

Human IL-7 ELISA Set (without plates) ab47357

1 图像

概述

产品名称	人IL-7 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	6.25 pg/ml - 200 pg/ml
检测时间	2h 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-7 Human <i>in vitro</i> ELISA Set is designed for the quantitative measurement of IL-7 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>This ELISA Set is intended to measure human IL-7 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-7 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-7.</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	15 x 96 tests	5 x 96 tests	10 x 96 tests
Chromogen TMB Substrate Solution		6 vials	2 vials	4 vials
Detection biotinylated anti-IL-7 antibody		3 x 0.55ml	1 x 0.55ml	2 x 0.55ml
IL-7 Capture antibody		3 x 0.25ml	1 x 0.25ml	2 x 0.25ml
IL-7 standard	Yellow	15 vials	5 vials	10 vials
Streptavidin-HRP		3 vials	1 vial	2 vials

相关性 Interleukin 7 (IL-7) is a lymphoid cell growth factor that affects pre-B, pro-B, and early T cells. IL-7 was previously known as pre-B cell growth factor and lymphopoietin 1. IL-7 supports the growth of early B cells from long-term lymphoid bone marrow cultures. It is mitogenic to thymocytes and enhances the response of cells to other stimuli such as polyhydroxyalkanoate (PHA) and concanavalin A (ConA). IL-7 stimulates the proliferation of CD4+/CD8+ cells. The proliferative response of thymocytes to IL-7 is not affected by antibodies to the T cell growth factors such as IL2, IL4 and IL6, suggesting that IL-7 is capable of stimulating T cell proliferation through a pathway independent of the known T cell growth factors. Mature T cells respond to IL-7 and Con A, but not to IL-7 alone. In myeloid lineage cells, IL-7 upregulates the production of pro-inflammatory cytokines and stimulates the tumoricidal activity of monocytes/macrophages. IL-7 is expressed by adherent stromal cells from various tissues.

细胞定位 Secreted

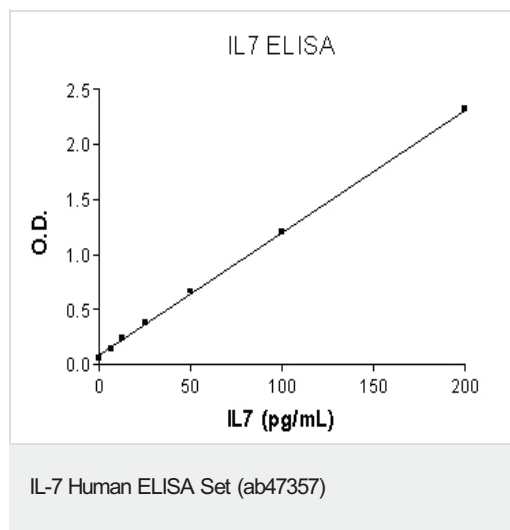
应用

The Abpromise guarantee **Abpromise™** 承诺保证使用ab47357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-7 Standard curve ranging from 6.25 to 200 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