

Human MMP-3 ELISA Kit

Enzyme-linked Immunosorbent Assay for quantitative detection of human MMP-3

Catalog Numbers BMS2014-3 and BMS2014-3TEN

Pub. No. MAN0016472 **Rev.** A.0 (30)

WARNING! Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. Safety Data Sheets (SDSs) are available from thermofisher.com/support.

Product description

The Human MMP-3 ELISA Kit is an enzyme-linked immunosorbent assay for the quantitative detection of human MMP-3.

Summary

The matrix metalloproteinases (MMPs) represent a family of more than 15 Zn-dependent endopeptidases with extracellular activity. They share a common conserved protease domain sequence in which three His residues form a complex with a catalytic Zn ion. These enzymes are responsible for the degradation and remodelling of extracellular matrix (ECM) components such as collagen, aggrecan, laminin, fibrilin, and fibronectin.

MMP-3, also known as matrix metalloproteinase-3 or stromelysin-1, is secreted as a zymogen by connective tissue cells that requires proteolytic cleavage of the attached pro-peptide domain for activation. Increased serum levels of MMP-3 are observed in inflammatory rheumatic diseases. MMP expression also correlates strongly with invasive behavior of tumor cells in culture and with metastasis in animal models. Overexpression of SL-1 can facilitate cell infiltration into surrounding tissue and degrade the ECM leading to induction of apoptosis.

For literature update refer to our website.

Principles of the test

An anti-human MMP-3 coating antibody is adsorbed onto microwells.

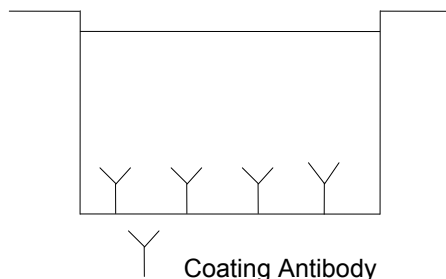


Fig. 1 Coated microwell.

Human MMP-3 present in the sample or standard binds to antibodies adsorbed to the microwells. A biotin-conjugated anti-human MMP-3 antibody is added and binds to human MMP-3 captured by the first antibody.

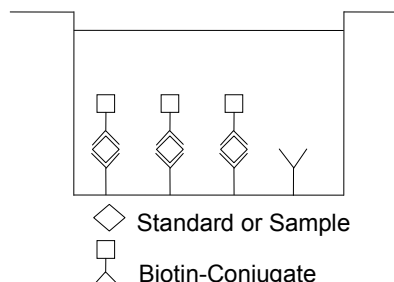


Fig. 2 First incubation.

Following incubation unbound biotin-conjugated anti-human MMP-3 antibody is removed during a wash step. Streptavidin-HRP is added and binds to the biotin-conjugated anti-human MMP-3 antibody.

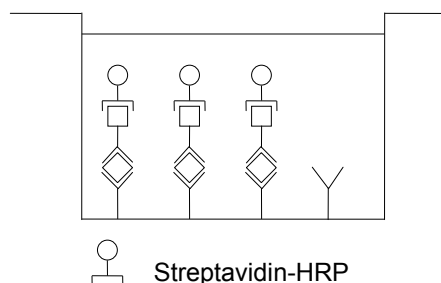


Fig. 3 Second incubation.

Following incubation unbound Streptavidin-HRP is removed during a wash step, and substrate solution reactive with HRP is added to the wells.

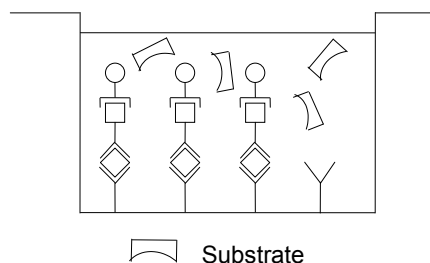


Fig. 4 Third incubation.

A colored product is formed in proportion to the amount of human MMP-3 present in the sample or standard. The reaction is terminated by addition of acid and absorbance is measured at 450 nm. A standard curve is prepared from 7 human MMP-3 standard dilutions and human MMP-3 sample concentration determined.

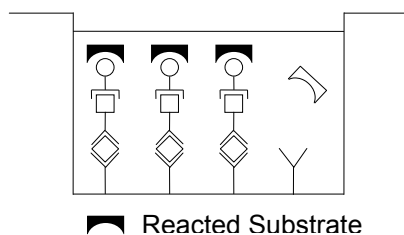


Fig. 5 Stop reaction.

Reagents provided

Reagents for human MMP-3 ELISA BMS2014-3 (96 tests)

1 aluminum pouch with a Microwell Plate (12 strips of 8 wells each) coated with polyclonal antibody to human MMP-3

1 vial (70 μ L) Biotin-Conjugate anti-human MMP-3 polyclonal antibody

1 vial (150 μ L) Streptavidin-HRP

2 vials human MMP-3 Standard lyophilized, 8 ng/mL upon reconstitution

1 vial (5 mL) Assay Buffer Concentrate 20x (PBS with 1% Tween[™] 20, 10% BSA)

1 bottle (50 mL) Wash Buffer Concentrate 20x (PBS with 1% Tween[™] 20)

1 bottle (12 mL) Sample Diluent

1 vial (15 mL) Substrate Solution (tetramethyl-benzidine)

1 vial (15 mL) Stop Solution (1M Phosphoric acid)

4 Adhesive Films

Reagents for human MMP-3 ELISA BMS2014-3TEN (10x96 tests)

10 aluminum pouches with a Microwell Plate (12 strips of 8 wells each) coated with polyclonal antibody to human MMP-3

10 vials (70 μ L) Biotin-Conjugate anti-human MMP-3 polyclonal antibody

10 vials (150 μ L) Streptavidin-HRP

10 vials human MMP-3 Standard lyophilized, 8 ng/mL upon reconstitution

2 vials (5 mL) Assay Buffer Concentrate 20x (PBS with 1% Tween[™] 20, 10% BSA)

6 bottles (50 mL) Wash Buffer Concentrate 20x 10 bottles (12 mL) Sample Diluent

10 vials (15 mL) Substrate Solution (tetramethyl-benzidine)

1 vial (100 mL) Stop Solution (1M Phosphoric acid)

20 Adhesive Films

Storage instructions – ELISA kit

Store kit reagents between 2°C and 8°C. Immediately after use remaining reagents should be returned to cold storage (2°C to 8°C). Expiry of the kit and reagents is stated on labels.

Expiry of the kit components can be guaranteed only if the components are stored properly, and if, in case of repeated use of one component, this reagent is not contaminated by the first handling.

Sample collection and storage instructions

Cell culture supernatant, serum, and plasma (citrate, heparin) were tested with this assay. Other biological samples might be suitable for use in the assay. Remove serum or plasma from the clot or cells as soon as possible after clotting and separation.

Pay attention to a possible *Hook Effect* due to high sample concentrations (see “Calculation of results” on page 4).

Samples containing a visible precipitate must be clarified prior to use in the assay. Do not use grossly hemolyzed or lipemic samples.

Samples should be aliquoted and must be stored frozen at –20°C to avoid loss of bioactive human MMP-3. If samples are to be run within 24 hours, they may be stored at 2°C to 8°C (refer to “Sample stability” on page 6).

Avoid repeated freeze-thaw cycles. Prior to assay, the frozen sample should be brought to room temperature slowly and mixed gently.

Materials required but not provided

- 5 mL and 10 mL graduated pipettes
- 5 μ L to 1000 μ L adjustable single channel micropipettes with disposable tips
- 50 μ L to 300 μ L adjustable multichannel micropipette with disposable tips
- Multichannel micropipette reservoir
- Beakers, flasks, and cylinders necessary for preparation of reagents
- Device for delivery of wash solution (multichannel wash bottle or automatic wash system)
- Microwell strip reader capable of reading at 450 nm (620 nm as optional reference wave length)
- Glass-distilled or deionized water
- Statistical calculator with program to perform regression analysis

Precautions for use

- All reagents should be considered as potentially hazardous. We therefore recommend that this product is handled only by those persons who have been trained in laboratory techniques and that it is used in accordance with the principles of good laboratory practice. Wear suitable protective clothing such as laboratory overalls, safety glasses and gloves. Care should be taken to avoid contact with skin or eyes. In the case of contact with skin or eyes wash immediately with water. See material safety data sheet(s) and/or safety statement(s) for specific advice.
- Reagents are intended for research use only and are not for use in diagnostic or therapeutic procedures.
- Do not mix or substitute reagents with those from other lots or other sources.
- Do not use kit reagents beyond expiration date on label.
- Do not expose kit reagents to strong light during storage or incubation.
- Do not pipet by mouth.
- Do not eat or smoke in areas where kit reagents or samples are handled.
- Avoid contact of skin or mucous membranes with kit reagents or samples.
- Rubber or disposable latex gloves should be worn while handling kit reagents or samples.
- Avoid contact of substrate solution with oxidizing agents and metal.
- Avoid splashing or generation of aerosols.
- To avoid microbial contamination or cross-contamination of reagents or samples that may invalidate the test, use disposable pipette tips and/or pipettes.
- Use clean, dedicated reagent trays for dispensing the conjugate and substrate reagent.
- Exposure to acid inactivates the conjugate.
- Glass-distilled water or deionized water must be used for reagent preparation.
- Substrate solution must be at room temperature prior to use.
- Decontaminate and dispose samples and all potentially contaminated materials as if they could contain infectious agents. The preferred method of decontamination is autoclaving for a minimum of 1 hour at 121.5°C.
- Liquid wastes not containing acid and neutralized waste may be mixed with sodium hypochlorite in volumes such that the final mixture contains 1.0% sodium hypochlorite. Allow 30 minutes for effective decontamination. Liquid waste containing acid must be neutralized prior to the addition of sodium hypochlorite.

Preparation of reagents

1. Buffer Concentrates should be brought to room temperature and should be diluted before starting the test procedure.
2. If crystals have formed in the Buffer Concentrates, warm them gently until they have completely dissolved.

Wash buffer (1x)

1. Pour entire contents (50 mL) of the Wash Buffer Concentrate (20x) into a clean 1,000 mL graduated cylinder. Bring to final volume of 1,000 mL with glass-distilled or deionized water. Mix gently to avoid foaming.
2. Transfer to a clean wash bottle and store at 2°C to 25°C. The Wash Buffer (1x) is stable for 30 days.
3. Wash Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Wash Buffer Concentrate (20x) (mL)	Distilled Water (mL)
1–6	25	475
1–12	50	950

Assay buffer (1x)

1. Pour the entire contents (5 mL) of the Assay Buffer Concentrate (20x) into a clean 100 mL graduated cylinder. Bring to final volume of 100 mL with distilled water. Mix gently to avoid foaming.
2. Store at 2°C to 8°C. The Assay Buffer (1x) is stable for 30 days.
3. Assay Buffer (1x) may also be prepared as needed according to the following table:

Number of Strips	Assay Buffer Concentrate (20x) (mL)	Distilled Water (mL)
1–6	2.5	47.5
1–12	5.0	95.0

Biotin-Conjugate

Note: The Biotin-Conjugate should be used within 30 minutes after dilution.

Make a 1:100 dilution of the concentrated Biotin-Conjugate solution with Assay Buffer (1x) in a clean plastic tube as needed according to the following table:

Number of Strips	Biotin-Conjugate (mL)	Assay Buffer (1x) (mL)
1–6	0.03	2.97
1–12	0.06	5.94

Streptavidin-HRP

Note: The Streptavidin-HRP should be used within 30 minutes after dilution.

Make a 1:100 dilution of the concentrated Streptavidin-HRP solution with Sample Diluent in a clean plastic tube as needed according to the following table:

Number of Strips	Streptavidin-HRP (mL)	Sample Diluent (mL)
1–6	0.06	5.94
1–12	0.12	11.88

Human MMP-3 standard

1. Reconstitute human MMP-3 Standard by addition of distilled water. Reconstitution volume is stated on the label of the standard vial. Swirl or mix gently to insure complete and homogeneous solubilization (concentration of reconstituted standard = 8 ng/mL).
2. Allow the standard to reconstitute for 10–30 minutes. Mix well prior to making dilutions.

The standard has to be used immediately after reconstitution and cannot be stored.

External standard dilution

1. Label 7 tubes, one for each standard point: S1, S2, S3, S4, S5, S6, S7.
2. Prepare 1:2 serial dilutions for the standard curve as follows: Pipette 225 µL of Assay Buffer (1x) into each tube.
3. Pipette 225 µL of reconstituted standard (concentration of standard = 8 ng/mL) into the first tube, labeled S1, and mix (concentration of standard 1 = 4 ng/mL).
4. Pipette 225 µL of this dilution into the second tube, labeled S2, and mix thoroughly before the next transfer.
5. Repeat serial dilutions 5 more times thus creating the points of the standard curve (see Figure 6).

Assay Buffer (1x) serves as blank.

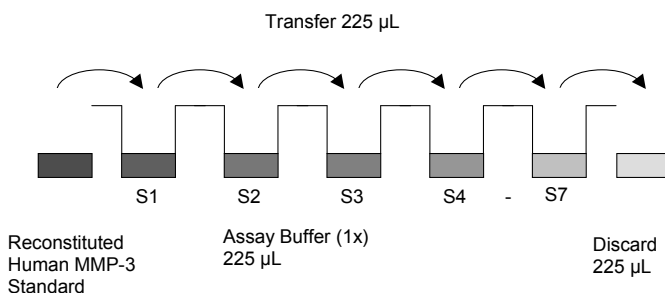


Fig. 6 Dilute standards - tubes.

Test protocol

Note: If instructions of this protocol have been followed, samples have been diluted 1:10 and the concentration read from the standard curve must be multiplied by the dilution factor (x 10).

Shaking is absolutely necessary for an optimal test performance.

1. Determine the number of microwell strips required to test the desired number of samples plus appropriate number of wells needed for running blanks and standards. Each sample, standard and blank should be assayed in duplicate. Remove extra microwell strips from holder and store in foil bag with the desiccant provided at 2°C to 8°C sealed tightly.
2. Wash the microwell strips twice with approximately 400 µL Wash Buffer per well with thorough aspiration of microwell contents between washes. Allow the Wash Buffer to sit in the wells for about 10–15 seconds before aspiration. Take care not to scratch the surface of the microwells.

After the last wash step, empty wells and tap microwell strips on absorbent pad or paper towel to remove excess Wash Buffer. Use the microwell strips immediately after washing. Alternatively, microwell strips can be placed upside down on a wet absorbent paper for not longer than 15 minutes. Do not allow wells to dry.

3. Standard dilution on the microwell plate (alternatively, the standard dilution can be prepared in tubes, see “External standard dilution” on page 3):

Add 100 µL of Assay Buffer (1x) in duplicate to all standard wells. Pipet 100 µL of prepared standard (see “Human MMP-3 standard” on page 3, concentration = 8 ng/mL) in duplicate into well A1 and A2 (see Table 1). Mix the contents of wells A1 and A2 by repeated aspiration and ejection (concentration of standard 1, S1 = 4 ng/mL), and transfer 100 µL to wells B1 and B2, respectively (see Figure 7). Take care not to scratch the inner surface of the microwells. Continue this procedure 5 times, creating two rows of human MMP-3 standard dilutions ranging from 4.0 to 0.06 ng/mL. Discard 100 µL of the contents from the last microwells (G1, G2) used.

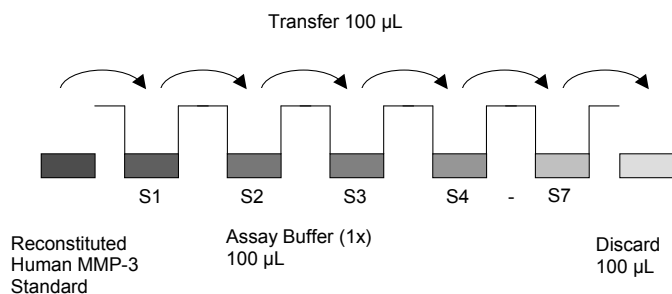


Fig. 7 Dilute standards - microwell plate.

In case of an external standard dilution (see “External standard dilution” on page 3), pipet 100 µL of these standard dilutions (S1–S7) in the standard wells according to Table 1.

Table 1 Example of the arrangement of blanks, standards, and samples in the microwell strips.

	1	2	3	4
A	Standard 1 4.0 ng/mL	Standard 1 4.0 ng/mL	Sample 1	Sample 1
B	Standard 2 2.0 ng/mL	Standard 2 2.0 ng/mL	Sample 2	Sample 2
C	Standard 3 1.0 ng/mL	Standard 3 1.0 ng/mL	Sample 3	Sample 3
D	Standard 4 0.50 ng/mL	Standard 4 0.50 ng/mL	Sample 4	Sample 4
E	Standard 5 0.25 ng/mL	Standard 5 0.25 ng/mL	Sample 5	Sample 5
F	Standard 6 0.13 ng/mL	Standard 6 0.13 ng/mL	Sample 6	Sample 6
G	Standard 7 0.06 ng/mL	Standard 7 0.06 ng/mL	Sample 7	Sample 7
H	Blank	Blank	Sample 8	Sample 8

4. Add 100 µL of Assay Buffer (1x) in duplicate to the blank wells.
5. Add 90 µL of Assay Buffer (1x) to the sample wells.
6. Add 10 µL of each sample in duplicate to the sample wells.
7. Prepare Biotin-Conjugate (see “Biotin-Conjugate” on page 3).
8. Add 50 µL of Biotin-Conjugate to all wells.
9. Cover with an adhesive film and incubate at room temperature (18°C to 25°C) for 2 hours on a microplate shaker.
10. Prepare Streptavidin-HRP (see “Streptavidin-HRP” on page 3).
11. Remove adhesive film and empty wells. Wash microwell strips 6 times according to step 2. Proceed immediately to the next step.
12. Add 100 µL of diluted Streptavidin-HRP to all wells, including the blank wells.
13. Cover with an adhesive film and incubate at room temperature (18°C to 25°C) for 1 hour on a microplate shaker.
14. Remove adhesive film and empty wells. Wash microwell strips 6 times according to step 2. Proceed immediately to the next step.
15. Pipet 100 µL of TMB Substrate Solution to all wells.
16. Incubate the microwell strips at room temperature (18°C to 25°C) for about 30 minutes. Avoid direct exposure to intense light.

The color development on the plate should be monitored and the substrate reaction stopped (see next step of this protocol) before positive wells are no longer properly recordable. Determination of the ideal time period for color development has to be done individually for each assay.

It is recommended to add the stop solution when the highest standard has developed a dark blue color. Alternatively the color development can be monitored by the ELISA reader at 620 nm. The substrate reaction should be stopped as soon as Standard 1 has reached an OD of 0.9–0.95.

17. Stop the enzyme reaction by quickly pipetting 100 µL of Stop Solution into each well. It is important that the Stop Solution is spread quickly and uniformly throughout the microwells to completely inactivate the enzyme. Results must be read immediately after the Stop Solution is added or within one hour if the microwell strips are stored at 2°C to 8°C in the dark.
18. Read absorbance of each microwell on a spectro-photometer using 450 nm as the primary wave length (optionally 620 nm as the reference wave length; 610 nm to 650 nm is acceptable). Blank the plate reader according to the manufacturer's instructions by using the blank wells. Determine the absorbance of both the samples and the standards.

Calculation of results

- Calculate the average absorbance values for each set of duplicate standards and samples. Duplicates should be within 20% of the mean value.
- Create a standard curve by plotting the mean absorbance for each standard concentration on the ordinate against the human MMP-3 concentration on the abscissa. Draw a best fit curve through the points of the graph (a 5-parameter curve fit is recommended).
- To determine the concentration of circulating human MMP-3 for each sample, first find the mean absorbance value on the ordinate and extend a horizontal line to the standard curve. At the point of intersection, extend a vertical line to the abscissa and read the corresponding human MMP-3 concentration.
- If instructions of this protocol have been followed, samples have been diluted 1:10 and the concentration read from the standard curve must be multiplied by the dilution factor (x 10).
- Calculation of samples with a concentration exceeding standard 1 will result in incorrect, low human MMP-3 levels (Hook Effect). Such samples require further external predilution according to expected human MMP-3 values with Assay Buffer (1x) in order to precisely quantitate the actual human MMP-3 level.
- It is suggested that each testing facility establishes a control sample of known human MMP-3 concentration and runs this additional control with each assay. If the values obtained are not within the expected range of the control, the assay results may be invalid.

- A representative standard curve is shown in Figure 8.
Note: Do not use this standard curve to derive test results. Each laboratory must prepare a standard curve for each group of microwell strips assayed.

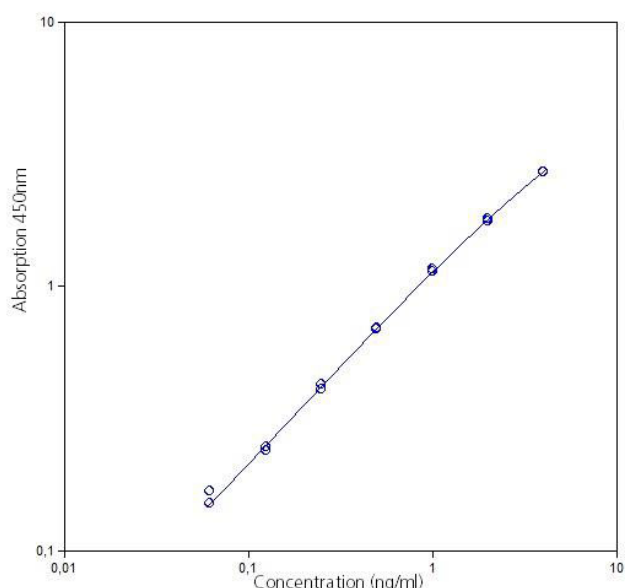


Fig. 8 Representative standard curve for Human MMP-3 ELISA Kit. Human MMP-3 was diluted in serial 2-fold steps in Assay Buffer (1x).

Table 2 Typical data using the Human MMP-3 ELISA Kit.

Measuring wavelength: 450 nm

Reference wavelength: 620 nm

Standard	Human MMP-3 concentration (ng/mL)	O.D. at 450 nm	Mean O.D. at 450 nm	C.V. (%)
1	4.00	2.697 2.689	2.647	0.1
2	2.00	1.785 1.752	1.722	0.9
3	1.00	1.150 1.132	1.094	0.8
4	0.50	0.690 0.681	0.639	0.7
5	0.25	0.421 0.404	0.366	2.0
6	0.13	0.246 0.237	0.195	1.9
7	0.06	0.150 0.167	0.112	5.5
Blank	0.00	0.047 0.046	0.47	1.4

The OD values of the standard curve may vary according to the conditions of assay performance (e.g., operator, pipetting technique, washing technique, or temperature effects). Furthermore, shelf life of the kit may affect enzymatic activity and thus color intensity. Values measured are still valid.

Limitations

- Because exact conditions may vary from assay to assay, a standard curve must be established for every run.
- Bacterial or fungal contamination of either screen samples or reagents or cross-contamination between reagents may cause erroneous results.

- Disposable pipette tips, flasks, or glassware are preferred, reusable glassware must be washed and thoroughly rinsed of all detergents before use.
- Improper or insufficient washing at any stage of the procedure will result in either false positive or false negative results. Empty wells completely before dispensing fresh wash solution, fill with Wash Buffer as indicated for each wash cycle, and do not allow wells to sit uncovered or dry for extended periods.

Performance characteristics

Sensitivity

The limit of detection of human MMP-3 defined as the analyte concentration resulting in an absorbance significantly higher than that of the dilution medium (mean plus 2 standard deviations) was determined to be 0.0005 ng/mL (mean of 3 independent assays).

Reproducibility

Intra-assay

Reproducibility within the assay was evaluated in 3 independent experiments. Each assay was carried out with 6 replicates of 8 serum samples containing different concentrations of human MMP-3. Two standard curves were run on each plate. Data below show the mean human MMP-3 concentration and the coefficient of variation for each sample. The calculated overall intra-assay coefficient of variation was 6.1%.

Table 3 The mean human MMP-3 concentration and the coefficient of variation for each sample.

Sample	Experiment	Mean concentration (ng/mL)	Coefficient of variation (%)
1	1	8.4	3.1
	2	8.8	6.2
	3	8.5	3.1
2	1	19.9	6.0
	2	18.5	6.9
	3	19.6	4.6
3	1	38.0	6.9
	2	38.1	5.8
	3	39.0	4.3
4	1	12.9	6.4
	2	11.9	2.9
	3	12.1	9.4
5	1	7.1	5.7
	2	6.4	6.4
	3	6.6	6.2
6	1	10.6	8.2
	2	10.6	9.3
	3	11.0	8.4
7	1	6.4	4.7
	2	5.7	6.4
	3	6.0	3.2
8	1	4.5	8.6
	2	4.1	10.0
	3	4.0	4.0

Inter-assay

Assay-to-assay reproducibility within one laboratory was evaluated in 3 independent experiments. Each assay was carried out with 6 replicates of 8 serum and plasma samples containing different concentrations of human MMP-3. Two standard curves were run on each plate. Data below show the mean human MMP-3 concentration and the coefficient of variation calculated on 18 determinations of each

sample. The calculated overall inter-assay coefficient of variation was 4.0%.

Table 4 The mean human MMP-3 concentration and the coefficient of variation of each sample.

Sample	Mean concentration (ng/mL)	Coefficient of variation (%)
1	8.6	2.1
2	19.3	3.7
3	38.3	1.4
4	12.3	4.3
5	6.7	5.5
6	10.7	2.1
7	6.1	6.0
8	4.2	7.1

Spike recovery

The spike recovery was evaluated by spiking 2 levels of human MMP-3 into serum, plasma (heparin, citrate), and cell culture supernatant. Recoveries were determined with 2 replicates each. The amount of endogenous human MMP-3 in unspiked samples was subtracted from the spike values.

Sample matrix	Spike medium mean (%)	Spike low mean (%)
Serum	78	93
Plasma (citrate)	96	93
Plasma (heparin)	78	99
Cell culture supernatant	115	95

Dilution parallelism

Serum and plasma (citrate, heparin) samples with different levels of human MMP-3 were analyzed at serial 2-fold dilutions with 4 replicates each.

Sample matrix	Recovery of exp. val.	
	Dilution	Mean (%)
Serum	1:20	103
	1:40	118
	1:80	90
Plasma (citrate)	1:20	104
	1:40	106
	1:80	109
Plasma (heparin)	1:20	122
	1:40	104
	1:80	105

Sample stability

Freeze-thaw stability

Aliquots of serum, plasma, and cell culture supernatant samples (spiked or unspiked) were stored at -20°C and thawed 3 times, and the human MMP-3 levels determined.

There was no significant loss of human MMP-3 immunoreactivity detected by freezing and thawing.

Storage stability

Aliquots of serum, plasma, and cell culture supernatant samples (spiked or unspiked) were stored at -20°C, 2°C to 8°C, and room temperature, and the human MMP-3 level determined after 24 hours.

There was no significant loss of human MMP-3 immunoreactivity detected during storage under above conditions.

Specificity

The cross-reactivity and interference of circulating factors of the immune system was evaluated by spiking these proteins at physiologically relevant concentrations into a human MMP-3 positive sample. No cross-reactivity or interference was detected.

Expected values

Forty serum and plasma (citrate, heparin) samples from apparently healthy donors were tested for human MMP-3. The detected human MMP-3 levels ranged between 2.0 and 28 ng/mL for serum, between 1.2 and 28.2 ng/mL for plasma (citrate), and between 4.5 and 29.5 ng/mL for plasma (heparin). Elevated human MMP-3 levels depend on the type of immunological disorder.

Reagent preparation summary

Wash buffer (1x)

Add Wash Buffer Concentrate 20x (50 mL) to 950 mL distilled water.

Number of Strips	Wash Buffer Concentrate (mL)	Distilled Water (mL)
1–6	25	475
1–12	50	950

Assay buffer (1x)

Add Assay Buffer Concentrate 20x (5 mL) to 95 mL distilled water.

Number of Strips	Assay Buffer Concentrate (mL)	Distilled Water (mL)
1–6	2.5	47.5
1–12	5.0	95.0

Biotin-Conjugate

Make a 1:100 dilution of Biotin-Conjugate in Assay Buffer (1X).

Number of Strips	Biotin-Conjugate (mL)	Assay Buffer (1x) (mL)
1–6	0.03	2.97
1–12	0.06	5.94

Streptavidin-HRP

Make a 1:100 dilution of Streptavidin-HRP in Sample Diluent.

Number of Strips	Streptavidin-HRP (mL)	Sample Diluent (mL)
1–6	0.06	5.94
1–12	0.12	11.88

Human MMP-3 standard

Reconstitute lyophilized human MMP-3 standard with distilled water. (Reconstitution volume is stated on the label of the standard vial.)

Test protocol summary

Note: If instructions of this protocol have been followed, samples have been diluted 1:10 and the concentration read from the standard curve must be multiplied by the dilution factor (x 10).

Shaking is absolutely necessary for an optimal test performance.

1. Determine the number of microwell strips required.
2. Wash microwell strips twice with Wash Buffer.
3. Standard dilution on the microwell plate: Add 100 µL Assay Buffer (1x), in duplicate, to all standard wells. Pipet 100 µL prepared standard into the first wells and create standard dilutions by transferring 100 µL from well to well. Discard 100 µL from the last wells.

Alternatively, external standard dilution in tubes (see “External standard dilution” on page 3): Pipet 100 µL of these standard dilutions in the microwell strips.

4. Add 100 µL Assay Buffer (1x), in duplicate, to the blank wells.
5. Add 90 µL Assay Buffer (1x) to sample wells.
6. Add 10 µL sample in duplicate, to designated sample wells.
7. Prepare Biotin-Conjugate.
8. Add 50 µL Biotin-Conjugate to all wells.
9. Cover microwell strips and incubate 2 hours at room temperature (18°C to 25°C) on a microplate shaker.
10. Prepare Streptavidin-HRP.
11. Empty and wash microwell strips 6 times with Wash Buffer.

12. Add 100 µL diluted Streptavidin-HRP to all wells.
13. Cover microwell strips and incubate 1 hour at room temperature (18°C to 25°C) on a microplate shaker.
14. Empty and wash microwell strips 6 times with Wash Buffer.
15. Add 100 µL of TMB Substrate Solution to all wells.
16. Incubate the microwell strips for about 30 minutes at room temperature (18°C to 25°C).
17. Add 100 µL Stop Solution to all wells.
18. Blank microwell reader and measure color intensity at 450 nm.

Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs

- Software, patches, and updates
- Training for many applications and instruments
- Order and web support
- Product documentation
 - 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 at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



Bender MedSystems GmbH | Campus Vienna Biocenter 2 | 1030 Vienna, Austria

For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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.

©2019 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 the property of their respective owners.