

Version 7c Last updated 21 April 2026

ab157718

Mouse IgE ELISA Kit

For the quantitative measurement of IgE in mouse serum and plasma samples.

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	5
9. Reagent Preparation	6
10. Standard Preparation	8
11. Sample Preparation	10
12. Plate Preparation	11
13. Assay Procedure	12
14. Calculations	13
15. Typical Data	14
16. Typical Sample Values	15
17. Assay Specificity	15
18. Troubleshooting	16
19. Notes	17

1. Overview

Abcam's IgE mouse ELISA kit is an in vitro enzyme-linked immunosorbent assay (ELISA) for the quantitative measurement of IgE in biological samples of mice.

In this assay the IgE present in samples reacts with the anti-IgE antibodies which have been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti-IgE antibodies conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound IgE. Following another washing step, the amount of enzyme bound in complex is measured by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies proportionately with the concentration of IgE in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of IgE in the test sample. The quantity of IgE in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

2. Protocol Summary

Remove appropriate number of antibody coated well strips.



Equilibrate all reagents to room temperature. Prepare all the reagents, samples, and standards as instructed.



Add standard or sample to each well used. Incubate at room temperature.



Aspirate and wash each well. Add prepared HRP labeled secondary detector antibody. Incubate at room temperature



Aspirate and wash each well. Add Chromogen Substrate Solution to each well.



Immediately begin recording the color development.

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- If applicable, please refer to the current Safety Data Sheet (SDS) provided with this product for safety, handling, and disposal information. The most up to date and current versions are available on our website <https://www.abcam.com/en-us>.

4. Storage and Stability

Store kit at +4°C immediately upon receipt.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

Item	Amount	Storage Condition (Before Preparation)
Mouse IgE ELISA Microplate	96 wells	4°C
Mouse IgE Calibrator (lyophilized)	1 vial	4°C
5X Diluent Concentrate	50 mL	4°C
20X Wash Buffer Concentrate	50 mL	4°C
Enzyme-Antibody Conjugate	150 µL	4°C
Chromogen Substrate Solution	12 mL	4°C
Stop Solution	12 mL	4°C

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Precision pipette (2 μ L to 200 μ L) for making and dispensing dilutions
- Test tubes
- Microtitre washer/aspirator
- Distilled or Deionized H₂O
- Microtitre Plate reader
- Assorted glassware for the preparation of reagents and buffer solutions
- Timer

8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers
- Avoid foaming or bubbles when mixing or reconstituting components
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions
- Ensure plates are properly sealed or covered during incubation steps
- Complete removal of all solutions and buffers during wash steps

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells.
- Prepare only as much reagent as is needed on the day of the experiment.

9.1 1X Diluent Solution

The diluent solution is supplied as 5X Diluent Concentrate and must be diluted 1/5 with distilled or deionized water (1 part buffer concentrate, 4 parts dH₂O). The 1X Diluent Solution is stable for at least one week from the date of preparation and should be stored at 4°C.

9.2 1X Wash Buffer

The wash solution is supplied as 20X Concentrate and must be diluted 1/20 with distilled or deionized water (1 part buffer concentrate, 19 parts dH₂O). Crystal formation in the concentrate is not uncommon when storage temperatures are low. Warming of the concentrate to 30-35°C before dilution can dissolve crystals. The 1X Wash Buffer is stable for at least one week from the date of preparation and can be stored at room temperature (16-25°C) or at 4°C.

9.3 **1X Enzyme-Antibody Conjugate**

Calculate the required amount of 1X Enzyme-Antibody Conjugate solution for each microtitre plate test strip by adding 10 μL Enzyme-Antibody Conjugate to 990 μL of 1X Diluent for each test strip to be used for testing. Mix uniformly, but gently. Avoid foaming. The working conjugate solution is stable for up to 1 hour when stored in the dark.

9.4 **Chromogen Substrate Solution**

Ready to use as supplied.

9.5 **Stop Solution**

Ready to use as supplied.

9.6 **Mouse IgE Calibrator**

Add 1.0 mL of distilled or de-ionized water to the Mouse IgE Calibrator and mix gently until dissolved. The amount of calibrator is shown on the vial, and after reconstitution will have a concentration of $X \mu\text{g}/\text{mL}$, where X is the amount on the vial (the reconstituted calibrator should be aliquoted and frozen if future use is intended). Mouse IgE standards need to be prepared immediately prior to use. Mix well between each step. Avoid foaming.

10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

- 10.1 The liquid mouse IgE Calibrator should be aliquoted and stored frozen. Avoid multiple freeze-thaw cycles. The calibrator is provided at the concentration stated on the vial.
- 10.2 Label tubes numbers 1-7.
- 10.3 Prepare **Standard #1** by adding the appropriate volume of 1X Diluent Solution (derived below) to tube #1. Add 300 μ L of stock mouse IgE Calibrator to obtain a concentration at 400 ng/mL.

*Example:

NOTE: This example is for demonstration purposes only. Please remember to check your calibrator vial for the actual concentration of calibrator provided.

C_S = Starting concentration of reconstituted mouse IgE Calibrator
(variable e.g. 693.3 ng/mL)

C_F = Final concentration of mouse IgE Calibrator for the assay
procedure (400 ng/mL)

V_A = Total volume of stock mouse IgE Calibrator to dilute (e.g. 300 μ L)

V_D = Total volume of 1X Diluent Solution required to dilute stock mouse
IgE Calibrator to prepare **Standard #1**

V_T = Total volume of **Standard #1**

D_F = Dilution factor

Calculate the dilution factor (D_F) between stock calibrator and the **Standard #1** final concentration:

$$C_S/C_F = D_F$$
$$693.3 / 400 = 1.733$$

Calculate the final volume V_D required to prepare the **Standard #1** at 400 ng/mL

$$V_A * D_F = V_T$$

$$V_D = V_T - V_A$$

$$300 * 1.733 = 520 \mu\text{L}$$

$$V_D = 520 - 300 = 220 \mu\text{L}$$

To tube #1, add 300 μL of mouse IgE Calibrator to 220 μL of 1X Diluent Solution to obtain a concentration of 400 ng/mL (**Standard #1**).

- 10.4 Add 250 μL 1X Diluent Solution into tube numbers 2-7.
- 10.5 Prepare **Standard #2** by adding 250 μL **Standard #1** to tube #2. Mix thoroughly and gently.
- 10.6 Prepare **Standard #3** by adding 250 μL from **Standard #2 to #3**. Mix thoroughly and gently.
- 10.7 Using the table below as a guide to prepare further serial dilutions.
- 10.8 1X Diluent Solution serves as the zero standard (0 ng/mL).

Standard Dilution Preparation Table

Standard #	Volume to Dilute (μL)	Diluent (μL)	Total Volume (μL)	Starting Conc. (ng/mL)	Final Conc. (ng/mL)
1	See step 10.3				400
2	250	250	500	400	200
3	250	250	500	200	100
4	250	250	500	100	50
5	250	250	500	50	25
6	250	250	500	25	12.5
7	250	250	500	12.5	6.25

11. Sample Preparation

11.1 Serum:

Blood should be collected by venipuncture. The serum should be separated from the cells after clot formation by centrifugation.

11.2 Plasma:

For plasma samples, blood should be collected into a container with an anticoagulant and then centrifuged. Care should be taken to minimize hemolysis, excessive hemolysis can impact your results.

Assay immediately or aliquot and store samples at -20°C . Avoid repeated freeze-thaw cycles.

Precautions

For any sample that might contain pathogens, care must be taken to prevent contact with open wounds.

Additives and Preservatives

No additives or preservatives are necessary to maintain the integrity of the specimen. Avoid azide contamination.

General Sample information:

The assay for quantification of IgE in samples requires that each test sample be diluted before use. For a single step determination a dilution of 1/50 is appropriate for most serum/plasma samples. For absolute quantification, samples that yield results outside the range of the standard curve, a lesser or greater dilution might be required. If unsure of sample level, a serial dilution with one or two representative samples before running the entire plate is highly recommended.

To prepare a 1/50 dilution of sample, transfer 5 μL of sample to 245 μL of 1X diluent. This gives you a 1/50 dilution. Mix thoroughly.

Refer to Dilution Guidelines for further instruction.

Guidelines for Dilutions of 100-fold or Greater <i>(for reference only; please follow the insert for specific dilution suggested)</i>	
100x	10000x
<p>4 μl sample + 396 μl buffer (100X) = 100-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 400 μl</i></p>	<p>A) 4 μl sample + 396 μl buffer (100X) B) 4 μl of A + 396 μl buffer (100X) = 10000-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 400 μl</i></p>
1000x	100000x
<p>A) 4 μl sample + 396 μl buffer (100X) B) 24 μl of A + 216 μl buffer (10X) = 1000-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 240 μl</i></p>	<p>A) 4 μl sample + 396 μl buffer (100X) B) 4 μl of A + 396 μl buffer (100X) C) 24 μl of A + 216 μl buffer (10X) = 100000-fold dilution</p> <p><i>Assuming the needed volume is less than or equal to 240 μl</i></p>

12. Plate Preparation

- The 96 well plate strips included with this kit is supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused well plate strips should be returned to the plate packet and stored at 4°C.
- For each assay performed, a minimum of 2 wells must be used as blanks, omitting primary antibody from well additions
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

13. Assay Procedure

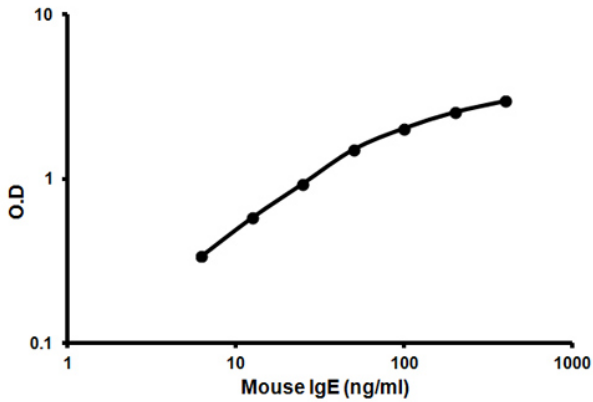
- Equilibrate all materials and prepared reagents to room temperature prior to use.
 - We recommend that you assay all standards, controls and samples in duplicate.
- 13.1 Pipette 100 μ L of each standard, including zero control, in duplicate, into pre designated wells.
 - 13.2 Pipette 100 μ L of sample (in duplicate) into pre designated wells.
 - 13.3 Incubate the microtiter plate at room temperature for thirty (30 ± 2) minutes. Keep plate covered and level during incubation.
 - 13.4 Following incubation, aspirate the contents of the wells.
 - 13.5 Completely fill each well with appropriately diluted 1X Wash Buffer and aspirate. Repeat three times, for a total of four washes. If washing manually: completely fill wells with wash buffer, invert the plate then pour/shake out the contents in a waste container. Follow this by gently striking the wells on absorbent paper to remove residual buffer. Repeat 3 times for a total of four washes.
 - 13.6 Pipette 100 μ L of appropriately 1X Enzyme-Antibody Conjugate to each well. Incubate at room temperature for thirty (30 ± 2) minutes. Keep plate covered in the dark and level during incubation.
 - 13.7 Wash and blot the wells as described in 13.4 - 13.5.
 - 13.8 Pipette 100 μ L of Chromogen Substrate Solution into each well.
 - 13.9 Incubate in the dark at room temperature for precisely ten (10) minutes.
 - 13.10 After ten minutes, add 100 μ L of Stop Solution to each well.
 - 13.11 Determine the absorbance (450 nm) of the contents of each well. Calibrate the plate reader to manufacturer's specifications.

14. Calculations

Average the duplicate standard reading for each standard, sample and control blank. Subtract the control blank from all mean readings. Plot the mean standard readings against their concentrations and draw the best smooth curve through these points to construct a standard curve. Most plate reader software or graphing software can plot these values and curve fit. A four-parameter algorithm (4PL) usually provides the best fit, though other equations can be examined to see which provides the most accurate (e.g. linear, semi-log, log/log, 4-parameter logistic). Extrapolate protein concentrations for unknown and control samples from the standard curve plotted. Samples producing signals greater than that of the highest standard should be further diluted in 1X Incubation Buffer and reanalyzed, then multiplying the concentration found by the appropriate dilution factor.

15. Typical Data

Typical standard curve – data provided **for demonstration purposes only**. A new standard curve must be generated for each assay performed.



Standard Curve Measurements	
Conc. (ng/mL)	Mean O.D.
6.25	0.341
12.5	0.583
25	0.939
50	1.513
100	2.036
200	2.551
400	2.981

Figure 1. Example of mouse IgE standard curve.

16. Typical Sample Values

SENSITIVITY –

Calculated minimum detectable dose = 1.83 ng/mL.

PRECISION –

	% CV
Inter-Assay	< 10%
Intra-Assay	< 10%

RECOVERY –

Control Serum Recovery = > 85%.

17. Assay Specificity

INTERFERENCES –

These chemicals or biologicals will cause interferences in this assay causing compromised results or complete failure.

Azide and thimerosal at concentrations higher than 0.1% inhibits the enzyme reaction.

Please contact our Technical Support team for more information.

18. Troubleshooting

Problem	Cause	Solution
Poor standard curve	Inaccurate pipetting	Check pipettes
	Improper standards dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times; change to overnight standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	Store the reconstituted protein at -80°C, all other assay components 4°C. Keep substrate solution protected from light.

19. Notes

Technical Support

Copyright © 2026 Abcam. All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

For all technical or commercial enquiries please go to:

<https://www.abcam.com/en-us/contact-us>

<https://www.abcam.cn/contact-us> (China)

<https://www.abcam.co.jp/contact-us> (Japan)