



Ngf (Rat) ELISA Kit

Catalog Number KA0401

96 assays

Version: 53

Intended for research use only

www.abnova.com

Table of Contents

Introduction	3
Intended Use	3
Background	3
Principle of the Assay	3
General Information	4
Materials Supplied	4
Storage Instruction	4
Materials Required but Not Supplied	4
Precautions for Use	5
Assay Protocol	6
Reagent Preparation	6
Sample Preparation	7
Assay Procedure	7
Data Analysis.....	9
Calculation of Results.....	9
Performance Characteristics	10
Resources.....	11
Plate Layout	11

Introduction

Intended Use

For the quantitation of Rat Ngf concentrations in cell culture supernates and serum.

Background

Nerve growth factor (NGF) is a polypeptide involved in the regulation of growth and differentiation of sympathetic and certain sensory neurons. NGF is thought to have a profound effect on the development and maintenance of sympathetic and embryonic sensory neurones. NGF activity isolated from the male mouse submaxillary gland (MSG) consists of three types of subunits, alpha, beta and gamma, which specifically interact to form a 7S, approximately 130,000-molecular weight (Mr) complex. The 7S complex contains two identical 118-amino acid beta-chains, which are solely responsible for the nerve growth-stimulating activity of NGF. NGF, which is expressed by inflammatory cells and effects changes that lead to increased neural responsiveness, could be a pivotal mediator in allergic rhinitis.

Principle of the Assay

The Ngf (Rat) ELISA Kit is a solid phase immunoassay specially designed to measure Rat Ngf with a 96-well strip plate that is pre-coated with antibody specific for Ngf. The detection antibody is a biotinylated antibody specific for Ngf. The capture antibody is monoclonal antibody from mouse, the detection antibody is polyclonal antibody from goat. The kit contains recombinant Rat Ngf with immunogen: Expression system for standard: NSO; Immunogen sequence: S122-G241. The kit is analytically validated with ready to use reagents.

To measure Rat Ngf, add standards and samples to the wells, then add the biotinylated detection antibody. Wash the wells with PBS or TBS buffer, and add Avidin-Biotin-Peroxidase Complex (ABC-HRP). Wash away the unbounded ABC-HRP with PBS or TBS buffer and add TMB. TMB is substrate to HRP and will be catalyzed to produce a blue color product, which changes into yellow after adding acidic stop solution. The density of the yellow product is linearly proportional to Rat Ngf in the sample. Read the density of the yellow product in each well using a plate reader, and benchmark the sample wells' readings against the standard curve to determine the concentration of Rat Ngf in the sample.

General Information

Materials Supplied

List of component

Component	Amount
Anti-Rat Ngf Pre-coated 96-well strip microplate	96 (8x12) wells
Rat Ngf Standard	10 ng/tube x 2
Rat Ngf Biotinylated antibody (100x)	130 µL
Avidin-Biotin-Peroxidase Complex (100x)	130 µL
Sample Diluent	30 mL
Antibody Diluent	12 mL
Avidin-Biotin-Peroxidase Diluent	12 mL
Color Developing Reagent (TMB)	10 mL
Stop Solution	10 mL
Plate Sealers	4 slides

Storage Instruction

Store at 4°C for 6 months, at -20°C for 12 months. Avoid multiple freeze-thaw cycles.

Materials Required but Not Supplied

- ✓ Microplate Reader capable of reading absorbance at 450 nm.
- ✓ Automated plate washer (optional).
- ✓ Pipettes and pipette tips capable of precisely dispensing 0.5 µL through 1 mL volumes of aqueous solutions.
- ✓ Multichannel pipettes are recommended for large amount of samples.
- ✓ Deionized or distilled water.
- ✓ 500 mL graduated cylinders.
- ✓ Test tubes for dilution.
- ✓ Washing buffer

Prepare standard 1X PBS as wash buffer. Wash buffer can be prepared in-house.

Preparation of wash buffer: Add 8.5 g NaCl, 1.4 g Na₂HPO₄ and 0.2 g NaH₂PO₄ to 1000 mL distilled water and adjust pH to 7.2-7.6.

Precautions for Use

This protocol must be read in its entirety before using this product. For research use only. Not for use in diagnostic procedures.

✓ Notice Before Application

Please read the following instructions before starting the experiment.

1. To inspect the validity of experiment operation and the appropriateness of sample dilution proportion, pilot experiment using standards and a small number of samples is recommended.
2. Before using the Kit, spin tubes and bring down all components to the bottom of tubes.
3. Don't let 96-well plate dry, for dry plate will inactivate active components on plate.
4. Don't reuse tips and tubes to avoid cross contamination.
5. Avoid using the reagents from different batches together.

Assay Protocol

Reagent Preparation

- ✓ Bring all reagents to 37°C prior to use. The assay can also be done at room temperature however we recommend doing it at 37°C for best consistency with our QC results. Also the TMB incubation time estimate (15-25 min) is based on 37°C.
 - Biotinylated Anti-Rat Ngf antibody
It is recommended to prepare this reagent immediately prior to use by diluting the Rat Ngf Biotinylated antibody (100x) 1:100 with Antibody Diluent. Prepare 100 µL by adding 1 µL of Biotinylated antibody (100x) to 99 µL of Antibody Diluent for each well. Mix gently and thoroughly and use within 2 hours of generation.
 - Avidin-Biotin-Peroxidase Complex
It is recommended to prepare this reagent immediately prior to use by diluting the Avidin-Biotin-Peroxidase Complex (100x) 1:100 with Avidin-Biotin-Peroxidase Diluent. Prepare 100 µL by adding 1 µL of Avidin-Biotin-Peroxidase Complex (100x) to 99 µL of Avidin-Biotin-Peroxidase Diluent for each well. Mix gently and thoroughly and use within 2 hours of generation.
 - Rat Ngf Standard
It is recommended that the standards be prepared no more than 2 hours prior to performing the experiment. Use one 10 ng of lyophilized Rat Ngf standard for each experiment. Gently spin the vial prior to use. Reconstitute the standard to a stock concentration of 10 ng/mL using 1 mL of sample diluent. Allow the standard to sit for a minimum of 10 minutes with gentle agitation prior to making dilutions.
 - Microplate
The included microplate is coated with capture antibodies and ready-to-use. It does not require additional washing or blocking. The unused well strips should be sealed and stored in the original packaging.
- ✓ Dilution of Rat Ngf Standard
 1. Number tubes 1-8. Final Concentrations to be Tube # 1 –1000 pg/mL, #2 –500 pg/mL, #3 – 250 pg/mL, #4 – 125 pg/mL, #5 – 62.5 pg/mL, #6 – 31.25 pg/mL, #7 – 15.625 pg/mL , #8 –Sample Diluent serves as the zero standard (0 pg/mL).
 2. To generate standard #1, add 100 µL of the reconstituted standard stock solution of 10 ng/mL and 900 µL of sample diluent to tube #1 for a final volume of 1000 µL. Mix thoroughly.
 3. Add 300 µL of sample diluent to tubes # 2-7.
 4. To generate standard #2, add 300 µL of standard #1 from tube #1 to tube #2 for a final volume of 600 µL. Mix thoroughly.
 5. To generate standard #3, add 300 µL of standard #2 from tube #2 to tube #3 for a final volume of 600 µL. Mix thoroughly.
 6. Continue the serial dilution for tube #4-7.

Sample Preparation

✓ Sample Preparation and Storage

These sample collection instructions and storage conditions are intended as a general guideline and the sample stability has not been evaluated.

- Cell culture supernatants: Clear sample of particulates by centrifugation, assay immediately or store samples at -20°C.
- Serum: Use a serum separator tube (SST) and allow serum to clot at room temperature for about four hours. Then, centrifuge for 15 min at approximately 1,000 x g. assay immediately or store samples at -20°C.

✓ Sample Dilution

The target protein concentration should be estimated and appropriate sample dilutions should be selected such that the final protein concentration lies near the middle of the linear dynamic range of the assay.

It is recommended to prepare 150 µL of sample for each replicate to be assayed. The samples should be diluted with sample diluent and mixed gently.

Assay Procedure

It is recommended that all reagents and materials be equilibrated to 37°C/room temperature prior to the experiment (see Reagent Preparation if you have missed this information).

1. Prepare all reagents and working standards as directed previously.
2. Remove excess microplate strips from the plate frame and seal and store them in the original packaging.
3. Add 100 µL of the standard, samples, or control per well. Add 100 µL of the sample diluent buffer into the zero well. At least two replicates of each standard, sample, or control is recommended.
4. Cover with the plate sealer provided and incubate for 120 minutes at RT (or 90 minutes at 37°C).
5. Remove the cover and discard the liquid in the wells into an appropriate waste receptacle. Invert the plate on the benchtop onto a paper towel and tap the plate to gently blot any remaining liquid. It is recommended that the wells are not allowed to completely dry at any time.
6. Add 100 µL of the prepared 1x Biotinylated Anti-Rat Ngf antibody to each well.
7. Cover with plate sealer and incubate for 90 minutes at RT (or 60 minutes at 37°C).
8. Wash the plate 3 times with the 1x wash buffer.
 - a. Discard the liquid in the wells into an appropriate waste receptacle. Then, invert the plate on the benchtop onto a paper towel and tap the plate to gently blot any remaining liquid. It is recommended that the wells are not allowed to completely dry at any time.
 - b. Add 300 µL of the 1x wash buffer to each assay well. (For cleaner background incubate for 60 seconds between each wash).

- c. Repeat steps a-b 2 additional times.
9. Add 100 μL of the prepared 1x Avidin-Biotin-Peroxidase Complex into each well. Cover with the plate sealer provided and incubate for 40 minutes at RT (or 30 minutes at 37°C).
10. Wash the plate 5 times with the 1x wash buffer.
 - a. Discard the liquid in the wells into an appropriate waste receptacle. Then, invert the plate on the benchtop onto a paper towel and tap the plate to gently blot any remaining liquid. It is recommended that the wells are not allowed to completely dry at any time.
 - b. Add 300 μL of the 1x wash buffer to each assay well. (For cleaner background incubate for 60 seconds between each wash).
 - c. Repeat steps a-b 4 additional times.
11. Add 90 μL of Color Developing Reagent to each well. Cover with the plate sealer provided and incubate in the dark for 30 minutes at RT (or 15-25 minutes at 37°C). (The optimal incubation time must be empirically determined. A guideline to look for is blue shading the top four standard wells, while the remaining standards remain clear.).
12. Add 100 μL of Stop Solution to each well. The color should immediately change to yellow.
13. Within 30 minutes of stopping the reaction, the O.D. absorbance should be read with a microplate reader at 450 nm.

Data Analysis

Calculation of Results

Average the duplicate readings for each standard, sample, and control. Subtract the average zero standard O.D. reading.

It is recommended that a standard curve be created using computer software to generate a four parameter logistic (4-PL) curve assay.

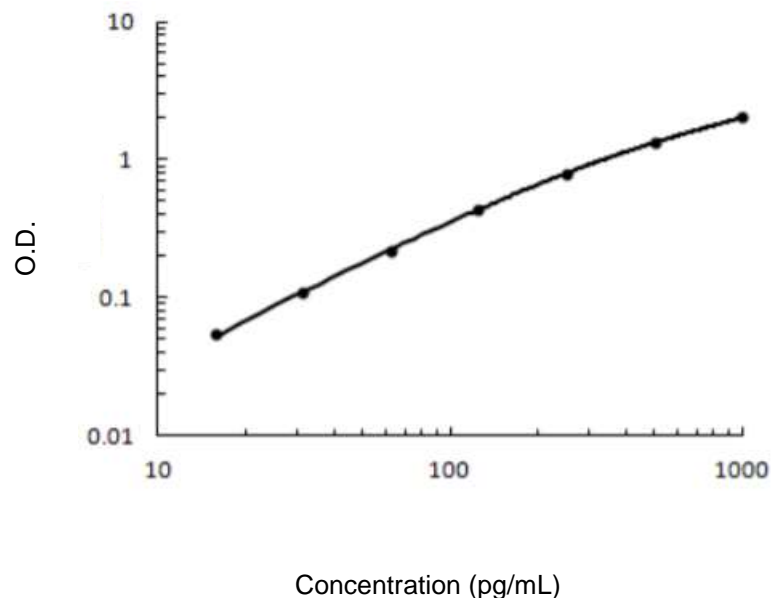
Alternatively, plot the mean absorbance for each standard against the concentration. The measured concentration in the sample can be interpolated by using linear regression of each average relative OD against the standard curve generated using curve fitting software. This will generate an adequate but less precise fit of the data.

For diluted samples, the concentration reading from the standard curve must be multiplied by the dilution factor.

✓ The Ngf (Rat) ELISA Kit Standard Curve Example

Highest O.D. value might be higher or lower than in the example. The experiment result is statistically significant if the highest O.D. value is no less than 1.0.

Concentration (pg/mL)	0	15.6	31.2	62.5	125	250	500	1000
O.D	0.079	0.133	0.190	0.296	0.521	0.876	1.414	2.125



A standard curve is provided for demonstration only. A standard curve should be generated for each set of samples assayed.

Performance Characteristics

- ✓ Detection Range: 15.6 pg/mL to 1000 pg/mL
- ✓ Sensitivity: < 1 pg/mL
 - * The sensitivity or the minimum detectable dose (MDD) is the lower limit of target protein that can be detected by the kit. It is determined by adding two standard deviations to the mean O.D. value of twenty (20) blank wells and calculating the corresponding concentration.
- ✓ Specificity: Natural and recombinant Rat Ngf
- ✓ Cross-reactivity: There is no detectable cross-reactivity with BDNF, GDNF, CNTF, NT3, and NT4.
- ✓ Intra/Inter Assay Variability
 - Intra-Assay Precision (Precision within an assay)
 - Three samples of known concentration were tested on one plate to assess intra-assay precision.
 - Inter-Assay Precision (Precision across assays)
 - Three samples of known concentration were tested in separate assays to assess inter-assay precision.

Sample	Intra-Assay Precision			Inter-Assay Precision		
	1	2	3	1	2	3
n	16	16	16	24	24	24
Mean (pg/mL)	33	191	480	29	201	491
Standard deviation	1.84	8.02	32.64	1.62	11.65	35.84
CV (%)	4.2%	4.2%	6.8%	5.6%	5.8%	7.3%

- ✓ Reproducibility
 - To assay reproducibility, three samples with differing target protein concentrations were assayed using four different lots.

Lots	Lot1 (pg/mL)	Lot2 (pg/mL)	Lot3 (pg/mL)	Lot4 (pg/mL)	Mean (pg/mL)	Standard Deviation	CV (%)
Sample 1	33	32	33	35	33	1.08	3.2%
Sample 2	191	172	185	192	185	7.96	4.3%
Sample 3	480	451	440	436	451	17.21	3.8%

*number of samples for each test n=16.

Resources

Plate Layout

12								
11								
10								
9								
8								
7								
6								
5								
4								
3								
2								
1								
	A	B	C	D	E	F	G	H