate contains snap-off strips. Unused wells must be stored at 2-8°C in the sealed foil pouch and used in the frame provided.
5. Pipetting of samples and reagents must be done as quickly as possible and in the same sequence for each step.
6. Use reservoirs only for single reagents. This especially applies to the substrate reservoirs. Using a reservoir for dispensing a substrate solution that had previously been used for the conjugate solution may turn solution colored. Do not pour reagents back into vials as reagent contamination may occur.
7. Mix the contents of the microplate wells thoroughly to ensure good test results. Do not reuse microwells.
8. Do not let wells dry during assay; add reagents immediately after completing the rinsing steps.
9. Allow the reagents to reach room temperature (21-26°C) before starting the test. Temperature will affect the absorbance readings of the assay. However, values for the samples will not be affected.
10. Never pipet by mouth and avoid contact of reagents and specimens with skin and mucous membranes.
11. Do not smoke, eat, drink, or apply cosmetics in areas where specimens or kit reagents are handled.
12. Wear disposable latex gloves when handling specimens and reagents. Microbial contamination of reagents or specimens may give false results.
13. Handling should be done in accordance with the procedures defined by an appropriate national biohazard safety guideline or regulation.
14. Do not use reagents beyond expiry date as shown on the kit labels.
15. All indicated volumes have to be performed according to the protocol. Optimal test results are only obtained when using calibrated pipettes and microtiterplate readers.

16. Do not mix or use components from kits with different lot numbers. It is advised not to exchange wells of different plates even of the same lot. The kits may have been shipped or stored under different conditions and the binding characteristic of the plates may result slightly different.
17. Avoid contact with Stop Solution containing 0.5 M H₂SO₄. It may cause skin irritation and burns.
18. Some reagents contain Proclin 300, BND and MIT as preservatives. In case of contact with eyes or skin, flush immediately with water.
19. TMB substrate has an irritant effect on skin and mucosa. In case of possible contact, wash eyes with an abundant volume of water and skin with soap and abundant water. Wash contaminated objects before reusing them. If inhaled, take the person to open air.
20. Chemicals and prepared or used reagents have to be treated as hazardous waste according to the national biohazard safety guideline or regulation.
21. For information on hazardous substances included in the kit please refer to Safety Data Sheet. Safety Data Sheet for this product is available upon request directly from the manufacturer.

✓ Disposal of the Kit

The disposal of the kit must be made according to the national regulations. Special Information for this product is given in the Safety Data Sheet.

✓ Damaged Test Kits

In case of any severe damage to the test kit or components, the manufacturer has to be informed in writing, at the latest, one week after receiving the kit. Severely damaged single components should not be used for a test run. They have to be stored until a final solution has been found. After this, they should be disposed according to the official regulations.

✓ Limitations of use

Reliable and reproducible results will be obtained when the assay procedure is performed with a complete understanding of the package insert instruction and with adherence to good laboratory practice. Any improper handling of samples or modification of this test might influence the results.

✓ Interfering Substances

The following substances have no influence on the assay results up to the below stated concentrations.

	Serum	Urine
	Concentration up to	
Haemoglobin	4 mg/mL	
Bilirubin	0.5 mg/mL	
Triglyceride	7.5 mg/mL	

✓ Drug interferences

Until today no substances (drugs) are known to us, which have an influence to the measurement of aldosterone in a sample.

✓ High-Dose-Hook effect

A High-Dose-Hook Effect is not detected up to 20,000 pg/mL of aldosterone.

Assay Protocol

Reagent Preparation

Bring all reagents and required number of strips to room temperature prior to use.

✓ Standards

Reconstitute the lyophilized contents of the standard vials with 1.0 mL deionized water and let stand for at least 10 minutes. Mix several times before use.

Note: The reconstituted standards are stable for 8 weeks at 2°C to 8°C.

For longer storage freeze - only once - at -20°C.

✓ Controls

Reconstitute the lyophilized content of the controls with 1.0 mL deionized water and let stand for at least 10 minutes. Mix several times before use.

Note: The reconstituted controls are stable for 8 weeks at 2°C to 8°C.

For longer storage freeze - only once - at -20°C.

✓ Wash Solution

Add deionized water to the 40X concentrated Wash Solution.

Dilute 30 mL of concentrated Wash Solution with 1170 mL deionized water to a final volume of 1200 mL.

The diluted Wash Solution is stable for 2 weeks at room temperature.

Sample Preparation

Serum or plasma (EDTA-, heparin- or citrate plasma) and urine can be used in this assay.

Do not use haemolytic, icteric or lipaemic specimens.

Please note: Samples containing sodium azide should not be used in the assay.

✓ Serum / Plasma Samples

- Specimen Collection

Serum:

Collect blood by venipuncture (e.g. Sarstedt Monovette for serum), allow to clot, and separate serum by centrifugation at room temperature. Do not centrifuge before complete clotting has occurred.

Plasma:

Whole blood should be collected into centrifuge tubes containing anti-coagulant (e.g. Sarstedt Monovette with the appropriate plasma preparation) and centrifuged immediately after collection.

- Specimen Storage and Preparation

Specimens should be capped and may be stored for up to 5 days at 2 °C to 8 °C prior to assaying.

Specimens held for a longer time (up to two months) should be frozen only once at -20 °C prior to assay. Thawed samples should be inverted several times prior to testing.

- Specimen Dilution

If in an initial assay, a specimen is found to contain more than the highest standard, the specimens can be diluted with Standard A and reassayed as described in Assay Procedure.

For the calculation of the concentrations this dilution factor has to be taken into account.

Example:

a) dilution 1:10: 10 μ L sample + 90 μ L Standard A (mix thoroughly)

b) dilution 1:100: 10 μ L dilution a) 1:10 + 90 μ L Standard A (mix thoroughly).

- ✓ Urine Samples

Aldosterone concentration can also be determined from urine samples. However, urine samples must be pretreated before analysis. This will need additional reagents that are not included in this kit.

- Sample Collection

First clean genital area with mild disinfectant to prevent contamination. Then collect clean-catch midstream urine in an appropriate sterile container. Directly after collection, the urine should be centrifuged for 5 - 10 minutes (e.g. at 2,000 g) to remove cellular debris. Use supernatant for analyte quantification. The supernatant may be stored for up to 8 hours at 2 °C to 8 °C prior to assaying. Specimens held for a longer time should be frozen at -20 °C. Thawed supernatant should be inverted several times prior to testing.

- Protocol for Urine Sample Pre-treatment

1. Secure the desired number of vials (e.g. 0.5 - 1.5 mL plastic tubes; not included in this kit).
2. Dispense 25 μ L of urine with new disposable tips into appropriate tubes.
3. Dispense 25 μ L Release Reagent into each tube. Thoroughly mix for 10 seconds. It is important to have a complete mixing in this step.
4. Incubate over night at 2 °C to 8 °C.
5. Add 25 μ L Neutralization Reagent to each tube and mix thoroughly.
6. Add 400 μ L Dilution Buffer to each tube and mix thoroughly.
(This pre-treatment leads to a 1:19 dilution. Therefore the dilution factor 19 has to be taken into account for calculation of the final concentration of the urine sample.)
7. Transfer 100 μ L of pre-treated and diluted urine samples directly to the microtiter well and continue with step 3 of Assay Procedure.

- Storage of pre-treated urine samples

Pre-treated and diluted urine samples should be capped and may be stored for up to 7 days at 2 °C to 8 °C prior to assaying.

Samples held for a longer time (up to two months) should be frozen only once at -20 °C prior assay. Thawed samples should be inverted several times prior to testing.

- Specimen Dilution

If in an initial assay, an urine sample is found to contain more than the highest standard, the pre-treated and diluted urine sample can be further diluted with Dilution Buffer and reassayed as described in Assay Procedure.

For the calculation of the concentrations this dilution factor has to be taken into account too.

Example:

a) dilution 1:10: 10 μ L pre-treated and diluted urine sample + 90 μ L Dilution Buffer (mix thoroughly)
(final dilution factor = $19 \times 10 = 190$)

Assay Procedure

- All reagents and specimens must be allowed to come to room temperature before use. All reagents must be mixed without foaming.
 - Once the test has been started, all steps should be completed without interruption.
 - Use new disposal plastic pipette tips for each standard, control or sample in order to avoid cross contamination.
 - Absorbance is a function of the incubation time and temperature. Before starting the assay, it is recommended that all reagents are ready, caps removed, all needed wells secured in holder, etc. This will ensure equal elapsed time for each pipetting step without interruption.
 - As a general rule the enzymatic reaction is linearly proportional to time and temperature.
 - Each run must include a standard curve.
-
1. Secure the desired number of Microtiter wells in the frame holder.
 2. Dispense 100 μ L of each Standard, Control and samples with new disposable tips into appropriate wells. For urine samples dispense 100 μ L of the pre-treated and diluted urine samples (see section Protocol for Urine Sample Pre-treatment, step 7).
 3. Incubate for 30 minutes at room temperature.
 4. Dispense 100 μ L Enzyme Conjugate into each well. Thoroughly mix for 10 seconds. It is important to have a complete mixing in this step.
 5. Incubate for 60 minutes at room temperature.
 6. Wash the wells as follows: Manually:
 - i. Briskly shake out the contents of the wells.
 - ii. Rinse the wells 3 times with 300 μ L diluted wash solution per well.
 7. If an automated plate washer is used:
 - i. Rinse the wells 3 times with 400 μ L diluted wash solution per well.
 8. At the end of the wash step, always strike the wells sharply on absorbent paper to remove residual droplets.
 9. Add 100 μ L of Substrate Solution to each well.
 10. Incubate for 30 minutes at room temperature.
 11. Stop the enzymatic reaction by adding 50 μ L of Stop Solution to each well.
 12. Determine the absorbance (OD) of each well at 450 ± 10 nm with a microtiter plate reader. It is recommended that the wells be read within 10 minutes after adding the Stop Solution.

Data Analysis

Calculation of Results

1. Calculate the average absorbance values for each set of standards, controls and samples.
2. Using scale paper or semi-logarithmic graph paper, construct a standard curve by plotting the mean absorbance obtained from each standard against its concentration with absorbance value on the vertical (Y) axis and concentration on the horizontal (X) axis.
3. Using the mean absorbance value for each sample determine the corresponding concentration from the standard curve.
4. Automated method: The results in the Instructions for Use have been calculated automatically using a 4 Parameter curve fit. (4 Parameter Rodbard or 4 Parameter Marquardt are the preferred methods.) Other data reduction functions may give slightly different results.
5. The concentration of the serum/plasma samples can be read directly from this standard curve.
For urine samples the concentration read from the standard curve, has to be multiplied with the dilution factor 19 (see section Protocol for Urine Sample Pre-treatment).
6. Samples with concentrations higher than that of the highest standard have to be further diluted or reported as > 1000 pg/mL. For the calculation of the concentrations this dilution factor has to be taken into account too.

✓ **Example of Typical Standard Curve**

The following data is for demonstration only and cannot be used in place of data generations at the time of assay.

Standard	Optical Units (450 nm)
Standard A (0 pg/mL)	2.040
Standard B (20 pg/mL)	1.634
Standard C (80 pg/mL)	0.966
Standard D (200 pg/mL)	0.516
Standard E (500 pg/mL)	0.223
Standard F (1000 pg/mL)	0.130

✓ **Final Calculation for Urine Samples**

Calculate the 24 hours excretion for each urine sample: $\mu\text{g}/24 \text{ h} = \mu\text{g}/\text{L} \times \text{L}/24 \text{ h}$

Example:

Concentration for urine sample read from the standard curve = 500 pg/mL

Result after correction with the dilution factor 19 = 9500 pg/mL

$9500 \text{ pg/L} / 1000 = 9.5 \text{ } \mu\text{g/L}$

Total volume of 24 h-urine = 1.3 L (example)

$9.5 \text{ } \mu\text{g/L} \times 1.3 \text{ L}/24 \text{ h} = 12.35 \text{ } \mu\text{g}/24 \text{ h}$

✓ Excepted Normal Values

It is strongly recommended that each laboratory should determine its own normal and abnormal values.

- Serum/Plasma

In a study conducted with EDTA plasma samples of apparently normal healthy adults, using the Aldosterone ELISA the following values are observed:

Health Adults	n	Mean (pg/mL)	Median (pg/mL)	2.5th - 97.5 th Percentile (pg/mL)	Range (min - max) (pg/mL)
Supine position	60	56.14	39.71	14.21 - 156.47	8.58 – 272.30
Upright position	60	77.48	58.00	13.37 – 233.55	12.87 – 358.50

These values are also valid for serum, heparin plasma and citrate plasma.

These results correspond well to published reference ranges ^{8,9}.

In a study conducted with apparently normal healthy adults, using the Aldosterone ELISA and the Renin ELISA the following Aldosterone-Renin Ratios were determined in plasma:

Ratio Aldosterone Renin (pg/mL / pg/mL)

	n	Mean	Median	2.5 th - 97.5 th Percentile
Healthy Adults	89	8.68	5.30	0.52 – 37.83

- Urine samples

In a study conducted with urine samples of apparently normal healthy adults, using the Aldosterone ELISA the following values are observed:

	n	Mean (µg/24 h)	Median (µg/24 h)	2.5th - 97.5 th Percentile (µg/24 h)	Range (min - max) (µg/24 h)
Health Adults	8	11.34	9.40	3.31 – 25.09	3.06 – 27.17

These results correspond well to published reference ranges ⁸.

✓ Quality Control

Good laboratory practice requires that controls be run with each standard curve. A statistically significant number of controls should be assayed to establish mean values and acceptable ranges to assure proper performance.

It is recommended to use control samples according to state and federal regulations. The use of control samples is advised to assure the day to day validity of results. Use controls at both normal and pathological levels.

The controls and the corresponding results of the QC-Laboratory are stated in the QC Report added to the kit. The values and ranges stated on the QC sheet always refer to the current kit lot and should be used for direct comparison of the results.

It is also recommended to make use of national or international Quality Assessment programs in order to ensure the accuracy of the results.

Employ appropriate statistical methods for analysing control values and trends. If the results of the assay

do not fit to the established acceptable ranges of control materials results should be considered invalid. In this case, please check the following technical areas: Pipetting and timing devices; photometer, expiration dates of reagents, storage and incubation conditions, aspiration and washing methods. After checking the above mentioned items without finding any error contact your distributor or the manufacturer directly.

Performance Characteristics

✓ **Assay Dynamic Range**

For serum and plasma the range of the assay is between 7.75 pg/mL – 1000 pg/mL.

For urine the range of the assay is between 9.81 pg/mL – 1000 pg/mL.

✓ **Specificity of Antibodies (Cross Reactivity)**

The following substances were tested for cross reactivity of the assay:

Substance	Conc. Range of spiked substance	Mean cross-reactivity (%)
11-Deoxy Cortisol	10 – 1000 ng/mL	0.01
17-OH progesterone	1.2 – 120 ng/mL	0.00
21-OH progesterone	3.5 – 350 ng/mL	0.04
Androstenedione	0.2 – 22 ng/mL	0.0
Androsterone	10 – 1000 ng/mL	0.00
BSA	1 – 100 mg/mL	0.00
Cholesterol	0.5 – 50 mg/mL	0.00
Corticosterone	0.5 – 50 ng/mL	0.15
Cortisol	8 – 800 ng/mL	0.0
Cortisone	16 – 1600 ng/mL	0.01
Creatinine	50 – 5000 ug/mL	0.00
DHEA	0.5 – 50 ng/mL	0.03
DHEA-S	300 -30000 ng/mL	0.00
Estradiol	0.02 – 2 ng/mL	0.01
Estriol	1.5 – 150 ng/mL	0.00
Estrone	0.01 – 1 ng/mL	0.05
Glucose	1 – 100 mg/mL	0.00
Prednisolone	35 – 3500 ng/mL	0.00
Prednisone	35 – 3500 ng/mL	0.00
Pregnenolone	35 – 3500 ng/mL	0.00
Progesterone	42.2 – 4220 ng/mL	0.00
Testosterone	0.01 – 1 ng/mL	0.00

✓ Sensitivity

	Serum(pg/mL)	Urine (pg/mL)
The Limit of Blank (LoB)	5.359	
The Limit of Detection (LoD)	7.374	8.902
The Limit of Quantification (LoQ)	10.647	15.665
Measuring range	7.374 – 1000	8.902 – 1000
Linear range	16.37 – 1000	23.73 - 1000

✓ Reproducibility

Intra Assay

The within assay variability is shown below:

Sample	n	Mean (pg/mL)	CV %
Serum 1	5	55.66	6.5
Serum 2	5	88.88	6.7
Serum 3	5	325.82	3.8
Serum 4	5	709.07	4.5
Urine 1	5	88.47	4.3
Urine 2	5	167.59	6.4
Urine 3	5	343.98	5.5
Urine 4	5	634.55	5.9

Inter Assay

The within assay variability is shown below:

Sample	n	Mean (pg/mL)	CV %
Serum 1	25	55.66	7.4
Serum 2	25	88.88	9.8
Serum 3	25	325.82	6.5
Serum 4	25	709.07	9.6
Urine 1	25	88.47	8.2
Urine 2	25	167.59	7.3
Urine 3	25	343.98	7.5
Urine 4	25	634.55	8.3

✓ Recovery

Samples have been spiked by adding aldosterone solutions with known concentrations in a 1:1 ratio. The % recovery has been calculated by multiplication of the ratio of the measurements and the expected values with 100 (expected value = (endogenous aldosterone + added aldosterone) / 2; because of a 1:2 dilution of serum with spike material).

	Serum 1	Serum 2	Serum 3	Serum 4	Urine 1	Urine 2	Urine 3	Urine 4
Concentration [pg/mL]	186.56	282.32	445.44	589.33	56.28	108.41	215.98	240.38
Average Recovery	87.7	90.5	93.5	94.5	108.4	105.4	96.9	111.6
Range of Recovery from	86	87.9	92.6	90.9	106.8	104.0	88.8	108.8
[%] to	89.5	94.2	95.5	98.2	110.1	108.0	104.6	113.8

✓ Linearity

	Serum 1	Serum 2	Serum 3	Serum 4	Urine 1	Urine 2	Urine 3	Urine 4
Concentration [pg/mL]	85.9	328.62	514.73	537.69	96.87	127	379.7	592.9
Average Recovery	111	106.2	96.8	98.5	100.3	87.6	90.5	108.9
Range of Recovery from	108.5	102.2	93.2	88	90.7	87.6	85.8	106.5
[%] to	112.3	108.9	101.8	107.7	110.8	87.7	93.5	110.2

Resources

References

1. Brown RD, Strott CA, and Liddle GW. Site of stimulation of Aldosterone biosynthesis by angiotensin and potassium. *J Clin Invest.* (1972), 51 (6), 1413–8.
2. Bauer JH, Gauntner WC. Effect of potassium chloride on plasma renin activity and plasma aldosterone during sodium restriction in normal man. *Kidney Int.* (1979), 15 (3): 286–93.
3. Williams GH, Dluhy RG. Aldosterone biosynthesis. Interrelationship of regulatory factors. *Am J Med.* (1972), 53 (5), 595–605.
4. Tiu SC et al. The use of aldosterone-renin ratio as a diagnostic test for primary hyperaldosteronism and its test characteristics under different conditions of blood sampling. *J Clin Endocrinol Metab.* (2005), 90 (1), 72-8.
5. Mulatero P et al. Confirmatory tests in the diagnosis of primary aldosteronism. *Horm Metab Res.* (2010), 42 (6), 406-10.
6. Quillo AR. Primary aldosteronism: results of adrenalectomy for nonsingle adenoma. *J Am Coll Surg.* (2011), 213 (1), 106-12.
7. Grossmann C and Gekle M. New aspects of rapid aldosterone signaling. *Mol Cell Endocrinology* (2009), 308 (1-2), 53-62.
8. Thomas L (editor). *Renin-Angiotensin-Aldosteron-System (RAAS). Labor und Diagnose* (2005); 1406-24.
9. Perschel FH et al. Rapid Screening test for primary hyperaldosteronism: ratio of plasma aldosterone to renin concentration determined by fully automated chemiluminescence immunoassays. *Clin. Chemistry* (2004); 50 (9), 1650-55.

Plate Layout

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