



PRODUCT INFORMATION & MANUAL

BCG Albumin Assay Kit (Colorimetric) *NBP3-24461*

For research use only.
Not for diagnostic or therapeutic
procedures.

www.novusbio.com - P: 303.730.1950 - P: 888.506.6887 - F: 303.730.1966 - technical@novusbio.com

Novus kits are guaranteed for 6 months from date of receipt

BCG Albumin Assay Kit (Colorimetric)

Catalog No: NBP3-24461

Method: Colorimetric method

Specification: 96T (Can detect 80 samples without duplication)

Instrument: Microplate reader

Sensitivity: 0.08 g/L

Detection range: 0.08-15 g/L

Average intra-assay CV (%): 1.5

Average inter-assay CV (%): 4.6

Average recovery rate (%): 95

- ▲ This kit is for research use only.
- ▲ Instructions should be followed strictly, changes of operation may result in unreliable results.
- ▲ Please kindly provide us the lot number (on the outside of the box) of the kit for more efficient service.

General information

▲ Intended use

This kit can be used to measure the albumin (ALB) content in serum, plasma, cell culture supernatant samples.

▲ Background

Albumin is the most abundant plasma protein, with a molecular weight of 66.5 kDa. Albumin is synthesized in the liver at a rate of 9 to 12 grams per day and regulated by insulin, amino acid intake and low colloidal osmotic pressure. Changes of albumin content in urine and serum are predictors of diabetic nephropathy, cardiovascular disease, liver disease and sepsis.

▲ Detection principle

Bromocresol green (BCG) is widely used as protein staining agent. BCG can combine the albumin in pH 4.0~4.2 to form an albumin-BCG complex. And the color changed from yellow to green. The depth of color is proportional to the concentration of albumin. The content of albumin in serum can be calculated indirectly by measuring the OD value at 630 nm.

▲ Kit components & storage

| Item | Component | Specification | Storage |
|---|----------------------------------|------------------|----------------------------------|
| Reagent 1 | Chromogenic Agent Stock Solution | 6 mL ×1 vial | 2-8°C , 12 months, shading light |
| Reagent 2 | 20 g/L Standard Solution | 1.2 mL × 2 vials | -20°C , 12 months |
| | Microplate | 96 wells | No requirement |
| | Plate Sealer | 2 pieces | |
| Note: The reagents must be stored strictly according to the preservation conditions in the above table. The reagents in different kits cannot be mixed with each other. | | | |

▲ Materials prepared by users

Instruments

Microplate reader(620-640 nm), Micropipettor, Vortex mixer

Reagents

Double distilled water, Normal saline (0.9% NaCl)

▲ Safety data

Some of the reagents in the kit contain dangerous substances. It should be avoided to touch the skin and clothing. Wash immediately with plenty of water if touching it carelessly. All the samples and waste material should be treated according to the relevant rules of laboratory's biosafety.

▲ Precautions

Before the experiment, please read the instructions carefully, and wear gloves and work clothes.

▲ The key points of the assay

1. Prevent the formulation of bubbles in the microplate.
2. Standard should be avoid repeated freezing and thawing.
3. Reagent 1 working solution should be stored with shading light.

Pre-assay preparation

▲ Reagent preparation

1. The preparation of reagent 1 working solution
Dilute the reagent 1 with double distilled water at a ratio of 1:4 and mix fully.
Prepare the fresh solution before use.
2. Take the reagent 2 from -20 °C and place on ice to thaw slowly. It is recommended to aliquot the reagent 2 to avoid repeated freezing and thawing.

▲ Sample preparation

The samples should be prepared as conventional methods. Also please refer to appendix II.

▲ Dilution of sample

It is recommended to take 2~3 samples with expected large difference to do pre-experiment before formal experiment and dilute the sample according to the result of the pre-experiment and the detection range (0.08-15 g/L).

The recommended dilution factor for different samples is as follows (for reference only)

| Sample type | Dilution factor |
|-------------------|-----------------|
| Human serum | 8-15 |
| Human plasma | 8-15 |
| HepG2 supernatant | 1 |
| Mouse plasma | 8-15 |
| Rat serum | 8-15 |

Note: The diluent is normal saline (0.9% NaCl) or PBS (0.01 M, pH 7.4).

Assay protocol

▲ Plate set up

| | 1 | 2 | 3 | 4 | 5 | 6 | 7 | 8 | 9 | 10 | 11 | 12 |
|---|---|---|----|-----|-----|-----|-----|-----|-----|-----|-----|-----|
| A | A | A | S1 | S9 | S17 | S25 | S33 | S41 | S49 | S57 | S65 | S73 |
| B | B | B | S2 | S10 | S18 | S26 | S34 | S42 | S50 | S58 | S66 | S74 |
| C | C | C | S3 | S11 | S19 | S27 | S35 | S43 | S51 | S59 | S67 | S75 |
| D | D | D | S4 | S12 | S20 | S28 | S36 | S44 | S52 | S60 | S68 | S76 |
| E | E | E | S5 | S13 | S21 | S29 | S37 | S45 | S53 | S61 | S69 | S77 |
| F | F | F | S6 | S14 | S22 | S30 | S38 | S46 | S54 | S62 | S70 | S78 |
| G | G | G | S7 | S15 | S23 | S31 | S39 | S47 | S55 | S63 | S71 | S79 |
| H | H | H | S8 | S16 | S24 | S32 | S40 | S48 | S56 | S64 | S72 | S80 |

Note: A, blank well; B-H, standard wells; S1-S80, sample wells.

▲ Detailed operation steps

The preparation of standard curve

Dilute 20 g/L standard solution with double distilled water to a serial concentration. The recommended dilution gradient is as follows: 0, 1, 2, 3.5, 5, 8, 12, 15 g/L. Reference is as follows:

| Number | Standard concentrations (g/L) | 20 g/L Standard solution (μL) | Double distilled water (μL) |
|--------|-------------------------------|-------------------------------|-----------------------------|
| A | 0 | 0 | 200 |
| B | 1 | 10 | 190 |
| C | 2 | 20 | 180 |
| D | 3.5 | 35 | 165 |
| E | 5 | 50 | 150 |
| F | 8 | 80 | 120 |
| G | 12 | 120 | 80 |
| H | 15 | 150 | 50 |

The measurement of samples

- (1) **Standard well:** add 10 μL standard with different concentration into the wells.
Sample tube: add 10 μL of sample into the wells.
- (2) Add 250 μL of the reagent 1 working solution to each well.
- (3) Stand for 10 min at room temperature.
- (4) Measure the OD value of each well at 630 nm with microplate reader.

▲ Summary operation table

| | Standard well | Sample well |
|---|---------------|-------------|
| BSA Standard solution with different concentrations (μL) | 10 | |
| Sample (μL) | | 10 |
| Reagent 1 working solution (μL) | 250 | 250 |
| Stand at room temperature for 10 min. Measure the OD values of each tube. | | |

▲ Calculation

Plot the standard curve by using OD value of standard and correspondent concentration as y-axis and x-axis respectively. Create the standard curve with graph software (or EXCEL). The concentration of the sample can be calculated according to the formula based on the OD value of sample. If the sample have been diluted, the concentration calculated from the standard curve must be multiplied by the dilution factor. The actual concentration is the calculated concentration multiplied by dilution factor.

The standard curve is: $y = ax + b$.

$$\text{ALB content (g/L)} = (\Delta A_{630} - b) \div a \times f$$

Note:

y: The absolute OD value of standard

x: The concentration of standard

a: The slope of standard curve

b: The intercept of standard curve

ΔA_{630} : $OD_{\text{Sample}} - OD_{\text{Blank}}$

f: Dilution factor of sample before test

Appendix I Data

▲ Example analysis

Dilute human serum with normal saline at a ratio of 1:9, take 10 μ L of diluted sample, carry the assay according to the operation table.

The results are as follows:

standard curve: $y = 0.0505x + 0.00779$, the average OD value of the sample is 0.364, the average OD value of the blank is 0.113, and the calculation result is:

ALB content (g/L) = $(0.364 - 0.113 - 0.0079) \div 0.0505 \times 10 = 48.75$ g/L

Appendix II Sample preparation

The following sample pretreatment methods are for reference only.

▲ Serum

Collect fresh blood and stand at 25°C for 30 min to clot the blood. Then centrifuge at 2000 g for 15 min at 4°C . Take the serum (which is the upper light yellow clarified liquid layer) to preserve it on ice for detection. If not detected on the same day, the serum can be stored at -80°C for a month.

▲ Plasma

Take fresh blood into the tube which has anticoagulant, centrifuge at 700-1000 g for 10 min at 4°C . Take the plasma (which is the upper light yellow clarified liquid layer, don't take white blood cells and platelets in the middle layer) to preserve it on ice for detection. If not detected on the same day, the plasma can be stored at -80°C for a month.

▲ Cell culture supernatant

Collect the cell culture supernatant, centrifuge at 10000 g for 10 min at 4°C ,and then take the supernatant to preserve it on ice for detection.

Appendix III References

1. Evans T W. Review article: albumin as a drug--biological effects of albumin unrelated to oncotic pressure. *Alimentary Pharmacology & Therapeutics*, 2002, 16(5): 6-11.
2. Arques S. Human serum albumin in cardiovascular diseases. *European Journal of Internal Medicine*, 2018, 52: 8-12.
3. Siotto M, Squitti R. Copper imbalance in Alzheimer's disease: Overview of the exchangeable copper component in plasma and the intriguing role albumin plays. *Coordination Chemistry Reviews*, 2018, 371: 86-95.
4. Hovind P, Tarnow L, Rossing P, et al. Predictors for the development of microalbuminuria and macroalbuminuria in patients with type 1 diabetes: inception cohort study. *Bmj British Medical Journal*, 2004, 328(7448): 1105-1108.