



# C3 (Dog) ELISA Kit

Catalog Number KA1923

96 assays

Version: 02

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Intended Use .....	3
Background .....	3
Principle of the Assay .....	3
<b>General Information .....</b>	<b>5</b>
Materials Supplied .....	5
Storage Instruction .....	5
Materials Required but Not Supplied .....	6
Precautions for Use .....	6
<b>Assay Protocol .....</b>	<b>7</b>
Reagent Preparation .....	7
Sample Preparation .....	8
Assay Procedure .....	8
<b>Data Analysis .....</b>	<b>10</b>
Calculation of Results .....	10
<b>Resources .....</b>	<b>11</b>
<b>Plate Layout .....</b>	<b>11</b>

## **Introduction**

### **Intended Use**

The C3 (Dog) ELISA Kit is a highly sensitive two-site enzyme linked immunoassay (ELISA) for measuring C3 in biological samples of dogs. For Research Use Only, Not for Diagnostic Purposes.

### **Background**

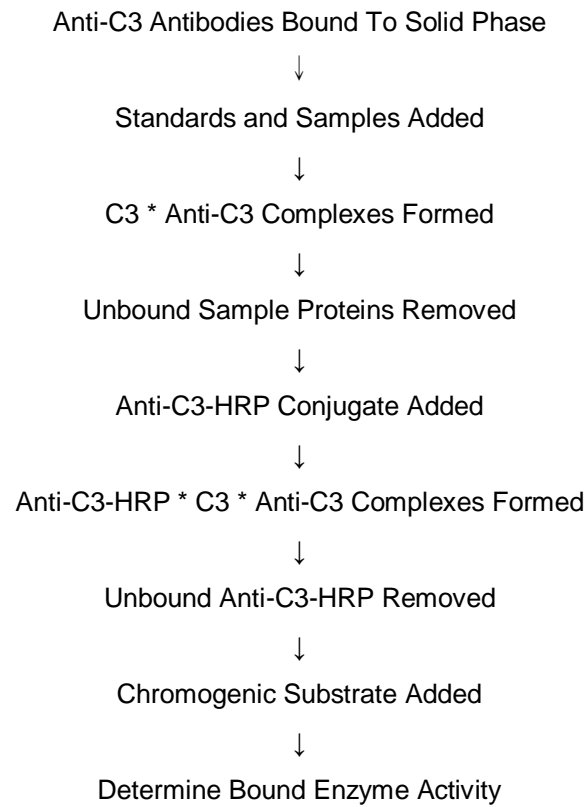
A number of serum proteins participate in acute inflammatory reactions. These include the complement, coagulation and kinin systems as well as a number of other proteins, known as acute phase proteins, that regulate acute inflammation.

The complement system is a complex set of up to 20 serum proteins. The most abundant and pivotal of the complement components is C3, which has a molecular weight of about 187 kD and consists of an alpha and beta chain.

### **Principle of the Assay**

The principle of the double antibody sandwich ELISA is represented in Figure 1. In this assay the C3 present in the sample reacts with the anti-C3 antibody which has been adsorbed to the surface of polystyrene microtiter wells. After the removal of unbound proteins by washing, anti-C3 antibody conjugated with horseradish peroxidase (HRP), are added. These enzyme-labeled antibodies form complexes with the previously bound C3. Following another washing step, the enzyme bound to the immunosorbent is assayed by the addition of a chromogenic substrate, 3,3',5,5'-tetramethylbenzidine (TMB). The quantity of bound enzyme varies directly with the concentration of C3 in the sample tested; thus, the absorbance, at 450 nm, is a measure of the concentration of C3 in the test sample. The quantity of C3 in the test sample can be interpolated from the standard curve constructed from the standards, and corrected for sample dilution.

Figure 1



## General Information

### Materials Supplied

List of component

Component	Amount
Diluent Concentrate (Running buffer): One bottle containing 5X concentrated diluent running buffer.	50 mL
Wash Solution Concentrate: One bottle containing 20X concentrated wash solution.	50 mL
Enzyme-Antibody Conjugate 100X: One vial containing affinity purified anti- Dog C3 antibody conjugated with horseradish peroxidase in a stabilizing buffer.	150 $\mu$ L
Chromogen-Substrate Solution: One vial containing 3,3',5,5'-tetramethylbenzidine (TMB) and hydrogen peroxide in citric acid buffer at pH 3.3.	12 mL
Stop Solution: One vial containing 0.3 M sulfuric acid. WARNING: Avoid contact with skin.	12 mL
Anti-Dog C3 ELISA Micro Plate: Twelve removable eight (8) well strips in well holder frame. Each well is coated with affinity purified anti-Dog C3.	96 (8x12) wells
Dog C3 Calibrator: One vial containing a lyophilized Dog C3 calibrator.	1 vial

### Storage Instruction

The expiration date for the package is stated on the box label.

- **Diluent**  
The 5X Diluent Concentrate is stable until the expiration date. The 1X working solution is stable for at least one week from the date of preparation. Both solutions should be stored at 4-8°C.
- **Wash Solution**  
The 20X Wash Solution Concentrate is stable until the expiration date. The 1X working solution is stable for at least one week from the date of preparation. Both solutions can be stored at room temperature (16-25°C) or at 4-8°C.
- **Enzyme-Antibody Conjugate**  
Undiluted horseradish peroxidase anti-C3 conjugate should be stored at 4-8°C and diluted immediately prior to use. The working conjugate solution is stable for up to 1 hour when stored in the dark.
- **Chromogen-Substrate Solution**  
The Substrate Solution should be stored at 4-8°C and is stable until the expiration date.
- **Stop Solution**  
The Stop Solution should be stored at 4-8°C and is stable until the expiration date.
- **Anti-Dog C3 ELISA Micro Plate**  
Anti-Dog C3 coated wells are stable until the expiration date and should be stored at 4-8°C in the sealed foil pouch with a desiccant pack

- Dog C3 Calibrator

The lyophilized Dog C3 calibrator should be stored at 4°C or frozen until reconstituted. The reconstituted calibrator should be aliquoted out and stored frozen (Avoid multiple freeze-thaw cycles). The working standard solutions should be prepared immediately prior to use and are stable for up to 8 hours.

### **Materials Required but Not Supplied**

- ✓ Precision pipette (2 µL to 200 µL) for making and dispensing dilutions
- ✓ Test tubes
- ✓ Microtiter washer/aspirator
- ✓ Distilled or Deionized H<sub>2</sub>O
- ✓ Microtiter reader
- ✓ Assorted glassware for the preparation of reagents and buffer solutions
- ✓ Timer

### **Precautions for Use**

- Precaution
  - ✓ For any sample that might contain pathogens, care must be taken to prevent contact with open wounds.
  - ✓ No additives or preservatives are necessary to maintain the integrity of the specimen. Avoid azide contamination.
  - ✓ Azide and thimerosal at concentrations higher than 0.1% inhibits the enzyme reaction.
- Limitation of the procedure
  - ✓ Reliable and reproducible results will be obtained when the assay procedure is carried out with a complete understanding of the instructions and with adherence to good laboratory practice.
  - ✓ Factors that might affect the performance of the assay include proper instrument function, cleanliness of glassware, quality of distilled or deionized water, and accuracy of reagent and sample pipettings.

## Assay Protocol

### Reagent Preparation

- Diluent Concentrate**  
 The Diluent solution supplied is a 5X concentrate and must be diluted 1/5 (1 part buffer concentrate, 4 parts dH<sub>2</sub>O) with distilled or deionized water.
- 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 is not uncommon 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 microtiter 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.
- Chromogen-Substrate Solution**  
 Ready to use as supplied.
- Stop Solution**  
 Ready to use as supplied.
- Anti-Dog C3 ELISA Micro Plate**  
 Ready to use as supplied. Unseal Microtiter 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. Take clear plastic tape and cover tops of strips to avoid particulates from contaminating wells along with desiccant.
- Dog C3 Calibrator**  
 Add 1.0 ml of distilled or de-ionized water to the Dog C3 calibrator and mix gently until dissolved. The calibrator is now at a concentration of 8.49 µg/ml (the reconstituted calibrator should be aliquoted and frozen if future use is intended). Dog C3 standards need to be prepared immediately prior to use (see chart below). Mix well between each step. Avoid foaming.

Standard	ng/ml	Volume added to 1x Diluent	Volume of 1x Diluent
6	200	20 µl Dog C3 Calibrator	829 µl
5	100	300 µl standard 6	300 µl
4	50	300 µl standard 5	300 µl
3	25	300 µl standard 4	300 µl
2	12.5	300 µl standard 3	300 µl
1	6.25	300 µl standard 2	300 µl
0	0		600 µl

## **Sample Preparation**

Blood should be collected by venipuncture and the serum separated from the cells, after clot formation, by centrifugation. Care should be taken to minimize hemolysis, excessive hemolysis can impact your results. Specimens may be shipped at room temperature and then stored refrigerated at 2-8°C if testing is to take place within one week after collection. If testing is to take place later than one week, specimens should be stored at -20°C. Avoid repeated freeze-thaw cycles.

## **Assay Procedure**

### 1. Dilution of Serum Samples

The assay for quantification of C3 in serum requires that each test sample be diluted before use. For a single step determination a dilution of serum at 1/40,000 is appropriate for most 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

2. To prepare a 1/40,000 dilution of sample, transfer 5 µL of sample to 995 µL of 1X diluent. This gives you a 1/200 dilution. Next, dilute the 1/200 samples by transferring 5 µL, to 995 µL of 1X diluent. You now have a 1/40,000 dilution of your sample. Mix thoroughly at each stage.

3. Bring all reagents to room temperature before use.

4. Pipette 100 µL of

Standard 0 (0.0 ng/mL) into duplicate

Standard 1 (6.25 ng/mL) into duplicate

Standard 2 (12.5 ng/mL) into duplicate

Standard 3 (25 ng/mL) into duplicate

Standard 4 (50 ng/mL) into duplicate

Standard 5 (100 ng/mL) into duplicate

Standard 6 (200 ng/mL) into duplicate

5. Pipette 100 µL of sample (in duplicate) into pre designated wells.

6. Incubate the microtiter Plate at room temperature for twenty (20 ± 2) minutes. Keep plate covered and level during incubation.

7. Following incubation, aspirate the contents of the wells.

8. 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 diluted Wash Solution, invert the plate and pour/shake out the contents in a waste container. Follow this by sharply striking the wells on absorbent paper to remove residual Wash Solution. Repeat 3 times for a total of four washes.

9. Pipette 100 µL of appropriately diluted Enzyme-Antibody Conjugate to each well. Incubate at room temperature for twenty (20 ± 2) minutes. Keep plate covered in the dark and level during incubation.

10. Wash and blot the wells as described in Steps 7/8.



11. Pipette 100  $\mu$ L of TMB Substrate Solution into each well.
  12. Incubate in the dark at room temperature for precisely ten (10) minutes.
  13. After ten minutes, add 100  $\mu$ L of Stop Solution to each well.
  14. Determine the absorbance (450 nm) of the contents of each well. Calibrate the plate reader to manufacturer's specifications.
- **Stability of the final reaction mixture**  
The absorbance of the final reaction mixture can be measured up to 2 hours after the addition of the Stop Solution. However, good laboratory practice dictates that the measurement be made as soon as possible.

## Data Analysis

### Calculation of Results

1. Subtract the average background value from the test values for each sample.
2. Using the results observed for the standards construct a Standard Curve. The appropriate curve fit is that of a four-parameter logistics curve. A second order polynomial (quadratic) or other curve fits may also be used.
3. Interpolate test sample values from the standard curve. Correct for sera dilution factor to arrive at C3 concentration in original samples.

**Resources****Plate Layout**

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