

Human IL-10 ELISA Kit ab46034

★★★★★ [1 Abreviews](#) [24 References](#) [2 图像](#)

概述

产品名称	人IL-10 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	sample	6			3.2%
	批次间				
	样品	n	Mean	SD	CV%
	sample	6			7.3%
样品类型	Cell culture supernatant, Serum, Plasma				
检测类型	Sandwich (quantitative)				
灵敏度	< 5 pg/ml				
范围	12.5 pg/ml - 400 pg/ml				
回收率	111 %				
检测时间	2h 40m				
实验步骤	Multiple steps standard assay				
种属反应性	与反应: Human				
产品概述	Abcam's IL-10 Human <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 in supernatants, buffered solutions, serum and plasma samples.				

A monoclonal antibody specific for IL-10 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-10 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-10 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-10 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-10.

Get results in 90 minutes with Human IL-10 ELISA Kit ([ab185986](#)) from our SimpleStep ELISA<sup>®</sup> range.

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

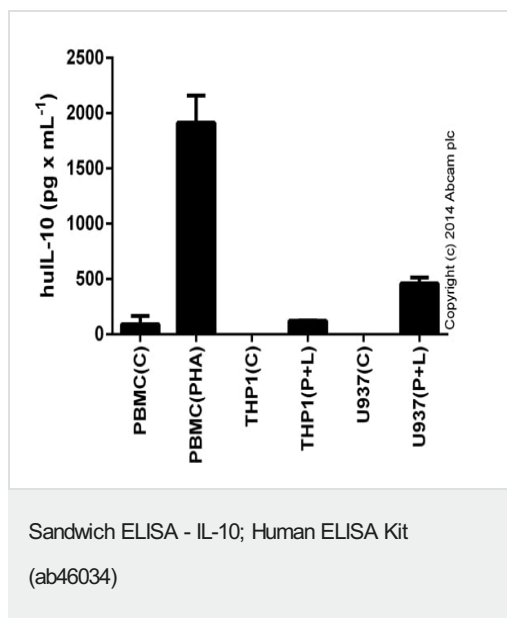
功能 Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

组织特异性 Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types.

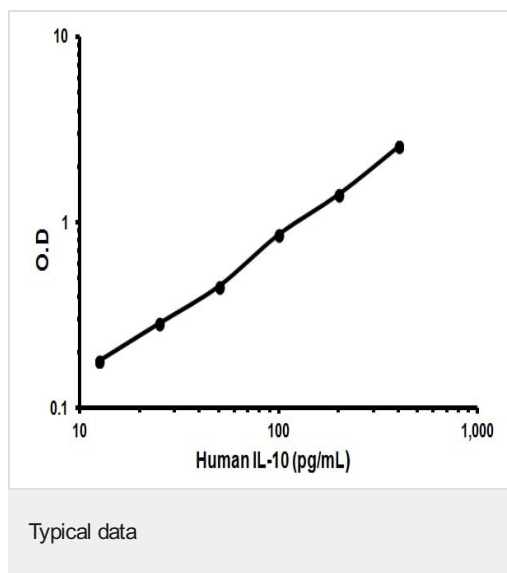
序列相似性 Belongs to the IL-10 family.

细胞定位 Secreted.

## 图片



IL-10 detected in supernatants from control cells (C) or cells stimulated for 24 hours with 50 ng x mL<sup>-1</sup> of PMA ([ab120297](#)), with the addition of 1 ug x mL<sup>-1</sup> of LPS (Sigma) (P/L) for the last 6 hours. PBMCs were stimulated for 48 hours with 2 % PHA-M (LifeTechnologies). Results shown after background signal was subtracted (duplicates +/- SD).



Representative standard curve using ab46034

**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