

Technical Bulletin

48i Multi-Point Calibration

Bulletin# 48i-MPC

Rev. 09/2014

The multi-point calibrations are divided into three calibration points consisting of:

- Cal point 1
- Cal point 2
- Cal point 3

The range is user definable and the following calibration points are recommended:

- Cal point 1: 80% of range
- Cal point 2: 50% of range
- Cal point 3: 20% of range

Note: If performing a multi-point calibration after a bad calibration or changing ranges, it is recommended to start off with default values for the Cal point coefficients.

Main Menu > Service Menu > Multi-Point Cal > Default Coef

To calibrate Cal points 1, 2, & 3:

1. Introduce a CO concentration of 80% of the upper range limit (URL) to the analyzer. The total flow must exceed the total demand of the analyzer connected to the output manifold to ensure that no ambient air is pulled into the manifold vent.
2. Allow the analyzer to sample the CO concentration until a stable response is obtained.
3. Main Menu > Service > Multi-Point Cal > Calibrate Point 1.
4. Using the arrow keys, enter the concentration value of Cal Gas 1 (third line).
5. Press the Menu button to back up a step in the Service Menu.
6. Introduce a CO concentration 50% of the URL.
7. Select Cal Point 2.
8. Using the arrow keys, enter the concentration value of Cal Gas 2 (third line).
9. Press the Menu button to back up a step in the Service Menu.
10. Introduce a CO concentration 20% of the URL.
11. Select Cal Point 3.
12. Using the arrow keys, enter the concentration value of Cal Gas 3 (third line).
13. Press the Menu button to back up a step in the Service Menu.
14. Select Coefficients.
15. Press the Enter button, the instrument will automatically calculate and save the new coefficients.