

Human IL-17α ELISA Kit ab83688

6 References 2 Images

Overview

Product name	Human IL-17a ELISA Kit				
Detection method	Colorimetric				
Precision	Intra-assay				
	Sample	n	Mean	SD	CV%
	Overall	6			3.3%
	Inter-assay				
	Sample	n	Mean	SD	CV%
	Overall	6			5.2%
Sample type	Cell culture supernatant, Serum, Plasma				
Assay type	Sandwich (quantitative)				
Sensitivity	< 3 pg/ml				
Range	3.125 pg/ml - 100 pg/ml				
Recovery	105 %				
	Sample specific recovery				
	Sample type	Average %		Range	
	Serum	121		113% - 130%	
Assay time	3h 45m				
Assay duration	Multiple steps standard assay				
Species reactivity	Reacts with: Human				
Product overview	Abcam's IL-17a Human <i>in vitro</i> 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.

## Platform

Microplate

## Properties

### Storage instructions

Store at +4°C. Please refer to protocols.

Components	Identifier	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

### Function

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

### Tissue specificity

Restricted to activated memory T-cells.

### Sequence similarities

Belongs to the IL-17 family.

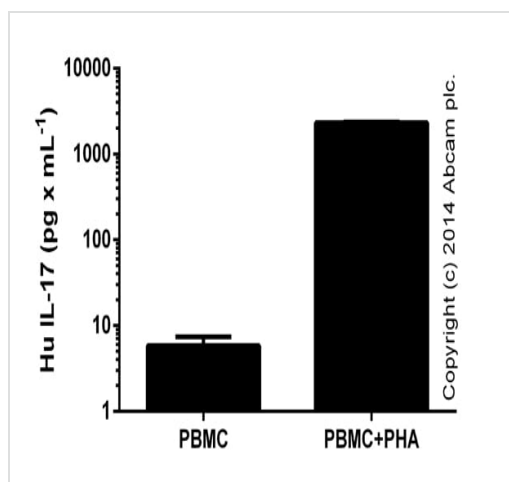
### Post-translational modifications

Found both in glycosylated and nonglycosylated forms.

### Cellular localization

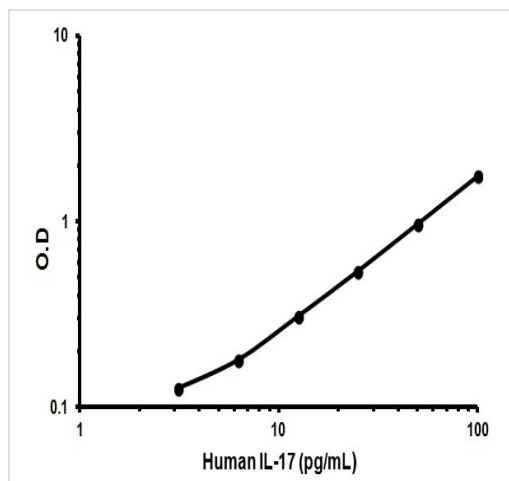
Secreted.

## Images



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.com/abpromise> or contact our technical team.

## Terms and conditions

---

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