

### Human IL-13 ELISA Set (without plates) ab47353

[1 References](#) [1 图像](#)

#### 概述

产品名称	人IL-13 ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 1.5 pg/ml
范围	3.125 pg/ml - 100 pg/ml
检测时间	2h 15m
实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	<p>Abcam's IL-13 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-13 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure human IL-13 in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>The IL-13 ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>Recognizes both natural and recombinant human IL-13.</p>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

#### 性能

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

组件	标识符	10 x 96 tests	5 x 96 tests	15 x 96 tests
IL-13 standard		—	—	

组件	标识符	10 x 96 tests	5 x 96 tests	15 x 96 tests
Chromogen TMB Substrate Solution		4 vials	2 vials	6 vials
Detection biotinylated anti-IL-13 antibody		2 x 0.55ml	1 x 0.55ml	3 x 0.55ml
IL-13 Capture antibody		2 x 0.5ml	1 x 0.5ml	3 x 0.5ml
5 vials				
Streptavidin-HRP		2 vials	1 vial	3 vials

**相关性** Human IL-13 was originally identified by differential screening of an anti CD28 activated human peripheral blood mononuclear cell cDNA library as an induction specific novel cytokine. It was also isolated from cDNA libraries of human T cell clones using the murine IL-13 (P600) cDNA as a probe. Human IL-13, a pleiotropic cytokine, is produced by activated Th0, Th1 like, Th2 like, and CD8 T cells. The gene for human IL-13 maps to chromosome 5 and is closely linked to the genes for IL3, IL4, IL5, and GMCSF. IL-13 inhibits proinflammatory cytokine production and stimulates antibody production. It induces proliferation in the human pre myeloid cell line TF1. IL-13 has multiple effects on the differentiation and functions of monocytes and macrophages. It suppresses cytotoxic functions and induces changes in the morphology of human monocytes and in the phenotype of human monocytes and B cells by upregulating MHC class II expression. IL-13 will also decrease the production of nitric oxide by activated murine macrophages, leading to impaired parasitocidal activity. Human and mouse interleukin 13 share approximately 58% amino acid sequence identity. Although human and mouse IL-13 are equally active on human cells, human IL-13 is much less active than mouse IL-13 on mouse cells. Human IL-13 and human IL-4 also share approximately 30% sequence homology and have similar biological functions.

**细胞定位** Secreted

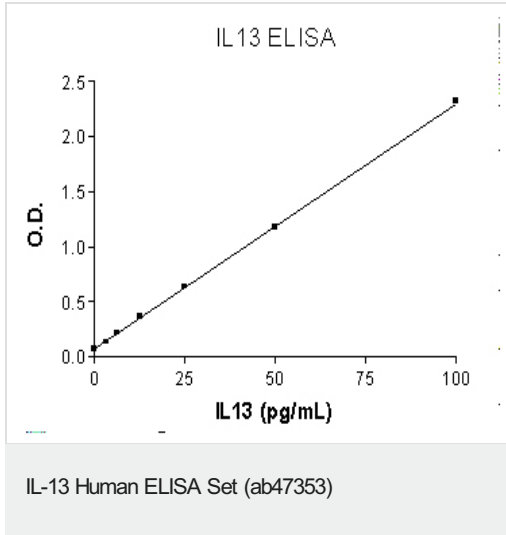
**应用**

**The Abpromise guarantee** **Abpromise™**承诺保证使用ab47353于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
<b>Sandwich ELISA</b>		Use at an assay dependent concentration.

**图片**



IL-13 Standard curve ranging from 3.12 to 100 pg/mL

**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