

Human IL-10 ELISA Kit High Sensitivity ab46059

5 References 1 图像

| 概述 | | | | | | | | | | | |
|--------------------------|---|-----------|------|------|--------------------------|-----|---------|-------|----|-----------|------|
| 产品名称 | 人IL-10 ELISA试剂盒High Sensitivity | | | | | | | | | | |
| 检测方法 | Colorimetric | | | | | | | | | | |
| 精确度 | 批次内 | | | | | | | | | | |
| | <table><tr><th>样品</th><th>n</th><th>Mean</th><th>SD</th><th>CV%</th></tr><tr><td>Overall</td><td>3</td><td></td><td></td><td>7.8%</td></tr></table> | 样品 | n | Mean | SD | CV% | Overall | 3 | | | 7.8% |
| | 样品 | n | Mean | SD | CV% | | | | | | |
| Overall | 3 | | | 7.8% | | | | | | | |
| 批次间 | | | | | | | | | | | |
| 样品类型 | Cell culture supernatant, Serum, Plasma | | | | | | | | | | |
| | 检测类型 | | | | | | | | | | |
| | Sandwich (quantitative) | | | | | | | | | | |
| 灵敏度 | < 1.3 pg/ml | | | | | | | | | | |
| 范围 | 1.56 pg/ml - 50 pg/ml | | | | | | | | | | |
| 回收率 | 87 % | | | | | | | | | | |
| | 特定样本回收率 | | | | | | | | | | |
| | <table><tr><th>样品类型</th><th>平均 %</th><th>范围</th></tr><tr><td>Cell culture supernatant</td><td>94</td><td>% - %</td></tr><tr><td>Serum</td><td>80</td><td>59% - 93%</td></tr></table> | 样品类型 | 平均 % | 范围 | Cell culture supernatant | 94 | % - % | Serum | 80 | 59% - 93% | |
| 样品类型 | 平均 % | 范围 | | | | | | | | | |
| Cell culture supernatant | 94 | % - % | | | | | | | | | |
| Serum | 80 | 59% - 93% | | | | | | | | | |
| 检测时间 | 3h 00m | | | | | | | | | | |
| 实验步骤 | Multiple steps standard assay | | | | | | | | | | |
| 种属反应性 | 与反应: Human | | | | | | | | | | |
| 产品概述 | Abcam's IL-10 Human High Sensitivity <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 in Human sera, plasmas, buffered solutions and cell culture media. | | | | | | | | | | |

A monoclonal antibody specific for IL-10 has been coated onto the wells of the microtiter strips provided. Samples, including standards of known IL-10 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-10 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-10 present in the samples.

This kit will recognize both endogenous and recombinant Human IL-10.

Get results in 90 minutes with Human IL-10 ELISA Kit ([**ab185986**](#)) from our SimpleStep ELISA[®] range.

平台 Microplate

性能

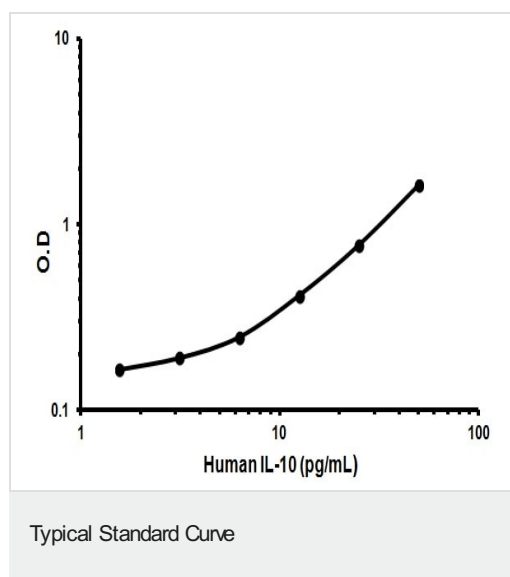
存放说明 Store at +4°C. Please refer to protocols.

| 组件 | 标识符 | 1 x 96 tests |
|---------------------------------------|-------------------|--------------|
| 10X Standard Diluent Buffer | Black | 1 x 15ml |
| 200X Wash Buffer | White | 1 x 10ml |
| Amplification Diluent | Brown & blue spot | 1 x 15ml |
| Amplifier | Yellow | 1 x 200µl |
| Biotinylated anti IL-10 | Red | 1 x 400µl |
| Biotinylated Antibody Diluent | Red | 1 x 7.5ml |
| Chromogen TMB Substrate Solution | | 1 x 11ml |
| HRP Diluent | Red | 1 x 23ml |
| IL-10 Microplate (12 x 8 well strips) | | 1 unit |
| IL-10 Standard (Lyophilized) | Yellow | 2 vials |
| Stop Reagent | Black | 1 x 11ml |
| | | |

| 组件 | 标识符 | 1 x 96 tests |
|------------------|-----|--------------|
| Streptavidin-HRP | | 2 x 5µl |

| | |
|--------------|---|
| 功能 | Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells. |
| 组织特异性 | Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types. |
| 序列相似性 | Belongs to the IL-10 family. |
| 细胞定位 | Secreted. |

图片



Representative Standard Curve using ab46059

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