



**EAGLE**  
BIOSCIENCES

## **Mouse FGF-21 ELISA Assay Kit**

Catalog Number: FGF11-K01

For the quantitative determination of mouse FGF-21 concentrations  
in serum, plasma and cell culture supernate samples.

This package insert must be read in its entirety before using this product.  
Use only the current version of product data sheet enclosed with the kit.

**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](http://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	2
PREPARATION OF STANDARDS AND SAMPLES ASSAY	3
PROCEDURE	4
CALCULATION	5
TYPICAL STANDARD CURVE	5
ASSAY CHARACTERISTICS	6
REFERENCES	7
SUMMARY OF ASSAY PROCEDURE	8
	9



## **ASSAY BACKGROUND**

Fibroblast growth factor 21 (FGF-21) is a novel protein that has been implicated in the regulation of lipid and glucose metabolism under fasting and ketotic conditions<sup>1,2</sup>. In murine models, FGF-21 is predominantly expressed in liver, but it also expressed in adipose tissue and pancreatic  $\beta$  cells<sup>3,4</sup>. FGF-21 stimulates glucose uptake in adipocytes. It also protects animals from diet-induced obesity when overexpressed in transgenic mice and lowers blood glucose and triglyceride levels when administered to diabetic rodents<sup>5</sup>. When administered daily for 6 weeks to diabetic rhesus monkeys, FGF-21 caused a dramatic decline in fasting plasma glucose, fructosamine, triglycerides, insulin, and glucagon<sup>6</sup>. Furthermore, elevated plasma FGF-21 concentrations in humans appear to be related to the presence of hepatic and peripheral insulin resistance<sup>7</sup>.

## **ASSAY PRINCIPLE**

This assay is a quantitative sandwich ELISA. The immunoplate is pre-coated with a rabbit polyclonal antibody specific for mouse FGF-21. Standards and samples are pipetted into the wells and any mouse FGF-21 present is bound by the immobilized antibody. After washing away any unbound substances, a biotin labelled polyclonal antibody specific for mouse FGF-21 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 FGF-21 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 FGF-21, the unknown sample concentration can be interpolated from a reference curve included in each assay.

## **INTENDED USE**

The Eagle Biosciences Mouse FGF-21 ELISA Assay Kit is designed for quantification of mouse FGF-21 in serum, plasma and cell culture samples. The Mouse FGF-21 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 FGF-21, 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 FGF-21, 0.12 ml.
5. Mouse FGF-21 standard-2000 pg of recombinant Mouse FGF-21 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 FGF-21 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.*

### **1. 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.

### **2. 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.

### **3. 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  $\mu$ 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.

### **4. 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  $\mu$ 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 FGF-21 standards:** Reconstitute the lyophilised standard with 1 ml of 1×Assay buffer to generate a standard stock solution of 2000 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
2000 pg/ml stock	-	2000 pg/ml
250 $\mu$ l of 2000 pg/ml std	250 $\mu$ l	1000 pg/ml
250 $\mu$ l of 1000 pg/ml std	250 $\mu$ l	500 pg/ml
250 $\mu$ l of 500 pg/ml std	250 $\mu$ l	250 pg/ml
250 $\mu$ l of 250 pg/ml std.	250 $\mu$ l	125 pg/ml
250 $\mu$ l of 125 pg/ml std.	250 $\mu$ l	62.5 pg/ml
250 $\mu$ l of 62.5 pg/ml std.	250 $\mu$ l	31.2 pg/ml

1x Assay buffer serves as the zero standard (0 ng/ml).

**Note:** The reconstituted standard stock should be aliquoted and frozen at -80°C for 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 2-fold dilution in 1×Assay buffer.



## ASSAY PROCEDURE

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

1. Add 100  $\mu$ l 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  $\mu$ 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  $\mu$ 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  $\mu$ 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  $\mu$ l of Substrate solution to each well, incubate at room temperature for 15 minutes. **Protect from light.**
8. Add 100  $\mu$ 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 FGF-21 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 FGF-21 concentration of samples from standard curve and multiply the value by the dilution factor.

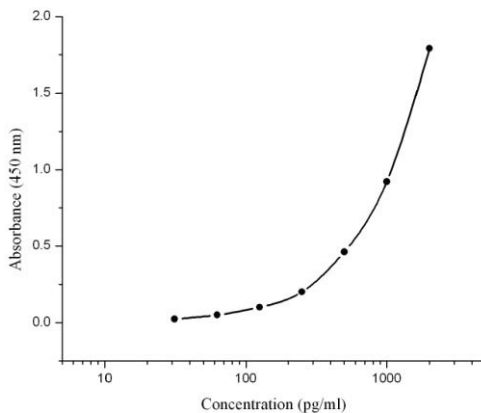


### TYPICAL STANDARD CURVE

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

FGF-21 (pg/ml)	Absorbance (450 nm)	Blanked Absorbance
0	0.109	0
31.2	0.132	0.023
62.5	0.158	0.049
125	0.209	0.1
250	0.309	0.2
500	0.571	0.462
1000	1.031	0.992
2000	1.9	1.791

Mouse FGF-21 standard curve (4-parameter)







## ASSAY CHARACTERISTICS

### A. Sensitivity:

The lowest level of mouse FGF-21 that can be detected by this assay is 31.2 pg/ml.

### B. Specificity:

Cross reactivity of recombinant proteins

Analyte	Cross Reactivity
Human FGF-21	Yes
Mouse Mup-1	No
Mouse LCN2	No
Mouse Adiponectin	No
Mouse ANGPL4	No

### C. Precision:

Intra-assay Precision (Precision within an assay) C.V <8%.

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

### D. Recovery:

The recovery of the assay was determined by adding various amounts FGF21 to a sample. The measured concentration of the spiked sample in the assay was compared to the expected concentration. The average recovery was 92%.

### E. Linearity:

To assess the linearity of the assay, a sample with high level of FGF-21 was serially diluted with the 1×Assay buffer to produce samples with values within the dynamic range of the assay.

Dilution	Measured (pg/ml)	Expected (pg/ml)	Recovery (%)
1/2	1360	1360	100
1/4	680	680	101
1/8	380	340	111
1/16	200	170	115



## REFERENCES:

- [1] Kharitononkov A, et al. (2005) *J Clin Invest*; 115: 1627– 1635.
- [2] Badman MK, et al. (2007) *Cell Metab*; 5: 426– 437.
- [3] Nishimura T, et al. (2000) *Biochim Biophys Acta*; 1492: 203– 206.
- [4] Kurosu H, et al. (2007) *J Biol Chem*; 282: 26687– 26695
- [5] Kharitononkov A, et al. (2005) *J. Clin. Invest.* 115: 1627–35.
- [6] Kharitononkov A, et al. (2007) *Endocrinology*;148:774-81.
- [7] Chavez AO, et al. (2009) *Diabetes Care*; 32:1542-6.



## SUMMARY OF ASSAY PROCEDURE

Add 100  $\mu$ l of Standard or sample to each well.



Incubate at room temperature for 1 hour.



Aspirate and wash each well three times.



Add 100  $\mu$ 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  $\mu$ 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  $\mu$ l of Substrate solution to each well.



Incubate at room temperature for 15 minutes.



Add 100  $\mu$ 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](mailto: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](http://www.EagleBio.com) • [info@eaglebio.com](mailto:info@eaglebio.com)



**EAGLE**  
BIOSCIENCES