Instruction Sheet

Thermo Scientific AquaSensors™ RDO® Pro Sensor Calibration Kit, Cat. No. RDK1

Kit Contents

- Calibration and storage chamber
- Calibration cap
- Storage cap
- Sodium sulfite

1 Point Calibration

Water-saturated Air

1. Remove the storage cap from the top of calibration chamber and replace it with the calibration cap (cap with the vent hole).





Calibration Cap

Storage Cap

- 2. Fill the calibration chamber to the lower fill line with approximately 10 mL of water.
 - **Note:** The water temperature should be similar to the sample temperature.
- 3. Gently dry the sensor and foil with a paper towel, making sure that no water is on the body of the sensor and that the sensing foil is free of water and fouling.
- 4. Place the sensor in the calibration chamber so that the sensing foil is about 1 inch above the surface of the water.

Note: Make sure that there are no water droplets on the surface of the foil when it is in the calibration chamber.

5. Allow 5 to 10 minutes for temperature stabilization prior to calibration. Do not leave the sensor in the calibration chamber for more than 30 minutes. This can allow condensation to form on the surface of the foil, giving false low readings after calibration. If condensation does occur, remove the sensor, dry the foil, place the sensor in the chamber and calibrate.

Lower fill line





2 Point Calibration

100% Calibration Point

Perform a water-saturated air calibration as described above.

0% Calibration Point

1. Place the calibration cap (cap with the vent hole) onto the top of the calibration chamber.



Calibration Cap

- 2. Fill the calibration chamber to the upper fill line with approximately 60 mL of fresh sodium sulfite solution.
- 3. Completely submerge the foil of the sensor, making sure that the temperature thermistor is also completely submerged under the surface of the solution. Do not allow the foil to rest completely on the bottom of the calibration chamber. Leave at least ½ inch between the surface of the foil and the bottom of the chamber.
- 4. Allow at least five minutes for the temperature to stabilize prior to calibration.
- 5. Once calibrated, remove the sensor and thoroughly rinse it to remove any excess sodium sulfite.

— Upper fill line



©2009 Thermo Fisher Scientific Inc. All rights reserved. RDO is a registered trademark of In-Situ Inc., Fort Collins, CO USA. All other trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.



