

## QuantiQuik™ Oxalate Quick Test Strips

Catalog Number: QQOXL10

### DESCRIPTION

Oxalate, or Oxalic Acid, is a metabolic breakdown product of the Krebs's Cycle in eukaryotes and the glyoxylate cycle in other microorganisms. Typically present in the urine of humans and other mammals, oxalate concentration can be used as a measure of kidney function. A high level of oxalate is an indicator for kidney stones, which are primarily made of the insoluble salt calcium oxalate. Measuring oxalate is more accurate than measuring calcium as a marker for kidney stones because calcium is excreted at high concentrations even in normal urine.

BioAssay Systems' QuantiQuik™ Oxalate Quick Test Strips are based on the enzymatic oxidation of oxalate coupled with a color reaction in one step. The intensity of product color is directly proportional to the oxalate concentration in the sample.

### Product Information

Catalog No: QQOXL10

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

Contents:

- Ten Test Strips
- Instruction Manual

Shipping/Storage: The kit is shipped at room temperature and should be stored in the refrigerator (4°C).

Expiry: 6 months upon receipt.

### Product Accessories

Urine samples require special treatment. Carbon Powder, EDTA, and a centrifuge or cutoff filters are required for treatment and can be purchased separately. We offer the following:

- Ten tubes containing 50 mg Carbon Powder, Cat. No. CARB-10
- 150 µL 0.5 M EDTA, Cat. No. EDTA-150
- Mini Centrifuge, Cat. No. CTFG-100
- Ten 5 kDa Cutoff Filters, Cat. No. CUTF-10

### SAMPLE PREPARATION

Samples: Whole blood samples are not compatible with the strips. Most serum and plasma samples can be tested undiluted and require no sample preparation. Urine samples require special treatment (please see below for more information). For other samples, please contact Technical Support at [info@bioassaysys.com](mailto:info@bioassaysys.com) for dilution recommendations.

### URINE

1. Take a tube containing 50 mg of carbon powder and use a pipetteman to add 240 µL of urine and 10 µL 0.5 M EDTA to the tube.
2. Place tube on a shaker for 5 min. (Note: If a shaker is not available the sample can be shaken by hand).
3. Using a pipetteman, carefully add the sample to a 5 kDa cutoff filter. Centrifuge for 2 minutes. Reserve the effluent for sample testing.

**Note:** If you do not wish to purchase cut off filters an alternative approach is available: Following step 2, centrifuge the tube for 10 minutes. Then, carefully transfer the supernatant to a clean tube. Centrifuge again for 5 min. Reserve the supernatant for sample testing.

### TEST PROCEDURE

1. Carefully add 10 µL of sample directly to the strip reaction pad.
2. Let color develop on strip for 5 minutes.
3. Compare the color of the reaction pad of the strip to the color on the provided Oxalate Chart shown on the test strip bag. If the sample was diluted, multiply the concentration on the chart by the dilution used (e.g. 1.04 for urine) to determine the oxalate concentration in the sample.