



Irisin ELISA Kit

(For Human, Mouse, Rat, Canine)

Competitive Enzyme Immunoassay for the quantification of Irisin in plasma (EDTA) samples

Catalog number: ARG81195

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

For research use only. Not for use in diagnostic procedures.

TABLE OF CONTENTS

SECTION	Page
INTRODUCTION	3
PRINCIPLE OF THE ASSAY	3
MATERIALS PROVIDED & STORAGE INFORMATION	4
MATERIALS REQUIRED BUT NOT PROVIDED	4
TECHNICAL HINTS AND PRECAUTIONS	5
SAMPLE COLLECTION & STORAGE INFORMATION	5
REAGENT PREPARATION.....	5
ASSAY PROCEDURE	7
CALCULATION OF RESULTS	9
EXAMPLE OF TYPICAL STANDARD CURVE	9
QUALITY ASSURANCE.....	10

MANUFACTURED BY:

Arigo Biolaboratories Corporation

Address: No. 22, Ln. 227, Gongyuan Rd., Hsinchu City 300, Taiwan

Phone: +886 (3) 562 1738

Fax: +886 (3) 561 3008

Email: info@arigobio.com

INTRODUCTION

This gene encodes a secreted protein that is released from muscle cells during exercise. The encoded protein may participate in the development of brown fat. Translation of the precursor protein initiates at a non-AUG start codon at a position that is conserved as an AUG start codon in other organisms. Alternative splicing results in multiple transcript variants. [provided by RefSeq, Jun 2013]

Irisin: mediates beneficial effects of muscular exercise. Induces browning of white adipose tissue by stimulating UCP1 expression, at least in part, via the nuclear receptor PPARA (By similarity) [UniProt]

PRINCIPLE OF THE ASSAY

This is a Competitive Enzyme Immunoassay for the quantification of Irisin in Serum and plasma samples

This assay employs the competitive enzyme immunoassay technique. A secondary antibody has been pre-coated onto a microtiter plate and non-specific binding sites were blocked. Fc regions of primary antibodies specific for target peptides can bind to the secondary antibodies on microtiter plate. The Fab regions of primary antibodies are competitively bound by biotinylated peptide and targeted peptides in samples or standards. The biotinylated peptide interacts with streptavidin-horseradish peroxidase (SA-HRP) which catalyzes the substrate solution. The reaction is monitored by a color change which is readable at OD of 450 nm±2 nm. Quantification of unknown samples is achieved by comparing their absorbance with a reference curve prepared

Irisin ELISA Kit ARG81195

with known standards. The intensity of color development is inversely proportional to the amount of Irisin in the samples.

MATERIALS PROVIDED & STORAGE INFORMATION

Store the unopened kit at 2-8 °C. Use the kit before expiration date.

Component	Quantity	Storage information
Microtiter Plate	12 x 8 wells	4°C
20X assay buffer concentrate	50 ml	4°C
Acetate Plate sealer (APS)	3 pieces	4°C
Primary antibody (Rabbit anti-peptide IgG)	1 vial	4°C
Standard peptide	1 vial	4°C
Biotinylated peptide	1 vial	4°C
Streptavidin-HRP	30 µl	4°C
Positive control	2 vials	4°C
Substrate Solution (TMB)	12 ml	4°C, ready for use
2N HCl	15 ml	4°C, ready for use

MATERIALS REQUIRED BUT NOT PROVIDED

- Microplate reader capable of measuring absorbance at 450nm
- Pipettes and pipette tips
- Multi-channel pipette capable of dispensing 50-100µl
- Solution reservoir
- Aprotinin (0.6TIU/ml of blood) (optional)
- Deionized or distilled water
- Orbital microplate shaker capable of 300-400rpm (Recommended)
- Automated microplate washer (optional)

TECHNICAL HINTS AND PRECAUTIONS

- Wear protective gloves, clothing, eye, and face protection especially while handling blood or body fluid samples.
- Store the kit at 4°C at all times.
- Briefly spin down the standards and solutions before use.
- Ensure complete reconstitution and dilution of reagents prior to use.
- It is highly recommended that the standards, samples and controls be assayed in duplicates.
- Change pipette tips between the addition of different reagent or samples.

SAMPLE COLLECTION & STORAGE INFORMATION

The sample collection and storage conditions listed below are intended as general guidelines.

Blood collection- Collect blood samples into tubes containing EDTA. Gently rock the tubes several times immediately to prevent coagulation. Transfer the blood from tubes to centrifuge tubes containing aprotinin (0.6 TIU/ml blood) and gently rock for a few times to inhibit activity of proteinases. Centrifuge the blood at 1600 x g for 15 minutes at 4°C and collect plasma. Plasma can be kept at - 80°C up to 1 month.

REAGENT PREPARATION

- **1X Assay Buffer:** Dilute 20X Assay Buffer into distilled water to yield 1X Assay Buffer (e.g. 50ml of 20X Assay Buffer + 950ml distilled water). Keep 1X Assay Buffer at 4°C. If crystals appear in 20X Assay Buffer, warm the

Irisin ELISA Kit ARG81195

buffer in warm water bath for 30 minutes or until crystals disappear. Mix well before use.

- **Primary antibody:** Reconstitute the Primary antibody vial with 5 ml of 1X assay buffer. Allow to sit for at least 5 minutes and mix well to dissolve completely before use. Keep rehydrated solution at 4°C and use as soon as possible.
- **Biotinylated peptide:** Reconstitute the Biotinylated peptide vial with 5 ml of 1X assay buffer. Allow to sit for at least 5 minutes and mix well to dissolve completely before use. Keep rehydrated solution at 4°C and use as soon as possible.
- **Positive control:** Reconstitute the Positive control vial with 200 µl of 1X assay buffer. Allow to sit for at least 5 minutes and mix well to dissolve completely before use. Keep rehydrated solution at 4°C and use as soon as possible.
- **Sample:** It is recommended each laboratory to do pre-test to determine the appropriate dilution factors for the samples to be measured to ensure that the samples are within the dynamic range of the standard curve. Samples can be diluted with 1x assay buffer if needed. For the calculation of the concentrations this dilution factor has to be taken into account.
- **Standard peptide:** Reconstitute the Standard peptide vial with 1 ml of 1X assay buffer. Vortex it. The concentration of this stock solution is 1000 ng/ml. Allow to sit for 10 minutes at room temperature (20-23°C) to dissolve completely. Mix well and spin down before use. Dilute standard solutions according to the table below and make serial dilutions of 1000

Irisin ELISA Kit ARG81195

ng/ml, 100 ng/ml, 10 ng/ml, 1 ng/ml and 0.1 ng/ml.

Tube No.	Standard volume	1X Assay Buffer	Concentrations (ng/ml)
Stock	1000 μ l	0 μ l	1000
1	100 μ l of stock	900 μ l	100
2	100 μ l of Tube 1	900 μ l	10
3	100 μ l of Tube 2	900 μ l	1
4	100 μ l of Tube 3	900 μ l	0.1

ASSAY PROCEDURE

Note: All materials should be equilibrated to room temperature (20-23°C) before use or opening. It is recommended that the solutions be used as soon as possible after rehydration. Samples and controls should be assayed in duplicates. Unused microplate strips should be placed in the foil pouch with a dessicant and stored at 4°C. Do not allow moisture to enter the wells.

1. Add 50 μ l of 1X assay buffer as Total Binding. Keep another Two Empty Wells as Blank.
2. Add 50 μ l of peptide standards (1000 ng/ml, 100 ng/ml, 10 ng/ml, 1 ng/ml and 0.1 ng/ml), 50 μ l positive controls or 50 μ l samples into corresponding wells. It is advisable to assay each condition in duplicates.
3. Add 25 μ l of primary antibody into each well except the Blank wells.
4. Add 25 μ l of Biotinylated peptide into each well except the Blank wells. It is not recommended to use a multi-channel pipette to load the primary antibody and biotinylated peptide.
5. Seal the microtiter plate with acetate plate sealer (APS). Incubate for 2 hours at room temperature (20-23°C). Orbital shaking at 300-400 rpm with a microplate shaker is recommended.

Irisin ELISA Kit ARG81195

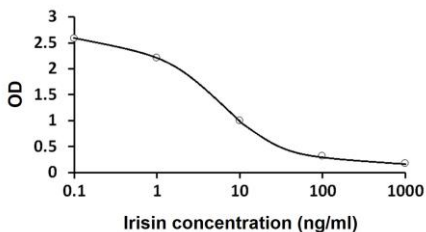
6. Centrifuge Streptavidin-HRP vial to spin down the solution in the vial and pipette 12 μ l of Streptavidin-HRP into 12ml of 1X Assay buffer. Vortex thoroughly. Prepare fresh.
7. Remove sealer from plate.
8. Aspirate each well and wash, repeating the process 3 times for a total 4 washes. Wash by filling each well with 1 \times Assay Buffer (350 μ l) using a squirt bottle, manifold dispenser, or autowasher. Complete removal of liquid at each is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating, decanting or blotting against clean paper towels.
9. Add 100 μ l diluted Streptavidin-HRP solution into each well.
10. Reseal the plate with sealer. Incubate for 1 hour at room temperature (20-23°C). Orbital shaking at 300-400 rpm with a microplate shaker is recommended.
11. Remove sealer from plate.
12. Wash as according to step 8.
13. Add 100 μ l TMB substrate solution into each well.
14. Reseal the plate with sealer. Incubate for 1 hour at RT. Orbital shaking at 300-400 rpm with a microplate shaker is recommended. (Protect from light)
15. Remove sealer from plate. Add 100 μ l 2N HCl into all wells to stop the reaction, gently tap the plate to ensure thorough mixing. The color in the well should change from blue to yellow. Go to the next step within 20 minutes.
16. Read the OD with a microplate reader at 450 nm immediately.

CALCULATION OF RESULTS

1. Calculate the average absorbance values for each set of standards, controls and patient samples.
2. Using semi-log graph paper, construct a standard curve by plotting the mean absorbance obtained from each standard against its concentration with absorbance value on the vertical (Y) axis and concentration on the horizontal (X) axis.
3. Using the mean absorbance value for each sample determine the corresponding concentration from the standard curve.
4. Automated method: The results in the IFU have been calculated automatically using a 4 PL (4 Parameter Logistics) curve fit. 4 Parameter Logistics is the preferred method. Other data reduction functions may give slightly different results.
5. If samples have been diluted prior to the assay, the measured concentration must be multiplied by their respective dilution factors.

EXAMPLE OF TYPICAL STANDARD CURVE

The following data is for demonstration only and cannot be used in place of data generations at the time of assay.



QUALITY ASSURANCE

Sensitivity

Standard Range: 0.1- 1000 ng/ml

Linear Range: 1.29- 27.5 ng/ml

Sensitivity: 1.29 ng/ml

Cross Reactivity

The cross-reactivity with the following factors were as below:

Irisin, Recombinant (Human, Rat, Mouse, Dog): 100%

Irisin (aa. 42-112) (Human, Rat, Mouse, Dog): 100%

FNDC5, isoform 4 recombinant (Human, Rat, Mouse, Dog): 9%

Irisin precursor, C-terminal 48mer / FNDC5 (aa. 165-212) (Human) /

FNDC5(aa. 162-209) (Rat, Mouse): 0%

Irisin (aa. 42-95) (Human, Rat, Mouse, Dog): 0%

FNDC4 (aa. 47-101) (Human) / FNDC4 (aa. 43-97) (Mouse): 0%

FNDC4 (aa. 103-158) (Human) / FNDC4 (aa. 99-154) (Mouse): 0%

FNDC4 / FRCP1 (aa. 190-234) (Human): 0%

Intra-assay and Inter-assay precision

Intra-assay variation: <10%

Inter-assay variation: <15%

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



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.

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.