



2-Methoxyestradiol ELISA Kit

Item No. 582261

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 | 19 | Preparation of Assay-Specific Reagents |
| | 22 | Plate Set Up |
| | 23 | Performing the Assay |
| ANALYSIS | 26 | Calculations |
| | 28 | 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 |
|---------------|------------------------------------|---------------------------|----------------------------|
| 482262 | 2-Methoxyestradiol ELISA Antiserum | 1 vial/100 dtn | 1 vial/500 dtn |
| 482260 | 2-Methoxyestradiol AChE Tracer | 1 vial/100 dtn | 1 vial/500 dtn |
| 482264 | 2-Methoxyestradiol 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 |
| 400004/400006 | 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 AChE ELISA Kits. 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
Fax: 734-971-3641
Email: techserv@caymanchem.com
Hours: M-F 8:00 AM to 5:30 PM EST

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 including glucuronidase used for Sample Preparation (see pages 12 and 14).

Background

2-Methoxyestradiol (2-ME2) is a natural metabolite of estradiol with potent antitumor and antiangiogenic properties.¹⁻³ Several clinical trials have demonstrated the antitumor effects of 2-ME2 in prostate cancer, breast cancer, renal cell carcinoma, ovarian cancer, glioblastoma, and multiple myeloma.⁴ Other studies indicate that 2-ME2 exhibits more potent cardioprotective effects than estradiol.^{4,5}

2-ME2 is produced by cytochrome P450-dependant hydroxylation of 17 β -estradiol at the 2-position, followed by O-methylation catalyzed by catechol-O-methyltransferase. Deficiency in catechol-O-methyltransferase and 2-ME2 is associated with pre-eclampsia.⁶ Therefore, 2-ME2 may be utilized as a plasma and urinary diagnostic marker for this disease, as well as a therapeutic supplement for its prevention.⁶ 2-ME2 has very low affinity for estrogen receptors compared to estradiol and other estradiol metabolites,⁷ suggesting a different mechanism of action for this compound.

Levels of 2-ME2 in plasma vary from less than 0.1 ng/ml in males and non-pregnant females to 3.0 ng/ml in pregnant women.⁶ Urinary excretion of 2-ME2 in non-pregnant women was reported to be 0.91-2.42 nanomoles per 24 hours,^{8,9} which increases dramatically during pregnancy.¹⁰

About This Assay

Cayman's 2-Methoxyestradiol ELISA Kit is a competitive assay that can be used for quantification of 2-ME2 in plasma, urine, and other sample matrices. The assay has a range from 13.1-8,000 pg/ml and a sensitivity (80% B/B₀) of approximately 40 pg/ml.

Description of AChE Competitive ELISAs^{11,12}

This assay is based on the competition between free 2-ME2 and a 2-Methoxyestradiol Tracer (2-ME2 linked to an acetylcholinesterase (AChE) molecule) for a limited number of 2-ME2-specific rabbit antiserum binding sites. The concentration of the 2-Methoxyestradiol Tracer is held constant while the concentration of free 2-ME2 (standard or sample) varies. Thus, the amount of 2-Methoxyestradiol Tracer that is able to bind to the rabbit antiserum will be inversely proportional to the concentration of free 2-ME2 in the well. This rabbit antiserum-2-ME2 (either free or tracer) complex binds to the mouse monoclonal 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 2-Methoxyestradiol Tracer bound to the well, which is inversely proportional to the amount of free 2-ME2 present in the well during the incubation; or

$$\text{Absorbance} \propto [\text{Bound 2-Methoxyestradiol Tracer}] \propto 1/[\text{2-ME2}]$$

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

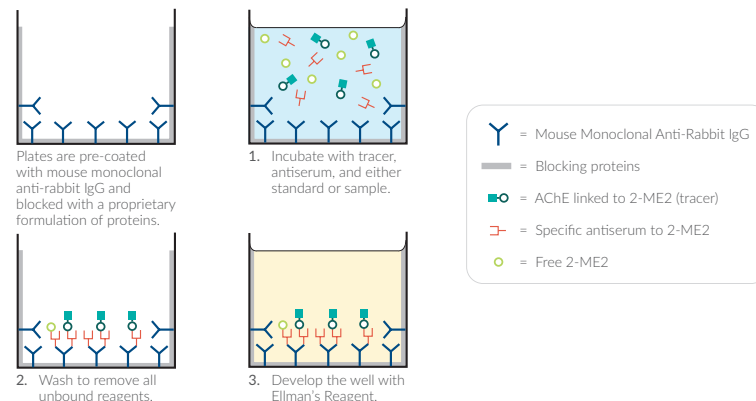


Figure 1. Schematic of the AChE ELISA

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.

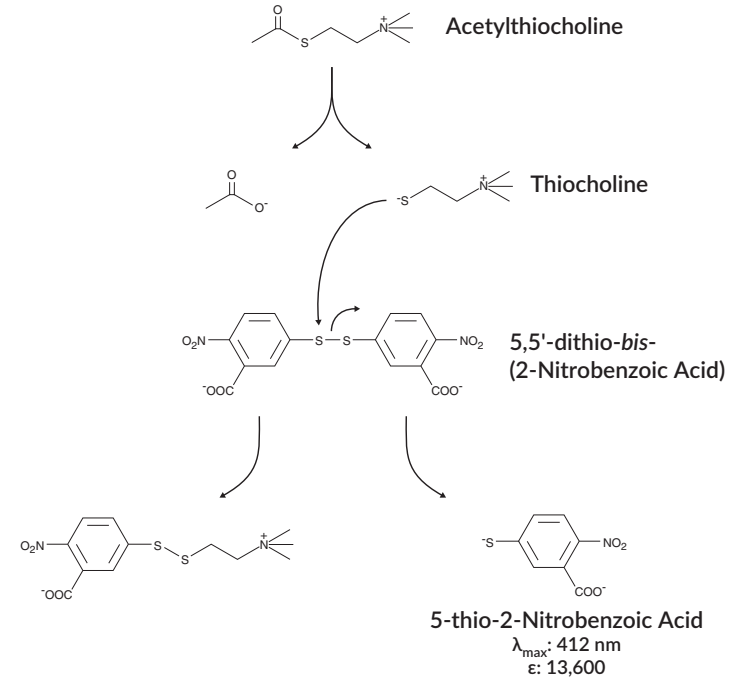


Figure 2. Reaction catalyzed by acetylcholinesterase

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 Concentrate (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 Concentrate (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

2-ME2 is present in plasma and urine mostly as glucuronide and sulfate conjugates. Treatment of samples with glucuronidase and sulfatase in our lab indicates that the antiserum in this kit reacts with only the unconjugated form of 2-ME2 (see Figure 3). Therefore, it is recommended that plasma and urine samples be hydrolyzed with glucuronidase and sulfatase prior to assay. *NOTE: b-glucuronidase from some sources also has sulfatase activity. If hydrolysis is performed with such a b-glucuronidase, it may not be necessary to add additional sulfatase.* The Hydrolysis protocol, on page 14, is one method that may be used for hydrolysis of either plasma or urine.

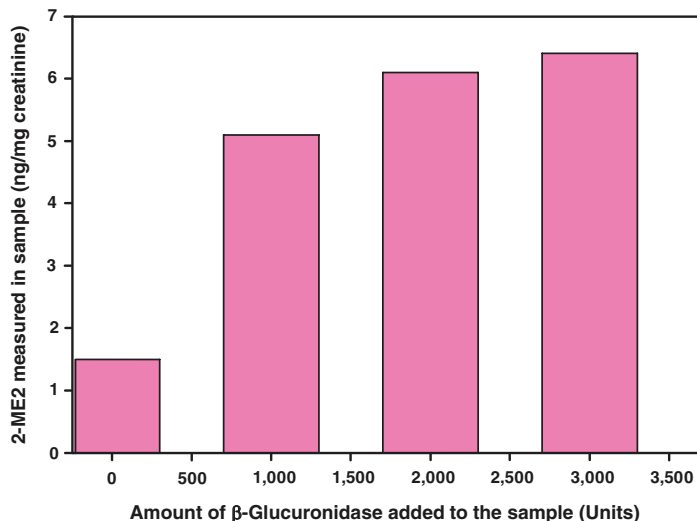


Figure 3. Treatment with glucuronidase increases measurable 2-ME2 in urine. Urine samples were incubated with increasing amounts of β -glucuronidase overnight at 37°C prior to analysis using the 2-Methoxyestradiol ELISA Kit.

Testing for Interference

Cayman strongly recommends that you test for interference before embarking on a large number of sample measurements. To test for interference, dilute one or two test samples to obtain at least two different dilutions of each sample between approximately 40-3,000 pg/ml (*i.e.*, between 20-80% B/B₀). If the two dilutions of the sample show good correlation (differ by 20% or less) in the final calculated 2-ME2 concentration, purification is not required. If you do not see good correlation of the different dilutions, purification is advised. The Extraction protocol found on page 15 may be used for this purpose.

General Precautions

- All samples must be free of organic solvents prior to assay.
- Samples should be assayed immediately after collection; samples that cannot be assayed immediately should be stored at -80°C.
- Samples of rabbit origin may contain antibodies which interfere with the assay by binding to the mouse anti-rabbit IgG-coated plate. We recommend that all rabbit samples be purified prior to use in this assay.

Plasma

Collect blood in tubes containing heparin or sodium citrate. To obtain plasma, spin samples at 1,000 x g for 15 minutes. Transfer the top yellow layer to a clean tube, being careful not to disturb the buffy layer. Samples should be assayed immediately after collection; samples that cannot be assayed immediately should be stored at -80°C.

Plasma or Urine Hydrolysis

1. Aliquot a known volume of each sample (500 µl is recommended) into a clean tube. If your samples need to be concentrated, a larger volume should be used (e.g., a 5 ml sample will be concentrated by a factor of 10, a 10 ml sample will be concentrated by a factor of 20, etc.)
2. Add an equal volume of 2 mg/ml β-glucuronidase from *Helix pomatia* (Sigma Item No. G0751) in 0.1 M acetate buffer, pH 5.0, vortex to mix, and incubate overnight at 37°C.

It may be possible to use hydrolyzed urine samples directly in the immunoassay without extraction. To determine if this is the case for your samples, test for interference following the recommendations on page 13. If extraction is required, the protocol described for plasma is one method that may be used (see page 13).

NOTE: Hydrolysis causes a two-fold dilution of samples. Make sure to account for this dilution in calculating the final concentration of 2-ME2 in your samples if you do not extract them.

Plasma Extraction

1. Add 3-6 volumes of ethanol to each sample, vortex to mix, and incubate at room temperature for 10 minutes.
2. Centrifuge samples at 1,250 x g for 10 minutes. Transfer supernatants into clean test tubes.
3. Evaporate ethanol by heating to 30°C under a gentle stream of nitrogen.
4. Reconstitute extracts in a volume of UltraPure water equal to the original sample volume.
5. Add 4 volumes of methylene chloride and mix thoroughly by vortexing. Allow the layers to separate. Transfer the methylene chloride (lower) layer to a clean tube using a transfer pipette. Repeat this extraction procedure three times.*
6. Evaporate the methylene chloride by heating to 30°C under a gentle stream of nitrogen.
7. Dissolve the extract in 500 µl of diluted ELISA Buffer.

NOTE: Purified plasma samples should be diluted an additional two-fold (e.g., 500 µl reconstituted sample + 500 µl of ELISA Buffer) to minimize interference.

*If it is necessary to stop during this purification, samples may be stored in the methylene chloride solution at -20°C or -80°C.

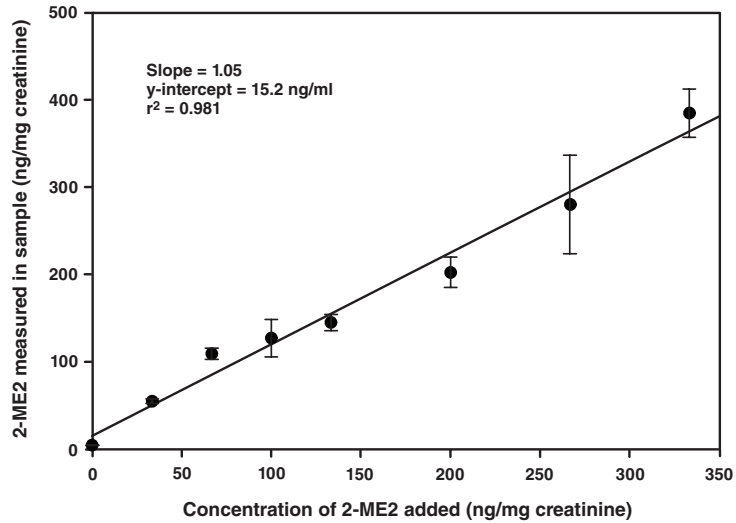


Figure 4. Recovery of 2-ME2 from urine

Urine samples were spiked with 2-ME2, hydrolyzed as described in the Sample Preparation section, and analyzed using the 2-Methoxyestradiol ELISA Kit. The y-intercept corresponds to the amount of 2-ME2 in unspiked urine. Error bars represent standard deviations obtained from multiple dilutions of each sample.

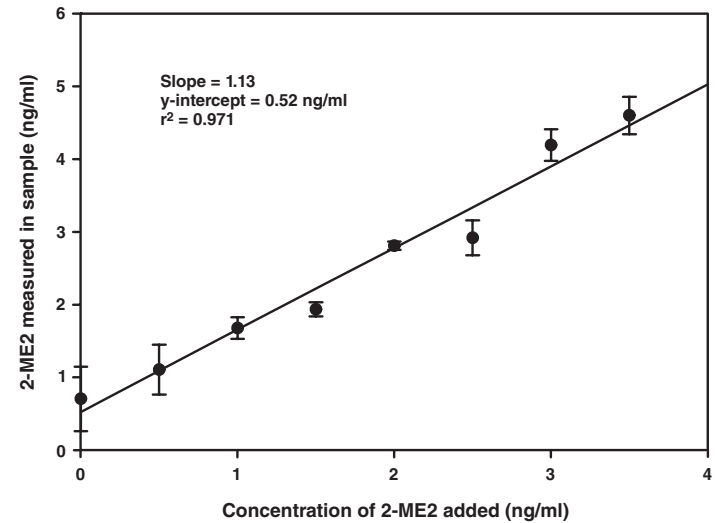


Figure 5. Recovery of 2-ME2 from plasma

Plasma samples were spiked with 2-ME2, hydrolyzed as described in the Sample Preparation section, and analyzed using the 2-Methoxyestradiol ELISA Kit. The y-intercept corresponds to the amount of 2-ME2 in unspiked plasma. Error bars represent standard deviations obtained from multiple dilutions of each sample.

| Level | %CV Intra-assay variation | Average (pg/ml) | %CV Inter-assay variation | Average (pg/ml) |
|--------|---------------------------|-----------------|---------------------------|-----------------|
| High | 11.49 | 3,037 | 7.10 | 3,080 |
| Medium | 16.56 | 1,832 | 9.65 | 1,921 |
| Low | 22.00 | 347 | 10.14 | 323 |

Table 1. Plasma sample validation

Plasma samples containing a high, medium, or low level of 2-ME2 were measured 60 times each using a single set of reagents. The calculated coefficient of variance (%CV) is reported as intra-assay variance. The same plasma samples containing a high, medium, or low level of 2-ME2 were measured four times each using eight independent sets of reagents. The calculated CV is reported as inter-assay variance.

ASSAY PROTOCOL

Preparation of Assay-Specific Reagents

2-Methoxyestradiol 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 2-Methoxyestradiol ELISA Standard (Item No. 482264) into a clean test tube, then dilute with 900 μ l UltraPure water. The concentration of this solution (the bulk standard) will be 80 ng/ml. Do not store this solution for more than 24 hours.

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 600 μ l ELISA Buffer to tubes #2-8. Transfer 100 μ l of the bulk standard (80 ng/ml) to tube #1 and mix thoroughly. Serially dilute the standard by removing 400 μ l from tube #1 and placing in tube #2; mix thoroughly. Next, remove 400 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.

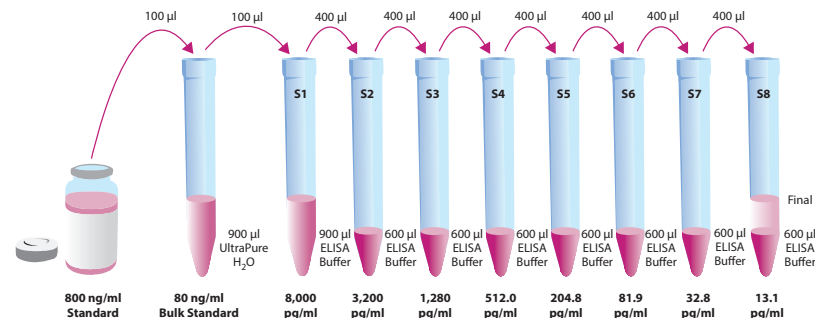


Figure 6. Preparation of the 2-ME2 standards

2-Methoxyestradiol AChE Tracer

Reconstitute the 2-Methoxyestradiol AChE Tracer as follows:

100 dtn 2-Methoxyestradiol AChE Tracer (96-well kit; Item No. 482260):
Reconstitute with 6 ml ELISA Buffer.

OR

500 dtn 2-Methoxyestradiol AChE Tracer (480-well kit; Item No. 482260):
Reconstitute with 30 ml ELISA Buffer.

Store the reconstituted 2-Methoxyestradiol AChE Tracer at 4°C (*do not freeze!*) and use within five 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). *NOTE: Do not store tracer with dye for more than 24 hours.*

2-Methoxyestradiol ELISA Antiserum

Reconstitute the 2-Methoxyestradiol ELISA Antiserum as follows:

100 dtn 2-Methoxyestradiol ELISA Antiserum (96-well kit; Item No. 482262): Reconstitute with 6 ml ELISA Buffer.

OR

500 dtn 2-Methoxyestradiol ELISA Antiserum (480-well kit; Item No. 482262): Reconstitute with 30 ml ELISA Buffer.

Store the reconstituted 2-Methoxyestradiol ELISA Antiserum at 4°C and use within five weeks. A 20% surplus of antiserum has been included to account for any incidental losses.

Antiserum Dye Instructions (optional)

This dye may be added to the antiserum, if desired, to aid in visualization of antiserum-containing wells. Add the dye to the reconstituted antiserum at a final dilution of 1:100 (add 60 µl of dye to 6 ml antiserum or add 300 µl of dye to 30 ml of antiserum). *NOTE: Do not store antiserum with dye for more than 24 hours.*

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 (B₀), 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 below in Figure 7. 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 **Analysis**, page 26, 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
 B₀ - 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 100 µl ELISA Buffer to NSB wells. Add 50 µl ELISA Buffer to B₀ wells. If culture medium was used to dilute the standard curve, substitute 50 µl of culture medium for ELISA Buffer in the NSB and B₀ wells (i.e., add 50 µl culture medium to NSB and B₀ wells and 50 µl ELISA Buffer to NSB wells).

2. 2-Methoxyestradiol ELISA Standard

Add 50 µl from tube #8 to both of the lowest standard wells (S8). Add 50 µl 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 50 µl 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. 2-Methoxyestradiol AChE Tracer

Add 50 µl to each well *except* the TA and the Blk wells.

5. 2-Methoxyestradiol ELISA Antiserum

Add 50 µl to each well *except* the TA, the NSB, and the Blk wells.

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

Table 2. Pipetting summary

Incubation of the Plate

Cover each plate with plastic film (Item No. 400012) and incubate overnight 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 be 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.*, B₀ 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 B₀ wells have reached a minimum of 0.3 A.U. (blank subtracted). The plate should be read when the absorbance of the B₀ wells in the range of 0.3-1.0 A.U. (blank subtracted). If the absorbance of the wells exceeds 2.0, wash the plate, add fresh Ellman's Reagent and let it develop again.

ANALYSIS

Many plate readers come with data reduction software that plots data automatically. Alternatively a spreadsheet program can be used. The data should be plotted as either %B/B₀ versus log concentration using a four-parameter logistic fit or as logit B/B₀ versus log concentration using a linear fit. *NOTE: Cayman Chemical 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 B₀ wells.
3. Subtract the NSB average from the B₀ average. This is the corrected B₀ or corrected maximum binding.
4. Calculate the B/B₀ (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 B₀ (from Step 3). Repeat for S2-S8 and all sample wells. (To obtain %B/B₀ 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 B₀ 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 28). 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 2-ME2 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. *NOTE: Remember to account for any concentration or dilution of the sample prior to the addition to the well. Samples that were hydrolyzed as described on page 14, but were not extracted, will be diluted two-fold from their original concentration. 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

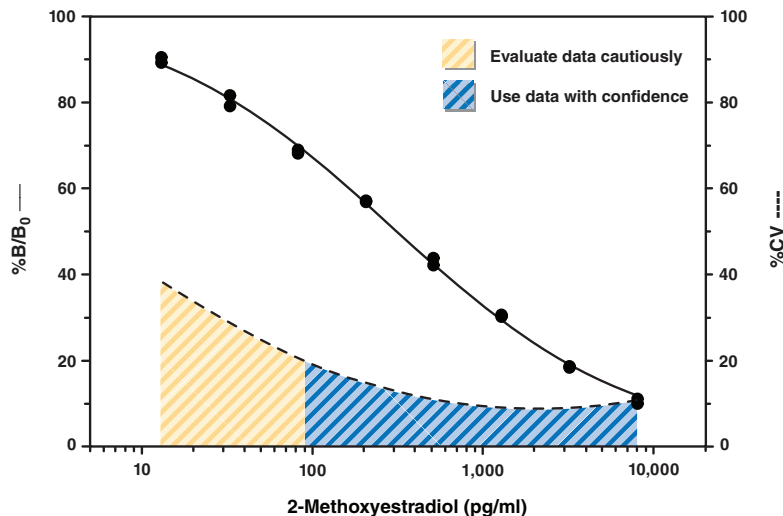
Sample Data

The standard curve presented here is an example of the data typically obtained 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 value of your samples. Your results could differ substantially.

| | Raw Data | | Average | Corrected |
|----------------|----------|-------|---------|-----------|
| Total Activity | 3.029 | 3.101 | 3.065 | |
| NSB | 0.003 | 0.003 | 0.003 | |
| B ₀ | 0.957 | 0.968 | | |
| | 0.965 | 1.030 | 0.980 | 0.977 |

| Dose (pg/ml) | Raw Data | | Corrected | | %B/B ₀ | |
|--------------|----------|-------|-----------|-------|-------------------|------|
| | | | | | | |
| 8,000 | 0.101 | 0.110 | 0.098 | 0.107 | 10.0 | 10.9 |
| 3,200 | 0.183 | 0.185 | 0.180 | 0.182 | 18.4 | 18.6 |
| 1,280 | 0.299 | 0.303 | 0.296 | 0.300 | 30.2 | 30.6 |
| 512.0 | 0.431 | 0.416 | 0.428 | 0.413 | 43.7 | 42.1 |
| 204.8 | 0.560 | 0.562 | 0.557 | 0.559 | 56.8 | 57.0 |
| 81.9 | 0.679 | 0.670 | 0.676 | 0.667 | 69.0 | 68.1 |
| 32.8 | 0.802 | 0.778 | 0.799 | 0.775 | 81.5 | 79.1 |
| 13.1 | 0.878 | 0.888 | 0.875 | 0.885 | 89.3 | 90.3 |

Table 3. Typical results



Assay Range = 13.1-8,000 pg/ml
Sensitivity (defined as 80% B/B₀) = 40 pg/ml
Mid-point (defined as 50% B/B₀) = 330-390 pg/ml

The sensitivity and mid-point were derived from the standard curve shown above. The standard was diluted with ELISA Buffer.

Figure 8. Typical standard curve

Precision:

The intra- and inter-assay CVs have been determined at multiple points on the standard curve. These data are summarized in the graph on page 29 and in the table below.

| Dose (pg/ml) | %CV* Intra-assay variation | %CV* Inter-assay variation |
|--------------|-------------------------------|-------------------------------|
| 8,000 | 11.3 | 18.6 |
| 3,200 | 7.0 | 9.8 |
| 1,280 | 11.1 | 11.3 |
| 512.0 | 11.8 | 11.3 |
| 204.8 | 13.4 | 9.8 |
| 81.9 | 18.6 | 13.2 |
| 32.8 | † | † |
| 13.1 | † | † |

Table 3. Intra- and inter-assay variation

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

†Outside of the recommended usable range of the assay.

Cross Reactivity:

| Compound | Cross Reactivity |
|------------------------|------------------|
| 2-Methoxyestradiol | 100% |
| Estradiol | 0.29% |
| Estrone | 0.04% |
| Progesterone | 0.04% |
| Testosterone | 0.02% |
| Androstenolone | 0.01% |
| Estriol | 0.01% |
| Cholesterol | <0.01% |
| 17 α -Estradiol | <0.01% |
| Pregnenolone | <0.01% |

Table 4. Cross Reactivity of the 2-Methoxyestradiol 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. Re-wash 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 ¹³ |
| 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. D'Amato, R.J., Lin, C.M., Flynn, E., et al. *Proc. Nat Acad. Sci. USA* **91**, 3964-3968 (1994).
2. Sweeney, C., Liu, G., Yiannoutsos, C., et al. *Clin. Cancer Res.* **11(18)**, 6625-6633 (2005).
3. Roswall, P., Bu, S., Rubin, K., et al. *Thyroid* **16(2)**, 143-150 (2006).
4. Mueck, A.O. and Seeger, H. *Steroids* **75**, 625-631 (2010).
5. Bourghardt, J., Bergström, G., Krettek, A., et al. *Endocrinology* **148(9)**, 4128-4132 (2007).
6. Kanasaki, K., Palmsten, K., Sugimoto, H., et al. *Nature* **453**, 1117-1123 (2008).
7. LaVallee, T.M., Zhan, X.H., Herbstritt, C.J., et al. *Cancer Res.* **62**, 3691-3697 (2002).
8. Xu, X., Duncan, A.M., Merz-Demlow, B.E., et al. *J. Clin. Endocrin. Metab.* **84(11)**, 3914-3918 (1999).
9. Xu, X., Veenstra, T.D., Fox, S.D., et al. *Anal. Chem.* **77**, 6646-6654 (2005).
10. Berg, F.D. and Kuss, E. *Arch. Gynecol. Obstet.* **251**, 17-27 (1992).
11. Maclouf, J., Grassi, J., and Pradelles, P. Chapter 5, in *Prostaglandin and Lipid Metabolism in Radiation Injury*. Walden, T.L., Jr. and Hughes, H.N., editors, Plenum Press, Rockville, 355-364 (1987).
12. Pradelles, P., Grassi, J. and Maclouf, J. *Anal. Chem.* **57**, 1170-1173 (1985).
13. Maxey, K.M., Maddipati, K.R., and Birkmeier, J. *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.

©08/18/2016, Cayman Chemical Company, Ann Arbor, MI, All rights reserved.
Printed in U.S.A.

