

Rat IL-10 ELISA Kit ab100764

[8 References](#) [2 图像](#)

概述

产品名称 大鼠IL-10 ELISA试剂盒

检测方法 Colorimetric

精确度 批次内

样品	n	Mean	SD	CV%
Overall				< 10%

批次间

样品	n	Mean	SD	CV%
Overall				< 12%

样品类型 Cell culture supernatant, Serum, Plasma

检测类型 Sandwich (quantitative)

灵敏度 < 10 pg/ml

范围 8.23 pg/ml - 6000 pg/ml

回收率 97 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	98.69	90% - 109%
Serum	96.75	84% - 104%
Plasma	95.97	82% - 102%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Rat

产品概述 Abcam's IL-10 Rat ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of rat IL-10 in serum, plasma, and cell culture supernatants.

This assay employs an antibody specific for rat IL-10 coated on a 96-well plate. Standards and

samples are pipetted into the wells and IL-10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-rat IL-10 antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of IL-10 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

平台 Microplate

性能

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

组件	1 x 96 tests
140X HRP-Streptavidin Concentrate	1 x 200µl
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
Biotinylated anti-Rat IL-10	2 vials
IL-10 Microplate (12 strips x 8 wells)	1 unit
Recombinant rat IL-10 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

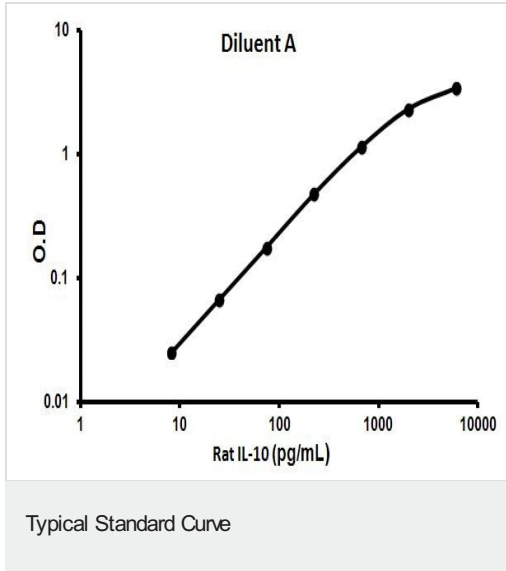
功能 Inhibits the synthesis of a number of cytokines, including IFN-gamma, IL-2, IL-3, TNF and GM-CSF produced by activated macrophages and by helper T-cells.

组织特异性 Produced by a variety of cell lines, including T-cells, macrophages, mast cells and other cell types.

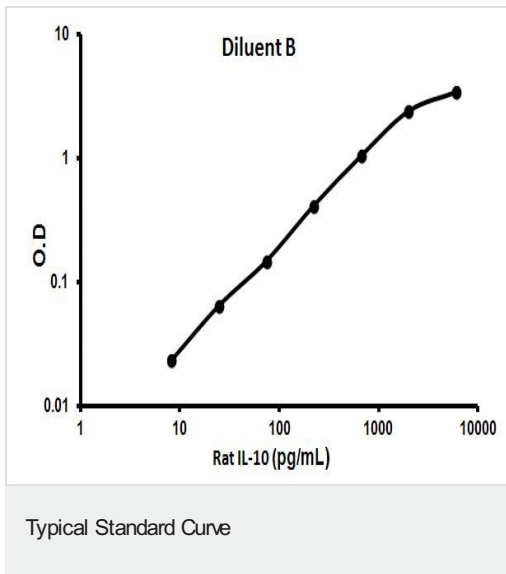
序列相似性 Belongs to the IL-10 family.

细胞定位 Secreted.

图片



Representative standard curve using ab100764.



Representative standard curve using ab100764.

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