

Product datasheet

Human MCP1 ELISA Kit (CCL2) ab179886

SimpleStep ELISA[®]

1 References 2 Images

Overview

Product name Human MCP1 ELISA Kit (CCL2)

Detection method Colorimetric

Precision

Intra-assay				
Sample	n	Mean	SD	CV%
Overall	8			4.3%

Inter-assay				
Sample	n	Mean	SD	CV%
Overall	3			1.6%

Sample type Cell culture supernatant, Serum, Plasma, Heparin Plasma, EDTA Plasma, Citrate Plasma

Assay type Sandwich (quantitative)

Sensitivity 1.1 pg/ml

Range 1.1 pg/ml - 800 pg/ml

Recovery

Sample specific recovery		
Sample type	Average %	Range
Serum	91	86% - 101%
Cell culture media	115	103% - 122%
Heparin Plasma	91	85% - 100%
EDTA Plasma	105	100% - 107%
Citrate Plasma	89	78% - 102%

Assay time 1h 30m

Assay duration One step assay

Species reactivity**Reacts with:** Human**Product overview**

Abcam's MCP1 (CCL2) *in vitro* SimpleStep ELISA® (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of MCP1 (CCL2) protein in Human serum, plasma, and cell culture supernatant.

The SimpleStep ELISA® employs an affinity tag labeled capture antibody and a reporter conjugated detector antibody which immunocapture the sample analyte in solution. This entire complex (capture antibody/analyte/detector antibody) is in turn immobilized via immunoaffinity of an anti-tag antibody coating the well. To perform the assay, samples or standards are added to the wells, followed by the antibody mix. After incubation, the wells are washed to remove unbound material. TMB substrate is added and during incubation is catalyzed by HRP, generating blue coloration. This reaction is then stopped by addition of Stop Solution completing any color change from blue to yellow. Signal is generated proportionally to the amount of bound analyte and the intensity is measured at 450 nm. Optionally, instead of the endpoint reading, development of TMB can be recorded kinetically at 600 nm.

Notes

MCP1 (CCL2) is a 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.

Tested applications**Suitable for:** Sandwich ELISA**Platform**

Microplate

Properties**Storage instructions** Store at +4°C. Please refer to protocols.

Components	1 x 96 tests
10X MCP1 Capture Antibody	1 x 600µl
10X MCP1 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
Antibody Diluent CPI - HAMA Blocker (ab193969)	1 x 6ml
MCP1 Human Lyophilized Recombinant Protein	2 vials
Plate Seal	1 unit
Sample Diluent NS (ab193972)	1 x 50ml
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Substrate	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.

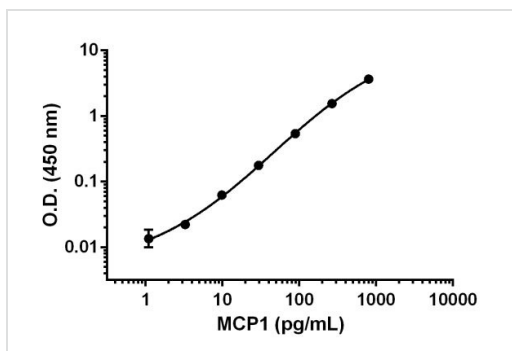
Applications

Our [Abpromise guarantee](#) covers the use of **ab179886** in the following tested applications.

The application notes include recommended starting dilutions; optimal dilutions/concentrations should be determined by the end user.

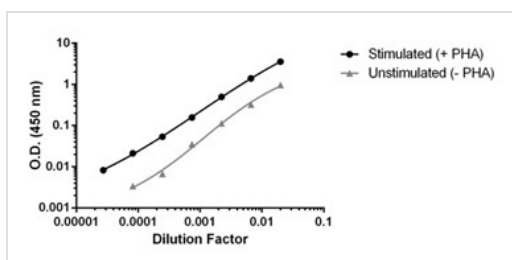
Application	Abreviews	Notes
Sandwich ELISA		Use at an assay dependent concentration.

Human MCP1 ELISA Kit (CCL2) images



Background-subtracted data values (mean +/- SD) are graphed.

Example of MCP1 (CCL2) standard curve.



Background subtracted data from triplicate measurements are plotted.

Titration of PBMC +/- PHA conditioned media within the working range of the assay.

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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 <http://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