

### Rabbit IL-13 ELISA Kit ab277391

[1 Image](#)

#### Overview

**Product name** Rabbit IL-13 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** Cell culture supernatant, Serum, Plasma

**Assay type** Sandwich (quantitative)

**Sensitivity** 0.042 ng/ml

**Range** 0.041 ng/ml - 10 ng/ml

**Recovery**

Sample specific recovery

Sample type	Average %	Range
Serum	89.69	81% - 97%
Plasma	75.91	71% - 80%
Cell culture media	95.33	87% - 102%

**Assay duration** Multiple steps standard assay

**Species reactivity** **Reacts with:** Rabbit

**Product overview** Rabbit IL-13 ELISA Kit (ab277391) is an in-vitro enzyme-linked immunosorbent assay for the quantitative measurement of Rabbit IL-13 in serum (Rabbit IL-13 concentration is low in normal serum/plasma, and may not be detectable in this assay), plasma and cell culture supernatants.

This assay employs an antibody specific for Rabbit IL-13 coated on a 96-well plate. Standards

and samples are pipetted into the wells and Rabbit IL-13 present in a sample is bound to the wells by the immobilized antibody. The wells are washed, and biotinylated anti-Rabbit IL-13 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 Rabbit IL-13 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

**Platform** Pre-coated microplate (12 x 8 well strips)

## 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 B	1 x 15ml
Assay Diluent C	1 x 30ml
Biotinylated Anti-Rabbit IL-13 Antibody	2 vials
Anti-Rabbit IL-13 coated Microplate	1 unit
Rabbit IL-13 standard protein (Lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

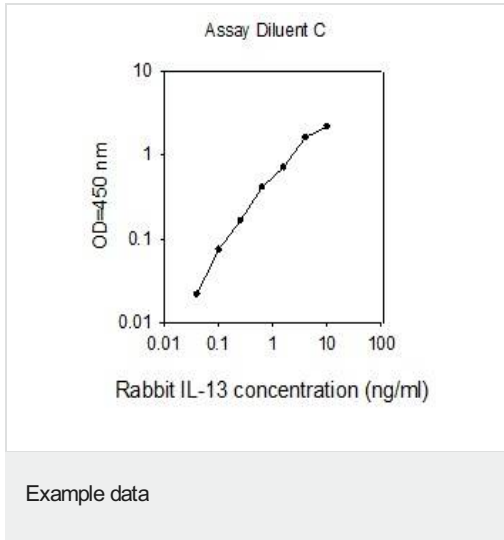
**Function** Cytokine (PubMed:8096327, PubMed:8097324). Inhibits inflammatory cytokine production (PubMed:8096327). Synergizes with IL2 in regulating interferon-gamma synthesis (PubMed:8096327). May be critical in regulating inflammatory and immune responses (PubMed:8096327, PubMed:8097324). Positively regulates IL31RA expression in macrophages.

**Involvement in disease** Allergic rhinitis

**Sequence similarities** Belongs to the IL-4/IL-13 family.

**Cellular localization** Secreted.

## Images



These standard curves are for demonstration only. A standard curve must be run with each assay.

**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