

Human IL-5 ELISA Set (without plates) ab47356

1 图像

概述

产品名称	人IL-5 ELISA试剂盒(without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 5 pg/ml
范围	7.8 pg/ml - 250 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-5 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-5 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure human IL-5 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>These reagents are intended for research use only, not for diagnostic purposes.</p> <p>The IL-5 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>Recognizes both natural and recombinant human IL-5.</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	标识符	10 x 96 tests	5 x 96 tests	15 x 96 tests
Chromogen TMB Substrate Solution		4 vials	2 vials	6 vials
Detection biotinylated anti-IL-5 antibody		2 x 0.55ml	1 x 0.55ml	3 x 0.55ml
IL-5 Capture antibody		2 x 0.5ml	1 x 0.5ml	3 x 0.5ml
IL-5 standard	Yellow	10 vials	5 vials	15 vials
Streptavidin-HRP		2 vials	1 vial	3 vials

相关性 Interleukin 5 (IL-5) is a cytokine produced primarily by activated T lymphocytes. It exists as an antiparallel disulfide linked homodimeric glycoprotein with 115 amino acid residues in each chain. Known also as EDF (eosinophil differentiating factor), functions predominantly as an eosinophilopoietic factor. Analysis of its crystal structure reveals a novel two domain structure, with each domain showing significant homology to the cytokine fold in GM-CSF, MCSF, IL2, IL4 and growth hormone. Human and mouse IL5 have 70% amino acid sequence homology.

细胞定位 Secreted

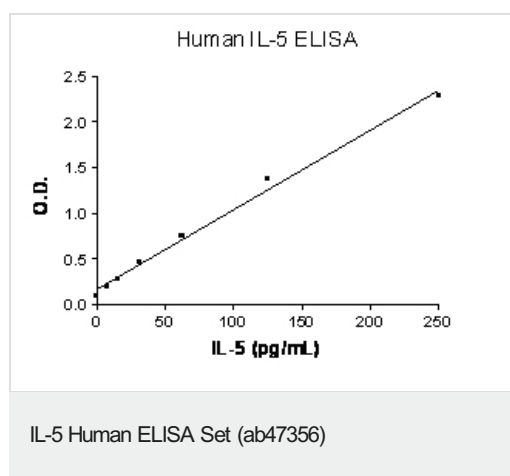
应用

The Abpromise guarantee **Abpromise™** 承诺保证使用 ab47356 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-5 Standard curve ranging from 7.8 to 250 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