

Mouse MCP1 ELISA Kit ab100721

★★★★★ [2 Abreviews](#) [15 References](#) [3 Images](#)

Overview

Product name	Mouse MCP1 ELISA Kit
Detection method	Colorimetric
Sample type	Cell culture supernatant, Serum, Plasma
Assay type	Sandwich (quantitative)
Sensitivity	< 3 pg/ml
Range	2.74 pg/ml - 2000 pg/ml
Recovery	94 %

Sample specific recovery

Sample type	Average %	Range
Cell culture supernatant	95.29	84% - 104%
Serum	94.28	83% - 103%
Plasma	92.79	82% - 102%

Assay duration Multiple steps standard assay

Species reactivity **Reacts with:** Mouse

Product overview 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 serum, plasma and cell culture supernatants.

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.

Platform Microplate

Properties

**Storage instructions**

Store at -20°C. Please refer to protocols.

Components	1 x 96 tests
20X Wash Buffer	1 x 25ml
400X HRP-Streptavidin Concentrate	1 x 200µl
5X Assay Diluent B	1 x 15ml
Assay Diluent A	1 x 30ml
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

**Function**

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.

**Sequence similarities**

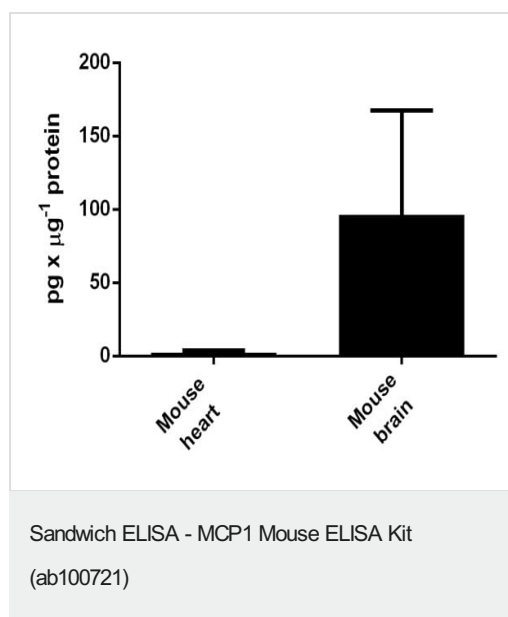
Belongs to the intercrine beta (chemokine CC) family.

**Post-translational modifications**

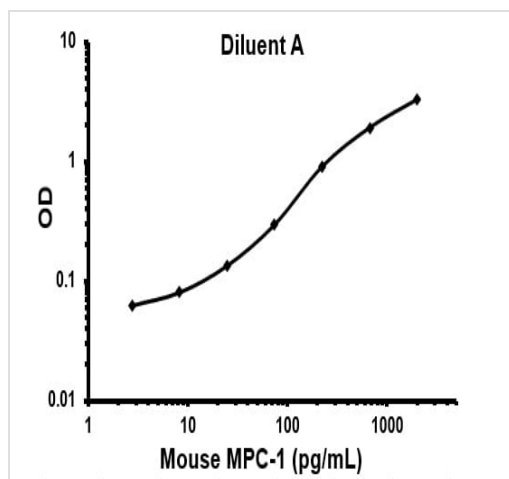
Processing at the N-terminus can regulate receptor and target cell selectivity. Deletion of the N-terminal residue converts it from an activator of basophil to an eosinophil chemoattractant.

**Cellular localization**

Secreted.

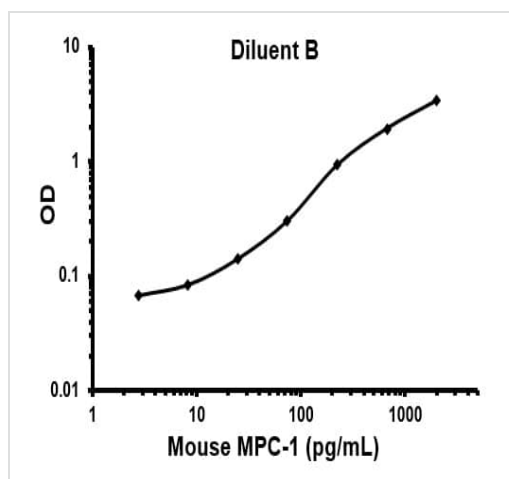
**Images**

MCP1 measured in mouse lysates showing quantity (pg) per microgram of total protein. Values retrieved from 1:100 and 1:1000 dilutions.



Representative standard curve using ab100721.

Typical Standard Curve



Representative standard curve using ab100721.

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.com/abpromise> or contact our technical team.

### Terms and conditions

- 
- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors