

Product datasheet

Human IL-7 ELISA Kit (Interleukin-7) ab46031

1 图像

概述

产品名称 人IL-7 ELISA试剂盒(Interleukin-7)

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Sample A	8	198.7pg/ml	1.667	0.83%
Sample B	8	28.69pg/ml	1.118	3.86%

批次间

样品	n	Mean	SD	CV%
Sample A	29	198.9pg/ml	3.772	1.89%
Sample B	29	25.68pg/ml	1.505	5.84%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 3 pg/ml

范围 6 pg/ml - 200 pg/ml

复检 97.6 %

检测时间 2h 40m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Human

产品概述 Abcam's IL-7 Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-7 in Human serum, plasma, buffered solutions and cell culture medium.

A monoclonal antibody specific for IL-7 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-7 concentrations, control specimens or unknowns are pipetted into these wells. During the first incubation, the standards or samples and a biotinylated monoclonal antibody specific for IL-7 are simultaneously incubated. After washing,

the enzyme Streptavidin-HRP, that binds the biotinylated antibody is added, incubated and washed. A TMB substrate solution is added which acts on the bound enzyme to induce a colored reaction product. The intensity of this colored product is directly proportional to the concentration of IL-7 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-7.

Get higher sensitivity in only 90 minutes with Human IL-7 ELISA Kit ([ab185984](#)) from our SimpleStep ELISA[®] range.

经测试应用

适用于: Sandwich ELISA

平台

Microplate

性能

存放说明

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

组件	标识符	1 x 96 tests	2 x 96 tests
10X Standard Diluent Buffer	Black	1 x 25ml	1 x 25ml
200X Wash Buffer	White	1 x 10ml	2 x 10ml
Biotinylated Antibody Diluent	Red	1 x 7.5ml	1 x 13ml
Biotinylated anti-IL-7	Red	1 x 400µl	2 x 400µl
Chromogen TMB Substrate Solution		1 x 11ml	1 x 24ml
IL-7 Control	Silver	2 vials	4 vials
HRP Diluent	Red	1 x 23ml	1 x 23ml
IL-7 Microplate (12 x 8 well strip)		1 unit	2 units
IL7 standard	Yellow	2 vials	4 vials
Standard Diluent (Human Serum)		1 x 7ml	2 x 7ml
Stop Reagent	Black	1 x 11ml	2 x 11ml
Streptavidin-HRP		2 x 5µl	4 x 5µl

相关性

IL7 is a cytokine important for B and T cell development. It and the hepatocyte growth factor form a heterodimer that functions as a pre-pro-B cell growth-stimulating factor. IL7 is found to be a cofactor for V(D)J rearrangement of the T cell receptor beta (TCRB) during early T cell development. It can be produced locally by intestinal epithelial and epithelial goblet cells, and may serve as a regulatory factor for intestinal mucosal lymphocytes. Knockout studies in mice suggested that IL7 plays an essential role in lymphoid cell survival.

细胞定位

Secreted

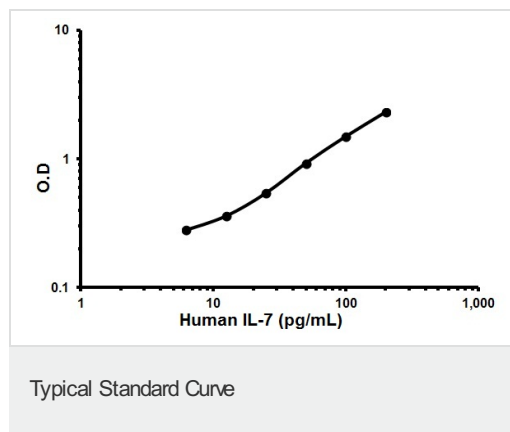
应用

Our [Abpromise guarantee](#) covers the use of **ab46031** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

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

图片



Representative Standard Curve using ab46031

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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