



Cayman Practice ELISA Kit

Item No. 10009658

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	About This Assay
	6	Principle of This Assay
	8	Biochemistry of AChE
	10	Definition of Key Terms
PRE-ASSAY PREPARATION	13	Buffer Preparation
ASSAY PROTOCOL	14	Preparation of Assay-Specific Reagents
	16	Plate Set Up
	17	Performing the Assay
ANALYSIS	20	Calculations
	22	Performance Characteristics
RESOURCES	24	Troubleshooting
	25	Plate Template
	26	Notes
	27	Warranty and Limitation of Remedy

GENERAL INFORMATION

Materials Supplied

Number	Item	96 wells Quantity/Size
400814	Practice ELISA Antiserum	1 vial/100 dtn
400815	Practice ELISA AChE Tracer	1 vial/100 dtn
400816	Practice ELISA Standard	1 vial
400060	ELISA Buffer Concentrate (10X)	1 vial/10 ml
400062	Wash Buffer Concentrate (400X)	1 vial/5 ml
400035	Polysorbate 20	1 vial/3 ml
400004	Mouse Anti-Rabbit IgG-Coated Plate	1 plate
400012	96-Well Cover Sheet	1 ea
400050	Ellman's Reagent	4 vials/100 dtn
400040	ELISA Tracer Dye	1 vial
400042	ELISA Antiserum Dye	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 Cayman Practice ELISA Kit. This kit may not perform as described if any reagent or procedure is replaced or modified.

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, with a resistivity of 18.2 MΩ·cm and total organic carbon (TOC) levels of <10 ppb, is recommended. Pure water - glass-distilled or deionized - may not be acceptable. *NOTE: UltraPure Water is available for purchase from Cayman (Item No. 400000).*
4. An orbital microplate shaker

INTRODUCTION

About This Assay

This assay has been developed for researchers that do not yet have experience performing enzyme-linked immunosorbent assays (ELISAs). It can be used as a practice tool allowing the user to become comfortable with running Cayman ELISAs. Practicing these assays can help decrease errors when samples, time, and costs are at risk.

This kit contains enough reagents to run at least four complete standard curves, including blank wells, non-specific binding wells, and maximum binding wells. For more information regarding the development and science behind the manufacturing of ELISAs please contact our technical service department.

Principle of This Assay

Competitive enzyme immunoassays are based on the competition between an unlabeled analyte and an enzyme-labeled analyte. In this particular assay, the enzyme labeled to the analyte is acetylcholinesterase (AChE tracer). These analytes compete for a limited number of analyte-specific rabbit antiserum binding sites. Because the concentration of the tracer is held constant while the concentration of free analyte varies, the amount of tracer that is able to bind to the rabbit antiserum will be inversely proportional to the concentration of free analyte in the well. This rabbit antiserum-analyte (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 for 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 tracer bound to the well, which is inversely proportional to the amount of free analyte present in the well during the incubation, as described in the equation:

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

A schematic of this process is shown in Figure 1 on page 7.

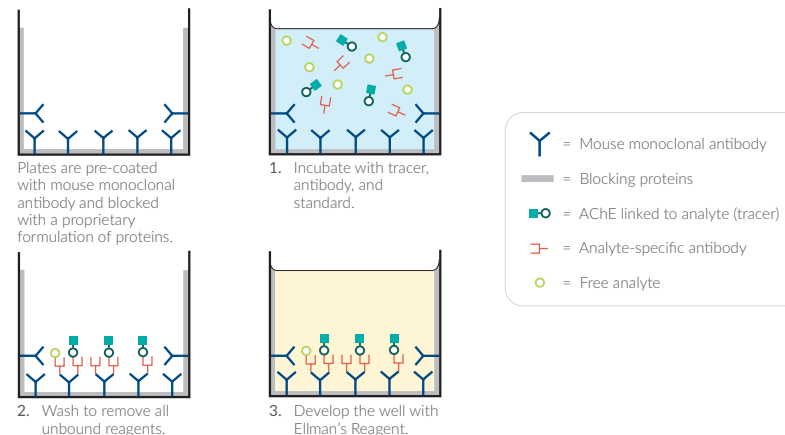


Figure 1. Schematic of the ELISA

Biochemistry of AChE

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.

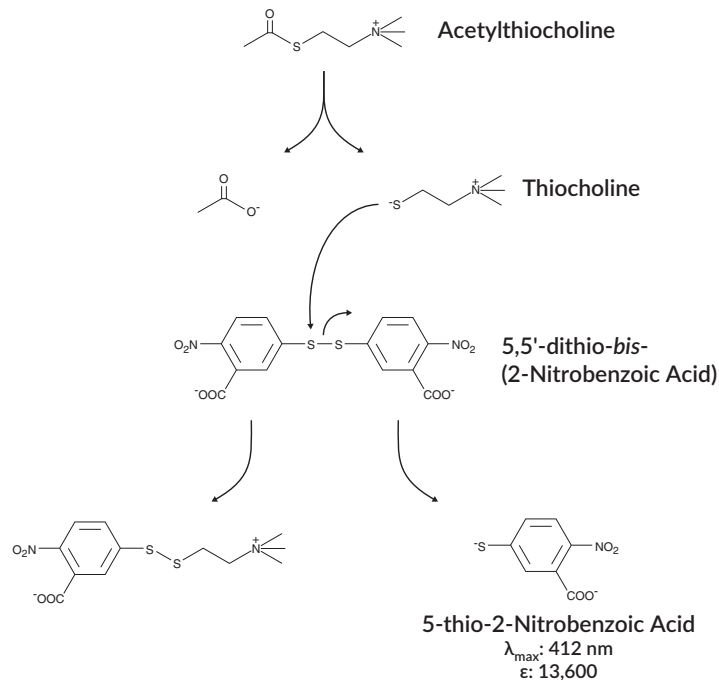


Figure 2. Reaction catalyzed by AChE

Definition of Key Terms

Blk (Blank): background absorbance caused by Ellman's Reagent. The Blk absorbance should be subtracted from the absorbance readings of all the other wells, including the non-specific binding (NSB) wells.

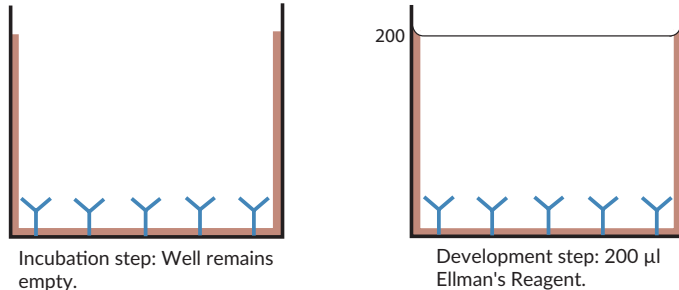


Figure 3. Blank well

TA (Total Activity): total enzymatic activity of the AChE-linked tracer.

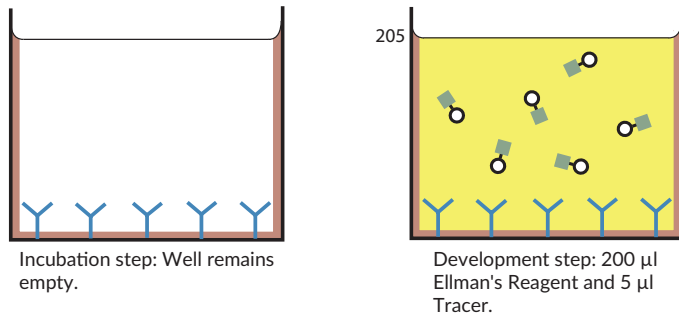


Figure 4. Total activity well

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 Blk absorbance values.

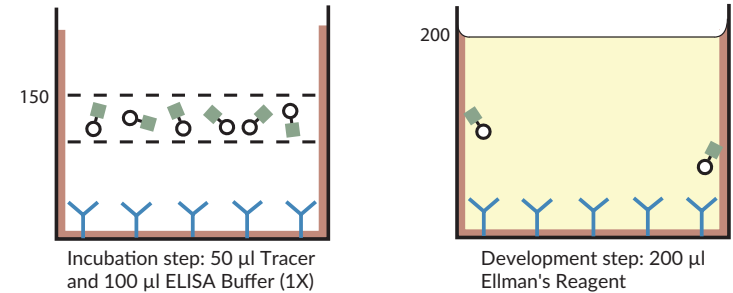


Figure 5. Non-specific binding well

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

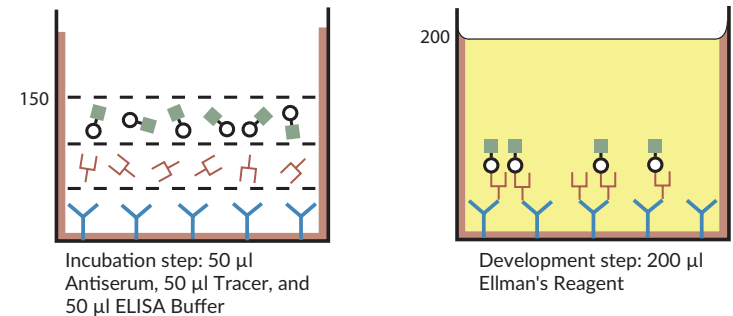


Figure 6. Maximum binding (B₀) well

Standard Curve: a plot of the $\%B/B_0$ values versus concentration of a series of wells containing various known amounts of analyte.

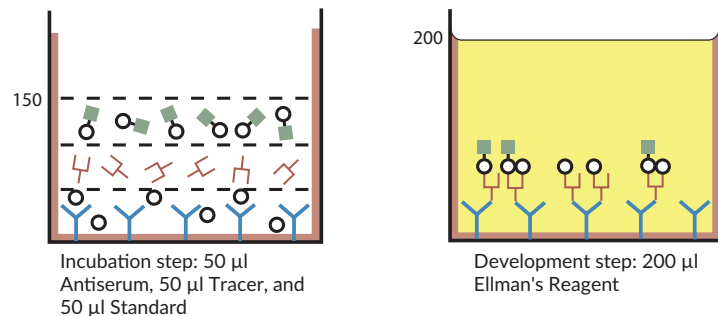


Figure 7. Standard curve well

PRE-ASSAY PREPARATION

Buffer Preparation

Store all diluted buffers at 4°C; they will be stable for approximately two months. *NOTE: It is normal for the concentrated buffers to contain crystalline salts. These will completely dissolve upon dilution with ultrapure water.*

1. ELISA Buffer (1X) 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.

2. Wash Buffer (1X) Preparation

Dilute the contents of the vial of Wash Buffer Concentrate (400X) (Item No. 400062) to a total volume of 2 L with ultrapure water and add 1 ml of Polysorbate 20 (Item No. 400035). Smaller volumes of Wash Buffer (1X) can be prepared by diluting the Wash Buffer Concentrate (400X) 1:400 and adding 0.5 ml of Polysorbate 20 per 1 L of Wash Buffer (1X). *NOTE: Polysorbate 20 is a viscous liquid and cannot be measured by a regular pipette. A positive displacement pipette or syringe should be used to deliver small quantities accurately.*

ASSAY PROTOCOL

Preparation of Assay-Specific Reagents

Practice 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 Practice ELISA Standard (Item No. 400816) into a clean test tube, then dilute with 900 μ l ultrapure water. The concentration of this solution (the bulk standard) will be 5 ng/ml.

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 (1X) to tube #1 and 500 μ l ELISA Buffer (1X) to tubes #2-8. Transfer 100 μ l of the bulk standard (5 ng/ml) to tube #1 and mix thoroughly. 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. The diluted standards may be stored at 4°C for no more than 24 hours.

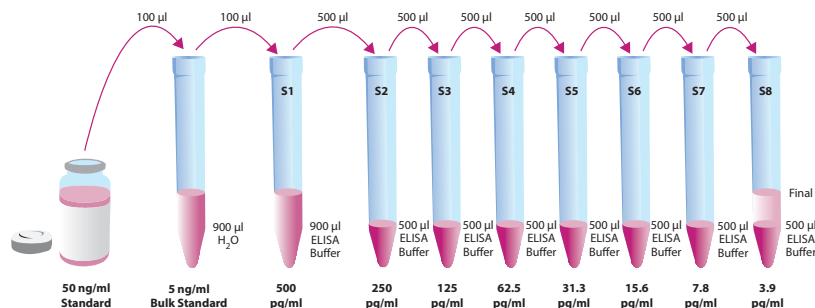


Figure 8. Preparation of the ELISA standards

Practice ELISA AChE Tracer

Reconstitute the Practice ELISA AChE Tracer (Item No. 400815) with 6 ml ELISA Buffer (1X). Store the reconstituted Practice ELISA AChE Tracer at 4°C (*do not freeze!*) and use within four 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 μ l of dye to 6 ml tracer). *NOTE: Do not store tracer with dye for more than 24 hours.*

Practice ELISA Antiserum

Reconstitute the Practice ELISA Antiserum (Item No. 400814) with 6 ml ELISA Buffer (1X). Store the reconstituted Practice ELISA Antiserum at 4°C. It will be stable for at least four 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 μ l of dye to 6 ml 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. *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 set of strips must contain a minimum of two Blk, two NSB, and three B₀ wells, and an eight-point standard curve run in duplicate. TA wells are recommended as a diagnostic tool. Low or no absorbance from a TA well could indicate a dysfunction in the enzyme-substrate system.

A suggested plate format is shown in Figure 9, below. The plate format provided below has been designed to allow for easy data analysis using a convenient spreadsheet offered by Cayman (see page 20, for more details). We suggest you record the contents of each well on the template sheet provided (see page 25).

	1	2	3	4	5	6	7	8	9	10	11	12
A	Blk	S1	S1	Blk	S1	S1	Blk	S1	S1	Blk	S1	S1
B	Blk	S2	S2	Blk	S2	S2	Blk	S2	S2	Blk	S2	S2
C	NSB	S3	S3	NSB	S3	S3	NSB	S3	S3	NSB	S3	S3
D	NSB	S4	S4	NSB	S4	S4	NSB	S4	S4	NSB	S4	S4
E	B ₀	S5	S5	B ₀	S5	S5	B ₀	S5	S5	B ₀	S5	S5
F	B ₀	S6	S6	B ₀	S6	S6	B ₀	S6	S6	B ₀	S6	S6
G	B ₀	S7	S7	B ₀	S7	S7	B ₀	S7	S7	B ₀	S7	S7
H	TA	S8	S8	TA	S8	S8	TA	S8	S8	TA	S8	S8

Blk - Blank
 TA - Total Activity
 NSB - Non-Specific Binding
 B₀ - Maximum Binding
 S1-S8 - Standards 1-8

Figure 9. 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 (1X)

Add 100 µl ELISA Buffer (1X) to NSB wells. Add 50 µl ELISA Buffer (1X) to B₀ wells.

2. Practice ELISA Standard

Add 50 µl from tube #8 to the lowest standard wells (S8). Add 50 µl from tube #7 to each of the next 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. Practice ELISA AChE Tracer

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

4. Practice ELISA Antiserum

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

Well	ELISA Buffer (1X)	Practice ELISA Standard	Practice ELISA AChE Tracer	Practice ELISA Antiserum
Blk	-	-	-	-
TA	-	-	5 µl (at devel. step)	-
NSB	100 µl	-	50 µl	-
B ₀	50 µl	-	50 µl	50 µl
Standard	-	50 µl	50 µl	50 µl

Table 1. Pipetting summary

Incubation of the Plate

Cover each plate with a 96-Well Cover Sheet (Item No. 400012) and incubate for two hours at room temperature on an orbital shaker.

Development of the Plate

1. Reconstitute Ellman's Reagent (Item No. 400050) immediately before use with 20 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 ~300 µl Wash Buffer (1X).
3. Add 200 µl of Ellman's Reagent to each well.
4. Add 5 µl of the reconstituted tracer to the TA wells.
5. Cover the plate with the 96-Well Cover Sheet. 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. (Blk subtracted)) in 60-90 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 (1X) 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. (Blk subtracted). The plate should be read when the absorbance of the B₀ wells is in the range of 0.3-1.0 A.U. (Blk 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 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 Blk 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 (Standard Bound/Maximum Bound) for the standard 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. (To obtain %B/B₀ for a logistic four-parameter fit, multiply these values by 100.)

$$\% B/B_0 = \left[\frac{(\text{Standard Absorbance} - \text{NSB})}{(\text{Average } B_0 \text{ Absorbance}) - \text{NSB}} \right] \times 100$$

Plot the Standard Curve

Plot %B/B₀ for standards S1-S8 versus analyte concentration using linear (y) and log (x) axes and perform a four-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.*

$$\text{logit } (B/B_0) = \ln [B/B_0 / (1 - B/B_0)]$$

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

Performance Characteristics

Sample Data

Analyte (pg/ml)	Blk-Subtracted Absorbance	NSB-Corrected Absorbance	%B/B ₀
NSB	0.000	--	--
B0	0.743	0.743	--
500	0.136	0.136	18.2
250	0.200	0.200	26.9
125	0.284	0.284	38.2
62.5	0.381	0.381	51.2
31.3	0.485	0.485	65.2
15.6	0.571	0.571	76.8
7.8	0.649	0.649	87.4
3.9	0.685	0.685	92.2
TA	0.483	--	--

Table 2. Typical results

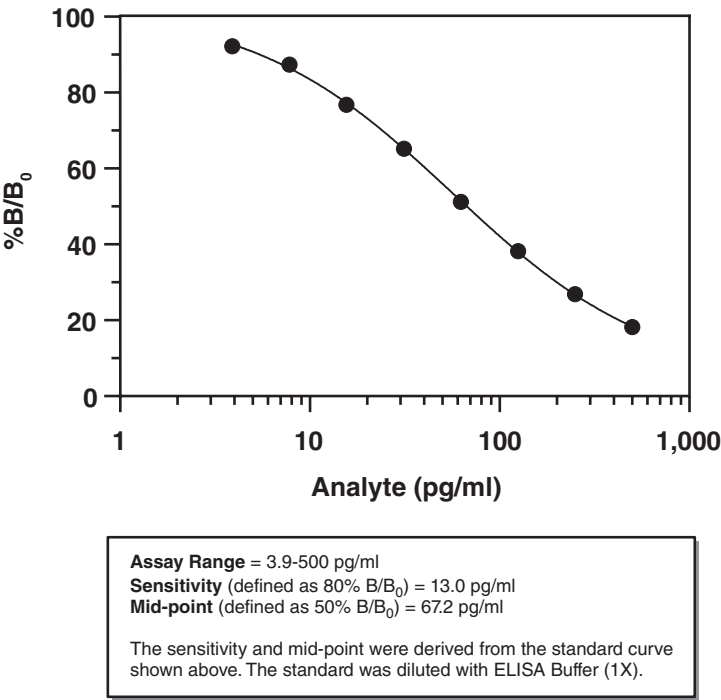


Figure 10. Typical standard curve

Troubleshooting

Problem	Possible Causes	Recommended Solutions
Erratic values; dispersion of replicates	A. Trace organic contaminants in the water 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 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
Only TA wells develop	Trace organic contaminants in the water	Replace activated carbon filter or change source of ultrapure water

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

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.

©01/28/2025, Cayman Chemical Company, Ann Arbor, MI, All rights reserved.
Printed in U.S.A.

