



EAGLE
BIOSCIENCES

Mouse GDF-15 ELISA Assay Kit

Catalogue Number: MHF11-K01

For the quantitative determination of mouse 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 polyclonal antibody specific for mouse GDF-15. Standards and samples are pipetted into the wells and any mouse GDF-15 present is bound by the immobilized antibody. After washing away any unbound substances, a biotin labelled polyclonal antibody specific for mouse 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 mouse 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 mouse GDF-15, the unknown sample concentration can be interpolated from a reference curve included in each assay.

INTENDED USE

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



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 polyclonal antibody against mouse 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 mouse GDF-15, 0.12 ml.
5. Mouse GDF-15 standard-250 pg of recombinant mouse 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 mouse 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

Mouse GDF-15 standards: Reconstitute the lyophilised standard with 1 ml of 1×Assay buffer to generate a standard stock solution of 250 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
250 pg/ml stock	-	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
250 l of 7.8 pg/ml	250 l	3.9 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 μ l of standard or sample per well, incubate at room temperature for 2 hours.
2. Discard the content and tap the plate on a clean paper towel to remove residual solution in each well. Add 300 μ l 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 μ l 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 μ l 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 μ l of Substrate solution to each well, incubate at room temperature for 15 minutes. Protect from light.
8. Add 100 μ l 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 mouse 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 mouse 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.

Mouse GDF-15 (pg/ml)	Absorbance (450 nm)	Blanked Absorbance
0	0.091	0
3.9	0.126	0.035
7.8	0.164	0.073
15.6	0.246	0.155
31.2	0.404	0.313
62.5	0.692	0.601
125	1.259	1.168
250	2.216	2.125

ASSAY CHARACTERISTICS

A. Sensitivity:

The lowest level of mouse GDF-15 that can be detected by this assay is 3.9 pg/ml.

B. Precision:

Intra-assay Precision (Precision within an assay) < 2.9%.

Inter-assay Precision (Precision between assays) < 6.1%.

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 2 hours.

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