

## Mouse CCL2 (MCP-1) Uncoated ELISA

**Catalog Number:** 88-7391

**Also known as:** Monocyte chemotactic protein 1

**RUO: For Research Use Only. Not for use in diagnostic procedures.**

### Product Information

**Contents:** Mouse CCL2 (MCP-1) Uncoated ELISA  
**Catalog Number:** 88-7391  
**Sensitivity:** 15 pg/mL  
**Standard Curve Range:** 15 - 2000 pg/mL

REF



**Temperature Limitation:** Store at 2-8°C except standard which should be stored at less than or equal to -70°C.

LOT



**Batch Code:** Refer to vial



**Use By:** Refer to vial

**Caution, contains Azide**

### Description

This Mouse CCL2 (MCP-1) Uncoated ELISA reagent set (with or without high-affinity binding microwell plates) contains the necessary reagents, buffers and diluents for performing quantitative enzyme linked immunosorbent assays (ELISA). This ELISA reagent set is specifically engineered for accurate and precise measurement of mouse CCL2 protein levels from samples including serum, plasma, and supernatants from cell cultures.

### Components

Capture Antibody: Pre-titrated, purified antibody  
 Detection Antibody: Pre-titrated, biotin-conjugated antibody  
 Standard: Recombinant protein for generating standard curve and calibrating samples  
 10X Coating Buffer : Buffer for plating the Capture Antibody  
 5X ELISA/ELISPOT Diluent : Buffer for blocking and diluting the Detection Antibody and Enzyme  
 Enzyme: Pre-titrated Avidin-HRP  
 Substrate: 1X TMB Solution  
 Certificate of Analysis: Lot-specific instructions for dilution of antibodies and standards  
 96 Well Plate: Corning Costar 9018 (included with product Cat. #'s ending in suffixes -22, -76, -86)

### Applications Tested

This assay has been validated for the detection of endogenous mouse CCL2 with supernatant collected from cultures of mouse thioglycolate-elicited peritoneal macrophages stimulated with LPS.

It is recommended that serum samples be diluted at least 2-fold with 1X ELISA Diluent prior to evaluation. Plasma samples have not been tested in this assay.

This assay was evaluated for specificity on a panel of 72 recombinant cytokines at 100 ng/mL. No significant crossreactivity was observed, including to human CCL2.

### References

Soehnlein O, Zernecke A, Eriksson EE, Rothfuchs AG, Pham CT, Herwald H, Bidzhekov K, Rottenberg ME, Weber C, Lindbom L. Neutrophil secretion products pave the way for inflammatory monocytes. *Blood*. 2008 Aug 15;112(4):1461-71. (RSG ELISA kit, **lavage**, PubMed)

Senaratne RH, Sidders B, Sequeira P, Saunders G, Dunphy K, Marjanovic O, Reader JR, Lima P, Chan S, Kendall S, McFadden J, Riley LW. Mycobacterium tuberculosis strains disrupted in mce3 and mce4 operons are attenuated in mice. *J Med Microbiol*. 2008 Feb;57(Pt 2):164-70. (RSG ELISA kit, **TC supernatant**, PubMed)

### Related Products

88-7399 Human CCL2 (MCP-1) Uncoated ELISA

Not for further distribution without written consent.

Copyright © 2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Tel: 888.999.1371 or 858.642.2058 • Fax: 858.642.2046 • [thermofisher.com/ebioscience](http://thermofisher.com/ebioscience) • [info@ebioscience.com](mailto:info@ebioscience.com)

# Enzyme Linked Immunosorbent Assay (ELISA)

## Protocol: Uncoated ELISA

The following protocol is a general guideline for the Uncoated Sets

### Materials Provided

- Please refer to the Certificate of Analysis (C of A) for components

### Other Materials Needed

- Buffers\*
  - Wash Buffer: 1x PBS, 0.05% Tween-20 (or Thermo Fisher ELISA Wash Buffer Powder, Cat. No. 00-0400)
  - Stop Solution: 1M H<sub>3</sub>PO<sub>4</sub> (recommended) or 2N H<sub>2</sub>SO<sub>4</sub>
- Pipettes
- Refrigerator & frost-free -20°C freezer
- 96-well plate (Corning Costar 9018 or NUNC Maxisorp™)

**NOTE:** The use of ELISA plates which are not high affinity protein binding plates will result in suboptimal performance, e.g., low signal or inconsistent data. Do not use tissue culture plates or low protein absorption plates. Use only the Corning Costar 9018 or NUNC Maxisorp (Cat. No. 44-2404) 96-well plates.

- 96-well ELISA plate reader (microplate spectrophotometer)
- ELISA plate washer (highly recommended)

**NOTE:** To ensure optimal results from this Uncoated ELISA Set, please only use the components included in the set. Exchanging of components is not recommended as a change in signal may occur.

### Time Requirements

- 1 overnight incubation
- 4½-hour incubations
- 1 hour washing and analyzing samples

### Experimental Procedure

1. Coat Corning Costar 9018 (or Nunc Maxisorp™) ELISA plate with 100 µL/well of capture antibody in 1X Coating Buffer (dilute as noted on C of A, which is included with the reagent set). Seal the plate and incubate overnight at 4°C.
2. Aspirate wells and wash 3 times with >250 µL/well Wash Buffer\*. Allowing time for soaking (~ 1 minute) during each wash step increases the effectiveness of the washes. Blot plate on absorbent paper to remove any residual buffer.
3. Dilute 1 part 5X ELISA/ELISPOT Diluent with 4 parts DI water\*. Block wells with 200 µL/well of 1X ELISA/ELISPOT Diluent. Incubate at room temperature for 1 hour.
4. Optional: Aspirate and wash at least once with Wash Buffer.
5. Using 1X ELISA/ELISPOT Diluent \*, dilute standards as noted on the C of A to prepare the top concentration of the standard. Add 100 µL/well of top standard concentration to the appropriate wells. Perform 2-fold serial dilutions of the top standards to make the standard curve for a total of 8 points. Add 100 µL/well of your samples to the appropriate wells. Seal the plate and incubate at room temperature for 2 hours (or overnight at 4°C for maximal sensitivity).

6. Aspirate/wash as in step 2. Repeat for a total of 3-5 washes\*\*.
7. Add 100  $\mu$ L/well of detection antibody diluted in 1X ELISA/ELISPOT Diluent \* (dilute as noted on C of A). Seal the plate and incubate at room temperature for 1 hour.
8. Aspirate/wash as in step 2. Repeat for a total of 3-5 washes\*\*.
9. Add 100  $\mu$ L/well of Avidin-HRP\* diluted in 1X ELISA/ELISPOT Diluent (dilute as noted on C of A). Seal the plate and incubate at room temperature for 30 minutes.
10. Aspirate and wash as in step 2. In this wash step, soak wells in Wash Buffer\* for 1 to 2 minutes prior to aspiration. Repeat for a total of 5-7 washes\*\*.
11. Add 100  $\mu$ L/well of 1X TMB Solution to each well. Incubate plate at room temperature for 15 minutes.
12. Add 50  $\mu$ L of Stop Solution to each well.
13. Read plate at 450 nm. If wavelength subtraction is available, subtract the values of 570 nm from those of 450 nm and analyze data.

**NOTES:**

**\* Be certain that no sodium azide is present in the solutions used in this assay, as this inhibits HRP enzyme activity.**

**\*\*The number of washes in the protocol was adapted to an automatic plate washer. This can be decreased when using other methods but should be tested empirically. Allowing time for soaking (~ 1 minute) during each wash step increases the effectiveness of the washes.**

## Quick Guide: Standard Calibration

The following table indicates the protein standard contained in the Uncoated kit is calibrated against NIBSC standards.

Table of Standard Calibration				
Cytokine	ng of eB standard	ng of NIBSC standard	U of NIBSC standard	NIBSC Lot #
hIL-2	1	1.1	14.6	86/564
hIL-4	1	2.2	22	88/656
hIL-5	1	2.2	22	90/586
hIL-6	1	1.7	170	89/548
hIL-8	1	1.8	180	89/520
hIL-10	1	0.8	4	93/722
hIL-12	1	0.8	8	95/544
hIL-17A	1	0.9	9000	01/420
hIFN-g	1	1.1	22	87/586
hTNF-a	1	0.9	36	87/650
mIL-2	1	3.1	310	93/566
mIL-4	1	3	30	91/656
mIL-6	1	8.5	850	93/730
mIFN-g*	1		4.5	Gg02-901-533
mTNF-a	1	1.7	340	88/532

\* Mouse IFN-g is calibrated using NIH standard (Lot Gg02-901-533) and is measured in Units (U)

ELISA Troubleshooting Guide		
Problem	Possibility	Solution
A. High background	1. Improper and inefficient washing	1. Improve efficiency of washing. Fill plates completely, soak for 1 minute per wash, as directed.
	2. Cross contamination from other specimens or positive control	2. Repeat ELISA being careful when washing and pipetting.
	3. Contaminated substrate	3. Substrate should be colorless. Replace.
	4. Incorrect dilutions, e.g., conjugate concentration was too high	4. Repeat using correct dilutions.
B. No signal	1. Improper, low protein binding capacity plates were used	1. Repeat ELISA using recommended high binding capacity plates.
	2. Wrong substrate was used	2. Repeat ELISA using the correct substrate
	3. Enzyme inhibitor present in buffers; e.g., sodium azide in the washing buffer and Assay Diluent inhibits peroxidase activity	3. Repeat ELISA making no enzyme inhibitor is present in any buffers.
	4. Coated capture antibody in ELISA/ELISPOT Diluent rather than Coating Buffer	4. Repeat ELISA using Coating Buffer contained in the set as the diluent for the capture antibody.

ELISA Troubleshooting Guide		
Problem	Possibility	Solution
C. Very weak signal	1. Improper and inefficient washing	1. Make sure washing procedure is done correctly, with a soak time.
	2. Incorrect dilutions of standard	2. Follow recommendations of standard preparation exactly as written on the C of A.
	3. Insufficient incubation time	3. Repeat ELISA following the protocol carefully for each step.
	4. Incorrect storage of reagents	4. Store reagents at the correct temperature as indicated on the Technical Data Sheet. Freezing certain components will severely impact results. Do not re-use the standards.
	5. Wrong filter in ELISA reader was used	5. Use the correct wavelength setting.
	6. Wrong plate used	6. Use the recommended Corning Costar 9018 or NUNC Maxisorp flat bottom 96 well plates.
D. Variation amongst replicates	1. Improper and inefficient washing	1. Make sure washing procedure is done correctly; see C of A. Edge effects can be avoided by moving samples and standards in from the edge of the plate.
	2. Poor mixing of samples	2. Mix samples and reagents gently and equilibrate to proper temperature.
	3. Plates not clean	3. Plates should be wiped on bottom before measuring absorbance
	4. Reagents have expired	4. Order a new Ready-Set-Go ELISA.

## Documentation and support

### Customer and technical support

Visit [thermofisher.com/support](http://thermofisher.com/support) for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
  - Product FAQs
  - Software, patches, and updates
- Order and web support
- Product documentation, including:
  - User guides, manuals, and protocols
  - Certificates of Analysis
  - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

### Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [thermofisher.com/support](http://thermofisher.com/support).

The information in this guide is subject to change without notice.

**DISCLAIMER:** TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

**Important Licensing Information:** These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

**Corporate entity:** Life Technologies | Carlsbad, CA 92008 USA | Toll Free in USA 1.800.955.6288

©2017 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are properties of their respective owners.

For support visit [thermofisher.com/support](http://thermofisher.com/support) or email [techsupport@lifetech.com](mailto:techsupport@lifetech.com)

[thermofisher.com](http://thermofisher.com)

23 January 2017

**ThermoFisher**  
SCIENTIFIC