



EAGLE
BIOSCIENCES

Active Ghrelin ELISA Assay Kit

Catalog Number:

GHA31-K01 (1 x 96 wells)

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

v. 1.0 (08 OCT 24)

Eagle Biosciences, Inc.

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

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

www.EagleBio.com



INTRENDED USE

The Active Ghrelin ELISA Assay Kit measures the active form of ghrelin in human, rat, or mouse plasma. The Active Ghrelin ELISA Assay Kit is for research use only and not to be used for diagnostic procedures.

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

ASSAY BACKGROUND

Ghrelin, a novel growth hormone releasing peptide is an acylated peptide that stimulates the release of growth hormone from pituitary. It was isolated from rat stomach and the structure was determined as a peptide consisting of 28-amino acid by Dr. Kenji Kangawa (National Cardiovascular Center in Japan). The Ser3 residue of Ghrelin is modified by n-octanoic acid, a modification necessary for hormone activity.

ASSAY PRINCIPLE

This Active Ghrelin ELISA kit measures the active form of Ghrelin based on the principle of 2 Site Sandwich enzyme-linked immunosorbent assay (ELISA). It can detect not only octanoylated human Ghrelin but also octanoylated rat/mouse Ghrelin (1-28). This kit is manufactured using the high specific antibody pairs generated by Dr. Kangawa and by following his protocol. (patent pending PCT WO 01/07475 A1).

Active Ghrelin ELISA Assay Kit	Contents
<ul style="list-style-type: none">♦ The assay can measure active ghrelin within the range of 2.9~184 fmol/mL♦ The assay is completed within 2 hr + 1 hr + 0.5 hr.♦ With one assay kit, 40 samples can be measured in duplicate.♦ Test Sample: human/rat/mouse plasma♦ Sample Volume: 50 uL♦ The 96-wells plate in kit is consisted by 8-wells strips, and the strips can be used separately.♦ Stability and Storage Store all of the components at 2-8°C The kit is stable under the condition for 12 months from the date of manufacturing The expiry date is stated on the label of the kit.	<ol style="list-style-type: none">1. Monoclonal antibody against active ghrelin coated plate2. Standard3. HRP labeled monoclonal antibody against ghrelin solution4. Substrate solution (TMB)5. HRP dilution buffer6. Stopping solution7. Assay buffer8. Washing buffer (concentrated)9. Transparent sheet



COMPONENTS

Component	Form	Quantity	Main Ingredient
1. Antibody coated plate	Microtiter plate	1 plate (96 wells)	Mouse anti-active ghrelin monoclonal antibody coated
2. Standard	Lyophilized	1 vial (184 fmol)	Synthetic human active ghrelin
3. HRP labeled antibody solution	Liquid	1 tube (0.15 mL)	HRP labeled mouse anti-ghrelin monoclonal antibody
4. Substrate Solution	Liquid	2 bottles (11 mLx2)	3,3',5,5'-Tetramethylbenzidine (TMB)
5. HRP dilution buffer	Liquid	1 bottle (22 mL)	Phosphate buffer
6. Stopping Solution	Liquid	1 bottle (6 mL)	1 mol/L H ₂ SO ₄
7. Assay Buffer	Liquid	1 bottle (22 mL)	Buffer containing a reaction accelerator
8. Washing Buffer Concentrate	Liquid	1 bottle (40 mL)	Concentrated phosphate buffer
9. Transparent sheet		3 pieces	

MATERIALS REQUIRED

1. Plate Washer
2. Plate Reader (450 nm measurement available)
3. Vortex Mixer

PREPARATION OF SAMPLE

Ghrelin is very unstable. Be careful to avoid any fragmentation or inactivation. All biological fluid should be treated with protease inhibitor such as aprotinin. It is also required to inhibit the esterase activity. Standard procedure of human blood sample preparation is as described below.

Collect into the bleeding tubes which contain 500 KIU (Kallikrein Inhibitor Unit) of aprotinin and 1.25 mg of EDTA-2Na per 1 mL of whole blood. Rock the tubes gently and then immediately centrifuge the blood sample (1500xg, 15 min at 4°C). Earned plasma should be immediately treated with 1/10 volume of 1 mol/L HCl. Sample must be kept below -40°C for long term storage.

REAGENT PREPARATION

1. Dilute the washing buffer concentrate with x20 volume of distilled water. Store the diluted washing buffer in refrigerator and use within 2 weeks.
2. Reconstitute the Standard (Lyophilized) with 1 mL of distilled water (→ Standard #1). **Keep still approximately 10 minutes and vortex well.**



Then dilute the standard as follows:

Standard No.	Std Vol.	Assay Buffer
#2	#1 → 500 μL	500 μL
#3	#2 → 500 μL	500 μL
#4	#3 → 500 μL	500 μL
#5	#4 → 500 μL	500 μL
#6	#5 → 500 μL	500 μL
#7	#6 → 500 μL	500 μL

The lyophilized standard contains 184 fmol of human active ghrelin

The dilution procedure described above is for 2.9~184 fmol/mL of the standard curve.

3. **Dilute only the required volume of the HRP labeled antibody solution with x200 volume of the HRP dilution buffer and vortex well.** (Prepare the diluted HRP labeled antibody solution more than one hour before using and use it within a day.)

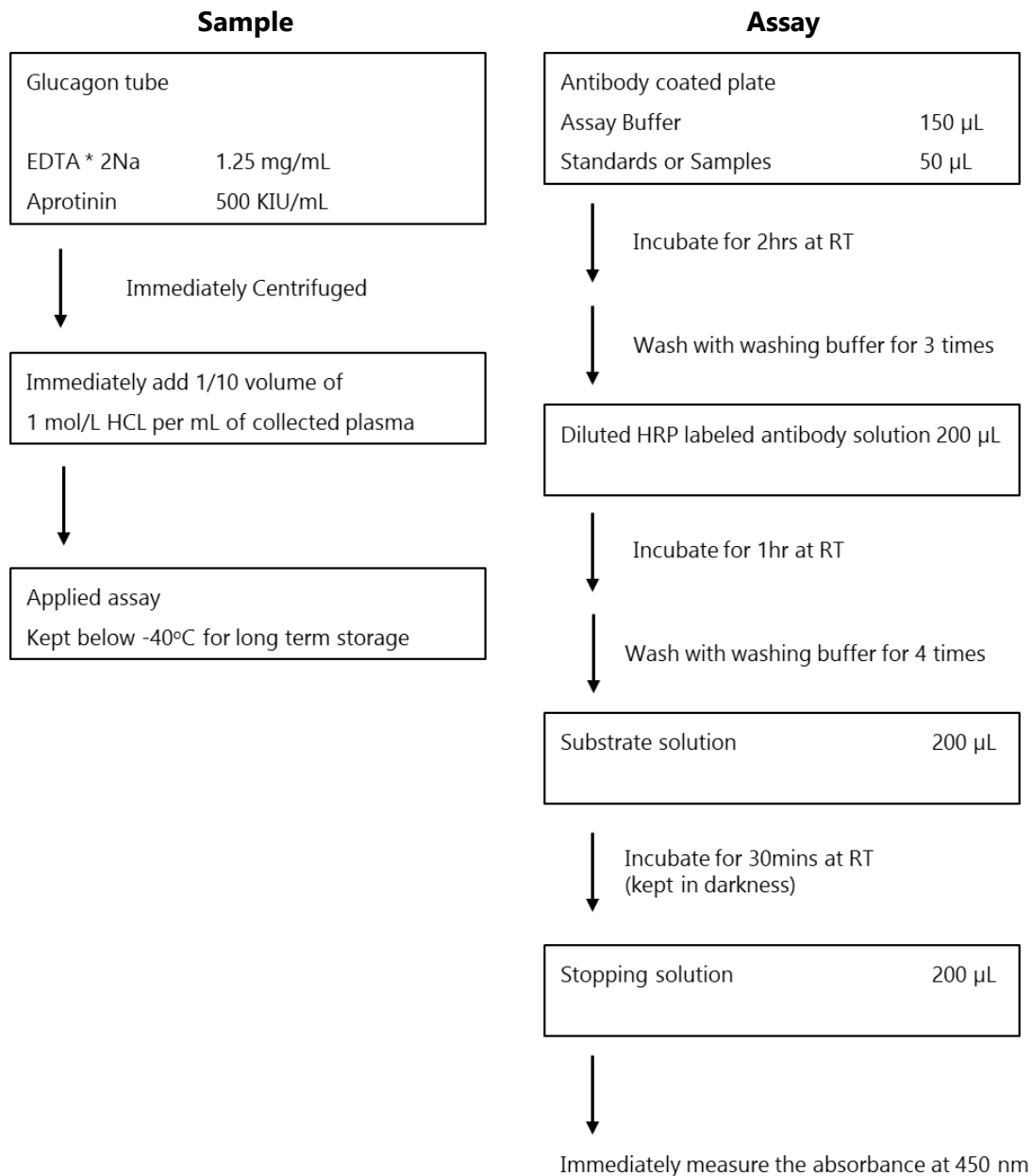
ASSAY PROCEDURE

Pre-warm all reagents to room temperature prior to setting up the assay. Do not dry up the wells during a measurement.

1. 150 μL of assay buffer is poured into the testing well. 50 μL of samples and standards are added into the each well. As a "Blank", 50 μL of assay buffer is added into the testing well. Manufacturer recommends to test duplicate for each sample. Plate is covered with transparent sheet and is incubated for 2 hours at RT.
2. Aspirate samples from wells and wash by washing buffer for 3 times. **Washing buffer volume: 350 μL.** Keep 1 min of interval before removing the washing buffer from wells. For removing the remnant completely, testing plate is tapped on a paper towel upside down. 200 μL of diluted HRP labeled antibody solution is poured into the wells. Testing plate is covered with transparent sheet and is incubated for 1 hour at RT.
3. Aspirate samples from wells and wash by washing buffer for 4 times. **Washing buffer volume: 350 μL.** Keep 1 min of interval before aspirating the washing buffer from wells. For removing the remnant completely, testing plate is tapped on a paper towel upside down. 200 μL of substrate solution is poured into each well and is incubated for 30 min at RT with shading. After the incubation, 50 μL of stopping solution is added to each well to stop reaction. Then shake the plate gently.
4. Measure the absorbance of each well at 450 nm immediately.
5. The results of unknown samples can be calculated with any computer program having a 4 (or 5)-parameter logistic function. Otherwise plot the standard concentration (X-axis) and its corresponding absorbance (Y-axis). The concentration of active ghrelin in unknown sample is determined by plotting the sample's absorbance on the standard curve. When HCl is added to the samples, multiply the results by 1.1 to offset the dilution.



ACTIVE GHRELIN ELISA PROTOCOL



PRECAUTION

1. This kit is for research purpose only. Not for diagnostic use.
2. Warning – potential biohazardous material. Samples should be handled at the Biosafety level 2 as recommended for any potentially infection human serum or blood samples in the Centers for Disease Control/National Institutes of Health manual "Biosafety in Microbiological and Biomedical Laboratories", 1984. In addition, handle and dispose of the Antibody coated plate as well as all material coming into contact with them or with the ample as if capable of transmitting infection.



3. Substrate solution is sensitive to contamination from a variety of oxidizing agents such as bacteria, dust, metals and commonly used laboratory glassware. Avoid contacting with any potential source of these contaminations. Substrate solution is also sensitive to light. Avoid unnecessary exposure to light.
4. Stopping solution contains 1M sulfuric acid solution. Sulfuric acid is corrosive and can cause eye and skin burns. Avoid contact with skin and eyes. To prevent any contact, wear protective equipment such as safety gloves, rubber gloves, as appropriate.
5. Reagents are stored between +2°C and +8°C. Before the measurement, all reagents must be equilibrated to room temperature. Put unused strips back in the aluminum pouch immediately, because strips are affected by humidity.
6. Do not mix the reagents from different kits unless they have the same lot numbers.
7. Do not use reagents after the expiration date printed on the label.
8. Occasionally, the assay buffer and the washing buffer concentrate generate some precipitates. However, they can be resolved by raising the temperature from room temperature to approximately 30°C. After then, you can use their buffers.
9. Dilute the high level samples with the assay buffer

STABILITY AND STORAGE

- Storage: Store all of the components at 2-8°C.
- Shelf Life: The kit is stable under the conditions for 12 months from the date of manufacturing. The expiry date is stated on the label of the kit.
- Package: For 96 tests per one kit including standards.

REFERENCES

1. Kojima, M. et al.: Nature, 402: 656, 1999
2. Inui, A.: Nat. Rev. Neurosci., 2: 551, 2001
3. Date, Y. et al. : Endocrinology, 141 : 4255, 2000
4. Gualillo, O. et al. : Endocrinology, 142: 788, 2001
5. Matsumoto, M. et al.: Biochem. Biophys. Res. Commun., 287: 142, 2001
6. Hosoda, H. et al.: Biochem. Biophys. Res. Commun., 279: 909, 2000



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.