

Version 2 Last updated 12 December 2016

ab215415

Human CDNF ELISA kit

For the quantitative determination of human CDNF in cell culture supernatant, serum and plasma

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1. Overview	1
2. Protocol Summary	2
3. Precautions	3
4. Storage and Stability	3
5. Limitations	4
6. Materials Supplied	4
7. Materials Required, Not Supplied	5
8. Technical Hints	6
9. Reagent Preparation	7
10. Sample Preparation	7
11. Plate Preparation	8
12. Assay Procedure	9
13. Calculations	10
14. Typical data	11
15. Typical sample values	12
16. Troubleshooting	14
17. Notes	15

1. Overview

The human CDNF ELISA kit (ab215415) provides a rapid and easy method for the quantitative determination of human CDNF in cell culture supernatant, serum and plasma. The kit includes ready-to-use reagents necessary to analyze up to 88 samples in 2 hours.

The human CDNF test is based on the quantitative sandwich enzyme immunoassay technique. Microplate wells are pre-coated with anti-human CDNF-specific monoclonal capture antibodies. Samples and standards are pipetted into microwells and human CDNF molecules present in the sample are bound by the capture antibodies. After incubation, unbound material is removed by washing the wells. Then, horseradish peroxidase (HRP) conjugated human CDNF-specific monoclonal detection antibodies bind to a different epitope of human CDNF molecules. After washing, the ready to use HRP substrate (TMB) is added to wells. The intensity of the color produced is directly proportional to the amount of human CDNF in the sample. Color development is then stopped by the addition of stop solution. Absorbance is measured at 450 nm.

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed.



Add 100 μ L of samples and standards into appropriate wells.



Incubate for 1 hour at RT. Discard the solution and wash the wells four times with 300 μ L of washing solution.



Add 100 μ L of enzyme conjugate into each well.



Incubate for 30 minutes at RT. Discard the solution and wash the wells 4 times with 300 μ L of washing solution.



Add 100 μ L of substrate solution into each well.



Incubate for 10 - 25 minutes (the precise incubation time comes with the kit) at RT.



Add 50 μ L of Stop solution into each well. Read the absorbance at 450 nm immediately.

3. Precautions

Please read these instructions carefully prior to beginning the ELISA assay.

- All kit components have been formulated and quality control tested to function successfully as a kit.
- We understand that, occasionally, experimental protocols might need to be modified to meet unique experimental circumstances. However, we cannot guarantee the performance of the product outside the conditions detailed in this protocol booklet.
- Reagents should be treated as possible mutagens and should be handled with care and disposed of properly. Please review the Safety Datasheet (SDS) provided with the product for information on the specific components.
- Observe good laboratory practices. Gloves, lab coat, and protective eyewear should always be worn. Never pipet by mouth. Do not eat, drink or smoke in the laboratory areas.
- All biological materials should be treated as potentially hazardous and handled as such. They should be disposed of in accordance with established safety procedures.

4. Storage and Stability

Store ELISA kit at 2-8°C immediately upon receipt. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Materials Supplied section.

Aliquot components in working volumes before storing at the recommended temperature.

5. Limitations

- ELISA kit intended for research use only. Not for use in diagnostic procedures.
- Do not mix or substitute reagents or materials from other kit lots or vendors. Kits are QC tested as a set of components and performance cannot be guaranteed if utilized separately or substituted.

6. Materials Supplied

Item	Quantity	Storage Condition (Before prep)	Storage Condition (After prep)
anti-Human CDNF coated microplate (12 x 8 wells)	96 well	+2-8°C	+2-8°C
Human CDNF sample diluent	25 mL	+2-8°C	+2-8°C
Human CDNF standard A (0 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard B (15 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard C (30 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard D (60 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard E (120 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard F (240 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard G (480 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF standard H (960 pg/mL)	1 mL	+2-8°C	+2-8°C
Human CDNF enzyme conjugate	12 mL	+2-8°C	+2-8°C
10X Wash concentrate	50 mL	+2-8°C	+2-8°C
TMB Substrate solution	12 mL	+2-8°C	+2-8°C
Stop solution	12 mL	+2-8°C	+2-8°C

7. Materials Required, Not Supplied

These materials are not included in the kit, but will be required to successfully perform this assay:

- Pipettes and tips (100-500 μ L)
- ELISA plate washer
- Microplate reader (450 nm)
- Lid or sealing tape for microwell plate
- Microwell plate shaker

8. Technical Hints

- Samples generating values higher than the highest standard should be further diluted in the appropriate sample dilution buffers.
- Avoid foaming or bubbles when mixing or reconstituting components.
- Avoid cross contamination of samples or reagents by changing tips between sample, standard and reagent additions.
- Ensure plates are properly sealed or covered during incubation steps.
- Complete removal of all solutions and buffers during wash steps is necessary to minimize background.
- All samples should be mixed thoroughly and gently.
- Avoid multiple freeze/thaw of samples.
- When generating positive control samples, it is advisable to change pipette tips after each step.
- **This kit is sold based on number of tests. A 'test' simply refers to a single assay well. The number of wells that contain sample, control or standard will vary by product. Review the protocol completely to confirm this kit meets your requirements. Please contact our Technical Support staff with any questions.**

9. Reagent Preparation

- Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells.
- Prepare only as much reagent as is needed on the day of the experiment.

9.1 Anti-Human CDNF coated microplate (12 x 8 wells)

96 well microtiter plate coated with anti-human CDNF mouse monoclonal antibodies. 96 tests. Ready to use. Store at +2-8°C.

9.2 Human CDNF standards

8 tubes of 1 mL. Ready to use. Store at +2-8°C
(0, 15, 30, 60, 120, 240, 480, 960 pg/mL)

9.3 Human CDNF sample diluent

25 mL. Ready to use. Store at +2-8°C.

9.4 Human CDNF enzyme conjugate

12 mL. Ready to use. Store at +2-8°C.

9.5 10X Wash concentrate

Dilute 50 mL of 10X Wash concentrate with 450 mL of distilled water to prepare 1X Washing solution. 50 mL. Store at +2-8°C.

9.6 TMB Substrate solution

12 mL. Ready to use. Store at +2-8°C.

9.7 Stop solution

12 mL. Ready to use. Store at +2-8°C.

10. Sample Preparation

10.1 Samples

Dilute the samples in sample diluent.

11. Plate Preparation

- The 96 well plate included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).
- Differences in well absorbance or “edge effects” have not been observed with this assay.

12. Assay Procedure

- Equilibrate all materials and prepared reagents to room temperature prior to use.
 - It is recommended to assay all standards, controls and samples in duplicate.
- 12.1** Dilute 50 mL of wash concentrate with 450 mL of distilled water to prepare washing solution.
 - 12.2** Perform dilutions of each sample in sample diluent
 - 12.3** Add 100 μ L of samples and standards into appropriate wells in duplicate.
 - 12.4** Incubate the covered microplate for 1 hour at RT on a microwell plate shaker (300 rpm).
 - 12.5** Discard the solution and wash the wells 4 times with 300 μ L of washing solution.
 - 12.6** Add 100 μ L of enzyme conjugate into each well.
 - 12.7** Incubate the covered microplate for 30 min at RT on a microwell plate shaker (300 rpm).
 - 12.8** Discard the solution and wash the wells 4 times with 300 μ L of washing solution.
 - 12.9** Add 100 μ L of substrate solution into each well.
 - 12.10** Incubate the covered microplate for 10 - 25 minutes (the precise incubation time comes with the kit) at RT on a microwell plate shaker (300 rpm).
 - 12.11** Stop the reaction by adding 50 μ L of Stop solution into each well in the same order and time as for TMB distribution.
 - 12.12** Read the absorbance at 450 nm immediately.

13. Calculations

- 13.1 Standard curve: Calculate the mean absorbance for each standard. Subtract the blank value (standard A) from the mean absorbance's. Plot the value (absorbance) of each standard on a log-log scale. The use of software to generate a cubic spline fit curve is recommended
- 13.2 When generating a linear regression fit curve instead of a cubic spline fit curve only minor differences occur in human CDNF concentration calculation.
- 13.3 Validation of the assay: The mean absorbance of the Standard A (blank) should be below 0.1 AU (absorbance unit). The mean absorbance of the Standard H is usually above 1.0 AU.

14. Typical data

Typical standard curve – Data provided for demonstration purposes only. A new standard curve must be generated for each assay performed.

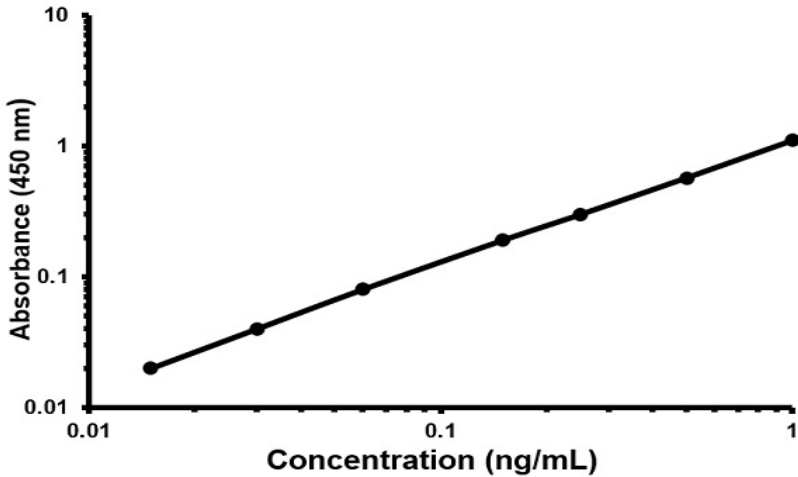


Figure 1. Cubic spline fit curve on a log-log scale

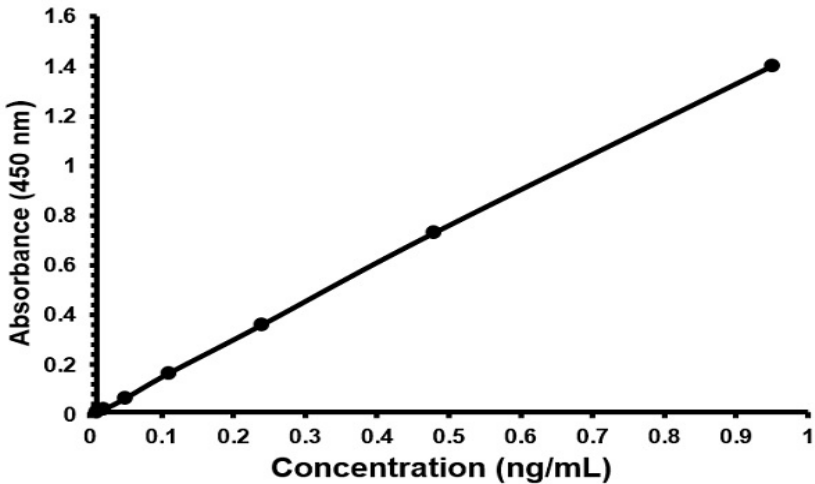


Figure 2. Linear regression (LLS) fit curve on a lin-lin scale

15. Typical sample values

SENSITIVITY –

The detection range is from 15 pg/mL to 960 pg/mL.

The detection limit is 1 pg/mL to 8 pg/mL, defined by the minimum human CDNF concentration deviating by 2 standard deviations (2SD) from that of the standard A. The test was performed by using 16 replicate determinations of standard A (blank) and standard B.

RECOVERY –

Human CDNF standards of 120, 480 and 960 pg/mL were added to equal volumes of three samples (serum or plasma) containing a low (50 pg/mL), a medium (220 pg/mL) and a high (730 pg/mL) concentration of human CDNF. The theoretical concentration and the recovered concentration were calculated.

Sample	Added conc. (pg/mL)	Expected conc. (pg/mL)	Obtained conc. (pg/mL)	Recovery %
Low	0	N/A	50	100
	120	85	70	82
	480	290	260	90
	960	510	510	100
Medium	0	N/A	220	100
	120	170	120	71
	480	350	330	94
	960	590	620	105
High	0	N/A	730	100
	120	430	310	72
	480	610	600	98
	960	850	880	103

LINEARITY OF DILUTION –

Three samples (plasma or serum) were diluted with sample diluent. The concentration of human CDNF in each diluted sample was measured. The results are shown as a change in percentage from the lowest dilution (corrected with the dilution factor).

Sample	Dilution	Conc. (pg/mL)	%
Serum	1:10	870	100
	1:20	950	109
	1:40	950	109
Plasma #1	1:10	1420	100
	1:20	1270	89
	1:40	1280	90
Plasma #2	1:03	120	100
	1:06	100	83
	1:12	130	108

PRECISION –

Intra-assay precision:

Sample	Number of measures	Mean (pg/mL)	CV%
1	15	677	3.7
2	15	272	2.8
3	15	167	2.5

Inter-assay precision:

Sample	Number of assays	Mean (pg/mL)	CV%
1	2	19	8.1
2	2	155	16.0
3	2	439	20.8

Problem	Cause	Solution
Poor standard curve	Inaccurate Pipetting	Check pipettes
	Improper standard dilution	Prior to opening, briefly spin the stock standard tube and dissolve the powder thoroughly by gentle mixing
Low Signal	Incubation times too brief	Ensure sufficient incubation times standard/sample incubation
	Inadequate reagent volumes or improper dilution	Check pipettes and ensure correct preparation
	Incubation times with TMB too brief	Ensure sufficient incubation time until blue color develops prior addition of Stop solution
Large CV	Plate is insufficiently washed	Review manual for proper wash technique. If using a plate washer, check all ports for obstructions.
	Contaminated wash buffer	Prepare fresh wash buffer
Low sensitivity	Improper storage of the ELISA kit	All components 4°C. Keep TMB substrate solution protected from light.

16. Troubleshooting

17. Notes

Technical Support

Copyright © 2016 Abcam, All Rights Reserved. The Abcam logo is a registered trademark. All information / detail is correct at time of going to print.

Austria

wissenschaftlicherdienst@abcam.com | 019-288-259

France

supportscientifique@abcam.com | 01.46.94.62.96

Germany

wissenschaftlicherdienst@abcam.com | 030-896-779-154

Spain

soportecientifico@abcam.com | 91-114-65-60

Switzerland

technical@abcam.com

Deutsch: 043-501-64-24 | Français: 061-500-05-30

UK, EU and ROW

technical@abcam.com | +44(0)1223-696000

Canada

ca.technical@abcam.com | 877-749-8807

US and Latin America

us.technical@abcam.com | 888-772-2226

Asia Pacific

hk.technical@abcam.com | (852) 2603-6823

China

cn.technical@abcam.com | 400 921 0189 / +86 21 2070 0500

Japan

technical@abcam.co.jp | +81-(0)3-6231-0940

Singapore

sg.technical@abcam.com | 800 188-5244

Australia

au.technical@abcam.com | +61-(0)3-8652-1450

New Zealand

nz.technical@abc.com | +64-(0)9-909-7829