

## ab291069- Mouse C1q ELISA Kit

Immunoperoxidase Assay for Determination of C1q in Mouse Samples.  
This product is for research use only and is not intended for diagnostic use.

For overview, typical data and additional information please visit:

<http://www.abcam.com/ab291069>

### Storage and Stability

On receipt, the entire assay kit should be stored at 4°C, protected from light.

### Materials Supplied

Item	Quantity	Storage Condition
20X Wash Buffer Concentrate	50 mL	4°C
5X Diluent Concentrate	50 mL	4°C
Calibrator	1 vial	4°C
Chromogen Substrate Solution	12 mL	4°C
ELISA Micro Plate, antibody coated	1 unit	4°C
Enzyme Conjugated Detection Antibody	150 µL	4°C
Stop Solution	12 mL	4°C

### Materials Required, Not Supplied

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

- Precision pipettes (2 µL to 100 µL) for making and dispensing dilutions
- Test tubes
- Squirt bottle or Microtitre washer/aspirator
- Distilled or Deionized H<sub>2</sub>O
- Microtitre Plate reader
- Assorted glassware for the preparation of reagents and buffer solutions
- Centrifuge for sample collection
- Anticoagulant for plasma collection
- Timer
- Microplate shaker

### Reagent Preparation

- Bring all reagents to room temperature (16°C to 25°C) before use.

**Diluent Concentrate:** The Diluent Solution supplied is a 5X Concentrate and must be diluted 1:5 with distilled or deionized water (1 part buffer concentrate, 4 parts dH<sub>2</sub>O).

**Wash Solution Concentrate:** The Wash Solution supplied is a 20X Concentrate and must be diluted 1:20 with distilled or deionized water (1 part buffer concentrate, 19 parts dH<sub>2</sub>O). Crystal formation in the concentrate may occur when storage temperatures are low. Warming of the concentrate to 30-35°C before dilution can dissolve crystals.

**Enzyme-Antibody Conjugate:** Calculate the required amount of working 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. Dilute immediately before use and protect from light. Mix uniformly, but gently. Avoid foaming.

**Pre-coated ELISA Micro Plate:** Ready to use as supplied. Unseal foil pouch and remove plate from pouch. Remove all strips and wells that will not be used in the assay and place back in pouch and re-seal along with desiccant.

**C1q Calibrator:** Prepare according to the lot specific vial.

Version 1b, Last updated February 9, 2023

### Sample Preparation

#### Sample Collection and Handling

- All blood components and biological materials should be handled as potentially hazardous. Follow universal precautions when handling and disposing.
- If blood samples are clotted, grossly haemolyzed, lipemic, or the integrity of the sample is of concern, make a note and interpret results with caution.
- The sample collection and storage conditions listed below are intended as general guidelines. Sample stability has not been evaluated.

**Serum:** Blood should be collected by venipuncture. The serum should be separated from the cells after clot formation by centrifugation. Remove serum and assay immediately or aliquot and store samples at -80°C (preferably) or -20°C. Avoid repeated freeze-thaw cycles.

**Plasma:** Blood should be collected into a container with an anticoagulant and then centrifuged. Assay immediately or aliquot and store samples at -80°C (preferably) or -20°C. Avoid repeated freeze-thaw cycles.

**Urine:** Collect mid-stream using sterile or clean urine collector. Centrifuge to remove cell debris. Assay immediately or aliquot and store samples at -80°C (preferably) or -20°C. Avoid repeated freeze-thaw cycles.

**Known interfering substances:** Azide and thimerosal at concentrations higher than 0.1% inhibits the enzyme reaction.

#### Sample Dilution

- The assay requires that each test sample be diluted before use. All samples should be assayed in duplicate each time the assay is performed. The recommended dilutions are only suggestions. Dilutions should be based on the expected concentration of the unknown sample such that the diluted sample falls within the dynamic range of the standard curve. If unsure of sample level, a serial dilution with one or two representative samples before running the entire plate is highly recommended.

**Serum:** Recommended starting dilution is 1:1000. To prepare a 1:1000 dilution of a sample, transfer 5 µL of sample to 495 µL of 1X diluent. This gives you a 1:100 dilution. Next, dilute the 1:100 by transferring 30 µL into 270 µL of 1X diluent. This gives you a 1:1000 dilution. Mix thoroughly each stage.

**Plasma:** Recommended starting dilution is 1:1000. To prepare a 1:1000 dilution of a sample, transfer 5 µL of sample to 495 µL of 1X diluent. This gives you a 1:100 dilution. Next, dilute the 1:100 by transferring 30 µL into 270 µL of 1X diluent. This gives you a 1:1000 dilution. Mix thoroughly each stage.

## Assay Protocol

- All samples and standards should be assayed in duplicates.
1. The Standards and the test sample(s) should be loaded into the ELISA wells as quickly as possible to avoid a shift in OD readings. Using a multichannel pipette would reduce this occurrence.
  2. Pipette 100  $\mu$ L of:

Standard 0	(0.0 ng/mL) in duplicate
Standard 1	(1.57 ng/mL) in duplicate
Standard 2	(3.13 ng/mL) in duplicate
Standard 3	(6.25 ng/mL) in duplicate
Standard 4	(12.50 ng/mL) in duplicate
Standard 5	(25 ng/mL) in duplicate
Standard 6	(50 ng/mL) in duplicate
Standard 7	(100 ng/mL) in duplicate

3. Pipette 100  $\mu$ L of sample (in duplicate) into pre designated wells.
4. Incubate the microtiter plate at room temperature for 60 ( $\pm$  2) minutes. Keep plate covered and level during incubation.
5. Following incubation, aspirate the contents of the wells.
6. Completely fill each well with appropriately diluted Wash Solution 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 sharply striking the wells on absorbent paper to remove residual buffer. Repeat 3 times for a total of four washes.
7. Pipette 100  $\mu$ L of appropriately diluted Enzyme-Antibody Conjugate to each well. Incubate at room temperature for 30 ( $\pm$  2) minutes. Keep plate covered in the dark and level during incubation.
8. Wash and blot the wells as described in Steps 5/6.
9. Pipette 100  $\mu$ L of TMB Substrate Solution into each well.
10. Incubate in the dark at room temperature for precisely 10 minutes. Keep plate covered in the dark and level during incubation.
11. After 10 minutes, add 100  $\mu$ L of Stop Solution to each well.
12. Determine the absorbance (450 nm) of the contents of each well within 30 minutes. Calibrate the plate reader to manufacturer's specifications.

## Calculation

- Subtract the average background value (Average absorbance reading of Standard zero) from the test values for each sample.
- Average the duplicate readings for each standard and use the results to construct a Standard Curve. Construct the standard curve by reducing the data using computer software capable of generating a four parameter logistic curve fit. A second order polynomial (quadratic) or other curve fits may also be used; however, they will be a less precise fit of the data.
- Interpolate test sample values from standard curve. Correct for sera dilution factor to arrive at the C1q concentration in original samples.

## Technical Support

Copyright © 2023 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:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)