

Mouse IL-10 ELISA kit ab108870

★★★★☆ [1 Abreviews](#) [34 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%
Overall				5.8%

批次间

样品	n	Mean	SD	CV%
Overall				10%

样品类型 Cell culture supernatant, Serum, Plasma, Tissue, Cell Lysate

检测类型 Sandwich (quantitative)

灵敏度 0.014 ng/ml

范围 0.125 ng/ml - 1 ng/ml

回收率 97 %

检测时间 5h 00m

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 plasma, serum, tissue extracts and cell culture supernatants.

An IL-10 specific antibody has been precoated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an IL-10 specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Conjugate is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of IL-10 captured in plate.

The entire kit may be stored at -20°C for long term storage before reconstitution - Avoid repeated freeze-thaw cycles.

平台 Microplate

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	1 x 96 tests
100X Streptavidin-Peroxidase Conjugate	1 x 80µl
10X Diluent N Concentrate	1 x 30ml
1X Standard Diluent	1 x 2ml
20X Wash Buffer Concentrate	2 x 30ml
30X Biotinylated Mouse IL-10 Antibody	1 x 180µl
Chromogen Substrate	1 x 7ml
IL-10 Microplate (12 x 8 well strips)	1 unit
IL-10 Standard	1 vial
Sealing Tapes	3 units
Stop Solution	1 x 11ml

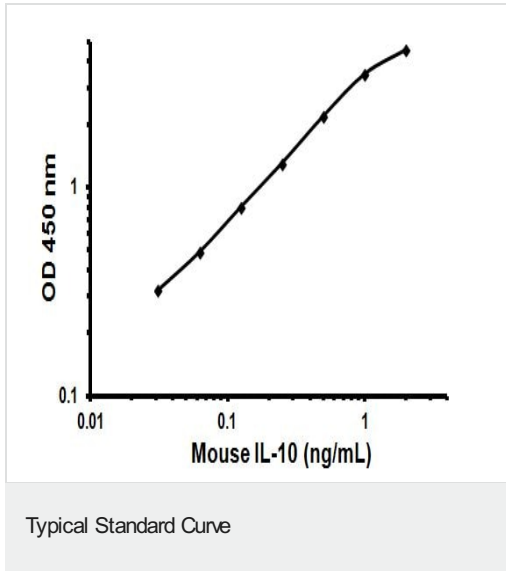
功能 Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

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 ab108870

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