

## Product datasheet

# Human IL-1R3 ELISA Kit ab267659

[1 Image](#)

### Overview

**Product name** Human IL-1R3 ELISA Kit

**Detection method** Colorimetric

**Precision**

Intra-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 10%

Inter-assay

Sample	n	Mean	SD	CV%
--------	---	------	----	-----

Overall < 12%

**Sample type** Serum, Plasma, Cell culture media

**Assay type** Sandwich (quantitative)

**Sensitivity** 18 pg/ml

**Range** 16.38 pg/ml - 4000 pg/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
-------------	-----------	-------

Serum 106 91% - 129%

Plasma 113 102% - 132%

Cell culture media 105 92% - 111%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Human

**Product overview** Human IL-1R3 ELISA Kit is designed for the quantitative determination of IL-1R3 in cell culture supernatants, plasma and serum samples.

This assay employs an antibody specific for human IL-1R3 coated on a 96-well plate. Standards

and samples are pipetted into the wells and IL-1R3 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-human IL-1R3 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IL-1R3 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Microplate

## Properties

**Storage instructions** Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
200X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent	1 x 15ml
Anti-Human IL-1R3 coated Microplate (12 x 8 wells)	1 unit
Biotinylated Anti-Human IL-1R3 Detection Antibody	2 vials
Human IL-1R3 Standard (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB Substrate Solution	1 x 12ml

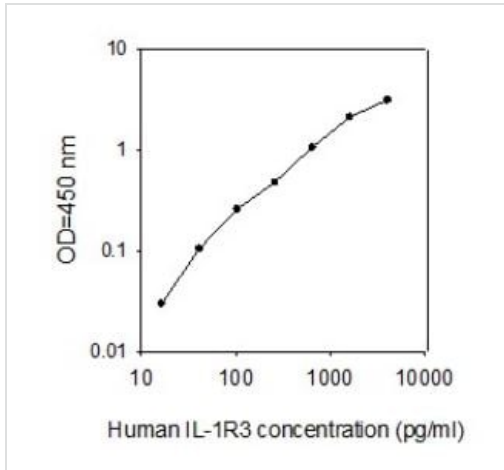
**Function** Coreceptor with IL1R1. Associates with IL1R1 bound to IL1B to form the high affinity interleukin-1 receptor complex which mediates interleukin-1-dependent activation of NF-kappa-B and other pathways. Signaling involves the recruitment of adapter molecules such as TOLLIP, MYD88, and IRAK1 or IRAK2 via the respective TIR domains of the receptor/coreceptor subunits. Recruits TOLLIP to the signaling complex. Does not bind to interleukin-1 alone; binding of IL1RN to IL1R1, prevents its association with IL1R1 to form a signaling complex. The cellular response is modulated through a non-signaling association with the membrane IL1R2 decoy receptor. Secreted forms (isoforms 2 and 3) associate with secreted ligand-bound IL1R2 and increase the affinity of secreted IL1R2 for IL1B; this complex formation may be the dominant mechanism for neutralization of IL1B by secreted/soluble receptors.

**Tissue specificity** Detected in liver, skin, placenta, thymus and lung.

**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 IL-1R3 ELISA kit (ab267659) Standard curve.

Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

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