

Human IL-1 RL1 ELISA Kit (ST2) ab213796

1 Image

Overview

Product name Human IL-1 RL1 ELISA Kit (ST2)

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	1	16	0.53ng/ml	0.028	5.2%
	2	16	2.05ng/ml	0.094	4.6%
	3	16	4.31ng/ml	0.142	3.3%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	1	24	0.69ng/ml	0.06	8.7%
	2	24	2.54ng/ml	0.157	6.2%
	3	24	4.68ng/ml	0.229	4.9%

Sample type Cell culture supernatant, Serum, Cell Lysate, Hep Plasma, EDTA Plasma, Tissue Homogenate

Assay type Sandwich (quantitative)

Sensitivity 10 pg/ml

Range 156 pg/ml - 10000 pg/ml

Assay time 3h 30m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview The Human IL-1 RL1 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ST2) (ab213796) is designed for the quantitative detection of Human IL1RL1 in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immunosorbent assay technology. A

monoclonal antibody from mouse specific for IL-1 RL1 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: K19-S328) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL-1 RL1 is added 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 IL1RL1 amount of sample captured in plate.

## Notes

Interleukin 1 receptor-like 1, also known as IL-1 RL1 or ST2, is a protein that in humans is encoded by the IL1RL1 gene. The protein encoded by this gene is a member of the interleukin 1 receptor family. Studies of the similar gene in mouse suggested that this receptor can be induced by pro-inflammatory stimuli and may be involved in the function of helper T cells. TL1RL1 is necessary for endotoxin tolerance and, by inhibiting TLR responses, enhances Th2 responses. This gene, interleukin 1 receptor, type I (IL-1 RL1), interleukin 1 receptor, type II (IL-1 R2) and interleukin 1 receptor like 2 (IL-1 RL2) form a cytokine receptor gene cluster in a region mapped to chromosome 2q12.

## Platform

Pre-coated microplate (12 x 8 well strips)

## Properties

### Storage instructions

Store at -20°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	1 x 96 tests
ABC Diluent Buffer	Blue Cap	1 x 12ml	1 x 12ml
Adhesive Plate Seal		4 units	4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml	1 x 12ml
Anti-Human IL1RL1 coated Microplate (12 x 8 wells)		1 unit	1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 100µl	1 x 100µl
Biotinylated anti-Human IL1RL1 antibody		1 x 100µl	1 x 100µl
Lyophilized recombinant Human IL1RL1 standard		2 vials	2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml	1 x 10ml
Wash Buffer (25X)		1 x 20ml	1 x 20ml

## Function

Receptor for interleukin-33 (IL-33), its stimulation recruits MYD88, IRAK1, IRAK4, and TRAF6, followed by phosphorylation of MAPK3/ERK1 and/or MAPK1/ERK2, MAPK14, and MAPK8. Possibly involved in helper T-cell function.

## Tissue specificity

Highly expressed in kidney, lung, placenta, stomach, skeletal muscle, colon and small intestine.

Isoform A is prevalently expressed in the lung, testis, placenta, stomach and colon. Isoform B is more abundant in the brain, kidney and the liver. Isoform C is not detected in brain, heart, liver, kidney and skeletal muscle.

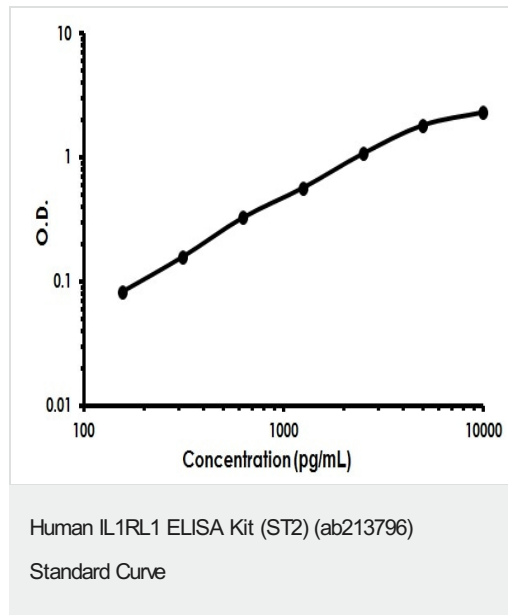
#### Sequence similarities

Belongs to the interleukin-1 receptor family.  
Contains 3 Ig-like C2-type (immunoglobulin-like) domains.  
Contains 1 TIR domain.

#### Cellular localization

Secreted and Cell membrane.

#### Images



Human IL1RL1 ELISA Kit (ST2) (ab213796) Standard Curve.

**Please note:** All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

#### Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.com/abpromise> or contact our technical team.

#### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors