

User Guide

MilliSentials™ Aliquoting Pipette Controller User Calibration

MISEPTCON

Introduction

If accuracy readings are determined to be outside of acceptable levels using the Gravimetric Verification Procedure, the MilliSentials™ Aliquoting Pipette Controller may be re-calibrated or reset to factory specification by the end user using the procedures outlined here.

Recalibrating Procedure

The Gravimetric Verification Procedure should be done prior to and after recalibrating or resetting to factory calibration.

NOTE: After entered, calibration mode may be exited (without saving changes) by quickly pressing and releasing the aliquot button. This will be followed by the Status LED flashing **RED** five (5) times to confirm exit.

Enter Calibration Mode:

1. Set the volume dial to BELOW 0.0mL.
2. Hold the MilliSentials™ Aliquoting Pipette Controller upside down. Maintain this orientation and hold the aliquot button for five (5) seconds. Three tones will sound to indicate entry into calibration mode and the Status LED will begin quickly blinking **PURPLE**. Release the trigger and return the unit to normal orientation. Watch for the Status LED to blink **PURPLE** before proceeding.

Enter accuracy error values:

1. Once the Status LED is flashing quickly, it is ready to accept the 0.5mL error number for Accuracy (%) calculated after the Gravimetric Verification Procedure.
2. Set the volume dial to the number obtained (for example, if the mean error is 1.5%, then set the aliquot volume dial to 1.5mL)
 - a. If the MilliSentials™ Aliquoting Pipette Controller dispensing was high, (for example, Accuracy value of +1.5%), then press the aspirate button to enter the number.
 - b. If the MilliSentials™ Aliquoting Pipette Controller dispensing was low, (for example, Accuracy value of -1.5%), then press the dispense button to enter the number.

3. The Battery and Status LEDs will alternate **PURPLE** five (5) times, accompanied by a tone, to confirm entry.
4. The Status LED will now be slowly blinking **PURPLE**, indicating that the unit is ready to accept the calculated error at 5 mL.
5. Repeat Step 2 to enter this value for 5.0 mL and confirm entry.
6. The Battery and Status LEDs will alternate **PURPLE** five (5) times, accompanied by a tone, to confirm entry.
7. If the Battery and Status LEDs blink **GREEN**, the unit has been successfully calibrated. The unit will now return to normal operation.

NOTE: A post-calibration Gravimetric Verification should always be performed to confirm the user calibration was successful and that the unit demonstrates acceptable performance.

Reset to Factory Calibration

The MilliSentials™ Aliquoting Pipette Controller has been calibrated from the factory to $\pm 3\%$ at 5.0 mL and $\pm 4\%$ for 0.5 mL. However, the MilliSentials™ Aliquoting Pipette Controller can easily be reset to the factory calibration at any time.

1. Set the volume dial to BELOW 0.0mL.
2. Hold the MilliSentials™ Aliquoting Pipette Controller upside down. Maintain this orientation and hold the aliquot button for five (5) seconds. The Status LED will now be slowly blinking **PURPLE**. Release the trigger and return the unit to normal orientation.
3. Press and hold the aliquot button for approximately five (5) seconds, until the Battery and Status LEDs begin to blink **GREEN**. Release the button.
4. The Battery and Status LEDs will blink **GREEN** five (5) times to confirm that a factory reset has been applied.

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page on our web site at SigmaAldrich.com/TechService.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/Terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/Offices.

Merck, Millipore, Millisentials, and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.
© 2020-2023 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

The life science business of Merck operates as MilliporeSigma in the U.S. and Canada.