

Human IL-17 ELISA Set (without plates) ab83707

2 References 1 图像

概述

产品名称	人IL-17 ELISA试剂盒(without plates)
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
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	<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>
经测试应用	适用于: Sandwich ELISA
平台	Reagents

性能

存放说明Store at +4°C. Please refer to protocols.

组件	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

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.

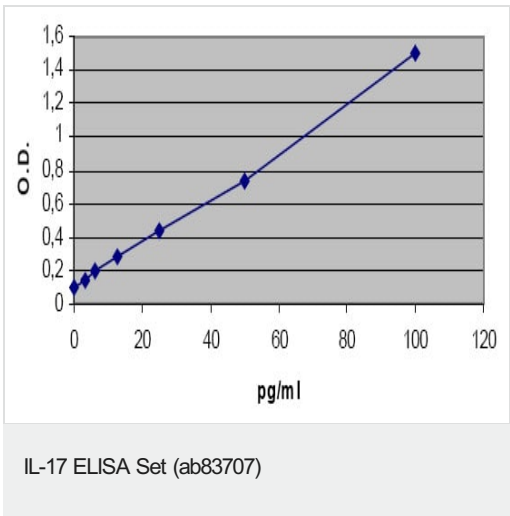
应用

The Abpromise guaranteeAbpromise™承诺保证使用ab83707于以下的经测试应用

“应用说明”部分 下显示的仅为推荐的起始稀释度;实际最佳的稀释度/浓度应由使用者检定。

应用	Ab评论	说明
Sandwich ELISA		Use at an assay dependent concentration.

图片



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.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