

## ab284005 – Human HDHD2 ELISA Kit

For the quantitative measurement of HDHD2 in Human CSF and cell lysate samples.  
For research use only - not intended for diagnostic use.

For overview, typical data and additional information please visit: [www.abcam.com/ab284005](http://www.abcam.com/ab284005)

**Storage and Stability:** Store kit at +4°C immediately upon receipt, apart from the Standard, Biotinylated Antibody and SP Conjugate, which should be stored at -20°C. Kit has a storage time of 1 year from receipt, providing components have not been reconstituted.

### Materials Supplied

Item	Quantity	Storage Condition
100X Streptavidin-Peroxidase Conjugate	80 µl	-20°C
10X Diluent M Concentrate	20 ml	4°C
1X Standard Diluent	2 ml	4°C
20X Wash Buffer Concentrate	2 x 30 ml	4°C
50X Biotinylated Human HDHD2 Antibody	1 vial	-20°C
Chromogen Substrate	7 mL	4°C
Human HDHD2 Microplate	1 unit	4°C
Human HDHD2 Standard	1 vial	-20°C
Sealing Tapes	3 units	4°C
Stop Solution	11 mL	4°C

### Materials Required, Not Supplied

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

Microplate reader capable of measuring absorbance at 450 nm  
Pipettes (1-20 µl, 20-200 µl, 200-1000 µl, and multiple channel)  
Deionized or distilled reagent grade water

### Reagents 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.

If crystals have formed in the concentrate, mix gently until the crystals have completely dissolved.

**Δ Note:** Concentration of the kit components are lot-specific, and the end user should always refer to the vial label.

**10X Diluent M Concentrate:** Dilute the Diluent M Concentrate 10-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved. Store for up to 30 days at 2-8°C.

**50X Biotinylated Human HDHD2 Antibody:** Spin down the antibody briefly and dilute the desired amount of the antibody 50-fold with Diluent M to produce a 1x solution. The undiluted antibody should be stored at -20°C.

**20X Wash Buffer Concentrate:** Dilute the Wash Buffer Concentrate 20-fold with reagent grade water to produce a 1x solution. When diluting the concentrate, make sure to rinse the bottle thoroughly to extract any precipitates left in the bottle. Mix the 1x solution gently until the crystals have completely dissolved.

**100X SP Conjugate:** Spin down the SP Conjugate briefly and dilute the desired amount of the conjugate 100-fold with Diluent M to produce a 1x solution. The undiluted conjugate should be stored at -20°C.

### Standard Preparation

Always prepare a fresh set of standards for every use.

Prepare serially diluted standards immediately prior to use.

Any remaining standard should be stored at -20°C after reconstitution and used within 30 days. The following section describes the preparation of a standard curve for duplicate measurements (recommended).

Reconstitution of the HDHD2 Standard vial to prepare a 4 ng/mL Stock Standard.

- First consult the HDHD2 Standard vial to determine the mass of protein in the vial.
- Calculate the appropriate volume of 1X Diluent M to add when resuspending the HDHD2 Standard vial to produce a 4 ng/mL Stock Standard by using the following equation:

- o  $C_s$  = Starting mass of HDHD2 Standard (see vial label) (ng)
- o  $C_f$  = The 4 ng/ml HDHD2 Stock Standard final required concentration
- o  $V_D$  = Required volume of 1X Diluent M for reconstitution (µL)
- o Calculate total required volume 1X Diluent M for resuspension:

$$(C_s / C_f) \times 1,000 = V_D$$

**Δ Note:** This example is for demonstration purposes only. Please remember to check your standard vial for the actual amount of standard provided.

$C_s$  = 1 ng of HDHD2 Standard in vial

$C_f$  = 2 ng/mL HDHD2 Standard #1 final concentration

$V_D$  = Required volume of 1X Diluent M for reconstitution (1 ng / 2 ng/mL) x 1,000 = 500 µL

- Reconstitute the HDHD2 Standard vial by adding the appropriate calculated amount VD of 1X Diluent M to the vial to generate the 2 ng/mL HDHD2 Standard #1. Mix gently and thoroughly.
- Allow the reconstituted 2 ng/mL HDHD2 Standard #1 to sit for 10 minutes with gentle agitation prior to making subsequent dilutions
- Label tubes #2 – 8.
- Add 120 µL of 1X Diluent M to tube #2 – 8.
- To prepare Standard #2, add 120 µL of the Standard #1 into tube #2 and mix gently.
- To prepare Standard #3, add 120 µL of the Standard #2 into tube #3 and mix gently
- Using the table below as a guide, prepare subsequent serial dilutions. 1X Diluent M serves as the zero standard (0 ng/mL).

Standard #	Volume to dilute (µL)	Volume 1X Diluent M (µL)	HDHD2 (ng/ml)
1	240 (Stock Standard)	-	2.0
2	120 (Standard #1)	120	1.0
3	120 (Standard #2)	120	0.5
4	120 (Standard #3)	120	0.25
5	120 (Standard #4)	120	0.125
6	120 (Standard #5)	120	0.063
7	120 (Standard #6)	120	0.031
8	-	120	0.0

## Sample Preparation

**CSF:** Collect cerebrospinal fluid (CSF) using sample pot. Centrifuge samples at 3000 x g for 10 minutes. The sample is suggested for use at 1x; however, user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -80°C for up to 3 months. Avoid repeated freeze-thaw cycles.

**Cell Lysate:** Rinse cell with cold PBS and then scrape the cell into a tube with 5 ml of cold PBS and 0.5 M EDTA. Centrifuge suspension at 1500 rpm for 10 minutes at 4°C and aspirate supernatant. Resuspend pellet in ice-cold Lysis Buffer (PBS, 1% Triton X-100, protease inhibitor cocktail). For every 1 x 10<sup>6</sup> cells, add approximately 100 µl of ice-cold Lysis Buffer. Incubate on ice for 60 minutes. Centrifuge at 13000 rpm for 30 minutes at 4°C and collect supernatant. If necessary, dilute samples into Diluent M; user should determine optimal dilution factor depending on application needs. The undiluted samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.

## Plate Preparation

The 96 well plate strips included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.

Unused plate strips should be immediately returned to the foil pouch containing the desiccant pack, resealed and stored at 4°C.

For each assay performed, a minimum of two wells must be used as the zero control.

For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).

## Assay Procedure

Equilibrate all materials and prepared reagents to room temperature prior to use. We recommend that you assay all standards, controls and samples in duplicate

1. Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccants inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
2. Add 50 µl of Human HDHD2 Standard or sample to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 2 hours. Start the timer after the last addition.
3. Wash the microplate manually or automatically using a microplate washer. Invert the plate and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If washing manually, wash five times with 200 µl of Wash Buffer per well. Invert the plate each time and decant the contents; hit 4-5 times on absorbent material to completely remove the liquid. If using a microplate washer, wash six times with 300 µl of Wash Buffer per well; invert the plate and hit 4-5 times on absorbent material to completely remove the liquid.
4. Add 50 µl of Biotinylated Human HDHD2 Antibody to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 1 hour.
5. Wash the microplate as described above.
6. Add 50 µl of SP Conjugate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Cover wells with a sealing tape and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.
7. Wash the microplate as described above.

8. Add 50 µl of Chromogen Substrate to each well. Gently tap plate to thoroughly coat the wells. Break any bubbles that may have formed. Incubate in ambient light for 15 minutes or until the optimal blue color density develops.
9. Add 50 µl of Stop Solution to each well. The color will change from blue to yellow. Gently tap plate to ensure thorough mixing. Break any bubbles that may have formed.
10. Read the absorbance on a microplate reader at a wavelength of 450 nm immediately. If wavelength correction is available, subtract readings at 570 nm from those at 450 nm to correct optical imperfections. Otherwise, read the plate at 450 nm only. Please note that some unstable black particles may be generated at high concentration points after stopping the reaction for about 10 minutes, which will reduce the readings.

## Ab284005 – Human HDHD2 ELISA Kit

### Additional information

#### CALCULATION

1. Calculate the mean value of the duplicate or triplicate readings for each standard and sample.
2. To generate a standard curve, plot the graph using the standard concentrations on the x-axis and the corresponding mean 450 nm absorbance (OD) on the y-axis. The best fit line can be determined by regression analysis using log-log or four-parameter logistic curve fit.
3. Determine the unknown sample concentration from the Standard Curve and multiply the value by the dilution factor.

#### TYPICAL DATA

Typical data provided for demonstration purposes only.

Standard #	ng/ml	Average OD
1	2.0	2.219
2	1.0	1.381
3	0.5	0.752
4	0.25	0.433
5	0.125	0.265
6	0.063	0.180
7	0.031	0.140
8	0.0	0.092

#### PERFORMANCE CHARACTERISTICS

1. This assay recognizes both natural and recombinant human HDHD2.
2. The minimum detectable dose of human HDHD2 as calculated by 2SD from the mean of a zero standard was established to be 21 pg/ml.
3. Intra-assay precision was determined by testing three reference control samples twenty times in one assay.
4. Inter-assay precision was determined by testing three reference control samples in twenty assays.

	Intra-Assay Precision	Inter-Assay Precision
CV (%)	5.8	9.3

#### RECOVERY

Standard Added Value	0.063 – 0.5 ng/ml
Recovery %	88 – 115
Average Recovery %	96

#### LINEARITY

Average Percentage of Expected Value (%)	
Sample Dilution	293T (Human Embryonic Kidney) Cell Lysate
1:5	89
1:10	100
1:20	111

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:

[www.abcam.com/protocols/the-complete-elisa-guide](http://www.abcam.com/protocols/the-complete-elisa-guide)

For all technical or commercial enquiries please go to:

[www.abcam.com/contactus](http://www.abcam.com/contactus)

[www.abcam.cn/contactus](http://www.abcam.cn/contactus) (China)

[www.abcam.co.jp/contactus](http://www.abcam.co.jp/contactus) (Japan)

Copyright © 2024 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.

Version 2a | Thursday, 14 November 2024