

Mouse MCP1 ELISA Kit ab100722

[4 References](#) [1 图像](#)

概述

产品名称	小鼠MCP1 ELISA试剂盒
检测方法	Colorimetric
样品类型	Cell culture extracts, Tissue Extracts
检测类型	Sandwich (quantitative)
灵敏度	< 3 pg/ml
范围	2.74 pg/ml - 2000 pg/ml
回收率	86 %

特定样本回收率

样品类型	平均%	范围
Cell culture extracts	87.56	80% - 102%
Tissue Extracts	85.89	78% - 101%

实验步骤 Multiple steps standard assay

种属反应性 与反应: Mouse

产品概述 Abcam's MCP1 Mouse ELISA (Enzyme-Linked Immunosorbent Assay) kit is an *in vitro* enzyme-linked immunosorbent assay for the quantitative measurement of mouse MCP1 in cell lysates and tissue lysates.

This assay employs an antibody specific for mouse MCP1 coated on a 96-well plate. Standards and samples are pipetted into the wells and MCP1 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-mouse MCP1 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 MCP1 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

说明 Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

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
2X Cell Lysis Buffer	1 x 5ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent	1 x 15ml
5X Sample Diluent Buffer	1 x 10ml
Biotinylated anti-Mouse MCP1	2 vials
MCP1 Microplate (12 x 8 wells)	1 unit
Recombinant Mouse MCP1 Standard (lyophilized)	2 vials
Stop Solution	1 x 8ml
TMB One-Step Substrate Reagent	1 x 12ml

功能

Chemotactic factor that attracts monocytes and basophils but not neutrophils or eosinophils. Augments monocyte anti-tumor activity. Has been implicated in the pathogenesis of diseases characterized by monocytic infiltrates, like psoriasis, rheumatoid arthritis or atherosclerosis. May be involved in the recruitment of monocytes into the arterial wall during the disease process of atherosclerosis.

序列相似性

Belongs to the intercrine beta (chemokine CC) family.

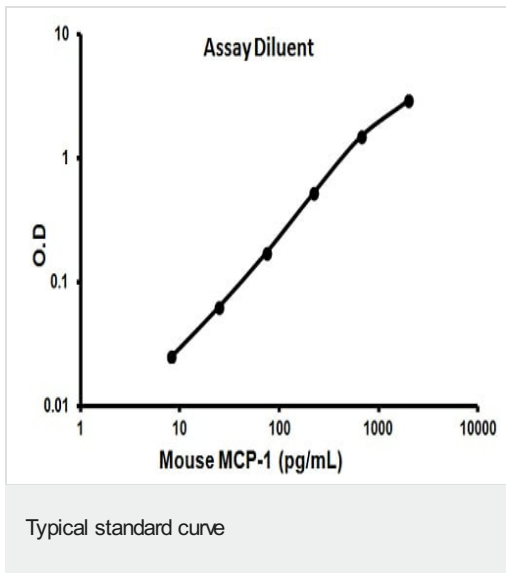
翻译后修饰

Processing at the N-terminus can regulate receptor and target cell selectivity. Deletion of the N-terminal residue converts it from an activator of basophil to an eosinophil chemoattractant.

细胞定位

Secreted.

图片



Representative standard curve using ab100722

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