



EAGLE  
BIOSCIENCES

# **Anti-IgE Receptor (FcεR1α) Autoantibody ELISA Kit**

Catalog Number:  
KTR-802 (1 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.  
v. 2.2 (30 SEP 24)

---

Eagle Biosciences, Inc.  
20A Northwest Blvd., Suite 112, Nashua, NH 03063  
Phone: 617-419-2019 Fax: 617-419-1110  
[www.EagleBio.com](http://www.EagleBio.com)



## **INTENDED USE**

The Eagle Biosciences Anti-IgE Receptor (FcεR1α) Autoantibody ELISA (enzyme-linked immunosorbent assay) Kit is produced for the quantitative determination of human anti-human IgE Receptor (FcεR1α) autoantibody levels in human serum or plasma samples. For research use only and not to be used in diagnostic procedures.

## **SUMMARY OF PHYSIOLOGY**

The presence of anti-IgE receptor (FcεR1α) antibodies in sample serum or plasma has been associated with chronic spontaneous urticaria (CSU), which is a common skin disorder affecting 0.5% to 1.8% of the general population. It is characterized by repeated occurrence of short-lived cutaneous wheals accompanied by redness and itching. Although the CSU symptom is very much similar to those of acute urticaria triggered by allergens, in most CSU cases, there is no definite identifiable direct external triggering factor. As of today the pathogenesis of CSU has not been fully elucidated, a proportion of people with CSU have been found to have functional autoantibodies. In CSU samples, circulation human anti-FcεR1α autoantibodies are seen in 30% to 60% and anti-IgE autoantibodies in 5% to 10%. These human autoantibodies are mainly IgG subtype.

## **PRINCIPLE OF THE ASSAY**

This ELISA is designed, developed and produced for the quantitative measurement of human anti-hIgE receptor (FcεR1α) autoantibodies in serum and plasma samples.

Assay calibrators, controls and diluted samples are directly added to wells of a microplate that is coated with recombinant human IgE receptor protein. After the first incubation period, anti-IgE receptor antibodies bind to the human IgE receptor protein on the wall of microtiter well and unbound proteins in each microtiter well are washed away. Highly purified Protein A labeled with horseradish peroxidase is then added to each microtiter well. After the second incubation period, a complex of coated human IgE receptor Anti- hIgE receptor antibody peroxidase-labeled Protein A is formed. The unbound protein is removed in the subsequent washing step. The wells are then incubated with a substrate solution in a timed reaction and subsequently measured in a spectrophotometric microplate reader. The enzymatic activity of the immunocomplex bound to the IgE receptor protein on the wall of the microtiter well is directly proportional to the amount of anti-IgE receptor antibody in the sample. A calibration curve is generated by plotting the absorbance versus the respective anti-IgE receptor concentration for each calibrator on a cubic spline curve fit. The concentration of IgE receptor antibody in test samples is determined directly from this calibration curve.

## **REAGENTS: Preparation and Storage**

This test kit must be stored at 2 – 8 °C upon receipt. For the expiration date of the kit refer to the label on the kit box. All components are stable until this expiration date.

**Prior to use, allow all reagents to come to room temperature.** Reagents from different kit lot numbers should not be combined or interchanged.



### **1. Human IgE Receptor Protein Coated Microplate**

One well-breakable microplate with 12 x eight strips (96 wells total) coated with purified recombinant human IgE receptor protein. The plate is framed and sealed in a zipper foil bag with a desiccant. This reagent should be stored at 2 – 8 °C and is stable until the expiration date on the kit box.

### **2. HRP Conjugated Protein A**

One bottle containing 0.6 mL of concentrated horseradish peroxidase conjugated Protein A. Before the use of it, this concentrated reagent should be diluted with HRP Conjugated Protein A Diluent. This reagent should be stored at 2 – 8 °C and is stable until the expiration date on the kit box

### **3. HRP Conjugated Protein A Diluent**

One bottle containing 12 mL of ready-to-use HRP Conjugated Protein A Diluent. This reagent is used to dilute the HRP Conjugated Protein A. It should be stored at 2 – 8°C and stable until the expiration date on the kit box.

### **4. ELISA Wash Concentrate**

One bottle contains 30 mL of 30-fold concentrated wash buffer. Before use, the contents must be diluted with 870 mL of distilled water and mixed well. Upon dilution this yields a working wash solution containing a surfactant in phosphate- buffered saline with a non-azide and non-mercury based preservative. The diluted wash buffer should be stored at room temperature and is stable until the expiration date on the kit box.

### **5. ELISA HRP Substrate**

One bottle contains 12 mL of ready-to-use tetramethylbenzidine (TMB) with stabilized hydrogen peroxide. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

### **6. ELISA Stop Solution**

One bottle contains 12 mL of ready-to-use 0.5 M sulfuric acid. This reagent should be stored at 2 – 8°C or room temperature and is stable until the expiration date on the kit box.

### **7. Anti-human IgE Receptor Calibrators**

Five vials each containing 1 mL of a different level of ready-to- use anti-human IgE Receptor antibodies in a liquid protein matrix with a non-azide based preservative. Refer to vials for exact concentration for each calibrator. These reagents should be stored at 2 – 8 °C and are stable until the expiration date on the kit box.

### **8. Anti-human IgE Receptor Controls**

Two vials each containing 1 mL of a different level of ready-to- use anti-human IgE Receptor antibodies in a liquid protein matrix with a non-azide based preservative. Refer to vials for exact concentration range for each control. Both controls should be stored at 2 – 8 °C and are stable until the expiration date on the kit box.

### **9. Anti-human IgE Receptor Sample Diluent**

One bottle containing 125 mL of ready-to-use anti-human IgE Receptor Sample Diluent. This reagent is used to dilute the samples. It should be stored at 2 – 8°C and is stable until the expiration date.



## SAFETY PRECAUTIONS

The reagents must be used for professional use only. Source material (e.g. highly purified bovine serum albumin) was derived in the contiguous 48 United States. It was obtained only from healthy donor animals maintained under veterinary supervision and found free of contagious diseases. Wear gloves while performing this assay and handle these reagents as if they are potentially infectious. Avoid contact with reagents containing TMB, hydrogen peroxide, or sulfuric acid. TMB may cause irritation to skin and mucous membranes and cause an allergic skin reaction. TMB is a suspected carcinogen. Sulfuric acid may cause severe irritation on contact with skin. Do not get in eyes, on skin, or on clothing. Do not ingest or inhale fumes. On contact, flush with copious amounts of water for at least 15 minutes. Use Good Laboratory Practices.

## MATERIALS REQUIRED BUT NOT PROVIDED

- Precision single channel pipettes capable of delivering 10  $\mu$ L, 100  $\mu$ L, and 1000  $\mu$ L etc.
- Disposable pipette tips suitable for above volume dispensing.
- Disposable 12 x 75 mm or 13 x 100 plastic tubes.
- Disposable plastic 100 mL and 1000 mL bottle with caps.
- Aluminum foil.
- Deionized or distilled water
- Clean test tubes.
- Plastic microtiter well cover or polyethylene film.
- ELISA multichannel wash bottle or automatic (semi-automatic) washing system.
- Spectrophotometric microplate reader capable of reading
- Absorbance at 450 nm.

## SPECIMEN COLLECTION

Only 10 $\mu$ L of human serum or plasma is required for anti-human IgE receptor antibody measurement in duplicate. No special preparation of individual is necessary prior to specimen collection. In the case of serum, whole blood should be collected and must be allowed to clot for a minimum of 30 minutes at room temperature before the serum is separated by centrifugation (850 – 1500 x g for 10 minutes). The serum should be separated from the clot within three hours of blood collection and transferred to a clean test tube. Serum or plasma samples should be stored at 2 – 8°C if the assay is to be performed within 72 hours. Otherwise, samples should be stored at -20°C or below until measurement. Avoid repeated (more than three times) freezing and thawing of specimen.

## ASSAY PROCEDURE

### 1. Sample Preparation

1. The undiluted serum and plasma can be stored at 2-8°C for up to 1 week before use, otherwise, samples need to be stored at < -20°C before use. Avoid more than 3x freeze/thaw cycle.
2. Before testing, each serum and plasma sample should be diluted 1:100 with Sample diluent. For example, add 1mL of sample diluent + 10  $\mu$ L sample into a clean test tube.



3. The unused extracted (1:100) samples should be sealed and stored at  $< -20^{\circ}\text{C}$  for future use. It is optional to store the extracted samples at  $2-8^{\circ}\text{C}$  and is stable for up to 1 week and/or room temperature ( $20\text{C} - 25^{\circ}\text{C}$ ) for up to 3 days. Otherwise, extracted samples need to be stored at  $< -20^{\circ}\text{C}$ . Avoid more than 3x freeze/thaw cycle.

## 2. Reagent Preparation

1. Prior to use allow all reagents to come to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.
2. ELISA Wash Concentrate must be diluted to working solution prior use. Please see REAGENTS section for details.
3. HRP Conjugated Protein A must be diluted 1:21 with HRP Conjugated Protein A Diluent before use. Refer to the table below:

Strip no.	Protein A HRP Diluent	Protein A HRP
1	1 mL	50 $\mu\text{L}$
2	2 mL	100 $\mu\text{L}$
3	3 mL	150 $\mu\text{L}$
4	4 mL	200 $\mu\text{L}$
5	5 mL	250 $\mu\text{L}$
6	6 mL	300 $\mu\text{L}$
7	7 mL	350 $\mu\text{L}$
8	8 mL	400 $\mu\text{L}$
9	9 mL	450 $\mu\text{L}$
10	10 mL	500 $\mu\text{L}$
11	11 mL	550 $\mu\text{L}$
12	12 mL	600 $\mu\text{L}$

## 3. Assay Procedure

1. Place a sufficient number of human IgE Receptor Protein coated microwell strips/wells in a holder to run Anti- IgE Receptor calibrators, controls and unknown diluted samples in duplicate.

### Test Configuration:

ROW	STRIP 1	STRIP 2	STRIP 3
A	Calibrator 1	Calibrator 5	SAMPLE 2
B	Calibrator 1	Calibrator 5	SAMPLE 2
C	Calibrator 2	C 1	SAMPLE 3
D	Calibrator 2	C 1	SAMPLE 3
E	Calibrator 3	C 2	SAMPLE 4
F	Calibrator 3	C 2	SAMPLE 4
G	Calibrator 4	SAMPLE 1	
H	Calibrator 4	SAMPLE 1	



2. Add 100  $\mu\text{L}$  of calibrators, controls and diluted (1:100) samples into the designated microwells.
3. Cover the plate with one plate sealer and foil and incubate plate at room temperature (20– 25°C), static for 1 hour.
4. Remove foil and plate sealer. Aspirate the contents of each well. Wash each well 5 times by dispensing 350  $\mu\text{L}$  of working wash solution into each well and then completely aspirating the contents. Alternatively, an automated microplate washer can be used.
5. Add 100  $\mu\text{L}$  of diluted HRP Conjugated Protein A to each of the wells.
6. Cover the plate with a plate sealer and aluminum foil and incubate plate at room temperature (20 – 25°C), static for 30 minutes.
7. Remove foil and plate sealer. Aspirate the contents of each well. Wash each well 5 times by dispensing 350  $\mu\text{L}$  of working wash solution into each well and then completely aspirating the contents. Alternatively, an automated microplate washer can be used.
8. Add 100  $\mu\text{L}$  of ELISA HRP Substrate into each of the wells.
9. Cover the plate with aluminum foil to avoid exposure to light.
10. Incubate plate at room temperature (20 – 25°C), static for 20 minutes.
11. Remove the aluminum foil and plate sealer. Add 100  $\mu\text{L}$  of ELISA Stop Solution into each of the wells. Mix gently.
12. Read the absorbance at 450 nm within 10 minutes in a microplate reader

## PROCEDURAL NOTES

1. It is recommended that all calibrators, controls and unknown samples be assayed in duplicate. The average absorbance reading of each duplicate should be used for data reduction and the calculation of results.
2. For samples with concentration higher than level 5 calibrator, it is recommended to measure diluted the specimen with sample diluent at 1:200, 1:400, etc. for a more accurate report. The result is then multiplied by 2, 4 etc. to obtain the corrected anti IgE receptor antibody concentration.
3. Keep light-sensitive reagents in the original amber bottles.
4. Store any unused human IgE receptor protein coated strips sealed in the foil bag with desiccant to protect from moisture.
5. Careful technique and use of properly calibrated pipeting devices are necessary to ensure reproducibility of the test.
6. Incubation times or temperatures other than those stated in this insert may affect the results.
7. Avoid introducing air bubbles into the microwell as this could result in lower binding efficiency and higher CV% of duplicate readings.
8. . All reagents should be mixed gently and thoroughly prior use. Avoid foaming.

## INTERPRETATION OF THE RESULTS

1. Calculate the average absorbance for each pair of duplicate test results.
2. Subtract the average absorbance of the calibrator 1 (0U/mL) from the average absorbance of all other readings to obtain corrected absorbance.
3. The calibration curve is generated by the corrected absorbances of all standard levels on the ordinate against the calibrator concentration on the abscissa using point-to- point or log-log paper. Appropriate computer assisted data reduction programs may also be used for the calculation of results. We recommend using a cubic spline curve fit.

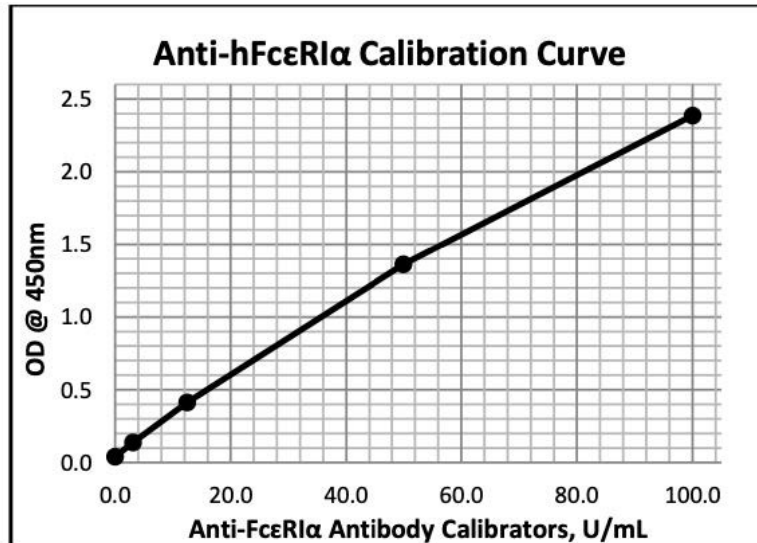


The Anti-IgE receptor antibody concentrations for the controls and samples are read directly from the standard curve using their respective corrected absorbance.

### EXAMPLE DATA AND STANDARD DATA

A typical absorbance data and the resulting calibration curve from this ELISA are represented. This curve should not be used in lieu of standard curve run with each assay.

Well I.D.	OD 450 nm Absorbance			Results U/mL
	Readings	Average	Corrected	
0 U/mL	0.040 0.041	0.041	0.000	
3.1 U/mL	0.140 0.137	0.139	0.098	
12.5 U/mL	0.411 0.415	0.413	0.373	
50 U/mL	1.384 1.342	1.363	1.323	
100 U/mL	2.373 2.400	2.387	2.346	
Control 1	0.290 0.275	0.283	0.242	7.95
Control 2	0.879 0.924	0.902	0.861	30.8





## **EXPECTED VALUES**

Serum samples from 60 normal donors (male: 13, Female: 47) with average age of 39 (range: 19 – 67) were measured with this test. The 95th percentile normal range is 0.8 U/mL – 7.5 U/mL, with a Median (P50) of 1.89U/mL, P25 of 1.32U/mL, and P75 of 2.51U/mL.

Epitope Diagnostics recommends that the normal cut off of this test is 7.5 U/mL. It is highly recommended that each laboratory establish its own normal cut-off level.

## **LIMITATION OF THE PROCEDURE**

1. Since there is no Gold Standard concentration or international standard available for anti-IgE measurement, the values of assay standards were established and validated by Epitope Diagnostics. Results obtained with different assay methods or kits cannot be used interchangeably.
2. Bacterial or fungal contamination of serum specimens or reagents, or cross-contamination between reagents may cause erroneous results.
3. Water deionized with polyester resins may reduce the activity the horseradish peroxidase enzyme.

## **QUALITY CONTROL**

To assure the validity of the results each assay should include adequate controls with known anti-IgE receptor IgG levels. We recommend that all assays include the laboratory's own anti-IgE receptor controls in addition to those provided with this kit.

## **PERFORMANCE CHARACTERISTICS**

### **Sensitivity**

The sensitivity of this Human Anti h-IgE Receptor ELISA as determined by the 95% confidence limit on 16 replicates determinations are the following:

- Limit of Blank (LoB) = 0.054 U/mL
- Limit of Detection (LoD) = 0.374 U/mL
- Limit of Quatification (LoQ) = 0.694 U/mL

### **Precision**

The intra-assay precision was validated by measuring three (3) samples in 16 replicates determinations.

The inter-assay precision was validated by measuring two samples in 12 separate assays.





	Inter-Assay		Intra-Assay		
	Sample 1	Sample 2	Sample 1	Sample 2	Sample 3
<b>Mean</b>	7.693	31.172	1.562	7.591	29.826
<b>Std Dev</b>	0.396	0.967	0.124	0.611	1.587
<b>%CV</b>	5.2%	3.1%	7.9%	8.1%	5.3%

### Linearity

Two samples were serially diluted with sample diluent and tested. The results of dilution recovery value are summarized as follows:

#	DILUTION	OBSERVED VALUE (U/mL)	EXPECTED VALUE (U/mL)	RECOVERY %
1	1:100	17.0	-	-
	1:200	9.3	8.5	110%
	1:400	4.5	4.2	105%
	1:800	2.2	2.1	103%
2	1:100	38.3	-	-
	1:200	20.8	19.2	109%
	1:400	9.4	9.6	98%
	1:800	3.9	4.8	82%

Three Standards (level 5, level 4, level 3) were serially diluted with sample diluent and tested. The results of dilution recovery value are summarized as follows:

#	DILUTION	OBSERVED VALUE (U/mL)	EXPECTED VALUE (U/mL)	RECOVERY %
1	Level 5	100.0	-	-
	1:2	48.1	50.0	96%
	1:4	22.2	25.0	89%
	1:8	11.3	12.5	90%
2	Level 4	50.0	-	-
	1:2	24.5	25.0	98%
	1:4	11.0	12.5	88%
	1:8	5.3	6.3	85%
2	Level 3	12.5	-	-
	1:2	6.1	6.3	97%
	1:4	3.0	3.1	96%
	1:8	1.4	1.6	87%



### Spiked Recovery

Two samples were spiked (50%-50%) with Calibrators 2-4 in equal volume and assayed. The results indicate below:

Sample	Expected	Observed	% Recovery
<b>A</b>	-	7.1	-
+ Level 2 :			
3.125U/mL	5.1	5.1	101%
+ Cal 3 :			
12.5U/mL	9.8	9.0	92%
+ Cal 4:			
50U/mL	28.6	28.4	99%
<b>B</b>	-	15.5	-
+ Level 2 :			
3.125U/mL	9.3	9.7	104%
+ Cal 3 :			
12.5U/mL	14.0	13.4	95%
+ Cal 4:			
50U/mL	32.8	29.3	89%

### High Dose "hook" effect

This assay has showed that it didn't exhibit any high dose "hook" effect up to 64,000 U/mL.

### Interference

One positive and one negative sample are added with 5% volume of interference materials to reach a final concentration shown in the table below. All samples are tested in an assay in duplicate

Interferant	Test Control (U/mL)	Interference Result (U/mL)	%Bias
Bilirubin 0.4mg/mL	1.20	1.17	-3%
	26.89	26.71	-1%
Bilirubin 2mg/mL	1.20	1.28	7%
	26.89	25.80	-4%
Bilirubin 2mg/mL	1.20	1.42	18%
	26.89	31.98	19%
Hemoglobin 0.4mg/mL	1.22	1.11	-9%
	29.42	36.49	24%
Hemoglobin 2mg/mL	1.22	1.18	-3%
	29.42	34.82	18%
Hemoglobin 2mg/mL	1.22	1.07	-12%
	29.42	32.41	10%
Lipid 8mg/mL	1.22	1.21	0%
	29.42	31.57	7%
Lipid 40mg/mL	1.22	1.87	54%
	29.42	33.34	13%
Lipid 200mg/mL	1.22	1.54	26%
	29.42	33.57	14%



## REFERENCES

1. Christine B Cho, et al. Autoantibodies in chronic idiopathicurticaria and nonurticarial systemic autoimmune disorders. *AnnAllergy Asthma Immunol.* 2013;110:29-33
2. Hide M,et al. Autoantibodies against the high-affinity IgE receptor as a cause of histamine release in chronic urticaria. *N Engl J Med.*1993;328:1599-1604.
3. Bastsetseg Ulambayer, et al. Detection of circulating IgG autoantibody to FcεR1α in sera from chronic spontaneous urticaria patients. *J Microbiology* 2017;
4. Marta Ferrer, et al. Progress and Challenges in the Understanding of Chronic Urticaria. *Allergy, Asthma, and Clinical Immunology* 2017;3:31-35

## 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.