

Surface Sampling Device

GENERAL DESCRIPTION

The LIGHTNING MVP ICON Surface Sampling Device is a self-contained device for use with the LIGHTNING MVP ICON™ or LIGHTNING MVP® systems for quality and HACCP-related monitoring of the environment. The MVP ICON and LIGHTNING MVP systems detect food and microbiological residues by measuring Adenosine triphosphate (ATP) using luciferase bioluminescence. ATP is found in most food residues and all bacteria, yeast and mold cells.

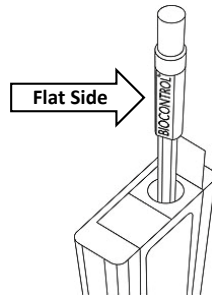
Part No. 64003-100 (100 tests)

KIT COMPONENTS

One box of 100 surface sampling devices

IMPORTANT USAGE GUIDELINES

It is critical to insert the sampling device into the MVP ICON with the two flat sides of the sampling device's base chamber facing left and right. Improper sampling device orientation may result in damage to the instrument.



Do not insert wet or contaminated sampling devices into the instrument. Wipe condensate from exterior of device before use with a clean, lint-free cloth. Avoid handling the device with wet or contaminated hands.

Always remove sampling devices from instrument after reading.

Do not store the MVP ICON or MVP instrument on its side or back. Avoid storing in humid areas where water leaks are possible.

TEST PROCEDURE

A. Collecting Sample

- Twist to remove sampling device from body tube. The swab tip must be moist. If fully saturated, any excess wetting agent should be blotted off by gently pressing the swab tip against the interior of the body tube. Be careful not to touch the swab or inside the sampling device with fingers.
- Thoroughly swab a standard 10 x 10 cm (4 x 4 in) area of interest for a typical flat surface. For irregularly shaped surfaces, ensure swabbing technique remains consistent each time you swab. Reinsert the swab into the body tube using a twisting motion to ensure it is sealed.

Note: After swabbing, the sampling device is stable for one hour prior to activation.

B. Activating Device and Reading Results

- To activate the sampling device, hold upright and push the plunger down until it is fully depressed. The reagents will be dispensed through the sampling device washing the sample off the swab tip and into the collection tube.
- Insert the sampling device into the MVP ICON or MVP and read the results. **When inserting into the MVP ICON, ensure the flat sides of the sampling device base chamber face left and right in the instrument.**

Note: Results should be read within two minutes of activation. Please refer to the MVP ICON User Guide or MVP System Guide for operating instructions.

INTERPRETING RESULTS

- When using default thresholds in the MVP ICON or MVP systems, readings of 2.5 Zones or less indicate the surface is considered clean.
- Readings between 2.6 - 3.0 Zones indicate a warning the surface is not adequately clean.
- If the reading is greater than 3.0 Zones, the surface is considered dirty and should be re-cleaned and re-tested.

CONTROLS

A. Positive Control

It is recommended LIGHTNING MVP ICON ATP Positive Controls (Part No. 64001-25) be run every other day to verify integrity of the reagent system.

- A positive control should yield a result from 3.1 to 4.3 Zones. See ATP Positive Control Directions for Use for additional information.

B. Negative Control

A negative control is an unused yet activated LIGHTNING MVP ICON Surface Sampling Device.

- Negative control results should read between 1.0 to 2.5 Zones.
- Readings above 2.5 may indicate contamination of the MVP ICON or MVP sampling chamber or sampling device.

See ATP Positive Control Directions for Use for additional information. Contact BioControl Technical Support at 800.245.0113 for more information.

PRECAUTIONS

The appearance of a yellow band on the swab tip is normal and will not affect performance.

If accidental activation prior to sampling occurs, do not use device.

Hold sampling device upright when activating.

Always insert the sampling device into the MVP ICON with the base chamber's two flat sides facing left-to-right.

Read sampling device within two minutes of activation.

When reading the sampling device, the MVP ICON or MVP must be kept upright (no greater than 45° angle). See MVP ICON User Guide or MVP System Guide for additional information.

STORAGE

Sampling devices should be stored at a temperature between 2 °C – 30 °C (36 °F – 86 °F). Protect from direct sunlight.

Unused devices must be resealed in the foil bags.

WARRANTY

BioControl Systems, Inc. (BCS) warrants this product to be free from defects in materials and workmanship, when stored under labeled conditions and used as intended until the expiration date stated on the package. BCS agrees during the applicable warranty period to replace all defective products after return to BCS. BCS shall not have obligation under this Limited Warranty to make replacements which result, in whole or in part, from negligence of the Buyer, or from improper use of the products, or use of the product in a manner for which it was not indicated. Buyer shall notify BCS of any products which it believes to be defective during the warranty period. At BCS option, such products shall be returned to BCS, transportation and insurance prepaid. BCS shall replace any such product found to be defective, at no charge. Should BCS examination not disclose any defect covered by the foregoing warranty, BCS shall so advise Buyers and dispose of the product in accordance with Buyer's instructions.

©BioControl Systems, Inc. 2013. Printed in USA. LIGHTNING MVP ICON is a trademark of BioControl, pending registration. LIGHTNING MVP is a registered trademark of BioControl. US patents No. 7,399,984; 7,544,961; 7,030,403; 6924498. Australian patents 2002357107; 2003231770. New Zealand patents 533051, 535819. Japanese patents 4,467,304; 4,431,034. Mexico patents 259897; 246269; 272356.

All rights reserved.

IF YOU REQUIRE MORE INFORMATION ABOUT THE LIGHTNING MVP ICON SURFACE SAMPLING DEVICE, ITS USE, OR OTHER BIOCONTROL PRODUCTS, PLEASE CONTACT: