

# User Guide

Cadmium  
Ion Selective  
Electrode



ROSS and the COIL trade dress are trademarks of Thermo Fisher Scientific Inc.

AQUAfast, Cahn, ionplus, KNipHE, No Cal, ORION, perpHect, PerpHecT, PerpHecTion, pHISA, pHuture, Pure Water, Sage, Sensing the Future, SensorLink, ROSS, ROSS Ultra, Sure-Flow, Titrator PLUS and TURBO2 are registered trademarks of Thermo Fisher.

1-888-pHAX-ION, A+, All in One, Aplus, AQUAsnap, AssuredAccuracy, AUTO-BAR, AUTO-CAL, AUTO DISPENSER, Auto-ID, AUTO-LOG, AUTO-READ, AUTO-STIR, Auto-Test, BOD AutoEZ, Cable-Free, CERTI-CAL, CISA, DataCOLLECT, DataPLUS, digital LogR, DirectCal, DuraProbe, Environmental Product Authority, Extra Easy/Extra Value, FAST QC, GAP, GLPcal, GLPcheck, GLPdoc, ISEasy, KAP, LabConnect, LogR, Low Maintenance Triode, Minimum Stir Requirement, MSR, NISS, One-Touch, One-Touch Calibration, One-Touch Measurement, Optimum Results, Orion Star, Pentrode, pHuture MMS, pHuture Pentrode, pHuture Quatrode, pHuture Triode, Quatrode, QuiKcheK, rf link, ROSS Resolution, SAOB, SMART AVERAGING, Smart CheK, SMART STABILITY, Stacked, Star Navigator 21, Stat Face, The Enhanced Lab, ThermaSense, Triode, TRIUMpH, Unbreakable pH, Universal Access are trademarks of Thermo Fisher.

Guaranteed Success and The Technical Edge are service marks of Thermo Fisher.

PerpHecT meters are protected by U.S. patent 6,168,707.

PerpHecT ROSS are protected by U.S. patent 6,168,707.

ORION Series A meters and 900A printer are protected by U.S. patents 5,198,093, D334,208 and D346,753.

ionplus electrodes and Optimum Results solutions are protected by US Patent 5,830,338.

ROSS Ultra electrodes are protected by US patents 6,793,787.

Orion ORP Standard is protected by US Patent 6,350,367.

Orion NoCal electrodes are protected by US Patent 7,276,142.

© 2008 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific Inc. and its subsidiaries.

The specifications, descriptions, drawings, ordering information and part numbers within this document are subject to change without notice.

This publication supersedes all previous publications on this subject.

# Table of Contents

<b>Introduction</b>	<b>1</b>
<b>Required Equipment</b>	<b>2</b>
<b>Serial Dilutions</b>	<b>3</b>
<b>Electrode Setup</b>	<b>4</b>
Electrode Preparation	4
Checking Electrode Operation (Slope)	6
Measurement Units	7
Sample Requirements	7
Measuring Hints	8
Electrode Storage	9
Electrode Maintenance	10
<b>Analytical Techniques</b>	<b>12</b>
Direct Calibration Technique	14
Small Volume Direct Calibration Technique	18
Low Level Calibration Technique	22
Known Addition Technique	24
Cadmium Titration Technique	32
<b>Electrode Characteristics</b>	<b>35</b>
Electrode Response	35
Reproducibility	35
Temperature Effects	36
Interferences	37
pH Effects	38
Complexation	39
Theory of Operation	40
<b>Troubleshooting</b>	<b>42</b>
Assistance	43
Warranty	43
Troubleshooting Checklist	44
<b>Ordering Information</b>	<b>45</b>
<b>Specifications</b>	<b>46</b>



# Introduction

This user guide contains information on the preparation, operation and maintenance for the cadmium ion selective electrode (ISE). General analytical procedures, electrode characteristics and electrode theory are also included in this user guide. Cadmium electrodes measure free cadmium ions in aqueous solutions quickly, simply, accurately and economically.

Technical Support Chemists can be consulted for assistance and troubleshooting advice. Within the United States call 1.800.225.1480 and outside the United States call 978.232.6000 or fax 978.232.6031. In Europe, the Middle East and Africa, contact your local authorized dealer. For the most current contact information, visit [www.thermo.com/contactwater](http://www.thermo.com/contactwater).

For the latest application and technical resources for Thermo Scientific Orion products, visit [www.thermo.com/waterapps](http://www.thermo.com/waterapps).

## **Cadmium ionplus® Sure-Flow® Solid State Combination ISE, Cat. No. 9648BNWP**

The cadmium combination electrode has the sensing and reference half-cells built into one electrode, which decreases the amount of required solutions and reduces waste. The built-in Sure-Flow reference junction prevents electrode clogging and provides fast and stable readings. The cadmium combination electrode is available with a waterproof BNC connector, Cat. No. 9648BNWP. Electrodes with a waterproof BNC connector can be used on any ISE or mV meter with a BNC connection.

## **Cadmium Solid State Half-Cell ISE, Cat. No. 9448BN and 9448SC**

The cadmium half-cell electrode must be used with the double junction reference electrode, Cat. No. 900200. The cadmium half-cell electrode is available with a BNC connector, Cat. No. 9448BN, and a screw cap connector, Cat. No. 9448SC. Electrodes with a screw cap connector require a separate cable.

# Required Equipment

1. Thermo Scientific Orion ISE meter, such as the 4-Star pH/ISE meter or 5-Star pH/ISE/DO/conductivity meter; equivalent ISE meter; or mV meter with a 0.1 mV resolution.

Cadmium electrodes can be used on any ISE or mV meter with a BNC connection. The electrodes can also be used on meters with a variety of inputs when an adapter cable is used. Visit [www.thermo.com/water](http://www.thermo.com/water) for details.

2. Thermo Scientific Orion cadmium electrode.

The 9448BN and 9448SC cadmium half-cell electrodes require a separate reference electrode, Cat. No. 900200.

3. Magnetic stirrer or Thermo Scientific Orion stirrer probe, Cat. No. 096019. The stirrer probe can be used with 3-Star, 4-Star and 5-Star benchtop meters.
4. Volumetric flasks, graduated cylinders and beakers. Plastic labware is required for low level cadmium analysis.
5. Distilled or deionized water.
6. Cadmium electrode filling solution.

Use Optimum Results™ A filling solution, Cat. No. 900061, for the 9648BNWP cadmium combination electrode.

Use inner chamber filling solution, Cat. No. 900002, and outer chamber filling solution, Cat. No. 900003, for the double junction reference electrode that is used with the 9448BN and 9448SC cadmium half-cell electrodes.

7. 0.1 M or 1000 ppm cadmium calibration standards.

0.1 M  $\text{Cd}(\text{NO}_3)_2 \cdot 4\text{H}_2\text{O}$  solution – In a 1 liter flask, place 30.8 grams of reagent-grade cadmium nitrate. Dissolve the solid and dilute to volume with distilled water.

1000 ppm cadmium solution – Weigh out 2.74 grams of reagent-grade cadmium nitrate and add to a 1 liter volumetric flask. Dissolve the solid and dilute to volume with distilled water.

8. Cadmium ionic strength adjuster (ISA), Cat. No. 940011. ISA provides a constant background ionic strength for samples and standards.

# Serial Dilutions

Serial dilution is the best method for the preparation of standards. Serial dilution means that an initial standard is diluted, using volumetric glassware, to prepare a second standard solution. The second standard is similarly diluted to prepare a third standard, and so on, until the desired range of standards has been prepared.

1. **To prepare a  $10^{-2}$  M standard (1124 ppm cadmium) –**  
Pipet 10 mL of the 0.1 M standard into a 100 mL volumetric flask. Dilute to the mark with deionized water and mix well.
2. **To prepare a  $10^{-3}$  M standard (112.4 ppm cadmium) –**  
Pipet 10 mL of the  $10^{-2}$  M standard into a 100 mL volumetric flask. Dilute to the mark with deionized water and mix well.
3. **To prepare a  $10^{-4}$  M standard (11.24 ppm cadmium) –**  
Pipet 10 mL of the  $10^{-3}$  M standard into a 100 mL volumetric flask. Dilute to the mark with deionized water and mix well.

To prepare standards with a different concentration use the following formula:

$$C_1 * V_1 = C_2 * V_2$$

$C_1$  = concentration of original standard

$V_1$  = volume of original standard

$C_2$  = concentration of standard after dilution

$V_2$  = volume of standard after dilution

For example, to prepare 100 mL of a 100 ppm cadmium standard from a 11240 ppm cadmium standard:

$$C_1 = 11240 \text{ ppm cadmium}$$

$$V_1 = \text{unknown}$$

$$C_2 = 100 \text{ ppm cadmium}$$

$$V_2 = 100 \text{ mL}$$

$$11240 \text{ ppm} * V_1 = 100 \text{ ppm} * 100 \text{ mL}$$

$$V_1 = (100 \text{ ppm} * 100 \text{ mL}) / 11240 \text{ ppm} = 0.9 \text{ mL}$$

# Electrode Setup

## Electrode Preparation

**9448BN and 9448SC Cadmium Half-Cell Electrode** – Remove the protective shipping cap from the sensing element and save the cap for storage.

**900200 Double Junction Reference Electrode** – Prepare the reference electrode according to the reference electrode user guide. Fill the reference electrode with inner chamber filling solution, Cat. No. 900002, and outer chamber filling solution, Cat. No. 900003.

**9648BNWP Cadmium Combination Electrode** – Remove the protective shipping cap from the sensing element and save the cap for storage. Fill the electrode with Optimum Results A filling solution, Cat. No. 900061.

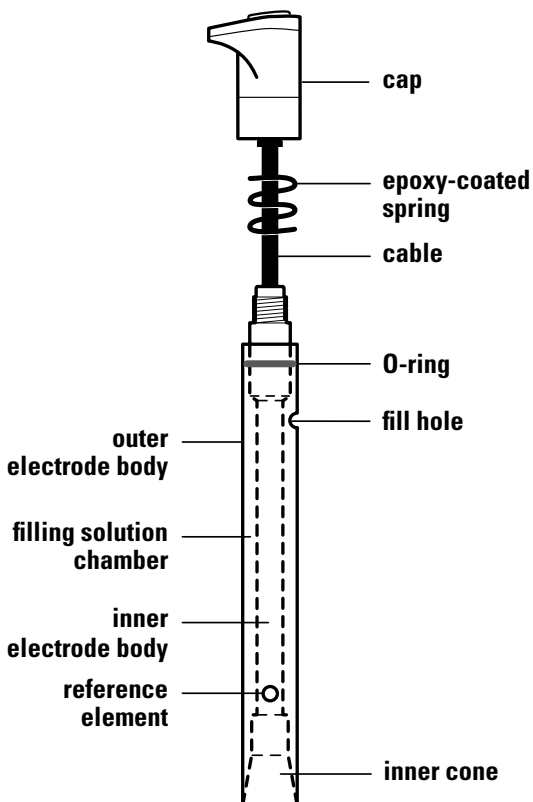
## 9648BNWP Cadmium Combination Electrode Filling Instructions

1. Lift the flip spout on the filling solution bottle to a vertical position.
2. Insert the spout into the filling hole on the outer body of the electrode and add a small amount of filling solution to the reference chamber. Invert the electrode to moisten the top O-ring and then return the electrode to the upright position.
3. Hold the electrode body with one hand and use your thumb to push down on the electrode cap to allow a few drops of filling solution to drain out of the electrode.
4. Release the electrode cap. If the sleeve does not return to its original position, check if the O-ring is moist and repeat steps 2 through 4 until the sleeve returns to the original position.
5. Add filling solution to the electrode up to the filling hole.

**Note:** Add filling solution each day before using the electrode. The filling solution level should be at least one inch above the level of sample in the beaker to ensure a proper flow rate. The fill hole should always be open when taking measurements.



**Figure 1**  
**9648BNWP Cadmium Combination Electrode**



## Checking Electrode Operation (Slope)

These are general instructions that can be used with most meters to check the electrode operation. Refer to the meter user guide for more specific information.

This procedure measures electrode slope. Slope is defined as the change in millivolts observed with every tenfold change in concentration. Obtaining the slope value provides the best means for checking electrode operation.

1. If the electrode has been stored dry, prepare the electrode as described in the **Electrode Preparation** section.
2. Connect the electrode to a meter with a mV mode. Set the meter to the mV mode.
3. Add 100 mL of distilled water and 2 mL of ISA into a 150 mL beaker. Stir the solution thoroughly.
4. Rinse the electrode with distilled water and place the electrode into the solution prepared in step 3.
5. Select either a 0.1 M or 1000 ppm cadmium standard. Pipet 1 mL of the standard into the beaker and stir the solution thoroughly. When a stable reading is displayed, record the electrode potential in millivolts.
6. Pipet 10 mL of the same standard into the same beaker and stir the solution thoroughly. When a stable reading is displayed, record the electrode potential in millivolts.
7. There should be a 25 to 30 mV difference between the two millivolt readings when the solution temperature is between 20 to 25 °C. If the millivolt potential is not within this range, refer to the **Troubleshooting** section.

# Measurement Units

Cadmium concentration can be measured in moles per liter (M), parts per million (ppm) or any convenient concentration unit.

**Table 1**  
**Concentration Unit Conversion Factors**

Moles/Liter (M)	ppm
1.0	112400
$10^{-1}$	11240
$10^{-2}$	1124
$8.9 \times 10^{-3}$	1000
$10^{-3}$	112.4
$10^{-4}$	11.24
$8.9 \times 10^{-6}$	1

## Sample Requirements

The epoxy body of the cadmium electrode is resistant to damage by aqueous solutions. The electrode may be used intermittently in solutions that contain methanol, benzene or acetone. Contact Technical Support for information on using the electrode for specific applications.

Samples and standards should be at the same temperature. A 1 °C difference in temperature for a  $10^{-3}$  M cadmium solution will give rise to about a 4% error.

The solution temperature must be less than 80 °C.

Cadmium samples must be below pH 7 to avoid precipitation of  $\text{Cd}(\text{OH})_2$ . Acidify cadmium samples with 1 M  $\text{HNO}_3$  if necessary, but the hydrogen ion will interfere if the pH is too low. See the **pH Effects** section to determine the optimum pH working range for your sample.

In all analytical procedures, ISA must be added to all samples and standards before measurements are taken.

## Measuring Hints

- Stir all standards and samples at a uniform, moderate rate. Place a piece of insulating material, such as Styrofoam or cardboard, between the magnetic stir plate and beaker to prevent measurement errors from the transfer of heat to the sample.
- Always use freshly prepared standards for calibration.
- Always rinse the electrode with distilled water between measurements and shake the electrode to remove the water and prevent sample carryover. Do not wipe or rub the electrode sensing element.
- Allow all standards and samples to reach the same temperature for precise measurements.
- Concentrated samples (greater than  $10^{-1}$  M cadmium) should be diluted before measurement.
- Verify the electrode calibration every two hours by placing the electrode in a fresh aliquot of the least concentrated standard used for calibration. If the value has changed by more than 2%, recalibrate the electrode.
- After immersing the electrode in a solution, check the electrode sensing surface for air bubbles and remove air bubbles by reimmersing the electrode in the solution and gently tapping it.
- For high ionic strength samples, prepare standards with a background composition similar to the sample.
- The fill hole cover must be open during measurements to ensure a uniform flow of filling solution.
- If the combination electrode is used and the electrode is used in dirty or viscous samples or the electrode response becomes sluggish, empty the electrode completely, hold the junction open and flush the junction with distilled water. Empty any water from the electrode and refill it with fresh filling solution. Press down on the electrode cap to let a few drops of the filling solution flow out of the electrode and then replenish any lost solution.

# Electrode Storage

## **Cadmium Half-Cell Electrode Storage, Cat. No. 9448BN and 9448SC**

The cadmium half-cell electrode should be rinsed thoroughly with distilled water and stored dry in the air at all times. When storing the electrode for long periods of time, cover the sensing element with the protective shipping cap.

## **Double Junction Reference Electrode Storage, Cat. No. 900200**

The double junction reference electrode may be stored in the outer chamber filling solution, Cat. No. 900003, between sample measurements and up to one week. The filling solution inside the electrode should not be allowed to evaporate, as crystallization will result.

For storage longer than one week, drain the reference electrode, flush the inside with distilled water and store the electrode dry.

## **Cadmium Combination Electrode Storage, Cat. No. 9648BNWP**

For storage between measurements and up to one week, store the electrode in a 4 M potassium chloride solution. Do not add ISA to the storage solution. The filling solution inside the electrode should not be allowed to evaporate, as crystallization will result.

For storage longer than one week, drain the electrode, flush the chamber with distilled water and store the electrode dry with the protective shipping cap covering the sensing element.

# Electrode Maintenance

## Polishing the Cadmium Combination Electrode and Cadmium Half-Cell Electrode

The sensing surface of solid state electrodes can wear over time, which causes drift, poor reproducibility and loss of response in low level samples. The electrode can be restored by polishing the sensing surface with a polishing strip, Cat. No. 948201. The polishing strip can also be used if the sensing surface has been etched or chemically poisoned.

1. Cut off about an inch of the polishing strip.
2. Hold the electrode with the sensing surface facing up.
3. Place a few drops of distilled water on the sensing surface.
4. With the frosted side of the polishing strip facing down, use light finger pressure to place the polishing strip on top of the sensing surface.
5. Rotate the electrode for about 30 seconds.
6. Rinse the electrode with distilled water and soak the electrode in a 1 ppm or  $10^{-5}$  M cadmium standard for ten minutes.

## Cadmium Combination Electrode and Double Junction Reference Electrode Flushing

If the area between the electrode sleeve and inner cone becomes clogged with sample or precipitate, flush the area with filling solution or distilled water.

1. Hold the electrode body with one hand and use your thumb to push down on the electrode cap to drain the electrode. Push down on the cap until all the filling solution is drained from the chamber.
2. Fill the electrode with distilled water and then push down on the cap until all the water is drained from the chamber.
3. Fill the electrode with fresh filling solution up to the fill hole. Push down on the cap to allow a few drops of filling solution to drain out of the electrode and replenish the lost filling solution.

## Disassembling and Reassembling the Cadmium Combination Electrode

**Note:** *Disassembly is usually not required and should not be done unless a thorough cleaning is required.*

1. Tip the electrode so the filling solution moistens the O-ring on the electrode body. Hold the electrode body with one hand and use your thumb to push down on the electrode cap to drain the electrode.
2. Unscrew the cap counterclockwise and then slide the cap and spring up the cable.
3. Hold the outer sleeve with one hand and firmly push down on the threaded portion with your thumb and forefinger to separate the inner body from the sleeve.
4. Grasp the inner cone with a clean, lint-free tissue and withdraw the body from the sleeve using a gentle twisting motion. Do not touch the pellet above the cone, as it will damage to the pellet. Rinse the outside of the electrode body and the entire sleeve with distilled water. Allow it to air dry.
5. Moisten the O-ring on the electrode body with a drop of filling solution. Insert the screw-thread end of the electrode body into the tapered, ground end of the sleeve.
6. Push the body into the sleeve using a gentle twisting motion until the bottom surface of the inner cone is flush with the tapered end of the sleeve.
7. Place the spring onto the electrode body and screw on the cap. Refill the electrode with filling solution.

# Analytical Techniques

A variety of analytical techniques are available to the analyst. The following is a description of these techniques.

**Direct Calibration** is a simple procedure for measuring a large number of samples. Only one meter reading is required for each sample. Calibration is performed using a series of standards. The concentration of the samples is determined by comparison to the standards. ISA is added to all solutions to ensure that samples and standards have similar ionic strength.

**Low Level Calibration** is similar to the direct calibration technique. This method is recommended when the expected sample concentration is less than 1 ppm or  $10^{-5}$  M cadmium. A minimum three point calibration is recommended to compensate for the electrode's non-linear response at these concentrations. A special calibration standard preparation procedure is the best means of preparing low level calibration standards.

**Incremental Techniques** provide a useful method for measuring samples, since a calibration is not required. The different incremental techniques are described below. They can be used to measure the total concentration of a specific ion in the presence of a large (50 to 100 times) excess of complexing agents. As in direct calibration, any convenient concentration unit can be used.

**Known Addition** is useful for measuring dilute samples, checking the results of direct calibration (when no complexing agents are present), or measuring the total concentration of an ion in the presence of an excess complexing agent. The electrode is immersed in the sample solution and an aliquot of a standard solution containing the measured species is added to the sample. From the change in potential before and after the addition, the original sample concentration is determined.



**Titration**s are quantitative analytical techniques for measuring the concentration of a species by incremental addition of a reagent (titrant) that reacts with the sample species. Sensing electrodes can be used for determination of the titration end point. Ion selective electrodes are useful as end point detectors, because they are unaffected by sample color or turbidity. Titrations are approximately 10 times more precise than direct calibration, but are more time-consuming.

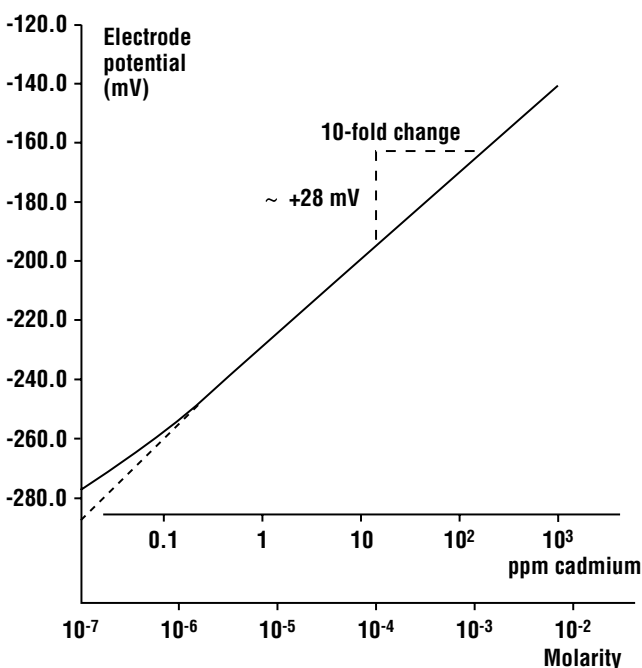
**Indicator Titration Method** is useful for measuring ionic species where an ion specific electrode does not exist. With this method the electrodes sense a reagent species that has been added to the sample before titration. The cadmium electrode may be used in indicator titrations for many different metal ions.

# Direct Calibration Technique

## Typical Direct Calibration Curve

In the direct calibration procedure, a calibration curve is constructed either in the meter memory or on semi-logarithmic paper. Electrode potentials of standard solutions are measured and plotted on the linear axis against their concentrations on the log axis. In the linear regions of the curves, only two standards are needed to determine a calibration curve. In non-linear regions, more points must be taken. These direct calibration procedures are given for concentrations in the region of linear electrode response. Low level measurement procedures are given in a following section for measurements in the non-linear electrode region.

**Figure 2**  
**Typical Direct Calibration Curve**



## Direct Calibration Overview

The following direct measurement procedures are recommended for moderate to high level measurements. Samples must be in the linear range of the electrode – greater than 1 ppm or  $10^{-5}$  M cadmium. A two point calibration is sufficient, although more points can be used. When using an ISE meter, sample concentrations can be read directly from the meter. When using a mV meter, a calibration curve can be prepared on semi-logarithmic graph paper, or a linear regression (against logarithmic concentration values) can be performed using a spreadsheet or graphing program.

### Calibration Hints

- Standard concentrations should bracket the expected sample concentrations.
- Always add 2 mL of ISA, Cat. No. 940011, per 100 mL of standard or sample.
- For high ionic strength samples that have an ionic strength of 0.1 M or greater, prepare standards with a background composition similar to that of the samples, or measure the samples using the known addition method.
- During calibration, measure the least concentrated standard first, and work up to the most concentrated standard.

## Direct Calibration Setup

1. Prepare the electrode as described in the **Electrode Preparation** section. If using the 9648BNWP combination cadmium electrode, fill the electrode with Cat. No. 900061. If using the 9448BN or 9448SC half-cell cadmium electrode with the 900200 reference electrode, fill the reference electrode with inner chamber filling solution, Cat. No. 900002, and outer chamber filling solution, Cat. No. 900003.
2. Connect the electrode to the meter.
3. Prepare at least two standards that bracket the expected sample range and differ in concentration by a factor of ten. Standards can be prepared in any concentration unit to suit the particular analysis requirement. See the **Serial Dilution** section for instructions on how to prepare standards. All standards should be at the same temperature as the samples. For details on temperature effects on electrode performance, refer to the **Temperature Effects** section.

## Direct Calibration Procedure Using a Meter with an ISE Mode

**Note:** See the meter user guide for more specific information.

1. Add 100 mL of the less concentrated standard and 2 mL of ISA to a 150 mL beaker and stir the solution thoroughly.
2. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the less concentrated standard. Wait for a stable reading and adjust the meter to display the value of the standard, as described in the meter user guide.
3. Add 100 mL of the more concentrated standard and 2 mL of ISA to a second 150 mL beaker and stir the solution thoroughly.
4. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the more concentrated standard. Wait for a stable reading and adjust the meter to display the value of the second standard, as described in the meter user guide.
5. Record the resulting slope value. The slope should be between 25 and 30 mV when the standards are between 20 and 25 °C.
6. Add 100 mL of sample and 2 mL of ISA to a clean 150 mL beaker and stir the solution thoroughly.
7. Rinse the electrode with distilled water, blot it dry and place it into the sample. The concentration of the sample will be displayed on the meter.

**Note:** Other solution volumes may be used, as long as the ratio of solution to ISA remains 50:1.

## Direct Calibration Procedure Using a Meter with a mV Mode

**Note:** See the meter user guide for more specific information.

1. Set the meter to the mV mode.
2. Add 100 mL of the less concentrated standard and 2 mL of ISA to a 150 mL beaker and stir the solution thoroughly.
3. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the less concentrated standard. When a stable reading is displayed, record the mV value and corresponding standard concentration.
4. Add 100 mL of the more concentrated standard and 2 mL of ISA to a second 150 mL beaker and stir the solution thoroughly.
5. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the more concentrated standard. When a stable reading is displayed, record the mV value and corresponding standard concentration.
6. Using semi-logarithmic graph paper, prepare a calibration curve by plotting the millivolt values on the linear axis and the standard concentration values on the logarithmic axis.
7. Add 100 mL of sample and 2 mL of ISA to a clean 150 mL beaker and stir the solution thoroughly.
8. Rinse the electrode with distilled water, blot it dry and place it into the beaker. When a stable reading is displayed, record the mV value.
9. Using the calibration curve prepared in step 6, determine the unknown concentration of the sample.

**Note:** Other solution volumes may be used, as long as the ratio of solution to ISA remains 50:1.

# Small Volume Direct Calibration Technique

Take advantage of special design features available with the 9648BNWP ionplus combination cadmium electrode to meet your measuring needs. Due to the Sure-Flow reference, this electrode is able to measure sample volumes as small as 5 mL using a modified direct measurement procedure. Because less solution volume is required, the chemical usage of cadmium standards and ISA is reduced. This method is also convenient when making field measurements, since the 9648BNWP combination cadmium electrode does not require a separate reference electrode. All samples should have a concentration greater than 1 ppm or  $10^{-5}$  M cadmium. A two point calibration is sufficient, although more points can be used. The following procedure recommends using 25 mL of sample. Smaller sample volumes can be used, as long as the final volume of solution is sufficient to cover the bottom of the electrode.

## Calibration Hints

- Use the 9648BNWP ionplus combination cadmium electrode.
- Standard concentrations should bracket the expected sample concentrations.
- Always keep the ratio of standard or sample to ISA at 50:1.
- For high ionic strength samples that have an ionic strength of 0.1 M or greater, prepare standards with a background composition similar to that of the samples, or measure the samples using the known addition method.
- During calibration, measure the least concentrated standard first, and work up to the most concentrated standard.
- Calibrate with the same volume of standard as the volume of sample that is available for analysis.

## Small Volume Direct Calibration Setup

1. Prepare the 9648BNWP combination cadmium electrode as described in the **Electrode Preparation** section and fill the electrode with Optimum Results A filling solution, Cat. No. 900061.
2. Connect the electrode to the meter.
3. Prepare at least two standards that bracket the expected sample range and differ in concentration by a factor of ten. Standards can be prepared in any concentration unit to suit the particular analysis requirement. See the **Serial Dilution** section for instructions on how to prepare standards. All standards should be at the same temperature as the samples. For details on temperature effects on electrode performance, refer to the **Temperature Effects** section.

## Small Volume Direct Calibration Procedure Using a Meter with an ISE Mode

**Note:** See the meter user guide for more specific information.

1. Add 25 mL of the less concentrated standard and 0.5 mL of ISA to a 50 mL beaker and swirl the solution to mix.
2. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the less concentrated standard. Wait for a stable reading and adjust the meter to display the value of the standard, as described in the meter user guide.
3. Add 25 mL of the more concentrated standard and 0.5 mL of ISA to a second 50 mL beaker and swirl the solution to mix.
4. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the more concentrated standard. Wait for a stable reading and adjust the meter to display the value of the second standard, as described in the meter user guide.
5. Record the resulting slope value. The slope should be between 25 and 30 mV when the standards are between 20 and 25 °C.
6. Add 25 mL of sample and 0.5 mL of ISA to a clean 50 mL beaker and swirl the solution to mix.
7. Rinse the electrode with distilled water, blot it dry and place it into the sample. The concentration of the sample will be displayed on the meter.

**Note:** Other solution volumes may be used, as long as the ratio of solution to ISA remains 50:1.



## Small Volume Direct Calibration Procedure Using a Meter with a mV Mode

**Note:** See the meter user guide for more specific information.

1. Set the meter to the mV mode.
2. Add 25 mL of the less concentrated standard and 0.5 mL of ISA to a 50 mL beaker and swirl the solution to mix.
3. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the less concentrated standard. When a stable reading is displayed, record the mV value and corresponding standard concentration.
4. Add 25 mL of the more concentrated standard and 0.5 mL of ISA to a second 50 mL beaker and swirl the solution to mix.
5. Rinse the electrode with distilled water, blot it dry and place it into the beaker with the more concentrated standard. When a stable reading is displayed, record the mV value and corresponding standard concentration.
6. Using semi-logarithmic graph paper, prepare a calibration curve by plotting the millivolt values on the linear axis and the standard concentration values on the logarithmic axis.
7. Add 25 mL of sample and 0.5 mL of ISA to a clean 50 mL beaker and swirl the solution to mix.
8. Rinse the electrode with distilled water, blot it dry and place it into the beaker. When a stable reading is displayed, record the mV value.
9. Using the calibration curve prepared in step 6, determine the unknown concentration of the sample.

**Note:** Other solution volumes may be used, as long as the ratio of solution to ISA remains 50:1.

# Low Level Calibration Technique

These procedures are for solutions that have a cadmium concentration of less than 1 ppm or  $10^{-5}$  M cadmium. For solutions low in cadmium but high in total ionic strength (greater than  $10^{-1}$  M), perform the same procedure by preparing a calibrating solution with a composition similar to the sample.

Accurate results require that the following conditions be met:

- Prepare at least three calibration standards that bracket the expected sample concentration.
- Always use low level ISA for standards and samples.
- Plastic labware must be used for all low level cadmium measurements.
- Adequate time must be allowed for electrode stabilization. Longer response time will be needed at low level measurements.
- Stir all standards and samples at a uniform rate.

## Low Level Setup

1. Prepare the electrode as described in the **Electrode Preparation** section.
2. Connect the electrode to the meter. Set the meter to the mV mode.
3. Prepare the low level ISA by pipetting 20 mL of the ISA, Cat. No. 940011, into a 100 mL volumetric flask and diluting to the mark with distilled water. Use low level ISA for low level measurements only.
4. Select a standard solution. Use either a 10 ppm cadmium standard or a  $10^{-4}$  M cadmium standard.

To prepare the 10 ppm cadmium standard, pipet 10 mL of the 1000 ppm cadmium standard into a 1 liter volumetric flask. Dilute to the mark with distilled water and mix the solution thoroughly.

To prepare the  $10^{-4}$  M cadmium standard, pipet 1 mL of the 0.1 M cadmium standard into a 1 liter volumetric flask. Dilute to the mark with distilled water and mix the solution thoroughly.

## Low Level Calibration and Measurement

1. Add 100 mL of distilled water and 1 mL of low level ISA to a 150 mL beaker.
2. Rinse the electrode with distilled water, blot it dry and place it into the beaker. Stir the solution thoroughly.
3. Add increments of the 10 ppm or  $10^{-4}$  M cadmium standard mixed with low level ISA to the beaker using the steps outlined in **Table 2**. Record the stable millivolt reading after each increment.
4. On semi-logarithmic paper, plot the concentration (log axis) against the millivolt potential (linear axis). Prepare a new calibration curve with fresh standards each day.
5. Measure 100 mL of sample and 1 mL of low level ISA and pour the solutions into a clean 150 mL beaker. Rinse the electrode with distilled water, blot it dry and place the electrode into the sample.
6. Stir the solution thoroughly. When a stable reading is displayed, record the mV value.
7. Determine the sample concentration corresponding to the measured potential from the low level calibration curve.

**Table 2**

### Calibration Curve For Low Level Calibrations

Additions of standard to 100 mL distilled water and 1 mL low level ISA solution.

Step	Pipet Size	Volume Added	Concentration ppm	Concentration M
1	1 mL	0.1 mL	0.01	$1.0 \times 10^{-7}$
2	1 mL	0.3 mL	0.04	$4.0 \times 10^{-7}$
3	1 mL	0.6 mL	0.10	$1.0 \times 10^{-6}$
4	2 mL	2.0 mL	0.30	$3.0 \times 10^{-6}$

# Known Addition Technique

Known addition is a convenient technique for measuring samples in the linear range of the electrode (greater than 0.2 ppm cadmium) because no calibration curve is required. It can be used to verify the results of a direct calibration or to measure the total concentration of an ion in the presence of a large excess of a complexing agent. The sample potential is measured before and after addition of a standard solution.

Accurate results require that the following conditions be met:

- Concentration should approximately double as a result of the addition.
- Sample concentration should be known to within a factor of three.
- Either no complexing agent or a large excess of the complexing agent may be present.
- The ratio of the uncomplexed ion to complexed ion must not be changed by addition of the standard.
- All samples and standards should be at the same temperature.
- With double or multiple known addition, the final addition should be 10 to 100 times the sample concentration.
- Add 2 mL of ISA to every 100 mL of sample before analysis.

## Known Addition Setup

1. Prepare the electrode as described in the **Electrode Preparation** section.
2. Connect the electrode to the meter.
3. Prepare a standard solution that will cause the cadmium concentration of the sample to double when added to the sample solution. Refer to **Table 3** for guidelines.
4. Determine the electrode slope by performing the procedure in the **Checking Electrode Operation (Slope)** section.
5. Rinse the electrode with distilled water.

**Table 3**  
**Guideline For Known Addition**

Volume of Addition	Concentration of Standard
1 mL	100 times sample concentration
5 mL	20 times sample concentration
10 mL*	10 times sample concentration

\* Most convenient volume to use

## Known Addition Using a Meter with a Known Addition Mode

**Note:** See the meter user guide for more specific information.

1. Set the meter to measure in the known addition mode.
2. Measure 100 mL of the sample and 2 mL of ISA and pour the solutions into a beaker. Rinse the electrode with distilled water and place it into the sample solution. Stir the solution thoroughly.
3. When a stable reading is displayed, set the meter as described in the meter user guide, if required.
4. Pipet the appropriate amount of the standard solution into the beaker. Stir the solution thoroughly.
5. When a stable reading is displayed, record the sample concentration.

## Known Addition Using a Meter with a Millivolt Mode

1. Set the meter to the relative millivolt mode. If a relative millivolt mode is not available, use the millivolt mode.
2. Measure 100 mL of sample and 2 mL of ISA and pour the solutions into a 150 mL beaker. Stir the solution thoroughly.
3. Rinse the electrode with distilled water, blot it dry and place the electrode into the beaker. When a stable reading is displayed, set the meter to read 0.0 mV. If the reading cannot be adjusted to 0.0 mV, record the actual mV value.
4. Pipet the appropriate amount of standard solution into the beaker. Stir the solution thoroughly.
5. When a stable reading is displayed, record the mV value. If the meter could not be set to 0.0 mV in step 3, subtract the first reading from the second reading to calculate  $\Delta E$ .
6. Use **Table 5** to find the Q value that corresponds to the change in potential,  $\Delta E$ . To determine the original sample concentration, multiply Q by the concentration of the added standard:

$$C_{\text{sample}} = Q * C_{\text{standard}}$$

$C_{\text{standard}}$  = standard concentration

$C_{\text{sample}}$  = sample concentration

Q = value from **Table 5**

The table of Q values is calculated for a 10% volume change. The equation for the calculation of Q for different slopes and volume changes is given below.

$$Q = (p * r) / \{[(1 + p) * 10^{\Delta E/S}] - 1\}$$

Q = value from **Table 5**

$\Delta E = E_2 - E_1$

S = slope of the electrode

p = volume of standard / volume of sample and ISA

r = volume of sample and ISA / volume of sample

# Calculating Known Addition for Samples using Lotus, Excel, or Quattro Spreadsheets

If it is more convenient, a simple spreadsheet can be set up to calculate the known addition results, using any ratio of sample to addition. A typical worksheet is shown in **Table 4**. The numbers shown are examples, but the formulas and their locations should be copied exactly.

**Table 4**  
**Known Addition Calculations using Lotus, Excel, or Quattro Spreadsheets**

A	B	C
1		Enter Value
2	Volume of sample and ISA (mL)	102
3	Volume of addition (mL)	10
4	Concentration of addition	10
5	Volume of sample	100
6	Initial mV reading	45.3
7	Final mV reading	63.7
8	Electrode slope	28.2
9		
10		Derived Values
11	Delta E	+C7 - C6
12	Solution volume ratio	+C3/C2
13	Antilog term	+10^ (C11/C8)
14	Sample volume ratio	+C2/C5
15	Q term	+C12*C14/ (((1+C12)*C13)-1)
16	Calculated initial concentration in same units as addition	+C15*C4

**Note:** For Excel, use = instead of + at start of formulas.



**Table 5**

**Q Values for a 10% volume change,  
slopes (in column heading) are in units of mV/decade**

<b>ΔE</b>	<b>Q Concentration Ratio</b>			
	<b>28.6</b>	<b>29.1</b>	<b>29.6</b>	<b>30.1</b>
2.5	0.2917	0.2957	0.2996	0.3035
2.6	0.2827	0.2867	0.2906	0.2944
2.7	0.2742	0.2781	0.2820	0.2858
2.8	0.2662	0.2700	0.2738	0.2775
2.9	0.2585	0.2623	0.2660	0.2697
3.0	0.2512	0.2550	0.2586	0.2623
3.1	0.2443	0.2480	0.2516	0.2552
3.2	0.2377	0.2413	0.2449	0.2484
3.3	0.2314	0.2349	0.2384	0.2419
3.4	0.2253	0.2288	0.2323	0.2357
3.5	0.2196	0.2230	0.2264	0.2298
3.6	0.2140	0.2174	0.2208	0.2241
3.7	0.2087	0.2121	0.2154	0.2187
3.8	0.2037	0.2070	0.2102	0.2135
3.9	0.1988	0.2020	0.2052	0.2084
4.0	0.1941	0.1973	0.2005	0.2036
4.1	0.1896	0.1927	0.1959	0.1990
4.2	0.1852	0.1884	0.1914	0.1945
4.3	0.1811	0.1841	0.1872	0.1902
4.4	0.1770	0.1801	0.1831	0.1861
4.5	0.1732	0.1762	0.1791	0.1821
4.6	0.1694	0.1724	0.1753	0.1782
4.7	0.1658	0.1687	0.1716	0.1745
4.8	0.1623	0.1652	0.1680	0.1709
4.9	0.1590	0.1618	0.1646	0.1674
5.0	0.1557	0.1585	0.1613	0.1640
5.1	0.1525	0.1553	0.1580	0.1608
5.2	0.1495	0.1522	0.1549	0.1576
5.3	0.1465	0.1492	0.1519	0.1546
5.4	0.1437	0.1463	0.1490	0.1516
5.5	0.1409	0.1435	0.1461	0.1487
5.6	0.1382	0.1408	0.1434	0.1459
5.7	0.1356	0.1382	0.1407	0.1432
5.8	0.1331	0.1356	0.1381	0.1406
5.9	0.1306	0.1331	0.1356	0.1381
6.0	0.1282	0.1307	0.1331	0.1356
6.1	0.1259	0.1283	0.1308	0.1332
6.2	0.1236	0.1260	0.1284	0.1308
6.3	0.1214	0.1238	0.1262	0.1285
6.4	0.1193	0.1217	0.1240	0.1263
6.5	0.1172	0.1195	0.1219	0.1242
6.6	0.1152	0.1175	0.1198	0.1221
6.7	0.1132	0.1155	0.1178	0.1200
6.8	0.1113	0.1136	0.1158	0.1180
6.9	0.1094	0.1117	0.1139	0.1161

$\Delta E$	Q Concentration Ratio			
	28.6	29.1	29.6	30.1
7.0	0.1076	0.1098	0.1120	0.1142
7.1	0.1058	0.1080	0.1102	0.1123
7.2	0.1041	0.1063	0.1084	0.1105
7.3	0.1024	0.1045	0.1067	0.1088
7.4	0.1008	0.1029	0.1050	0.1071
7.5	0.0992	0.1012	0.1033	0.1054
7.8	0.0946	0.0966	0.0986	0.1006
8.0	0.0917	0.0936	0.0956	0.0976
8.3	0.0876	0.0895	0.0914	0.0933
8.5	0.0850	0.0869	0.0887	0.0906
8.8	0.0813	0.0831	0.0849	0.0868
9.0	0.0790	0.0808	0.0825	0.0843
9.3	0.0757	0.0774	0.0791	0.0809
9.5	0.0736	0.0753	0.0770	0.0787
9.8	0.0706	0.0722	0.0739	0.0755
10.0	0.0687	0.0703	0.0719	0.0735
10.3	0.0660	0.0675	0.0691	0.0707
10.5	0.0642	0.0658	0.0673	0.0689
10.8	0.0617	0.0633	0.0648	0.0663
11.0	0.0602	0.0617	0.0631	0.0646
11.3	0.0579	0.0593	0.0608	0.0623
11.5	0.0564	0.0579	0.0593	0.0607
11.8	0.0544	0.0558	0.0572	0.0585
12.0	0.0530	0.0544	0.0558	0.0572
12.3	0.0511	0.0525	0.0538	0.0551
12.5	0.0499	0.0512	0.0525	0.0539
12.8	0.0481	0.0494	0.0507	0.0520
13.0	0.0470	0.0483	0.0495	0.0508
13.3	0.0454	0.0466	0.0478	0.0491
13.5	0.0443	0.0455	0.0468	0.0480
13.8	0.0428	0.0440	0.0452	0.0464
14.0	0.0419	0.0430	0.0442	0.0454
14.3	0.0404	0.0416	0.0427	0.0439
14.5	0.0395	0.0407	0.0418	0.0429
14.8	0.0382	0.0393	0.0404	0.0416
15.0	0.0374	0.0385	0.0396	0.0407
15.5	0.0354	0.0365	0.0375	0.0386
16.0	0.0335	0.0345	0.0356	0.0366
16.5	0.0318	0.0328	0.0337	0.0347
17.0	0.0302	0.0311	0.0320	0.0330
17.5	0.0286	0.0295	0.0305	0.0314
18.0	0.0272	0.0281	0.0290	0.0298
18.5	0.0258	0.0267	0.0275	0.0284
19.0	0.0246	0.0254	0.0262	0.0270
19.5	0.0234	0.0242	0.0250	0.0258
20.0	0.0223	0.0230	0.0238	0.0246
20.5	0.0212	0.0219	0.0227	0.0234
21.0	0.0202	0.0209	0.0216	0.0224

<b>ΔE</b>	<b>Q Concentration Ratio</b>			
	<b>28.6</b>	<b>29.1</b>	<b>29.6</b>	<b>30.1</b>
21.5	0.0192	0.0199	0.0206	0.0213
22.0	0.0183	0.0190	0.0197	0.0204
22.5	0.0175	0.0181	0.0188	0.0195
23.0	0.0167	0.0173	0.0179	0.0186
23.5	0.0159	0.0165	0.0171	0.0178
24.0	0.0152	0.0158	0.0164	0.0170
24.5	0.0145	0.0151	0.0157	0.0162
25.0	0.0139	0.0144	0.0150	0.0155
25.5	0.0132	0.0138	0.0143	0.0149
26.0	0.0126	0.0132	0.0137	0.0142
26.5	0.0121	0.0126	0.0131	0.0136
27.0	0.0116	0.0120	0.0125	0.0131
27.5	0.0110	0.0115	0.0120	0.0125
28.0	0.0106	0.0110	0.0115	0.0120
28.5	0.0101	0.0106	0.0110	0.0115
29.0	0.0097	0.0101	0.0105	0.0110
29.5	0.0093	0.0097	0.0101	0.0105
30.5	0.0085	0.0089	0.0093	0.0097
31.5	0.0078	0.0081	0.0085	0.0089
32.0	0.0074	0.0078	0.0082	0.0085
32.5	0.0071	0.0075	0.0078	0.0082
33.0	0.0068	0.0072	0.0075	0.0079
33.5	0.0065	0.0069	0.0072	0.0076
34.0	0.0063	0.0066	0.0069	0.0072
34.5	0.0060	0.0063	0.0066	0.0070
35.0	0.0058	0.0061	0.0064	0.0067
35.5	0.0055	0.0058	0.0061	0.0064
36.0	0.0053	0.0056	0.0059	0.0062
36.5	0.0051	0.0053	0.0056	0.0059
37.0	0.0049	0.0051	0.0054	0.0057
37.5	0.0047	0.0049	0.0052	0.0055
38.0	0.0045	0.0047	0.0050	0.0052
38.5	0.0043	0.0045	0.0048	0.0050
39.0	0.0041	0.0043	0.0046	0.0048
39.5	0.0039	0.0042	0.0044	0.0046
40.0	0.0038	0.0040	0.0042	0.0045
40.5	0.0036	0.0038	0.0041	0.0043
41.0	0.0035	0.0037	0.0039	0.0041
41.5	0.0033	0.0035	0.0037	0.0040
42.0	0.0032	0.0034	0.0036	0.0038
42.5	0.0031	0.0033	0.0035	0.0037
43.0	0.0029	0.0031	0.0033	0.0035
43.5	0.0028	0.0030	0.0032	0.0034
44.0	0.0027	0.0029	0.0031	0.0032
44.5	0.0026	0.0028	0.0029	0.0031
45.0	0.0025	0.0027	0.0028	0.0030
45.5	0.0024	0.0026	0.0027	0.0029
46.0	0.0023	0.0024	0.0026	0.0028

# Cadmium Titration Technique

The cadmium electrode makes a highly sensitive endpoint detector for titration with EDTA of cadmium samples. Titrations are more time consuming than direct electrode measurement, but results are more accurate and reproducible. With careful technique, titrations accurate to  $\pm 0.1\%$  of the total cadmium ion concentration of the sample can be performed.

EDTA complexes other cations besides cadmium ion. Interferences from other ions, whose EDTA complexes are stable only at low pH, can be eliminated by performing the titration for cadmium ion at a high pH, about pH 10 (adjusted with ammonia). In many cases, other interferences can be eliminated by a suitable choice of sample pH and the addition of masking agents to the sample solution. A comprehensive list of methods is given in the Handbook of Analytical Chemistry, L. Meites, (ed.) McGraw Hill Book Co., New York, (1st edit.), pp. 3-76, 3-225.

## Cadmium Titration Setup

1. Prepare the electrode as described in the **Electrode Preparation** section. If using the 9648BNWP combination cadmium electrode, fill the electrode with Cat. No. 900061. If using the 9448BN or 9448SC half-cell cadmium electrode with the 900200 reference electrode, fill the reference electrode with inner chamber filling solution, Cat. No. 900002, and outer chamber filling solution, Cat. No. 900003.
2. Connect the electrode to the meter.
3. Prepare a 1 M EDTA stock solution by adding 38.0 grams of reagent-grade  $\text{Na}_4\text{EDTA}$  to a 100 mL volumetric flask. Dissolve the solids with about 75 mL of distilled water and then dilute to the mark with distilled water.
4. Prepare an EDTA titrant solution 10 to 20 times as concentrated as the sample by dilution of the 1 M EDTA stock solution. For a good endpoint break, the sample concentration should be at least  $10^{-4}$  M in total cadmium.

## Cadmium Titration Procedure

1. Place 100 mL of sample into a 150 mL beaker and adjust the pH of the sample to 10 using  $\text{NH}_4\text{OH}$ . Place the electrode in the sample and stir the solution thoroughly.
2. Using a 10 mL burette, add increments of titrant and plot the electrode potential against mL of titrant added. The endpoint is the point of greatest slope (inflection point). See **Figure 3**.

3. Calculate the sample concentration before dilution:

$$C_{\text{sample}} = C_t (V_t / V_{\text{sample}})$$

$C_{\text{sample}}$  = sample concentration

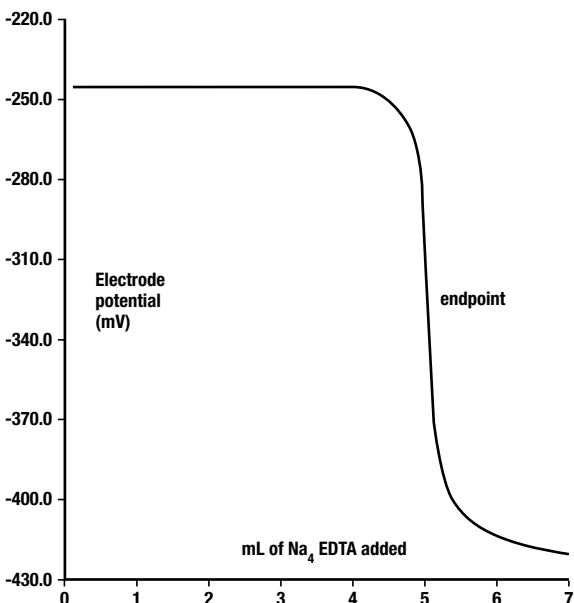
$C_t$  = titrant concentration

$V_{\text{sample}}$  = sample volume

$V_t$  = titrant volume added at endpoint.

**Figure 3**

**Typical Titration of 100 mL of  $5 \times 10^{-3}$  M  $\text{Cd}(\text{NO}_3)_2$  with 0.1 M  $\text{Na}_4\text{EDTA}$  when pH was adjusted to 10 with ammonia**



## Indicator Titrations

The cadmium electrode can be substituted for the cupric electrode to detect the endpoint in titrations of other metal ions. A small amount of cadmium complex is added to the sample, and a complexometric titration is done. The endpoint volume of titrant is used to calculate the sample concentration. The minimum level of sample ion that can be determined by indicator titration is above  $10^{-4}$  M. Titrations of barium, calcium, cobalt (+2), magnesium, manganese (+2), nickel and strontium are possible. For more information refer to Chelometric Indicator Titrations with the Solid State Cupric Ion Selective Electrode, Ross J.W., and Frant, M.S.; Anal. Chem., 1969 41(13), 1900.

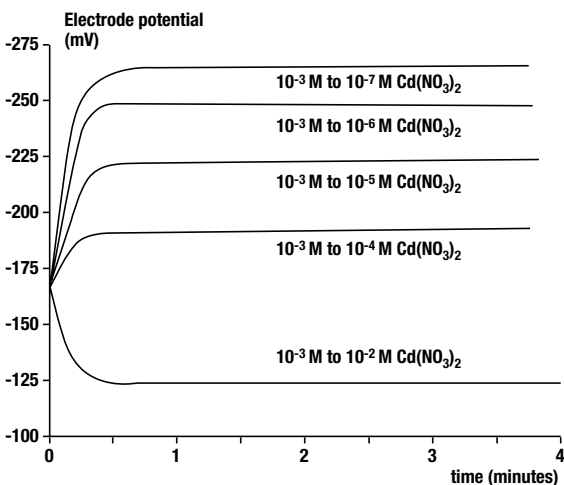
# Electrode Characteristics

## Electrode Response

The electrode potential plotted against concentration on semi-logarithmic paper results in a straight line with a slope of about 25 to 30 mV per decade change in concentration.

The time response of the electrode (the time required to reach 99% of the stable potential reading) varies from several seconds in concentrated solutions to several minutes near the limit of detection.

**Figure 4**  
**Typical Electrode Response to Step Changes in  $\text{Cd}(\text{NO}_3)_2$  Concentration**



## Reproducibility

Reproducibility is limited by factors such as temperature fluctuations, drift and noise. Within the operating range of the electrode, reproducibility is independent of concentration. With hourly calibrations, direct electrode measurements reproducible to  $\pm 4\%$  can be obtained.

## Temperature Effects

Since electrode potentials are affected by changes in temperature, samples and standard solutions should be within  $\pm 1\text{ }^{\circ}\text{C}$  ( $\pm 2\text{ }^{\circ}\text{F}$ ) of each other. At the  $10^{-3}\text{ M}$  level, a  $1\text{ }^{\circ}\text{C}$  difference in temperature results in errors greater than 4 %. The absolute potential of the reference electrode changes slowly with temperature because of the solubility equilibria on which the electrode depends. The slope of the electrode also varies with temperature, as indicated by the factor  $S$  in the Nernst equation. Theoretical values of the slope at different temperatures are given in **Table 6**. If the temperature changes, the meter and electrode should be recalibrated.

The electrode can be used at temperatures from 0 to  $80\text{ }^{\circ}\text{C}$ , provided that temperature equilibrium has occurred. For use at temperatures substantially different from room temperature, calibration standards should be at the same temperature as samples. The electrode must be used only intermittently at solution temperatures above  $80\text{ }^{\circ}\text{C}$ .

**Table 6**  
**Theoretical Slope vs. Temperature Values**

Temperature ( $^{\circ}\text{C}$ )	Slope (mV)
0	27.1
10	28.1
20	29.1
25	29.6
30	30.1
40	31.1
50	32.1

If sample temperatures vary, use of the 9648BNWP combination cadmium electrode is recommended. The Optimum Results A filling solution that is included with the electrode will minimize junction potentials and provide optimum temperature and time response. Optimum Results A filling solution produces an isopotential point of  $1.7 \times 10^{-3}\text{ M}$  cadmium. The isopotential point is the concentration at which the potential of the electrode does not vary with temperature. Since the isopotential point of this electrode is known, the combination cadmium electrode may be used on meters that allow automatic temperature compensation for ISE measurements. By programming in the isopotential point and placing an ATC probe into the sample, any time the temperature changes the meter will automatically adjust the slope of the calibration curve, resulting in more accurate measurement results.



## Interferences

Mercury and silver ions poison the cadmium electrode sensing element and must be absent from the sample solution. Exposure to either of these species at levels greater than  $10^{-7}$  M will require polishing of the electrode sensing surface. Ferric ions affect the sensing element only if the ferric ion level is greater than one tenth of cadmium ion level (ferric ion can be eliminated from the sample by adding sodium fluoride and adjusting the sample to pH 4 to 6). Lead ions affect the membrane surface if the level of lead ion exceeds the level of cadmium ion present in sample. Copper ions may also be an interference.

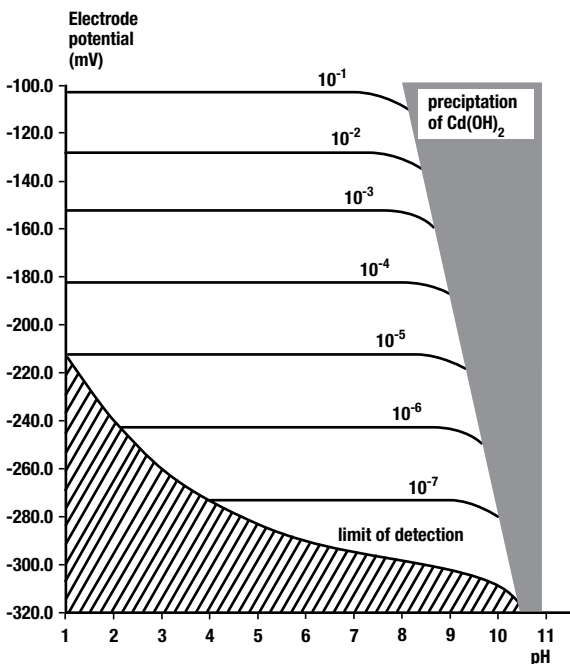
If the electrode is exposed to high levels of interfering ions, it may become unstable and sluggish in response. When this happens, restore normal electrode performance by polishing it. Refer to the **Electrode Maintenance** section.

## pH Effects

The electrode response to cadmium ions in solutions at various pH values is shown in **Figure 5**. Although the electrode can be used over a wide pH range, hydrogen ions interfere with measurements of low levels of cadmium ion. The edge of the shaded area to the left in **Figure 5** indicates the minimum pH at which low level cadmium measurements can be made without hydrogen ion interference.

At high solution pH, sufficient hydroxide ions are present to form a precipitate with a portion of the cadmium ions, reducing the level of free cadmium ions in the sample. As shown in **Figure 5**,  $\text{Cd}(\text{OH})_2$  forms at a higher pH in dilute solutions than in concentrated solutions. Since the electrode responds only to free, unbound cadmium ion, it does not detect that portion of the cadmium precipitated by hydroxide ion. Precipitation can be avoided by adjusting the pH of sample and standards to below 7.

**Figure 5**  
**Electrode Potential Behavior vs, Solution pH in Pure  $\text{Cd}(\text{NO}_3)_2$  Solutions at 25 °C**



## Complexation

Cadmium forms complexes with a wide variety of species including acetate, ammonia, bromide, chloride, citrate, cyanide and EDTA. The extent of complexation depends on the concentration of cadmium and complexing agent and the pH. Since the electrode only responds to free cadmium ions, complexation reduces the measured concentration. In a large excess (50 to 100 times) of a complexing agent, the total cadmium concentration can be measured by known addition.

Soluble cadmium salts are precipitated by sulfide, carbonate, oxalate, phosphate, hydroxide and other ions. The formation of a precipitate depends on the level of cadmium ion, the level of the precipitating ion in the sample solution and the solution pH.

## Theory of Operation

The cadmium electrode consists of a sensing element bonded into an epoxy body. When the sensing element is in contact with a solution containing cadmium ions, an electrode potential develops across the sensing element. This potential, which depends on the level of free cadmium ion in solution, is measured against a constant reference potential with a digital pH/mV meter or ISE (concentration) meter. The measured potential corresponding to the level of cadmium ion in solution is described by the Nernst equation.

$$E = E_o + S * \log (A)$$

E = measured electrode potential

E<sub>o</sub> = reference potential (a constant)

A = cadmium ion activity level in solution

S = electrode slope (about 28 mV per decade)

$S = (2.3 RT) / nF$

R and F are constants, T = temperature in degrees K and n = ionic charge

The level of cadmium ions, A, is the activity or “effective concentration” of free cadmium ions in solution. The cadmium ion activity is related to free cadmium ion concentration, C<sub>f</sub>, by the activity coefficient, γ.

$$A = \gamma * C_f$$

Ionic activity coefficients are variable and largely depend on total ionic strength. The ionic strength of a solution is determined by all of the ions present. It is calculated by multiplying the concentration of each individual ion by the square of its charge, adding all these values up and then dividing by two.

$$\text{Ionic strength} = 1/2 \sum (C_i Z_i^2)$$

C<sub>i</sub> = concentration of ion i

Z<sub>i</sub> = charge of ion i

Σ symbolizes the sum of all the types of ions in solutions

If background ionic strength is high and constant relative to the sensed ion concentration, the activity coefficient is constant and activity is directly proportional to concentration. Ionic strength adjustor (ISA) is added to all cadmium standards and samples so that the background ionic strength is high and constant relative to variable concentrations of cadmium. For cadmium, the recommended ISA is 5 M NaNO<sub>3</sub>. Other solutions can be used as long as they do not contain ions that would interfere with the electrode response to cadmium.

If samples have a high ionic strength (above 0.1 M), standards should be prepared with a composition similar to the samples.

Reference electrode conditions must also be considered. Liquid junction potentials arise any time when two solutions of different composition are brought into contact. The potential results from the interdiffusion of ions in the two solutions. Since ions diffuse at different rates, the electrode charge will be carried unequally across the solution boundary resulting in a potential difference between the two solutions. In making electrode measurements, it is important that this potential is the same when the reference is in the standardizing solution as well as in the same solution; otherwise, the change in liquid junction potential will appear as an error in the measured specific ion electrode potential.

The most important variable that analysts have under their control is the composition of the liquid junction filling solution. The filling solution should be equitransferent. That is, the speed with which the positive and negative ions in the filling solution diffuse into the sample should be nearly as equal as possible. If the rate at which positive and negative charge is carried into the sample solution is equal, then no junction potential can result. Optimum Results filling solutions are specifically designed to meet all reference electrode conditions.

# Troubleshooting

Follow a systematic procedure to isolate the problem. The measuring system can be divided into four components for ease in troubleshooting: meter, electrode, sample/application and technique.

## Meter

The meter is the easiest component to eliminate as a possible cause of error. Thermo Scientific Orion meters include an instrument checkout procedure and shorting cap for convenience in troubleshooting. Consult the meter user guide for directions.

## Electrode

1. Rinse the electrode thoroughly with distilled water.
2. Verify the electrode performance by performing the procedure in the **Checking Electrode Operation (Slope)** section.
3. If the electrode fails this procedure, review the **Measuring Hints** section. Clean the electrode thoroughly as directed in the **Electrode Maintenance** section. Drain and refill the electrode with fresh filling solution.
4. Repeat the procedure in the **Checking Electrode Operation (Slope)** section.
5. If the electrode fails this procedure again and the half-cell cadmium electrode is being used, determine whether the cadmium or reference electrode is at fault. To do this, substitute a known working electrode for the electrode in question and repeat the procedure in the **Checking Electrode Operation (Slope)** section.
6. If the electrode passes the procedure, but measurement problems persist, the sample may contain interferences or complexing agents, or the technique may be in error.
7. Before replacing a faulty electrode, review this user guide and be sure to thoroughly clean the electrode; correctly prepare the electrode; use the proper filling solution, ISA, and standards; correctly measure the samples and review the **Troubleshooting Checklist** section.

## Sample/Application

The quality of results depends greatly upon the quality of the standards. Always prepare fresh standards when problems arise, it could save hours of frustrating troubleshooting! Errors may result from contamination of prepared standards, accuracy of dilution, quality of distilled water, or a mathematical error in calculating the concentrations.

The best method for preparation of standards is serial dilution. Refer to the **Serial Dilution** section. The electrode and meter may operate with standards, but not with the sample. In this case, check the sample composition for interferences, incompatibilities or temperature effects. Refer to the **Sample Requirements, Temperature Effects, Interferences, pH Effects** and **Complexation** sections.

## Technique

If trouble persists, review operating procedures. Review calibration and measurement sections to be sure proper technique has been followed. Verify that the expected concentration of the ion of interest is within the limit of detection of the electrode.

Check the method of analysis for compatibility with your sample. Direct measurement may not always be the method of choice. If a large amount of complexing agents are present, known addition may be the best method. If working with low level samples, follow the procedure in the **Low Level Calibration** section.

## Assistance

After troubleshooting all components of your measurement system, contact Technical Support. Within the United States call 1.800.225.1480 and outside the United States call 978.232.6000 or fax 978.232.6031. In Europe, the Middle East and Africa, contact your local authorized dealer. For the most current contact information, visit [www.thermo.com/contactwater](http://www.thermo.com/contactwater).

For the latest application and technical resources for Thermo Scientific Orion products, visit [www.thermo.com/waterapps](http://www.thermo.com/waterapps).

## Warranty

For the most current warranty information, visit [www.thermo.com/water](http://www.thermo.com/water).

# Troubleshooting Checklist

- No electrode filling solution added –  
Fill the electrode with filling solution up to the fill hole. Refer to the **Electrode Preparation** section for details.
- Incorrect electrode filling solution used –  
Refer to the **Electrode Preparation** section to verify the correct electrode filling solution.
- Electrode junction is dry –  
Push down on the electrode cap to allow a few drops of filling solution to drain out of the electrode.
- No reference electrode present –  
The 9448BN and 9448SC cadmium half-cell electrodes require a separate reference electrode, Cat. No. 900200.
- Electrode is clogged or dirty –  
Refer to the **Electrode Maintenance** section for cleaning instructions.
- Sensing element is dirty or etched –  
Refer to the **Electrode Maintenance** section for cleaning instructions.
- Standards are contaminated or made incorrectly –  
Prepare fresh standards. Refer to the **Measurement Hints** and **Analytical Techniques** sections.
- ISA not used or incorrect ISA used –  
ISA must be added to all standards and samples. Refer to the **Required Equipment** section for information on the ISA.
- Samples and standards at different temperatures –  
Allow solutions to reach the same temperature.
- Air bubble on sensing element –  
Remove air bubble by reimmersing the electrode in solution.
- Electrode not properly connected to meter –  
Unplug and reconnect the electrode to the meter.
- Meter or stir plate not properly grounded –  
Check the meter and stir plate for proper grounding.
- Static electricity present –  
Wipe plastic parts on the meter with a detergent solution.
- Defective meter –  
Check the meter performance. See the meter user guide.



# Ordering Information

Cat. No.	Description
9648BNWP	Cadmium ionplus Sure-Flow combination electrode, waterproof BNC connector
900061	Optimum Results A electrode filling solution, 5 x 60 mL bottles
9448BN	Cadmium half-cell electrode, BNC connector (requires separate reference electrode)
9448SC	Cadmium half-cell electrode, screw cap connector (requires separate reference electrode)
900200	Double junction reference electrode, pin tip connector
900002	Inner chamber filling solution for the double junction reference electrode, 5 x 60 mL bottles
900003	Outer chamber filling solution for the double junction reference electrode, 5 x 60 mL bottles
940011	ISA for cadmium measurements, 475 mL bottle
984201	Polishing strips

# Specifications

## Concentration Range

$10^{-7}$  M to 0.1 M (0.01 ppm to 11,200 ppm)

## pH Range

2 to 12

## Temperature Range

0 to 80 °C continuous use, 80 to 100 °C intermittent use

## Electrode Resistance

Less than 1 megohms

## Reproducibility

± 4%

## Minimum Sample Size (9648BNWP)

5 mL in a 50 mL beaker

## Size— 9648BNWP

Body Diameter: 13 mm

Body Length: 110 mm

Cap Diameter: 16 mm

Cable Length: 1 meter

## Size— 9448BN and 9448SC

Body Diameter: 12 mm

Body Length: 110 mm

Cap Diameter: 16 mm

Cable Length: 1 meter (9448BN only)

*\* Specifications are subject to change without notice*

258500-001 Rev. A 12-08

**Thermo Fisher Scientific**  
**Environmental Instruments**  
Water Analysis Instruments



**North America**

166 Cummings Center  
Beverly, MA 01915 USA  
Toll Free: 1-800-225-1480  
Tel: 1-978-232-6000  
Dom. Fax: 1-978-232-6015  
Int'l Fax: 978-232-6031

**Europe**

P.O. Box 254, 3860 AG Nijkerk  
Wallerstraat 125K, 3862 BN  
Nijkerk, Netherlands  
Tel: (31) 033-2463887  
Fax: (31) 033-2460832

**Asia Pacific**

Blk 55, Ayer Rajah Crescent  
#04-16/24, Singapore 139949  
Tel: 65-6778-6876  
Fax: 65-6773-0836

[www.thermo.com/water](http://www.thermo.com/water)

**Thermo**  
S C I E N T I F I C