

Mouse IP10 ELISA Kit ab100675

1 References 2 图像

概述

产品名称	小鼠IP10 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture supernatant, Serum, Plasma
检测类型	Sandwich (quantitative)
灵敏度	< 15 pg/ml
范围	8.2 pg/ml - 2000 pg/ml
回收率	97 %

特定样本回收率

样品类型	平均%	范围
Cell culture supernatant	98.86	88% - 107%
Serum	97.68	86% - 106%
Plasma	96.42	85% - 104%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's IP10 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of Mouse IP10 in serum, plasma and cell culture supernatants.

This assay employs an antibody specific for mouse IP10 coated on a 96- well plate. Standards and samples are pipetted into the wells and IP10 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse IP10 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 IP10 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Get results in 90 minutes with Mouse IP-10 ELISA Kit (CXCL10) (**ab214563**) from our SimpleStep ELISA® range.

平台Microplate

性能

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

组件	1 x 96 tests
20X Wash Buffer	1 x 25ml
5X Assay Diluent B	1 x 15ml
80x HRP-Streptavidin Concentrate	1 x 200μl
Assay Diluent A	1 x 30ml
Biotinylated anti-Mouse IP10	2 vials
IP10 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse IP10 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

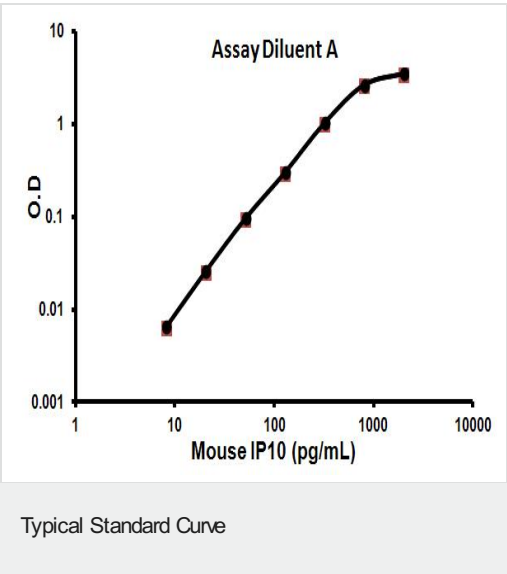
功能Chemotactic for monocytes and T-lymphocytes. Binds to CXCR3.

序列相似性Belongs to the intercrine alpha (chemokine CxC) family.

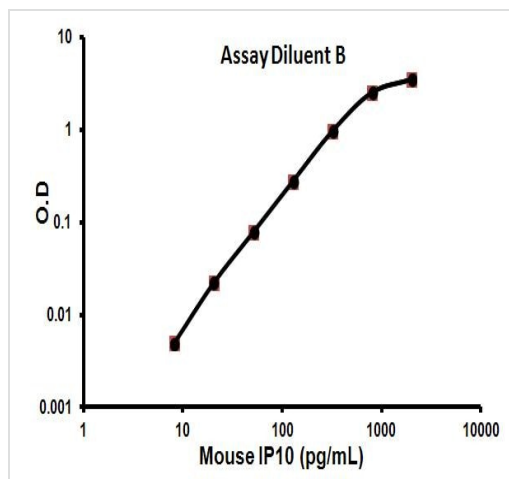
翻译后修饰CXCL10(1-73) is produced by proteolytic cleavage after secretion from keratinocytes.

细胞定位Secreted.

图片



Representative Standard Curve using ab100675



Representative Standard Curve using ab100675

Typical Standard Curve

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