

Version 16 Last updated 12 April 2018

ab108788 – Albumin Human ELISA Kit

For the quantitative measurement of Human Albumin in urine, saliva, milk, plasma, serum, cerebrospinal fluid, cell culture supernatants and cell lysate

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. Standard Preparation	10
11. Sample Preparation	12
12. Plate Preparation	13
13. Assay Procedure	13
14. Calculations	14
15. Typical Data	15
16. Typical Sample Values	16
17. Assay Specificity	16
18. Species Reactivity	17
19. Troubleshooting	18
20. Notes	20

1. Overview

Abcam's Albumin Human *in vitro* ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the quantitative measurement of albumin levels in urine, saliva, milk, cerebrospinal fluid, and cell culture supernatants.

An albumin specific antibody has been pre-coated onto 96-well plates and blocked. Standards or test samples are added to the wells and subsequently an albumin specific biotinylated detection antibody is added and then followed by washing with wash buffer. Streptavidin-Peroxidase Complex is added and unbound conjugates are washed away with wash buffer. TMB is then used to visualize Streptavidin-Peroxidase enzymatic reaction. TMB is catalyzed by Streptavidin-Peroxidase to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the amount of albumin captured in plate.

Albumin, a serum hepatic protein, is the most abundant protein in serum. It contributes to the maintenance of oncotic pressure as well as the transport of hydrophobic molecules. Serum albumin level has been linked in clinical practice to several diseases. Low albumin levels can suggest liver disease, kidney disease, inflammation, shock, and malnutrition. On the other hand, high albumin levels usually reflect dehydration.

2. Protocol Summary

Prepare all reagents, samples, and standards as instructed



Add standard or sample to appropriate wells.

Incubate at room temperature.



Wash and add prepared biotin antibody to each well. Incubate at room temperature.



Wash and add prepared Streptavidin-Peroxidase Conjugate. Incubate at room temperature.



Add Chromogen Substrate to each well. Incubate at room temperature



Add Stop Solution to each well. Read immediately.

3. Precautions

Please read these instructions carefully prior to beginning the 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 kit at +4°C immediately upon receipt, apart from the SP Conjugate & Biotinylated Antibody, which should be stored at -20°C. 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.

5. Limitations

- Assay 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
Albumin Microplate (12 x 8 wells)	96 wells	4°C
Albumin Standard	1 vial	4°C
10X Diluent N Concentrate	30 mL	4°C
Biotinylated Human Albumin Antibody	1 vial	-20°C
100X Streptavidin-Peroxidase Conjugate (SP Conjugate)	80 µL	-20°C
Chromogen Substrate	8 mL	4°C
Stop Solution	12 mL	4°C
20X Wash Buffer Concentrate	2 x 30 mL	4°C
Sealing Tapes	3	N/A

7. Materials Required, Not Supplied

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

- 1 Microplate reader capable of measuring absorbance at 450 nm.
- Precision pipettes to deliver 1 μ L to 1 mL volumes.
- Adjustable 1-25 mL pipettes for reagent preparation.
- 100 mL and 1 liter graduated cylinders.
- Absorbent paper.
- Distilled or deionized water.
- Log-log graph paper or computer and software for ELISA data analysis.
- 8 tubes to prepare standard or sample dilutions.

8. Technical Hints

- 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.
- Selected components in this kit are supplied in surplus amount to account for additional dilutions, evaporation, or instrumentation settings where higher volumes are required. They should be disposed of in accordance with established safety procedures.
- Make sure all buffers and solutions are at room temperature before starting the experiment.
- 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.
- Make sure you have the right type of plate for your detection method of choice.
- Make sure the heat block/water bath and microplate reader are switched on before starting the experiment.

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 1X Diluent N

Dilute the 10X Diluent N Concentrate 1:10 with reagent grade water. Mix gently and thoroughly. *Store for up to 1 month at 4°C.*

9.2 1X Wash Buffer

Dilute the 20X Wash Buffer Concentrate 1:20 with reagent grade water. Mix gently and thoroughly.

9.3 1X Biotinylated Albumin Detector Antibody

9.3.1 The stock Biotinylated Albumin Antibody must be diluted with 1X Diluent N according to the label concentration to prepare 1X Biotinylated Albumin Antibody for use in the assay procedure. Observe the label for the “X” concentration on the vial of Biotinylated Albumin Antibody.

9.3.2 Calculate the necessary amount of 1X Diluent N to dilute the Biotinylated Albumin Antibody to prepare a 1X Biotinylated Albumin Antibody solution for use in the assay procedure according to how many wells you wish to use and the following calculation:

Number of Wells Strips	Number of Wells	(V _T) Total Volume of 1X Biotinylated Detector Antibody (µL)
4	32	1,760
6	48	2,640
8	64	3,520
10	80	4,400
12	96	5,280

Any remaining solution should be frozen at -20°C.

Where:

C_S = Starting concentration (X) of stock Biotinylated Albumin Antibody (variable)

C_F = Final concentration (always = 1X) of 1X Biotinylated Albumin Detector Antibody solution for the assay procedure

V_T = Total required volume of 1X Biotinylated Albumin Detector Antibody solution for the assay procedure

V_A = Total volume of (X) stock Biotinylated Albumin Antibody

V_D = Total volume of 1X Diluent N required to dilute (X) stock Biotinylated Albumin Antibody to prepare 1X Biotinylated Detector Antibody solution for assay procedures

Calculate the volume of (X) stock Biotinylated Antibody required for the given number of desired wells:

$$(C_F / C_S) \times V_T = V_A$$

Calculate the final volume of 1X Diluent N required to prepare the 1X Biotinylated Albumin Detector Antibody:

$$V_T - V_A = V_D$$

Example:

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

C_S = 50X Biotinylated Albumin Antibody stock

C_F = 1X Biotinylated Albumin Detector Antibody solution for use in the assay procedure

V_T = 3,520 μ L (8 well strips or 64 wells)

$$(1X/50X) \times 3,520 \mu\text{L} = 70.4 \mu\text{L}$$

$$3,520 \mu\text{L} - 70.4 \mu\text{L} = 3,449.6 \mu\text{L}$$

V_A = 70.4 μ L total volume of (X) stock Biotinylated Albumin Antibody required

V_D = 3,449.6 μ L total volume of 1X Diluent N required to dilute the 50X stock Biotinylated Antibody to prepare 1X Biotinylated Albumin Detector Antibody solution for assay procedures.

- 9.3.3 First spin the Biotinylated Albumin Antibody vial to collect the contents at the bottom.
- 9.3.4 Add calculated amount V_A of stock Biotinylated Albumin Antibody to the calculated amount V_D of 1X Assay Diluent N. Mix gently and thoroughly.

9.4 1X SP Conjugate

Spin down the 100X Streptavidin-Peroxidase Conjugate (SP Conjugate) briefly and dilute the desired amount of the conjugate 1:100 with 1X Diluent N.

Any remaining solution should be frozen at -20°C.

10. Standard Preparation

- Always prepare a fresh set of standards for every use.
- Discard working standard dilutions after use as they do not store well.
- The following section describes the preparation of a standard curve for duplicate measurements (recommended).

10.1 Reconstitute the Albumin Stock to generate a 200 ng/mL Standard #1.

- 10.1.1 First consult the Albumin Standard vial to determine the mass of protein in the vial.
- 10.1.2 Calculate the appropriate volume of 1X Diluent N to add when resuspending the Albumin Standard vial to produce a 200 ng/mL Albumin Standard stock by using the following equation:

C_S = Starting mass of Albumin Standard (see vial label) (ng)

C_F = The 200 ng/mL Albumin **Standard #1** final required concentration

V_D = Required volume of 1X Diluent N for reconstitution (μ L)

Calculate total required volume 1X Diluent N for resuspension:

$$(C_S / C_F) \times 1,000 = V_D$$

Example:

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 = 800 ng of Albumin Standard in vial

C_F = 200 ng/mL Albumin **Standard #1** final concentration

V_D = Required volume of 1X Diluent N for reconstitution

$$(800 \text{ ng} / 200 \text{ ng/mL}) \times 1,000 = 4,000 \mu\text{L}$$

- 10.1.3 Reconstitute the Albumin Standard vial by adding the appropriate calculated amount V_D of 1X Diluent N to the vial to generate the 200 ng/mL Albumin **Standard #1**. Mix gently and thoroughly.
- 10.2 Allow the reconstituted 200 ng/mL Albumin **Standard #1** to sit for 10 minutes with gentle agitation prior to making subsequent dilutions
- 10.3 Label five tubes #2 – 8.
- 10.4 Add 120 μ L of 1X Diluent N to tube #2 – 8.
- 10.5 To prepare **Standard #2**, add 120 μ L of the **Standard #1** into tube #2 and mix gently.
- 10.6 To prepare **Standard #3**, add 120 μ L of the **Standard #2** into tube #3 and mix gently.
- 10.7 Using the table below as a guide, prepare subsequent serial dilutions. 1X Diluent N serves as the zero standard (0 ng/mL).

Standard #	Volume to Dilute (μ L)	Volume Diluent N (μ L)	Total Volume (μ L)	Starting Conc. (ng/mL)	Final Conc. (ng/mL)
1	Step 10.1				200.0
2	120	120	240	200.0	100.0
3	120	120	240	100.0	50.00
4	120	120	240	50.00	25.00
5	120	120	240	25.00	12.50
6	120	120	240	12.50	6.250
7	120	120	240	6.250	3.125
8	-	120	120	-	0

11. Sample Preparation

11.1 Cell Culture Supernatants:

Centrifuge cell culture media at 1500 rpm for 10 minutes at 4°C to remove debris and collect supernatant. Samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.

11.2 Urine:

Collect urine using sample pot. Centrifuge samples at 800 x *g* for 10 minutes. Dilute urine samples 1:100 - 1:500 into 1X Diluent N and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.3 Saliva:

Collect saliva using sample tube. Centrifuge samples at 800 x *g* for 10 minutes. Dilute saliva samples 1:200 – 1:2000 into 1X Diluent N and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.4 Milk:

Collect milk using sample tube. Centrifuge samples at 800 x *g* for 10 minutes. Dilute milk samples 1:2000 - 1:10,000 into 1X Diluent N and assay. The undiluted samples can be stored at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.5 Cerebrospinal fluid:

Collect cerebrospinal fluid (CSF) using sample tube. Centrifuge samples at 3,000 x *g* for 10 minutes. Dilute CSF samples 1:2000 - 1:10 000 into 1X Diluent N and assay. The undiluted samples can be stored at -80°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

11.6 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 (10 mM Tris, pH 8.0, 130 mM NaCl, 1% Triton X-100, protease inhibitor cocktail). For every 1×10^6 cells, add approximately 100 μ L of ice-cold Lysis Buffer. Incubate on ice for 60 minutes. Centrifuge at 13000 rpm for 30 minutes at 4°C and collect supernatant. Samples can be stored at -80°C. Avoid repeated freeze-thaw cycles.

Applicable samples may also include biofluids, cell culture, and tissue homogenates. If necessary, user should determine optimal dilution factor depending on application needs.

12. 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 well plate strips should be returned to the plate packet and stored at 4°C.
- For statistical reasons, we recommend each sample should be assayed with a minimum of two replicates (duplicates).

13. 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.

- 13.1** Prepare all reagents, working standards and samples as instructed. Equilibrate reagents to room temperature before use. The assay is performed at room temperature (18-25°C).
- 13.2** Remove excess microplate strips from the plate frame and return them immediately to the foil pouch with desiccant inside. Reseal the pouch securely to minimize exposure to water vapor and store in a vacuum desiccator.
- 13.3** Add 50 µL of Albumin Standard or sample per well. Cover wells with a sealing tape and incubate for 1 hour. Start the timer after the last addition.
- 13.4** Wash five times with 200 µL of 1X Wash Buffer manually. Invert the plate each time and decant the contents; tap it 4-5 times on absorbent paper towel to completely remove the liquid. If using a machine wash six times with 300 µL of 1X Wash Buffer and then invert the plate, decant the contents; tap it 4-5 times on absorbent paper towel to completely remove the liquid.
- 13.5** Add 50 µL of 1X Biotinylated Albumin Antibody to each well and incubate for 30 minutes.
- 13.6** Wash microplate as described above.
- 13.7** Add 50 µL of 1X SP Conjugate to each well and incubate for 30 minutes. Turn on the microplate reader and set up the program in advance.

- 13.8** Wash the microplate as described above.
- 13.9** Add 50 μL of Chromogen Substrate per well and incubate for about 25 minutes or till the optimal blue colour density develops. Gently tap plate to ensure thorough mixing and break the bubbles in the well with pipette tip.
- 13.10** Add 50 μL of Stop Solution to each well. The color will change from blue to yellow.
- 13.11** 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.

14. Calculations

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

15. Typical Data

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

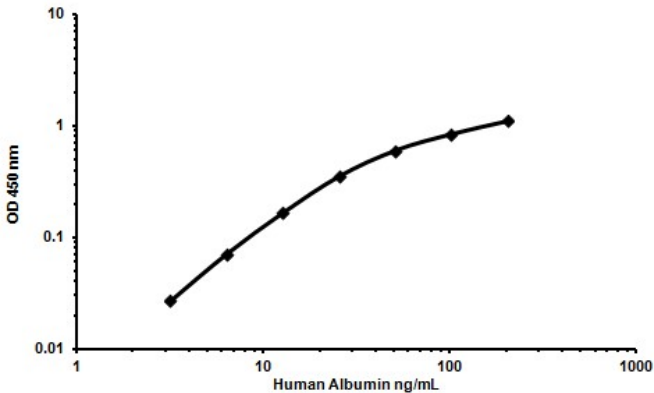


Figure 1. Example of Albumin standard curve. The standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean +/- SD) are graphed.

16. Typical Sample Values

SENSITIVITY –

The minimum detectable dose (MDD) of Albumin is typically 1.9 ng/ml.

PRECISION –

	Intra-assay Precision	Inter-Assay Precision
CV (%)	4.4	9

RECOVERY –

Standard Added Value	12.5 – 100 ng/ml
Recovery (%)	91-114 %
Average Recovery (%)	97 %

Linearity of Dilution

Linearity of dilution is determined based on interpolated values from the standard curve. Linearity of dilution defines a sample concentration interval in which interpolated target concentrations are directly proportional to sample dilution.

Average Percentage of Expected Value (%)	
Dilution Factor	Milk
1:3,000	92
1:6,000	98
1:12,000	105

17. Assay Specificity

This kit recognizes Albumin in urine, saliva, milk, plasma, serum, cerebrospinal fluid and cell culture supernatants

18. Species Reactivity

This kit recognizes Albumin.

Species	Cross Reactivity (%)
Canine	None
Mouse	None
Monkey	None
Bovine	None
Rat	None
Swine	None
Rabbit	None

Please contact our Technical Support team for more information.

19. Troubleshooting

Problem	Reason	Solution
Poor standard curve	Improper standard dilution	Confirm dilutions made correctly
	Standard improperly reconstituted (if applicable)	Briefly spin vial before opening; thoroughly re-suspend powder (if applicable)
	Standard degraded	Store sample as recommended
	Curve doesn't fit scale	Try plotting using different scale
Low signal	Incubation time too short	Try overnight incubation at 4°C
	Target present below detection limits of assay	Decrease dilution factor; concentrate samples
	Precipitate can form in wells upon substrate addition when concentration of target is too high	Increase dilution factor of sample
	Using incompatible sample type (e.g. serum vs. cell extract)	Detection may be reduced or absent in untested sample types
	Sample prepared incorrectly	Ensure proper sample preparation/dilution

Large CV	Bubbles in wells	Ensure no bubbles present prior to reading plate
	All wells not washed equally/thoroughly	Check that all ports of plate washer are unobstructed wash wells as recommended
	Incomplete reagent mixing	Ensure all reagents/master mixes are mixed thoroughly
	Inconsistent pipetting	Use calibrated pipettes and ensure accurate pipetting
	Inconsistent sample preparation or storage	Ensure consistent sample preparation and optimal sample storage conditions (eg. minimize freeze/thaws cycles)
High background/ Low sensitivity	Wells are insufficiently washed	Wash wells as per protocol recommendations
	Contaminated wash buffer	Make fresh wash buffer
	Waiting too long to read plate after adding STOP solution	Read plate immediately after adding STOP solution
	Improper storage of ELISA kit	Store all reagents as recommended. Please note all reagents may not have identical storage requirements.
	Using incompatible sample type (e.g. Serum vs. cell extract)	Detection may be reduced or absent in untested sample types

20. Notes

Technical Support

Copyright © 2017 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 | +86-21-5110-5938 | 400-628-6880

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@abcam.com | +64-(0)9-909-7829