ab100610 Human MMP9 ELISA Kit

For the quantitative measurement of Human MMP9 pro and active forms in serum, plasma, and cell culture supernatants. (Collect plasma using heparin as an anticoagulant. EDTA and Citrate are not recommended).

This product is for research use only and is not intended for diagnostic use.

Table of Contents

| 1. | Overview | 1 |
|-----|----------------------------------|----|
| 2. | Protocol Summary | 2 |
| 3. | Precautions | 3 |
| 4. | Storage and Stability | 3 |
| 5. | Limitations | 4 |
| 6. | Materials Supplied | 4 |
| 7. | Materials Required, Not Supplied | 5 |
| 8. | Technical Hints | 6 |
| 9. | Reagent Preparation | 7 |
| 10. | Standard Preparation | 8 |
| 11. | Sample Preparation | 9 |
| 12. | Plate Preparation | 11 |
| 13. | Assay Procedure | 12 |
| 14. | Calculations | 13 |
| 15. | Typical Data | 14 |
| 16. | Typical Sample Values | 15 |
| 17. | Assay Specificity | 16 |
| 18. | Troubleshooting | 17 |
| 19. | Notes | 18 |

1. Overview

Abcam's MMP9 Human ELISA (Enzyme-Linked Immunosorbent Assay) kit is an in vitro enzyme-linked immunosorbent assay for the quantitative measurement of Human MMP9 pro and active forms in serum, plasma (Collect plasma using heparin as an anticoagulant. EDTA and Citrate are not recommended), and cell culture supernatants.

This assay employs an antibody specific for Human MMP9 coated on a 96-well plate. Standards and samples are pipetted into the wells and MMP9 present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Human MMP9 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 MMP9 bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

ab100610 Human MMP9 ELISA Kit

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed

Add standard or sample to each well used. Incubate at room temperature

Add prepared biotin antibody to each well. Incubate at room temperature.

Add prepared Streptavidin solution. Incubate at room temperature.

Add TMB One-Step Development Solution to each well. Incubate at room temperature

Add Stop Solution to each well. Read at 450nm immediately.

3. Precautions

Please read these instructions carefully prior to beginning the assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances.
 However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth.
 Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

4. Storage and Stability

Store kit at -20°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

5. Limitations

- Assay kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

| Item | Amount | Storage Condition (Before Preparation) |
|-----------------------------------|----------|---|
| MMP9 Microplate (12 x 8 wells) | 96 wells | -20°C |
| 20X Wash Buffer Concentrate | 25 mL | -20°C |
| Recombinant Human MMP9 Standard | 2 vials | -20°C |
| 5X Assay Diluent | 15 mL | -20°C |
| Biotinylated anti-Human MMP9 | 2 vials | -20°C |
| 400X HRP-Streptavidin Concentrate | 200 µL | -20°C |
| TMB One-Step Substrate Reagent | 12 mL | -20°C |
| Stop Solution | 8 mL | -20°C |

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 2 µL to 1 mL volumes.
- Adjustable 1-25 mL pipettes for reagent preparation.
- 100 mL and 1 liter graduated cylinders.
- Absorbent paper.
- Distilled or deionized water.
- Log-log graph paper or computer and software for ELISA data analysis.
- Tubes to prepare standard or sample dilutions.

8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps.
- When preparing your standards, it is very critical to briefly spin down the vial first. The powder may drop off from the cap when opening it if you do not spin down. Be sure to dissolve the powder thoroughly when reconstituting. After adding Assay Diluent to the vial, we recommend inverting the tube a few times, then flick the tube a few times, and then spin it down; repeat this procedure 3-4 times. This is a technique we find very effective for thoroughly mixing the standard without too much mechanical force.
- Do not vortex the standard during reconstitution, as this will destabilize the protein.
- Once your standard has been reconstituted, it should be used right away or else frozen for later use.
- Keep the standard dilutions on ice while during preparation, but the ELISA procedure should be done at room temperature.
- Be sure to discard the working standard dilutions after use they do not store well.
- This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use.
 The kit contains enough reagents for 96 wells.
- Prepare only as much reagent as is needed on the day of the experiment.

9.1 1X Assay Diluent

5X Assay Diluent should be diluted 5-fold with deionized or distilled water before use.

9.2 1X Wash Solution

If the 20X Wash Concentrate contains visible crystals, equilibrate to room temperature and mix gently until dissolved. Dilute 20 mL of 20X Wash Buffer Concentrate into deionized or distilled water to yield 400 mL of 1X Wash Solution.

9.3 1X Biotinylated MMP9 Detection Antibody

Briefly spin the Biotinylated anti-Human MMP9 vial before use. Add 100 µL of 1X Assay Diluent into the vial to prepare a detection antibody concentrate. Pipette up and down to mix gently (the concentrate can either be stored at 4°C for 5 days or aliquoted and frozen at -20°C for 2 months). The detection antibody concentrate must be diluted 80-fold with 1X Assay Diluent prior to use in the Assay Procedure.

9.4 1X HRP-Streptavidin Solution

Briefly spin the 400X HRP-Streptavidin concentrate vial and pipette up and down to mix gently before use. HRP-Streptavidin concentrate must be diluted 400-fold with 1X Assay Diluent prior to use in the Assay Procedure.

For example: Briefly spin the vial and pipette up and down to mix gently. Add 30 µL of 400X HRP-Streptavidin concentrate into a tube with 12 mL 1X Assay Diluent to prepare a 400-fold diluted 1X HRP-Streptavidin solution (don't store the diluted solution for next day use). Mix well.

10. Standard Preparation

- Prepare serially diluted standards immediately prior to use. Always prepare a fresh set of standards for every use.
- Standard (recombinant protein) should be stored at -20°C or 80°C (recommended at -80°C) after reconstitution.
- 10.1 Reconstitute Briefly spin the vial of MMP9 Standard. Prepare the 50 ng/mL MMP9 **Stock Standard** by adding 400 μ L 1X Assay Diluent into the vial (see table below).
- 10.2 Dissolve the powder thoroughly by a gentle mix.
- **10.3** Label tubes #1-7.
- 10.4 Prepare Standard #1 by adding 80 µL 50 ng/mL Stock Standard to 586.7 µL Assay Diluent into tube #1. Mix thoroughly and gently.
- 10.5 Pipette 400 µL 1X Assay Diluent into remaining tubes.
- 10.6 Prepare **Standard #2** by transferring 200 μ L from tube #1 to #2, mix thoroughly.
- 10.7 Prepare **Standard #3** by transferring 200 μ L from tube #2 to #3, mix thoroughly.
- 10.8 Using the table below as a guide, prepare further serial dilutions
- 10.9 1X Assay Diluent serves as the zero standard (0 pg/mL).

Standard Dilution Preparation Table

| Standard # | Volume to Dilute (µL) | Diluent (µL) | Total Volume (µL) | Starting Conc. (pg/mL) | Final Conc. (pg/mL) |
|---------------|-----------------------------|-----------------|-------------------------|------------------------------|---------------------------|
| 1 | 80 | 586.7 | 666.7 | 50,000 | 6,000 |
| 2 | 200 | 400 | 600 | 6,000 | 2,000 |
| 3 | 200 | 400 | 600 | 2,000 | 666.7 |
| 4 | 200 | 400 | 600 | 666.7 | 222.2 |
| 5 | 200 | 400 | 600 | 222.2 | 74.07 |
| 6 | 200 | 400 | 600 | 74.07 | 24.69 |
| 7 | 200 | 400 | 600 | 24.69 | 8.23 |
| 8 | 0 | 400 | 400 | 0 | 0 |

11. Sample Preparation

General Sample Information:

- We recommend preparing serum-free or low-serum medium samples, as serum tends to contain cytokines which may produce significant background signals. If it is necessary to test serum containing medium, we recommend also running an uncultured media blank to assess baseline signals. This baseline can then be subtracted from the cultured media sample data.
- If your samples need to be diluted, 1X Assay Diluent should be used for dilution of serum, plasma and culture supernatants.
- Suggested dilution for normal serum/plasma: 300-3,000 fold.
- Please note that levels of the target protein may vary between different specimens. Optimal dilution factors for each sample must be determined by the investigator.

Cell culture supernatant preparation:

- 11.1 On day 0, seed ~1 million cells in 100 mm tissue culture plate with complete medium.*
- 11.2 On day 3, remove medium and replace medium with 6-8 mL of serum-free or low serum containing medium (e.g. medium containing 0.2% calf serum).
- 11.3 On day 5, collect medium into 15 mL tube. Centrifuge at 2,000 rpm in centrifuge at 4°C for 10 minutes. Save the supernatant. Transfer the supernatant into 1.5 mL Eppendorf tubes. Store supernatant at -80°C until experiment. Most samples can be stored this way for at least a year.

^{*}The optimal number of seeded cells varies from one cell type to another and may need to be empirically determined.

Refer to Dilution Guidelines for further instruction.

| Guidelines for Dilutions of 100-fold or Greater (for reference only; please follow the insert for specific dilution suggested) | | |
|--|--|--|
| 100x | 10000x | |
| 4 μl sample + 396 μl buffer (100X) = 100-fold dilution | A) 4 µl sample + 396 µl buffer (100X) B) 4 µl of A + 396 µl buffer (100X) = 10000-fold dilution | |
| Assuming the needed volume is less than or equal to 400 µl | Assuming the needed volume is less than or equal to 400 µl | |
| 1000x | 100000x | |
| A) 4 µl sample + 396 µl buffer (100X) B) 24 µl of A + 216 µl buffer (10X) = 1000-fold dilution | A) 4 µl sample + 396 µl buffer (100X) B) 4 µl of A + 396 µl buffer (100X) C) 24 µl of A + 216 µl buffer (10X) = 100000-fold dilution | |
| Assuming the needed volume is less than or equal to 240 µl | Assuming the needed volume is less than or equal to 240 µl | |

12. Plate Preparation

- The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- Unused well strips should be returned to the plate packet and stored at 4°C.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Well effects have not been observed with this assay. Contents of each well can be recorded on the template sheet included in the Resources section.

13. Assay Procedure

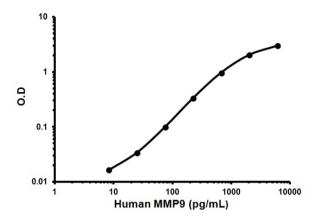
- Equilibrate all materials and prepared reagents to room temperature prior to use.
- We recommend that you assay all standards, controls and samples in duplicate.
- 13.1 Add 100 µL of each standard (see Standard Preparation section 10) and sample into appropriate wells. Cover well and incubate for 2.5 hours at room temperature or overnight at 4°C with gentle shaking.
- 13.2 Discard the solution and wash 4 times with 1X Wash Solution. Wash by filling each well with 1X Wash Solution (300 µL) using a multi-channel Pipette or auto washer. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining liquid by aspirating or decanting. Invert the plate and blot it against clean paper towels.
- 13.3 Add 100 µL of 1X Biotinylated MMP9 Detection Antibody (Standard Preparation section 10) to each well. Incubate for 1 hour at room temperature with gentle shaking.
- **13.4** Discard the solution. Repeat the wash as in step 13.2.
- 13.5 Add 100 μ L of 1X HRP-Streptavidin solution (see Standard Preparation section 10) to each well. Incubate for 45 minutes at room temperature with gentle shaking.
- **13.6** Discard the solution. Repeat the wash as in step 13.2.
- 13.7 Add 100 µL of TMB One-Step Substrate Reagent to each well.
- **13.8** Incubate for 30 minutes at room temperature in the dark with gentle shaking.
- 13.9 Add 50 μ L of Stop Solution to each well. Read at 450 nm immediately.

14. Calculations

Calculate the mean absorbance for each set of duplicate standards, controls and samples, and subtract the average zero standard optical density. Plot the standard curve on log-log graph paper, with standard concentration on the x-axis and absorbance on the y-axis. Draw the best-fit straight line through the standard points.

15. Typical Data

Typical standard curve – data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.



| | O.D. |
|------------------|------------------|
| Conc. (pg/mL) | Assay Diluent |
| 8.23 | 0.017 |
| 24.69 | 0.034 |
| 74.07 | 0.100 |
| 222.2 | 0.336 |
| 666.7 | 0.975 |
| 2,000 | 2.054 |
| 6,000 | 3.000 |

Figure 1. Example of human MMP standard curve.

16. Typical Sample Values

SENSITIVITY -

The minimum detectable dose of MMP9 is typically less than 10 pg/mL.

PRECISION -

| | Intra-Assay | Inter-Assay |
|--------|-------------|-------------|
| CV (%) | <10% | <12% |

RECOVERY -

Recovery was determined by spiking Human MMP9 into normal Human serum, plasma and cell culture media. Mean recoveries are as follows:

| Sample Type | Average % Recovery | Range (%) |
|--------------------|-----------------------|-----------|
| Serum | 96.23 | 84-103 |
| Plasma | 94.64 | 83-102 |
| Cell Culture Media | 95.38 | 84-104 |

Linearity of Dilution

| Serum Dilution | Average % Expected Value | Range (%) |
|----------------|--------------------------|-----------|
| 1:2 | 95 | 84-103 |
| 1:4 | 96 | 85-104 |

| Plasma Dilution | Average % Expected Value | Range (%) |
|-----------------|--------------------------|-----------|
| 1:2 | 93 | 83-102 |
| 1:4 | 97 | 86-105 |

| Cell Culture Media Dilution | Average % Expected Value | Range (%) |
|-----------------------------|--------------------------|-----------|
| 1:2 | 94 | 85-104 |
| 1:4 | 96 | 83-103 |

17. Assay Specificity

Cross Reactivity: This ELISA kit shows no cross-reactivity with the following cytokines tested: Human Angiogenin, BDNF, BLC, ENA-78, FGF-4, IL-1a, IL-1 β , IL-2, IL-3, IL-4, IL-5, IL-6, IL-7, IL-8, IL-9, IL-10, IL-11, IL-12 p70, IL-12 p40, IL-13, IL-15, IL-309, IP-10, G-CSF, GM-CSF, IFN- γ , Leptin (OB), MCP-1, MCP-3, MDC, MIP-1a, MIP-1 β , MIP-1 δ , MMP-1, -2, -3, -10, PARC, RANTES, SCF, TARC, TGF- β , TIMP-1, TIMP-2, TNF-a, TNF- β , TPO, VEGF).

Please contact our Technical Support team for more information.

18. Troubleshooting

| Problem | Cause | Solution |
|-------------------|---|---|
| Poor | Inaccurate pipetting | Check pipettes |
| standard curve | Improper standards dilution | Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing |
| Low Signal | Incubation times too brief | Ensure sufficient incubation times; change to overnight standard/sample incubation |
| Low Signal | Inadequate reagent volumes or improper dilution | Check pipettes and ensure correct preparation |
| Large CV | Plate is insufficiently washed | Review manual for proper wash technique. If using a plate washer, check all ports for obstructions |
| | Contaminated wash buffer | Prepare fresh wash buffer |
| Low sensitivity | Improper storage of the ELISA kit | Store the reconstituted protein at -80°C, all other assay components 4°C. Keep substrate solution protected from light. |

19. Notes

Technical Support

Copyright © 2017 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 / +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abcam.com | +64-(0)9-909-7829