

Human IL-13 Receptor alpha ELISA Kit ab46066

1 图像

概述

产品名称 人IL-13 Receptor alpha ELISA试剂盒

检测方法 Colorimetric

批次内				
样品	n	Mean	SD	CV%
A	14	3024pg/ml	93.67	3.1%
B	14	1054pg/ml	29.69	2.82%

批次间				
样品	n	Mean	SD	CV%
A	8	2744pg/ml	90.1	3.28%
B	8	725pg/ml	43.9	6.06%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 150 pg/ml

范围 156 pg/ml - 5000 pg/ml

特定样本回收率		
样品类型	平均%	范围
Serum	90	156pg/ml - 5000pg/ml

检测时间 3h 45m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 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.

**平台** Microplate

## 性能

**存放说明** Store at +4°C. Please refer to protocols.

组件	标识符	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

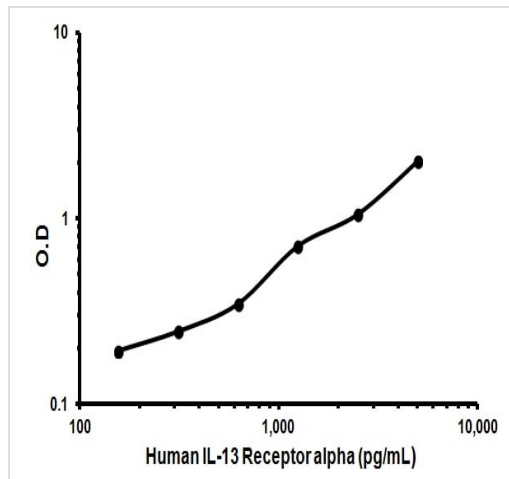
**功能** Binds as a monomer with high affinity to interleukin-13 (IL13), but not to interleukin-4 (IL4).

**序列相似性** Belongs to the type I cytokine receptor family. Type 5 subfamily.  
Contains 1 fibronectin type-III domain.

**结构域** 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.

**细胞定位** Membrane.

## 图片



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.cn/abpromise> or contact our technical team.

### Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors