

Mouse IL-10 ELISA Set (without plates) ab47599

1 图像

概述	
产品名称	小鼠IL-10 ELISA试剂盒 (without plates)
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 20 pg/ml
范围	31.25 pg/ml - 1000 pg/ml
实验步骤	Multiple steps standard assay
种属反应性	与反应: Mouse
产品概述	<p>Abcam's IL-10 Murine <i>in vitro</i> ELISA Set is designed for the quantitative measurement of Murine IL-10 in cell culture supernatants, buffered solutions, serum, plasma and other body fluids.</p> <p>The murine IL-10 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>This ELISA Set is intended to measure murine IL-10 in cell -culture supernatants, as well as in serum 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>Recognizes both natural and recombinant Murine IL-10</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	5 x 96 tests	15 x 96 tests	10 x 96 tests
Chromogen TMB Substrate Solution	2 vials	6 vials	4 vials
Detection biotinylated anti-murine IL-10 antibody	1 x 0.55ml	3 x 0.55ml	2 x 0.55ml

组件	5 x 96 tests	15 x 96 tests	10 x 96 tests
Murine IL-10 Capture antibody	1 x 0.5ml	3 x 0.5ml	2 x 0.5ml
Murine IL-10 standard	5 vials	15 vials	10 vials
Streptavidin-HRP	1 vial	3 vials	2 vials

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.

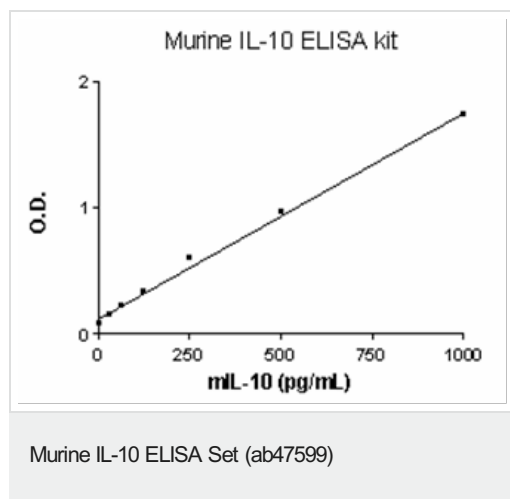
应用

The Abpromise guarantee Abpromise™承诺保证使用ab47599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



Murine IL-10 standard curve ranging from 31.25 to 1000 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