

Temperature Probe

GENERAL DESCRIPTION

The MVP ICON® Temperature Probe is a thermistor housed in a stainless steel probe designed for general purpose temperature measurement of liquids or semi-solids ranging from -55 to 125 °C. The MVP ICON Temperature Probe is designed for use with the MVP ICON instrument only. Any other combination might cause loss of performance or irreversible damage to both probe and instrument.

Part No. 78087-00

KIT COMPONENTS

Each kit contains one MVP ICON Temperature Probe.

TEST PROCEDURE

A. Using the MVP ICON Temperature Probe

- Ensure the Temperature Probe is securely connected to the MVP ICON instrument by inserting the probe's connector with the arrows facing toward the back of the instrument and the white circular sticker facing the front. Insert probe into the sample. Follow instructions on the MVP ICON to complete the reading. Results will display in either °C or °F based on instrument settings.

Note: To disconnect the probe, squeeze both sides of the connector while unplugging from the MVP ICON.

- When using the MVP ICON Temperature Probe with semi-solids, insert probe to desired depth then rotate left to right several times and tilt to ensure sample contact.
- If the sample temperature is fluctuating, the MVP ICON Temperature Probe will need to stabilize before a reading is given.

B. Calibration

The MVP ICON Temperature Probe features a single-point calibration that can be done at any known temperature within the operating range.

Note: Calibration should be performed prior to first initial use. The calibration is most accurate when using a known temperature close to the sample temperature range.

- From MVP ICON, enter **TEMPERATURE** mode, press **Device** menu, then press **Calibration**.
- Insert probe into a liquid of known temperature, enter value into the **Reference Value** field, then press **Calibrate**.

- The date of last calibration will appear on the MVP ICON Home screen.

STORAGE

Clean probe thoroughly with water and a mild detergent solution. Rinse with deionized water. Store probe in original packaging when not in use.

PRECAUTIONS

- The MVP ICON Temperature Probe is rugged and durable and requires little maintenance. To ensure lasting performance read and follow all operating guidelines.
- Avoid prolonged exposure within samples at either end of the specified temperature range. Do not use the probe outside the specified temperature range (see **SPECIFICATIONS**) as this might result in probe failure or irreversible damage.
- Samples must be liquids or semi-solids.

SPECIFICATIONS

Sensor:	Thermistor
Temperature range	-55 to 125 °C
Resolution	0.1 °C
Accuracy	± 0.5 °C

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, for 6 months from date of purchase. 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 product which it believes to be defective during the warranty period. At BCS option, such product 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. 2016. Printed in USA.

LIGHTNING MVP and MVP ICON are trademarks of BioControl Systems, Inc., US patents No. 7,399,984; 7,544,961; 7,030,403; 6,924,498. 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.