

Human IL-17 ELISA Set (without plates) ab83707

2 References [画像数 1](#)

製品の概要

製品名	Human IL-17 ELISA Set (without plates)
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
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	<p>Abcam's IL-17 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-17 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The IL-17A ELISA Set includes both the capture and the detection antibodies as well as Horse-Radish peroxidase conjugated Streptavidin (HRP-Strep) and TMB which are the reagents used in the colour development step.</p> <p>This ELISA Set is intended to measure human IL-17A in cell -culture supernatants, as well as in serum and plasma samples. The enclosed procedures provide general guidelines for ELISA preparation. This procedure may need some optimization depending on the plastic, matrix, standard range.</p> <p>SPECIFICITY: Recognizes both natural and recombinant human IL-17A</p> <p>RANGE: 3.12pg/ml - 100pg/ml</p> <p>SENSITIVITY: <2.3pg/ml</p>
アプリケーション	適用あり: Sandwich ELISA
試験プラットフォーム	Reagents

製品の特性

保存方法

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

内容	5 x 96 tests	10 x 96 tests
Detection biotinylated anti-IL-17 antibody	1 x 0.55ml	2 x 0.55ml
IL-17 Capture Antibody	1 x 0.5ml	2 x 0.5ml
IL-17 standard	5 vials	10 vials
Ready-to-use TMB	2 vials	4 vials
Streptavidin-HRP	1 vial	2 vials

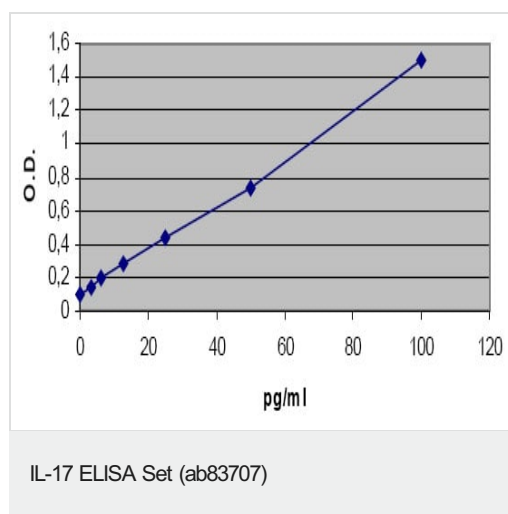
機能	Induces stromal cells to produce proinflammatory and hematopoietic cytokines. Enhances the surface expression of the intracellular adhesion molecule-1 (ICAM-1) in fibroblasts.
組織特異性	Restricted to activated memory T-cells.
配列類似性	Belongs to the IL-17 family.
翻訳後修飾	Found both in glycosylated and nonglycosylated forms.
細胞内局在	Secreted.

アプリケーション

The Abpromise guarantee Abpromise保証は、 次のテスト済みアプリケーションにおけるab83707の使用に適用されます
アプリケーションノートには、推奨の開始希釈率がありますが、適切な希釈率につきましてはご検討ください。

アプリケーション	Abreviews	特記事項
Sandwich ELISA		Use at an assay dependent concentration.

画像



IL-17A Standard curve ranging from 3.125 to 100 pg/mL

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