



**EAGLE**  
BIOSCIENCES

# Human C5 ELISA Assay Kit

Catalog Number:

C5H31-K01 (1 x 96 wells)

C5H31-K02 (2 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.

v. 04.19 (*effective 05May23*)

---

**Eagle Biosciences, Inc.**

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

[www.EagleBio.co](http://www.EagleBio.co)



## INTENDED USE

The Human C5 ELISA Assay Kit is to be used for the in vitro quantitative determination of C5 in plasma and serum samples. This Human C5 ELISA Assay Kit is intended for laboratory research use only and is not for use in diagnostic or therapeutic procedures. The analysis should be performed by trained laboratory professionals.

## INTRODUCTION

Complement C5 (molecular weight app. 190kDa) is a key component of the complement system. This system plays important roles in both innate and adaptive immune response and can produce an inflammatory and protective reaction to challenges from pathogens before an adaptive response can occur. The complement system consist of a complex family of proteins and receptors which are found in the circulation, in tissues and other body-fluids. Excessive and uncontrolled activation/regulation of the complement system have been implicated in various diseases such as asthma, lupus erythematosus, glomerulonephritis, various forms of arthritis, autoimmune heart disease, multiple sclerosis, inflammatory bowel disease, paroxysmal nocturnal hemoglobinuria, atypical hemolytic uremic syndrome and ischemia-reperfusion injuries, and rejection of transplanted organs. Three major pathways within the complement system have been described, namely the classical, alternative and lectin pathways which are activated by different stimuli. The classical pathway is initiated by Immune complexes; the lectin pathway by surface bound lectins; and the AP by all the surfaces that are not specifically protected against. All the 3 pathways converge on Complement C3. Cleavage of C3 induces, in turn, the cleavage of C5, successively leading to the production of the membrane attack complex (MAC). This protein complex is the final common effector responsible for lysis of susceptible microorganisms or damage cells. Mutations in C5 gene cause complement component 5 deficiency, a disease characterized by recurrent bacterial infections.

## KIT FEATURES

- Working time of 3½ hours.
- Minimum concentration which can be measured is 3.1 ng/ml.
- Measurable concentration range of 3.1 – 200 ng/ml.
- Working volume of 100 µL/well.

## Cross-reactivity

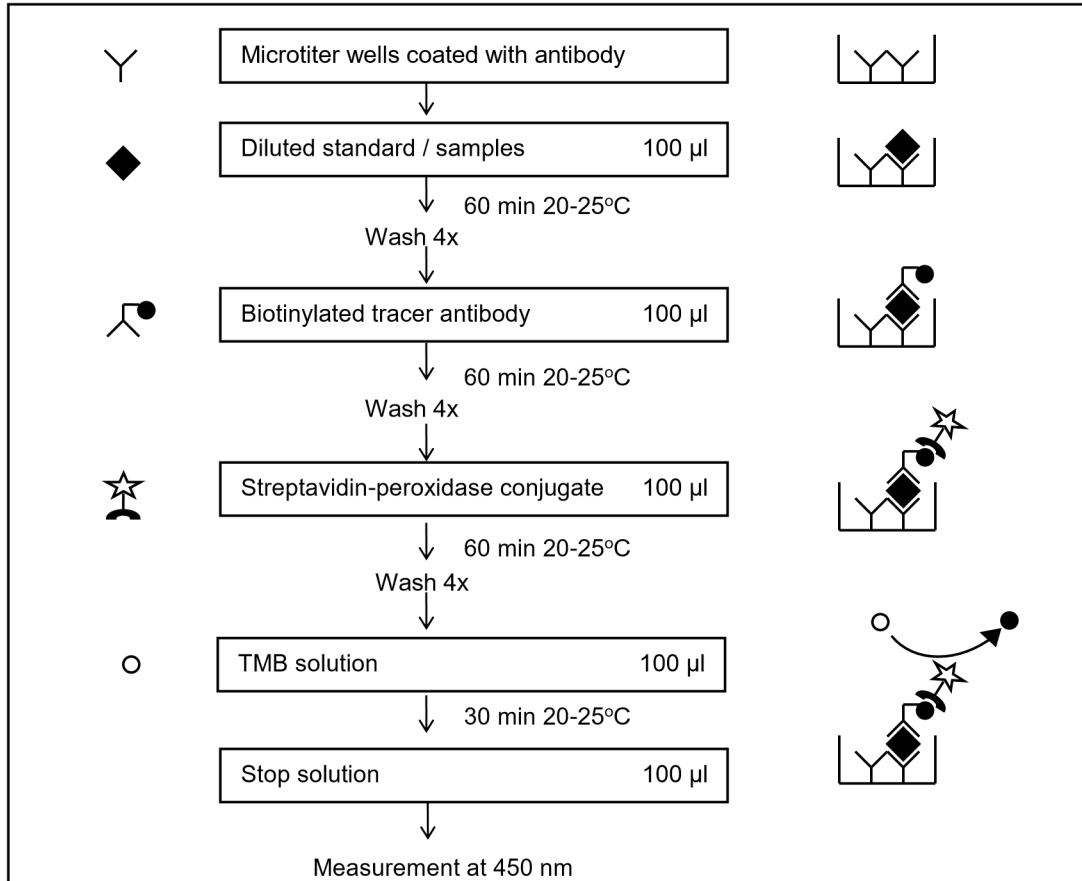
Potential cross-reacting proteins detected in the human C5 ELISA:

Cross reactant	Reactivity
Human C5a	No
Human C5a desArg	No
Mouse C5	No
Pig C5	No
Horse C5	No
Rat C5	No

Table 1 Cross-reactivity for other proteins has not been tested.



## PROTOCOL OVERVIEW



- The Human C5 ELISA is a ready-to-use solid-phase enzyme-linked immunosorbent assay based on the sandwich principle with a working time of 3½ hours.
- The efficient format of a plate with twelve disposable 8-well strips allows free choice of batch size for the assay.
- Samples and standards are incubated in microtiter wells coated with antibodies recognizing Human C5.
- Biotinylated tracer antibody will bind to the captured Human C5.
- Streptavidin-peroxidase conjugate will bind to the biotinylated tracer antibody.
- Streptavidin-peroxidase conjugate will react with the substrate, tetramethylbenzidine (TMB).
- The enzyme reaction is stopped by the addition of oxalic acid.
- The absorbance at 450 nm is measured with a spectrophotometer. A standard curve is obtained by plotting the absorbance (linear) versus the corresponding concentrations of the Human C5 standards (log).
- The Human C5 concentration of samples, which are run concurrently with the standards, can be determined from the standard curve.

## KIT COMPONENTS AND STORAGE INSTRUCTIONS

Kit component	Quantity C5H31-K01	Quantity C5H31-K02	Color code
Wash buffer 20x	1 vial (60 ml)	1 vial (60 ml)	Colorless
Dilution buffer 10x	1 vial (60 ml)	1 vial (60 ml)	Green
Standard	2 vials, lyophilized	4 vials, lyophilized	White
Tracer, biotinylated	1 vial, 1 ml lyophilized	2 vials, 1 ml lyophilized	White



Streptavidin-peroxidase	1 tube, 0.25 ml in solution	1 tube, 0.25 ml in solution	Brown
TMB substrate	1 vial (22 ml)	1 vial (22 ml)	Brown
Stop solution	1 vial (22 ml)	1 vial (22 ml)	Red
12 Microtiter strips, pre-coated	1 plate	2 plates	
Certificate of Analysis	1	1	
Manual	1	1	
Data collection sheet	2	2	

Table 2

- Upon receipt, store individual components at 2 - 8°C. Do not freeze.
- Do not use components beyond the expiration date printed on the kit label.
- The standard and tracer in lyophilized form and the streptavidin-peroxidase in concentrated solution are stable until the expiration date indicated on the kit label, if stored at 2 - 8°C.
- The exact amount of the standard is indicated on the label of the vial and the Certificate of Analysis.
- The standard is single use. After reconstitution the standard cannot be stored.
- Once reconstituted the tracer is stable for 1 month if stored at 2 - 8°C.
- The streptavidin-peroxidase can only be stored in concentrated solution and is not stable when stored diluted.
- Upon receipt, foil pouch around the plate should be vacuum-sealed and unpunctured. Any irregularities to aforementioned conditions may influence plate performance in the assay.
- Return unused strips immediately to the foil pouch containing the desiccant pack and reseal along the entire edge of the zip-seal. Quality guaranteed for one month if stored at 2 - 8°C.

#### Materials required but not provided

- Calibrated micropipettes and disposable tips.
- Distilled or de-ionized water.
- Plate washer: automatic or manual.
- Polypropylene tubes.
- Calibrated ELISA plate reader capable of measuring absorbance at 450 nm.
- Adhesive covers can be ordered separately. Please contact your local distributor.
- Centrifuge for 1 ml tubes.

#### WARNINGS AND PRECAUTIONS

##### For research use only, not for diagnostic or therapeutic use.

- This kit should only be used by qualified laboratory staff.
- Do not under any circumstances add sodium azide as preservative to any of the components.
- Do not use kit components beyond the expiration date.
- Do not mix reagents from different kits and lots. The reagents have been standardized as a unit for a given lot. Use only the reagents supplied by manufacturer.
- The assay has been optimized for the indicated standard range. Do not change the standard range.
- Open vials carefully: vials are under vacuum.
- It is advised to spin down streptavidin-peroxidase tubes before use.
- Do not ingest any of the kit components.
- Kit reagents contain 2-chloroacetamide as a preservative. 2-Chloroacetamide is harmful in contact with skin and toxic if swallowed. In case of accident or if you feel unwell, seek medical advice immediately.



- The TMB substrate is light sensitive, keep away from bright light. The solution should be colorless until use.
- The stop solution contains 2% oxalic acid and can cause irritation or burns to respiratory system, skin and eyes. Direct contact with skin and eyes should be strictly avoided. If contact occurs, rinse immediately with plenty of water and seek medical advice.
- Incubation times, incubation temperature and pipetting volumes other than those specified may give erroneous results.
- Do not reuse microwells or pour reagents back into their bottles once dispensed.
- Handle all biological samples as potentially hazardous and capable of transmitting diseases.
- Hemolyzed, hyperlipemic, heat-treated or contaminated samples may give erroneous results.
- Use polypropylene tubes for preparation of standard and samples. Do not use polystyrene tubes or sample plates.
- The standard is of human origin. It was tested for various viruses and found negative. Since no test method can offer complete assurance that infectious agents are absent, this reagent should be handled as any potentially infectious human serum or blood specimen. Handle all materials in contact with this reagent according to guide-lines for prevention of transmission of blood-borne infections.

## **SAMPLE PREPARATION**

### **Collection and handling**

#### **Serum or plasma**

Collect blood using normal aseptic techniques. Blood samples should be kept on ice. If serum is used, separate serum from blood after clotting at room temperature within one hour by centrifugation (1500xg at 4°C for 15 min). Transfer the serum to a fresh polypropylene tube.

If plasma is used, separate plasma from blood within 20 minutes after blood sampling by centrifugation (1500xg at 4°C for 15 min). Transfer the plasma to a fresh propylene tube.

#### **Storage**

Store samples below -20°C, preferably at -70°C in polypropylene tubes. Storage at -20°C can affect recovery of human C5. Use samples within 24 hours after thawing. Avoid multiple freeze-thaw cycles which may cause loss of human C5 and give erroneous results.

Do not use hemolyzed, hyperlipemic, heat-treated or contaminated samples. Before performing the assay, samples should be brought to room temperature (18 - 25°C) and mixed gently. Prepare all samples (controls and test samples) prior to starting the assay procedure. Avoid foaming.

### **Dilution procedures**

#### **Plasma and Serum samples**

Human C5 can be measured accurately if serum or plasma samples of healthy donors are diluted at least 500x with supplied dilution buffer in propylene tubes. Optimal dilution is dependent on sample quality and expected C5 quantity and should be determined for the specific sample set/study.

#### **Comment regarding recommended sample dilution**

The mentioned dilution for samples is a minimum dilution and should be used as a guideline. The recovery of human C5 from an undiluted sample is not 100% and may vary from sample to sample. When testing less diluted samples it is advisable to run recovery experiments to determine the influence of the matrix on the detection of human C5.

Do not use polystyrene tubes or sample plates for preparation or dilution of the samples.



## Guideline for dilution of samples

Please see the table 3 for recommended sample dilutions. Volumes are based on a total volume of at least 230  $\mu$ L of diluted sample, which is sufficient for one sample in duplicate in the ELISA. For dilution of samples we recommend to use at least 10  $\mu$ l of sample.

Dilution	Pre-dilution	Amount of sample or pre-dilution required	Amount of Dilution buffer required
1. 10x	Not necessary	25 $\mu$ l (sample)	225 $\mu$ l
2. 20x	Not necessary	15 $\mu$ l (sample)	285 $\mu$ l
3. 50x	Not necessary	10 $\mu$ l (sample)	490 $\mu$ l
4. 100x	Not necessary	10 $\mu$ l (sample)	990 $\mu$ l
5. 500x	Recommended: 10x (see nr.1)	10 $\mu$ l (pre-dilution)	490 $\mu$ l
6. 1000x	Recommended: 10x (see nr.1)	10 $\mu$ l (pre-dilution)	990 $\mu$ l
7. 5000x	Recommended: 50x (see nr.2)	10 $\mu$ l (pre-dilution)	990 $\mu$ l
8. 20000x	Recommended: 100x (see nr.3)	10 $\mu$ l (pre-dilution)	1990 $\mu$ l

Table 3

## REAGENT PREPARATION

Allow all the reagents to equilibrate to room temperature (20 – 25°C) prior to use. Return to proper storage conditions immediately after use.

### Wash buffer

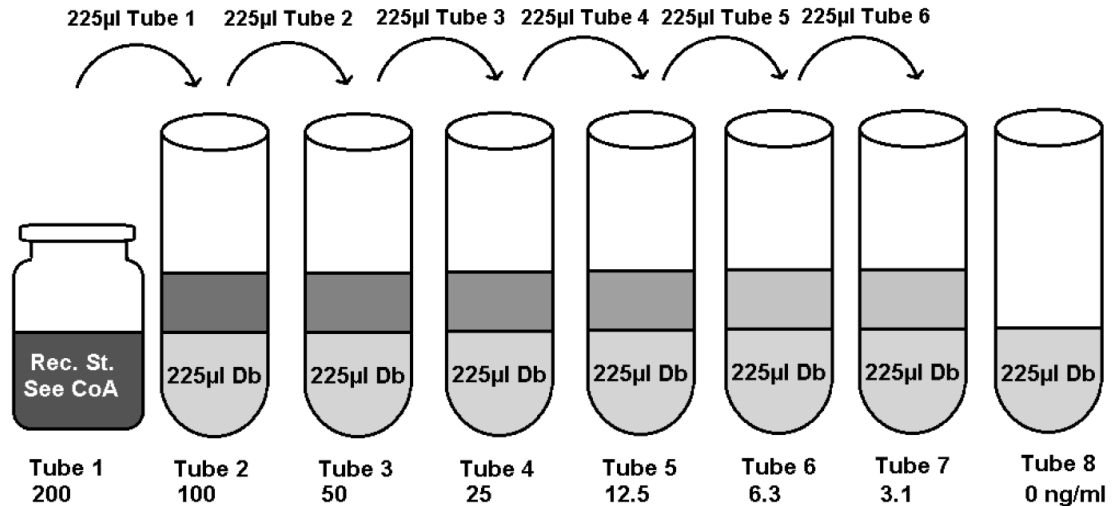
Prepare wash buffer by mixing 60 ml of 20x wash buffer with 1140 ml of distilled or de-ionized water, which is sufficient for 2 x 96 tests. In case less volume is required, prepare the desired volume of wash buffer by diluting 1 part of the 20x wash buffer with 19 parts of distilled or deionized water.

### Dilution buffer

Prepare dilution buffer by mixing 60 mL of the 10x dilution buffer with 540 mL of distilled or de-ionized water, which is sufficient for 2 x 96 tests. In case less volume is required, prepare the desired volume of dilution buffered by dilution 1 part of the 10x dilution buffer with 9 parts of distilled or de-ionized water. Concentrated dilution buffer may contain crystals. In case of crystals do not disappear at room temperature within one hour, concentrated dilution buffer can be warmed up to 37°C. Do not shake the solution.

### Standard solution

The standard is reconstituted by pipetting the amount of dilution buffer mentioned on the CoA in the standard vial. Use the standard vial as Tube 1 in Figure 1. Prepare each human C5 standard in polypropylene tubes by serial dilution of the reconstituted standard with dilution buffer as shown in Figure 1\*. After reconstitution the standard cannot be stored for repeated use.



\*) CoA: Certificate of Analysis; Rec.St: Reconstituted Standard; Db: dilution buffer

Figure 1

### Tracer solution

The tracer is reconstituted by pipetting 1 ml distilled or de-ionized water. Dilute the reconstituted 1 ml tracer with 11 ml dilution buffer, which is sufficient for 1 x 96 tests. Where less volume is required, prepare the desired volume of tracer by diluting 1 part of the reconstituted tracer with 11 parts of dilution buffer.

### Streptavidin-peroxidase solution

It is advised to spin down streptavidin-peroxidase tubes before use. Prepare the streptavidin-peroxidase solution by mixing 0.25 ml of the 100x streptavidin-peroxidase solution with 24.75 ml dilution buffer, which is sufficient for 2 x 96 tests. In case less volume is required, prepare the desired volume of streptavidin-peroxidase solution by diluting 1 part of the 100x streptavidin-peroxidase solution with 99 parts of dilution buffer.

### ELISA PROTOCOL

Bring all reagents to room temperature (20 - 25°C) before use.

1. Determine the number of test wells required, put the necessary microwell strips into the supplied frame, and fill out the data collection sheet. Return the unused strips to the storage bag with desiccant, seal and store at 2 - 8°C.
2. Transfer 100 µl in duplicate of standard, samples, or controls into appropriate wells. Do not touch the side or bottom of the wells.
3. Cover the tray and tap the tray to eliminate any air bubbles. Be careful not to splash liquid onto the cover.
4. Incubate the strips or plate for 1 hour at room temperature.
5. Wash the plates 4 times with wash buffer as follows\*:
  - a. Carefully remove the cover, avoid splashing.
  - b. Empty the plate by inverting plate and shaking contents out over the sink, keep inverted and tap dry on a thick layer of tissues.
  - c. Add 200 µl of wash buffer to each well, wait 20 seconds, empty the plate as described in 5b.
  - d. Repeat the washing procedure 5b/5c three times.
  - e. Empty the plate and gently tap on thick layer of tissues.
6. Add 100 µl of diluted tracer to each well using the same pipetting order as applied in step



2. Do not touch the side or bottom of the wells.
7. Cover the tray and incubate the tray for 1 hour at room temperature.
8. Repeat the wash procedure described in step 5a-e.
9. Add 100  $\mu$ l of diluted streptavidin-peroxidase to each well, using the same pipetting order as applied in step 2. Do not touch the side or bottom of the wells.
10. Cover the tray and incubate the tray for 1 hour at room temperature.
11. Repeat the wash procedure described in step 5a-e.
12. Add 100  $\mu$ l of TMB substrate to each well, using the same pipetting order as applied in step 2. Do not touch the side or bottom of the wells.
13. Cover the tray and incubate the tray for 30 minutes at room temperature. It is advised to control the reaction on the plate regularly. In case of strong development the TMB reaction can be stopped sooner. Avoid exposing the microwell strips to direct sunlight. Covering the plate with aluminum foil is recommended.
14. Stop the reaction by adding 100  $\mu$ l of stop solution with the same sequence and timing as used in step 12. Mix solutions in the wells thoroughly by gently swirling the plate. Gently tap the tray to eliminate any air bubbles trapped in the wells.
15. Read the plate within 30 minutes after addition of stop solution at 450 nm using a plate reader, following the instructions provided by the instrument's manufacturer.

\*) In case plate washer is used, please note: use of a plate washer can result in higher background and decrease in sensitivity. We advise validation of the plate washer with the manual procedure. Make sure the plate washer is used as specified for the manual method.

## INTERPRETATION OF RESULTS

- Calculate the mean absorbance for each set of duplicate standards, control and samples.
- If individual absorbance values differ by more than 15% from the corresponding mean value, the result is considered suspect and the sample should be retested.
- The mean absorbance of the zero standard should be less than 0.3.
- Create a standard curve using computer software capable of generating a good curve fit. The mean absorbance for each standard concentration is plotted on the vertical (Y) axis versus the corresponding concentration on the horizontal (X) axis (logarithmic scale).
- If samples have been diluted, the concentration read from the standard curve must be multiplied by the dilution factor.
- Samples that give a mean absorbance above the absorbance for the highest standard concentration are out of range of the assay. These samples should be retested at a higher dilution.

## TECHNICAL HINTS

- User should be trained and familiar with ELISA assays and test procedure.
- If you are not familiar with the ELISA technique it is recommended to perform a pilot assay prior to evaluation of your samples. Perform the assay with a standard curve only following the instructions.
- Improper or insufficient washing at any stage of the procedure will result in either false positive or false negative results. Completely empty wells before dispensing wash buffer, fill with wash buffer as indicated for each cycle and do not allow wells to sit uncovered or dry for extended periods.
- Since exact conditions may vary from assay to assay, a standard curve must be established for every run. Sample should be referred to the standard curve prepared on the same plate.
- Do not mix reagents from different batches, or other reagents and strips. Remainders should not be mixed with contents of freshly opened vials.





- Each time the kit is used, fresh dilutions of standard, sample, tracer, streptavidin-peroxidase and buffers should be made.
- Caps and vials are not interchangeable. Caps should be replaced on the corresponding vials.
- To avoid cross-contamination, change pipette tips between reagent additions of each standard, between sample additions, and between reagent additions. Also, use separate reservoirs for each reagent.
- The waste disposal should be performed according to your laboratory regulations.

### Technical support

Do not hesitate to contact our technical support team at [info@eaglebio.com](mailto:info@eaglebio.com) for inquiries and technical support regarding the human C5 ELISA.

### QUALITY CONTROL

The Certificate of Analysis included in this kit is lot-specific and is to be used to verify results obtained by your laboratory. The absorption values provided on the Certificate of Analysis are to be used as a guideline only. The results obtained by your laboratory may differ.

This assay is designed to eliminate interference by soluble receptors, binding proteins, and other factors present in biological samples. Until all factors have been tested in the Eagle Biosciences, Inc. immunoassay, the possibility of interference cannot be excluded. For optimal performance of this kit, it is advised to work according to good laboratory practice.

### PERFORMANCE CHARACTERISTICS

#### Recovery

Normal human blood samples (plasma) containing baseline levels of human C5, were spiked with human C5, in concentrations of 100 and 12.5 ng/mL. Samples with and without human C5, were incubated for 30 minutes at room temperature. Samples were measure using the ELISA. Values for human C5, ranged between 96% and 109%.

#### Precision and Reproducibility

The intra-assay was tested by testing 4 EDTA plasma samples in triplo

Sample	Average Amount (µg)	CV%
Sample 1	18.6	4.1
Sample 2	15.5	3.5
Sample 3	19.3	4.6
Sample 4	21.3	1.9

The inter-assay was tested by testing 4 EDTA plasma samples by 2 operators.

Sample	Average Amount (µg)	CV%
Sample 1	18.4	6.3
Sample 2	15.2	5.1
Sample 3	18.9	6.3
Sample 4	20.7	7.1

### TROUBLESHOOTING

Warranty claims and complaints in respect of deficiencies must be logged before expiry date of the product. A written complaint containing lot number of the product and experimental data should be sent to [info@eaglebio.com](mailto:info@eaglebio.com).



Suggestions summarized below in Table 4 can be used as a guideline in the case of unexpected assay results.

Low absorbance	High absorbance	Poor duplicates	All wells positive	All wells negative	Possible cause
<input type="checkbox"/>	<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>	Kit materials or reagents are contaminated or expired
<input type="checkbox"/>					Incorrect reagents used
<input type="checkbox"/>		<input type="checkbox"/>	<input type="checkbox"/>		Lyophilized reagents are not properly reconstituted
<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	Incorrect dilutions or pipetting errors
<input type="checkbox"/>		<input type="checkbox"/>			Improper plastics used for preparation of standard and/or samples
<input type="checkbox"/>	<input type="checkbox"/>				Improper incubation times or temperature
		<input type="checkbox"/>			Especially in case of 37°C incubation: plates are not incubated uniformly
<input type="checkbox"/>					Assay performed before reagents were brought to room temperature
<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>	Procedure not followed correctly
				<input type="checkbox"/>	Omission of a reagent or a step
		<input type="checkbox"/>			Poor mixing of samples
	<input type="checkbox"/>		<input type="checkbox"/>		Low purity of water
	<input type="checkbox"/>	<input type="checkbox"/>			Strips were kept dry for too long during/after washing
	<input type="checkbox"/>	<input type="checkbox"/>	<input type="checkbox"/>		Inefficient washing
	<input type="checkbox"/>	<input type="checkbox"/>			Cross-contamination from other samples or positive control
		<input type="checkbox"/>	<input type="checkbox"/>		TMB solution is not clear or colorless
<input type="checkbox"/>	<input type="checkbox"/>				Wrong filter in the microtiter reader
	<input type="checkbox"/>	<input type="checkbox"/>			Airbubbles
		<input type="checkbox"/>			Imprecise sealing of the plate after use
<input type="checkbox"/>					Wrong storage conditions
<input type="checkbox"/>					Lamp in microplate reader is not functioning optimally

Table 4



## Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit, please contact the Technical Service Team at Eagle Biosciences, Inc. at [info@eaglebio.com](mailto:info@eaglebio.com) or at 866-411-8023.