

Human IL-13 ELISA Kit ab100553

3 References 2 Images

Overview

Product name	Human IL-13 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 0.15 pg/ml
Range	0.16 pg/ml - 40 pg/ml
Recovery	100 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	100.65	88% - 107%
Serum	96.35	84% - 103%
Plasma	94.39	82% - 102%

Assay duration Multiple steps standard assay

Species reactivity Reacts with: Human

Product overview Abcam's IL-13 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Human IL-13 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for Human IL-13 coated on a 96-well plate. Standards and samples are pipetted into the wells and IL-13 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human 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 IL-13 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Human IL-13 ELISA Kit ([ab178014](#)) from our SimpleStep ELISA® range.

<b>Notes</b>	Optimization may be required with urine samples
<b>Platform</b>	Microplate

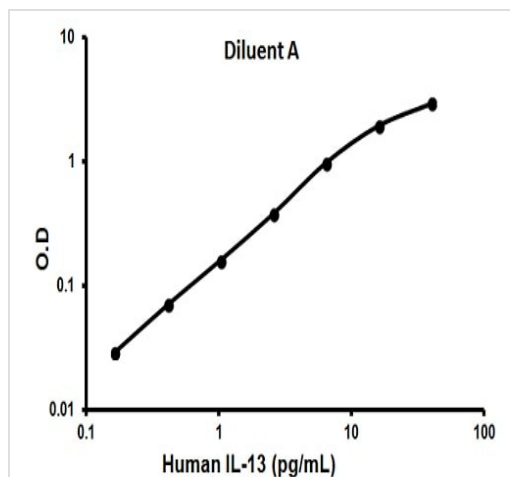
## Properties

<b>Storage instructions</b>	Store at -20°C. Please refer to protocols.
-----------------------------	--

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
120X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Human IL-13	2 vials
IL-13 Microplate (12 x 8 wells)	1 unit
Recombinant Human IL-13 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

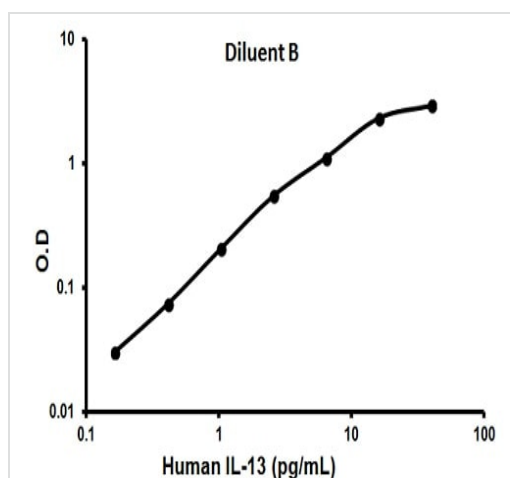
<b>Function</b>	Cytokine (PubMed:8096327, PubMed:8097324). Inhibits inflammatory cytokine production (PubMed:8096327). Synergizes with IL2 in regulating interferon-gamma synthesis (PubMed:8096327). May be critical in regulating inflammatory and immune responses (PubMed:8096327, PubMed:8097324). Positively regulates IL31RA expression in macrophages.
<b>Involvement in disease</b>	Allergic rhinitis
<b>Sequence similarities</b>	Belongs to the IL-4/IL-13 family.
<b>Cellular localization</b>	Secreted.

## Images



Representative standard curve using ab100553

Typical standard curve



Representative standard curve using ab100553

Typical 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