

# Human GDF15 ELISA Kit

## Enzyme-linked Immunosorbent Assay for quantitative detection of human GDF15

Catalog Numbers BMS2258 and BMS2258TEN

Pub. No. MAN0017159 Rev. D (33)



**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](https://www.thermofisher.com/support).

### Product description

The Human GDF15 ELISA Kit is a research use enzyme-linked immunosorbent assay for the quantitative detection of human GDF15.

### Summary

GDF15 (Growth/differentiation factor 15) is a protein belonging to the transforming growth factor beta superfamily that has a role in regulating inflammatory and apoptotic pathways in injured cells. This gene encodes a secreted ligand of the TGF-beta (transforming growth factor-beta) superfamily of proteins. Ligands of this family bind various TGF-beta receptors leading to recruitment and activation of SMAD family transcription factors regulating gene expression. The encoded preproprotein is proteolytically processed to generate each subunit of the disulfide-linked homodimer.

GDF15 is also known as TGF-PL, MIC-1, PDF, PLAB, and PTGFB. The protein is expressed in a broad range of cell types, acts as a pleiotropic cytokine, and is involved in the stress response program of cells after cellular injury.

For literature update visit our website.

### Principles of the test

An anti-human GDF15 coating antibody is adsorbed onto microwells.

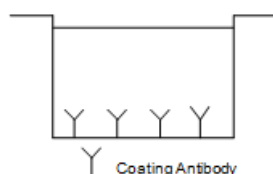


Figure 1 Coated microwell.

Human GDF15 present in the sample or standard binds to antibodies adsorbed to the microwells. A biotin-conjugated anti-human GDF15 antibody is added and binds to human GDF15 captured by the first antibody.

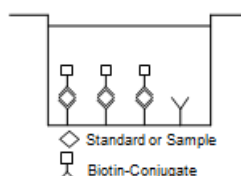


Figure 2 First incubation.

Following incubation, unbound biotin-conjugated anti-human GDF15 antibody is removed during a wash step. Streptavidin-HRP is added and binds to the biotin-conjugated anti-human GDF15 antibody.

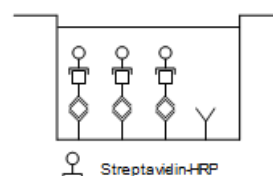


Figure 3 Second incubation.

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

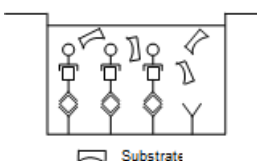
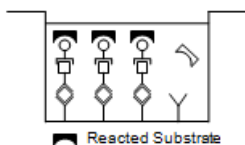


Figure 4 Third incubation.

A colored product is formed in proportion to the amount of human GDF15 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 GDF15 standard dilutions, and human GDF15 sample concentration is determined.



**Figure 5** Fourth incubation.

## Reagents provided

### Reagents for human GDF15 ELISA BMS2258 (96 tests)

1 aluminum pouch with a Microwell Plate coated with monoclonal antibody to human GDF15

1 vial (70  $\mu$ L) Biotin-Conjugate anti-human GDF15 polyclonal antibody

1 vial (150  $\mu$ L) Streptavidin-HRP

2 vials human GDF15 Standard lyophilized, 250 pg/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 (50 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 GDF15 ELISA BMS2258TEN (10x96 tests)

10 aluminum pouches with a Microwell Plate coated with monoclonal antibody to human GDF15

10 vials (70  $\mu$ L) Biotin-Conjugate anti-human GDF15 polyclonal antibody

10 vials (150  $\mu$ L) Streptavidin-HRP

10 vials human GDF15 Standard lyophilized, 250 pg/ml upon reconstitution

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

5 bottles (50 ml) Wash Buffer Concentrate 20x

5 bottles (50 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° and 8°C . Immediately after use remaining reagents should be returned to cold storage (2–8°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, 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 GDF15. Avoid repeated freeze-thaw cycles. Prior to assay, the frozen sample should be brought to room temperature slowly and mixed gently.

## Required materials not supplied

- 5 mL and 10 mL graduated pipettes
- 5  $\mu$ L to 1,000  $\mu$ L adjustable single channel micropipettes with disposable tips
- 50  $\mu$ L to 300  $\mu$ 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)
- Microplate shaker
- Microwell strip reader capable of reading at 450 nm (620 nm as optional reference wavelength)
- 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 of 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.

## Before you begin

- Equilibrate the buffer concentrates to room temperature (18–25°C), then dilute before use.
- If crystals have formed in the buffer concentrates, warm gently to dissolve the crystals.

## Prepare 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 2–25°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

## Prepare 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

## Prepare 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

## Prepare Streptavidin-HRP

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

Make a 1:200 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	0.03	5.97
1–12	0.06	11.94

## Prepare Human GDF15 Standard

1. Reconstitute Human GDF15 Standard by adding distilled water. Reconstitution volume is stated on the label of the standard vial. Swirl or mix gently to ensure complete and homogeneous solubilization (concentration of reconstituted standard = 250 pg/mL).
2. Allow the standard to reconstitute for 10–30 minutes. Mix well prior to making dilutions.
3. Remaining standard cannot be stored after usage and must be discarded.
4. Standard dilutions can be prepared directly on the microwell plate (see step 3) or alternatively in tubes (see “External standard dilution” on page 4).

## External standard dilution

1. Label 7 tubes, one for each standard point: S1, S2, S3, S4, S5, S6, and S7.
2. Prepare 1:2 serial dilutions for the standard curve as follows:  
Pipet 225 µL of Sample Diluent into each tube.
3. Pipet 225 µL of reconstituted standard (concentration of standard = 250 pg/mL) into the first tube, labeled S1, and mix (concentration of standard 1 = 125 pg/mL).
4. Pipet 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.

Sample Diluent serves as blank.

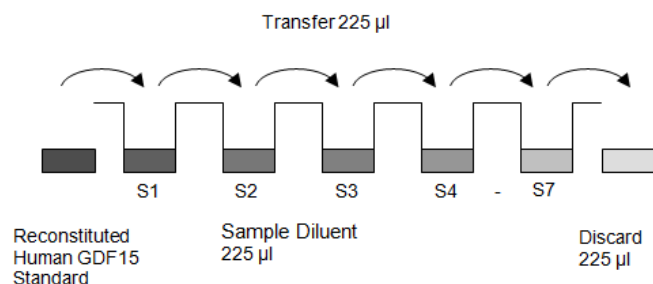


Figure 6 Dilute standards - tubes.

## Test protocol

**Note:** If instructions of this protocol have been followed, samples have been diluted 1:40 (1:20 external predilution: 10 µL Sample + 190 µL Sample Diluent, on the plate 1:2), and the concentration read from the standard curve must be multiplied by the dilution factor (x 40).

Shaking is absolutely necessary for an optimal test performance.

1. Predilute your samples before starting with the test procedure. Dilute serum, plasma, and cell culture samples 1:20 with Sample Diluent according to the following scheme:  
10 µL Sample + 190 µL Sample Diluent
2. 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 2–8°C sealed tightly.
3. Wash the microwell strips twice with approximately 400 µL Wash Buffer per well. Thoroughly aspirate 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.

4. Prepare standard dilution on the microwell plate (alternatively, the standard dilution can be prepared in tubes, see “External standard dilution” on page 4):

Add 100 µL of Sample Diluent in duplicate to all standard wells. Pipet 100 µL of prepared standard (see “Prepare Human GDF15 Standard” on page 3, concentration = 250 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 = 125 pg/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 GDF15 standard dilutions ranging from 125 to 2.0 pg/mL. Discard 100 µL of the contents from the last microwells (S7) used.

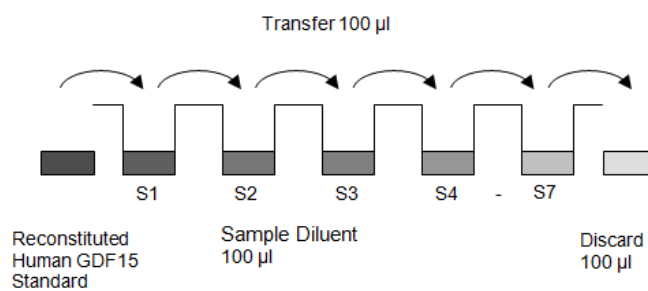


Figure 7 Dilute standards, microwell plate.

In case of an external standard dilution (see “External standard dilution” on page 4), 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 1 125 pg/mL	Standard 1 125 pg/mL	Sample 1	Sample 1
B	Standard 2 62.5 pg/mL	Standard 2 62.5 pg/mL	Sample 2	Sample 2
C	Standard 3 31.25 pg/mL	Standard 3 31.25 pg/mL	Sample 3	Sample 3
D	Standard 4 15.63 pg/mL	Standard 4 15.63 pg/mL	Sample 4	Sample 4
E	Standard 5 7.81 pg/mL	Standard 5 0.25 ng/mL	Sample 5	Sample 5
F	Standard 6 3.91 pg/mL	Standard 6 3.91 pg/mL	Sample 6	Sample 6
G	Standard 7 2.0 pg/mL	Standard 7 2.0 pg/mL	Sample 7	Sample 7
H	Blank	Blank	Sample 8	Sample 8

5. Add 100 µL of Sample Diluent in duplicate to the blank wells.
6. Add 50 µL of Sample Diluent to the sample wells.
7. Add 50 µL of each prediluted sample in duplicate to the sample wells.

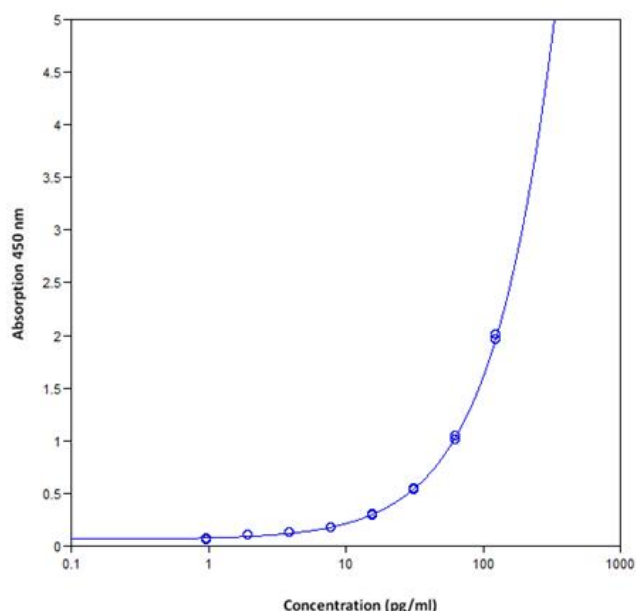
8. Prepare Biotin-Conjugate (see “Prepare Biotin-Conjugate” on page 3).
9. Add 50 µL of Biotin-Conjugate to all wells.
10. Cover with an adhesive film and incubate at room temperature for 2 hours on a microplate shaker set at 400 rpm.  
Shaking is absolutely necessary for an optimal test performance.
11. Prepare Streptavidin-HRP (see “Prepare Streptavidin-HRP” on page 3).
12. Remove adhesive film and empty wells. Wash microwell strips 4 times according to step 2. Proceed immediately to the next step.
13. Add 100 µL of diluted Streptavidin-HRP to all wells, including the blank wells.
14. Cover with an adhesive film and incubate at room temperature for 1 hour on a microplate shaker set at 400 rpm.  
Shaking is absolutely necessary for an optimal test performance.
15. Remove adhesive film and empty wells. Wash microwell strips 4 times according to step 2. Proceed immediately to the next step.
16. Pipet 100 µL of TMB Substrate Solution to all wells.
17. Incubate the microwell strips at room temperature 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) 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.
18. 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.
19. Read absorbance of each microwell on a spectrophotometer 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 GDF15 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 GDF15 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 GDF15 concentration.
- If instructions of this protocol have been followed, samples have been diluted 1:40 (1:20 external predilution: 10 µl Sample + 190 µl Sample Diluent, on the plate 1:2). The concentration read from the standard curve must be multiplied by the dilution factor (x40).
- Calculation of samples with a concentration exceeding standard 1 will result in incorrect, low human GDF15 levels. Such samples require further external predilution according to expected human GDF15 values with Sample Diluent in order to precisely quantify the actual human GDF15 level.
- It is suggested that each testing facility establishes a control sample of known human GDF15 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.



**Figure 8** Representative standard curve for Human GDF15 ELISA Kit. Human GDF15 was diluted in serial 2-fold steps in Sample Diluent.

**Table 2** Typical data using the Human GDF15 ELISA Kit (measuring wavelength of 450 nm, reference wavelength of 620 nm).

Standard	Human GDF15 concentration (pg/mL)	O.D. at 450 nm	Mean O.D. at 450 nm	C.V. (%)
1	125.00	2.006 1.966	1.986	1.0
2	62.50	1.041 1.011	1.026	1.5
3	31.25	0.550 0.531	0.541	1.8
4	15.63	0.293 0.300	0.297	1.2
5	7.81	0.178 0.171	0.175	2.0
6	3.91	0.126 0.127	0.127	0.4
7	2.0	0.110 0.100	0.105	4.8
Blank	0	0.064 0.057	0.061	5.8

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

- 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 may 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 GDF15 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 0.589 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 8 serum samples containing different concentrations of human GDF15. Two standard curves were run on each plate. Data below show the mean human GDF15 concentration and the coefficient of variation for each sample. The calculated overall intra-assay coefficient of variation was 4.2%.

**Table 3 The mean human GDF15 concentration and the coefficient of variation for each sample.**

Sample	Experiment	Mean human GDF15 concentration (pg/mL)	Coefficient of variation (%)
1	1	1,209.5	4.0
	2	1,222.0	5.1
	3	1,203.9	4.4
2	1	818.0	4.2
	2	800.5	7.0
	3	828.6	3.0
3	1	513.5	3.1
	2	524.8	4.3
	3	521.9	2.6
4	1	345.2	8.0
	2	404.5	8.7
	3	367.3	6.9
5	1	546.8	2.0
	2	555.9	4.6
	3	547.0	3.6
6	1	823.4	3.8
	2	834.0	3.5
	3	839.8	3.2
7	1	474.3	3.2
	2	504.0	5.8
	3	523.1	3.6
8	1	743.1	1.7
	2	745.2	3.5
	3	801.1	2.1

#### 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 samples containing different concentrations of human GDF15. Two standard curves were run on each plate. Data below show the mean human GDF15 concentration and the coefficient of variation calculated on 18 determinations of each sample. The calculated overall inter-assay coefficient of variation was 2.9%.

**Table 4 The mean human GDF15 concentration and the coefficient of variation of each sample.**

Sample	Mean human GDF15 concentration (pg/mL)	Coefficient of variation (%)
1	1211.8	0.8
2	815.7	1.7
3	520.1	1.1
4	372.3	8.1
5	549.9	0.9
6	832.4	1.0
7	500.4	4.9
8	763.2	4.3

### Spike recovery

The spike recovery was evaluated by spiking 3 levels of human GDF15 into serum, plasma (EDTA, heparin, citrate), and cell culture supernatant. Recoveries were determined with 2 replicates each. The amount of endogenous human GDF15 in unspiked samples was subtracted from the spike values.

Sample matrix	Spike high (%)		Spike medium (%)		Spike low (%)	
	Mean	Range	Mean	Range	Mean	Range
Serum	103	93–129	97	79–119	110	98–28
Plasma (EDTA)	83	77–94	102	89–117	98	93–106
Plasma (citrate)	98	84–120	94	73–112	103	93–126
Plasma (heparin)	84	74–100	98	89–112	96	88–108
Cell culture supernatant	96	93–99	99	98–101	108	96–121

## Dilution parallelism

Serum, plasma (EDTA, citrate, heparin), and cell culture supernatant samples with different levels of human GDF15 were analysed at serial 2-fold dilutions with 3 replicates each.

Sample matrix	Recovery of exp. val.	
	Dilution	Mean %
Serum	80	108
	160	113
	320	110
Plasma (EDTA)	80	112
	160	120
	320	117
Plasma (citrate)	80	118
	160	116
	320	98
Plasma (heparin)	80	96
	160	101
	320	74
Cell culture supernatant	80	87
	160	86
	320	88

## Sample stability

### Freeze-thaw stability

Aliquots of serum, plasma, and cell culture supernatant samples (spiked or unspiked) were stored at  $-20^{\circ}\text{C}$  and thawed 3 times, and the human GDF15 levels determined. There was no significant loss of human GDF15 immunoreactivity detected by freezing and thawing.

### Storage stability

Aliquots of serum, plasma, and cell culture supernatant samples (spiked or unspiked) were stored at  $-20^{\circ}\text{C}$ ,  $2-8^{\circ}\text{C}$ , room temperature, and at  $37^{\circ}\text{C}$ , and the human GDF15 level determined after 24 hours. No significant loss of human GDF15 immunoreactivity was detected under above conditions .

## Specificity

The cross-reactivity and interference of circulating factors of the immune system were evaluated by spiking these proteins at physiologically relevant concentrations into a human GDF15 positive sample. No cross-reactivity or interference was detected.

## Expected values

Panels of 40 serum samples, as well as plasma samples (EDTA, citrate, heparin), from randomly selected healthy donors (males and females) were tested for human GDF15.

Table 5

Sample matrix	Number of samples evaluated	Mean (pg/mL)	Range (pg/mL)	Standard deviation (pg/mL)	Positive samples (%)
Serum	40	680	300–1,401	278	40
Plasma (EDTA)	40	706	355–1,951	335	40
Plasma (citrate)	40	542	197–1,194	251	40
Plasma (heparin)	40	481	275–1,011	278	40

## 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 1:200 dilution of Streptavidin-HRP in Assay Buffer (1x) .

Number of Strips	Streptavidin-HRP (mL)	Assay Buffer (1x) (mL)
1–6	0.03	5.97
1–12	0.06	11.94



## Human GDF15 standard

Reconstitute lyophilized human GDF15 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 1:40 (1:20 external predilution: 10 µL Sample + 190 µL Sample Diluent, on the plate 1:2) and the concentration read from the standard curve must be multiplied by the dilution factor (x 40).

Shaking is absolutely necessary for an optimal test performance.

1. Predilute your samples before starting with the test procedure. Dilute serum, plasma, and cell culture samples 1:20 with 10 µL Sample + 190 µL Sample Diluent
2. Determine the number of microwell strips required.
3. Wash microwell strips twice with Wash Buffer.
4. Standard dilution on the microwell plate: Add 100 µL Sample Diluent, 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 4): Pipet 100 µL of these standard dilutions in the microwell strips.
5. Add 100 µL Sample Diluent, in duplicate, to the blank wells.
6. Add 50 µL Sample Diluent to sample wells.
7. Add 50 µL of each pre-diluted sample in duplicate, to designated sample wells.

8. Prepare Biotin-Conjugate.
9. Add 50 µL Biotin-Conjugate to all wells.
10. Cover microwell strips and incubate 2 hours at room temperature (18–25°C) on a microplate shaker.
11. Prepare Streptavidin-HRP.
12. Empty and wash microwell strips 4 times with Wash Buffer.
13. Add 100 µL diluted Streptavidin-HRP to all wells.
14. Cover microwell strips and incubate 1 hour at room temperature (18–25°C).
15. Empty and wash microwell strips 4 times with Wash Buffer.
16. Add 100 µL of TMB Substrate Solution to all wells.
17. Incubate the microwell strips for about 30 minutes at room temperature (18–25°C).
18. Add 100 µL Stop Solution to all wells.
19. Blank microwell reader and measure color intensity at 450 nm.

## Limited product warranty

Life Technologies Corporation and its affiliates 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](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have questions, contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).



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

**Revision history:** Pub. No. MAN0017159 D (33)

Revision	Date	Description
D (33)	28 January 2025	Minor updates throughout the document. Version format was changed in conformance with internal document control procedures.
C.0 (32)	30 November 2022	Heading of table column changed
B.0 (31)	1 January 2018	Revision not specified

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.

©2018-2025 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.