## MCP-1 monoclonal antibody (E10052) (biotin conjugate)

MCP-1 is one of several cytokine genes clustered on the q-arm of chromosome 17. Cytokines are a family of secreted proteins involved in immunoregulatory and inflammatory processes. The protein encoded by this gene is structurally related to the CXC subfamily of cytokines. Members of this subfamily are characterized by two cysteines separated by a single amino acid. This cytokine displays chemotactic activity for monocytes and basophils but not for neutrophils or eosinophils. It has been implicated in the pathogenesis of diseases characterized by monocytic infiltrates, like psoriasis, rheumatoid arthritis and atherosclerosis. It binds to chemokine receptors CCR2 and CCR4.

This antibody is covered by our Worry-Free Guarantee.

**Ordering Information** 

Order Online »

ENZ-ABS281-0100

100µg

Manuals, SDS & CofA

**View Online** »

## **Handling & Storage**

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## **Product Details**

Alternative Name CCL2, C-C motif chemokine 2, Monocyte chemoattractant

protein 1, Small-inducible cytokine A2

**Application** ELISA

**Clone** E10052

**Formulation** Liquid. In PBS containing no preservative.

**Host** Mouse

**Immunogen** Full-length recombinant human MCP-1.

lsotype lgG1

Purity Detail Protein G affinity purified.

Species Reactivity Human

UniProt ID P13500

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

.

Last modified: May 29, 2024

