



# BCP Albumin Assay Kit

Catalog Number KA1613

250 assays

Version: 03

Intended for research use only

[www.abnova.com](http://www.abnova.com)

## Table of Contents

<b>Introduction .....</b>	<b>3</b>
Intended Use .....	3
Background .....	3
<b>General Information .....</b>	<b>4</b>
Materials Supplied .....	4
Storage Instruction .....	4
Materials Required but Not Supplied .....	4
Precautions for Use .....	4
<b>Assay Protocol .....</b>	<b>5</b>
Assay Procedure .....	5
<b>Data Analysis.....</b>	<b>6</b>
Calculation of Results.....	6
<b>Resources.....</b>	<b>7</b>
References .....	7

## Introduction

### Intended Use

Applications:

- ✓ Direct Assays: albumin in serum, urine, biological preparations.
- ✓ Drug Discovery/Pharmacology: effects of drugs on albumin metabolism.

Features:

- ✓ Sensitive and accurate: Use as little as 20  $\mu$ L samples. Detection range 0.3 g/dL (45 $\mu$ M) to 5 g/dL (750 $\mu$ M) albumin in 96-well plate assay.
- ✓ Simple and high-throughput: The procedure involves addition of a single working reagent and incubation for 5 min. Can be readily automated as a high-throughput assay for thousands of samples per day.
- ✓ Improved reagent stability and versatility: The optimized formulation has greatly enhanced reagent and signal stability. Cuvet or 96-well plate assay.
- ✓ Low interference in biological samples. No pretreatments are needed. Assays can be directly performed on raw biological samples i.e., in the presence of lipid and protein.

### Background

Albumin is the most abundant plasma protein in human. It accounts for about 60% of the total serum protein. Albumin plays important physiological roles, including maintenance of colloid osmotic pressure, binding of key substances such as long-chain fatty acids, bile acids, bilirubin, haematin, calcium, magnesium. It has anti-oxidant and anticoagulant effects, and also acts as a carrier for nutritional factors and drugs, as an effective plasma pH buffer. Serum albumin is a reliable prognostic indicator for morbidity and mortality, liver disease, nephritic syndrome, malnutrition and protein-losing enteropathies. High levels are associated with dehydration.

Simple, direct and automation-ready procedures for measuring albumin concentration in biological samples are becoming popular in Research and Drug Discovery. BCP Albumin Assay Kit is designed to measure albumin directly in biological samples without any pretreatment. The improved method utilizes bromcresol purple that forms a colored complex specifically with albumin. The intensity of the color, measured at 610nm, is directly proportional to the albumin concentration in the sample. The optimized formulation substantially reduces interference by substances in the raw samples.

## General Information

### **Materials Supplied**

List of component

Component	Amount
Reagent	50 mL
Albumin standard	2 mL 5 g/dL BSA

### **Storage Instruction**

Store Reagent and standard at 4°C and -20°C, respectively. Shelf life of at least 6 months (see expiry dates on labels).

### **Materials Required but Not Supplied**

Pipeting devices and accessories (e.g. 5 µL).

- Procedure using 96-well plate:

Clear bottom 96-well plates (e.g. Corning Costar) and plate reader.

- Procedure using cuvette:

Spectrophotometer and cuvetts for measuring OD at 610nm.

### **Precautions for Use**

Precautions:

Reagents are for research use only. Normal precautions for laboratory reagents should be exercised while using the reagents.

## Assay Protocol

### Assay Procedure

- Reagent Preparation:

Important: bring reagent to room temperature and shake before use.

- Procedure using 96-well plate:

1. Dilute standards in distilled water as follows. Dilute serum samples 2 fold. Transfer 20  $\mu$ L diluted standards and diluted samples to wells of a clear bottom plate. Store diluted standards at -20 °C for future use.

No	STD + H <sub>2</sub> O	Vol ( $\mu$ L)	BSA (g/dL)
1	100 $\mu$ L + 0 $\mu$ L	100	5.0
2	80 $\mu$ L + 20 $\mu$ L	100	4.0
3	60 $\mu$ L + 40 $\mu$ L	100	3.0
4	40 $\mu$ L + 60 $\mu$ L	100	2.0
5	30 $\mu$ L + 70 $\mu$ L	100	1.5
6	20 $\mu$ L + 80 $\mu$ L	100	1.0
7	10 $\mu$ L + 90 $\mu$ L	100	0.5
8	0 $\mu$ L + 100 $\mu$ L	100	0

2. Add 200  $\mu$ L working reagent and tap lightly to mix. Avoid bubble formation!
3. Incubate 5 min at room temperature and read optical density at 590-630nm (peak absorbance at 610nm).

- Procedure using cuvette:

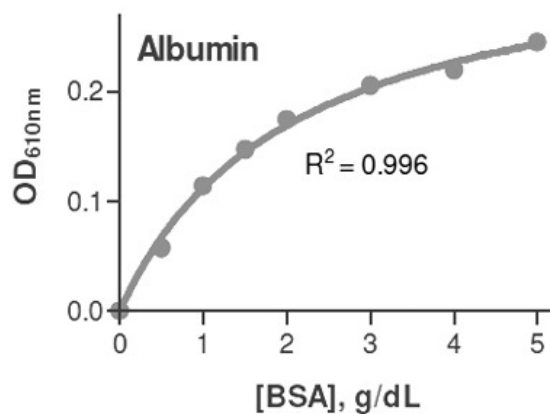
1. Transfer 60  $\mu$ L Blank, Standards and samples to appropriately labeled tubes. Add 1000  $\mu$ L working reagent and tap lightly to mix. Incubate 5 min at room temperature.
2. Transfer to cuvet and read optical density at 610nm.

Important: if sample OD is higher than the OD for standard, dilute samples in distilled water and repeat the assay.

## Data Analysis

### Calculation of Results

Albumin was assayed in duplicate using the 96-well assay protocol. The albumin content (g/dL) was  $3.6 \pm 0.0$ ,  $2.8 \pm 0.1$  and  $4.3 \pm 0.0$  in rat serum, fetal bovine serum and goat serum, respectively. The albumin content in a fresh human urine sample was below the detection limit (0.3 g/dL).



Standard Curve in 96-well plate assay

## Resources

### References

1. Maier SM et al (2007) Proteinuria of Nonautoimmune Origin in Wild-type FVB/NJ Mice. *Comp Med.* 57(3):255-66.
2. Sharifuzzaman SM and Austin B (2009). Kocuria SM1 controls vibriosis in rainbow trout (*Oncorhynchus mykiss*, Walbaum). *J Appl Microbiol.* 108(6): 2162 – 2170.
3. Shin SY et al (2009). Immunological investigation in the adenoid tissues from children with chronic rhinosinusitis. *Otolaryngol Head Neck Surg.* 141(1):91-6.