

QuantiQuik™ Malic Acid Quick Test Strips

Catalog Number: QQMALA10

DESCRIPTION

L-MALIC ACID, or L-malate, is a dicarboxylic acid that is made by all living organisms and plays an important role in the Calvin and Krebs Cycle. Malic acid is frequently used in food and beverage industries as an additive in products such as wine, beer, candies, etc. It can be used as an acidity regulator, preservative, or simply added for flavor. In the wine industry, decreasing levels of L-malic acid and increasing levels of L-Lactic Acid are monitored during malolactic fermentation. As L-malic acid is converted to L-lactic acid, the overall acidity of the wine is reduced which can lead to the improvement of the flavor of the wine.

BioAssay Systems' QuantiQuik™ L-Malic Acid Test Strips are based on L-malate dehydrogenase catalyzed oxidation of L-malic acid in which the formed NADH reduces a chromogenic reagent. The intensity of the product color is directly proportional to the L-malic acid concentration in the sample.

Product Information

Catalog No: QQMALA10

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

Contents:

- Ten Test Strips
- Ten Sample Development Tubes (400 μ L Development Reagent per tube)
- Instruction Manual

Shipping/Storage: The kit is shipped and stored at room temperature. Keep strips dry and out of direct sunlight. For long term storage (> 30 days), we recommend keeping the Sample Development Tubes at 4°C or below.

Expiry: 6 months upon receipt.

Product Accessories

Most samples require either a 2 \times , 5 \times , or 21 \times dilution. These dilutions can be performed either with a pipetteman if available or single use transfer pipettes can be purchased separately. We offer the following:

- Ten 20 μ L Transfer Pipettes (for 21 \times sample dilutions), Cat. No. TP20
- Ten 100 μ L Transfer Pipettes (for 5 \times sample dilutions), Cat. No. TP100
- Ten 400 μ L Transfer Pipettes (for 2 \times sample dilutions), Cat. No. TP400

TEST PROCEDURE

Samples: For white wine samples, especially those that have not gone through malolactic fermentation, we strongly recommend diluting samples 21 \times . Red wine and other acidic samples (fruit juice, beer, etc.) should be diluted 5 \times . Homogenized milk should be diluted 2 \times . Other samples such as yogurt require some extra treatment. Please see our website, www.bioassaysys.com, for the treatment instructions for these samples.

1. Unscrew the cap of one of the Sample Development tubes.
2. For samples requiring a 21 \times dilution, use a 20 μ L transfer pipette (a pipetteman can also be used if available), and carefully transfer 20 μ L of sample to the Sample Development Tube. For samples requiring a 5 \times dilution, use a 100 μ L transfer pipette and carefully transfer 100 μ L of sample to the Sample Development Tube. For samples requiring a 2 \times dilution, use a 400 μ L transfer pipette and carefully transfer 400 μ L of sample to the Sample Development Tube. (To use the transfer pipette: Squeeze top bulb of pipette and dip into sample and release bulb to take up sample. Next, place pipette tip into the Development tube and squeeze bulb again to release sample. *Important:* remove pipette from Sample Development Tube before releasing bulb).
3. Replace cap, securely close the vial and invert the vial a couple of times to mix diluted sample.
4. Unscrew cap and dip in one of the test strips making sure to fully submerge the yellow reaction pad at the end of the strip. Leave submerged for 5 seconds and then take out and shake a couple times to remove any drops clinging to strip.
5. Let color develop on strip for 5 minutes.
6. Compare the color of the reaction pad of the strip with the provided Malic Acid Chart. Multiply the Concentration in the chart by the dilution used (i.e. 2, 5, or 21).