RayBio[®] Rat BDNF ELISA Kit

Catalog #: ELR-BDNF

User Manual Last revised June 14, 2021

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



RayBio[®] Rat BDNF ELISA Kit Protocol

Table of Contents

Section				
Ι.	Introduction			
II.	Storage	4		
III.	Reagents	4		
IV.	Additional Materials Required			
V.	Reagent Preparation	5		
VI.	Assay Procedure			
VII.	Assay Procedure Summary			
VIII.	Calculation of Results A. Typical Data B. Sensitivity C. Spiking & Recovery D. Linearity E. Reproducibility	8 8 9 9 10		
IX.	Specificity	10		
Χ.	Troubleshooting Guide	11		

Please read the entire manual carefully before starting your experiment

I. INTRODUCTION

The RayBio[®] Rat BDNF ELISA kit is an in-vitro enzyme-linked immunosorbent assay for the quantitative measurement of Rat BDNF in serum, plasma and cell culture supernatants. This assay employs an antibody specific for BDNF coated on a 96-well plate. Standards and samples are pipetted into the wells and BDNF present in a sample is bound to the wells by the immobilized antibody. The wells are washed and biotinylated anti-Rat BDNF antibody is added. After washing away unbound biotinylated antibody, HRP-conjugated streptavidin is pipetted to the wells. The wells are again washed, a TMB substrate solution is added to the wells and color develops in proportion to the amount of BDNF bound. The Stop Solution changes the color from blue to yellow, and the intensity of the color is measured at 450 nm.

Need your results faster? Try RayBiotech's *SpeedELISA* platform for quantitative detection in just three hours.

https://www.raybiotech.com/speedelisa-kits/

Short on sample, or need higher sensitivity? Check out the IQELISA[™] Immuno-PCR platform. https://www.raybiotech.com/immuno-pcr/

II. STORAGE

The entire kit may be stored at -20°C for up to 1 year from the date of shipment. Avoid repeated freeze-thaw cycles. The kit may be stored at 4°C for up to 6 months. For extended storage, it is recommended to store at -80°C. For prepared reagent storage, see table below.

III. REAGENTS

Component	Size / Description	Storage / Stability After Preparation	
BDNF Microplate (Item A)	96 wells (12 strips x 8 wells) coated with anti-Rat BDNF.	1 month at 4°C*	
Wash Buffer Concentrate (20X) (Item B)	25 ml of 20X concentrated solution.	1 month at 4°C	
Standard Protein (Item C)	2 vials of Rat BDNF. 1 vial is enough to run each standard in duplicate.	1 week at -80°C	
Detection Antibody BDNF (Item F)	2 vials of biotinylated anti-Rat BDNF. Each vial is enough to assay half the microplate.	5 days at 4°C	
HRP-Streptavidin Concentrate (Item G)	200 µl 200X concentrated HRP-conjugated streptavidin.	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	
Assay Diluent C (Item L)	30 ml of diluent buffer.	N/A	
Assay Diluent B (Item E)	15 ml of 5X concentrated buffer.	1 month at 4°C	

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

IV. ADDITIONAL MATERIALS REQUIRED

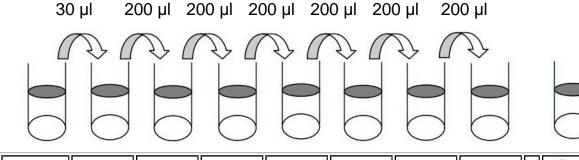
- 1. Microplate reader capable of measuring absorbance at 450 nm.
- 2. Precision pipettes to deliver 2 µl to 1 ml volumes.
- 3. Adjustable 1-25 ml pipettes for reagent preparation.
- 4. 100 ml and 1 liter graduated cylinders.
- 5. Absorbent paper.
- 6. Distilled or deionized water.
- 7. Log-log graph paper or computer and software for ELISA data analysis.
- 8. Tubes to prepare standard or sample dilutions.

V. REAGENT PREPARATION

- 1. Bring all reagents and samples to room temperature (18 25°C) before use.
- Assay Diluent B (Item E) should be diluted 5-fold with deionized or distilled water before use.
- Sample dilution: Assay Diluent C (Item L) should be used for dilution of serum, plasma, and cell culture supernatant samples. The suggested dilution for normal serum/plasma is 10 fold.

Note: Levels of BDNF may vary between different samples. Optimal dilution factors for each sample must be determined by the investigator.

4. Preparation of standard: Briefly spin a vial of Item C. Add 400 µl of Assay Diluent C into Item C vial to prepare a 50 ng/ml standard. Dissolve the powder thoroughly by a gentle mix. Add 30 µl BDNF standard from Item C into a tube with 470 µl Assay Diluent C to prepare a 3,000 pg/ml standard solution. Pipette 300 µl of Assay Diluent C into each tube. Use the 3,000 pg/ml standard solution to produce a dilution series (shown below). Mix each tube thoroughly before the next transfer. Assay Diluent C serves as the zero standard (0 pg/ml).



		Std1	Std2	Std3	Std4	Std5	Std6	Std7	Zero Standard
Diluent volume	Item C + 400 μΙ	470 µl	300 µl	300 µl	300 µl	300 µl	300 µl	300 µl	300 µl
Conc.	50 ng/ml	3000 pg/ml	1200 pg/ml	480 pg/ml	192 pg/ml	76.8 pg/ml	30.72 pg/ml	12.29 pg/ml	0 pg/ml

- 5. If the Wash Concentrate (20X) (Item B) contains visible crystals, warm to room temperature and mix gently until dissolved. Dilute 20 ml of Wash Buffer Concentrate into deionized or distilled water to yield 400 ml of 1X Wash Buffer.
- 6. Briefly spin the Detection Antibody vial (Item F) before use. Add 100 µl of 1X Assay Diluent B (Item E) into the vial to prepare a detection antibody concentrate. Pipette up and down to mix gently (the concentrate can be stored at 4°C for 5 days). The detection antibody concentrate should be diluted 80-fold with 1X Assay Diluent B (Item E) and used in step 5 of Part VI Assay Procedure.
- Briefly spin the HRP-Streptavidin concentrate vial (Item G) and pipette up and down to mix gently before use, as precipitates may form during storage. HRP-Streptavidin concentrate should be diluted 200-fold with 1X Assay Diluent B (Item E).

For example: Briefly spin the vial (Item G) and pipette up and down to mix gently. Add 50 µl of HRP-Streptavidin concentrate into a tube with 10 ml 1X Assay Diluent B to prepare a 200-fold diluted HRP-Streptavidin solution (do not store the diluted solution for next day use). Mix well.

VI. ASSAY PROCEDURE

- 1. Bring all reagents and samples to room temperature (18 25°C) before use. It is recommended that all standards and samples be run at least in duplicate.
- 2. Label removable 8-well strips as appropriate for your experiment.
- 3. Add 100 µl of each standard (see Reagent Preparation step 3) and sample into appropriate wells. Cover wells and incubate for 2.5 hours at room temperature with gentle shaking.
- 4. Discard the solution and wash 4 times with 1X Wash Solution. Wash by filling each well with Wash Buffer (300 μl) using a multi-channel Pipette or autowasher. Complete removal of liquid at each step is essential to good performance. After the last wash, remove any remaining Wash Buffer by aspirating or decanting. Invert the plate and blot it against clean paper towels.
- Add 100 µl of 1X prepared biotinylated antibody (Reagent Preparation step 6) to each well. Incubate for 1 hour at room temperature with gentle shaking.
- 6. Discard the solution. Repeat the wash as in step 4.
- Add 100 µl of prepared Streptavidin solution (see Reagent Preparation step 7) to each well. Incubate for 45 minutes at room temperature with gentle shaking.

- 8. Discard the solution. Repeat the wash as in step 4.
- 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.
- 10. Add 50 µl of Stop Solution (Item I) to each well. Read at 450 nm immediately.

VII. ASSAY PROCEDURE SUMMARY

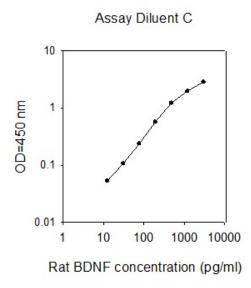
- 1. Prepare all reagents, samples and standards as instructed.
- 2. Add 100 µl standard or sample to each well. Incubate 2.5 hours at room temperature.
- 3. Add 100 µl prepared biotin antibody to each well. Incubate 1 hour at room temperature.
- 4. Add 100 µl prepared Streptavidin solution. Incubate 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.

VIII. CALCULATION OF RESULTS

Calculate the mean absorbance for each set of duplicate standards, controls and samples, and subtract the average zero standard optical density. Plot the standard curve on log-log graph paper or using Sigma plot software, with standard concentration on the x-axis and absorbance on the y-axis. Draw the best-fit straight line through the standard points.

A. TYPICAL DATA

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



B. SENSITIVITY

The minimum detectable dose of Rat BDNF was determined to be 12 pg/ml.

Minimum detectable dose is defined as the analyte concentration resulting in an absorbance that is 2 standard deviations higher than that of the blank (diluent buffer).

C. SPIKING & RECOVERY

Recovery was determined by spiking various levels of Rat BDNF into the sample types listed below. Mean recoveries are as follows:

Sample Type	Average % Recovery	Range (%)
Serum	124.2	102-148
Plasma	116.9	82-149
Cell culture media	113.7	94-131

D. LINEARITY

Sample Type		Serum	Plasma	Cell Culture Media		
1:2	Average % of Expected	110.7	109.1	109.6		
	Range (%)	106-115	94-122	101-118		
1:4	Average % of Expected	106.3	84.62	111.3		
	Range (%)	99-113	72-102	103-119		

E. REPRODUCIBILITY

Intra-Assay CV%: <10%

Inter-Assay CV%: <12%

IX. SPECIFICITY

This ELISA antibody pair detects rat and human BDNF. Other species not determined.

X. 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 3 may be done overnight at 4°C with gentle shaking (note: may increase overall signals including background). Check pipettes and ensure correct preparation 			
Large CV	 Inaccurate pipetting Air bubbles in wells 	Check pipettesRemove 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 	 Store your standard at <-70°C after reconstitution, others at 4°C. Keep substrate solution protected from light. Add stop solution to each well before reading plate 			

RayBio[®] ELISA Kits

Over 4,000 ELISA kits available, visit www.RayBiotech.com/ELISA-Kits for details.

This product is for research use only.



©2020 RayBiotech, Inc