

PDM3700 Quick Start Guide

Start a Sample

From WinPDM

1. Connect the PDM3700 to the charger. Connect the charger to your PC.
2. Start WinPDM.
3. Select the “Program Shift” button.
4. Enter the start date and time of the sample run.
5. Enter the sample run duration.
6. Select the sample type.
7. Determine the expected shift start and shift end ambient temperatures. Select the ambient temperature range from the WinPDM Shift Programming window so that the shift start and end temperatures fall within this programmed range.
8. Make sure all available fields are entered.
9. Select “OK.” The PDM3700 unit will warm up for 35 minutes before the sample run is scheduled to begin.

From the PDM3700 Unit

1. Press the “A” (“WAKE UP”) button to activate the LED display.
2. Press the “A” button and the “B” (“SCROLL”) button simultaneously to display the Start Sampling screen.
3. Press the “A” button to display the Start Sample screen.
4. Press the “B” button to adjust the sample time duration.
5. Press the “A” button to begin an immediate sample run. The Warming screen will now display. After the 35 minute warmup period has passed, the PDM3700 unit will begin the sample run, and the First Sample Screen #1 will display.

Stop a Sample

From WinPDM

1. Connect the PDM3700 to the charger. Connect the charger to your PC.
2. Start WinPDM.
3. Select the “Clear Program” button.

From the PDM3700 Unit

1. Press the “A” button and the “B” (“SCROLL”) button simultaneously to display the Reset Sample 2 screen.
2. Press the “B” button to display the Stop Sampling screen.
3. Press the “A” button to stop the immediate sample run and display the Sampling Complete Screen #1. This action will stop both the first and second immediate sample run (if a second immediate sample run has been started).

Note If starting a sample run from WinPDM, the programmed sample run can be stopped only by using the WinPDM software program. ▲

Note Samples collected for MSHA compliance must be programmed and started using WinPDM. ▲

PDM3700 Operating Notes

Programming a sample for MSHA reporting:

WinPDM must be used to program a sample to meet MSHA sampling and reporting requirements. Samples may be programmed to run from the instrument keypad, but may not meet MSHA requirements.

WinPDM version 7.52 or higher must be used for compliance with MSHA requirements for sampling and data submission. While earlier versions of WinPDM may operate with the PDM3700, they do not meet MSHA requirements. Always verify that the latest approved version of WinPDM is being used.

Sample Tubing Length

The PDM3700 is supplied with tubing of extra length (a 38 ± 0.5" tube or 50' roll) and must be trimmed before first use by the operator to meet agency requirements (35–37"). The tubing must be checked prior to use each day and may be trimmed due to breakage or fraying, but must always be between 35–37" during use. If the tubing is shorter than 35" it must be replaced. To measure the tubing, remove from the instrument and lay on a flat surface. Straighten the tubing and measure the total length without stretching the overall tubing length.

Installing the Tubing onto the PDM3700 Sample Inlet

When installing the tubing onto the inlet of the PDM3700, do not push the tubing completely to the shoulder of the inlet. Leave a small space to allow the use of the tubing removal tool. Care must be taken when installing the sample tube onto the inlet to prevent cracking, gouging, or otherwise damaging the tubing.

Removing the Tubing from the Sample Inlet

Create a small gap between the tubing end and inlet shoulder, if one isn't already present. Slip the tubing removal tool onto the sample inlet just above the sample tubing. Using a straight alignment, pull the tool away from the inlet bracket and pull the tubing off the inlet. Do NOT use the tubing removal tool on the instrument cyclone, since this may damage the plastic cyclone.

