

Mouse IL-10 ELISA Kit ab46103

12 References 1 图像

概述

产品名称 小鼠IL-10 ELISA试剂盒

检测方法 Colorimetric

| 精确度 | 批次内 | | | | |
|-----|-----|---|------------|------|------|
| | 样品 | n | Mean | SD | CV% |
| | A | 8 | 556.8pg/ml | 11.6 | 2% |
| | B | 8 | 137.4pg/ml | 6.8 | 4.9% |

| | 批次间 | | | | |
|--|-----|---|------------|------|------|
| | 样品 | n | Mean | SD | CV% |
| | A | 8 | 548.2pg/ml | 12.1 | 2.2% |
| | B | 8 | 133.6pg/ml | 5.4 | 4% |

样品类型 Cell culture supernatant, Serum

检测类型 Sandwich (quantitative)

灵敏度 < 20 pg/ml

范围 31.25 pg/ml - 1000 pg/ml

检测时间 3h 45m

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's IL-10 mouse *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of IL-10 in mouse serum, buffered solutions and cell culture medium.

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 mouse IL-10.

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 15ml | 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-10 | Red | 1 x 400µl | 2 x 400µl |
| Chromogen TMB Substrate Solution | | 1 x 11ml | 1 x 24ml |
| HRP Diluent | Red | 1 x 12ml | 1 x 23ml |
| IL-10 Microplate (12 x 8 well strips) | | 1 unit | 2 units |
| Murine IL-10 standard | | 2 vials | 4 vials |
| Stop Reagent | Black | 1 x 11ml | 2 x 11ml |
| Streptavidin-HRP | | 2 x 5µl | 4 x 5µl |

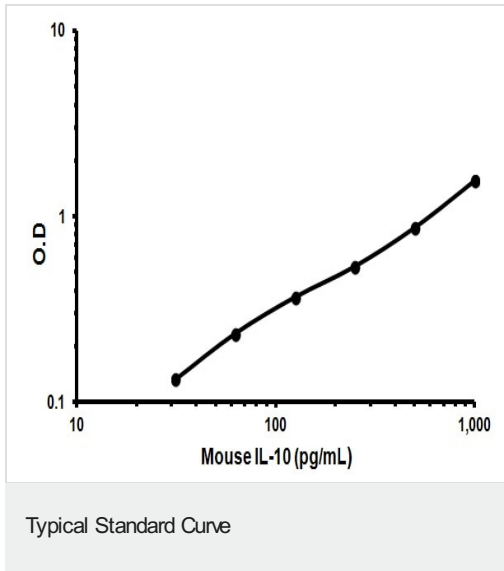
功能 Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

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 ab46103

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