

Human IL-17a ELISA Kit ab83688

6 References [画像数 2](#)

製品の概要

製品名 Human IL-17a ELISA Kit

検出方法 Colorimetric

再現性 Intra-Assay (同時再現性)

サンプル	N	平均値	SD	CV%
Overall	6			3.3%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall	6			5.2%

サンプルの種類 Cell culture supernatant, Serum, Plasma

アッセイタイプ Sandwich (quantitative)

検出感度 < 3 pg/ml

検出範囲 3.125 pg/ml - 100 pg/ml

添加回収試験 105 %

特定サンプルでの回収試験

サンプルの種類	平均 %	測定範囲
Serum	121	113% - 130%

全工程の試験時間 3h 45m

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 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.

試験プラットフォーム

Microplate

製品の特性

保存方法

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

内容	ラベル	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

機能

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

組織特異性

Restricted to activated memory T-cells.

配列類似性

Belongs to the IL-17 family.

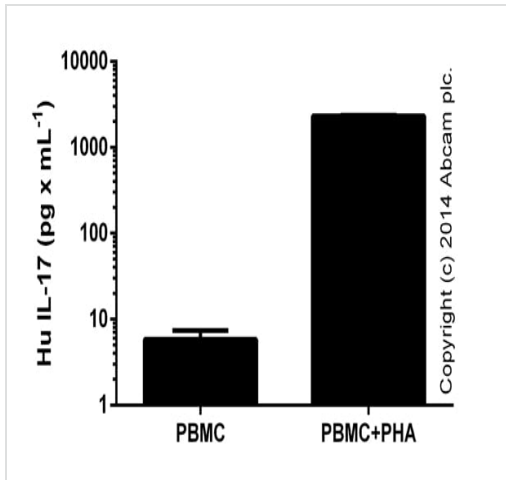
翻訳後修飾

Found both in glycosylated and nonglycosylated forms.

細胞内局在

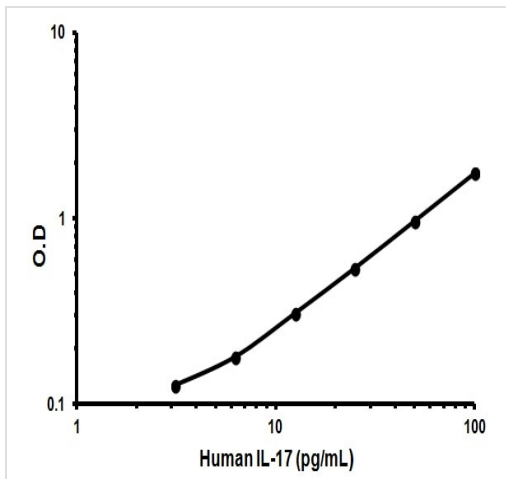
Secreted.

画像



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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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