



Aldosterone ELISA Kit

Item No. 501090

www.caymanchem.com

Customer Service 800.364.9897

Technical Support 888.526.5351

1180 E. Ellsworth Rd • Ann Arbor, MI • USA

TABLE OF CONTENTS

| | | |
|------------------------------|----|--|
| GENERAL INFORMATION | 3 | Materials Supplied |
| | 4 | Safety Data |
| | 4 | Precautions |
| | 5 | If You Have Problems |
| | 5 | Storage and Stability |
| | 5 | Materials Needed but Not Supplied |
| INTRODUCTION | 6 | Background |
| | 6 | About This Assay |
| | 7 | Description of AChE Competitive ELISAs |
| | 8 | Biochemistry of Acetylcholinesterase |
| | 10 | Definition of Key Terms |
| PRE-ASSAY PREPARATION | 11 | Buffer Preparation |
| | 12 | Sample Preparation |
| ASSAY PROTOCOL | 17 | Preparation of Assay-Specific Reagents |
| | 21 | Plate Set Up |
| | 22 | Performing the Assay |
| ANALYSIS | 25 | Calculations |
| | 27 | Performance Characteristics |
| RESOURCES | 32 | Troubleshooting |
| | 33 | References |
| | 34 | Plate Template |
| | 35 | Notes |
| | 35 | Warranty and Limitation of Remedy |

GENERAL INFORMATION

Materials Supplied

| Item Number | Item | 96 wells Quantity/Size | 480 wells Quantity/Size |
|-------------|------------------------------------|------------------------|-------------------------|
| 401092 | Aldosterone Polyclonal Antiserum | 1 vial/100 dtn | 1 vial/500 dtn |
| 401090 | Aldosterone AChE Tracer | 1 vial/100 dtn | 1 vial/500 dtn |
| 401094 | Aldosterone ELISA Standard | 1 vial | 1 vial |
| 400060 | ELISA Buffer Concentrate (10X) | 2 vials/10 ml | 4 vials/10 ml |
| 400062 | Wash Buffer Concentrate (400X) | 1 vial/5 ml | 1 vial/12.5 ml |
| 400035 | Polysorbate 20 | 1 vial/3 ml | 1 vial/3 ml |
| 400004 | Mouse Anti-Rabbit IgG Coated Plate | 1 plate | 5 plates |
| 400012 | 96-Well Cover Sheet | 1 cover | 5 covers |
| 400050 | Ellman's Reagent | 3 vials/100 dtn | 6 vials/250 dtn |
| 400040 | ELISA Tracer Dye | 1 vial | 1 vial |
| 400042 | ELISA Antiserum Dye | 1 vial | 1 vial |

If any of the items listed above are damaged or missing, please contact our Customer Service department at (800) 364-9897 or (734) 971-3335. We cannot accept any returns without prior authorization.



WARNING: THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

Safety Data

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent *via* email to your institution.

Precautions

Please read these instructions carefully before beginning this assay.

The reagents in this kit have been tested and formulated to work exclusively with Cayman Chemical's Aldosterone ELISA Kit. This kit may not perform as described if any reagent or procedure is replaced or modified.

When compared to quantification by LC/MS or GC/MS, it is not uncommon for immunoassays to report higher analyte concentrations. While LC/MS or GC/MS analyses typically measure only a single compound, antibodies used in immunoassays sometimes recognize not only the target molecule, but also structurally related molecules, including biologically relevant metabolites. In many cases, measurement of both the parent molecule and metabolites is more representative of the overall biological response than is the measurement of a short-lived parent molecule. It is the responsibility of the researcher to understand the limits of both assay systems and to interpret their data accordingly.

If You Have Problems

Technical Service Contact Information

Phone: 888-526-5351 (USA and Canada only) or 734-975-3888

Email: techserv@caymanchem.com

In order for our staff to assist you quickly and efficiently, please be ready to supply the lot number of the kit (found on the outside of the box).

Storage and Stability

This kit will perform as specified if stored as directed at -20°C and used before the expiration date indicated on the outside of the box.

Materials Needed But Not Supplied

1. A plate reader capable of measuring absorbance between 405-420 nm.
2. Adjustable pipettes and a repeating pipettor.
3. A source of 'UltraPure' water. Water used to prepare all ELISA reagents and buffers must be deionized and free of trace organic contaminants ('UltraPure'). Use activated carbon filter cartridges or other organic scavengers. Glass distilled water (even if double distilled), HPLC-grade water, and sterile water (for injections) are not adequate for ELISA. *NOTE: UltraPure water is available for purchase from Cayman (Item No. 400000).*
4. Materials used for **Sample Preparation** (see page 12).

Background

Aldosterone, a steroid hormone secreted by the adrenal cortex, is the principle mineralocorticoid controlling sodium and potassium balance.^{1,2} The primary role of aldosterone is to promote unidirectional salt reabsorption in a variety of epithelial tissues, the salivary gland, intestine, sweat glands, and the kidney. Aldosterone is synthesized from cholesterol in the zona glomerulosa of the adrenal cortex. Secretion of aldosterone is complex, being regulated by both hormones and electrolytes. However, the renin-angiotensin system (RAS) is the primary regulator of aldosterone secretion.³ Angiotensin II and potassium stimulate secretion of aldosterone by increasing the rate of synthesis of the hormone. Serum aldosterone levels vary depending on age and body posture, *i.e.*, supine *versus* upright. In normal upright adult individuals, serum aldosterone concentrations are generally less than 300 pg/ml. Only a fraction of urinary aldosterone is excreted intact, with larger percentages excreted as aldosterone-18-glucuronide and tetrahydroaldosterone-3-glucuronide.⁴ Urinary excretion of aldosterone in normal adults, assessed as a combination of free aldosterone and aldosterone-18-glucuronide, is generally less than 20 µg/24 hr.^{5,6} The average reported level of tetrahydroaldosterone in urine is 103 nmol/L (37 ng/ml) or less than 70 µg/24 hr.^{5,6}

About This Assay

Cayman's Aldosterone ELISA Kit is a competitive assay that can be used for quantification of aldosterone in serum, plasma, urine, and other sample matrices. The assay has a range from 15.6-2,000 pg/ml and a sensitivity (80% B/B₀) of approximately 30 pg/ml.

Description of AChE Competitive ELISAs

This assay is based on the competition between aldosterone and a aldosterone-acetylcholinesterase (AChE) conjugate (Aldosterone Tracer) for a limited amount of Aldosterone Polyclonal Antiserum. Because the concentration of the Aldosterone Tracer is held constant while the concentration of aldosterone varies, the amount of Aldosterone Tracer that is able to bind to the Aldosterone Polyclonal Antiserum will be inversely proportional to the concentration of aldosterone in the well. This antiserum-aldosterone complex binds to mouse anti-rabbit IgG that has been previously attached to the well. The plate is washed to remove any unbound reagents and then Ellman's Reagent (which contains the substrate to AChE) is added to the well. The product of this enzymatic reaction has a distinct yellow color and absorbs strongly at 412 nm. The intensity of this color, determined spectrophotometrically, is proportional to the amount of Aldosterone Tracer bound to the well, which is inversely proportional to the amount of free aldosterone present in the well during the incubation; or

$$\text{Absorbance} \propto [\text{Bound Aldosterone Tracer}] \propto 1/[\text{Aldosterone}]$$

A schematic of this process is shown in Figure 1, below.

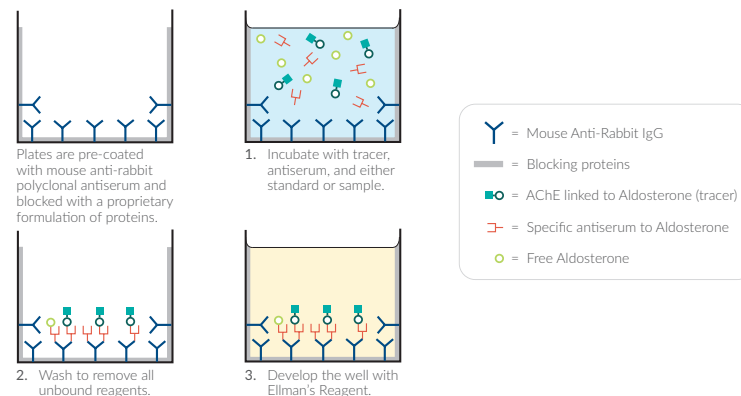


Figure 1. Schematic of the AChE ELISA

Biochemistry of Acetylcholinesterase

The electric organ of the electric eel, *E. electricus*, contains an avid AChE capable of massive catalytic turnover during the generation of its electrochemical discharges. The electric eel AChE has a clover leaf-shaped tertiary structure consisting of a triad of tetramers attached to a collagen-like structural fibril. This stable enzyme is capable of high turnover ($64,000 \text{ s}^{-1}$) for the hydrolysis of acetylthiocholine.

A molecule of the analyte covalently attached to a molecule of AChE serves as the tracer in AChE enzyme immunoassays. Quantification of the tracer is achieved by measuring its AChE activity with Ellman's Reagent. This reagent consists of acetylthiocholine and 5,5'-dithio-bis-(2-nitrobenzoic acid). Hydrolysis of acetylthiocholine by AChE produces thiocholine (see Figure 2, on page 9). The non-enzymatic reaction of thiocholine with 5,5'-dithio-bis-(2-nitrobenzoic acid) produces 5-thio-2-nitrobenzoic acid, which has a strong absorbance at 412 nm ($\epsilon = 13,600$).

AChE has several advantages over other enzymes commonly used for enzyme immunoassays. Unlike horseradish peroxidase, AChE does not self-inactivate during turnover. This property of AChE also allows redevelopment of the assay if it is accidentally splashed or spilled. In addition, the enzyme is highly stable under the assay conditions, has a wide pH range (pH 5-10), and is not inhibited by common buffer salts or preservatives. Since AChE is stable during the development step, it is unnecessary to use a 'stop' reagent, and the plate may be read whenever it is convenient.

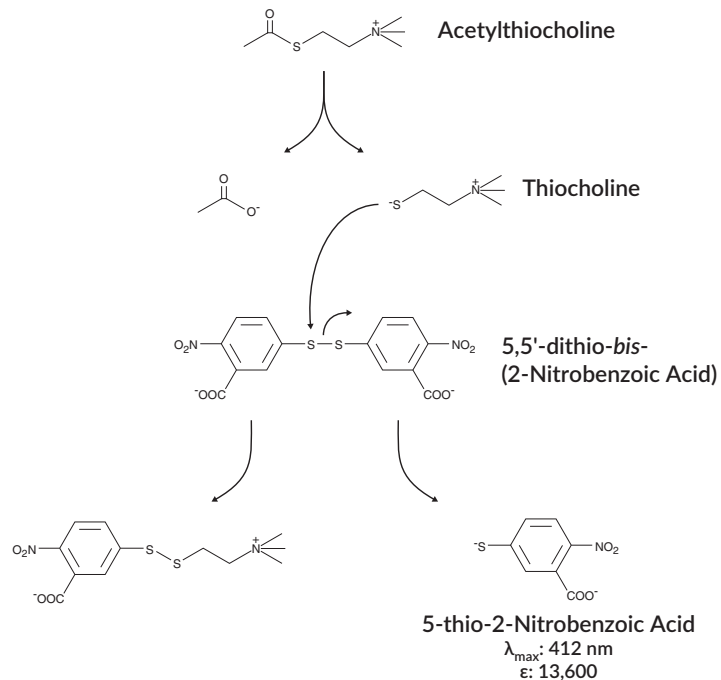


Figure 2. Reaction catalyzed by acetylcholinesterase

Definition of Key Terms

Blank: background absorbance caused by Ellman's Reagent. The blank absorbance should be subtracted from the absorbance readings of all the other wells, including NSB wells.

Total Activity: total enzymatic activity of the AChE-linked tracer. This is analogous to the specific activity of a radioactive tracer.

NSB (Non-Specific Binding): non-immunological binding of the tracer to the well. Even in the absence of specific antibody a very small amount of tracer still binds to the well; the NSB is a measure of this low binding. Do not forget to subtract the Blank absorbance values.

B₀ (Maximum Binding): maximum amount of the tracer that the antibody can bind in the absence of free analyte.

%B/B₀ (%Bound/Maximum Bound): ratio of the absorbance of a particular sample or standard well to that of the maximum binding (B₀) well.

Standard Curve: a plot of the %B/B₀ values *versus* concentration of a series of wells containing various known amounts of analyte.

Dtn: determination, where one dtn is the amount of reagent used per well.

Cross Reactivity: numerical representation of the relative reactivity of this assay towards structurally related molecules as compared to the primary analyte of interest. Biomolecules that possess similar epitopes to the analyte can compete with the assay tracer for binding to the primary antibody. Substances that are superior to the analyte in displacing the tracer result in a cross reactivity that is greater than 100%. Substances that are inferior to the primary analyte in displacing the tracer result in a cross reactivity that is less than 100%. Cross reactivity is calculated by comparing the mid-point (50% B/B₀) value of the tested molecule to the mid-point (50% B/B₀) value of the primary analyte when each is measured in assay buffer using the following formula:

$$\% \text{ Cross Reactivity} = \left[\frac{50\% \text{ B/B}_0 \text{ value for the primary analyte}}{50\% \text{ B/B}_0 \text{ value for the potential cross reactant}} \right] \times 100\%$$

PRE-ASSAY PREPARATION

NOTE: Water used to prepare all ELISA reagents and buffers must be deionized and free of trace organic contaminants ('UltraPure'). Use activated carbon filter cartridges or other organic scavengers. Glass distilled water (even if double distilled), HPLC-grade water, and sterile water (for injections) are not adequate for ELISA. UltraPure water may be purchased from Cayman (Item No. 400000).

Buffer Preparation

Store all diluted buffers at 4°C; they will be stable for about two months.

1. ELISA Buffer Preparation

Dilute the contents of one vial of ELISA Buffer Concentrate (10X) (Item No. 400060) with 90 ml of UltraPure water. Be certain to rinse the vial to remove any salts that may have precipitated. *NOTE: It is normal for the concentrated buffer to contain crystalline salts after thawing. These will completely dissolve upon dilution with water.*

2. Wash Buffer Preparation

5 ml vial Wash Buffer Concentration (400X) (96-well kit; Item No. 400062): Dilute to a total volume of 2 liters with UltraPure water and add 1 ml of Polysorbate 20 (Item No. 400035).

OR

12.5 ml vial Wash Buffer Concentration (400X) (480-well kit; Item No. 400062): Dilute to a total volume of 5 liters with UltraPure water and add 2.5 ml of Polysorbate 20 (Item No. 400035).

Smaller volumes of Wash Buffer can be prepared by diluting the Wash Buffer Concentrate 1:400 and adding Polysorbate 20 (0.5 ml/liter of Wash Buffer).

NOTE: Polysorbate 20 is a viscous liquid and cannot be measured by a regular pipette. A positive displacement pipette or a syringe should be used to deliver small quantities accurately.

Sample Preparation

General Precautions

- All samples must be free of organic solvents prior to assay.
- AEBSF (Pefabloc SC[®]) and PMSF inhibit acetylcholinesterase. Samples containing these protease inhibitors should not be used in this assay.
- Samples should be assayed immediately after collection; samples that cannot be assayed immediately should be stored at -20 or -80°C.

Materials Needed

1. Chloroform (for plasma and serum samples)
2. 3.2 N Hydrochloric Acid and 1M Sodium Hydroxide for hydrolysis of the urine samples

To prepare 3.2 N HCl add 10 ml of concentrated HCl (12N) to 27 ml of purified water. To prepare 1M NaOH dissolve 4 g of dry sodium hydroxide in 50-80 ml of purified water, then add water to the total volume of 100 ml.

Serum and Plasma

Serum and plasma contain a number of steroids that may interfere with this assay.⁷ It is essential that these samples be extracted in order to obtain accurate results. Failure to extract samples may result in spurious data.

Extraction Protocol for Serum and Plasma

The following protocol is recommended for extraction of serum and plasma samples.

NOTE: We do not recommend the use of plastic vials, caps, or pipettes for this procedure. Chloroform may extract interfering compounds from the plastic.

1. Aliquot a known amount of each sample into a clean test tube (600 μ l is recommended).
2. Add 4X the sample volume of chloroform and mix thoroughly with a vortexer. Allow layers to separate. Centrifugation of the samples at 30 x g for 5-10 minutes will ensure a better separation of the layers and also lower variability in the assay. Using a Pasteur pipette, transfer 90% of the chloroform (lower layer) into a clean test tube. Repeat this extraction procedure three times.*
3. Evaporate the combined chloroform extracts by heating to 30°C under a gentle stream of nitrogen.
4. Dissolve the extract in 0.6 ml of ELISA Buffer.

*If it is necessary to stop during this purification, samples may be stored in the chloroform solution at -20°C.

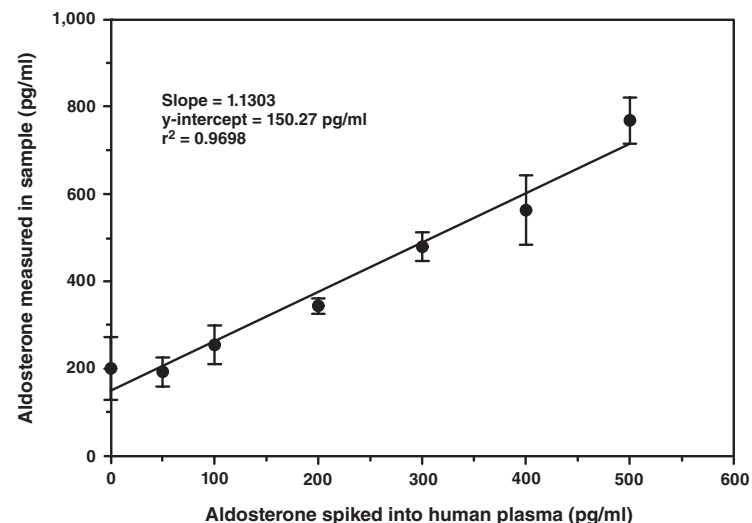


Figure 3. Recovery of aldosterone from plasma

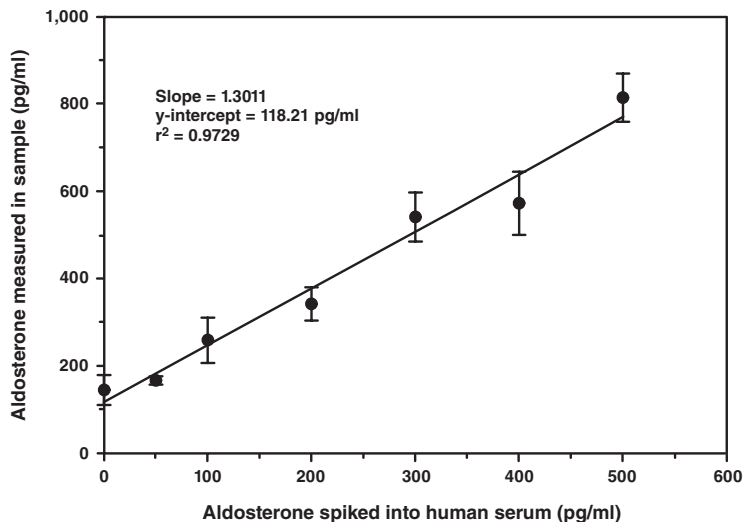


Figure 4. Recovery of aldosterone from serum

Urine

Urinary aldosterone is found in three forms: free aldosterone, aldosterone-18-glucuronide, and tetrahydroaldosterone-glucuronide. The antibody in this kit recognizes free aldosterone only; it does not recognize aldosterone-18-glucuronide or tetrahydroaldosterone (see **Specificity** on page 31). Treatment of the urine with hydrochloric acid (acid hydrolysis) converts the aldosterone-18-glucuronide into free aldosterone, allowing it to be detected in this assay.

Acid Hydrolysis Protocol for Urine

The following protocol is recommended for measurement of free aldosterone in urine samples.

1. Aliquot a known amount of each sample into a clean test tube (1,000 μ l is recommended).
2. Add 100 μ l of 3.2 N HCl, mix well and incubate overnight at room temperature.
3. Neutralize each sample with 225 μ l of 1 M NaOH.
4. Dilute at least 10 times with ELISA Buffer prior to the assay.

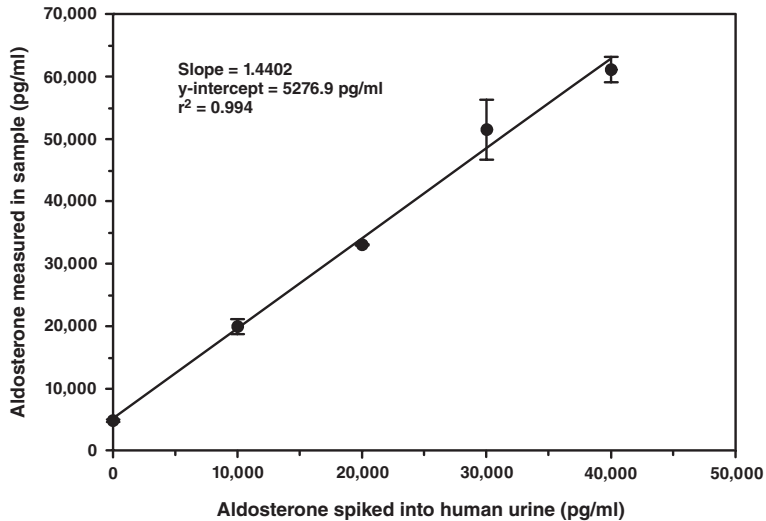


Figure 5. Recovery of aldosterone from urine

ASSAY PROTOCOL

Preparation of Assay-Specific Reagents

Aldosterone ELISA Standard

Equilibrate a pipette tip in ethanol by repeatedly filling and expelling the tip with ethanol several times. Using the equilibrated pipette tip, transfer 100 μ l of the Aldosterone ELISA Standard (Item No. 401094) into a clean test tube, then dilute with 900 μ l UltraPure water. The concentration of this solution (the bulk standard) will be 20 ng/ml.

NOTE: If assaying culture medium samples that have not been diluted with ELISA Buffer, culture medium should be used in place of ELISA Buffer for dilution of the standard curve.

To prepare the standard for use in ELISA: Obtain eight clean test tubes and number them #1 through #8. Aliquot 900 μ l ELISA Buffer to tube #1 and 500 μ l ELISA Buffer to tubes #2-8. Transfer 100 μ l of the bulk standard (20 ng/ml) to tube #1 and mix thoroughly. The concentration of this standard, the first point on the standard curve, will be 2 ng/ml (2,000 pg/ml). Serially dilute the standard by removing 500 μ l from tube #1 and placing in tube #2; mix thoroughly. Next, remove 500 μ l from tube #2 and place it into tube #3; mix thoroughly. Repeat this process for tubes #4-8. These diluted standards should not be stored for more than 24 hours.

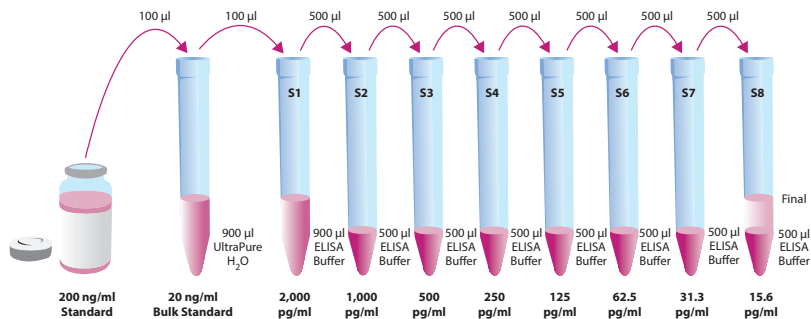


Figure 6. Preparation of the Aldosterone standards

Aldosterone AChE Tracer

Reconstitute the Aldosterone AChE Tracer as follows:

100 dtn Aldosterone AChE Tracer (96-well kit; Item No. 401090):
Reconstitute with 6 ml ELISA Buffer.

OR

500 dtn Aldosterone AChE Tracer (480-well kit; Item No. 401090):
Reconstitute with 30 ml ELISA Buffer.

Store the reconstituted Aldosterone AChE Tracer at 4°C (*do not freeze!*) and use within eight weeks. A 20% surplus of tracer has been included to account for any incidental losses.

Tracer Dye Instructions (optional)

This dye may be added to the tracer, if desired, to aid in visualization of tracer-containing wells. Add the dye to the reconstituted tracer at a final dilution of 1:100 (add 60 µl of dye to 6 ml tracer or add 300 µl of dye to 30 ml of tracer). Tracer with dye will be stable for at least five weeks.

Aldosterone Polyclonal Antiserum

Reconstitute the Aldosterone Polyclonal Antiserum as follows:

100 dtn Aldosterone Polyclonal Antiserum (96-well kit; Item No. 401092):

Reconstitute with 6 ml ELISA Buffer.

OR

500 dtn Aldosterone Polyclonal Antiserum (480-well kit; Item No. 401092):

Reconstitute with 30 ml ELISA Buffer.

Store the reconstituted Aldosterone Polyclonal Antiserum at 4°C. It will be stable for at least eight weeks. A 20% surplus of antibody has been included to account for any incidental losses.

Antiserum Dye Instructions (optional)

This dye may be added to the antibody, if desired, to aid in visualization of antibody-containing wells. Add the dye to the reconstituted antibody at a final dilution of 1:100 (add 60 µl of dye to 6 ml antibody or add 300 µl of dye to 30 ml of antibody). Antiserum with dye will be stable for at least five weeks.

Plate Set Up

The 96-well plate(s) included with this kit is supplied ready to use. It is not necessary to rinse the plate(s) prior to adding the reagents. *NOTE: If you do not need to use all the strips at once, place the unused strips back in the plate packet and store at 4°C. Be sure the packet is sealed with the desiccant inside.*

Each plate or set of strips must contain a minimum of two blanks (Blk), two non-specific binding wells (NSB), two maximum binding wells (B_0), and an eight point standard curve run in duplicate. *NOTE: Each assay must contain this minimum configuration in order to ensure accurate and reproducible results.* Each sample should be assayed at two dilutions and each dilution should be assayed in duplicate. For statistical purposes, we recommend assaying samples in triplicate.

A suggested plate format is shown in Figure 7, below. The user may vary the location and type of wells present as necessary for each particular experiment. The plate format provided below has been designed to allow for easy data analysis using a convenient spreadsheet offered by Cayman (see page 25, for more details). We suggest you record the contents of each well on the template sheet provided (see page 34).

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|-------|----|----|---|---|---|----|----|----|----|----|----|
| A | Blk | S1 | S1 | 1 | 1 | 1 | 9 | 9 | 9 | 17 | 17 | 17 |
| B | Blk | S2 | S2 | 2 | 2 | 2 | 10 | 10 | 10 | 18 | 18 | 18 |
| C | NSB | S3 | S3 | 3 | 3 | 3 | 11 | 11 | 11 | 19 | 19 | 19 |
| D | NSB | S4 | S4 | 4 | 4 | 4 | 12 | 12 | 12 | 20 | 20 | 20 |
| E | B_0 | S5 | S5 | 5 | 5 | 5 | 13 | 13 | 13 | 21 | 21 | 21 |
| F | B_0 | S6 | S6 | 6 | 6 | 6 | 14 | 14 | 14 | 22 | 22 | 22 |
| G | B_0 | S7 | S7 | 7 | 7 | 7 | 15 | 15 | 15 | 23 | 23 | 23 |
| H | TA | S8 | S8 | 8 | 8 | 8 | 16 | 16 | 16 | 24 | 24 | 24 |

Blk - Blank
TA - Total Activity
NSB - Non-Specific Binding
 B_0 - Maximum Binding
S1-S8 - Standards 1-8
1-24 - Samples

Figure 7. Sample plate format

Performing the Assay

Pipetting Hints

- Use different tips to pipette each reagent.
- Before pipetting each reagent, equilibrate the pipette tip in that reagent (*i.e.*, slowly fill the tip and gently expel the contents, repeat several times).
- Do not expose the pipette tip to the reagent(s) already in the well.

Addition of the Reagents

1. ELISA Buffer

Add 150 μ l ELISA Buffer to NSB wells. Add 100 μ l ELISA Buffer to B₀ wells. If culture medium was used to dilute the standard curve, substitute 100 μ l of culture medium for ELISA Buffer in the NSB and B₀ wells (*i.e.*, add 100 μ l culture medium to NSB and B₀ wells and 50 μ l ELISA Buffer to NSB wells).

2. Aldosterone ELISA Standard

Add 100 μ l from tube #8 to both of the lowest standard wells (S8). Add 100 μ l from tube #7 to each of the next two standard wells (S7). Continue with this procedure until all the standards are aliquoted. The same pipette tip should be used to aliquot all the standards. Before pipetting each standard, be sure to equilibrate the pipette tip in that standard.

3. Samples

Add 100 μ l of sample per well. Each sample should be assayed at a minimum of two dilutions. Each dilution should be assayed in duplicate (triplicate recommended).

4. Aldosterone AChE Tracer

Add 50 μ l to each well except the TA and the Blk wells.

5. Aldosterone Polyclonal Antiserum

Add 50 μ l to each well except the TA, the NSB, and the Blk wells.

| Well | ELISA Buffer | Standard/Sample | Tracer | Antiserum |
|----------------|--------------|-----------------|----------------------------|------------|
| Blk | - | - | - | - |
| TA | - | - | 5 μ l (at devel. step) | - |
| NSB | 150 μ l | - | 50 μ l | - |
| B ₀ | 100 μ l | - | 50 μ l | 50 μ l |
| Std/Sample | - | 100 μ l | 50 μ l | 50 μ l |

Table 1. Pipetting summary

Incubation of the Plate

Cover each plate with plastic film (Item No. 400012) and incubate 18 hours at 4°C.

Development of the Plate

1. Reconstitute Ellman's Reagent immediately before use (20 ml of reagent is sufficient to develop 100 wells):

100 dtn vial Ellman's Reagent (96-well kit; Item No. 400050): Reconstitute with 20 ml of UltraPure water.

OR

250 dtn vial Ellman's Reagent (480-well kit; Item No. 400050): Reconstitute with 50 ml of UltraPure water.

NOTE: Reconstituted Ellman's Reagent is unstable and should be used the same day it is prepared; protect the Ellman's Reagent from light when not in use. Extra vials of the reagent have been provided should a plate need to be re-developed or multiple assays run on different days.

2. Empty the wells and rinse five times with Wash Buffer.
3. Add 200 μl of Ellman's Reagent to each well.
4. Add 5 μl of tracer to the TA wells.
5. Cover the plate with plastic film. Optimum development is obtained by using an orbital shaker equipped with a large, flat cover to allow the plate(s) to develop in the dark. This assay typically develops (i.e., B_0 wells ≥ 0.3 A.U. (blank subtracted)) in 90-120 minutes.

Reading the Plate

1. Wipe the bottom of the plate with a clean tissue to remove fingerprints, dirt, etc.
2. Remove the plate cover being careful to keep Ellman's Reagent from splashing on the cover. *NOTE: Any loss of Ellman's Reagent will affect the absorbance readings. If Ellman's Reagent is present on the cover, use a pipette to transfer the Ellman's Reagent into the well. If too much Ellman's Reagent has splashed on the cover to easily redistribute back into the wells, wash the plate three times with wash buffer and repeat the development with fresh Ellman's Reagent.*
3. Read the plate at a wavelength between 405 and 420 nm. The absorbance may be checked periodically until the B_0 wells have reached a minimum of 0.3 A.U. (blank subtracted). The plate should be read when the absorbance of the B_0 wells are in the range of 0.3-1.0 A.U. (blank subtracted). If the absorbance of the wells exceeds 1.5, wash the plate, add fresh Ellman's Reagent and let it develop again.

ANALYSIS

Many plate readers come with data reduction software that plot data automatically. Alternatively a spreadsheet program can be used. The data should be plotted as either $\%B/B_0$ versus log concentration using a four-parameter logistic fit or as logit B/B_0 versus log concentration using a linear fit. *NOTE: Cayman has a computer spreadsheet available for data analysis. Please contact Technical Service or visit our website (www.caymanchem.com/analysis/elisa) to obtain a free copy of this convenient data analysis tool.*

Calculations

Preparation of the Data

The following procedure is recommended for preparation of the data prior to graphical analysis.

NOTE: If the plate reader has not subtracted the absorbance readings of the blank wells from the absorbance readings of the rest of the plate, be sure to do that now.

1. Average the absorbance readings from the NSB wells.
2. Average the absorbance readings from the B_0 wells.
3. Subtract the NSB average from the B_0 average. This is the corrected B_0 or corrected maximum binding.
4. Calculate the B/B_0 (Sample or Standard Bound/Maximum Bound) for the remaining wells. To do this, subtract the average NSB absorbance from the S1 absorbance and divide by the corrected B_0 (from Step 3). Repeat for S2-S8 and all sample wells. (To obtain $\%B/B_0$ for a logistic four-parameter fit, multiply these values by 100.)

NOTE: The TA values are not used in the standard curve calculations. Rather, they are used as a diagnostic tool; the corrected B_0 divided by the actual TA (10X measured absorbance) will give the %Bound. This value should closely approximate the %Bound that can be calculated from the Sample Data (see page 27). Erratic absorbance values and a low (or no) %Bound could indicate the presence of organic solvents in the buffer or other technical problems (see page 32 for Troubleshooting).

Plot the Standard Curve

Plot %B/B₀ for standards S1-S8 versus aldosterone concentration using linear (y) and log (x) axes and perform a 4-parameter logistic fit.

Alternative Plot - The data can also be linearized using a logit transformation. The equation for this conversion is shown below. *NOTE: Do not use %B/B₀ in this calculation.*

$$\text{logit (B/B}_0\text{)} = \ln [\text{B/B}_0\text{}/(1 - \text{B/B}_0\text{)}]$$

Plot the data as logit (B/B₀) versus log concentrations and perform a linear regression fit.

Determine the Sample Concentration

Calculate the B/B₀ (or %B/B₀) value for each sample. Determine the concentration of each sample using the equation obtained from the standard curve plot. *NOTE: Remember to account for any concentration or dilution of the sample prior to the addition to the well. Samples with %B/B₀ values greater than 80% or less than 20% should be re-assayed as they generally fall out of the linear range of the standard curve. A 20% or greater disparity between the apparent concentration of two different dilutions of the same sample indicates interference which could be eliminated by purification.*

Performance Characteristics

Sample Data

The standard curve presented here is an example of the data typically produced with this kit; however, your results will not be identical to these. You **must** run a new standard curve. Do not use the data below to determine the values of your samples. Your results could differ substantially.

| | Raw Data | | Average | Corrected |
|-----------------------|----------|-------|---------|-----------|
| Total Activity | 0.499 | 0.533 | 0.516 | |
| NSB | 0.007 | 0.007 | 0.007 | |
| B₀ | 1.224 | 1.254 | | |
| | 1.258 | 1.179 | 1.237 | 1.230 |

| Dose (pg/ml) | Raw Data | | Corrected | | %B/B ₀ | |
|--------------|----------|-------|-----------|-------|-------------------|------|
| | | | | | | |
| 2,000 | 0.212 | 0.221 | 0.205 | 0.214 | 16.7 | 17.4 |
| 1,000 | 0.288 | 0.291 | 0.281 | 0.284 | 22.8 | 23.1 |
| 500 | 0.379 | 0.377 | 0.372 | 0.370 | 30.2 | 30.1 |
| 250 | 0.485 | 0.487 | 0.478 | 0.480 | 38.9 | 39.0 |
| 125 | 0.647 | 0.656 | 0.640 | 0.649 | 52.0 | 52.8 |
| 62.5 | 0.869 | 0.907 | 0.862 | 0.900 | 70.1 | 73.2 |
| 31.3 | 0.972 | 0.939 | 0.965 | 0.932 | 78.5 | 75.8 |
| 15.6 | 1.078 | 1.014 | 1.071 | 1.014 | 87.1 | 82.4 |

Table 2. Typical results

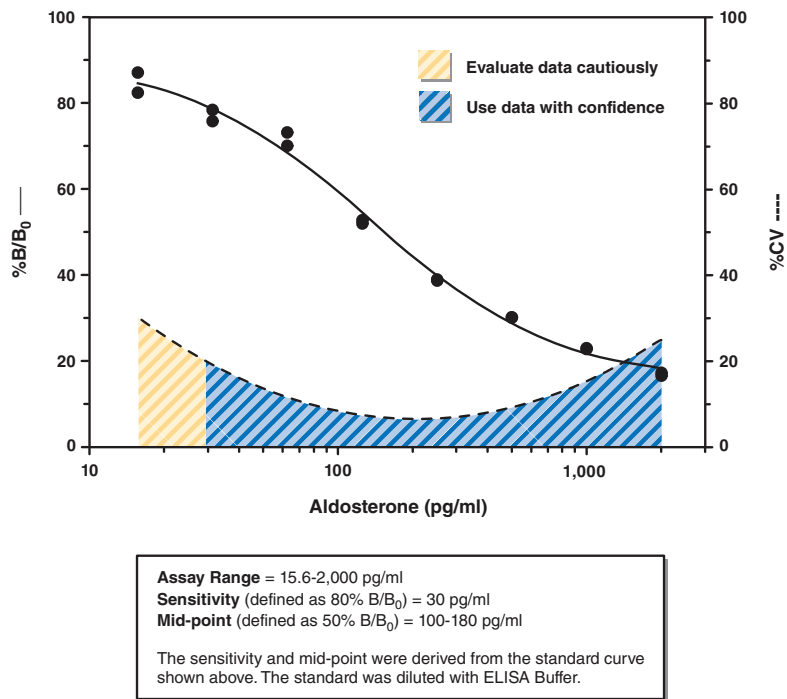


Figure 8. Typical standard curve

Precision:

The intra-assay CVs have been determined at multiple points on the standard curve. These data are summarized in the graph on page 28 and in the table below.

| Dose (pg/ml) | %CV* Intra-assay variation |
|--------------|-------------------------------|
| 2,000 | 25.9 |
| 1,000 | 13.8 |
| 500 | 8.4 |
| 250 | 8.6 |
| 125 | 8.6 |
| 62.5 | 12.0 |
| 31.3 | 15.6 |
| 15.6 | 32.2 |

Table 3. Intra-assay variation

*%CV represents the variation in concentration (not absorbance) as determined using a reference standard curve.

| Matrix | Level | Average (pg/ml) | %CV Inter-assay variation |
|--------|--------|-----------------|---------------------------|
| Plasma | High | 198.3 | 6.9 |
| | Medium | 128.3 | 9.0 |
| | Low | 71.3 | 17.4 |
| Serum | High | 255.0 | 11.5 |
| | Medium | 126.8 | 16.5 |
| | Low | 57.6 | 21.3 |
| Urine | High | 36,221.4 | 6.6 |
| | Medium | 5,107.3 | 9.2 |
| | Low | 3,016.7 | 11.1 |

Table 4. Sample Inter-assay variation for plasma, serum, and urine

Plasma, serum and urine samples from different individuals containing high, medium, or low level of aldosterone were measured eight times on four different days using a different set of reagents each time. The calculated %CV is reported as inter-assay variance.

Cross Reactivity:

| Compound | Cross Reactivity | Compound | Cross Reactivity |
|--|------------------|---------------------------------|------------------|
| Aldosterone | 100% | 11-Deoxycortisol | 0.0008% |
| 3 β ,5 β -Tetrahydroaldosterone | 41% | DHEA | 0.0008% |
| 3 α ,5 β -Tetrahydroaldosterone | 0.5457% | Cortisol | 0.0006% |
| Corticosterone | 0.1051% | Cortisone | 0.0006% |
| Progesterone | 0.0360% | DHEA Sulfate | 0.0005% |
| Pregnenolone | 0.0243% | 5 α -Dihydrotestosterone | 0.0004% |
| Androstenedione | 0.0212% | Estrone | 0.0001% |
| 11-Deoxycorticosterone | 0.0166% | Dexamethasone | <0.01% |
| Testosterone | 0.0072% | β -Estradiol | <0.01% |
| 17-Hydroxypregnenolone | 0.0027% | Estriol | <0.01% |
| 17-Hydroxyprogesterone | 0.0013% | Estrone Sulfate | <0.01% |

Table 5. Cross Reactivity of the Aldosterone ELISA

RESOURCES

Troubleshooting

| Problem | Possible Causes | Recommended Solutions |
|--|---|---|
| Erratic values; dispersion of duplicates | A. Trace organic contaminants in the water source B. Poor pipetting/technique | A. Replace activated carbon filter or change source of UltraPure water |
| High NSB (>10% of B ₀) | A. Poor washing B. Exposure of NSB wells to specific antibody | A. Rewash plate and redevelop |
| Very low B ₀ | A. Trace organic contaminants in the water source B. Plate requires additional development time C. Dilution error in preparing reagents | A. Replace activated carbon filter or change source of UltraPure water B. Return plate to shaker and re-read later |
| Low sensitivity (shift in dose response curve) | Standard is degraded | Replace standard |
| Analyses of two dilutions of a biological sample do not agree (i.e., more than 20% difference) | Interfering substances are present | Purify sample prior to analysis by ELISA ⁶ |
| Only Total Activity (TA) wells develop | Trace organic contaminants in the water source | Replace activated carbon filter or change source of UltraPure water |

References

1. Rogerson, F.M. and Fuller, P.J. Mineralocorticoid action. *Steroids* **65**, 61-73 (2000).
2. Agarwal, M.K. and Mirshahi, M. General overview of mineralocorticoid hormone action. *Pharmacol. Toxicol.* **84**, 273-326 (1999).
3. Lumbers, E.R. Angiotensin and aldosterone. *Regul. Pept.* **80**, 91-100 (1999).
4. Cartledge, S. and Lawson, N. Aldosterone and renin measurements. *Ann. Clin. Biochem.* **37**, 262-278 (2000).
5. Koopman, B.J., Lokerse, I.J.G., Verweij, H., *et al.* Improved gas chromatographic-mass fragmentographic assay for tetrahydroaldosterone and aldosterone in urine. *J. Chromatogr.* **378(2)**, 283-292 (1986).
6. Abdelhamid, S., Blomer, R., Hommel, G., *et al.* Urinary tetrahydroaldosterone as a screening method for primary aldosteronism: A comparative study. *Am. J. Hypertens.* **16(7)**, 522-530 (2003).
7. Maxey, K.M., Maddipati, K.R. and Birkmeier, J. Interference in enzyme immunoassays. *J. Clin. Immunoassay* **15**, 116-120 (1992).

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

NOTES

Warranty and Limitation of Remedy

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

This document is copyrighted. All rights are reserved. This document may not, in whole or part, be copied, photocopied, reproduced, translated, or reduced to any electronic medium or machine-readable form without prior consent, in writing, from Cayman Chemical Company.

©10/06/2022, Cayman Chemical Company, Ann Arbor, MI, All rights reserved. Printed in U.S.A.

