



EAGLE
BIOSCIENCES

GDF-15 ELISA Assay Kit

Catalog Number: GDF31-K01

For the quantitative determination of human GDF-15 concentrations
in serum, plasma and cell culture supernate samples.

This package insert must be read in its entirety before using this product.

1 x 96 Wells

For Research Use Only (RUO). Not for use in clinical, diagnostic or therapeutic procedures.

EAGLE BIOSCIENCES, INC.

20A Northwest Blvd., Suite 112, Nashua, NH 03063

Phone: 617-419-2019 Fax: 617-419-1110

www.EagleBio.com



TABLE OF CONTENTS

Contents	Page
ASSAY BACKGROUND	1
ASSAY PRINCIPLE	1
INTENDED USE	1
REAGENTS SUPPLIED	2
OTHER MATERIALS REQUIRED, BUT NOT PROVIDED	2
STORAGE	2
PREPARATION OF REAGENTS	3
PREPARATION OF STANDARDS AND SAMPLES	4
ASSAY PROCEDURE	5
CALCULATION	5
TYPICAL STANDARD CURVE	6
ASSAY CHARACTERISTICS	6
REFERENCES	6
SUMMARY OF ASSAY PROCEDURE	7



ASSAY BACKGROUND

Growth differentiation factor 15 (GDF-15) is a member of the transforming growth factor β cytokine superfamily, the protein is secreted as a 25 kDa disulfide linked dimer [1]. Circulating GDF-15 levels are associated with cancers, cardiovascular and kidney diseases [2-4]. GDF-15 regulates food intake, energy expenditure and body weight in response to metabolic and toxin-induced stresses [5].

ASSAY PRINCIPLE

This assay is a quantitative sandwich ELISA. The immunoplate is pre-coated with a monoclonal antibody specific for human GDF-15. Standards and samples are pipetted into the wells and any human GDF-15 present is bound by the immobilized antibody. After washing away any unbound substances, a biotin labelled polyclonal antibody specific for human GDF-15 is added to the wells. After wash step to remove any unbound reagents, streptavidin-HRP conjugate (STP-HRP) is added. After the last wash step, an HRP substrate solution is added and colour develops in proportion to the amount of human GDF-15 bound initially. The assay is stopped and the optical density of the wells determined using a microplate reader. Since the increases in absorbance are directly proportional to the amount of captured human GDF-15, the unknown sample concentration can be interpolated from a reference curve included in each assay.

INTENDED USE

The Eagle Biosciences GDF-15 ELISA Assay Kit is designed for quantification of human GDF-15 in serum, plasma and cell culture supernate samples. The GDF-15 ELISA Assay Kit is for research use only and not to be used for diagnostic procedures.



REAGENTS SUPPLIED

Each kit is sufficient for one 96-well plate and contains the following components:

1. Micro-titre Strips (96 wells)-Coated with a monoclonal antibody against human GDF-15, sealed.
2. 10×Wash buffer-50 ml.
3. 5×Assay buffer-20 ml.
4. 100×Detection antibody solution-A biotin labelled polyclonal antibody against human GDF-15, 0.12 ml.
5. Human GDF-15 standard-500 pg of recombinant human GDF-15 in a buffered protein base, lyophilised.
6. 200×STP-HRP solution- 0.06 ml.
7. Substrate solution- 12 ml, ready for use.
8. Stop solution- 12 ml, ready for use.

OTHER MATERIALS REQUIRED, BUT NOT PROVIDED

1. Pipettes and pipette tips.
2. 96-well plate or manual strip washer.
3. Buffer and reagent reservoirs.
4. Paper towels or absorbent paper.
5. Plate reader capable of reading absorbency at 450 nm.
6. Distilled water or deionized water.

STORAGE

The kit should be stored at 2-8°C upon receipt, and all reagents should be equilibrated to room temperature before use. Remove any unused antibody-coated strips from the human GDF-15 microplate, return them to the foil pouch and re-seal. Once opened, the strips may be stored at 2-8°C for up to one month.



PREPARATION OF REAGENTS

Bring all reagents and materials to room temperature before assay.

A. 1×Assay buffer.

Prepare 1×Assay buffer by mixing the 5×Assay buffer (20 ml) with 80 ml of distilled water or deionized water. If precipitates are observed in the 5×Assay buffer bottle, warm the bottle in a 37°C water bath until the precipitates disappear. The 1×Assay buffer may be stored at 2-8°C for up to one month.

B. 1×Wash buffer.

Prepare 1×Wash buffer by mixing the 10×Wash buffer (50 ml) with 450 ml of distilled water or deionized water. If precipitates are observed in the 10×Wash buffer bottle, warm the bottle in a 37°C water bath until the precipitates disappear. The 1×Wash buffer may be stored at 2-8°C for up to one month.

C. 1×Detection antibody solution.

Spin down the 100×Detection antibody solution briefly and dilute the desired amount of the antibody 1:100 with 1×Assay buffer, 100 l of the 1×Detection antibody solution is required per well. Prepare only as much 1×Detection antibody solution as needed. Return the 100×Detection antibody solution to 2-8°C immediately after the necessary volume is removed.

D. 1×STP-HRP solution.

Spin down the 200×STP-HRP solution briefly and dilute the desired amount of the 200×STP-HRP solution 1:200 with 1×Assay buffer, 100 l of the 1×STP-HRP solution is required per well. Prepare only as much 1×STP-HRP solution as needed. Return the 200×STP-HRP solution to 2-8°C immediately after the necessary volume is removed.



PREPARATION OF STANDARDS AND SAMPLES

Human GDF-15 standards: Reconstitute the lyophilised standard with 1 ml of 1×Assay buffer to generate a standard stock solution of 500 pg/ml. Allow the standard to sit for 10 minutes with gentle agitation prior to making dilutions. Prepare serially diluted standards using 1×Assay buffer as follows:

Standard volume	Volume of 1×Assay buffer	Concentration
500 pg/ml stock	-	500 pg/ml
250 l of 500 pg/ml	250 l	250 pg/ml
250 l of 250 pg/ml	250 l	125 pg/ml
250 l of 125 pg/ml	250 l	62.5 pg/ml
250 l of 62.5 pg/ml	250 l	31.2 pg/ml
250 l of 31.2 pg/ml	250 l	15.6 pg/ml
250 l of 15.6 pg/ml	250 l	7.8 pg/ml

1×Assay buffer serves as the zero standard (0 pg/ml).

Note: The reconstituted standard stock should be aliquoted and stored at -80°C for up to one month. Avoid repeating freezing/thawing cycles. Please do not store the diluted standard solutions.

Sample preparation

Serum or plasma sample is generally required a 10-fold dilution in the 1×Assay buffer.



Assay Procedure

It is recommended that all standards and samples should be assayed in duplicate.

1. Add 100 μl of standard or sample per well, incubate at room temperature for 1 hour.
2. Discard the content and tap the plate on a clean paper towel to remove residual solution in each well. Add 300 μl of 1 \times Wash buffer to each well and incubate for 1 minute. Discard the 1 \times Wash buffer and tap the plate on a clean paper towel to remove residual wash buffer. Repeat the wash step for a total 3 washes.
3. Add 100 μl of 1 \times Detection antibody solution to each well, incubate at room temperature for 1 hour.
4. Wash each well 3 times as in step 2.
5. Add 100 μl of 1 \times STP-HRP solution to each well, incubate at room temperature for 20 minutes.
6. Wash each well 4 times as described in step 2.
7. Add 100 μl of Substrate solution to each well, incubate at room temperature for 15 minutes. Protect from light.
8. Add 100 μl of Stop solution to each well, gently tap the plate frame for a few seconds to ensure thorough mixing.
9. Measure absorbance of each well at 450 nm immediately.

Calculation

1. Subtract the absorbance of the blank from that of standards and samples.
2. Generate a standard curve by plotting the absorbance obtained (y-axis) against human GDF-15 concentrations (x-axis). The best fit line can be generated with any curve-fitting software by regression analysis. Any curve of 4-parameter or log-log curve fitting can be used for calculation.
3. Determine human GDF-15 concentration of samples from standard curve and multiply the value by the dilution factor.



Typical standard curve

The following standard curve is provided for demonstration only. A standard curve should be generated for each set of sample assay.

Human GDF-15 (pg/ml)	Absorbance (450 nm)	Blanked Absorbance
0	0.074	0
7.8	0.13	0.056
15.6	0.197	0.123
31.2	0.338	0.264
62.5	0.577	0.503
125	1.017	0.943
250	1.719	1.645
500	2.626	2.552

ASSAY CHARACTERISTICS

A. Sensitivity:

The lowest level of human GDF-15 that can be detected by this assay is 7.8 pg/ml.

B. Precision:

Intra-assay Precision (Precision within an assay) <10%.

Inter-assay Precision (Precision between assays) <10%.

References:

- [1] Bootcov M. R., et al. (1997) *PNAS*; 94:11514-11519.
- [2] Welsh J.B., et al. (2003) *PNAS*;100:3410-3415.
- [3] Wang T. J, et al. (2012) *Circulation* 126; 1596– 1604.
- [4] Nair V., et al. (2017) *J.Am. Soc. Nephrol.* 28 2223-2240.
- [5] Hsu J-Y, et al. (2017) *Nature*; 550: 255-259.



Summary of Assay Procedure

Add 100 μ l of Standard or sample to each well.

Incubate at room temperature for 1 hour.

Aspirate and wash each well three times.

Add 100 μ l of 1 \times Detection antibody solution to each well.

Incubate at room temperature for 1 hour.

Aspirate and wash each well three times.

Add 100 μ l of 1 \times STP-HRP solution to each well.

Incubate at room temperature for 20 minutes.

Aspirate and wash each well four times.

Add 100 μ l of Substrate solution to each well.

Incubate at room temperature for 15 minutes.

Add 100 μ l of Stop solution to each well.

Measure absorbance of each well at 450 nm.

Calculation

Warranty Information

Eagle Biosciences, Inc. warrants its Product(s) to operate or perform substantially in conformance with its specifications, as set forth in the accompanying package insert. This warranty is expressly limited to the refund of the price of any defective Product or the replacement of any defective Product with new Product. This warranty applies only when the Buyer gives written notice to the Eagle Biosciences within the expiration period of the Product(s) by the Buyer. In addition, Eagle Biosciences has no obligation to replace Product(s) as result of a) Buyer negligence, fault, or misuse, b) improper use, c) improper storage and handling, d) intentional damage, or e) event of force majeure, acts of God, or accident.

Eagle Biosciences makes no warranties, either expressed or implied, except as provided herein, including without limitation thereof, warranties as to marketability, merchantability, fitness for a particular purpose or use, or non-infringement of any intellectual property rights. In no event shall the company be liable for any indirect, incidental, or consequential damages of any nature, or losses or expenses resulting from any defective product or the use of any product. Product(s) may not be resold, modified, or altered for resale without prior written approval from Eagle Biosciences, Inc.

For further information about this kit, its application or the procedures in this kit insert, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.

distributed in the US/Canada by:

EAGLE BIOSCIENCES, INC.

20A NW Blvd, Suite 112 Nashua, NH 03063

Phone: 617-419-2019 FAX: 617-419-1110

www.EagleBio.com • info@eaglebio.com



EAGLE
BIOSCIENCES