

## ab179887 – Human Albumin SimpleStep ELISA® Kit

For the quantitative measurement of Albumin in human serum, plasma (citrate), plasma (EDTA), plasma (heparin), and cell culture supernatant.  
For research use only - not intended for diagnostic use.

This kit is available in a 384-well plate format. This plate utilises smaller volumes of standards and samples per well. Directions for using this format can be found on pg. 9.

**Storage and Stability:** Store kit at 2-8°C immediately upon receipt. Refer to list of materials supplied for storage conditions of individual components. Observe the storage conditions for individual prepared components in the Standard Preparation and Reagent preparation sections.

### Materials Supplied

| Item                                       | Quantity<br>(1x96 tests) | Quantity<br>(10x96 tests) | Storage<br>Condition |
|--------------------------------------------|--------------------------|---------------------------|----------------------|
| 10X Human Albumin Capture Antibody         | 600 µL                   | 10 x 600 µL               | +4°C                 |
| 10X Human Albumin Detector Antibody        | 600 µL                   | 10x600 µL                 | +4°C                 |
| Human Albumin Lyophilized Purified Protein | 2 Vials                  | 20 Vials                  | +4°C                 |
| Antibody Diluent 4BI                       | 6 mL                     | 10 x 6 mL                 | +4°C                 |
| Sample Diluent NS                          | 2 x 50 mL                | 4 x 250 mL                | +4°C                 |
| 10X Wash Buffer PT                         | 20 mL                    | 200 mL                    | +4°C                 |
| TMB Development Solution                   | 12 mL                    | 120 mL                    | +4°C                 |
| Stop Solution                              | 12 mL                    | 120 mL                    | +4°C                 |
| SimpleStep Pre-Coated 96-Well Microplate   | 96 wells                 | 10x96 wells               | +4°C                 |
| Plate Seal                                 | 1                        | 10                        | +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 or 600 nm.  
Deionized water.  
Multi- and single-channel pipettes.  
Tubes for standard dilution.  
Plate shaker for all incubation steps.  
Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).

### Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for 96 wells. The sample volumes below are sufficient for 48 wells (6 x 8-well strips); adjust volumes as needed for the number of strips in your experiment.

Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

**1X Wash Buffer PT:** Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

**Antibody Cocktail:** Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 4BI. To make 3 mL of the Antibody Cocktail combine 300 µL 10X Capture Antibody and 300 µL 10X Detector Antibody with 2.4 mL Antibody Diluent 4BI. Mix thoroughly and gently.

### Standard Preparation

Always prepare a fresh set of standards for every use. Discard working standard dilutions after use as they do not store well. Excess Stock Standard can be aliquoted into single use sized aliquots and stored at - 80°C. Avoid repeated freeze-thaw cycles.

The following section describes the preparation of a standard curve for duplicate measurements (recommended).

1. Reconstitute the Albumin standard sample by adding the volume of Sample Diluent NS indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 100 ng/mL **Stock Standard** Solution.
2. Label eight tubes, Standards 1–8.
3. Add 220 µL of Sample Diluent NS into tube number 1 and 150 µL of Sample Diluent NS into numbers 2-8.
4. Use the **Stock Standard** to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:

| Standard # | Dilution<br>Sample    | Volume to<br>Dilute (µL) | Volume of<br>Diluent (µL) | Starting<br>Conc.<br>(ng/mL) | Final<br>Conc.<br>(ng/mL) |
|------------|-----------------------|--------------------------|---------------------------|------------------------------|---------------------------|
| 1          | <b>Stock Standard</b> | 180                      | 220                       | 100                          | 45                        |
| 2          | Standard#1            | 150                      | 150                       | 45                           | 22.5                      |
| 3          | Standard#2            | 150                      | 150                       | 22.5                         | 11.25                     |
| 4          | Standard#3            | 150                      | 150                       | 11.25                        | 5.63                      |
| 5          | Standard#4            | 150                      | 150                       | 5.63                         | 2.81                      |
| 6          | Standard#5            | 150                      | 150                       | 2.81                         | 1.41                      |
| 7          | Standard#6            | 150                      | 150                       | 1.41                         | 0.70                      |
| 8          | Blank Control         | 0                        | 150                       | 0                            | 0                         |

Sample Preparation

| Typical Sample Dynamic Range       |                             |
|------------------------------------|-----------------------------|
| Sample Type                        | Range                       |
| Human Serum                        | 1:64,000,000 – 1:4,000,000  |
| Human Plasma – EDTA                | 1:128,000,000 – 1:4,000,000 |
| Human Plasma – Citrate             | 1:64,000,000 – 1:8,000,000  |
| Human Plasma – Heparin             | 1:128,000,000 – 1:8,000,000 |
| Cell Culture Media (RPMI+10% FBS)* | 0.32% – 10%                 |

\*Based on spiked data

**Serum** Samples should be collected into a serum separator tube. After clot formation, centrifuge samples at 2,000 x g for 10 minutes and collect serum. Dilute samples at least 1:4,000,000 into Sample Diluent NS and assay. Store un-diluted serum at -20°C or below. Avoid repeated freeze-thaw cycles.

**Plasma** Collect plasma using citrate, EDTA or heparin. Centrifuge samples at 2,000 x g for 10 minutes. Dilute plasma – EDTA samples at least 1:4,000,000 into Sample Diluent NS and assay. Dilute plasma – Citrate and plasma – Heparin samples at least 1:8,000,000 into Sample Diluent NS and assay. Store un-diluted plasma samples at -20°C or below for up to 3 months. Avoid repeated freeze-thaw cycles.

**Cell Culture Supernatants** Centrifuge cell culture media at 2,000 x g for 10 minutes to remove debris. Collect supernatants. Dilute samples at least 1:10 into Sample Diluent NS and assay. Store un-diluted samples at -20°C or below. 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). Differences in well absorbance or “edge effects” have not been observed with this assay.

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. Prepare all reagents, working standards, and samples as directed in the previous sections.
2. Remove excess microplate strips from the plate frame, return them to the foil pouch containing the desiccant pack, reseal and return to 4°C storage.
3. Add 50 µL of all sample or standard to appropriate wells.
4. Add 50 µL of the Antibody Cocktail to each well.
5. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 400 rpm.
6. Wash each well with 3 x 350 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 350 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
7. Add 100 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 400 rpm.  
*Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.*  
Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.
8. Add 100 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading.
9. Alternative to 7 – 8: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

| Mode        | Kinetic                |
|-------------|------------------------|
| Wavelength: | 600 nm                 |
| Time:       | up to 20 min           |
| Interval:   | 20 sec - 1 min         |
| Shaking:    | Shake between readings |

**Note** that an endpoint reading can also be recorded at the completion of the kinetic read by adding 100 µL Stop Solution to each well and recording the OD at 450 nm.

## ab179887 – Human Albumin SimpleStep ELISA® Kit

### Additional information

#### ASSAY SPECIFICITY

This kit is designed for the quantification of human Albumin.

The standard protein in this kit is natively purified full length human Albumin.

Native signal was detected in serum, plasma (citrate), plasma (EDTA), and plasma (heparin) sample types.

Spiked protein experiments were used to validate cell culture supernatant sample types.

Saliva, urine, milk, CSF, cell extract, and tissue extract samples have not been tested with this kit.

#### SPECIES REACTIVITY

Other species reactivity was determined by measuring 1:4,000,000 serum samples of various species, interpolating the protein concentrations from the human standard curve, and expressing the interpolated concentrations as a percentage of the protein concentration in human serum assayed at the same dilution.

This assay recognizes monkey Albumin.

Reactivity < 3% was determined for the following species: Mouse, Rat, Cow

Other species reactivity not determined.

#### CALCULATION

- Calculate the average absorbance value for the blank control (zero) standards. Subtract the average blank control standard absorbance value from all other absorbance values.
- Create a standard curve by plotting the average blank control subtracted absorbance value for each standard concentration (y-axis) against the target protein concentration (x-axis) of the standard. Use graphing software to draw the best smooth curve through these points to construct the standard curve.  
 $\Delta$  Note: Most microplate reader software or graphing software will plot these values and fit a curve to the data. A four parameter curve fit (4PL) is often the best choice; however, other algorithms (e.g. linear, semi-log, log/log, 4 parameter logistic) can also be tested to determine if it provides a better curve fit to the standard values.
- Determine the concentration of the target protein in the sample by interpolating the blank control subtracted absorbance values against the standard curve. Multiply the resulting value by the appropriate sample dilution factor, if used, to obtain the concentration of target protein in the sample.
- Samples generating absorbance values greater than that of the highest standard should be further diluted and reanalyzed. Similarly, samples which measure at absorbance values less than that of the lowest standard should be retested in a less dilute form.

#### TYPICAL DATA

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

| Standard Curve Measurements |            |       |             |
|-----------------------------|------------|-------|-------------|
| Concentration<br>(ng/mL)    | O.D 450 nm |       | Mean<br>O.D |
|                             | 1          | 2     |             |
| 0                           | 0.102      | 0.102 | 0.102       |
| 0.70                        | 0.135      | 0.133 | 0.134       |
| 1.41                        | 0.169      | 0.161 | 0.165       |
| 2.81                        | 0.255      | 0.234 | 0.245       |
| 5.63                        | 0.414      | 0.389 | 0.402       |
| 11.25                       | 0.779      | 0.728 | 0.753       |
| 22.5                        | 1.348      | 1.369 | 1.359       |
| 45                          | 2.537      | 2.439 | 2.488       |

Table 1. Example of human Albumin standard curve in Sample Diluent NS. The Albumin standard curve was prepared as described in the Standard Preparation section. The table shows raw data values.

#### TYPICAL SAMPLE VALUES

##### Sensitivity:

The calculated minimal detectable dose (MDD) is 0.1 ng/mL. The MDD was determined by calculating the mean of zero standard replicates (n=16) and adding 2 standard deviations then extrapolating the corresponding concentration.

##### Recovery

Three concentrations of Albumin were spiked in duplicate to the indicated biological matrix to evaluate signal recovery in the working range of the assay.

| Sample Type                   | Average % Recovery | Range (%) |
|-------------------------------|--------------------|-----------|
| 1:16,000,000 Serum            | 100                | 92 – 110  |
| 1:4,000,000 Plasma – EDTA     | 104                | 97 – 109  |
| 1:16,000,000 Plasma – Citrate | 101                | 100 – 103 |
| 1:16,000,000 Plasma – Heparin | 98                 | 89 - 105  |
| 7% Cell Culture Media*        | 104                | 104 - 105 |

\*Media is DMEM containing 10% fetal bovine serum.

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.

Native Albumin was measured in the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

| Dilution Factor | Interpolated value | 1:4,000,000 Human Serum | 1:8,000,000 Human Plasma (Citrate) | 1:4,000,000 Human Plasma (EDTA) | 1:8,000,000 Human Plasma (Heparin) |
|-----------------|--------------------|-------------------------|------------------------------------|---------------------------------|------------------------------------|
| Undiluted       | ng/mL              | 17.0                    | 9.3                                | 21.6                            | 10.7                               |
|                 | % Expected value   | 100                     | 100                                | 100                             | 100                                |
| 2               | ng/mL              | 7.9                     | 4.3                                | 9.4                             | 5.0                                |
|                 | % Expected value   | 93                      | 93                                 | 87                              | 93                                 |
| 4               | ng/mL              | 3.5                     | 2.1                                | 4.5                             | 2.5                                |
|                 | % Expected value   | 82                      | 90                                 | 83                              | 92                                 |
| 8               | ng/mL              | 1.9                     | 1.2                                | 2.2                             | 1.3                                |
|                 | % Expected value   | 89                      | 103                                | 81                              | 96                                 |
| 16              | ng/mL              | 1.1                     | NL                                 | 1.2                             | 0.7                                |
|                 | % Expected value   | 104                     | NL                                 | 89                              | 105                                |

NL – Non-linear

Recombinant Albumin was spiked into the following biological samples in a 2-fold dilution series. Sample dilutions are made in Sample Diluent NS.

| Dilution Factor | Interpolated value | 10% RPMI 1640 + 10% FBS |
|-----------------|--------------------|-------------------------|
| Undiluted       | ng/mL              | 21.4                    |
|                 | % Expected value   | 100                     |
| 2               | ng/mL              | 10.1                    |
|                 | % Expected value   | 94                      |
| 4               | ng/mL              | 5.1                     |
|                 | % Expected value   | 95                      |
| 8               | ng/mL              | 2.7                     |
|                 | % Expected value   | 98                      |
| 16              | ng/mL              | 1.3                     |
|                 | % Expected value   | 97                      |

Precision

Mean coefficient of variations of interpolated values of Albumin from three concentrations of human serum within the working range of the assay.

|        | Intra-assay | Inter-assay |
|--------|-------------|-------------|
| N=     | 8           | 3           |
| CV (%) | 5.2         | 7.6         |

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips: <https://www.abcam.com/en-us/technical-resources/guides/elisa-guide>

For technical support contact information, visit:  
<https://www.abcam.com/en-us/contact-us>  
<https://www.abcam.cn/contact-us> (China)  
<https://www.abcam.co.jp/contact-us> (Japan)

Copyright © 2026 Abcam, All Rights Reserved. All information / detail is correct at time of going to print.  
Version 5a | 2026-06-25

DIRECTIONS FOR 384-WELL PLATE FORMAT:

Materials Supplied for 384-well Format

| Item                                       | Quantity  | Storage Condition |
|--------------------------------------------|-----------|-------------------|
| 10X Human Albumin Capture Antibody         | 600 µL    | +4°C              |
| 10X Human Albumin Detector Antibody        | 600 µL    | +4°C              |
| Human Albumin Lyophilized Purified Protein | 2 Vials   | +4°C              |
| Antibody Diluent 4BI                       | 6 mL      | +4°C              |
| Sample Diluent NS                          | 250 mL    | +4°C              |
| Wash Buffer PT 10X                         | 2 x 20 mL | +4°C              |
| TMB Development Solution                   | 2 x12 mL  | +4°C              |
| Stop Solution                              | 2 x12 mL  | +4°C              |
| SimpleStep Pre-Coated 384-Well Microplate  | 384 wells | +4°C              |
| Plate Seal                                 | 1         | +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 or 600 nm in a 384-well plate.  
Deionized water.  
Multi- and single-channel pipettes.  
Tubes for standard dilution.  
Plate shaker for all incubation steps.  
Optional: Phenylmethylsulfonyl Fluoride (PMSF) (or other protease inhibitors).  
Optional: Automated liquid handler.

Reagent Preparation

Equilibrate all reagents to room temperature (18-25°C) prior to use. The kit contains enough reagents for one full plate. The sample volumes below are sufficient for running all 384 wells; adjust volumes as needed for the number of samples and dilution scheme for your experiment.

Prepare only as much reagent as is needed on the day of the experiment. Capture and Detector Antibodies have only been tested for stability in the provided 10X formulations.

**1X Wash Buffer PT:** Prepare 1X Wash Buffer PT by diluting Wash Buffer PT 10X with deionized water. To make 50 mL 1X Wash Buffer PT combine 5 mL Wash Buffer PT 10X with 45 mL deionized water. Mix thoroughly and gently.

**Antibody Cocktail:** Prepare Antibody Cocktail by diluting the capture and detector antibodies in Antibody Diluent 4BI. To make 6 mL of the Antibody Cocktail combine 600 µL 10X Capture Antibody and 600 µL 10X Detector Antibody with 4.8 mL Antibody Diluent 4BI. Mix thoroughly and gently.

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).

1. Reconstitute the Albumin standard sample by adding the volume of Sample Diluent NS indicated on the protein vial label. Hold at room temperature for 10 minutes. Mix thoroughly and gently. This is the 100 ng/mL **Stock Standard** Solution.
2. Label eight tubes, Standards 1– 8.
3. Add 110 µL of Sample Diluent NS into tube number 1 and 75 µL of Sample Diluent NS into numbers 2-8.
4. Use the **Stock Standard** to prepare the following dilution series. Standard #8 contains no protein and is the Blank control:

| Standard # | Dilution Sample | Volume to Dilute (µL) | Volume of Diluent (µL) | Starting Conc. (ng/mL) | Final Conc. (ng/mL) |
|------------|-----------------|-----------------------|------------------------|------------------------|---------------------|
| 1          | Stock Standard  | 90                    | 110                    | 100                    | 45                  |
| 2          | Standard#1      | 75                    | 75                     | 45                     | 22.5                |
| 3          | Standard#2      | 75                    | 75                     | 22.5                   | 11.25               |
| 4          | Standard#3      | 75                    | 75                     | 11.25                  | 5.63                |
| 5          | Standard#4      | 75                    | 75                     | 5.63                   | 2.81                |
| 6          | Standard#5      | 75                    | 75                     | 2.81                   | 1.41                |
| 7          | Standard#6      | 75                    | 75                     | 1.41                   | 0.70                |
| 8          | Blank Control   | 0                     | 75                     | 0                      | 0                   |

Plate Preparation

The 384-well plate included with this kit are supplied ready to use. It is not necessary to rinse the plate prior to adding reagents.  
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).  
Differences in well absorbance or “edge effects” have not been observed with this assay.

Assay Procedure for 384-well plate format

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. Prepare all reagents, working standards, and samples as directed in the previous sections.
- 2. Add 12.5 µL of all sample or standard to appropriate wells.
- 3. Add 12.5 µL of the Antibody Cocktail to each well.
- 4. Seal the plate and incubate for 1 hour at room temperature on a plate shaker set to 700 rpm.
- 5. Wash each well with 3 x 100 µL 1X Wash Buffer PT. Wash by aspirating or decanting from wells then dispensing 100 µL 1X Wash Buffer PT into each well. Wash Buffer PT should remain in wells for at least 10 seconds. Complete removal of liquid at each step is essential for good performance. After the last wash invert the plate and tap gently against clean paper towels to remove excess liquid.
- 6. Add 25 µL of TMB Development Solution to each well and incubate for 10 minutes in the dark on a plate shaker set to 700 rpm.  
*Given variability in laboratory environmental conditions, optimal incubation time may vary between 5 and 20 minutes.*  
Note: The addition of Stop Solution will change the color from blue to yellow and enhance the signal intensity about 3X. To avoid signal saturation, proceed to the next step before the high concentration of the standard reaches a blue color of O.D.600 equal to 1.0.
- 7. Add 25 µL of Stop Solution to each well. Shake plate on a plate shaker for 1 minute to mix. Record the OD at 450 nm. This is an endpoint reading. Proper mixing of the Stop Solution is required for proper measurement.
- 8. Alternative to 6 – 7: Instead of the endpoint reading at 450 nm, record the development of TMB Substrate kinetically. Immediately after addition of TMB Development Solution begin recording the blue color development with elapsed time in the microplate reader prepared with the following settings:

| Mode        | Kinetic                |
|-------------|------------------------|
| Wavelength: | 600 nm                 |
| Time:       | up to 20 min           |
| Interval:   | 20 sec - 1 min         |
| Shaking:    | Shake between readings |

**Note** that an endpoint reading can also be recorded at the completion of the kinetic read by adding 25 µL Stop Solution to each well and recording the OD at 450 nm.

Download our ELISA guide for technical hints, results, calculation, and troubleshooting tips:  
<https://www.abcam.com/en-us/technical-resources/guides/elisa-guide>

For technical support contact information, visit:  
<https://www.abcam.com/en-us/contact-us>  
<https://www.abcam.cn/contact-us> (China)  
<https://www.abcam.co.jp/contact-us> (Japan)