RayBio[®] Human/Mouse/Rat Urocortin Enzyme Immunoassay Kit

Catalog #: EIA-URO, EIAM-URO, EIAR-URO

User Manual Last revised August 7, 2017

Caution: Extraordinarily useful information enclosed



ISO 13485 Certified

3607 Parkway Lane, Suite 100 Norcross, GA 30092 Tel: 1-888-494-8555 (Toll Free) or 770-729-2992, Fax:770-206-2393 Web: www.RayBiotech.com, Email: info@raybiotech.com

Table of Contents

Sectio	n	Page #			
1.	Introduction				
II.	General Description				
III.	How It Works				
IV.	Storage				
V.	Reagents				
VI.	Additional Materials Required				
VII.	Reagent Preparation A. Preparation of Plate and Anti-Urocortin Antibody B. Preparation of Biotinylated Peptide (Item F) C. Preparation of Standards D. Preparation of Positive Control E. Preparation of Samples F. Preparation of Wash Buffer and HRP-Strep				
VIII.	Assay Procedure				
IX.	Assay Procedure Summary	11			
Х.	Calculation of Results A. Typical Data B. Sensitivity C. Standard Curve Range D. Reproducibility E. Assay Diagram	12 12 12 12 12 12 13			
XI.	Specificity				
XII.	Select Publications				
XIII.	Troubleshooting Guide				

Please read the entire manual carefully before starting your experiment

I. Introduction

Urocortin is a member of the sauvagine/corticotropin-releasing factor/urotensin I family. It is an endogenous ligand for corticotropin-releasing factor (CRF) receptors. Urocortin, peptide of 40 amino acids, acts in vitro to stimulate the secretion of adrenocorticotropic hormone (ACTH). It binds with high affinity to CRF receptor types 1, 2-alpha, and 2-beta. Urocortin plays a role in the establishment of normal hearing thresholds. It reduces food intake and regulates ghrelin levels in gastric body and plasma.

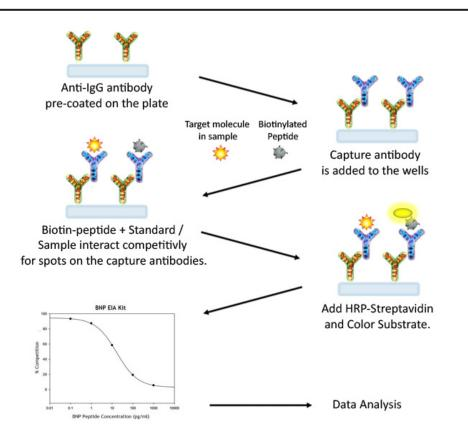
Urocortin is a potent anorexigenic peptide that induces fed-like motor activity when administered centrally or peripherally in fasted animals. Urocortin is also a potent and long-lasting hypotensive agent and increases coronary blood flow.

II. General Description

The RayBio[®] Urocortin Enzyme Immunoassay (EIA) Kit is an in vitro quantitative assay for detecting Urocortin peptide based on the competitive enzyme immunoassay principle.

In this assay, a biotinylated Urocortin peptide is spiked into the samples and standards. The samples and standards are then added to the plate, where the biotinylated Urocortin peptide competes with endogenous (unlabeled) Urocortin for binding to the anti-Urocortin antibody. After a wash step, any bound biotinylated Urocortin then interacts with horseradish peroxidase (HRP)-streptavidin, which catalyzes a color development reaction. The intensity of the colorimetric signal is directly proportional to the amount of captured biotinylated Urocortin peptide and inversely proportional to the amount of endogenous Urocortin in the standard or samples. A standard curve of known concentration of Urocortin peptide can be established and the concentration of Urocortin peptide in the samples can be calculated accordingly.

III. How It Works



IV. Storage

The entire kit may be stored at -20°C to -80°C for up to 6 months from the date of shipment. For extended storage, it is recommended to store at -80°C. **Avoid repeated freeze-thaw cycles.** For prepared reagent storage, see table below.

V. Reagents

Component	Size / Description	Storage / Stability After Preparation		
EIA Microplate (Item A)	96 wells (12 strips x 8 wells) coated with secondary antibody.	1 month at 4°C*		
Wash Buffer Concentrate (20X) (Item B)	25 ml of 20X concentrated solution.	1 month at 4°C		
Standard Urocortin Peptide (Item C)	2 vials of Urocortin Peptide. 1 vial is enough to run each standard in duplicate.	The first standard: 2-3 days at 4°C Additional dilutions: Do not store		
Anti-Urocortin Polyclonal Antibody (Item N)	2 vials of anti-Urocortin.	1 month at 4°C		
5X Assay Diluent B (Item E)	15 ml of 5X concentrated buffer. Diluent for both standards and samples including serum, plasma, cell culture media or other sample types.	1 month at 4°C		
Biotinylated Urocortin Peptide (Item F)	2 vials of Biotinylated Urocortin Peptide, 1 vial is enough to assay the whole plate.	2-3 days at 4°C		
HRP-Streptavidin Concentrate (Item G)	600 μl 100X concentrated HRP-conjugated streptavidin.	Do not store and reuse		
Positive Control (Item M)	1 vial of Positive Control.	2-3 days at 4°C		
TMB One-Step Substrate Reagent (Item H)	12 ml of 3,3,5,5'-tetramethylbenzidine (TMB) in buffer solution.	N/A		
Stop Solution (Item I)	8 ml of 0.2 M sulfuric acid.	N/A		

*Return unused wells to the pouch containing desiccant pack, reseal along entire edge.

VI. Additional Materials Required

- 1. Microplate reader capable of measuring absorbance at 450 nm
- 2. Precision pipettes to deliver 2 µl to 1 ml volumes
- 3. Adjustable 1-25 ml pipettes for reagent preparation
- 4. 100 ml and 1 liter graduated cylinders
- 5. Absorbent paper
- 6. Distilled or deionized water
- 7. SigmaPlot software (or other software which can perform four-parameter logistic regression models)
- 8. Tubes to prepare standard or sample dilutions
- 9. Orbital shaker
- 10. Aluminum foil
- 11. Plastic wrap

VII. Reagent Preparation

Keep kit reagents on ice during reagent preparation steps.

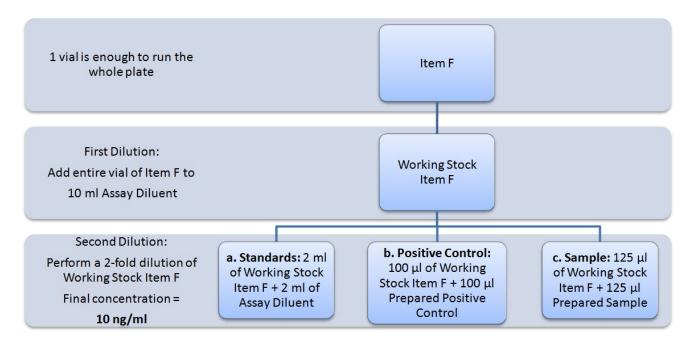
A. Preparation of Plate and Anti-Urocortin Antibody

- 1. Equilibrate plate to room temperature before opening the sealed pouch.
- 2. Label removable 8-well strips as appropriate for your experiment.
- 3. 5X Assay Diluent B (Item E) should be diluted 5-fold with deionized or distilled water.
- Briefly centrifuge the anti-Urocortin antibody vial (Item N). Then add 50 µl of 1X Assay Diluent B to the vial to prepare the antibody concentrate. Pipette up and down to mix gently.
- The antibody concentrate should then be diluted 100-fold with 1X Assay Diluent B. This is your anti-Urocortin antibody working solution, which will be used in step 2 of Assay Procedure (Section VIII).

Note: The following steps may be done during the antibody incubation procedure (step 2 of Assay Procedure)

B. Preparation of Biotinylated Urocortin (Item F)

- 6. Briefly centrifuge the vial of Biotinylated Urocortin (Item F) before use.
- 7. See the image below for proper preparation of Item F. Transfer the entire contents of the Item F vial into a tube containing 10 ml of 1X Assay Diluent B. This is your Working Stock of Item F. Pipette up and down to mix gently. *The final concentration of biotinylated Urocortin will be* **20 ng/ml**.
 - a. Second Dilution of Item F for Standards: Add 2 ml of Working Stock Item F to 2 ml of 1X Assay Diluent B. The final concentration of biotinylated Urocortin will be 10 ng/ml.
 - b. Second Dilution of Item F for Positive Control: Add 100 µl of Working Stock Item F to 100 µl of the prepared Positive Control (Item M). (See section D for Positive Control preparation) The final concentration of biotinylated Urocortin will be **10 ng/ml**.
 - c. Second Dilution of Item F for samples: Add 125 μl of Working Stock Item F to 125 μl of prepared sample (see section E for sample preparation). This is a 2-fold dilution of your sample. The final concentration of biotinylated Urocortin will be **10 ng/ml**.

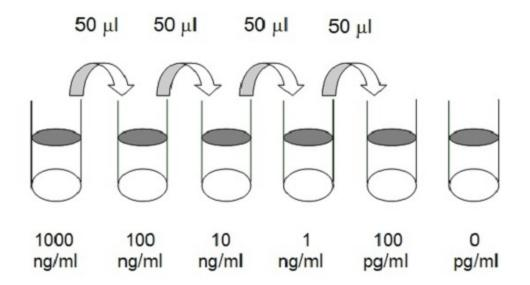


C. Preparation of Standards

Label 6 microtubes with the following concentrations: 1,000 ng/ml, 100 ng/ml, 10ng/ml, 1 ng/ml, 100 pg/ml and 0 pg/ml. Pipette 450 µl of biotinylated Urocortin Item F working solution (prepared in step 7a) into each tube, except the 1,000 ng/ml (leave this one empty).

It is very important to make sure the concentration of biotinylated Urocortin is 10 ng/ml in all standards.

- Briefly centrifuge the vial of Urocortin Standard (Item C). Pipette 8 µl of Item C and 792 µl of 10 ng/ml biotinylated Urocortin working solution (prepared in step 7a) into the tube labeled 1000 ng/ml. Mix thoroughly. This solution serves as the first standard (1,000 ng/ml Urocortin standard, 10 ng/ml biotinylated Urocortin).
- 10. To make the 100 ng/ml standard, pipette 50 μl of the 1000 ng/ml Urocortin standard into the tube labeled 100 ng/ml. Mix thoroughly.
- 11. Repeat this step with each successive concentration, preparing a dilution series as shown in the illustration below. Each time, use 450 µl of biotinylated Urocortin and 50 µl of the prior concentration until the 100 pg/ml is reached. Mix each tube thoroughly before the next transfer.



D. Positive Control Preparation

- 12. Briefly centrifuge the Positive Control vial (Item M).
- 13. Refer to step 7b. This is a 2-fold dilution of the Positive Control. The final concentration of biotinylated Urocortin should still be 10 ng/ml.

The Positive Control is a mouse serum sample that serves as a system control to verify that the kit components are working. The resulting OD will not be used in any calculations; if no positive competition is observed please contact RayBiotech Technical Support. The Positive Control may be diluted further if desired, but be sure the final concentration of biotinylated Urocortin is 10 ng/ml.

E. Sample Preparation

If you wish to perform a 2-fold dilution of your sample, proceed to step 7c. If you wish to perform a higher dilution of your sample, dilute your sample with 1X Assay Diluent B before performing step 7c.

EXAMPLE (to make a 4-fold dilution of sample):

- a. Dilute sample 2-fold (62.5 µl of sample + 62.5 µl of 1X Assay Diluent B.).
- b. Perform step 7c (125 μl of working solution Item F + 125 μl of sample prepared above).

The total volume is 250 μ l, enough for duplicate wells on the microplate. It is very important to make sure the final concentration of the biotinylated Urocortin is **10 ng/ml**.

Note: Optimal sample dilution factors should be determined empirically, however you may contact technical support (888-494-8555;

techsupport@raybiotech.com) to obtain recommended dilution factors for serum.

F. Preparation of Wash Buffer and HRP

- 15. If Item B (20X Wash Concentrate) contains visible crystals, warm to room temperature and mix gently until dissolved.
- 16. Dilute 20 ml of Wash Buffer Concentrate into deionized or distilled water to yield 400 ml of 1X Wash Buffer.
- 17. Briefly centrifuge the HRP-Streptavidin vial (Item G) before use.

18. Dilute the HRP-Streptavidin concentrate 100-fold with 1X Assay Diluent B. **VIII. Assay Procedure**

- 1. Keep kit reagents on ice during reagent preparation steps. It is recommended that all standards and samples be run at least in duplicate.
- Add 100 µl of Anti-Urocortin Antibody (Item N) (See Reagent Preparation step 5) to each well. Incubate for 1.5 hours at room temperature with gentle shaking (1-2 cycle/sec). You may also incubate overnight at 4°C.
- 3. Discard the solution and wash wells 4 times with 1X Wash Solution Buffer (200-300 µl each). Washing may be done with a multichannel pipette or an automated plate washer. Complete removal of liquid at each step is essential to good assay performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
- 4. Add 100 µl of each standard (see Reagent Preparation Section C), Positive Control (see Reagent Preparation Section D) and sample (see Reagent Preparation Section E) to appropriate wells. Be sure to include a blank well (Assay Diluent only). Cover wells and incubate for 2.5 hours at room temperature with gentle shaking (1-2 cycles/sec) overnight or at 4°C.
- 5. Discard the solution and wash 4 times as directed in Step 3.
- Add 100 µl of prepared HRP-Streptavidin solution (see Reagent Preparation step 18) to each well. Incubate for 45 minutes at room temperature with gentle shaking. It is recommended that incubation time should not be shorter or longer than 45 minutes.

- 7. Discard the solution and wash 4 times as directed in Step 3.
- Add 100 µl of TMB One-Step Substrate Reagent (Item H) to each well. Incubate for 30 minutes at room temperature in the dark with gentle shaking (1-2 cycles/sec).
- 9. Add 50 µl of Stop Solution (Item I) to each well. Read at 450 nm immediately.

IX. Assay Procedure Summary

- 1. Prepare all reagents, samples and standards as instructed.
- 2. Add 100 µl anti-Urocortin to each well. Incubate 1.5 hours at room temperature or overnight at 4°C.
- 3. Add 100 µl standard or sample to each well. Incubate 2.5 hours at room temperature or overnight at 4°C.
- 4. Add 100 μI prepared Streptavidin solution. Incubate 45 minutes at room temperature.
- 5. Add 100 µl TMB One-Step Substrate Reagent to each well. Incubate 30 minutes at room temperature.
- 6. Add 50 µl Stop Solution to each well. Read at 450 nm immediately.

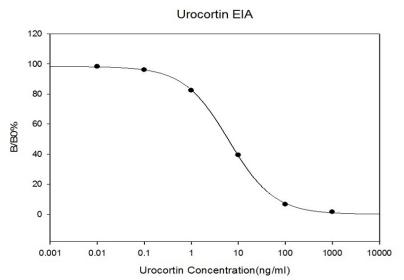
X. Calculation of Results

Calculate the mean absorbance for each set of duplicate stands, controls, and samples and subtract the blank optical density. Plot the standard curve using SigmaPlot software (or other software which can perform four-parameter logistic regression models), with standard concentration on the x-axis and percentage of absorbance (see calculation below) on the y-axis. Draw the best-fit curve through the standard points.

Percentage absorbance = $(B-blank OD)/(B_0-blank OD)$ where B = OD of sample or standard and B₀ = OD of zero standard (total binding)

A. Typical Data

These standard curves are for demonstration only. A standard curve must be run with each assay.



B. Sensitivity

The minimum detectable concentrations of Urocortin is 1.2 ng/ml.

C. Standard Curve Range

0.1-1,000 ng/ml

D. Reproducibility

Intra-Assay: CV<10% Inter-Assay: CV<15%

E. Assay Diagram

Recommended Plate Layout:

Blank	Blank	SA1	SA1	SA9	SA9	SA17	SA17	SA25	SA25	SA33	SA33
Total Binding	Total Binding	SA2	SA2	SA10	SA10	SA18	SA18	SA26	SA26	SA34	SA34
Standard1	Standard1	SA3	SA3	SA11	SA11	SA19	SA19	SA27	SA27	SA35	SA35
Standard2	Standard2	SA4	SA4	SA12	SA12	SA20	SA20	SA28	SA28	SA36	SA36
Standard3	Standard3	SA5	SA5	SA13	SA13	SA21	SA21	SA29	SA29	SA37	SA37
Standard4	Standard4	SA6	SA6	SA14	SA14	SA22	SA22	SA30	SA30	SA38	SA38
Standard5	Standard5	SA7	SA7	SA15	SA15	SA23	SA23	SA31	SA31	SA39	SA39
Pos Control	Pos Control	SA8	SA8	SA16	SA16	SA24	SA24	SA32	SA32	SA40	SA40

Key:

Blank = Buffer Only

Total Binding = Biotin- Urocortin only

Standard 1 = 1000 ng/ml

Standard 2 = 100 ng/ml

Standard 3 = 10 ng/ml

Standard 4 = 1 ng/ml

Standard 5 = 100 pg/ml

Pos Control = Biotin with Item M

XI. Specificity

This EIA kit is designed to detect human, mouse, and rat active Urocortin.

XIV. Select EIA Publications

- Plum L, Lin HV, Dutia R, Tanaka J, Aizawa KS, et al. The Obesity Susceptibility Gene Carboxypeptidase E Links FoxO1 Signaling in Hypothalamic Pro-opiomelanocortin Neurons with Regulation of Food Intake. Nature Med. 2009;15(10):1195-1201. (Ghrelin EIA, EIA-GHR-1)
- Hug C, Lodish HF. Visfatin: a new adipokine. Science. 2005; 307(5708):366-7.
- 3. Kim MK. Crystal structure of visfatin/pre-B cell colony-enhancing factor 1/nicotinamide phosphoribosyltransferase, free and in complex with the anti-cancer agent FK-866. J Mol Biol. 2006; 362(1):66-77.
- 4. Revollo, J.R., et al. The NAD biosynthesis pathway mediated by nicotinamide phosphoribosyltransferase regulates Sir2 activity in mammalian cells. J. Biol. Chem. 2004; 279: 50754-50763.
- 5. Oh-I S, Shimizu H, Satoh T, et al. Identification of nesfatin-1 as a satiety molecule in the hypothalamus. Nature 2006; 443 (7112): 709-12.
- 6. Zhang J, Ren P, Avsian-Kretchmer O, Luo C, Rauch R, Klein C, Hsueh A. Obestatin, a peptide encoded by the ghrelin gene, opposes ghrelin's effects on food intake. Science 2005; 310 (5750): 996-9.
- Cummings D, Weigle D, Frayo R, Breen P, Ma M, Dellinger E, Purnell J. Plasma ghrelin levels after diet-induced weight loss or gastric bypass surgery. N Engl J Med 2002; 346 (21): 1623-30.
- Tschop M, Smiley DL, Heiman ML. Ghrelin induces adiposity in rodents. Nature 2002; 407 (6806): 908-913.9. Kojima M, Hosoda H, Date Y, Nakazato M, Matsuo H, Kangawa K. Ghrelin is a growth-hormone-releasing acylated peptide from stomach. Nature 1999; 402 (6762): 656-60.

XIII. Troubleshooting Guide

Problem	Cause	Solution			
Poor standard curve	 Inaccurate pipetting Improper standard dilution 	 Check pipettes Briefly centrifuge Item C and dissolve the powder thoroughly by gently mixing 			
Low signal	 Improper preparation of standard and/or biotinylated antibody Too brief incubation times Inadequate reagent volumes or improper dilution 	 Briefly spin down vials before opening. Dissolve the powder thoroughly. Ensure sufficient incubation time; assay procedure step 2 may be done overnight Check pipettes and ensure correct preparation 			
Large CV	 Inaccurate pipetting Air bubbles in wells 	 Check pipettes Remove bubbles in wells 			
High background	 Plate is insufficiently washed Contaminated wash buffer 	 Review the manual for proper wash. If using a plate washer, ensure that all ports are unobstructed. Make fresh wash buffer 			
Low sensitivity	 Improper storage of the ELISA kit Stop solution 	 Follow storage recomendations in sections IV and V. Keep substrate solution protected from light. Add stop solution to each well before reading plate 			

Over 3,000 ELISA kits available, visit www.RayBiotech.com/ELISA-Kits.html for details.

This product is for research use only.



©2018 RayBiotech, Inc