



EAGLE
BIOSCIENCES

Human Intact Fibroblast Growth Factor (FGF-21) ELISA

Catalog Number:

F2131-K01 (1 x 96 wells)

For Research Use Only. Not for use in diagnostic procedures.

v. 1.0

EAGLE BIOSCIENCES, INC.

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

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

WWW.EAGLEBIO.COM



INTENDED USE

The Eagle Biosciences Human Intact Fibroblast Growth Factor 21 (FGF-21) ELISA Assay Kit (enzyme-linked immunoassay kit) is intended for the quantitative determination of human Intact Fibroblast Growth Factor 21 (FGF-21) levels in serum. The Eagle Biosciences Human Intact Fibroblast Growth Factor 21 (FGF-21) ELISA Assay Kit is for research use only and not to be used in diagnostic procedures.

INTRODUCTION

Fibroblast Growth Factor 21 (FGF-21) belongs to the FGF-19 subfamily, which includes FGF-19, FGF-21 and FGF-23. The FGF-19 family members are potent endocrine hormones in the regulation of a diverse physiological homeostasis.

The intact FGF-21 is a small protein comprising 181 amino acids. Administration of recombinant FGF21 lowered plasma glucose and insulin levels, reduced hepatic and circulating triglycerides and cholesterol levels, and improved insulin sensitivity, energy expenditure, hepatic steatosis and obesity in a range of insulin resistant animal models. The physiological functions of FGF-21 are relied on the intact molecular structure and amino acid sequence in its N-terminal and C-terminal region. An N-terminal truncated FGF-21 (7-181) is a potent inhibitor that competitively inhibits the biological activity of intact FGF-21 (1-181). Therefore, it is important to measure the circulation intact FGF-21 level in the assessment of the physiological and pathophysiological condition. An assay that determines the fragment of the FGF-21 might overestimate the biological activity of the protein in test sample.

Circulation FGF-21 is a biomarker and its levels is increased in patient with nonalcoholic fatty liver disease (NAFLD), type 2 diabetes, gestational diabetes and obesity. An increase of circulating FGF-21 is also found in patient with Cushing's syndrome, patient with lipodystrophy induced by HIV-1 and patient with chronic renal disease or end-stage renal disease (ESRD).

PRINCIPLE OF THE ASSAY

The Eagle Biosciences Human Intact Fibroblast Growth Factor 21 (FGF-21) ELISA Assay Kit is designed, developed and produced for the quantitative measurement of human intact FGF-21 in serum and EDTA-plasma sample. The assay utilizes the two-site "sandwich" technique with two selected antibodies that bind to different epitopes of human intact FGF-21. One of the antibodies is specifically binds to the N-terminal human FGF-21 (1-7) and the other is specifically to the C-terminal human FGF-21 (175-181).

Assay standards, controls and patient samples are added directly to wells of microplate that is coated with an anti-human FGF-21 (1-7) specific antibody. Simultaneously, a horseradish peroxidase conjugated anti-human FGF-21 (175-181) specific antibody is added to each well. After the first incubation period, the antibody on the wall of microtiter well captures human FGF-21 in the sample and an unbound protein in each microtiter well is washed away. A "sandwich" of "anti-FGF-21 antibody --- human intact FGF-21 --- HRP conjugated tracer antibody" is formed. The unbound tracer antibody is removed in the subsequent washing step. For the detection of this immunocomplex, the well is then incubated with a substrate solution in a timed reaction and then measured in a spectrophotometric microplate reader. The enzymatic activity of the immunocomplex bound to human intact FGF-21 on the wall of the microtiter well is directly proportional to the amount of intact FGF-21 in the sample. A standard curve is generated by plotting the absorbance versus the respective human intact FGF-21 concentration



for each standard on point-to-point or 4 parameter curve fit. The concentration of human intact FGF-21 in test samples is determined directly from this standard curve.

REAGENTS: Preparation and Storage

This Intact Fibroblast Growth Factor 21 (FGF-21) ELISA Assay Kit must be stored at 2 – 8°C upon receipt. For the expiration date of the kit refer to the label on the kit box. All components are stable until the expiration date.

Allow all reagents to come to room temperature prior to use. Regents from different kit lot numbers should not be combined or interchanged.

1. Anti-Human FGF-21 Antibody Coated Microplate

One microplate with 12 x 8 well-breakable strips (96 wells total) coated with antibody to human FGF-21. The plate is framed and sealed in a foil zipper bag with a desiccant. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

2. Human FGF-21 Tracer Antibody

One vial containing 0.4 mL concentrated HRP labeled anti-human FGF-21 antibody in a stabilized protein matrix. This reagent must be diluted with FGF-21 Tracer Antibody Diluent before use. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

3. FGF-21 Tracer Antibody Diluent

One vial containing 8 mL ready to use buffer. It should be only used for tracer antibody dilution according to the assay procedures. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

4. ELISA Wash Concentrate

One bottle contains 30 mL of 30 fold concentrate. Before use the contents must be diluted with 870 mL of demineralized water and mix well. Upon dilution this yields a working wash solution containing a surfactant in phosphate buffered saline with a non-azide, non-mercury preservative. The diluted should be stored at room temperature and is stable until the expiration date on the kit box.

5. ELISA HRP Substrate

One bottle contains 12 mL of tetramethylbenzidine (TMB) with stabilized hydrogen peroxide. This reagent should be stored at 2 – 8°C and is stable until the expiration date on the kit box.

6. ELISA Stop Solution

One bottle contains 12 mL of 0.5 M sulfuric acid. This reagent should be stored at 2 – 8°C or room temperature and is stable until the expiration date on the kit box.

7. Human FGF-21 Standards

Six vials each contain different concentration of human FGF-21 in a lyophilized bovine serum based matrix with a non-azide, non-mercury preservative. **Refer to vial for exact concentration for each standard.** The standards are ready to use. These reagents should be stored at 2 – 8°C and are stable until the expiration date on the kit box.



8. Human FGF-21 Controls

Two vials each contain different concentration of human FGF-21 in a lyophilized bovine serum based matrix with a non-azide, non-mercury preservative. Refer to vials for exact concentration range for each control. The controls are ready to use. Both controls should be stored at 2 – 8°C and are stable until the expiration date on the kit box.

SAFETY PRECAUTIONS

The Human Intact Fibroblast Growth Factor 21 (FGF-21) Assay Kit reagents must be used in a professional laboratory environment and is for Research Use Only and is not to be used in diagnostic procedures. Only source material from which reagents of bovine serum was derived in the contiguous 48 United States. It was obtained only from donor health animals maintained under veterinary supervision and found free of contagious diseases. Wear gloves while performing this assay and handle these reagents as if they were potentially infectious. Avoid contact with reagents containing TMB, hydrogen peroxide, or sulfuric acid. TMB may cause irritation to skin and mucous membranes and cause an allergic skin reaction. TMB is a suspected carcinogen. Sulfuric acid may cause severe irritation on contact with skin. Do not get in eyes, on skin, or on clothing. Do not ingest or inhale fumes. On contact, flush with copious amounts of water for at least 15 minutes. Use Good Laboratory Practices.

MATERIALS REQUIRED BUT NOT PROVIDED

- Precision single channel pipettes capable of delivering 25 µL, 50 µL, 100 µL, and 1000 µL
- Repeating dispenser suitable for delivering 100 µL.
- Disposable pipette tips suitable for above volume dispensing.
- Disposable 12 x 75 mm or 13 x 100 glass or plastic tubes.
- Disposable plastic 100 mL and 1000 mL bottle with caps.
- Aluminum foil.
- Deionized or distilled water.
- Plastic microtiter well cover or polyethylene film.
- ELISA multichannel wash bottle or automatic (semi-automatic) washing system.
- Spectrophotometric microplate reader capable of reading absorbance at 450 nm.

SPECIMEN COLLECTION

Only 50 µL of human EDTA-plasma is required for human FGF-21 measurement in singlet. No special preparation of individual is necessary prior to specimen collection. Whole blood should be collected with lavender-top Vacutainer and separate the plasma from cells by centrifugation (850 – 1500xg for 10 minutes). The plasma should be separated from the cells right after collection or at least within one hour of blood collection. The plasma should be transferred to a clean test tube right after centrifugation. Plasma samples should be stored at – 20°C if the assay is not to be performed within 48 hours. Avoid more than three times freeze-thaw cycles of specimen.

Serum sample can also be used for FGF-21 measurement. Serum sample collection should perform as suggested by manufacturer of the sample collection tubes.

SPECIMEN SHIPMENT

Collected EDTA-plasma or serum samples should be shipped to designated laboratory in frozen condition with dry ice. In case frozen condition is not available, samples should be shipped at



room temperature in an insulated container for maximum 48 hour delivery. Samples must not be shipped refrigerated, such as, with blue ice pack.

REAGENT PREPARATION

1. Prior to use allow all reagents to come to room temperature. Reagents from different kit lot numbers should not be combined or interchanged.
2. ELISA Wash Concentrate must be diluted to working solution prior use. Please see REAGENTS section for details.
3. Reconstitute kit standards and controls by adding 0.5 mL distilled water into each vial. Gently mix and dissolve the entire particle before use. The reconstituted standards and controls should be stored at -20°C right after use.
4. Prepare working human FGF-21 tracer antibody (Cat# 30620) by 1:21 fold dilution of the conjugation antibody with the FGF-21 Tracer Antibody Diluent (Cat# 30600). Following is a table that outlines the relationship of strips used and antibody mix prepared.

Strip no.	FGF-21 Tracer Antibody Diluent	FGF-21 Tracer Antibody
1	500 µL	25 µL
2	1000 µL	50 µL
3	1500 µL	75 µL
4	2000 µL	100 µL
5	2500 µL	125 µL
6	3000 µL	150 µL
7	3500 µL	175 µL
8	4000 µL	200 µL
9	4500 µL	225 µL
10	5000 µL	250 µL
11	5500 µL	275 µL
12	6000 µL	300 µL

Note: this antibody mixture must be freshly prepared right before testing.

Assay Procedure

1. Place a sufficient number of murine IgG coated microwell strips/wells in a holder to run human FGF-21 standards, controls and unknown samples in duplicate.
2. Test Configuration

ROW	STRIP 1	STRIP 2	STRIP 3
A	STD 1	STD 5	SAMPLE 1
B	STD 1	STD 5	SAMPLE 1
C	STD 2	STD 6	SAMPLE 2
D	STD 2	STD 6	SAMPLE 2
E	STD 3	C 1	SAMPLE 3
F	STD 3	C 1	SAMPLE 3



G	STD 4	C 2	
H	STD 4	C 2	

3. Add **50 µL** of standards, controls and patient plasma/serum samples into the designated microwell.
4. Add **50 µL** of 1:21 diluted tracer antibody to each well
5. Cover the plate with one plate sealer and incubate plate with orbital shaking 170 rpm at room temperature for **2 hours**.
6. Remove plate sealer. Aspirate the contents of each well. Wash each well 5 times by dispensing 350 µL of working wash solution into each well and then completely aspirating the contents. Alternatively, an automated microplate washer can be used.
7. Add **100 µL** of ELISA HRP Substrate into each of the wells.
8. Cover the plate with one plate sealer and also with aluminum foil to avoid exposure to light. Incubate plate at room temperature for **20 minutes**.
9. Remove the aluminum foil and plate sealer. Add **100 µL** of ELISA Stop Solution into each of the wells. Mix gently.
10. Read the absorbance at 450/650 nm within 10 minutes in a microplate reader

PROCEDURAL NOTES

1. It is recommended that all standards, controls and unknown samples be assayed in duplicate. The average absorbance reading of each duplicate should be used for data reduction and the calculation of results.
2. Keep light sensitive reagents in the original amber bottles.
3. Store any unused antibody coated strips in the foil zipper bag with desiccant to protect from moisture.
4. Careful technique and use of properly calibrated pipetting devices are necessary to ensure reproducibility of the test.
5. Incubation times or temperatures other than those stated in this insert may affect the results.
6. Avoid air bubbles in the microwell as this could result in lower binding efficiency and higher CV% of duplicate reading
7. All reagents should be mix gently and thoroughly prior use. Avoid foaming.

INTERPRETATION OF RESULTS

1. Calculate the average absorbance for each pair of duplicate test results.
2. Subtract the average absorbance of the STD 1 (0 ng/mL) from the average absorbance of all other readings to obtain corrected absorbance.
3. The standard curve is generated by the absorbance of all standards. Appropriate computer assisted data reduction programs may also be used for the calculation of results.

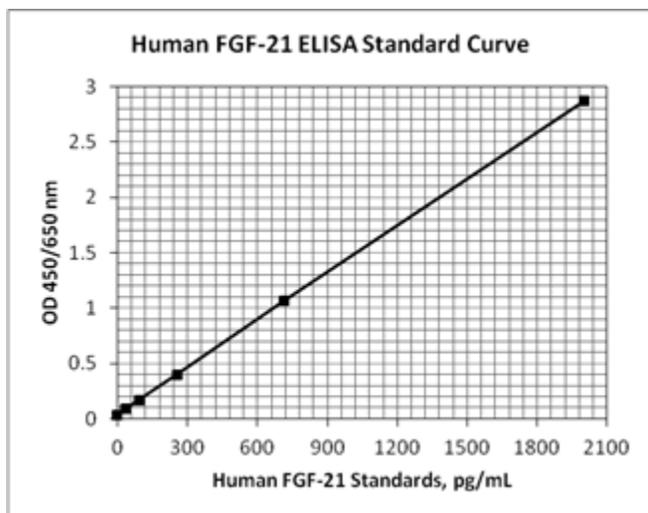
The human intact FGF-21 concentrations for the controls and patient samples are read directly from the standard curve using their respective corrected absorbance.



EXAMPLE DATA AND STANDARD CURVE

A typical absorbance data and the resulting standard curve from Human Intact Fibroblast Growth Factor 21 (FGF-21) ELISA Assay Kit are represented. This curve should not be used in lieu of standard curve run with each assay.

Well I.D.	OD 450/650 nm Absorbance			Results pg/mL
	Readings	Average	Corrected	
0 pg/mL	0.037 0.036	0.037	0.000	
32.5 pg/mL	0.087 0.086	0.087	0.050	
91 pg/mL	0.172 0.169	0.170	0.133	
255 pg/mL	0.398 0.399	0.399	0.302	
714 pg/mL	1.067 1.069	1.068	1.031	
2000 pg/mL	2.835 2.903	2.869	2.946	
Control 1	0.126 0.129	0.127	0.371	60.83
Control 2	0.736 0.721	0.729	1.200	481.29



EXPECTED VALUES

Thirty two normal adult plasma samples were measured with this human intact FGF-21 ELISA. The normal range was found to be less than 200 pg/mL. It is strongly recommended that each laboratory should establish its own normal range based on normal donor EDTA-plasma or serum samples.

LIMITATION OF THE PROCEDURE

1. Since there is no Gold Standard concentration available for human intact FGF-21 measurement, the values of assay standards were established by correlation to a highly purified FGF-21 standard.
2. For sample values reading greater than highest standard, it is recommend to re-assay samples with dilution.
3. Bacterial or fungal contamination of plasma specimens or reagents, or cross contamination between reagents may cause erroneous results.
4. Water deionized with polyester resins may inactive the horseradish peroxidase enzyme.

QUALITY CONTROL

To assure the validity of the results each assay should include adequate controls with known human intact FGF-21 levels. We recommend that all assays include the laboratory's own FGF-21 controls in addition to those provided with this kit.

PERFORMANCE CHARACTERISTICS

Sensitivity (LoD)

The sensitivity (lowest of of this human intact FGF-21 ELISA as determined by the corresponding OD value of 2 fold standard deviation above the mean on 20 duplicate determination of zero standard is 1.7 pg/mL.

High Dose "hook" effect

This assay has showed that it did not have any high dose "hook" effect up to 20,000 pg/mL.

Precision

The intra-assay precision is validated by measuring three donor EDTA-plasma samples in a single assay with 16-replicate determinations.



Mean Human Intact FGF-21 Value (pg/mL)	CV (%)
63.2	5.7
171	4.2
480	5.4

The inter-assay precision is validated by measuring three control samples in duplicate in 12 individual assays.

Mean Human Intact FGF-21 Value (pg/mL)	CV (%)
69.8	6.9
181	3.0
486	1.9

Linearity

Two human EDTA-plasma samples were diluted with 0.01M PBS, pH 7.4 and assayed. The results in the value of pg/mL are as follows:

#	DILUTION	OBSERVED VALUE	EXPECTED VALUE	RECOVERY %
1	Neat	286	-	-
	1:2	138	143	96
	1:4	75	72	104
	1:8	37.9	36	105
	1:16	19.5	18	108
2	Neat	61.8	-	-
	1:2	32.1	30.9	104
	1:4	15.9	15.5	103
	1:8	7.2	7.7	94

Spike Recovery

Two patient samples were spiked with various amounts of human intact FGF-21 (1 vol. + 1 vol. mixture) and assayed. The results in the value of ng/mL are as follows:

#	Orig. Value	Amount Spiked	Observed Value	Expected Value	Recovery %
1	45.9 (serum)	91	64.9	68.5	95
		255	150	151	100
		714	388	380	102
2	40.4 (plasma)	91	71.2	65.7	108
		255	148	148	100



REFERENCES

1. Yie J, et al. FGF21 N- and C-termini play different roles in receptor interaction and activation. FEBS Lett. 2009 Jan 5;583:19-24.
2. Micanovic R, et al. Different roles of N- and C- termini in the functional activity of FGF21. J Cell Physiol. 2009 May;219(2):227-34.
3. Yusuke Murata, et al. FGF21 as an Endocrine Regulator in Lipid Metabolism: From Molecular Evolution to Physiology and Pathophysiology. Journal of Nutrition and Metabolism, Vol 2011, Article ID 981315, 8 pages

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, please contact the Technical Service Team at Eagle Biosciences, Inc. at info@eaglebio.com or at 866-411-8023.