

Human IL-17a ELISA Kit ab83688

[6 References](#) [2 图像](#)

概述

产品名称 人IL-17a ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall	6			3.3%

批次间

样品	n	Mean	SD	CV%
Overall	6			5.2%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 3 pg/ml

范围 3.125 pg/ml - 100 pg/ml

回收率 105 %

特定样本回收率

样品类型	平均%	范围
Serum	121	113% - 130%

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-17a Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative of IL-17a in supernatants, buffered solutions, serum, plasma and other body fluids.

A monoclonal antibody specific for IL-17a has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-17a 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-17a 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-17a present in the samples.

This kit will recognize both endogenous and recombinant Human IL-17a.

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-IL17a	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-17a Microplate (12 x 8 well strips)		1 unit	2 units
IL-17a Standard (Lyophilized)	Orange	2 vials	4 vials
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

功能 Induces stromal cells to produce proinflammatory and hematopoietic cytokines. Enhances the surface expression of the intracellular adhesion molecule-1 (ICAM-1) in fibroblasts.

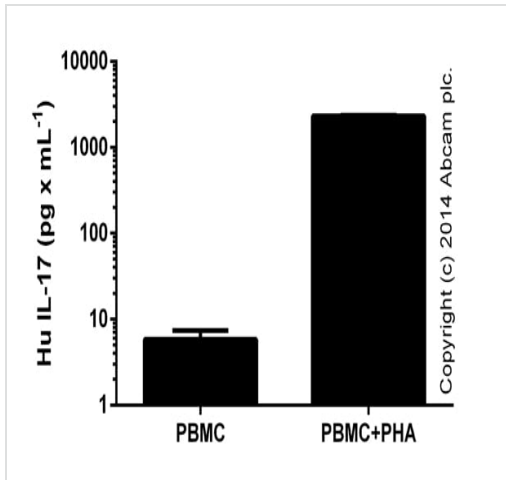
组织特异性 Restricted to activated memory T-cells.

序列相似性 Belongs to the IL-17 family.

翻译后修饰 Found both in glycosylated and nonglycosylated forms.

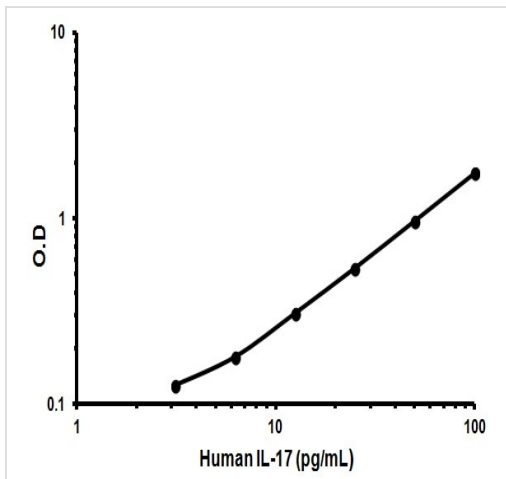
细胞定位 Secreted.

图片



IL-17 measured in cell culture supernatants (dilution range 1/1-1/100) from control (PBMC) or treated ((PBMC+PHA) 48 hours; 2% PHA-M; LifeTechnologies) cells (duplicates +/- SD).

ELISA: IL-17 Human ELISA Kit (ab83688)



Representative Standard Curve using ab83688

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