

Human IL-13 Receptor alpha ELISA Kit ab46066

1 Image

Overview

Product name Human IL-13 Receptor alpha ELISA Kit

Detection method Colorimetric

Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	A	14	3024pg/ml	93.67	3.1%
	B	14	1054pg/ml	29.69	2.82%

	Inter-assay				
	Sample	n	Mean	SD	CV%
	A	8	2744pg/ml	90.1	3.28%
	B	8	725pg/ml	43.9	6.06%

Sample type Cell culture supernatant, Serum, Plasma

Assay type Sandwich (quantitative)

Sensitivity < 150 pg/ml

Range 156 pg/ml - 5000 pg/ml

Recovery	Sample specific recovery		
	Sample type	Average %	Range
	Serum	90	156pg/ml - 5000pg/ml

Assay time 3h 45m

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Human

Product overview Abcam's IL-13 Receptor alpha Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of Interleukin 13 Receptor (IL-13 Receptor alpha, also known as CD213 alpha 2) in Human serum, plasma, buffered solutions and cell culture medium.

A monoclonal antibody specific for IL-13 Receptor alpha has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-13 Receptor alpha concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-13 Receptor alpha are simultaneously incubated. After washing, the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-13 Receptor alpha present in the samples.

This kit will recognize both endogenous and recombinant Human IL-13 Receptor alpha.

## Platform

Microplate

## Properties

### Storage instructions

Store at +4°C. Please refer to protocols.

Components	Identifier	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 15ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7ml	1 x 13ml
Biotinylated anti-IL-13 Receptor alpha	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
HRP Diluent	Red	1 x 12ml	1 x 23ml
IL-13 Receptor alpha Microplate (12 x 8 well strips)		1 unit	2 units
IL-13 Receptor alpha Standard (Lyophilized)	Yellow	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

## Function

Binds as a monomer with high affinity to interleukin-13 (IL13), but not to interleukin-4 (IL4).

## Sequence similarities

Belongs to the type I cytokine receptor family. Type 5 subfamily.

Contains 1 fibronectin type-III domain.

## Domain

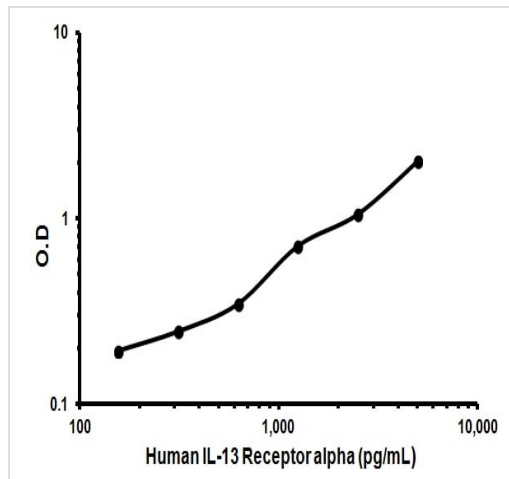
The WSXWS motif appears to be necessary for proper protein folding and thereby efficient intracellular transport and cell-surface receptor binding.

The box 1 motif is required for JAK interaction and/or activation.

## Cellular localization

Membrane.

## Images



Representative Standard Curve using ab46066

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