

Human MICA ELISA

Enzyme-linked Immunosorbent Assay for quantitative detection of human MICA

Catalog Numbers BMS2302 and BMS2302TEN

Pub. No. MAN0017640 **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 MICA ELISA is an enzyme-linked immunosorbent assay for the quantitative detection of human MICA. Cell lysates, cell culture supernatant, serum and plasma (citrate, heparin, EDTA) were tested with this assay. Other biological samples might be suitable for use in the assay.

Major histocompatibility complex (MHC) class I proteins are ubiquitously expressed and mediate the recognition of intracellular antigens by cytotoxic T cells. A related family, termed the MHC class I chain-related (MIC) proteins are recognized by NKG2D, a receptor on NK and T cells, and promote anti-tumor activity. MICA, a member of the MIC family, is widely expressed on many tumors, and it is the MICA/NKG2D interaction that is thought to stimulate the anti-tumor reactivity by T lymphocytes. Both MICA and MICB mRNA are widely expressed in normal tissues, with MICA being present in virtually every tissue except the nervous system, suggesting that MIC protein expression may only be one component of the anti-tumor activity of the immune system.

Principles of the test

An anti-human MICA coating antibody is adsorbed onto microwells.

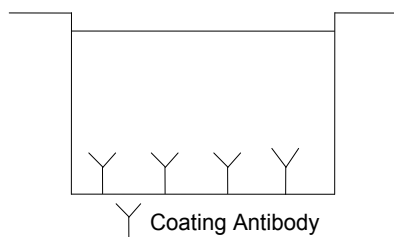


Fig. 1 Coated microwell

Human MICA present in the sample or standard binds to antibodies adsorbed to the microwells.

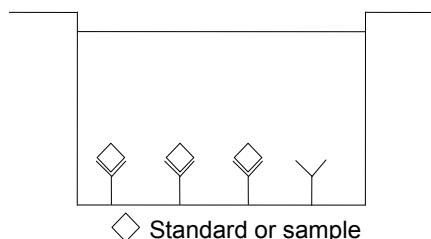


Fig. 2 First incubation

Following incubation, unbound standard or sample is removed during a wash step. A biotin-conjugated anti-human MICA antibody is added and binds to MICA captured by the first antibody.

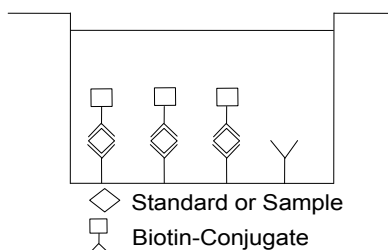


Fig. 3 Second incubation

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

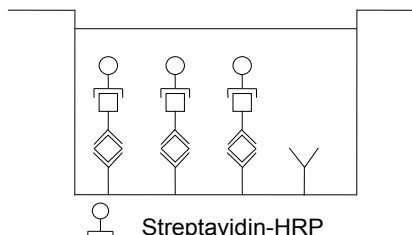


Fig. 4 Third incubation

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

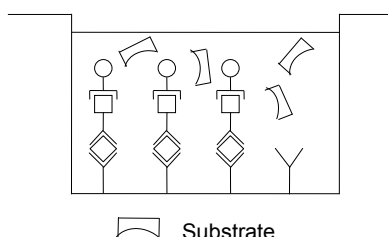


Fig. 5 Fourth incubation

A coloured product is formed in proportion to the amount of MICA 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 MICA standard dilutions and MICA sample concentration determined.

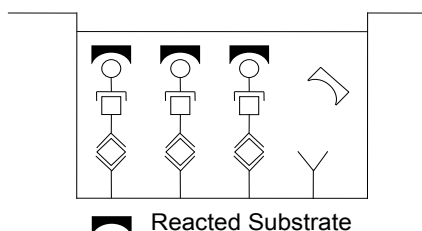


Fig. 6 Stop reaction

Reagents provided

Reagents for human MICA ELISA BMS2302 (96 tests)

1 aluminum pouch with a Microwell Plate coated with monoclonal antibody to human MICA

1 vial (120 µL) Biotin-Conjugate anti-human MICA monoclonal antibody

1 vial (150 µL) Streptavidin-HRP

2 vials human MICA Standard lyophilized, 200 pg/ml upon reconstitution

1 bottle (12 mL) Sample Diluent

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

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

6 Adhesive Films

Reagents for human MICA ELISA BMS2302TEN (10 x 96 tests)

10 aluminum pouches with a Microwell Plate coated with monoclonal antibody to human MICA

10 vials (120 µL) Biotin-Conjugate anti-human MICA monoclonal antibody to human MICA

10 vials (150 µL) Streptavidin-HRP

10 vials human MICA Standard lyophilized, 200 pg/mL upon reconstitution

10 vials (12 mL) Sample Diluent

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

6 bottles (50 mL) Wash Buffer Concentrate 20x (PBS with 1% Tween™ 20)

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

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

30 Adhesive Films

Storage instructions – ELISA kit

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

Expiry of the kit components can only be guaranteed 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, plasma (citrate, heparin, EDTA) were tested with this assay. Other biological samples might be suitable for use in the assay.

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 MICA. If samples are to be run within 24 hours, they may be stored at 2–8°C

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 µL to 1000 µL adjustable single channel micropipettes with disposable tips
- 50 µL to 300 µL adjustable multichannel micropipette with disposable tips
- Multichannel micropipette reservoir
- Beakers, flasks, 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 chemicals 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 1000 mL graduated cylinder. Bring to final volume of 1000 mL with glass-distilled or deionized water. Mix gently to avoid foaming.
2. Transfer to a clean wash bottle and store at 2° to 25°C. Please note that 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 2–8°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	0.06	5.94
1 - 12	0.12	11.88

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 Assay Buffer (1x) in a clean plastic tube as needed according to the following table:

Number of Strips	Streptavidin-HRP (mL)	Assay Buffer (1x) (mL)
1 - 6	0.06	5.94
1 - 12	0.12	11.88

Human MICA standard

1. Reconstitute human MICA standard by addition of distilled water.
2. 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 = 200 pg/mL).
3. Allow the standard to reconstitute for 10-30 minutes. Mix well prior to making dilutions.
4. After usage remaining standard cannot be stored and has to be discarded.
5. Standard dilutions can be prepared directly on the microwell plate (see "Test protocol" on page 4) or alternatively in tubes (see "External standard dilution" on page 3).

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 Sample Diluent into each tube.
3. Pipette 225 µL of reconstituted standard (concentration = 200 pg/mL) into the first tube, labeled S1, and mix (concentration of standard 1 = 100 pg/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 7).

Sample Diluent serves as blank.

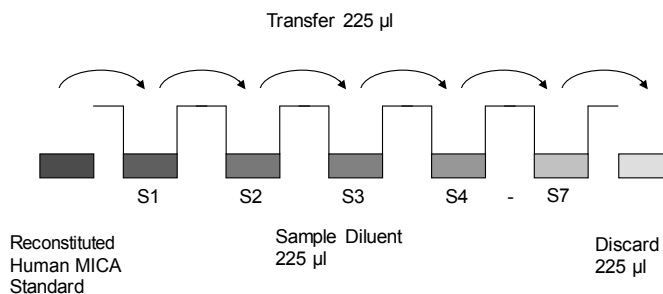


Fig. 7 Dilute standards - tubes

Test protocol

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, blank and optional control sample should be assayed in duplicate. Remove extra microwell strips from holder and store in foil bag with the desiccant provided at 2°-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 Sample Diluent in duplicate to all standard wells. Pipette 100 µL of prepared standard (see "Human MICA standard" on page 3, concentration = 200 pg/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 = 100 pg/mL), and transfer 100 µL to wells B1 and B2, respectively (see Figure 8). Take care not to scratch the inner surface of the microwells. Continue this procedure 5 times, creating two rows of human MICA standard dilutions ranging from 100 to 2 pg/mL. Discard 100 µL of the contents from the last microwells (S7) used.

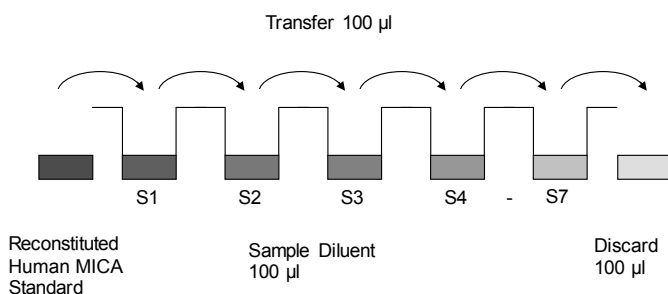


Fig. 8 Dilute standards - microwell plate

Table 1 Table depicting an example of the arrangement of blanks, standards and samples in the microwell strips:

	1	2	3	4
A	Standard 1 (100.0 pg/mL)	Standard 1 (100.0 pg/mL)	Sample 1	Sample 1
B	Standard 2 (50.0 pg/mL)	Standard 2 (50.0 pg/mL)	Sample 2	Sample 2
C	Standard 3 (25.0 pg/mL)	Standard 3 (25.0 pg/mL)	Sample 3	Sample 3
D	Standard 4 (12.5 pg/mL)	Standard 4 (12.5 pg/mL)	Sample 4	Sample 4
E	Standard 5 (6.3 pg/mL)	Standard 5 (6.3 pg/mL)	Sample 5	Sample 5
F	Standard 6 (3.1 pg/mL)	Standard 6 (3.1 pg/mL)	Sample 6	Sample 6
G	Standard 7 (1.6 pg/mL)	Standard 7 (1.6 pg/mL)	Sample 7	Sample 7
H	Blank	Blank	Sample 8	Sample 8

In case of an **external standard dilution** (see "External standard dilution" on page 3), pipette 100 µL of these standard dilutions (S1 - S7) in the standard wells according to Table 1.

4. Add 100 µL of Sample Diluent in duplicate to the blank wells.
5. Add 80 µL of Sample Diluent to the sample wells.
6. Add 20 µL of each sample in duplicate to the sample wells.

7. Cover with an adhesive film and incubate at room temperature (18 to 25°C) for 2 hours, on a microplate shaker.
8. Prepare Biotin-Conjugate (see "Biotin-Conjugate" on page 3).
9. Remove adhesive film and empty wells. Wash microwell strips 4 times according to point 2 of the test protocol.
10. Add 100 µL of Biotin-Conjugate to all wells, including blank wells.
11. Cover with an adhesive film and incubate at room temperature (18 to 25°C) for 1 hour, if available on a microplate shaker.
12. Prepare Streptavidin-HRP (see "Streptavidin-HRP" on page 3).
13. Remove adhesive film and empty wells. Wash microwell strips 4 times according to point 2 of the test protocol.
14. Pipette 100 µL of diluted Streptavidin-HRP to all wells, including blanks.
15. Cover with an adhesive film and incubate at room temperature (18 to 25°C) for 30 minutes on a microplate shaker.
16. Remove adhesive film and empty wells. Wash microwell strips 4 times according to point 2 of the test protocol.
17. Pipette 100 µL of TMB Substrate Solution to all wells.
18. Incubate the microwell strips at room temperature (18° to 25°C) for about 30 min. Avoid direct exposure to intense light.

The color development on the plate should be monitored and the substrate reaction stopped (see next point 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.

19. 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 - 8°C in the dark.
20. 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.

Note: If instructions of this protocol have been followed samples have been diluted 1:5, the concentration read from the standard curve must be multiplied by the dilution factor (x5).

Calculation of results

- Calculate the average absorbance values for each set of duplicate standards and samples. Duplicates should be within 20 percent of the mean value.
- Create a standard curve by plotting the mean absorbance for each standard concentration on the ordinate against the human MICA 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 MICA 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 MICA concentration.
- If instructions in this protocol have been followed, samples have been diluted 1:5 (20 µL sample + 80 µL Sample Diluent), the concentration read from the standard curve must be multiplied by the dilution factor (x 5).

- Calculation of samples with a concentration exceeding standard 1 will result in incorrect, low human MICA levels (Hook Effect). Such samples require further external predilution according to expected human MICA values with Sample Diluent in order to precisely quantitate the actual human MICA level.
- It is suggested that each testing facility establishes a control sample of known human MICA 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 9.

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.

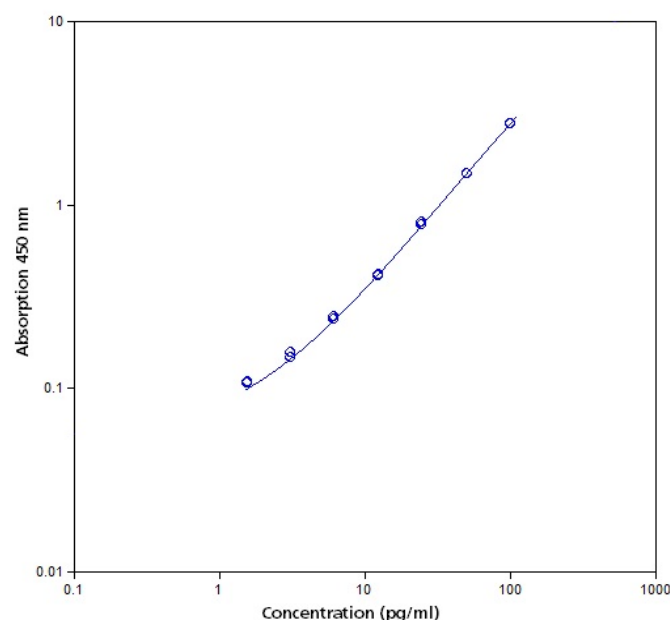


Fig. 9 Representative standard curve for human MICA ELISA. Human MICA was diluted in serial 2-fold steps in Sample Diluent.

Table 2 Typical data using the human MICA ELISA
Measuring wavelength: 450 nm, Reference wavelength: 620 nm

Standard	Human MICA Concentration (pg/mL)	O.D. at 450 nm	Mean O.D. at 450 nm	C.V. (%)
1	100	3.519 3.539	3.529	0.3
2	50	2.172 2.261	2.216	2.0
3	25.5	1.220 1.252	1.236	1.3
4	12.5	0.739 0.752	0.746	0.9
5	6.3	0.460 0.467	0.463	0.6
6	3.1	0.323 0.309	0.316	2.3
7	1.6	0.230 0.240	0.235	2.1
Blank	0	0.143 0.151	0.147	2.8

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

- Since 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 MICA 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 1.09 pg/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 serum and plasma samples containing different concentrations of human MICA. 2 standard curves were run on each plate. Data below show the mean human MICA concentration and the coefficient of variation for each sample (see Table 3). The calculated overall intra-assay coefficient of variation was 5.1%

Table 3 The mean human MICA concentration and the coefficient of variation for each sample

Sample	Experiment	Mean Human MICA Concentration (pg/mL)	Coefficient of Variation (%)
1	1	367.9	2.3
	2	374.1	2.7
	3	396.7	2.5
2	1	77.8	3.1
	2	76.2	2.8
	3	82.8	2.2
3	1	26.4	6.3
	2	25.8	5.3
	3	25.1	14.4
4	1	26.1	7.4
	2	26.0	4.4
	3	28.2	4.0
5	1	64.1	4.4
	2	62.4	2.8
	3	71.5	3.0
6	1	54.2	3.7
	2	50.5	2.8
	3	57.6	7.3
7	1	83.1	5.4
	2	56.4	3.4
	3	83.4	10.9
8	1	70.1	2.6
	2	67.2	6.6
	3	70.4	11.4

Inter-assay

Assay to assay reproducibility within one laboratory was evaluated in 3 independent experiments. Each assay was carried out with 6 replicates of serum and plasma samples containing different concentrations of human MICA. 2 standard curves were run on each plate. Data below show the mean human MICA concentration and the coefficient of variation calculated on 18 determinations of each sample (see Table 4). The calculated overall inter-assay coefficient of variation was 6.6%.

Table 4 The mean human MICA concentration and the coefficient of variation of each sample

Sample	Mean Human MICA Concentration (pg/mL)	Coefficient of Variation (%)
1	379.6	4.0
2	78.9	4.3
3	25.8	2.4
4	26.7	4.6
5	66.0	7.3
6	54.1	6.6
7	74.3	20.8
8	69.3	2.6

Spike recovery

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

Table 5

Sample matrix	Spike high		Spike medium	
	Mean (%)	Range (%)	Mean (%)	Range (%)
Serum	113	91 – 126	108	102 – 116
Plasma (EDTA)	88	64 – 113	87	64 – 100
Plasma (citrate)	103	80 – 127	82	68 – 96
Plasma (heparin)	100	91 – 114	101	99 – 102
Cell culture supernatant	101	100 – 102	107	102 – 112

Dilution parallelism

Serum, plasma (EDTA, citrate, heparin), cell culture supernatant samples with different levels of human MICA were analysed at serial 2 fold dilutions with 4 replicates each.

Sample matrix	Dilution	Expected Values	
		Mean (%)	Range (%)
Serum	10	97	94 - 100
	20	99	95 - 101
	40	105	95 - 113
Plasma (EDTA)	10	99	93 - 103
	20	101	93 - 111
	40	115	100 - 131
Plasma (citrate)	10	100	93 - 107
	20	105	95 - 116
	40	115	104 - 136
Plasma (heparin)	10	104	97 - 114
	20	110	102 - 116
	40	120	108 - 129
Cell culture supernatant	4	96	94 - 99
	8	95	90 - 100
	16	100	96 - 105

Sample stability

Freeze-Thaw stability

Aliquots of serum were stored at -20°C and thawed 3 times, and the human MICA levels determined. There was no significant loss of human MICA immunoreactivity detected by freezing and thawing.

Storage stability

Aliquots of serum were stored at -20°C, 2-8°C, room temperature (RT) and at 37°C, and the human MICA level determined after 24 h. There was no significant loss of human MICA immunoreactivity detected during storage under above conditions.

Specificity

The assay detects both natural and recombinant human MICA. There was no cross reactivity or interference detected.

Expected values

Panels of 40 serum as well as plasma samples (EDTA, citrate, heparin), from randomly selected healthy donors (males and females) and 9 Cell culture supernatants and lysates samples were tested for MICA.

Sample matrix	Number of samples evaluated	Mean (pg/mL)	Range (pg/mL)	Standard deviation (pg/mL)
Serum	40	74.2	0.0 - 205.1	51.8
Plasma (EDTA)	40	34.8	0.5 - 119.1	32.4
Plasma (citrate)	40	52.3	0.0 - 148.8	36.9
Plasma (heparin)	40	32.5	2.2 - 72.7	19.1

Note: The levels measured may vary with the sample collection used.

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 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	0.06	5.94
1 - 12	0.12	11.88

Streptavidin-HRP

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

Number of Strips	Streptavidin-HRP (mL)	Assay Buffer (1x) (mL)
1 - 6	0.06	5.94
1 - 12	0.12	11.88

Human MICA standard

Reconstitute human MICA standard with distilled water.
(Reconstitution volume is stated on the label of the standard vial.)

Test protocol summary

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 Sample Diluent, in duplicate, to all standard wells. Pipette 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): Pipette 100 µL of these standard dilutions in the microwell strips.
4. Add 100 µL Sample Diluent, in duplicate, to the blank wells.
5. Add 80 µL Sample Diluent to sample wells.
6. Add 20 µL sample in duplicate, to designated sample wells.
7. Cover microwell strips and incubate 2 hours at room temperature (18°-25°C) on a microplate shaker.
8. Prepare Biotin-Conjugate.
9. Empty and wash microwell strips 4 times with Wash Buffer.
10. Add 100 µL diluted Biotin-Conjugate to all wells.
11. Cover microwell strips and incubate 1 hour at room temperature (18° to 25°C).
12. Prepare Streptavidin-HRP.
13. Empty and wash microwell strips 4 times with Wash Buffer.
14. Add 100 µL diluted Streptavidin-HRP to all wells.
15. Cover microwell strips and incubate for about 30 minutes at room temperature (18° to 25°C) on a microplate shaker.
16. Empty and wash microwell strips 4 times with Wash Buffer.
17. Add 100 µL of TMB Substrate Solution to all wells.
18. Incubate the microwell strips for about 30 minutes at room temperature (18°C to 25°C).

19. Add 100 µL Stop Solution to all wells.

20. Blank microwell reader and measure colour intensity at 450 nm.

Note: If instructions in this protocol have been followed, samples have been diluted 1:5 (20 µL sample + 80 µL Sample Diluent) and the concentration read from the standard curve must be multiplied by the dilution factor (x 5).

Customer and technical support

Visit thermofisher.com/support for the latest in services and support, including:

- Worldwide contact telephone numbers
- Product support, including:
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- 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. If you have any questions, please contact Life Technologies at www.thermofisher.com/support.



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

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.

©2018 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.