



Csf1 (Mouse) ELISA Kit

Catalog Number KA0385

96 assays

Version: 03

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	5
Assay Protocol	6
Reagent Preparation	6
Sample Preparation	6
Assay Procedure	7
Data Analysis	9
Calculation of Results	9
Performance Characteristics	9
Resources	11
References	11
Plate Layout	12

Introduction

Intended Use

For quantitative detection of mouse M-CSF in sera, plasma, body fluids, tissue lysates or cell culture supernates.

Background

M-CSF, also called CSF1, consists of a 14-amino acid peptide, longer than the usual 8-to-11 mer recognized by most CTLs. The M-CSF gene is mapped to 1p21-p13 and contains 10 exons and 9 introns spanning 20 kb¹. Although it is a single-copy gene, its expression results in the synthesis of several mRNAs, ranging in size from about 1.5 to 4.5 kb². There are 2 forms of M-CSF, with 224 and 522 amino acids, resulting from alternative splicing³. Furthermore, M-CSF has a role in development of the placenta. Uterine M-CSF concentration is regulated by a synergistic action of estradiol and progesterone. M-CSF is produced by uterine glandular epithelial cells. It had been found that FMS, the M-CSF receptor, is expressed in placenta and choriocarcinoma cell lines⁴

Principle of the Assay

The mouse M-CSF ELISA Kit was based on standard sandwich enzyme-linked immune-sorbent assay technology. Mouse M-CSF specific polyclonal antibodies were precoated onto 96-well plates. The mouse specific detection polyclonal antibodies were biotinylated. The test samples and biotinylated detection antibodies were added to the wells subsequently and then followed by washing with PBS or TBS buffer. Avidin-Biotin-Peroxidase Complex was added and unbound conjugates were washed away with PBS or TBS buffer. HRP substrate TMB was used to visualize HRP enzymatic reaction. TMB was catalyzed by HRP to produce a blue color product that changed into yellow after adding acidic stop solution. The density of yellow is proportional to the mouse M-CSF amount of sample captured in plate.

General Information

Materials Supplied

List of component

Component	Amount
Lyophilized recombinant mouse M-CSF standard	10 ng/tube×2
One 96-well plate precoated with anti-mouse M-CSF antibody	1
Sample diluent buffer	30 ml
Biotinylated anti-mouse M-CSF antibody, dilution 1:100	130 µl
Antibody diluent buffer	12 ml
Avidin-Biotin-Peroxidase Complex (ABC), dilution 1:100	130 µl
ABC diluent buffer	12 ml
TMB color developing agent	10 ml
TMB stop solution	10 ml

Storage Instruction

Store at 4°C for frequent use, at -20°C for infrequent use. Avoid multiple freeze-thaw cycles (Shipped with wet ice.) Expiration: Four months at 4°C and eight months at -20°C.

Materials Required but Not Supplied

- ✓ Microplate reader in standard size.
- ✓ Automated plate washer.
- ✓ Adjustable pipettes and pipette tips. Multichannel pipettes are recommended in the condition of large amount of samples in the detection.
- ✓ Clean tubes and Eppendorf tubes.
- ✓ Washing buffer (neutral PBS or TBS).

Preparation of 0.01M TBS:

Add 1.2g Tris, 8.5g NaCl; 450µl of purified acetic acid or 700µl of concentrated hydrochloric acid to 1000ml H₂O and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

Preparation of 0.01 M PBS:

Add 8.5g sodium chloride, 1.4g Na₂HPO₄ and 0.2g NaH₂PO₄ to 1000ml distilled water and adjust pH to 7.2-7.6. Finally, adjust the total volume to 1L.

Precautions for Use

- Precautions
 - ✓ Before using Kit, spin tubes and bring down all components to bottom of tube.
 - ✓ Duplicate well assay was recommended for both standard and sample testing.
 - ✓ Don't let 96-well plate dry, dry plate will inactivate active components on plate.
 - ✓ In order to avoid marginal effect of plate incubation due to temperature difference (reaction may be stronger in the marginal wells), it is suggested that the diluted ABC and TMB solution will be pre-warmed in 37°C for 30 min before using.

Assay Protocol

Reagent Preparation

- Reconstitution of the mouse M-CSF standard: M-CSF standard solution should be prepared no more than 2 hours prior to the experiment. Two tubes of M-CSF standard (10ng per tube) are included in each kit. Use one tube for each experiment.
 - a. 10,000pg/ml of mouse M-CSF standard solution: Add 1 ml sample diluent buffer into one tube, keep the tube at room temperature for 10 min and mix thoroughly.
 - b. 1000pg/ml of mouse M-CSF standard solution: Add 0.1 ml of the above 10ng/ml M-CSF standard solution into 0.9 ml sample diluent buffer and mix thoroughly.
 - c. 500pg/ml→15.6pg/ml of mouse M-CSF standard solutions: Label 6 Eppendorf tubes with 500pg/ml, 250pg/ml, 125pg/ml, 62.5pg/ml, 31.3pg/ml, 15.6pg/ml, respectively. Aliquot 0.3 ml of the sample diluent buffer into each tube. Add 0.3 ml of the above 1000pg/ml M-CSF standard solution into 1st tube and mix. Transfer 0.3 ml from 1st tube to 2nd tube and mix. Transfer 0.3 ml from 2nd tube to 3rd tube and mix, and so on.

Note: The standard solutions are best used within 2 hours. The 10ng/ml standard solution may be stored at 4 °C for up to 12 hours, or at -20 °C for up to 48 hours. Avoid repeated freeze-thaw cycles.

- Preparation of biotinylated anti-mouse M-CSF antibody working solution: The solution should be prepared no more than 2 hours prior to the experiment.
 - a. The total volume should be: 0.1ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume)
 - b. Biotinylated anti-mouse M-CSF antibody should be diluted in 1:100 with the antibody diluent buffer and mixed thoroughly.
- Preparation of Avidin-Biotin-Peroxidase Complex (ABC) working solution: The solution should be prepared no more than 1 hour prior to the experiment.
 - a. The total volume should be: 0.1ml/well x (the number of wells). (Allowing 0.1-0.2 ml more than total volume)
 - b. Avidin- Biotin-Peroxidase Complex (ABC) should be diluted in 1:100 with the ABC dilution buffer and mixed thoroughly.

Sample Preparation

Store samples to be assayed within 24 hours at 2-8 °C. For long-term storage, aliquot and freeze samples at -20 °C. Avoid repeated freeze-thaw cycles.

- Cell culture supernate, tissue lysate or body fluids: Remove particulates by centrifugation, analyze immediately or aliquot and store at -20 °C
- Serum: Allow the serum to clot in a serum separator tube (about 2 hours) at room temperature. Centrifuge at approximately 2000 X g for 15 min. Analyze the serum immediately or aliquot and store

frozen at -20°C.

- Plasma: Use EDTA as an anticoagulant. Centrifuge for 15 min at 2000 x g within 30 min of collection. Analyze immediately or aliquot and store frozen at -20°C.

Sample Dilution Guideline

The user needs to estimate the concentration of the target protein in the sample and select a proper dilution factor so that the diluted target protein concentration falls near the middle of the linear regime in the standard curve. Dilute the sample using the provided diluent buffer. The following is a guideline for sample dilution. Several trials may be necessary in practice. The sample must be well mixed with the diluents buffer.

- High target protein concentration (10-100ng/ml). The working dilution is 1:100. i.e. Add 1 µl sample into 99 µl sample diluent buffer.
- Medium target protein concentration (1-10ng/ml). The working dilution is 1:10. i.e. Add 10 µl sample into 90 µl sample diluent buffer.
- Low target protein concentration (15.6-1000pg/ml). The working dilution is 1:2. i.e. Add 50 µl sample to 50 µl sample diluent buffer.
- Very Low target protein concentration (≤15.6pg/ml). No dilution necessary, or the working dilution is 1:2.

Assay Procedure

The ABC working solution and TMB color developing agent must be kept warm at 37°C for 30 min before use. When diluting samples and reagents, they must be mixed completely and evenly. Standard M-CSF detection curve should be prepared for each experiment. The user will decide sample dilution fold by crude estimation of M-CSF amount in samples.

1. Aliquot 0.1ml per well of the 1000pg/ml, 500pg/ml, 250pg/ml, 125pg/ml, 62.5pg/ml, 31.3pg/ml, 15.6pg/ml mouse M-CSF standard solutions into the precoated 96-well plate. Add 0.1ml of the sample diluent buffer into the control well (Zero well). Add 0.1ml of each properly diluted sample of mouse sera, plasma, body fluids, tissue lysates or cell culture supernatants to each empty well. See "Sample Dilution Guideline" above for details. We recommend that each mouse M-CSF standard solution and each sample is measured in duplicate.
2. Seal the plate with the cover and incubate at 37°C for 90 min.
3. Remove the cover, discard plate content, and blot the plate onto paper towels or other absorbent material. Do NOT let the wells completely dry at any time.
4. Add 0.1ml of biotinylated anti-mouse M-CSF antibody working solution into each well and incubate the plate at 37°C for 60 min.
5. Wash the plate three times with 0.01M TBS or 0.01M PBS, and each time let washing buffer stay in the wells for 1 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material.
6. Add 0.1ml of prepared ABC working solution into each well and incubate the plate at 37°C for 30 min.

7. Wash plate 5 times with 0.01M TBS or 0.01M PBS, and each time let washing buffer stay in the wells for 1-2 min. Discard the washing buffer and blot the plate onto paper towels or other absorbent material.
8. Add 90 µl of prepared TMB color developing agent into each well and incubate plate at 37°C in dark for 20-25 min (shades of blue can be seen in the wells with the four most concentrated mouse M-CSF standard solutions; the other wells show no obvious color).
9. Add 0.1ml of prepared TMB stop solution into each well. The color changes into yellow immediately.
10. Read the O.D. absorbance at 450nm in a microplate reader within 30 min after adding the stop solution.

Summary

1. Add samples and standards and incubate the plate at 37°C for 90 min. Do not wash.
2. Add biotinylated antibodies and incubate the plate at 37°C for 60 min. Wash plate 3 times with 0.01M TBS.
3. Add ABC working solution and incubate the plate at 37°C for 30 min. Wash plate 5 times with 0.01M TBS.
4. Add TMB color developing agent and incubate the plate at 37°C in dark for 25-30 min.
5. Add TMB stop solution and read.

Data Analysis

Calculation of Results

For calculation, (the relative O.D.450) = (the O.D.450 of each well) – (the O.D.450 of Zero well). The standard curve can be plotted as the relative O.D.450 of each standard solution (Y) vs. the respective concentration of the standard solution (X). The mouse M-CSF concentration of the samples can be interpolated from the standard curve.

Note: if the samples measured were diluted, multiply the dilution factor to the concentrations from interpolation to obtain the concentration before dilution.

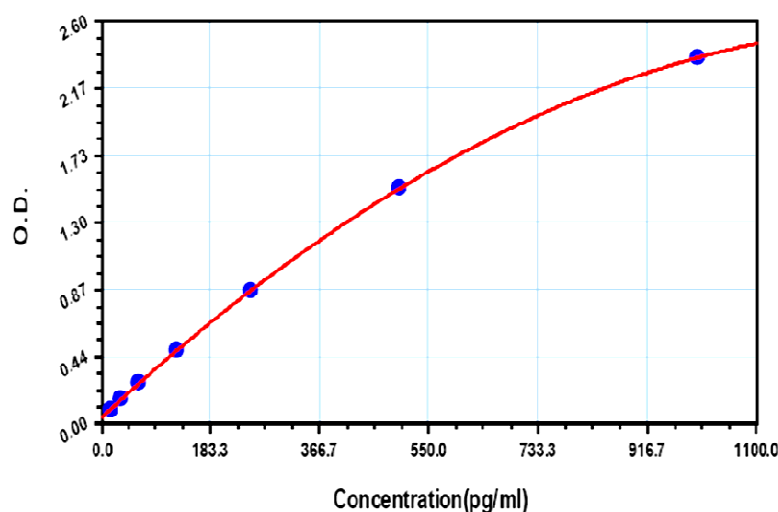
Typical result

Typical Data Obtained from Mouse M-CSF

Concentration	0.0pg/ml	15.6pg/ml	31.3pg/ml	62.5pg/ml	125pg/ml	250pg/ml	500pg/ml	1000pg/ml
O.D.	0.034	0.095	0.167	0.263	0.480	0.855	1.520	2.365

(TMB reaction incubate at 37 °C for 20 min)

Mouse M-CSF ELISA Kit



This standard curve was generated for demonstration purpose only. A standard curve must be run with each assay.

Performance Characteristics

1. Range
15.6pg/ml-1000pg/ml
2. Sensitivity
< 1 pg/ml

3. Specificity

No detectable cross-reactivity with any other cytokine.

Resources

References

1. Ladner, M. B.; Martin, G. A.; Noble, J. A.; Nikoloff, D. M.; Tal, R.; Kawasaki, E. S.; White, T. J. : Human CSF-1: gene structure and alternative splicing of mRNA precursors. EMBO J. 6: 2693-2698, 1987.
2. Kawasaki, E. S.; Ladner, M. B.; Wang, A. M.; Van Arsdell, J.; Warren, M. K.; Coyne, M. Y.; Schweickart, V. L.; Lee, M.-T.; Wilson, K. J.; Boosman, A.; Stanley, E. R.; Ralph, P.; Mark, D. F. : Molecular cloning of a complementary DNA encoding human macrophage-specific colony-stimulating factor (CSF-1). Science 230: 291-296, 1985.
3. Ladner, M. B.; Martin, G. A.; Noble, J. A.; Nikoloff, D. M.; Tal, R.; Kawasaki, E. S.; White, T. J. : Human CSF-1: gene structure and alternative splicing of mRNA precursors. EMBO J. 6: 2693-2698, 1987.
4. Pollard, J. W.; Bartocci, A.; Arceci, R.; Orlofsky, A.; Ladner, M. B.; Stanley, E. R. : Apparent role of the macrophage growth factor, CSF-1, in placental development. Nature 330: 484-486, 1987.

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H