

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

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

User Manual
Last revised September 6, 2023

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

Section		Page #
I.	Introduction	3
II.	General Description	4
III.	How It Works	4
IV.	Storage	5
V.	Reagents	5
VI.	Additional Materials Required	6
VII.	Reagent Preparation	6
	A. Preparation of Plate and Anti-PYY Antibody	6
	B. Preparation of Biotinylated Peptide (Item F)	7
	C. Preparation of Standards	8
	D. Preparation of Positive Control	9
	E. Preparation of Samples	9
	F. Preparation of Wash Buffer and HRP-Strep	10
VIII.	Assay Procedure	10
IX.	Assay Procedure Summary	11
X.	Calculation of Results	12
	A. Typical Data	12
	B. Sensitivity	12
	C. Standard Curve Range	12
	D. Reproducibility	12
	E. Assay Diagram	13
XI.	Specificity	14
XII.	Select Publications	14
XIII.	Troubleshooting Guide	15

Please read the entire manual carefully before starting your experiment

I. Introduction

Peptide YY is a 36 amino acid peptide released by cells in the ileum and colon in response to feeding. It is also known as PYY, Peptide Tyrosine Tyrosine, or Pancreatic Peptide YY3-36.

There are two major forms of Peptide YY: PYY1-36 and PYY3-36 which is the most common form of circulating PYY. Peptide YY3-36 (PYY) is a linear polypeptide consisting of 36 amino acids with structural homology to NPY and pancreatic polypeptide. Circulating PYY concentration increases postprandially and decreases by fasting.

PYY exerts its action through NPY receptors, inhibits gastric motility and increases water and electrolyte absorption in the colon. PYY may also suppress pancreatic secretion. It is secreted by the neuroendocrine cells in the ileum and colon in response to a meal, and has been shown to reduce appetite. PYY works by slowing the gastric emptying; hence, it increases efficiency of digestion and nutrient absorption after meal.

PYY has been shown to play an important role in obesity. Animal studies have shown that acute peripheral administration of PYY3-36 inhibits feeding of rodents and primates. Studies on Y2R-knockout mice have revealed that there is no anorectic effect on Y2R-knockout mice (Y2R is the receptor for PYY). These findings indicate that PYY3-36 has anorectic effect which is suggested to be mediated by Y2R. Studies on PYY-knockout mice have shown that they have higher fat mass and lower glucose tolerance when compared to control mice, indicating that PYY also plays very important role in energy homeostasis by balancing the food intake.

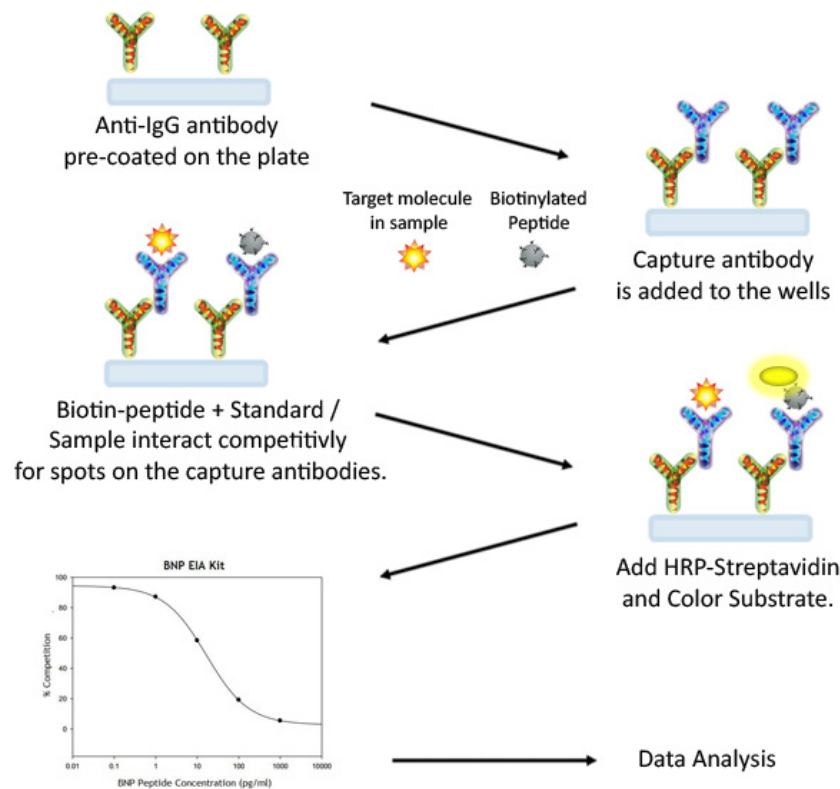
Studies have also shown that obese people secrete less PYY than non-obese people. The anorectic effect of PYY represents a possible anti-obesity therapy in the future.

II. General Description

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

In this assay, a biotinylated PYY peptide is spiked into the samples and standards. The samples and standards are then added to the plate, where the biotinylated PYY peptide competes with endogenous (unlabeled) PYY for binding to the anti-PYY antibody. After a wash step, any bound biotinylated PYY 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 PYY peptide and inversely proportional to the amount of endogenous PYY in the standard or samples. A standard curve of known concentration of PYY peptide can be established and the concentration of PYY 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 PYY Peptide (Item C)	2 vials of Lyophilized PYY Peptide. 1 vial is enough to run each standard in duplicate.	Do not store and reuse
Anti-PYY Polyclonal Antibody (Item N)	2 vials of Lyophilized anti-PYY.	Do not store and reuse
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 PYY Peptide (Item F)	2 vials of Lyophilized Biotinylated PYY Peptide, 1 vial is enough to assay the whole plate.	Do not store and reuse
HRP-Streptavidin Concentrate (Item G)	600 µl 100x concentrated HRP-conjugated streptavidin.	Do not store and reuse
Positive Control (Item M)	1 vial of Lyophilized Positive Control.	Do not store and reuse
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 μ l 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-PYY 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.
4. Briefly centrifuge the anti-PYY antibody vial (Item N) and reconstitute with 55 μ l of 1X Assay Diluent B to prepare the antibody concentrate. Pipette up and down to mix gently.
5. The antibody concentrate should then be diluted 100-fold with 1X Assay Diluent B. This is your anti-PYY 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)

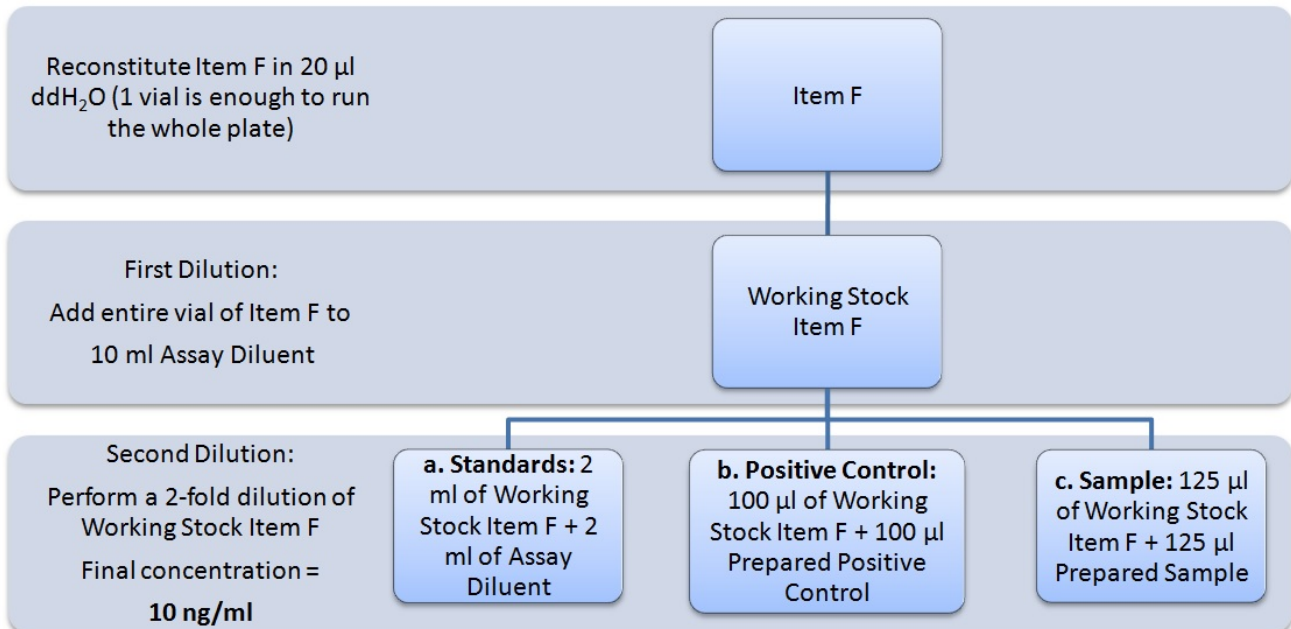
B. Preparation of Biotinylated PYY (Item F)

6. Briefly centrifuge the vial of Biotinylated PYY (Item F) and reconstitute with 20 μl of ddH₂O 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 PYY will be 20 $\mu\text{g/ml}$.*
 - 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 PYY will be **10 $\mu\text{g/ml}$** .

 - 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 PYY will be **10 $\mu\text{g/ml}$** .

 - 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 PYY will be **10 $\mu\text{g/ml}$** .

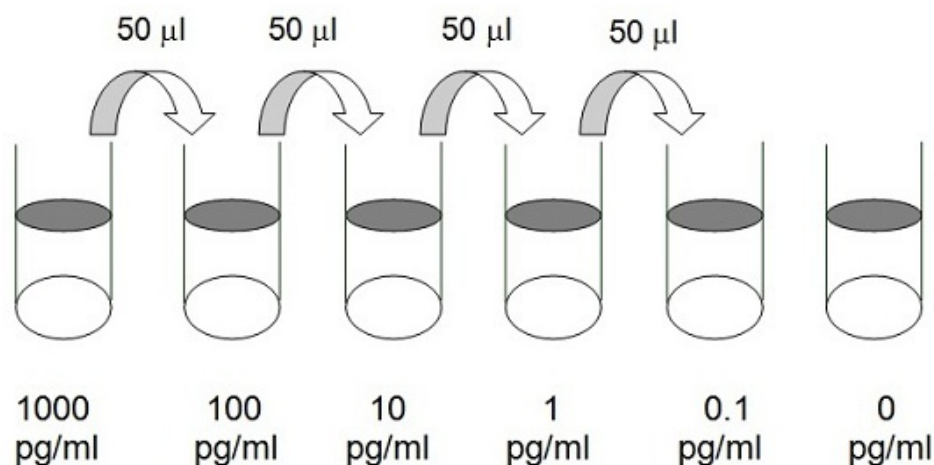


C. Preparation of Standards

- Label 6 microtubes with the following concentrations: 1,000 pg/ml, 100 pg/ml, 10 pg/ml, 1 pg/ml, 0.1 pg/ml and 0 pg/ml. Pipette 450 μ l of biotinylated PYY Item F working solution (prepared in step 7a) into each tube, except the 1,000 pg/ml (leave this one empty).

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

- Briefly centrifuge the vial of PYY Standard (Item C). Reconstitute with 10 μ l of ddH₂O and briefly vortex if desired. Pipette 8 μ l of Item C and 792 μ l of 10 pg/ml biotinylated PYY working solution (prepared in step 7a) into the tube labeled 1000 pg/ml. Mix thoroughly. This solution serves as the first standard (1000 pg/ml PYY standard, 10 pg/ml biotinylated PYY).
- To make the 100 pg/ml standard, pipette 50 μ l of the 1000 pg/ml PYY standard into the tube labeled 100 pg/ml. Mix thoroughly.
- Repeat this step with each successive concentration, preparing a dilution series as shown in the illustration below. Each time, use 450 μ l of biotinylated PYY and 50 μ l of the prior concentration until the 0.1 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) and reconstitute with 100 μl of ddH₂O.

13. Refer to step 7b. This is a 2-fold dilution of the Positive Control. The final concentration of biotinylated PYY should still be 10 pg/ml.

The Positive Control is a mouse serum sample 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 PYY is 10 pg/ml.

E. Sample Preparation

14. 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 μl , enough for duplicate wells on the microplate. It is very important to make sure the final concentration of the biotinylated PYY is **10 pg/ml**.

Note: Optimal sample dilution factors should be determined empirically, however you may reference below for recommended dilution factors for serum: Human=2x Mouse=2x Rat=2x .

If you have any questions regarding the recommended dilutions you may contact technical support at 888-494-8555 or techsupport@raybiotech.com.

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.
2. Add 100 μ l of Anti-PYY 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 μ l 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 μ l 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.

6. Add 100 μ l 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.
8. Add 100 μ l 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 μ l 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 μ l anti-PYY to each well. Incubate 1.5 hours at room temperature or overnight at 4°C.
3. Add 100 μ l standard or sample to each well. Incubate 2.5 hours at room temperature or overnight at 4°C.
4. Add 100 μ l prepared Streptavidin solution. Incubate 45 minutes at room temperature.
5. Add 100 μ l TMB One-Step Substrate Reagent to each well. Incubate 30 minutes at room temperature.
6. Add 50 μ l Stop Solution to each well. Read at 450 nm immediately.

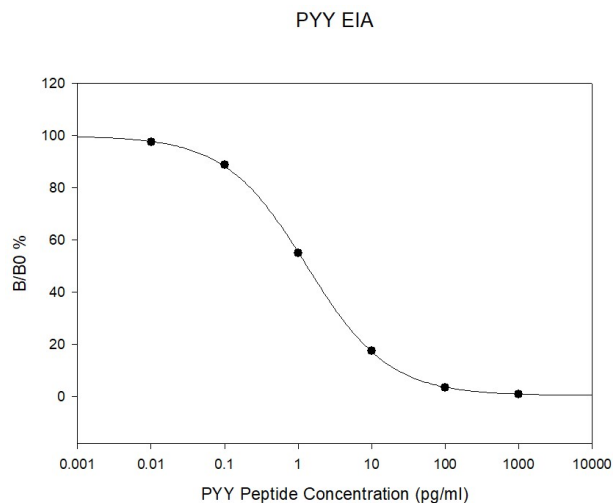
X. Calculation of Results

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 = $(B - \text{blank OD}) / (B_0 - \text{blank OD})$ where
B = OD of sample or standard and
 B_0 = 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 PYY is 5.6 pg/ml.

C. Standard Curve Range

0.1-1,000 pg/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-PYY only

Standard 1 = 1000 pg/ml

Standard 2 = 100 pg/ml

Standard 3 = 10 pg/ml

Standard 4 = 1 pg/ml

Standard 5 = 0.1 pg/ml

Pos Control = Biotin with Item M

XI. Specificity

Cross Reactivity: This EIA kit shows no cross-reactivity with any of the cytokines tested: Ghrelin, Nesfatin, Angiotensin II, NPY and APC.

This kit detects the 1-36 form PYY.

XIV. Publications Citing This Product

1. Bar F., et al. Carboxypeptidase E Modulates Intestinal Immune Homeostasis and Protects against Experimental Colitis in Mice. PLOS One. Published: July 22, 2014. DOI: 10.1371/journal.pone.0102347
Species: Mouse
Sample Type: Conditioned Media
2. Hand K., et al. Hormone profiling in a novel enteroendocrine cell line pGIP/neo: STC-1. METABOLISM CLINICAL AND EXPERIMENTAL 61 (2012) 1683–1686
Species: Mouse
Sample Type: Conditioned Media
3. Lin N, Cai DL, Jin D, Chen Y, Shi JJ. Ginseng Panaxoside Rb1 Reduces Body Weight in Diet-Induced Obese Mice. Cell Biochem Biophys. Cell Biochem Biophys (2014) 68:189–194 DOI 10.1007/s12013-013-9688-3
Species: Mouse
Sample Type: Serum

XIII. Troubleshooting Guide

Problem	Cause	Solution
Poor standard curve	<ul style="list-style-type: none"> ○ Inaccurate pipetting ○ Improper standard dilution 	<ul style="list-style-type: none"> ○ Check pipettes ○ Briefly centrifuge Item C and dissolve the powder thoroughly by gently mixing
Low signal	<ul style="list-style-type: none"> ○ Improper preparation of standard and/or biotinylated antibody ○ Too brief incubation times ○ Inadequate reagent volumes or improper dilution 	<ul style="list-style-type: none"> ○ 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	<ul style="list-style-type: none"> ○ Inaccurate pipetting ○ Air bubbles in wells 	<ul style="list-style-type: none"> ○ Check pipettes ○ Remove bubbles in wells
High background	<ul style="list-style-type: none"> ○ Plate is insufficiently washed ○ Contaminated wash buffer 	<ul style="list-style-type: none"> ○ Review the manual for proper wash. If using a plate washer, ensure that all ports are unobstructed. ○ Make fresh wash buffer
Low sensitivity	<ul style="list-style-type: none"> ○ Improper storage of the ELISA kit ○ Stop solution 	<ul style="list-style-type: none"> ○ Follow storage recommendations in sections IV and V. Keep substrate solution protected from light. ○ Add stop solution to each well before reading plate

RayBio[®] ELISA Kits

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