

Human IL-17 ELISA Kit ab119535

7 References [画像数 1](#)

製品の概要

製品名 Human IL-17 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				7.1%

Inter-Assay (日差再現性)

サンプル	N	平均値	SD	CV%
Overall				9.1%

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

アッセイタイプ Sandwich (quantitative)

検出感度 0.5 pg/ml

検出範囲 1.6 pg/ml - 100 pg/ml

添加回収試験 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
Cell culture supernatant	113.6	% - %
Serum	107	% - %

ステップ Multiple steps standard assay

種交差性 交差種: Human

製品の概要 Abcam's IL-17 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human IL-17 concentrations in Cell culture supernatant, serum and plasma (citrate, heparin).

Human IL-17 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated Human IL-17 detection antibody then incubated at room temperature. Following washing, a Streptavidin-HRP conjugate is added

to each well, incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after addition of an acidic stop solution. The density of yellow coloration is directly proportional to the amount of Human IL-17 captured on the plate.

試験プラットフォーム

Microplate

製品の特性

保存方法

Please refer to protocols.

内容	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-Human IL-17 monoclonal antibody	1 x 70µl
Control high (80 - 220 pg/ml)	1 vial
Control low (5 - 20 pg/ml)	1 vial
Human IL-17 Standard Lyophilized (200 pg/mL on reconstitution)	2 vials
Microplate coated with monoclonal antibody to Human IL-17 (12 x 8 wells)	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml

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.

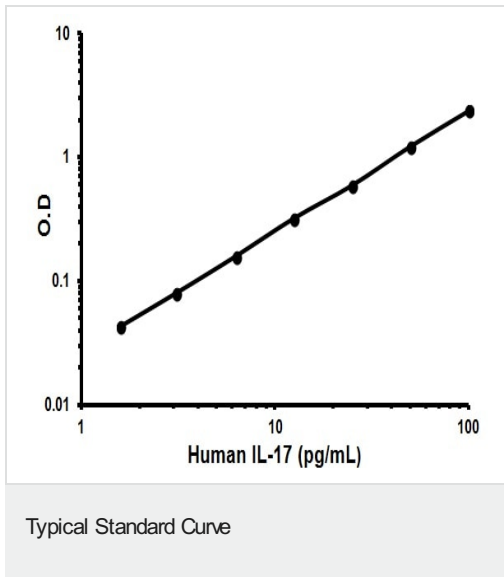
翻訳後修飾

Found both in glycosylated and nonglycosylated forms.

細胞内局在

Secreted.

画像



Representative Standard Curve using ab119535.

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