

Human IL-27 ELISA Kit ab83695

★★★★★ [1 Abreviews](#) [2 References](#) [1 图像](#)

概述

产品名称	人IL-27 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	A	9	343pg/ml	12.9	4.8%
	B	9	127pg/ml	8.1	5.1%
	C	9	427pg/ml	15.8	3.8%
	批次间				
	样品	n	Mean	SD	CV%
	A	9	343pg/ml	29	8.5%
	B	9	157pg/ml	17	10.8%
	C	9	427pg/ml	38	9%
样品类型	Cell culture supernatant, Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 12.8 pg/ml				
范围	31.25 pg/ml - 1000 pg/ml				
回收率	95 %				
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-27 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-27 in Human serum, plasma, buffered solutions and cell culture medium.				

A monoclonal antibody specific IL-27 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-27 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-27 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-27 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-27-A and IL-27-B.

平台 Microplate

## 性能

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

组件	标识符	2 x 96 tests	1 x 96 tests
1X Standard Diluent Buffer		2 x 25ml	1 x 25ml
200X Wash Buffer	White	2 x 10ml	1 x 10ml
Biotinylated anti IL-27		2 x 400µl	1 x 400µl
Biotinylated Antibody Diluent	Red	1 x 13ml	1 x 7ml
Chromogen TMB Substrate Solution		1 x 24ml	1 x 11ml
HRP Diluent	Red	1 x 23ml	1 x 12ml
IL-27 Microplate (12 x 8 well strips)		2 units	1 unit
IL-27 Standard (Lyophilized)		4 vials	2 vials
Plastic Plate Covers		4 units	2 units
Stop Reagent	Black	2 x 11ml	1 x 11ml
Streptavidin-HRP		4 x 5µl	2 x 5µl

功能 Cytokine with pro- and anti-inflammatory properties, that can regulate T helper cell development, suppress T-cell proliferation, stimulate cytotoxic T cell activity, induce isotype switching in B-cells, and that has diverse effects on innate immune cells. Among its target cells are CD4 T helper cells which can differentiate in type 1 effector cells (TH1), type 2 effector cells (TH2) and IL17 producing helper T-cells (TH17). It drives rapid clonal expansion of naive but not memory CD4 T-cells. It also strongly synergizes with IL-12 to trigger interferon-gamma/IFN-gamma production of naive CD4 T-cells, binds to the cytokine receptor WSX-1/TCCR which appears to be required but not sufficient for IL-27-mediated signal transduction. IL-27 potentiate the early phase of TH1 response and suppress TH2 and TH17 differentiation. It induces the differentiation of TH1 cells via two distinct pathways, p38 MAPK/TBX21- and ICAM1/ITGAL/ERK-dependent pathways. It also induces STAT1, STAT3, STAT4 and STAT5 phosphorylation and activates TBX21/T-Bet via STAT1 with

resulting IL12RB2 up-regulation, an event crucial to TH1 cell commitment. It suppresses the expression of GATA3, the inhibitor TH1 cells development. In CD8 T-cells, it activates STATs as well as GZMB. IL-27 reveals to be a potent inhibitor of TH17 cell development and of IL-17 production. Indeed IL-27 subunit p28 alone is also able to inhibit the production of IL17 by CD4 and CD8 T-cells. While IL-27 suppressed the development of proinflammatory Th17 cells via STAT1, it inhibits the development of anti-inflammatory inducible regulatory T-cells, iTreg, independently of STAT1. IL-27 has also an effect on cytokine production, it suppresses proinflammatory cytokine production such as IL2, IL4, IL5 and IL6 and activates suppressors of cytokine signaling such as SOCS1 and SOCS3. Apart from suppression of cytokine production, IL-27 also antagonizes the effects of some cytokines such as IL6 through direct effects on T cells. Another important role of IL-27 is its antitumor activity as well as its antiangiogenic activity with activation of production of antiangiogenic chemokines such as IP-10/CXCL10 and MIG/CXCL9. In vein endothelial cells, it induces IRF1/interferon regulatory factor 1 and increase the expression of MHC class II transactivator/CIITA with resulting up-regulation of major histocompatibility complex class II. IL-27 also demonstrates antiviral activity with inhibitory properties on HIV-1 replivation.

#### 组织特异性

Expressed in monocytes and in placenta.

#### 序列相似性

Belongs to the IL-6 superfamily.

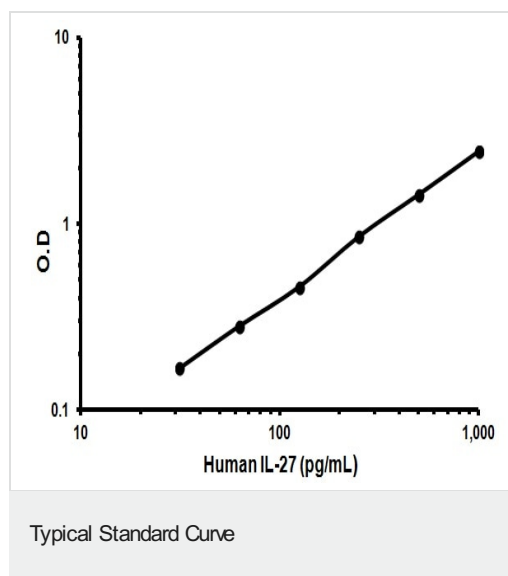
#### 翻译后修饰

O-glycosylated.

#### 细胞定位

Secreted. Does not seem to be secreted without coexpression of EBI3.

#### 图片



Representative Standard Curve using ab83695

**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