



# IL1RN (Human) ELISA Kit

Catalog Number KA1066

96 assays

Version: 03

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>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	5
<b>Assay Protocol .....</b>	<b>6</b>
Reagent Preparation .....	6
Sample Preparation .....	7
Assay Procedure .....	7
<b>Data Analysis .....</b>	<b>9</b>
Calculation of Results .....	9
Performance Characteristics .....	10
<b>Resources .....</b>	<b>11</b>
References .....	11
Plate Layout .....	12

## **Introduction**

### **Intended Use**

For quantitative detection of human IL-1RA in tissue lysates or cell culture supernates. The significance of detection in sera, plasma, body fluids is undetermined.

### **Background**

The interleukin-1 receptor antagonist (IL-1RA) is a protein that in humans is encoded by the *IL1RN* gene.<sup>1,2</sup> The human IL1RN gene was localized to the long arm of chromosome 2 at band 2q14.2 by fluorescence in situ hybridization.<sup>3</sup> IL-1RA is a member of the interleukin 1 cytokine family. This protein inhibits the activities of interleukin 1, alpha (IL1A) and interleukin 1, beta (IL1B), and modulates a variety of interleukin 1 related immune and inflammatory responses.<sup>4</sup>

### **Principle of the Assay**

IL1RN (Human) ELISA Kit was based on standard sandwich enzyme-linked immune-sorbent assay technology. Human IL-1RA specific polyclonal antibodies were precoated onto 96-well plates. The human specific detection polyclonal antibodies were biotinylated. The test samples and biotinylated detection antibodies were added to the wells subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the human IL-1RA amount of sample captured in plate.

## General Information

### Materials Supplied

List of component

Component	Amount
Lyophilized recombinant Human IL-1RA standard	10 ng/tube x 2
One 96-well plate precoated with anti-Human IL-1RA antibody	1
Sample diluent buffer	30 ml
Biotinylated anti- Human IL-1RA antibody, dilution 1:100	130 µl
Antibody diluent buffer	12 ml
Avidin-Biotin-Peroxidase Complex (ABC), dilution 1:100	130 µl
ABC diluent buffer	12 ml
TMB color developing agent	10 ml
TMB stop solution	10 ml

### Storage Instruction

Store at 4°C for frequent use, at -20°C for infrequent use. Avoid multiple freeze-thaw cycles. Expiration: Four months at 4°C and eight months at -20°C.

### Materials Required but Not Supplied

- ✓ Microplate reader in standard size.
- ✓ Automated plate washer.
- ✓ Adjustable pipettes and pipette tips. Multichannel pipettes are recommended in the condition of large amount of samples in the detection.
- ✓ Clean tubes and Eppendorf tubes.
- ✓ Washing buffer (neutral PBS or TBS).

#### Preparation of 0.01M TBS:

Add 1.2g Tris, 8.5g NaCl; 450µl of purified acetic acid or 700µl of concentrated hydrochloric acid to 1000ml H<sub>2</sub>O and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

#### Preparation of 0.01 M PBS:

Add 8.5g sodium chloride, 1.4g Na<sub>2</sub>HPO<sub>4</sub> and 0.2g NaH<sub>2</sub>PO<sub>4</sub> to 1000ml distilled water and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

**Precautions for Use**

- Precautions
  - ✓ Before using Kit, spin tubes and bring down all components to bottom of tube.
  - ✓ Duplicate well assay was recommended for both standard and sample testing.
  - ✓ Don't let 96-well plate dry, dry plate will inactivate active components on plate.
  - ✓ In order to avoid marginal effect of plate incubation due to temperature difference (reaction may be stronger in the marginal wells), it is suggested that the diluted ABC and TMB solution will be pre-warmed in 37°C for 30 min before using.

## Assay Protocol

### Reagent Preparation

- Reconstitution of the human IL-1RA standard: IL-1RA standard solution should be prepared no more than 2 hours prior to the experiment. Two tubes of IL-1RA standard (10ng per tube) are included in each kit. Use one tube for each experiment.
  - a. 10,000 pg/ml of human IL-1RA standard solution: Add 1 ml sample diluent buffer into one tube, keep the tube at room temperature for 10 min and mix thoroughly.
  - b. 2000 pg/ml of human IL-1RA standard solution: Add 0.2 ml of the above 10 ng/ml IL-1RA standard solution into 0.8 ml sample diluent buffer and mix thoroughly.
  - c. 1000 pg/ml→31.2 pg/ml of human IL-1RA standard solutions: Label 6 Eppendorf tubes with 1000 pg/ml, 500 pg/ml, 250 pg/ml, 125 pg/ml, 62.5 pg/ml, 31.2 pg/ml, respectively. Aliquot 0.3 ml of the sample diluent buffer into each tube. Add 0.3 ml of the above 2000 pg/ml IL-1RA standard solution into 1st tube and mix. Transfer 0.3 ml from 1st tube to 2nd tube and mix. Transfer 0.3 ml from 2nd tube to 3rd tube and mix, and so on.

*Note: The standard solutions are best used within 2 hours. The 10 ng/ml standard solution may be stored at 4 °C for up to 12 hours, or at -20 °C for up to 48 hours. Avoid repeated freeze-thaw cycles.*

- Preparation of biotinylated anti-human IL-1RA antibody working solution: The solution should be prepared no more than 2 hours prior to the experiment.
  - a. The total volume should be: 0.1 ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume)
  - b. Biotinylated anti-human IL-1RA antibody should be diluted in 1:99 with the antibody diluent buffer and mixed thoroughly.
- Preparation of Avidin-Biotin-Peroxidase Complex (ABC) working solution: The solution should be prepared no more than 1 hour prior to the experiment.
  - a. The total volume should be: 0.1 ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume)
  - b. Avidin- Biotin-Peroxidase Complex (ABC) should be diluted in 1:99 with the ABC dilution buffer and mixed thoroughly.

### **Sample Preparation**

Store samples to be assayed within 24 hours at 2-8°C. For long-term storage, aliquot and freeze samples at -20°C. Avoid repeated freeze-thaw cycles.

- Cell culture supernate, tissue lysate or body fluids: Remove particulates by centrifugation, analyze immediately or aliquot and store at -20°C
- Serum: Allow the serum to clot in a serum separator tube (about 30 min) at room temperature. Centrifuge at approximately 1000 X g for 15 min. Analyze the serum immediately or aliquot and store frozen at -20°C.
- Plasma: Collect plasma using EDTA as an anticoagulant. Centrifuge for 15 min at 2-8°C at 1500 x g within 30 min of collection. For eliminating the platelet effect, suggesting that further centrifugation for 10 min at 2-8°C at 10000 x g. Analyze immediately or aliquot and store frozen at -20°C.

### **Sample Dilution Guideline**

The user needs to estimate the concentration of the target protein in the sample and select a proper dilution factor so that the diluted target protein concentration falls near the middle of the linear regime in the standard curve. Dilute the sample using the provided diluent buffer. The following is a guideline for sample dilution. Several trials may be necessary in practice. The sample must be well mixed with the diluents buffer.

- High target protein concentration (20-200 ng/ml). The working dilution is 1:100. i.e. Add 3 µl sample into 297 µl sample diluent buffer.
- Medium target protein concentration (2-20 ng/ml). The working dilution is 1:10. i.e. Add 25 µl sample into 225 µl sample diluent buffer.
- Low target protein concentration (31.2-2000 pg/ml). The working dilution is 1:2. i.e. Add 100 µl sample to 100 µl sample diluent buffer.
- Very Low target protein concentration ( $\leq 31.2$  pg/ml). No dilution necessary, or the working dilution is 1:2.

### **Assay Procedure**

The ABC working solution and TMB color developing agent must be kept warm at 37°C for 30 min before use. When diluting samples and reagents, they must be mixed completely and evenly. Standard IL-1RA detection curve should be prepared for each experiment. The user will decide sample dilution fold by crude estimation of IL-1RA amount in samples.

1. Aliquot 0.1 ml per well of the 2000 pg/ml, 1000 pg/ml, 500 pg/ml, 250 pg/ml, 125 pg/ml, 62.5 pg/ml, 31.2 pg/ml human IL-1RA standard solutions into the precoated 96-well plate. Add 0.1 ml of the sample diluent buffer into the control well (Zero well). Add 0.1 ml of each properly diluted sample of human sera, plasma, body fluids, tissue lysates or cell culture supernatants to each empty well. See "Sample Dilution Guideline" above for details. We recommend that each human IL-1RA standard solution and each sample is measured in duplicate.

2. Seal the plate with the cover and incubate at 37°C for 90 min.
  3. Remove the cover, discard plate content, and blot the plate onto paper towels or other absorbent material. Do NOT let the wells completely dry at any time.
  4. Add 0.1 ml of biotinylated anti-human IL-1RA antibody working solution into each well and incubate the plate at 37°C for 60 min.
  5. Wash the plate three times with 0.01 M TBS or 0.01 M PBS, and each time let washing buffer stay in the wells for 1 min. (Plate Washing Method: Discard the solution in the plate without touching the side walls. Blot the plate onto paper towels or other absorbent material. Soak each well with at least 0.3 ml PBS or TBS buffer for 1~2 minutes. Repeat this process two additional times for a total of THREE washes. *Note: For automated washing, aspirate all wells and wash THREE times with PBS or TBS buffer, overfilling wells with PBS or TBS buffer. Blot the plate onto paper towels or other absorbent material.*)
  6. Add 0.1 ml of prepared ABC working solution into each well and incubate the plate at 37°C for 30 min.
  7. Wash plate 5 times with 0.01 M TBS or 0.01 M PBS, and each time let washing buffer stay in the wells for 1-2 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material.
  8. Add 90 µl of prepared TMB color developing agent into each well and incubate plate at 37°C in dark for 25-30 min (*Note: For reference only, the optimal incubation time should be determined by end user. And the shades of blue can be seen in the wells with the four most concentrated human IL-1RA standard solutions; the other wells show no obvious color*).
  9. Add 0.1 ml of prepared TMB stop solution into each well. The color changes into yellow immediately.
  10. Read the O.D. absorbance at 450nm in a microplate reader within 30 min after adding the stop solution.
- Summary
    1. Add samples and standards and incubate the plate at 37°C for 90 min. Do not wash.
    2. Add biotinylated antibodies and incubate the plate at 37°C for 60 min. Wash plate 3 times with 0.01M TBS.
    3. Add ABC working solution and incubate the plate at 37°C for 30 min. Wash plate 5 times with 0.01M TBS.
    4. Add TMB color developing agent and incubate the plate at 37°C in dark for 25-30 min.
    5. Add TMB stop solution and read.

## Data Analysis

### Calculation of Results

For calculation, (the relative O.D.<sub>450</sub>) = (the O.D.<sub>450</sub> of each well) – (the O.D.<sub>450</sub> of Zero well). The standard curve can be plotted as the relative O.D.<sub>450</sub> of each standard solution (Y) vs. the respective concentration of the standard solution (X). The human IL-1RA concentration of the samples can be interpolated from the standard curve.

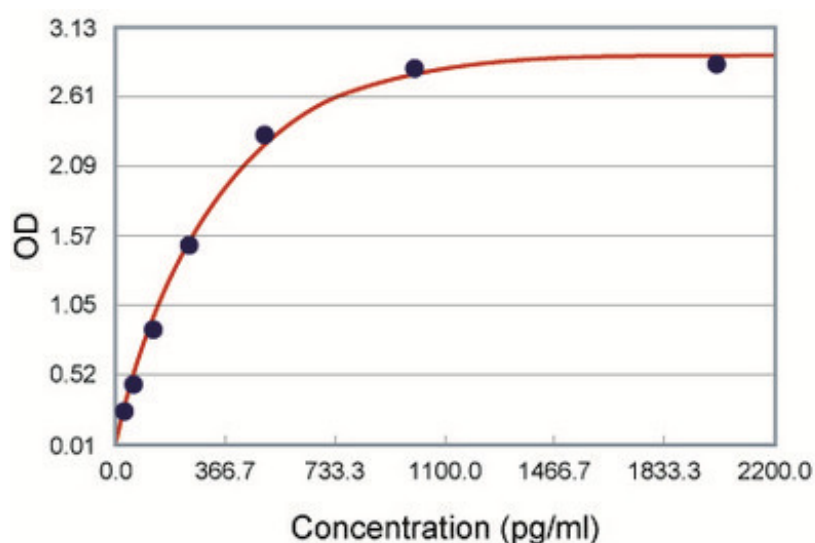
*Note: if the samples measured were diluted, multiply the dilution factor to the concentrations from interpolation to obtain the concentration before dilution.*

### Typical result

Typical Data Obtained from Human IL-1RA

(TMB reaction incubate at 37°C for 25 min)

Concentration pg/ml	0.0	31.2	62.5	125	250	500	1000	2000
O.D.	0.028	0.244	0.456	0.861	1.497	2.323	2.804	2.852



This standard curve was generated for demonstration purpose only. A standard curve must be run with each assay.

**Performance Characteristics**

- Range  
31.2 pg/ml-2000 pg/ml
- Sensitivity  
< 2 pg/ml
- Specificity  
No detectable cross-reactivity with any other protein.

## Resources

### References

1. Steinkasserer A, Spurr NK, Cox S, Jeggo P, Sim RB (July 1992). "The human IL-1 receptor antagonist gene (IL1RN) maps to chromosome 2q14-q21, in the region of the IL-1 alpha and IL-1 beta loci". Genomics 13 (3): 654–7.
2. Patterson D, Jones C, Hart I, Bleskan J, Berger R, Geyer D, Eisenberg SP, Smith MF, Arend WP (January 1993). "The human interleukin-1 receptor antagonist (IL1RN) gene is located in the chromosome 2q14 region". Genomics 15 (1): 173–6.
3. Patterson, D., Jones, C., Hart, I., Bleskan, J., Berger, R., Geyer, D., Eisenberg, S. P., Smith, M. F., Jr., Arend, W. P. The human interleukin-1 receptor antagonist (IL1RN) gene is located in the chromosome 2q14 region. Genomics 15: 173-176, 1993.
4. "Entrez Gene: IL1RN interleukin 1 receptor antagonist"

**Plate Layout**

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