

QuantiQuik™ Albumin Quick Test Strips

Catalog Number: QQALBM10

DESCRIPTION

Albumin is the most abundant plasma protein in humans. Albumin plays many important physiological roles, including maintenance of colloid osmotic pressure and binding of key substances such as long-chain fatty acids, bile acids, bilirubin, haematin, calcium, and magnesium. Serum albumin is a reliable prognostic indicator for liver disease, nephritic syndrome, malnutrition, protein-losing enteropathies, as well as morbidity and mortality. In addition, elevated levels of serum albumin are associated with dehydration. Detection of albumin in urine is also crucial for diagnosis of patients with albuminuria disease.

BioAssay Systems' QuantiQuik™ Albumin Test Strips are based on the immobilization of an albumin sensitive chromagen onto a pad. Albumin will react with the chromagen and produce a colored product. The intensity of the product color is directly proportional to the albumin concentration in the sample. The QuantiQuik™ Albumin Test strips are a rapid and inexpensive method for semi-quantitative determination of albumin that does not require sophisticated laboratory instruments.

Product Information

Catalog No: QQALBM10

Number of Tests: 10 per package (other sizes available upon request).

Contents:

- Ten Test Strips
- Ten Dilution Tubes
- Instruction Manual

Shipping/Storage: The kit is shipped and stored at room temperature. Keep strips dry and out of direct sunlight.

Expiry: 6 months upon receipt.

Product Accessories

Some samples require a 100× dilution. Deionized or distilled water is needed for samples that require a dilution and can also be purchased separately. We offer the following:

- Ten Sample Dilution Tubes (containing 495 µL H₂O for 100× sample dilutions), Cat No. DT495

TEST PROCEDURE

Samples: For most serum or plasma samples we recommend diluting samples 100×. Most urine samples can be tested undiluted. For other samples, please contact Technical Support at info@bioassaysys.com for dilution recommendations.

1. For samples that do not require a dilution, carefully transfer 10 µL of sample directly onto the strip reaction pad. Skip ahead to step 4.
For samples requiring a 100× dilution, carefully transfer 5 µL of sample to an empty Dilution Tube. Then add 495 µL of dH₂O to the Dilution Tube.
2. Replace cap, securely close the vial, and invert the vial a couple of times to mix diluted samples. Alternatively, pipette up and down to mix.
3. Unscrew cap and transfer 10 µL of the diluted sample from the Dilution Tube directly onto the strip reaction pad.
4. Allow color to develop on strip for 30 seconds.
5. Compare the color of the reaction pad of the strip to the color on the provided Albumin Chart shown on the test strip bag. Multiply the concentration on the chart by the dilution used (i.e. 1 or 100) to determine the albumin concentration in the sample.