

## Product datasheet

# Mouse IL-1 RL1 ELISA Kit (ST2) ab213871

1 Image

### Overview

**Product name** Mouse IL-1 RL1 ELISA Kit (ST2)

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
1	16	0.57ng/ml	0.029	5.1%
2	16	2.33ng/ml	0.086	3.7%
3	16	4.25ng/ml	0.136	3.2%

Inter-assay

Sample	n	Mean	SD	CV%
1	24	0.68ng/ml	0.058	8.6%
2	24	2.74ng/ml	0.17	6.2%
3	24	4.61ng/ml	0.244	5.3%

**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:** Mouse

**Product overview**

The Mouse IL1RL1 Enzyme-Linked Immunosorbent Assay (ELISA) kit (ab213871) is designed for the quantitative measurement of Mouse IL1RL1 in cell culture supernatants, cell lysates, tissue homogenates, serum and plasma (heparin, EDTA).

The ELISA kit is based on standard sandwich enzyme-linked immune-sorbent assay technology.

A monoclonal antibody from rat specific for IL1RL1 has been pre-coated onto 96-well plates. Standards (Expression system for standard: NSO; Immunogen sequence: S27-R332) and test samples are added to the wells, a biotinylated detection polyclonal antibody from goat specific for IL1RL1 is added subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex is added and unbound conjugates are washed away with PBS or TBS buffer. HRP substrate TMB is used to visualize HRP enzymatic reaction. TMB is 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 Mouse IL1RL1 amount of sample captured in plate.

**Notes**

Interleukin 1 receptor-like 1, also known as IL1RL1 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 proinflammatory stimuli, and may be involved in the function of helper T cells. IL1RL1 is necessary for endotoxin tolerance and, by inhibiting TLR responses, enhances Th2 responses. This gene, interleukin 1 receptor, type I (IL1R1), interleukin 1 receptor, type II (IL1R2) and interleukin 1 receptor-like 2 (IL1RL2) 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
ABC Diluent Buffer	Blue Cap	1 x 12ml
Adhesive Plate Seal		4 units
Antibody Diluent Buffer	Green Cap	1 x 12ml
Anti-Mouse IL1RL1 coated Microplate (12 x 8 wells)		1 unit
Avidin-Biotin-Peroxidase Complex (ABC)		1 x 130µl
Biotinylated anti- Mouse IL1RL1 antibody		1 x 130µl
Lyophilized recombinant Mouse IL1RL1 standard		2 vials
Sample Diluent Buffer	Green Cap	1 x 30ml
TMB Color Developing Agent	Black Cap	1 x 10ml
TMB Stop Solution	Yellow Cap	1 x 10ml

**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.

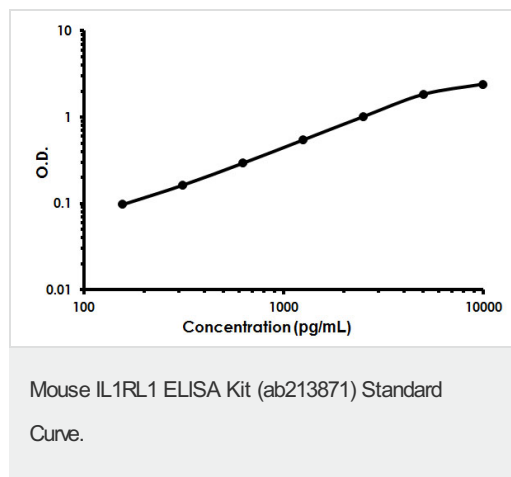
## Sequence similarities

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



Mouse IL1RL1 ELISA Kit (ab213871) 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