

# Thermo Scientific Orion Dual Star

## pH/ISE Dual Channel Benchtop Meter

255100-001 • Revision C • June 2015





### **Important Note**

Please read this user guide thoroughly before using your meter. Any use outside of these instructions may invalidate your warranty and cause permanent damage to the meter.

### **Contact Information**

For assistance with Thermo Scientific<sup>™</sup> Orion<sup>™</sup> products, contact Technical Support by email at <u>wai.techservbev@thermofisher.com</u> or by phone – within the United States call 1-800-225-1480 and outside the United States call +1-978-232-6000 or fax +1-978-232-6031.

For additional product information, contact your local authorized dealer, local Thermo Scientific Orion technical sales representative or contact us using the Water and Laboratory Products (WLP) information on the page back of this manual.

### **Applications and Technical Resources**

Visit <u>www.thermoscientific.com/water</u> to view Thermo Scientific Orion products and download product literature, user guides and manuals, software updates, and the latest application and technical resources including the Thermo Scientific<sup>™</sup> Orion<sup>™</sup> Dual Star<sup>™</sup> meter technical notes for Ammonia in Water and Wastewater, Fluoride in Water and Wastewater, Nitrate in Water and Wastewater, and pH of Water and Wastewater. These technical notes provide step-by-step illustrated instructions for performing common measurements on the Orion Dual Star meter.

# **Table of Contents**

Section 1	6
Meter Introduction	6
Meter Overview	6
Redesigned Orion Dual Star Meters	7
Meter Packing List	7
Intended Use	7
Section 2	8
Meter Basics	8
Using the Universal Power Adapter	8
Attaching the Electrode Stand and Holder	9
Using the Electrode Holder	10
Powering the Meter On and Off	
Meter Maintenance	
Meter Components	
Meter Connections	12
Channel 1 and Channel 2 Inputs	13 12
Connecting Electrodes and Cables to the Meter	13 14
Meter Reypau	14 15
Alphanumenc Reypau Matar Display	10 16
	10
Section 3	18
Meter Setup	
EZ Startup Menu	
Setup Menu	22
Navigating the Setup Menu	22
Channel 1 and Channel 2 Setup Menus	23
Instrument Parameters Setup Menu	28
Autosampler Setup Menu	33
Incremental Techniques Setup Menu	33
GLP Setup Menu	
EZ Startup Setup Menu	
Section 4	34
Calibration and Measurement	
pH Calibration and Measurement	
pH Calibration Hints	
·	
pH Calibration Procedure (1 to 6 points)	35
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure	35 36
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure ISE Calibration and Measurement	35 36 37
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure ISE Calibration and Measurement ISE Calibration Hints	35 36 37 37
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure ISE Calibration and Measurement ISE Calibration Hints ISE Calibration Procedure (2 to 6 points)	
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure ISE Calibration and Measurement ISE Calibration Hints ISE Calibration Procedure (2 to 6 points) ISE Measurement Procedure	35 36 37 37 37 37 39
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure ISE Calibration and Measurement ISE Calibration Hints ISE Calibration Procedure (2 to 6 points) ISE Measurement Procedure mV Measurement and RmV and ORP Calibration and Measurement	35 36 37 37 37 39 40
pH Calibration Procedure (1 to 6 points) pH Measurement Procedure ISE Calibration and Measurement ISE Calibration Hints ISE Calibration Procedure (2 to 6 points) ISE Measurement Procedure mV Measurement and RmV and ORP Calibration and Measurement mV Measurement Procedure	35 36 37 37 37 39 40 40

RmV Calibration Procedure	41
RmV Measurement Procedure	42
ORP Calibration Hints	42
ORP Calibration Procedure	43
ORP Measurement Procedure	44
Temperature Calibration and Measurement	45
Temperature Calibration Procedure	45
Section 5	46
Methods	46
Saving the Current Meter Settings as a New Method	47
Loading an Existing Method	47
Editing or Printing an Existing Method	48
Coping an Existing Method	48
Deleting an Existing Method	49
Section 6	50
Incremental Techniques	50
Electrode ID and Slope for Incremental Techniques	
Single Known Addition	
Single Known Addition Procedure	52
Double Known Addition	54
Double Known Addition Procedure	54
Single Known Subtraction	56
Single Known Subtraction Procedure	56
Double Known Subtraction	58
Double Known Subtraction Procedure	58
Section 7	58 60
Section 7 Data Storage and Retrieval	58 60 60
Section 7 Data Storage and Retrieval Data Storage Settings	58 60 60
Section 7         Data Storage and Retrieval         Data Storage Settings         Viewing the Calibration Log and Data Log	58 60 60 62
Section 7         Data Storage and Retrieval         Data Storage Settings         Viewing the Calibration Log and Data Log         Viewing the Calibration Log	58 60 60 62 62
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.	58 60 60 62 62 62 62
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Viewing the Data Log.         Interfacing a Printer	58 60 60 60 62 62 62 62 62
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Mater Seftware Upgrade	58 60 60 62 62 62 62 64 64
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade.	58 60 60 62 62 62 62 64 65
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade         Section 8	58 60 60 62 62 62 64 65 66
Section 7 Data Storage and Retrieval Data Storage Settings. Viewing the Calibration Log and Data Log. Viewing the Calibration Log. Viewing the Data Log. Interfacing a Printer Interfacing a Computer Meter Software Upgrade Section 8 Autosampler Operation	58 60 60 62 62 62 64 65 66
Section 7 Data Storage and Retrieval Data Storage Settings Viewing the Calibration Log and Data Log Viewing the Calibration Log Viewing the Data Log Viewing the Data Log Interfacing a Printer Interfacing a Computer Meter Software Upgrade Section 8 Autosampler Operation Interfacing the Meter and Autosampler	58 60 60 62 62 62 62 64 65 66
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade.         Section 8         Autosampler Operation         Interfacing the Meter and Autosampler         Autosampler Setup Menu	
Section 7 Data Storage and Retrieval Data Storage Settings Viewing the Calibration Log and Data Log Viewing the Calibration Log Viewing the Data Log Viewing the Data Log Interfacing a Printer Interfacing a Computer Meter Software Upgrade Section 8 Autosampler Operation Interfacing the Meter and Autosampler Autosampler Setup Menu Running the Autosampler	
Section 7 Data Storage and Retrieval Data Storage Settings. Viewing the Calibration Log and Data Log. Viewing the Calibration Log. Viewing the Data Log. Interfacing a Printer Interfacing a Computer Meter Software Upgrade. Section 8 Autosampler Operation. Interfacing the Meter and Autosampler. Autosampler Setup Menu Running the Autosampler Autosampler Error Messages.	
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade.         Section 8         Autosampler Operation.         Interfacing the Meter and Autosampler.         Autosampler Setup Menu         Running the Autosampler Error Messages.         Section 9	
Double Known Subtraction Procedure         Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade.         Section 8         Autosampler Operation         Interfacing the Meter and Autosampler         Autosampler Setup Menu         Running the Autosampler         Autosampler Error Messages         Section 9	
Section 7         Data Storage and Retrieval         Data Storage Settings         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade.         Section 8         Autosampler Operation         Interfacing the Meter and Autosampler         Autosampler Setup Menu         Running the Autosampler Irror Messages         Section 9         Customer Services         Troubleshooting Guide	
Section 7         Data Storage and Retrieval         Data Storage Settings.         Viewing the Calibration Log and Data Log.         Viewing the Calibration Log.         Viewing the Data Log.         Interfacing a Printer         Interfacing a Computer         Meter Software Upgrade.         Section 8         Autosampler Operation         Interfacing the Meter and Autosampler         Autosampler Setup Menu         Running the Autosampler Error Messages         Section 9         Customer Services         Troubleshooting Guide         Electrode Stability Test	58 60 60 62 62 62 64 64 64 65 66 66 
Section 7 Data Storage and Retrieval Data Storage Settings. Viewing the Calibration Log and Data Log Viewing the Calibration Log. Viewing the Data Log. Interfacing a Printer Interfacing a Computer Meter Software Upgrade. Section 8 Autosampler Operation Interfacing the Meter and Autosampler. Autosampler Setup Menu Running the Autosampler Autosampler Error Messages. Section 9 Customer Services Troubleshooting Guide Electrode Stability Test. Meter Self Test and Checkout.	

User Reset	74
Factory Configuration	74
About Meter	74
Electrode Condition Icon	75
Assistance	75
Warranty	75
Meter Error Codes	76
Over Range Measurements	76
Notice of Compliance	78
WEEE Compliance	78
Declaration of Conformity	79
Real Time Clock Battery Maintenance	80
Replacing and Disposing of the Battery	80
Ordering Information	81
Meter Specifications	84
Section 10	
Advanced Meter Features	
Multipoint Calibration Curve Feature	
Automatic Blank Correction Feature	
Error Code E306	
Isopotential Point Feature	
Determining an Isopotential Point	89
Automatic pH Buffer Recognition Feature	90
Testing a pH Electrode for Automatic Buffer Recognition	90
5	

# SECTION 1 Meter Introduction

# Meter Overview

This user manual contains information on the preparation, operation and maintenance of the Thermo Scientific Orion Dual Star pH/ISE benchtop meter. The Orion Dual Star meter has two measuring channels and is capable of measuring pH, ion concentration (ISE), raw millivolts (mV), relative millivolts (RmV), oxidation-reduction potential (ORP) and temperature.

### **Display Features**

The large, backlit LCD display on the Orion Dual Star meter shows both measuring channels simultaneously or each measuring channel separately. Step-by-step instructions and prompts are shown on the meter to assist operators with the setup, calibration and autosampler menus. A calibration graph, which includes the electrode slope between points and the average slope, can be viewed in the calibration mode.

### **Data Features**

The Orion Dual Star meter can save up to 10 calibrations per channel and 1000 data points in the meter data log. A graph of the calibration points with slope can be viewed on the meter display and printed. Calibration and measurement data can be exported from the meter to a printer or computer. The meter saves up to 10 methods per channel, so switching the meter to different testing methods is fast and simple.

### **Troubleshooting Features**

Troubleshooting procedures are available in the meter setup menu and include a meter stability test, meter self-test and meter reset functions.

### **Redesigned Orion Dual Star Meters**

Redesigned Orion Dual Star pH/ISE benchtop meters will be available beginning in late June 2015. These Orion Dual Star meters offer a redesigned housing that matches the appearance of our Thermo Scientific<sup>™</sup> Orion Star<sup>™</sup> A series benchtop meters and Thermo Scientific<sup>™</sup> Orion<sup>™</sup> Versa Star<sup>™</sup> benchtop meters.

Additional benefits of redesigned Orion Dual Star meters include the new housing that is compatible with the meter-attachable electrode stand (Cat. No. STARA-BEA) and this stand will be included with all redesigned Orion Dual Star meters and new Dual Star Com communication software.

The Orion Dual Star meter operation workflows, functions, user interface, specifications and catalog numbers are the same between current and redesigned Orion Dual Star meters.

# Meter Packing List

Thermo Scientific Orion Dual Star pH/ISE benchtop meters include:

- Orion Dual Star meter with two BNC shorting caps (Cat. No. 090045)
- Meter-attachable electrode stand and holder, Cat. No. STARA-BEA
- Universal power adapter, Cat. No. 1010003
- Literature CD
- Printed quick start guide
- Computer interface cable
- Meter test certificate

Each Orion Dual Star meter is fully tested and certified by Thermo Fisher Scientific and a certificate of calibration is included with each meter.

For specific meter and kit contents, refer to the <u>Ordering Information</u> section of this user manual and select from meter kits that include Thermo Scientific<sup>™</sup> Orion<sup>™</sup> ROSS Ultra<sup>™</sup> pH electrodes, Thermo Scientific<sup>™</sup> Orion<sup>™</sup> ROSS Ultra<sup>™</sup> Triode<sup>™</sup> pH/ATC electrodes Thermo Scientific<sup>™</sup> Orion<sup>™</sup> ROSS<sup>™</sup> Sure-Flow<sup>™</sup> pH electrodes and Thermo Scientific<sup>™</sup> Orion<sup>™</sup> ROSS<sup>™</sup> solutions.

To download free Orion Dual Star Com data transfer computer software and the Dual Star USB driver, please visit our website at <u>www.thermoscientific.com/OrionMeters</u>.

# Intended Use

Please read this reference guide thoroughly. Any use outside of these instructions may invalidate the meter warranty and cause permanent damage to the meter.



# SECTION 2 Meter Basics

# Using the Universal Power Adapter

A universal power adapter (Catalog Number 1010003) with US, EU, UK and China plug plates is included with the Orion Dual Star pH/ISE benchtop meters. This universal power adapter is specifically for use with Orion Dual Star benchtop meters. Use of other power adapters can damage the meter and void the warranty.

- 1. Select the appropriate plug plate for the power outlet that will be used.
- 2. Remove the clear plastic cover from the groove on the back of the power adapter.
- 3. Slide the appropriate plug plate into the groove on the back of the power adapter.



4. Connect the assembled power adapter to a power outlet and the meter input labeled Power. A surge protector or uninterrupted power supply (UPS) is also recommended.

# Attaching the Electrode Stand and Holder

Redesigned Orion Dual Star meters are compatible with the meter-attachable electrode stand (Cat. No. STARA-BEA) shown below. The electrode stand can be attached to either side of the meter and up to two stands can be attached to each meter. A weighted base (Catalog Number STARA-HB) is also available to support the stand without attachment to the meter. Older style Orion Dual Star meters can be used with this electrode stand, but require additional purchase of the weighted base (Catalog Number STARA-HB).

1. Open the box containing the electrode stand. The box will include a base plate, electrode arm, pivot pin, electrode holder, ATC spacer and RDO clip.



- 2. Turn the meter over, with the meter display facing down, on a clean dry surface.
- 3. Identify the side of the meter that the stand will be installed on and remove the screw between the circles on that side of the meter.
- 4. Align the base plate of the stand with the circles on the meter.
- 5. Replace the screw from step 3 to attach the base plate to the meter.
- 6. Turn the meter over, with the meter display facing up.
- 7. Insert the electrode arm into the metal post on the base plate.
- Connect the electrode holder to the electrode arm using the pivot pin.



### **Using the Electrode Holder**

Place electrodes in the stand for easy movement in and out of containers during calibration, sample measurement and storage. Recommended electrode positions are shown below.



# Powering the Meter On and Off

- 1. If the meter is not connected to a wall outlet via the universal power adapter, select the appropriate wall outlet plug and slide the plug plate it into the groove on the power adapter. Connect the power adapter to the meter and wall outlet.
- 2. Press and hold the *power* key for about three seconds to turn on the meter.
- 3. When the meter is on, press and release the *power* key to turn the display backlight off or to vary the intensity of the backlight. There are five levels of intensity that the *power* key can be used to select.
- To turn off the meter, press and hold the *power* key for about three seconds, until the meter display is blank.

# Meter Maintenance

For routine meter maintenance, dust and wipe the meter with a damp cloth. If necessary, warm water or a mild water-based detergent can be used. Meter maintenance can be performed on a daily, weekly or monthly basis, as required by the environment in which the meter is operated.

Immediately remove any spilled substance from the meter using the proper cleaning procedure for the type of spill.

# Meter Components



### **Meter Connections**



**BNC** – Input for pH electrode, ion selective electrode (ISE) or ORP/redox electrode with BNC or waterproof BNC connectors.

Ref. – Input for reference electrodes with standard 2.5 mm pin-tip connectors.

ATC - Input for ATC probes with 8 pin MiniDIN connectors.

**Stirrer** – Input for the Thermo Scientific<sup>™</sup> Orion Star<sup>™</sup> series stirrer probe, Catalog Number 096019. The stirrer probe can be purchased separately as a meter accessory.

USB - Input for the USB cable, used to connect the meter to a computer.

RS232 – Input for the RS232 cable, used to connect the meter to a printer or computer.

Ground – Input for a ground wire with a standard pin-tip connector.

Power - Input for the universal power adapter that is included with the meter.

### **Channel 1 and Channel 2 Inputs**

The **BNC**, **Ref.**, and **ATC** connections are labeled as Channel 1 or Channel 2 on the ridge above the connections. Using the meter orientation shown below, the channel 2 connections are on the left and the channel 1 connections are on the right.



### **Connecting Electrodes and Cables to the Meter**

- Connect the pH electrode, ion selective electrode (ISE) and/or ORP/redox electrodes to the BNC inputs on the meter. Attach a BNC connector to the meter input by sliding the connector onto the input, pushing down on the connector and turning the connector clockwise to lock it into position.
- Connect the reference electrodes to the reference inputs on the meter.
   Note: If a combination electrode is used, a reference electrode is not needed.
- 3. Connect the ATC probes to the 8 pin MiniDIN inputs on the meter. The ATC connector has a clip that locks the connection and prevents it from detaching from the meter.
- 4. Connect the stirrer probe, Catalog Number 096019, to the stirrer input on the meter.
- 5. To connect a printer to the meter, connect an RS232 printer cable to the RS232 inputs on the meter and printer. The Orion Star series printer, Catalog Number 1010006, is recommended for use with this meter and includes the appropriate RS232 printer cable.
- To connect a computer to the meter, connect the USB computer cable to the USB inputs on the meter and computer or connect the RS232 computer cable, Catalog Number 1010053, to the RS232 inputs on the meter and computer.

### **Meter Keypad**



Press the shortcut keys in the measurement mode to move quickly to different meter menus. Press the  $\blacktriangle$  or  $\checkmark$  key to scroll through a list of options on the display. The  $\blacktriangle$  or  $\checkmark$  key will loop through a list, so the meter returns to the first item on the list after scrolling past the last item on the list. Use the numeric keypad, **decimal** key and +/- key when entering a number.

Кеу	Description
	To turn the meter on, press and hold down the <i>power</i> key for about three seconds until the meter turns on.
¢	When the meter is on, press and quickly release the <i>power</i> key to turn the display backlight off or to vary the intensity of the backlight.
	To turn the meter off, press and hold down the <i>power</i> key for about three seconds until the meter turns off.
h 12 f3	Press the <i>f1</i> , <i>f2</i> and <i>f3</i> function keys to perform the action shown above each key on the display. The <i>f1</i> , <i>f2</i> and <i>f3</i> keys have menudependent functions.
channel	Press the <i>channel</i> key to display channel 1 only, channel 2 only or a split screen with channel 1 (top) and channel 2 (bottom).
(neasure) (esc)	Press the <i>measure (esc)</i> key to return to the meter to the measurement mode. The <i>measure (esc)</i> key can be used as an escape key in most meter menus.
$\smile$	Press to enter a value of one (1) when using the numeric keypad.
2	Press the $\blacktriangle$ key to scroll up through lists of items on the display. The $\blacktriangle$ key will to loop through a list, so the meter will go to the last item on the list after scrolling past the first item on the list.
	Press to enter a value of two (2) when using the numeric keypad.
3 ▼	Press the $\checkmark$ key to scroll down through lists of items on the display. The $\checkmark$ key will loop through a list, so the meter will go to the first item on the list after scrolling past the last item on the list.
$\smile$	Press to enter a value of three (3) when using the numeric keypad.

(log/print)	Press the <i>log/print</i> key to manually log, print or log and print a measurement. Press to enter a decimal place (.) when using the numeric keypad.
4	Press the <i>cal</i> key to enter the calibration mode.
cal	Press to enter a value of four ( <i>4</i> ) when using the numeric keypad.
5	Press the <i>mode</i> key to change the measurement mode for channel 1 or channel 2.
mode	Press to enter a value of five ( <i>5</i> ) when using the numeric keypad.
6 method	Press the <i>method</i> key to access the methods list for each channel, existing methods can be selected and run from the list. Press to enter a value of six (6) when using the numeric keypad.
+/- log view	Press the <i>log view</i> key to view the data log and calibration log. Press to make a value positive or negative (+/-) when using the numeric keypad.
(7	Press the <i>setup</i> key to enter the setup menu.
setup	Press to enter a value of seven (7) when using the numeric keypad.
8	Press the <i>res</i> key to change the displayed measurement resolution for channel 1 or channel 2.
res	Press to enter a value of eight ( <i>8</i> ) when using the numeric keypad.
9	Press the <i>auto</i> key to access the autosampler setup menu.
auto	Press to enter a value of nine (9) when using the numeric keypad.
0	Press the <i>stirrer</i> key to turn on and off the stirrer probe.
stirrer	Press to enter a value of zero (0) when using the numeric keypad.

### **Alphanumeric Keypad**

Use the alphanumeric keypad when entering the electrode serial number, sample ID number and passwords. The alphanumeric keypad works in the same way as a cell phone keypad. For example, press the 2 key to enter 2, A, B or C. Enter the letter B by pressing the 2 key three times  $(2 \rightarrow A \rightarrow B)$ . Uppercase or lowercase letters can be entered by pressing the same key. For example, enter the letter a by pressing the 2 key five times  $(2 \rightarrow A \rightarrow B \rightarrow C \rightarrow a)$ . The meter will automatically advance to the next digit when no keys are pressed for three seconds.



### **Meter Display**

**Dual Channel Meter Display** 



**Single Channel Meter Display** 



Display Icon	Description
<b>Ch.2</b> Channel Number – Indicates that channel number that corresponds with the displayed measurement. Press the <i>channel</i> key to scroll through a dual channel display for channel 1 and 2, single display for channel 1 and a single display for channel 2.	
F"	Electrode ID – Specifies that type of electrode that was selected by the operator. To change the electrode ID, refer to the <u>Setup Menu</u> section.
s/n:ABCDE1234	Electrode Serial Number (Single Channel Display Only) – Shows the operator assigned electrode serial number. To change the electrode serial number, refer to the <u>Setup Menu</u> section.
1.00	Measurement Value – Displays the current electrode reading based on the last saved calibration.

Display Icon	Description	
	Calibration Warning – Indicates that the electrode is calibrated. If the icon is flashing, it indicates that the electrode needs to be calibrated. Calibrate the electrode at least every twelve hours to prevent the icon from flashing.	
fL.	Electrode Condition – Represents the performance of the electrode, based on the last saved calibration and electrode measurement stability. If the electrode icon h two lines, the electrode condition is good. If the electrode icon has one line, the electrode condition is fair. If the electrode icon has a slash through it, the electrode condition is bad. Refer to the <u>Troubleshooting Guide</u> section for more information	
<b>M</b> 201	Method Number – Shows the method that is currently being used on that channel. To change the method number, refer to the <u>Methods</u> section.	
AR	Auto-Read ( <b>AR</b> ) Read Type Indicator – If Auto-Read was selected as the read type, the <b>AR</b> icon will appear when the reading is stable and the display has been locked. To change the read type, refer to the <u>Setup Menu</u> section.	
ready	Stabilizing/Ready Read Type Indicator – Specifies the measurement stability of the electrode. To change the read type, refer to the <u>Setup Menu</u> section.	
ppm	Measurement Mode – Indicates the unit of measurement for the measurement value and determines the type of calibration that will be performed. To change the measurement mode, refer to the <u>Setup Menu</u> section. If the ISE measurement mode is changed, the electrode must be re-calibrated with new standards that have the same measurement unit as the measurement mode.	
<b>25.0</b> ℃	Temperature Measurement – Displays the current temperature based on the selected ATC probe reading or entered temperature value. To change the temperature display to Celsius or Fahrenheit, refer to the <u>Setup Menu</u> section.	
ATC2	Temperature Source – Shows the origin of the current temperature measurement. To change the source of the temperature measurement, refer to the <i>Setup Menu</i> section.	
<b>0.0</b> mV	mV Measurement – Shows the raw millivolt reading of the electrode.	
ē	Printer Interface – Indicates if a printer is interfaced with the meter.	
USB	Computer Interface – Indicates if a computer is interfaced with the meter. The USB icon will only be shown when the meter is connected to a computer via USB input.	
$\square$	Timed Interval Read Type Indicator – Indicates when a timed reading is being taken by the meter. To change the read type, refer to the <u>Setup Menu</u> section.	
	Autosampler Interface – Indicates if the meter is running the autosampler.	
15:45 05/25/2009	Time and Date (Single Channel Display Only) – Displays the current time and date settings. To change the time and date, refer to the <u>Setup Menu</u> section.	
回 12345ABCD	Sample ID (Single Channel Display Only) – Shows the operator assigned sample ID number. To change the sample ID number, refer to the <u>Setup Menu</u> section.	
log/print	f1 Function – Displays the action that will be done when the <b>f1</b> key is pressed.	
cal	f2 Function – Displays the action that will be done when the <b>f2</b> key is pressed.	
setup	f3 Function – Displays the action that will be done when the <b>f3</b> key is pressed.	



# SECTION 3 Meter Setup

# EZ Startup Menu

It is highly recommended that the EZ Startup menu be completed the first time that the meter is used. The EZ Startup menu sets important meter parameters, such as the displayed language, date and time, measurement mode and read type for each channel and data output settings. To access the EZ Startup menu from the measurement mode, press the *setup* key, press the  $\land / \blacksquare$  keys to highlight EZ Startup and press the *f2 (select)* key.

- Press the ▲ /▼ keys to highlight English, Español, Deutsch, Français or Italiano as the language that will be displayed on the meter. Press the f2 (select) key to save the setting and proceed to the next EZ Startup menu.
- 2. Set the time, date and date format.
  - a. To set the time, press the ▲ /▼ keys to highlight Time and press the *f*2 (*select*) key. Use the numeric keypad to set the time in 24 hour format and press the *f*2 (*accept*) key.
  - b. To set the date, press the ▲ / ▼ keys to highlight Date and press the *f*2 (select) key. Use the numeric keypad to set the date (the date format is shown below the date setting) and press the *f*2 (accept) key.
  - c. To set the date format, press the ▲ /▼ keys to highlight Date Format and press the f2 (select) key. Press the ▲ /▼ keys to highlight MM/DD/YYYY to display the date as month/day/year or DD/MM/YYYY to display the date as day/month/year and press the f2 (accept) key.
  - d. Once all of the time and date settings are entered, press the *f*3 (*next*) key to save the settings and proceed to the next EZ Startup menu.

- 3. Press the ▲ /▼ keys to highlight pH, ISE (ion concentration), mV (raw millivolts), RmV (relative millivolts) or ORP (oxidation-reduction potential) as the measurement mode for channel 1 and press the *f2 (accept)* key. The meter will request information based on the selected measurement mode. If channel 1 will not be used, press the *f3 (no)* key to skip to the next EZ Startup menu.
  - a. ISE only: Press the ▲ /▼ keys to highlight the displayed ISE units as ppm (parts per million), mg/L (milligrams per liter, % (percent), M (moles per liter) or unitless (no units) and press the f2 (accept) key.
  - b. ISE, RmV and mV only: Press the ▲ /▼ keys to highlight the electrode ID and press the *f*2 (accept) key. The electrode ID can be set as pH, ORP, fluoride (F<sup>-</sup>), ammonia (NH<sub>3</sub>), ammonium (NH<sub>4</sub><sup>+</sup>), nitrate (NO<sub>3</sub><sup>-</sup>), nitrite (NO<sub>2</sub><sup>-</sup>), sulfide (S<sup>-2</sup>), chloride (CI<sup>-</sup>), chlorine (Cl<sub>2</sub>), bromide (Br<sup>-</sup>), iodide (I<sup>-</sup>), cyanide (CN<sup>-</sup>), sodium (Na<sup>+</sup>), potassium (K<sup>+</sup>), calcium (Ca<sup>+2</sup>), silver (Ag<sup>+</sup>), copper (Cu+2), lead (Pb+2), cadmium (Cd+2), perchlorate (ClO4<sup>-</sup>), fluoroborate (BF<sub>4</sub><sup>-</sup>), thiocyanate (SCN<sup>-</sup>), nitrogen oxide (NO<sub>x</sub>), carbon dioxide (CO<sub>2</sub>), oxygen (O<sub>2</sub>), monovalent cation (X<sup>+</sup>), divalent cation (X<sup>+2</sup>), monovalent anion (X<sup>-</sup>) or divalent anion (X<sup>-2</sup>). Some of the electrode ID options may not be accessible, depending on the measurement mode that was selected for the channel.
  - c. Press the *f2 (yes)* key to enter an electrode serial number or press the *f3 (no)* key to skip this step. To enter an electrode serial number, use the alphanumeric keypad to enter letters and/or numbers. The meter will automatically advance to the next digit when no keys are pressed for three seconds. Press the *f3 (clear)* key to delete any of the digits. Once the serial number is entered, press the *f2 (accept)* key. Refer to the meter pullout guide for the location of the letter keys on the numeric keypad.
  - d. **pH and ISE only:** Press the ▲ / ▼ keys to highlight the measurement resolution and press the *f*2 (*accept*) key to save the setting.
  - e. To change the displayed temperature unit to degree Celsius (°C) or degree Fahrenheit (°F), press the ▲ /▼ keys to highlight Temperature Unit and press the *f2 (accept)* key. Press the ▲ /▼ keys to highlight the desired temperature units and press the *f2 (accept)* key.
  - f. Press the ▲ /▼ keys to highlight ATC1, ATC2 or Manual as the source of the temperature compensation that will be used for measurements and press the *f2 (accept)* key to save the setting. Automatic temperature compensation can be performed by connecting an ATC probe to ATC input 1 (ATC1) or ATC input 2 (ATC2) on the meter. If manual temperature compensation is selected, enter the temperature using the numeric keypad and the *decimal* key. Once the temperature value has been entered, press the *f2 (accept)* key.
  - g. Press the ▲ /▼ keys to highlight Auto-Read, On Ready, At Time Intervals, Continuous, Value Change In Measurement or Timed Reading as the read type and press the *f*2 (accept) key to save the setting.

- i. Auto-Read The meter will display the measurement as it stabilizes and lock and hold the measurement when it is stable. Once a measurement is locked, press the *measure* key to take a new measurement. Once the measurement is locked, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.
- ii. On Ready The meter will show stabilizing on the display until the measurement is stable. When the measurement is stable, the meter will show ready on the display. The display will automatically update if the measurement changes and the measurement will not lock. Once ready is shown on the display, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.
- iii. At Time Intervals The meter will read and display a measurement at the set time interval that is programmed by the operator. At the set time interval, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. If At Time Intervals is selected, enter the time interval value in a hours : minutes : seconds format using the numeric keypad and press the f2 (accept) key.
- iv. Continuous The meter will continuously measure and update the display. No indicator will be shown on the display. This read type is useful when performing an experiment that requires continuous measurements to be taken, regardless of the measurement stability. Press the *log/print* key to export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.
- v. Value Change In Measurement The meter will display the measurement when the reading reaches or exceeds the set high or low value that is programmed by the operator. When the reading reaches or exceeds the set high or low value, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. If Value Change In Measurement is selected, press the ▲ /▼ keys to highlight Min Value, enter the minimum measurement value using the numeric keypad and the *decimal* key, press the ▲ /▼ keys to highlight Max Value, enter the maximum measurement value using the numeric keypad and the *decimal* key.
- vi. Timed Reading The meter will display a measurement after a set time delay and lock and hold the measurement after the time delay is reached. To start another time delay cycle, press the *measure* key. Once the measurement is locked, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is

connected to the meter and enabled in the setup menu. If Timed Reading is selected, enter the time delay value in a minutes : seconds format using the numeric keypad and press the *f*2 (*accept*) key.

- 4. Repeat step 3 for channel 2.
- Press the ▲ /▼ keys to highlight Printer, Computer or None as where the meter will export the measurement and calibration data and press the *f2 (accept)* key. If Printer or Computer is selected, the meter will prompt the operator to set the baud rate. Press the ▲ /▼ keys to highlight 1200, 2400, 4800, 9600, 19200 or 38400 as the baud rate and press the *f2 (accept)* key.
- 6. Press the *f*2 (*yes*) key to enable the exporting of the measurement data to the data log or press the *f*3 (*no*) key to disable the exporting of the measurement data to the data log.
- 7. Press the *f*2 (*measure*) key to save the EZ Startup menu settings and proceed to the measurement mode or press the *f*1 (*back*) key to edit any of the EZ Startup settings.

# Setup Menu



The setup menu can be accessed in the measurement mode by pressing the *setup* key or by pressing the *f3 (setup)* key in the dual channel display mode.

The setup menu contains seven submenus: Channel 1, Channel 2, Instrument Parameters, Autosampler, Incremental Techniques, GLP and EZ Startup. Settings for each menu can be selected and saved in the setup menu.

### **Navigating the Setup Menu**

- 1. In the setup menu, press the ▲ / ▼ keys to highlight the desired submenu and press the *f*2 (*select*) key to view the menu options.
- Press the ▲ / ▼ keys to highlight the desired menu option and press the f2 (select) key to view the current setting and change the selected setting, if required.
- If a new setting must be selected from a list of options, press the ▲ / ▼ keys to highlight the desired setting and press the *f2 (accept)* key. If a new numeric value must be entered, use the numeric keypad, *decimal* key and +/- key as required and press the *f2 (accept)* key.
- 4. Press the f1 (back) key to return to the list of setup submenus.
- 5. Press the *f1* (back) key to return to the measurement mode.

### **Channel 1 and Channel 2 Setup Menus**

**Note:** Channel 1 and Channel 2 have the same options. However, the options must be set separately for each channel.

Channel 1 Setur	)	Channel 2 Setu	ρ
Measure Mode	pН	Measure Mode	ppm
Electrode ID	pH	Electrode ID	F-
<ul> <li>Electrode Serial Number</li> </ul>	ABCDE1234	<ul> <li>Electrode Serial Number</li> </ul>	ABCDE1234
Resolution	0.001	Resolution	1.00
<ul> <li>Temperature Calibration</li> </ul>		<ul> <li>Temperature Calibration</li> </ul>	
<ul> <li>Temperature Input</li> </ul>	ATC1	<ul> <li>Temperature Input</li> </ul>	ATC2
<ul> <li>Calibration Setup</li> </ul>		<ul> <li>Calibration Setup</li> </ul>	
Read Type	Auto-Read	Read Type	Auto-Read
• Alarm	Off	Alarm	Off
Sample ID	Manual	Sample ID	Manual
Method	M100	Method	M200
back select		back select	

In the measurement mode, press the setup key to enter the setup menu. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Channel 1 or Channel 2 and press the *f2 (select)* key. Perform the following actions, as required.

### **Measure Mode**

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Measure Mode and press the *f2 (select)* key. Press the  $\land$  /  $\checkmark$  keys to highlight pH, ISE, mV, RmV or ORP as the measurement mode and press the *f2 (accept)* key. If ISE is selected, press the  $\land$  /  $\checkmark$  keys to highlight ppm (parts per million), mg/L (milligrams per liter, % (percent), M (moles per liter) or unitless (no units) as the ISE measurement unit that will be displayed and press the *f2 (accept)* key.

### Electrode ID

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Electrode ID and press the *f2 (select)* key. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight the type of electrode and press the *f2 (accept)* key. The type of electrode can be set as pH, ORP, fluoride (F<sup>-</sup>), ammonia (NH<sub>3</sub>), ammonium (NH<sub>4</sub><sup>+</sup>), nitrate (NO<sub>3</sub><sup>-</sup>), nitrite (NO<sub>2</sub><sup>-</sup>), sulfide (S<sup>-2</sup>), chloride (Cl<sup>-</sup>), chlorine (Cl<sub>2</sub>), bromide (Br<sup>-</sup>), iodide (l<sup>-</sup>), cyanide (CN<sup>-</sup>), sodium (Na<sup>+</sup>), potassium (K<sup>+</sup>), calcium (Ca<sup>+2</sup>), silver (Ag<sup>+</sup>), copper (Cu<sup>+2</sup>), lead (Pb<sup>+2</sup>), cadmium (Cd<sup>+2</sup>), perchlorate (ClO<sub>4</sub><sup>-</sup>), fluoroborate (BF<sub>4</sub><sup>-</sup>), thiocyanate (SCN<sup>-</sup>), nitrogen oxide (NO<sub>x</sub>), carbon dioxide (CO<sub>2</sub>), oxygen (O<sub>2</sub>), monovalent cation (X<sup>+</sup>), divalent cation (X<sup>+2</sup>), monovalent anion (X<sup>-</sup>) or divalent anion (X<sup>-2</sup>). Some or all of the electrode ID options may not be accessible, depending on the selected measurement mode for the channel.

### **Electrode Serial Number**

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Electrode Serial Number and press the *f2 (select)* key. Use the alphanumeric keys to enter a unique number and letter combination for the electrode. The meter will automatically advance to the next digit when no keys are pressed for three seconds. Press the *f3 (clear)* key to delete any of the numbers or letters. Once the serial number is entered, press the *f2 (accept)* key. Refer to the meter pullout guide for the location of the letter keys on the numeric keypad.

### Resolution

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Resolution and press the *f2 (select)* key. Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight the desired measurement display resolution and press the *f2 (accept)* key. This menu may not be accessible, depending on the measurement mode that was selected for the channel.

### **Temperature Calibration**

**Note:** The meter ATC temperature display has a relative accuracy of  $\pm 0.1$  °C. ATC probes have varying temperature accuracies, usually  $\pm 0.5$  °C to  $\pm 2$  °C. Use this function only if it is necessary to calibrate the temperature readings. Since the temperature offset calculated during the calibration is applied to all future temperature measurements, recalibrate if a different ATC probe is used.

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Temperature Calibration and press the *f2 (select)* key. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight ATC1, ATC2 or ATC1 & ATC2 to select the ATC probe(s) that will be calibrated and press the *f2 (select)* key. Place the ATC probe(s) and thermometers into a solution with a known, stable temperature. It is recommended that two NIST traceable thermometers be used to measure and verify the temperature of the solution. Wait for the readings of the ATC probe(s) and thermometers to stabilize (about five to ten minutes); use the numeric keypad, *decimal* key and +/- key to enter the temperature value read by the thermometer in the Reference Temperature: \_\_\_\_\_ °C field and press the *f2 (accept)* key. The meter will display the original temperature read by the ATC probe (i.e. ATC1: 24.2°C) and the temperature offset value that will be applied to all future temperature readings (i.e. ATC1 Offset: -0.8°C). Press the *f3 (log/print)* key to save and end the calibration, export the data to the calibration log, export the data to a printer or computer if one is connected to the meter and enabled in the setup menu, and proceed to the measurement mode or press the *f1 (escape)* key to return to the measurement mode without saving the calibration.

### Temperature Input

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Temperature Input and press the *f2 (select)* key. Select the source of the temperature measurement and the display units of the temperature measurement.

- ATC1, ATC2 or Manual To select the temperature source, press the ▲ /▼ keys to highlight ATC1, ATC2 or Manual and press the *f2 (accept)* key. Channel 1 and 2 are able to receive temperature readings from an ATC probe connected to the channel 1 input (ATC1) or the channel 2 input (ATC2). If Manual is selected, enter the temperature using the numeric keypad, *decimal* key and +/- key and press the *f2 (accept)* key.
- Temperature Unit To select the temperature units as Celsius or Fahrenheit, press the ▲ /▼ keys to highlight Temperature Unit, press the *f2 (accept)* key, press the ▲ /▼ keys to highlight Celsius or Fahrenheit and press the *f2 (accept)* key.

Once all of the temperature settings are entered, press the f1 (back) key.

### **Calibration Setup**

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Calibration Setup and press the *f2 (select)* key. View and change the pH or ISE calibration settings, depending on the measurement mode that was selected for the channel.

- pH Calibration Settings Entering the Calibration Setup menu will access the Buffer Set and ISO settings.
- Buffer Set Press the ▲ /▼ keys to highlight Buffer Set and press the *f*2 (*select*) key. Press the ▲ /▼ keys to select US or EURO as the calibration buffer set and press the *f*2 (*accept*) key. The meter will automatically recognize pH 1.68, 4.01, 7.00, 10.01 and 12.46 buffers if US is selected and pH 1.68, 4.01, 6.86 and 9.18 buffers if EURO is selected.
- ISO Press the ▲ /▼ keys to highlight ISO and press the *f*2 (*select*) key. To enter the pH isopotential point, use the numeric keypad and decimal key to enter the value and press the *f*2 (*accept*) key.

Once all of the pH calibration settings are entered, press the f1 (back) key.

- ISE calibration settings Entering the Calibration Setup menu will access the Autoblank, Low Level Stability and Isopotential settings.
  - Autoblank Press the ▲ /▼ keys to highlight Autoblank and press the *f*2 (select) key. Press the ▲ /▼ keys to highlight Off or On and press the *f*2 (accept) key.
  - Low Level Stability Press the ▲ /▼ keys to highlight Low Level Stability and press the *f2 (select)* key. Press the ▲ /▼ keys to highlight Off or On and press the *f2 (accept)* key.
  - Isopotential Press the ▲ /▼ keys to highlight Isopotential and press the f2 (select) key.
  - To turn off or on temperature compensation for ISE measurements, press the ▲ /▼ keys to highlight ISO On/Off and press the *f*2 (*select*) key. Press the ▲ /▼ keys to highlight Off or On and press the *f*2 (*accept*) key.
  - o To enter the ISE isopotential point, press the ▲ /▼ keys to highlight ISO Value and press the *f2* (*select*) key. Use the numeric keypad and decimal key to enter the value and press the *f2* (*accept*) key.
  - Once all of the ISE isopotential settings are entered, press the f1 (back) key.

Once all of the ISE calibration settings are entered, press the f1 (back) key.

### Read Type

Press the ▲ /▼ keys to highlight Read Type and press the *f*2 (*select*) key. Press the ▲ /▼ keys to highlight Auto-Read, On Ready, At Time Intervals, Continuous, Value Change In

Measurement or Timed Reading as the read type that will be used to display the measurements for the selected channel and press the *f*2 (*select*) key.

- Auto-Read The meter will display the measurement as it stabilizes and lock and hold the
  measurement when it is stable. Once a measurement is locked, press the measure key to
  take a new measurement. Once the measurement is locked, the meter will automatically
  export the measurement to the data log, if the data log is enabled in the setup menu, and
  to a printer or computer, if a printer or computer is connected to the meter and enabled in
  the setup menu.
- On Ready The meter will show stabilizing on the display until the measurement is stable. When the measurement is stable, the meter will show ready on the display. The display will automatically update if the measurement changes and the measurement will not lock. Once ready is shown on the display, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.
- At Time Intervals The meter will read and display a measurement at the set time interval that is programmed by the operator. At the set time interval, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. If At Time Intervals is selected, enter the time interval value in a hours : minutes : seconds format using the numeric keypad and press the *f2 (accept)* key.
- Continuous The meter will continuously measure and update the display. No indicator will
  be shown on the display. This read type is useful when performing an experiment that
  requires continuous measurements to be taken, regardless of the measurement stability.
  Press the log/print key to export the measurement to the data log, if the data log is enabled
  in the setup menu, and to a printer or computer, if a printer or computer is connected to the
  meter and enabled in the setup menu.
- Value Change In Measurement The meter will display the measurement when the reading reaches or exceeds the set high or low value that is programmed by the operator. When the reading reaches or exceeds the set high or low value, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. If Value Change In Measurement is selected, press the ▲ /▼ keys to highlight Min Value, enter the minimum measurement value using the numeric keypad and the *decimal* key, press the ▲ /▼ keys to highlight Max Value, enter the maximum measurement value using the numeric keypad and the *decimal* key.
- Timed Reading The meter will display a measurement after a set time delay and lock and hold the measurement after the time delay is reached. To start another time delay cycle, press the measure key. Once the measurement is locked, the meter will automatically export the measurement to the data log, if the data log is enabled in the setup menu, and

to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. If Timed Reading is selected, enter the time delay value in a minutes : seconds format using the numeric keypad and press the *f2* (accept) key.

### Alarm

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Alarm and press the *f2 (select)* key. View and change the alarm settings, if required.

- Alarm Press the ▲ /▼ keys to highlight Alarm and press the *f*2 (*select*) key. Press the ▲ /▼ keys to highlight Off, Low Limit, High Limit or Both Low/High and press the *f*2 (*select*) key.
- Parameter Press the ▲ /▼ keys to highlight Parameter and press the *f2 (select)* key.
   Press the ▲ /▼ keys to highlight pH, ISE-ppm, ISE-mg/L, ISE-%, ISE-M, ISE-unitless, mV, RmV, ORP or °C (°F will be shown instead of °C if it is the selected temperature unit) and press the *f2 (accept)* key.
- Low Limit Press the ▲ /▼ keys to highlight Low Limit and press the f2 (select) key. To
  enter a low limit alarm value, use the numeric keypad and decimal key to enter the low
  value and press the f2 (accept) key.
- High Limit Press the ▲ /▼ keys to highlight High Limit and press the *f2 (select)* key. To enter a high limit alarm value, use the numeric keypad and decimal key to enter the high value and press the *f2 (accept)* key.

**Note:** The read type must be set to continuous to use the alarm settings.

### Sample ID

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Sample ID and press the *f2* (select) key. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Off, Manual or Auto Incremental and press the *f2* (accept) key. If Auto Incremental is selected, use the alphanumeric keypad to enter the initial sample ID and press the *f2* (accept) key. The meter will save the entered sample ID and add a sequential number after it.

### Method

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Method and press the *f*2 (*select*) key. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight the desired method number, press the *f*2 (*load*) key and press the *f*2 (*yes*) key. To view a method, edit a method, copy an existing method to a new method, delete a method, print a method or save the current meter settings as a new method, refer to *Chapter 5 Methods*.

### **Instrument Parameters Setup Menu**



In the measurement mode, press the setup key to enter the setup menu. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Instrument Parameters and press the *f2 (select)* key. Perform the following actions, as required.

### Export Data (PC/Printer/Log)

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Export Data (PC/Printer/Log) and press the *f2* (select) key. View and change the data exporting settings, if required.

- Export Trigger Press the ▲ /▼ keys to highlight Export Trigger and press the f2 (select) key. Press the ▲ /▼ keys to highlight Channel 1, Channel 2 or Channel 1 & 2 and press the f2 (accept) key. The read type of the selected channel(s) will determine when the meter sends the measurement data to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.
- PC Press the ▲ /▼ keys to highlight PC and press the *f*2 (*select*) key. View and change the PC settings, if required.
  - PC On/Off Press the ▲ /▼ keys to highlight PC On/Off and press the *f*2 (*select*) key. To enable communication to a computer, press the ▲ /▼ keys to highlight On and press the *f*2 (*accept*) key. The meter will send the data to a computer in a comma delimited format. To disable communication to a computer, press the ▲ /▼ keys to highlight Off and press the *f*2 (*select*) key.
  - Baud Rate Press the ▲ /▼ keys to highlight Baud Rate and press the *f*2 (select) key. Press the ▲ /▼ keys to highlight a baud rate of 1200, 2400, 4800, 9600, 19200 or 38400 and press the *f*2 (accept) key.
  - Once all of the PC settings are entered, press the *f1 (back)* key.
- Printer Press the ▲ /▼ keys to highlight Printer and press the f2 (select) key. View and change the printer settings, if required.
  - Printer On/Off Press the ▲ /▼ keys to highlight Printer On/Off and press the f2 (select) key. To enable communication to a printer, press the ▲ /▼ keys to

highlight On and press the *f2 (accept)* key. To disable communication to a printer, press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Off and press the *f2 (accept)* key.

- Baud Rate Press the ▲ /▼ keys to highlight Baud Rate and press the *f*2 (select) key. Press the ▲ /▼ keys to highlight a baud rate of 1200, 2400, 4800, 9600, 19200 or 38400 and press the *f*2 (accept) key.
- Once all of the printer settings are entered, press the *f1 (back)* key.
- Data Log Press the ▲ /▼ keys to highlight Data Log and press the f2 (select) key. View and change the data log settings, if required.
  - Data Log On/Off Press the ▲ /▼ keys to highlight Data Log On/Off and press the f2 (select) key. To enable the data log, press the ▲ /▼ keys to highlight On and press the f2 (accept) key. To disable the data log, press the ▲ /▼ keys to highlight Off and press the f2 (accept) key.
  - Ring Buffer Press the ▲ /▼ keys to highlight Ring Buffer and press the *f2 (select)* key. Press the ▲ /▼ keys to highlight Off or On and press the *f2 (accept)* key. If Off is selected, an error message will be displayed when all 1000 data log points are filled and the data log must be emptied to clear the error message. If On is selected, the oldest data log point will be overwritten with the new data log point when all 1000 data log points are filled.
  - Once all of the data log settings are entered, press the *f1* (back) key.
- Once all of the exporting settings are entered, press the f1 (back) key.

### Stirrer Speed

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Stirrer Speed and press the *f*2 (*select*) key. Press the  $\blacktriangle$  /  $\checkmark$  keys to set the default stirrer speed to 1 (slowest speed), 2, 3, 4 or 5 (fastest speed) and press the *f*2 (*accept*) key. The default stirrer speed will be used whenever the stirrer key is pressed in the measurement mode.

### Log View

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Log View and press the *f2 (select)* key. Select Data Log or Calibration Log to view, export or clear the selected log.

Data Log – Press the ▲ /▼ keys to highlight Data Log and press the *f2 (accept)* key. The meter will display a list of the data points. The list shows the sequential number of the data point (D0001), the date that the data point was saved (07/05/2008) and the time that the data point was saved (14:05). To view the measurement information for an individual data point, press the ▲ /▼ keys to highlight the data point and press the *f2 (select)* key. Press the *f2 (print)* key to print the data point or press the *f1 (back)* key to return to the list of data points. To export the data log to a computer or printer or to clear the data log, press the *f3 (options)* key and select Log Export or Log Clear.

- Log Export Press the ▲ /▼ keys to highlight Log Export and press the *f2 (accept)* key. Press the ▲ /▼ keys to highlight Today, Last 10, Range or All as the data log points to be exported and press the *f2 (accept)* key. If Today, Last 10 or All is selected, press the *f2 (yes)* key to export the data. If Range is selected, use the numeric keypad to enter the number of the first data point to be exported, press the *f2 (accept)* key, use the numeric keypad to enter the number of the last data point to be exported, press the *log/print* key and press the *f2 (yes)* key to export the data.
- Log Clear Press the ▲ /▼ keys to highlight Log Clear and press the *f2* (accept) key. Press the *f2* (yes) key to clear the entire data log. Enter the password using the alphanumeric keypad and press the *f2* (accept) key. The default password is 111111111. Press the *f1* (back) key to return to the list of data points.
- Press the *f1 (back)* key to return to the log view settings.
- Calibration Log Press the ▲ /▼ keys to highlight Calibration Log and press the *f2 (accept)* key. Press the ▲ /▼ keys to highlight Channel 1 Cal or Channel 2 Cal and press the *f2 (select)* key. Press the ▲ /▼ keys to highlight pH, ISE, RmV, ORP or Temperature as the calibration type and press the *f2 (select)* key. The meter will display a list of calibrations for the selected channel and calibration type. The list shows the sequential number of the calibration (Cal\_101), the date that the point was saved (07/05/2008) and the time that the point was saved (12:05). To view the calibration data, press the ▲ /▼ keys to highlight the calibration point and press the *f2 (select)* key. Press the *f2 (print)* key to print the calibration point, press the *f3 (graph)* key to view the calibration graph or press the *f1 (back)* key to return to the list of calibration points.
- Press the *f1 (back)* key to return to the log view settings.
- Once all of the log view settings are entered, press the *f1* (back) key.

**Note:** When a data point is exported to a computer or printer, the time stamp will include the hour, minutes and seconds.

### Display

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Display and press the *f2 (select)* key. View and change the Display View, Backlight and Contrast settings, if required.

- Display View Press the ▲ /▼ keys to highlight Display View and press the *f2* (select) key. Press the ▲ /▼ keys to highlight Channel 1, Channel 2 or Channel 1 & 2 and press the *f2* (accept) key.
- Backlight Press the ▲ /▼ keys to highlight Backlight and press the f2 (select) key.
   Press the ▲ /▼ keys to set the backlight to Off, 1 (lowest setting), 2, 3, 4 or 5 (highest setting) and press the f2 (accept) key.

- Contrast Press the ▲ /▼ keys to highlight Contrast and press the *f2 (select)* key. Press the ▲ /▼ keys to set the contrast to 1 (least contrast), 2, 3, 4 or 5 (most contrast) and press the *f2 (accept)* key.
- Once all of the display settings are entered, press the f1 (back) key.

### Calendar

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Calendar and press the *f2 (select)* key. View and change the Time, Date and Date Format settings, if required.

- Time Press the ▲ /▼ keys to highlight Time and press the *f2* (*select*) key. Use the numeric keypad to set the time in 24 hour format and press the *f2* (*accept*) key.
- Date Press the ▲ /▼ keys to highlight Date and press the *f*2 (*select*) key. Use the numeric keypad to set the date (the date format is shown below the date setting) and press the *f*2 (*accept*) key.
- Date Format Press the ▲ /▼ keys to highlight Date Format and press the f2 (select) key. Press the ▲ /▼ keys to highlight MM/DD/YYYY to display the date as month/day/year or DD/MM/YYYY to display the date as day/month/year and press the f2 (accept) key.
- Once all of the date settings are entered, press the *f1 (back)* key.

### Troubleshooting

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Troubleshooting and press the *f2 (select)* key to access the stability test, self test, user reset, factory configuration or general meter information. Refer to the Troubleshooting Guide section for details.

- Stability Test Press the ▲ /▼ keys to highlight Stability Test and press the f2 (accept) key. Press the ▲ /▼ keys to highlight Channel 1 or Channel 2 and press the f2 (accept) key. Make sure that the electrode or probe to be tested is connected to the channel that was selected in the previous step and press the f2 (yes) key. Place the electrode or probe in an appropriate solution with a stable temperature (a pH 7 buffer is recommended for a pH electrode, a calibration standard with a moderate concentration of the ion of interest is recommended for an ISE and an ORP standard is recommended for an ORP electrode) and press the f2 (yes) key. The meter will perform the stability test and show the mV per minute drift and noise on the display. Press the f1 (escape) key to return to the measurement mode.
- Self Test Press the ▲ /▼ keys to highlight Self Test and press the *f2 (accept)* key. Disconnect all of the electrodes and probes from the meter, attach both BNC shorting caps to the two BNC inputs and press the *f2 (yes)* key. Press the *f2 (yes)* key to start the self test. The meter will perform an accuracy test. When the meter displays Accuracy Test Passed, press the *f2 (next)* key. The meter will perform an EPROM test. When the meter displays E2Prom Test Passed, press the *f2 (next)* key. The meter will perform a keypad test. Press each of the blinking keys shown on the display in the order that they are shown.

When the meter displays Keypad Test Passed, press the *f*2 (*next*) key. The meter will display Self Test Passed. Press the *f*1 (*escape*) key to return to the measurement mode.

- User Reset Press the ▲ /▼ keys to highlight User Reset and press the *f2 (accept)* key. If the user reset is performed, all of the current meter settings will be deleted and reset to the default settings. The data log, calibration log and methods will not be deleted with the user reset. Press the *f2 (yes)* key to perform the reset or press the *f3 (no)* key to abort the reset. The meter will enter the EZ Startup menu and then proceed to the measurement mode.
- Factory Config Press the ▲ /▼ keys to highlight Factory Config and press the f2 (accept) key. If the factory configuration is performed, all of the current meter settings will be deleted and reset to the default settings and the data log, calibration log and methods will be deleted. Press the f2 (yes) key to perform the reset or press the f3 (no) key to abort the reset. Use the alphanumeric keypad to enter the PIN number and press the f2 (accept) key to perform the reset to the default PIN number is 11111111. Press the f1 (measure) key, the meter will enter the EZ Startup menu and then proceed to the measurement mode.
- About Meter Press the ▲ /▼ keys to highlight About Meter and press the f2 (accept) key. The meter serial number and software version will be displayed. Press the f1 (back) key to return to the troubleshooting settings and press the f1 (back) key to return to the instrument parameter settings.

### **Beeper Volume**

Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Beeper Volume and press the *f2 (select)* key. Press the  $\land$  /  $\checkmark$  keys to highlight Off, Low or High and press the *f2 (accept)* key.

### Language

Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight Language and press the *f*2 (*select*) key. Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight English, Español, Deutsch, Français or Italiano and press the *f*2 (*accept*) key.

### **Autosampler Setup Menu**

Autosampler Setup		
Autosampler	On	
Ch.1 Method	M101	
Ch.2 Method	M201	
Calibration	On	
Temperature	ATC1	
Tray Size	28	
Rinse Beakers	3	
Rinse Time	5	
<ul> <li>Stirrer Speed</li> </ul>	1	
Samples		
back select		

The autosampler menu contains a list of meter parameters and settings that are used when the meter runs the autosampler and the command to turn the autosampler on or off. For instructions on navigating this menu, refer to Section 8 *Autosampler Operation*.

### **Incremental Techniques Setup Menu**



The incremental techniques menu allows the operator to access the single known addition, single known subtraction, double known addition and double known subtraction meter programs. For instructions on navigating this menu, refer to Chapter 6 *Incremental Techniques*.

### **GLP Setup Menu**

- 1. In the measurement mode, press the setup key to enter the setup menu.
- 2. Press the ▲ / ▼ keys to highlight GLP and press the *f*2 (*select*) key.
- 3. The meter will prompt for the old password. Enter the current password and press the *f2 (accept)* key. The default password is 111111111.
- 4. The meter will prompt for the new password. Enter a new password using the alphanumeric keypad and press the *f2* (accept) key.

### **EZ Startup Setup Menu**

For instructions on navigating this menu, refer to the EZ Startup Menu section.

# 

# **SECTION 4** Calibration and Measurement

The Orion Dual Star meter can perform a one to six point pH calibration, a two to six point ISE (concentration) calibration, a one point ORP calibration and a one point temperature calibration.

Step-by-step instructions and prompts are displayed on the meter to assist operators with the calibration menus. A calibration graph, which includes the electrode slope between points and the average slope, can be viewed in the calibration mode for pH and ISE (concentration) calibrations. The meter can save up to 10 calibrations per channel in the log and calibration data can be exported from the meter to a printer or computer.

# pH Calibration and Measurement

### pH Calibration Hints

- The measurement mode (displayed to the right of the measurement value) determines the type of calibration that the meter will perform. Set the measurement mode to pH for the channel that will be calibrated. Refer to the <u>Setup Menu</u> section for instructions on setting the measurement mode.
- Prepare and condition the pH electrode per the instructions in the electrode user guide.
- Make sure that the pH electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the channel that will be calibrated.
- Always use fresh pH buffers for calibration. Choose buffers that are 1 to 4 pH units apart.
- The meter will automatically recognize the US buffer set (pH 1.68, 4.01, 7.00, 10.01 and 12.46 buffers) or the EURO buffer set (pH 1.68, 4.01, 6.86 and 9.18 buffers). The default meter setting is the US buffer set.

### pH Calibration Procedure (1 to 6 points)

**Note:** In most calibration screens, press the *f1 (escape)* key to return to the measurement mode without saving the calibration.

- 1. In the measurement mode, press the *cal* key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel to be calibrated and press the *f*2 (*accept*) key.
- 3. Rinse the pH electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the buffer.
- 4. When the electrode and buffer are ready, press the f3 (start) key.
- 5. Wait for the pH value on the meter to stop flashing and perform one of the following actions:
  - a. Press the f2 (accept) key to accept the pH value.
  - b. Use the numeric keypad and *decimal* key to enter the value of the buffer at the measured temperature. Press the *f3 (clear)* key to delete any of the digits. Once the buffer value is correct, press the *f2 (accept)* key.
- 6. Press the *f*2 (*next*) key to proceed to the next buffer and repeat steps 3 through 5 or press the *f*3 (*cal done*) key to end the calibration. If a one point calibration is performed, enter the electrode slope using the numeric keypad and *decimal* key and press the *f*2 (*accept*) key.
- 7. Once the *f*3 (*cal done*) key is pressed, the meter will display a calibration summary including the average slope. Perform one of the following actions:



**Important Note:** The *f*2 (*log/print*) key must be pressed to save the calibration. Only pressing the *f*3 (*cal done*) key without pressing the *f*2 (*log/print*) key will not save the calibration.

- a. Press the f2 (log/print) key to save and end the calibration, export the calibration data to the calibration log and export the calibration data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. The meter will automatically proceed to the measurement mode.
- b. Press the f3 (graph) key to view a point-by-point graph of the calibration. After viewing the graph, press the f1 (back) key to return to the calibration summary, press the f3 (log/print) key to save and end the calibration, export the calibration data to the calibration log, export the calibration data to a printer or computer (if applicable) and automatically proceed to the measurement mode or press the f2 (escape) key to return to the measurement mode without saving the calibration.
- c. Press the *f1* (*cal edit*) key to edit the calibration points. Press the ▲ / ▼ keys to highlight the calibration point to be edited and press the *f2* (*select*) key. Press the ▲ / ▼ keys to highlight Remeasure, Edit or Delete and press the *f2* (*select*) key.

- If Remeasure is selected, repeat steps 3 through 5. If a one point calibration is performed, enter the electrode slope using the numeric keypad and decimal key and press the *f2 (accept)* key. Press the ▲ / ▼ keys to highlight another calibration point to be edited and press the *f2 (select)* key or press the *f1 (back)* key to return to the calibration summary.
- ii. If Edit is selected, use the numeric keypad and decimal key to enter a new buffer value and press the *f*2 (accept) key. Press the ▲ / ▼ keys to highlight another calibration point to be edited and press the *f*2 (select) key or press the *f*1 (back) key to return to the calibration summary.
- iii. If Delete is selected, press the *f*2 (*yes*) key to delete the selected calibration point. If deleting the calibration point changes the calibration to a one point calibration, enter the electrode slope using the numeric keypad and decimal key and press the *f*2 (*accept*) key. Press the ▲ /▼ keys to highlight another calibration point to be edited and press the *f*2 (*select*) key or press the *f*1 (*back*) key to return to the calibration summary.
- 8. Proceed to the pH Measurement Procedure section.

### **pH Measurement Procedure**

- 1. Rinse the pH electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample.
- If the stirrer probe is in use, press the *stirrer* key to turn on the stirrer probe. Press the
   ▲ /▼ keys to adjust the speed of the stirrer probe.
- 3. Wait for the measurement to stabilize and record the pH and temperature of the sample when the meter indicates that the measurement is stable. The read type selected in the EZ Startup menu or setup menu will determine how the pH measurements are displayed by the meter. The available read types are Auto-Read, on ready, at time intervals, continuous, value change in measurement and timed reading. To review or change the available read type, refer to the <u>Setup Menu</u> section.
- 4. If the stirrer probe is in use, press the *stirrer* key to turn off the stirrer probe.
- 5. Remove the electrode(s) from the sample, rinse with distilled water, blot dry, place the electrode(s) into the next sample.
- 6. Repeat steps 2 through 5 for all of the samples.
- When all of the samples have been measured, store the electrode(s) according to the instructions in the electrode user guides.
## **ISE** Calibration and Measurement

#### **ISE Calibration Hints**

- The measurement mode (displayed to the right of the measurement value) determines the type of calibration that the meter will perform. Set the measurement mode to ISE for the channel that will be calibrated and set the displayed ISE units as ppm, mg/L, %, M, or unitless. Refer to the <u>Setup Menu</u> section for instructions on setting the measurement mode and displayed units.
- Prepare and condition the ISE electrode per the instructions in the electrode user guide.
- Make sure that the ISE electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the channel that will be calibrated.
- Prepare at least two standards that bracket the expected sample range and differ in concentration by a factor of ten.
- Always use freshly prepared standards for calibration.
- Allow all standards and samples to reach the same temperature for precise measurements.
- 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 rinse the electrode with distilled water between measurements and blot or shake the electrode to remove the water and prevent sample carryover. Do not wipe or rub the electrode sensing element.

#### **ISE Calibration Procedure (2 to 6 points)**

**Note:** In most calibration screens, press the *f1 (escape)* key to return to the measurement mode without saving the calibration.

- 1. In the measurement mode, press the *cal* key.
- Dual channel display only: Press the ▲ / ▼ keys to select the channel to be calibrated and press the *f*2 (accept) key.
- Rinse the ISE electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the standard.
- 4. When the electrode and standard are ready, press the f3 (start) key.
- 5. Wait for the concentration value to stop flashing. Use the numeric keypad and *decimal* key to enter the concentration of the standard at the measured temperature. Press the *f3* (*clear*) key to delete any digits. Once the concentration value is correct, press the *f2* (*accept*) key.

- 6. Press the *f*2 (*next*) key to proceed to the next standard and repeat steps 3 through 5 or press the *f*3 (*cal done*) key to end the calibration.
- 7. Once the *f3 (cal done)* key is pressed, the meter will display a calibration summary including the average slope. Perform one of the following actions:



**Important Note:** The *f*2 (*log/print*) key must be pressed to save the calibration. Only pressing the *f*3 (*cal done*) key without pressing the *f*2 (*log/print*) key will not save the calibration.

- a. Press the f2 (log/print) key to save and end the calibration, export the calibration data to the calibration log and export the calibration data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. The meter will automatically proceed to the measurement mode.
- b. Press the f3 (graph) key to view a point-by-point graph of the calibration. After viewing the graph, press the f1 (back) key to return to the calibration summary, press the f3 (log/print) key to save and end the calibration, export the calibration data to the calibration log, export the calibration data to a printer or computer (if applicable) and automatically proceed to the measurement mode or press the f2 (escape) key to return to the measurement mode without saving the calibration.
- c. Press the *f1 (cal edit)* key to edit the calibration points. Press the ▲ /▼ keys to highlight the calibration point to be edited and press the *f2 (select)* key. Press the ▲ /▼ keys to highlight Remeasure, Edit or Delete and press the *f2 (select)* key.
  - i. If Remeasure is selected, repeat steps 3 through 5. Press the ▲ / ▼ keys to highlight another calibration point to be edited and press the *f*2 (*select*) key or press the *f*1 (*back*) key to return to the calibration summary.
  - If Edit is selected, use the numeric keypad and decimal key to enter a new concentration value and press the *f*2 (*accept*) key. Press the ▲ /▼ keys to highlight another calibration point to be edited and press the *f*2 (*select*) key or press the *f*1 (*back*) key to return to the calibration summary.
  - If Delete is selected, press the *f*2 (*yes*) key to delete the selected calibration point. If deleting the calibration point changes the calibration to a one point calibration, enter the electrode slope using the numeric keypad and decimal key and press the *f*2 (*accept*) key. Press the ▲ /▼ keys to highlight another calibration point to be edited and press the *f*2 (*select*) key or press the *f*1 (*back*) key to return to the calibration summary.
- 8. Proceed to the ISE Measurement Procedure section.

#### **ISE Measurement Procedure**

- 1. Rinse the ISE electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water and place the electrode(s) into the sample.
- If the stirrer probe is in use, press the *stirrer* key to turn on the stirrer probe. Press the
   ▲ /▼ keys to adjust the speed of the stirrer probe.
- 3. Wait for the measurement to stabilize and record the concentration and temperature of the sample when the meter indicates that the measurement is stable. The read type selected in the EZ Startup menu or setup menu will determine how the ISE measurements are displayed by the meter. The available read types are Auto-Read, on ready, at time intervals, continuous, value change in measurement and timed reading. To review or change the available read type, refer to the <u>Setup Menu</u> section.
- 4. If the stirrer probe is in use, press the stirrer key to turn off the stirrer probe.
- 5. Remove the electrode(s) from the sample, rinse with distilled water, blot dry, place the electrode(s) into the next sample.
- 6. Repeat steps 2 through 5 for all of the samples.
- 7. When all of the samples have been measured, store the electrode(s) according to the instructions in the electrode user guides.

# mV Measurement and RmV and ORP Calibration and Measurement

#### **mV Measurement Procedure**

Measure the raw millivolt (mV) values of an ORP, ISE or pH electrode in the mV mode. Oxidation-reduction potential (ORP) measurements are commonly reported as raw millivolts. The mV readings of an ISE electrode in standards can be used to draw a calibration curve on semi-logarithmic graph paper, which can be helpful for low level ISE measurements in the nonlinear range of the electrode. The millivolt readings of a pH electrode can be used to troubleshoot the electrode and verify it is reading the correct mV range of common pH buffers.

- Set the measurement mode to mV in the setup menu. Prepare the electrode according to the instructions in the electrode user guide. Ensure that the electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the channel that will be measured.
- 2. Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample.
- 3. If the stirrer probe is in use, press the stirrer key to turn on the stirrer probe.
- 4. Wait for the measurement to stabilize and record the mV and temperature of the sample when the meter indicates that the measurement is stable. The read type selected in the EZ Startup menu or setup menu will determine how the mV measurements are displayed by the meter. The available read types are Auto-Read, on ready, at time intervals, continuous, value change in measurement and timed reading. To review or change the available read type, refer to the <u>Setup Menu</u> section.
- 5. If the stirrer probe is in use, press the *stirrer* key to turn off the stirrer probe.
- 6. Remove the electrode(s) from the sample.
- 7. Repeat steps 2 through 6 for all of the samples.
- 8. When all of the samples have been measured, store the electrode(s) according to the instructions in the electrode user guides.

#### **RmV Calibration Hints**

- The measurement mode (displayed to the right of the measurement value) determines the type of calibration that the meter will perform. Set the measurement mode to RmV for the channel that will be calibrated. Refer to the <u>Setup Menu</u> section for instructions on setting the measurement mode.
- Prepare and condition the electrode per the instructions in the electrode user guide.
- Ensure that the electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the channel that will be calibrated.
- Always use a freshly prepared standard for calibration. Allow the standard and samples to reach the same temperature for precise measurements.

#### **RmV Calibration Procedure**

**Note:** In most calibration screens, press the *f1 (escape)* key to return to the measurement mode without saving the calibration.

- 1. In the measurement mode, press the *cal* key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel to be calibrated and press the *f*2 (*accept*) key.
- 3. Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the standard.
- 4. When the electrode(s) and standard are ready, press the f3 (start) key.
- 5. Wait for the value to stabilize and the meter to display ready next to the raw mV value. The raw mV value is the actual mV reading of the electrode. To enter a relative mV value, use the numeric keypad, *decimal* key and +/- key to enter the mV value of the standard. Press the *f3 (clear)* key to delete any digits. Once the value is correct, press the *f2 (accept)* key.
- 6. The meter will display the mV offset. Perform one of the following actions:



**Important Note:** The *f*2 (*log/print*) key must be pressed to save the calibration. Only pressing the *f*3 (*cal done*) key without pressing the *f*2 (*log/print*) key will not save the calibration.

- a. Press the *f2 (log/print)* key to save and end the calibration, export the calibration data to the calibration log and export the calibration data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. The meter will automatically proceed to the measurement mode.
- b. Press the *f1* (*escape*) key to return to the measurement mode without saving the calibration.
- 7. Proceed to the RmV Measurement Procedure section.

#### **RmV Measurement Procedure**

- 1. Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry and place the electrode into the sample.
- If the stirrer probe is in use, press the *stirrer* key to turn on the stirrer probe. Press the ▲ /▼ keys to adjust the speed of the stirrer probe.
- 3. Wait for the measurement to stabilize and record the relative mV and temperature of the sample when the meter indicates that the measurement is stable. The read type selected in the EZ Startup menu or setup menu will determine how the relative mV measurements are displayed by the meter. The available read types are Auto-Read, on ready, at time intervals, continuous, value change in measurement and timed reading. To review or change the available read type, refer to the Setup Menu section.
- 4. If the stirrer probe is in use, press the stirrer key to turn off the stirrer probe.
- 5. Remove the electrode(s) from the sample, rinse with distilled water, blot dry, place the electrode(s) into the next sample.
- 6. Repeat steps 2 through 5 for all of the samples.
- 7. When all of the samples have been measured, store the electrode(s) according to the instructions in the electrode user guides.

#### **ORP Calibration Hints**

The millivolt reading of an ORP standard is dependent on the reference system, reference material and filling solution of the ORP electrode and the temperature of the ORP standard. The ORP electrode can be automatically calibrated to read 420 mV relative to the Standard Hydrogen Electrode ( $E_H$ ) at 25 °C when the Thermo Scientific Orion ORP standard (Cat. No. 967901 or Cat. No. 967961), Thermo Scientific Orion ORP electrode (Cat. No. 9678BNWP, 9180BNMD or 9179BNMD) and 4 M KCl filling solution (Cat. No. 900011) or 4 M KCl gel (in Cat. No. 9179BNMD) are used. The raw mV reading of an electrode with the same conditions is 220 mV at 25 °C.

The Orion ORP standard is stable and will not change over time, so it is an ideal standard for any ORP measurement system. The Orion Dual Star meter offers a simple, automatic calibration to the E<sub>H</sub> value in the ORP mode when using Orion ORP electrodes and standard. Visit <u>www.thermoscientific.com/water</u> for ORP application notes that include a table of ORP standard values at different temperatures.

- The measurement mode (displayed to the right of the measurement value) determines the type of calibration that the meter will perform. Set the measurement mode to ORP for the channel that will be calibrated. Refer to the <u>Setup Menu</u> section for instructions on setting the measurement mode.
- Prepare and condition the ORP electrode per the instructions in the electrode user guide.

- Make sure that the ORP electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the channel that will be calibrated.
- The meter will automatically recognize the Orion ORP standard during an ORP calibration. Always use a freshly prepared ORP standard for calibration. Allow the standard and samples to come to the same temperature for precise measurements.

#### **ORP Calibration Procedure**

**Note:** In most calibration screens, press the *f1 (escape)* key to return to the measurement mode without saving the calibration.

- 1. In the measurement mode, press the *cal* key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel to be calibrated and press the *f*2 (*accept*) key.
- Rinse the ORP electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint free tissue and place the electrode(s) into the standard.
- 4. When the electrode and standard are ready, press the f3 (start) key.
- 5. Wait for the value to stabilize and the meter to display ready next to the  $E_H$  value. If the raw mV reading of the electrode is 220 mV ± 60 mV, the meter will automatically calculate and display the  $E_H$  value for the electrode at the measured temperature. Press the *f*2 (accept) key to accept the  $E_H$  value.
- 6. The meter will display the mV offset. Perform one of the following actions:



**Important Note:** The *f*2 (*log/print*) key must be pressed to save the calibration. Only pressing the *f*3 (*cal done*) key without pressing the *f*2 (*log/print*) key will not save the calibration.

- a. Press the *f*2 (*log/print*) key to save and end the calibration, export the calibration data to the calibration log and export the calibration data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. The meter will automatically proceed to the measurement mode.
- b. Press the *f1* (*escape*) key to return to the measurement mode without saving the calibration.
- 7. Proceed to the ORP Measurement Procedure section.

#### **ORP Measurement Procedure**

- 1. Rinse the ORP electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample.
- If the stirrer probe is in use, press the *stirrer* key to turn on the stirrer probe. Press the
   ▲ /▼ keys to adjust the speed of the stirrer probe.
- 3. Wait for the measurement to stabilize and record the ORP and temperature of the sample when the meter indicates that the measurement is stable. The read type selected in the EZ Startup menu or setup menu will determine how the ORP measurements are displayed by the meter. The available read types are Auto-Read, on ready, at time intervals, continuous, value change in measurement and timed reading. To review or change the available read type, refer to the <u>Setup Menu</u> section.
- 4. If the stirrer probe is in use, press the stirrer key to turn off the stirrer probe.
- 5. Remove the electrode(s) from the sample, rinse with distilled water, blot dry, place the electrode(s) into the next sample.
- 6. Repeat steps 2 through 5 for all of the samples.
- 7. When all of the samples have been measured, store the electrode(s) according to the instructions in the electrode user guides.

## Temperature Calibration and Measurement

#### **Temperature Calibration Procedure**

**Note:** The meter ATC temperature display has a relative accuracy of  $\pm 0.1$  °C. ATC probes have varying temperature accuracies, usually  $\pm 0.5$  °C to  $\pm 2$  °C. Use this function only if it is necessary to calibrate the temperature readings. Since the temperature offset calculated during the calibration is applied to all future temperature measurements, recalibrate if a different ATC probe is used.

- 1. In the measurement mode, press the setup key.
- 2. Press the ▲ / ▼ keys to highlight Channel 1 or Channel 2 and press the f2 (select) key.
- 3. Press the ▲ / ▼ keys to highlight Temperature Calibration and press the f2 (select) key.
- Press the ▲ /▼ keys to highlight ATC1, ATC2 or ATC1 & ATC2 to select the ATC probe(s) that will be calibrated and press the *f*2 (*select*) key.
- Place the ATC probe(s) and thermometers into a solution with a known, stable temperature. It is recommended that two NIST traceable thermometers be used to measure and verify the temperature of the solution.
- 6. Wait for the readings of the ATC probe(s) and thermometers to stabilize (about five to ten minutes); use the numeric keypad, *decimal* key and +/- key to enter the temperature value read by the thermometer in the Reference Temperature: \_\_\_\_\_°C field and press the *f2 (accept)* key.
- The meter will display the original temperature read by the ATC probe (i.e. ATC1: 24.2°C) and the temperature offset value that will be applied to all future temperature readings (i.e. ATC1 Offset: -0.8°C).
- 8. Press the f3 (log/print) key to save and end the calibration, export the data to the calibration log, export the data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu and proceed to the measurement mode or press the f1 (escape) key to return to the measurement mode without saving the calibration.

**Note:** The Temperature Calibration menu is for the calibration of ATC probes only. If manual temperature compensation is being used and the display shows MAN as the temperature source icon, refer to the Temperature Input menu for instructions on how to set the manual temperature value. Do not use the Temperature Calibration menu.



## SECTION 5 Methods

The Orion Dual Star meter can save up to ten methods on each channel and the corresponding calibration data for each method, for the fast and easy recollection of different meter setup parameters.

The default meter setting is that no method will be used on either channel. The meter will not display a method number if no method is currently be used on the channel. The current meter settings can be saved by pressing the method key to access the methods list. Existing methods can be selected from the methods list and run on the meter. Methods can be viewed, edited, copied to a new method, deleted, printed or the current meter settings can be saved as a new method.

# Saving the Current Meter Settings as a New Method

- 1. In the measurement mode, press the method key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel that the method will be saved on and press the *f*2 (accept) key.
- 3. The meter will display Current Settings and a list of the saved methods. The list will show the sequential number of the methods (M101) and if any methods were previously saved, the list will show the date that the method was last saved (10/05/2008) and the electrode ID used in the method (pH).
- 4. Press the ▲ / ▼ keys to highlight Current Settings and press the f3 (save) key.
- 5. Press the ▲ / ▼ keys to highlight the method number that the current meter settings will be saved as and press the *f*2 (*accept*) key.
- 6. Enter the current password (default password is 111111111) and press the f2 (accept) key.
- 7. The meter will automatically proceed to the measurement mode.

## Loading an Existing Method

- 1. In the measurement mode, press the method key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel that the method will be saved on and press the *f*2 (accept) key.
- 3. The meter will display Current Settings and a list of the saved methods. The list will show the sequential number of the methods (M101) and if any methods were previously saved, the list will show the date that the method was last saved (10/05/2008) and the electrode ID used in the method (pH).
- 4. Press the ▲ /▼ keys to highlight the method that will be loaded on the channel and press the *f2* (*load*) key.
- 5. Press the *f*2 (*yes*) key to load the method or press the *f*3 (*no*) key to retain the previous settings.
- 6. The meter will automatically proceed to the measurement mode. In the measurement mode, the meter will display the method number that was loaded (i.e. M102).

## Editing or Printing an Existing Method

- 1. In the measurement mode, press the *method* key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel that the method will be saved on and press the *f*2 (accept) key.
- 3. The meter will display Current Settings and a list of the saved methods. The list will show the sequential number of the methods (M101) and if any methods were previously saved, the list will show the date that the method was last saved (10/05/2008) and the electrode ID used in the method (pH).
- 4. Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight the method that will be edited press the f3 (options) key.
- 5. Enter the current password (default password is 111111111) and press the f2 (accept) key.
- 6. Press the  $\blacktriangle$  /  $\triangledown$  keys to highlight Edit and press the *f*2 (accept) key.
- 7. To print the method, press the f3 (print) key.
- To edit the method, press the ▲ /▼ keys to highlight Measure Mode, Electrode ID, Electrode Serial Number, Resolution, Temperature Calibration, Calibration Setup, Temperature Input, Read Type, Alarm or Sample ID and press the *f2 (select)* key.
- Review and change any of the method settings listed above. To select a setting from a list of options, press the ▲ /▼ keys to highlight the desired option and press the *f2 (accept)* key. To enter a numeric value, use the numeric keypad, *decimal* key and +/- key as required and press the *f2 (accept)* key. See the <u>Setup Menu</u> section for details on each method setting.
- 10. Press the *f1* (*back*) key to return to the list of methods. Press the *f1* (*back*) key to return to the measurement mode.

## Coping an Existing Method

- 1. In the measurement mode, press the method key.
- 2. Dual channel display only: Press the ▲ / ▼ keys to select the channel that the method will be saved on and press the *f*2 (accept) key.
- 3. The meter will display Current Settings and a list of the saved methods. The list will show the sequential number of the methods (M101) and if any methods were previously saved, the list will show the date that the method was last saved (10/05/2008) and the electrode ID used in the method (pH).
- 4. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight the method that will be copied press the f3 (options) key.
- 5. Enter the current password (default password is 111111111) and press the f2 (accept) key.
- 6. Press the ▲ / ▼ keys to highlight Copy and press the *f*2 (accept) key.

- 7. Press the ▲ / ▼ keys to highlight the method number that the existing method will be copied to and press the *f*2 (*accept*) key.
- 8. The meter will automatically proceed to the measurement mode.

### Deleting an Existing Method

- 1. In the measurement mode, press the *method* key.
- Dual channel display only: Press the ▲ / ▼ keys to select the channel that the method will be saved on and press the *f*2 (accept) key.
- 3. The meter will display Current Settings and a list of the saved methods. The list will show the sequential number of the methods (M101) and if any methods were previously saved, the list will show the date that the method was last saved (10/05/2008) and the electrode ID used in the method (pH).
- 4. Press the  $\blacktriangle$  /  $\blacktriangledown$  keys to highlight the method that will be deleted press the f3 (options) key.
- 5. Enter the current password (default password is 111111111) and press the *f2 (accept)* key.
- 6. Press the  $\blacktriangle$  /  $\triangledown$  keys to highlight Delete and press the *f*2 (accept) key.
- 7. Press the *f*2 (*yes*) key to delete the method or press the *f*3 (*no*) key to retain the previous settings.
- 8. The meter will automatically proceed to the measurement mode.



# **SECTION 6** Incremental Techniques

The Orion Dual Star meter has preset methods for single known addition, single known subtraction, double known addition and double known subtraction.

- When performing incremental techniques, the sample concentration must be known within an order of magnitude so the increment may be correctly chosen and the sample and standard must be volumetrically determined.
- It is very important that any complexing agent is present in excess (at least 50 to 100 times) or not at all. The ratio of free to complexed ions must remain constant over the addition. An indication that a complexing agent is present, but not in great enough quantity to maintain a constant free ion to complexed ion ratio, is an abnormally high or low slope. If this is a suspected problem, add an excess amount of complexing or decomplexing agent to the original solution in the beaker.
- Electrode interferences should be at a minimum, since the effect of the interference might change as the concentration of the ion of interest changes. This situation may also result in high or low electrode slopes. Consult the electrode user guide for information on reducing or eliminating some electrode interferences.
- Known addition is a useful method for measuring dilute samples, occasional samples or samples that contain an excess of complexing agent. It can also be used to verify the results of a direct measurement.
- Known subtraction is useful when measuring ions for which stable standards do not exist. It can also be used as a quick version of a titration. When performing a known subtraction, it is necessary to know the stoichiometric ratio between the standard and sample.
- Double incremental techniques calculate the electrode slope and sample concentration simultaneously. The electrode slope is determined directly in the sample, which results in greater accuracy for samples with complex matrices and greater analysis speed.

# Electrode ID and Slope for Incremental Techniques

The meter uses the electrode ID to determine the valence of the ion of interest and the theoretical slope of the electrode. The following table shows the electrode ID options and their corresponding valence and theoretical slope values.

Electrode ID	Valence Number	Theoretical Slope
рН	+1	+59.16
Fluoride (F <sup>-</sup> )	-1	-59.16
Ammonia (NH <sub>3</sub> )	-1	-59.16
Ammonium (NH₄+)	+1	+59.16
Nitrate (NO <sub>3</sub> -)	-1	-59.16
Nitrite (NO2 <sup>-</sup> )	-1	-59.16
Sulfide (S <sup>-2</sup> )	-2	-29.58
Chloride (Cl <sup>-</sup> )	-1	-59.16
Chlorine (Cl <sub>2</sub> )	+2	+29.58
Bromide (Br)	-1	-59.16
lodide (l <sup>-</sup> )	-1	-59.16
Cyanide (CN <sup>-</sup> )	-1	-59.16
Sodium (Na⁺)	+1	+59.16
Potassium (K⁺)	+1	+59.16

Electrode ID	Valence Number	Theoretical Slope
Calcium (Ca <sup>+2</sup> )	+2	+29.58
Silver (Ag⁺)	+1	+59.16
Copper (Cu <sup>+2</sup> )	+2	+29.58
Lead (Pb <sup>+2</sup> )	+2	+29.58
Cadmium (Cd <sup>+2</sup> )	+2	+29.58
Perchlorate (ClO4 <sup>-</sup> )	-1	-59.16
Fluoroborate (BF4 <sup></sup> )	-1	-59.16
Thiocyanate (SCN <sup>-</sup> )	-1	-59.16
Nitrogen Oxide (NO <sub>X</sub> )	+1	+59.16
Carbon Dioxide (CO <sub>2</sub> )	+1	+59.16
Monovalent Cation (X⁺)	+1	+59.16
Divalent Cation (X <sup>+2</sup> )	+2	+29.58
Monovalent Anion (X <sup>-</sup> )	-1	-59.16
Divalent Anion (X <sup>-2</sup> )	-2	-29.58

## Single Known Addition

A standard of known concentration is added to a sample of unknown concentration. The total concentration should approximately double upon the addition of the standard, therefore sample concentration should be known within an order of magnitude. Refer to the following table to choose the correct standard concentration and volume of addition.

Amount of standard to be added per 100 mL of sample	Standard concentration compared to expected sample concentration
1 mL	100 times more concentrated
5 mL	20 times more concentrated
10 mL	10 times more concentrated

- Refer to the appropriate electrode user guide for instructions for preparing the electrode, required solutions (ionic strength adjusters, standards, etc) and any special requirements. Determine the electrode slope as directed in the electrode user guide.
- Ensure that the electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the selected channel.

#### **Single Known Addition Procedure**

- 1. Prepare the sample as directed in the electrode user guide.
- 2. In the measurement mode, press the *setup* key. Press the ▲ /▼ keys to highlight Incremental Techniques and press the *f*2 (*select*) key.
- 3. Dual channel display only: Press the ▲ / ▼ keys to select the channel to perform the single known addition on and press the *f*2 (*accept*) key.
- 4. Press the ▲ /▼ keys to highlight Single Known Addition and press the *f*2 (*select*) key.
- 5. Use the numeric keypad, *decimal* key and +/- key to enter the electrode slope and press the *f2 (accept)* key. Determine the electrode slope as directed in the electrode user guide or use the theoretical slope from the <u>Electrode ID and Slope for Incremental Techniques</u> section.
- 6. Use the numeric keypad and *decimal* key to enter the sample volume, in mL, and press the *f2 (accept)* key.
- 7. Use the numeric keypad and *decimal* key to enter the total volume, in mL, and press the *f2 (accept)* key. The total volume includes the sample volume, ISA volume and diluting agent volume, but not the standard volume.
- 8. Use the numeric keypad and *decimal* key to enter the concentration of the standard and press the *f*2 (*accept*) key.

- Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample. When the electrode and sample are ready, press the *f2 (next)* key.
- 10. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 11. Use the numeric keypad and decimal key to enter the volume of the standard, in mL, to be added to the sample and press the *f2* (*accept*) key.
- 12. Add the standard to the sample and press the f2 (yes) key.
- 13. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 14. The meter will display the result of the single known addition. Press the *f3 (log/print)* key to save and end the single known addition, export the data to the data log and export the data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.

**Note:** Press the *f1 (escape)* key to return to the measurement mode without saving the single known addition.

- 15. Press the f2 (next) key.
- 16. To perform a single known addition using the values that were entered for the previous single known addition, press the *f2 (yes)* key. To perform a single known addition and enter new values, press the *f3 (no)* key. To end the incremental technique and return to the measurement mode, press the *f1 (escape)* key.

## Double Known Addition

- For most double known addition analysis, it is recommended that:
  - The sample volume should be 100 mL.
  - The concentration of the standard added should be 100 times the expected sample concentration.
  - The volume of the first addition should be 1 mL and the volume of the second addition should be 10 mL.
- Refer to the appropriate electrode user guide for instructions for preparing the electrode, required solutions (ionic strength adjusters, standards, etc) and any special requirements.
- Ensure that the electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the selected channel.

#### **Double Known Addition Procedure**

- 1. Prepare the sample as directed in the electrode user guide.
- 2. In the measurement mode, press the *setup* key. Press the ▲ /▼ keys to highlight Incremental Techniques and press the *f*2 (*select*) key.
- 3. Dual channel display only: Press the ▲ / ▼ keys to select the channel to perform the double known addition on and press the *f*2 (*accept*) key.
- 4. Press the ▲ /▼ keys to highlight Double Known Addition and press the *f*2 (select) key.
- 5. Use the numeric keypad and *decimal* key to enter the sample volume, in mL, and press the *f2 (accept)* key.
- 6. Use the numeric keypad and *decimal* key to enter the total volume, in mL, and press the *f2 (accept)* key. The total volume includes the sample volume, ISA volume and diluting agent volume, but not the standard volume.
- 7. Use the numeric keypad and *decimal* key to enter the concentration of the standard and press the *f2* (*accept*) key.
- Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample. When the electrode and sample are ready, press the *f*2 (*next*) key.
- 9. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 10. Use the numeric keypad and *decimal* key to enter the volume of the first increment of standard, in mL, to be added to the sample and press the *f2 (accept)* key.
- 11. Add the first increment of standard to the sample and press the f2 (yes) key.
- 12. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.

- 13. Use the numeric keypad and *decimal* key to enter the volume of the second increment of standard, in mL, to be added to the sample and press the *f*2 (*accept*) key.
- 14. Add the second increment of standard to the sample and press the *f*2 (yes) key.
- 15. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 16. The meter will display the result of the double known addition. Press the *f3 (log/print)* key to save and end the double known addition, export the data to the data log and export the data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.

**Note:** Press the *f1 (escape)* key to return to the measurement mode without saving the double known addition.

- 17. Press the *f*2 (*next*) key.
- 18. To perform a double known addition using the values that were entered for the previous double known addition, press the *f*2 (*yes*) key. To perform a double known addition and enter new values, press the *f*3 (*no*) key. To end the incremental technique and return to the measurement mode, press the *f*1 (*escape*) key.

## Single Known Subtraction

In known subtraction, each addition of standard subtracts an increment of the species to be measured. Enter the correct reaction ratio of the species to be measured with the standard to be added. For example, a sulfide (S<sup>-2</sup>) sample is to be subtracted by the addition of silver (Ag<sup>+</sup>). Two silver ions are required to subtract each sulfide ion and the reaction ratio is 0.5.

Sample concentration should be known within an order of magnitude and the total sample concentration should be approximately halved upon addition of the standard. Refer to the following table to choose the correct standard concentration and volume of addition.

Amount of standard to be added per 100 mL of sample	Standard concentration compared to expected sample concentration
1 mL	50 times more concentrated
5 mL	10 times more concentrated
10 mL	5 times more concentrated

- Refer to the appropriate electrode user guide for instructions for preparing the electrode, required solutions (ionic strength adjusters, standards, etc) and any special requirements. Determine the electrode slope as directed in the electrode user guide.
- Ensure that the electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the selected channel.

#### **Single Known Subtraction Procedure**

- 1. Prepare the sample as directed in the electrode user guide.
- 2. In the measurement mode, press the *setup* key. Press the ▲ / ▼ keys to highlight Incremental Techniques and press the *f*2 (*select*) key.
- 3. Dual channel display only: Press the ▲ / ▼ keys to select the channel to perform the single known subtraction on and press the *f*2 (accept) key.
- 4. Press the ▲ /▼ keys to highlight Single Known Subtraction and press the *f*2 (select) key.
- Use the numeric keypad, *decimal* key and +/- key to enter the electrode slope and press the f2 (accept) key. Determine the electrode slope as directed in the electrode user guide or use the theoretical slope from the table in the <u>Electrode ID and Slope for Incremental</u> <u>Techniques</u> section.
- 6. Use the numeric keypad and *decimal* key to enter the sample volume, in mL, and press the *f2 (accept)* key.
- 7. Use the numeric keypad and *decimal* key to enter the total volume, in mL, and press the *f2 (accept)* key. The total volume includes the sample volume, ISA volume and diluting agent volume, but not the standard volume.

- 8. Use the numeric keypad and *decimal* key to enter the concentration of the standard and press the *f2* (*accept*) key.
- 9. Use the numeric keypad and *decimal* key to enter the reaction ratio of sample to standard and press the *f2 (accept)* key. For example, if a sulfide (S<sup>-2</sup>) sample is to be subtracted by the addition of silver (Ag<sup>+</sup>), two silver ions are required to subtract each sulfide ion and the reaction ratio is 0.5.
- 10. Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample. When the electrode and sample are ready, press the *f*2 (*next*) key.
- 11. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 12. Use the numeric keypad and decimal key to enter the volume of the standard, in mL, to be added to the sample and press the *f2 (accept)* key.
- 13. Add the standard to the sample and press the f2 (yes) key.
- 14. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 15. The meter will display the result of the single known subtraction. Press the *f*3 (*log/print*) key to save and end the single known subtraction, export the data to the data log and export the data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.

**Note:** Press the f1 (escape) key to return to the measurement mode without saving the single known subtraction.

- 16. Press the f2 (next) key.
- 17. To perform a single known subtraction using the values that were entered for the previous single known subtraction, press the *f*2 (*yes*) key. To perform a single known subtraction and enter new values, press the *f*3 (*no*) key. To end the incremental technique and return to the measurement mode, press the *f*1 (*escape*) key.

## **Double Known Subtraction**

- Ensure that the subtraction ability of the chosen standard concentration and volume added does not exceed the concentration of sample.
  - o For most double known subtraction analysis, it is recommended that:
  - The sample volume should be 100 mL.
  - The concentration of the standard added should be 50 times the expected sample concentration.
- The volume of the first addition of standard should be 1 mL and the volume of the second addition of standard should be 10 mL.
- Refer to the appropriate electrode user guide for instructions for preparing the electrode, required solutions (ionic strength adjusters, standards, etc) and any special requirements.
- Ensure that the electrode, ATC probe and reference electrode (if applicable) are connected to the meter inputs that correspond with the selected channel.

#### **Double Known Subtraction Procedure**

- 1. Prepare the sample as directed in the electrode user guide.
- 2. In the measurement mode, press the *setup* key. Press the ▲ / ▼ keys to highlight Incremental Techniques and press the *f2* (*select*) key.
- 3. Dual channel display only: Press the ▲ / ▼ keys to select the channel to perform the double known subtraction on and press the *f*2 (*accept*) key.
- 4. Press the ▲ /▼ keys to highlight Double Known Subtraction and press the f2 (select) key.
- 5. Use the numeric keypad and *decimal* key to enter the sample volume, in mL, and press the *f2 (accept)* key.
- 6. Use the numeric keypad and *decimal* key to enter the total volume, in mL, and press the *f2 (accept)* key. The total volume includes the sample volume, ISA volume and diluting agent volume, but not the standard volume.
- 7. Use the numeric keypad and *decimal* key to enter the concentration of the standard and press the *f*2 (*accept*) key.
- 8. Use the numeric keypad and *decimal* key to enter the reaction ratio of sample to standard and press the *f2* (*accept*) key. For example, if a sulfide (S<sup>-2</sup>) sample is to be subtracted by the addition of silver (Ag<sup>+</sup>), two silver ions are required to subtract each sulfide ion and the reaction ratio is 0.5.
- Rinse the electrode (ATC probe, stirrer probe and reference electrode, if applicable) with distilled water, blot dry with a lint-free tissue and place the electrode(s) into the sample. When the electrode and sample are ready, press the *f2 (next)* key.

- 10. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 11. Use the numeric keypad and *decimal* key to enter the volume of the first increment of standard, in mL, to be added to the sample and press the *f2* (*accept*) key.
- 12. Add the first increment of standard to the sample and press the *f*2 (yes) key.
- 13. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 14. Use the numeric keypad and *decimal* key to enter the volume of the second increment of standard, in mL, to be added to the sample and press the *f*2 (*accept*) key.
- 15. Add the second increment of standard to the sample and press the *f*2 (yes) key.
- 16. Wait for the reading to stabilize. When the reading is stable, press the f2 (yes) key.
- 17. The meter will display the result of the double known subtraction. Press the *f*3 (*log/print*) key to save and end the double known subtraction, export the data to the data log and export the data to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu.

**Note:** Press the *f1 (escape)* key to return to the measurement mode without saving the double known subtraction.

- 18. Press the f2 (next) key.
- 19. To perform a double known subtraction using the values that were entered for the previous double known subtraction, press the *f2 (yes)* key. To perform a double known subtraction and enter new values, press the *f3 (no)* key. To end the incremental technique and return to the measurement mode, press the *f1 (escape)* key.



# **SECTION 7** Data Storage and Retrieval

## Data Storage Settings

The data storage settings include the export trigger, computer, printer and data log settings. The export trigger uses the read type of the selected channel or channels to determine when the meter sends the measurements to the data log, if the data log is enabled in the setup menu, and to a printer or computer, if a printer or computer is connected to the meter and enabled in the setup menu. The computer setting enables communication from the meter to a computer. The printer setting enables communication from the meter to a printer. The data log setting enables or disables the data log.

- 1. In the measurement mode, press the setup key.
- 2. Press the ▲ / ▼ keys to highlight Instrument Parameters and press the f2 (select) key.
- Press the ▲ / ▼ keys to highlight Export Data (PC/Printer/Log) and press the f2 (select) key.
- 4. Press the ▲ /▼ keys to highlight Export Trigger and press the *f2 (select)* key. Press the ▲ /▼ keys to highlight Channel 1, Channel 2 or Channel 1 & 2 as the measurement export trigger and press the *f2 (accept)* key.
- 5. The Orion Dual Star meter can be connected to a computer or a printer, but not both at the same time. If communication to a printer is enabled in the setup menu, communication to a computer will automatically be disabled. If communication to a computer is enabled in the in the setup menu, communication to a printer will automatically be disabled.
  - a. To enable communication to a computer, press the ▲ / ▼ keys to highlight PC and press the *f*2 (*select*) key. Press the ▲ / ▼ keys to highlight PC On/Off and press the

*f2 (select)* key. Press the ▲ / ▼ keys to highlight On and press the *f2 (accept)* key. Press the ▲ / ▼ keys to highlight Baud Rate and press the *f2 (select)* key. Press the ▲ / ▼ keys to highlight 1200, 2400, 4800, 9600, 19200 or 38400 as the baud rate and press the *f2 (accept)* key. Once computer settings are entered, press the *f1 (back)* key.

- b. To enable communication to a printer, press the ▲ /▼ keys to highlight Printer and press the f2 (select) key. Press the ▲ /▼ keys to highlight Printer On/Off and press the f2 (select) key. Press the ▲ /▼ keys to highlight On and press the f2 (accept) key. Press the ▲ /▼ keys to highlight Baud Rate and press the f2 (select) key. Press the ▲ /▼ keys to highlight 1200, 2400, 4800, 9600, 19200 or 38400 as the baud rate and press the f2 (accept) key. Once the printer settings are entered, press the f1 (back) key.
- c. To disable communication to a computer, press the ▲ / ▼ keys to highlight PC and press the *f2* (*select*) key. Press the ▲ / ▼ keys to highlight PC On/Off and press the *f2* (*select*) key. Press the ▲ / ▼ keys to highlight Off and press the *f2* (*accept*) key. Press the *f1* (*back*) key.
- d. To disable communication to a printer, press the ▲ /▼ keys to highlight Printer and press the f2 (select) key. Press the ▲ /▼ keys to highlight Printer On/Off and press the f2 (select) key. Press the ▲ /▼ keys to highlight Off and press the f2 (accept) key. Press the f1 (back) key.
- 6. Press the ▲ /▼ keys to highlight Data Log and press the f2 (select) key. Press the ▲ /▼ keys to highlight Data Log On/Off and press the f2 (select) key. Press the ▲ /▼ keys to highlight On to enable the data log and press the f2 (accept) key or press the ▲ /▼ keys to highlight Off to disable the data log and press the f2 (accept) key. Press the ▲ /▼ keys to highlight Ring Buffer and press the f2 (select) key. Press the ▲ /▼ keys to highlight Ring Buffer and press the f2 (select) key. Press the ▲ /▼ keys to highlight Ring Buffer and press the f2 (select) key. Press the ▲ /▼ keys to highlight Off or On and press the f2 (accept) key. If Off is selected, an error message will be displayed when all 1000 data log points are filled and the data log must be emptied to clear the error message. If On is selected, the oldest data log point will be overwritten with the new data log point when all 1000 data log points are filled. Once all of the data log settings are entered, press the f1 (back) key.
- 7. Once all of the exporting settings are entered, press the *f1 (back)* key until the meter returns to the measurement mode.

## Viewing the Calibration Log and Data Log

#### Viewing the Calibration Log

The Orion Dual Star meter saves up to ten of the most recent calibrations on each channel. Each of the ten calibrations per channel is saved according to the calibration type (pH, ISE, RmV or ORP), so the meter will use the last calibration performed for the calibration type (as long as it is one of the last ten most recent calibrations) and pH, ISE, RmV and ORP electrodes can be interchanged without recalibrating every time the electrodes are switched.

For example: 1. The channel one measurement mode is set to pH, a pH electrode is connected to the channel one input and a pH calibration is performed. 2. The channel one measurement mode is set to ORP, an ORP electrode is connected to the channel one input, an ORP calibration is performed. 3. The channel one measurement mode is set back to pH and the pH electrode is reconnected to the channel one input. In this scenario, the meter will use the pH calibration performed in step 1 for the pH measurements displayed after step 3.

- 1. In the measurement mode, press the log view key.
- 2. Press the ▲ / ▼ keys to highlight Calibration Log and press the *f*2 (accept) key.
- 3. Press the ▲ / ▼ keys to highlight Channel 1 Cal or Channel 2 Cal and press the *f*2 (*select*) *key*.
- 4. Press the ▲ /▼ keys to highlight pH, ISE, RmV, ORP or Temperature as the calibration type and press the *f*2 (*select*) key.
- 5. The meter will display a list of calibrations for the selected channel and calibration type. The list shows the sequential number of the calibration (Cal\_101), the date that the point was saved (07/05/2008) and the time that the point was saved (12:05).
- 6. To view the calibration data, press the ▲ /▼ keys to highlight the calibration point and press the *f*2 (*select*) key. Press the *f*2 (*print*) key to print the calibration point, press the *f*3 (*graph*) key to view the calibration graph or press the *f*1 (*back*) key to return to the list of calibration points.
- 7. Press the f1 (back) key twice to return to the measurement mode.

#### **Viewing the Data Log**

The Orion Dual Star meter saves up to 1000 points of data and assigns a log number to each data point. The read type of the displayed channel or channels determines when measurements are exported to the data log. When the meter is in the dual channel display mode, the measurements from both channels will be saved as two separate data points. When a data point is exported to a computer or printer, the time stamp will include the hour, minutes and seconds.

- 1. In the measurement mode, press the *log view* key.
- Press the ▲ /▼ keys to highlight Data Log and press the *f2 (accept)* key. The meter will display a list of the data points. The list shows the sequential number of the data point (D0001), the date that the data point was saved (07/05/2008) and the time that the data point was saved (14:05).
- 3. To view the measurement information for an individual data point, press the ▲ /▼ keys to highlight the data point and press the *f2* (*select*) key. Press the *f2* (*print*) key to print the data point or press the *f1* (*back*) key to return to the list of data points.
- 4. To export the data log to a computer or printer, press the f3 (options) key, press the ▲ /▼ keys to highlight Log Export and press the f2 (accept) key. Press the ▲ /▼ keys to highlight Today, Last 10, Range or All as the data log points to be exported and press the f2 (accept) key.
  - a. If Today, Last 10 or All is selected, press the *f*2 (*yes*) key to export the data. After exporting the data, the meter will automatically return to the measurement mode.
  - b. If Range is selected, use the numeric keypad to enter the number of the first data point to be exported, press the *f2 (accept)* key, use the numeric keypad to enter the number of the last data point to be exported, press the log/print key and press the *f2 (yes)* key to export the data. After exporting the data, the meter will automatically return to the measurement mode.
  - c. To clear the data log, press the f3 (options) key, press the ▲ / ▼ keys to highlight Log Clear and press the f2 (accept) key. Press the f2 (yes) key to clear the entire data log. Enter the password using the alphanumeric keypad and press the f2 (accept) key. The default password is 11111111. Press the f1 (back) key to return to the list of data points.
- 5. Press the f1 (back) key twice to return to the measurement mode.

## Interfacing a Printer

The Orion Star series printer, Cat. No. 1010006, is recommended for use with the Orion Dual Star meter. The Orion Star series printer has a 9600 baud rate and it is packaged with the RS232 printer cable, Cat. No. 250302-001, which is required to interface the Orion Dual Star meter to a printer. The Orion Star printer is also designed to print the calibration graphs and will show the data in the graph clearly with the best picture and detail. Other printers may print the calibration graphs as a bunch of information that looks like random text and numbers; however, the raw calibration data should print fine.

The baud rate of the Orion Dual Star meter can be set to 1200, 2400, 4800, 9600, 19200 or 38400 for communication to different printers. The fixed meter settings are number of bits = 8, stop bits = 1 and parity = none. The flow control for the printer should be set to XON/XOFF.

The meter will send measurements and calibration data to the printer if the printer option is turned on in the setup menu. Alternatively, measurements and calibration data can be sent to the data log and calibration log. From the data log and calibration log, the operator can chose to print a single data log point, a range of data log points or the entire data log.

## Interfacing a Computer

The Orion Dual Star meter can send measurements and calibration data to a computer in a comma delimited format that is easy to parse in computer programs like Microsoft<sup>®</sup> Excel<sup>®</sup>. Connect the USB computer cable to the USB inputs on the meter and computer and install the Dual Star USB driver, downloadable on our website at <u>www.thermoscientific.com/OrionMeters</u>. Alternatively, connect the RS232 computer cable, Catalog Number 1010053, to the RS232 inputs on the meter and computer.

The baud rate of the Orion Dual Star meter can be set to 1200, 2400, 4800, 9600, 19200 or 38400. The fixed meter settings are number of bits = 8, stop bits = 1 and parity = none. The flow control for the computer should be set to XON/XOFF.

The meter will send measurements and calibration data to the computer if the computer option is turned on in the setup menu. Alternatively, measurements and calibration data can be sent to the data log and calibration log. From the data log and calibration log, the operator can chose to send a single data log point, a range of data log points or the entire data log to the computer.

## Meter Software Upgrade

The Orion Dual Star meter can be updated with the latest software, so the meter programs are always up-to-date.

- 1. Disconnect the power adapter from the meter.
- 2. Connect the USB cable that is packaged with the Orion Dual Star meter to the USB input on a computer.
- 3. Connect the other end of the USB cable to the USB input on the Orion Dual Star meter.
- 4. Press and hold down the power key and the f1 key on the meter keypad.
- 5. Connect the power adapter to the meter.
- 6. Release the power key and the *f*1 key when the meter powers up and beeps twice.
- 7. A window will open on the computer for the USB drive that the meter is connected to. Delete the software file in this window.
- 8. Go to our website at <u>www.thermoscientific.com/OrionMeters</u> and select the link for upgrading the Orion Dual Star meter software. Download the new software file, unzip/extract the file and then copy and paste the new software file into the window that used to contain the old software file.
- 9. Wait for the software file to paste completely into the window.
- 10. Disconnect the power adapter from the meter and then reconnect the power adapter to the meter. Wait for the new software to load.

**Note:** After a new revision of meter software is loaded to the meter, all of the meter settings will be reset to the factory default settings and the data log, calibration log and methods will be cleared.



# **SECTION 8** Autosampler Operation

## Interfacing the Meter and Autosampler



- 1. Select a location for the system and unpack the autosampler and Orion Dual Star meter components. Prepare the autosampler according to the autosampler user guide.
- Connect the meter to the autosampler using the RS232 interface cable. Connect the 2.5 mm stereo jack end of the cable to the RS232 input on the meter. Connect the 9 pin DIN end of the cable to the COM1 RS232 port on the autosampler.
- 3. If a printer will be used, use the autosampler to printer interface cable to connect the COM2 port on the autosampler to the 25 pin RS232 port on the printer.

**Note:** The printer must have a 9600 baud rate to be connected to the autosampler. The Orion Star series printer, Catalog Number 1010006, is recommended for use with the Orion Dual Star meter and autosampler.

4. Prepare and install the electrodes. Prepare the electrodes according to the electrode user guides. Connect the electrodes to the meter using the extension cables and cable management system provided with the autosampler. Insert the electrodes into the electrode holder on the autosampler. If a stirrer probe will be used, connect the stirrer probe to the stirrer jack on the back of the meter and insert the stirrer probe into the electrode holder on the autosampler.

**Note:** The minimum cable length required to connect electrodes to the autosampler is 6.25 feet (1.9 meters). Extension cables or electrodes with an appropriate cable length are needed.

5. Connect the meter, autosampler and printer (if applicable) to a power supply and turn on the power. The default Orion Dual Star meter baud rate is 9600 and the default autosampler baud rate for the COM 1 port is 1200, so the meter baud rate needs to be set to 1200 or the autosampler dip switches need to be adjusted to 9600 for the COM 1 port.

Autosampler Setup		
<ul> <li>Autosampler</li> </ul>	On	
Ch.1 Method	M101	
Ch.2 Method	M201	
<ul> <li>Calibration</li> </ul>	On	
Temperature	ATC1	
Tray Size	28	
Rinse Beakers	3	
Rinse Time	5	
<ul> <li>Stirrer Speed</li> </ul>	1	
Samples		

## Autosampler Setup Menu

The autosampler setup menu can be accessed from the setup menu. The autosampler menu contains ten autosampler parameters: Autosampler, Ch.1 Method, Ch.2 Method, Calibration, Temperature, Tray Size, Rinse Beakers, Rinse Time, Stirrer Speed and Samples.

In the measurement mode, press the setup key. Press the  $\blacktriangle$  /  $\checkmark$  keys to highlight Autosampler and press the *f2* (*select*) key.

- Autosampler Press the ▲ /▼ keys to highlight Autosampler and press the f2 (select) key. Press the ▲ /▼ keys to highlight On or Off to turn communication between the meter and the autosampler on or off and press the f2 (accept) key.
- Ch.1 Method Press the ▲ /▼ keys to highlight Ch.1 Method and press the *f*2 (select) key. To set the method that the autosampler will use for channel 1, press the ▲ /▼ keys to highlight the method number and press the *f*2 (accept) key.

**Note:** A method can be selected when the autosampler setup parameters are entered, but any method editing must be done from the methods list. Calibration values from the last calibration performed in the method are stored to the method, so calibrations can be performed on the autosampler.

Ch.2 Method – Press the ▲ /▼ keys to highlight Ch.2 Method and press the *f*2 (*select*) key. To set the method that the autosampler will use for channel 2, press the ▲ /▼ keys to highlight the method number and press the *f*2 (*accept*) key.

**Note:** A method can be selected when the autosampler setup parameters are entered, but any method editing must be done from the method list. Calibration values from the last calibration performed in the method are stored to the method, so calibrations can be performed on the autosampler.

Calibration – Press the ▲ /▼ keys to highlight Calibration and press the *f2 (select)* key.
 Press the ▲ /▼ keys to highlight On or Off to turn the calibration function for the autosampler on or off and press the *f2 (accept)* key.

- Temperature Press the ▲ /▼ keys to highlight Temperature and press the f2 (select) key. Press the ▲ /▼ keys to highlight ATC1 or ATC2 as the temperature source of the ATC probe that will be used on the autosampler and press the f2 (accept) key.
- Tray Size Press the ▲ /▼ keys to highlight Tray Size and press the *f2 (select)* key.
   Press the ▲ /▼ keys to highlight 28 or 48 as the size of the tray that will be used on the autosampler and press the *f2 (accept)* key.
- Rinse Beakers Press the ▲ /▼ keys to highlight Rinse Beakers and press the *f2 (select)* key. Press the ▲ /▼ keys to highlight 0, 1, 2 or 3 as the number of the rinse beakers that will be used on the autosampler and press the *f2 (accept)* key.
- Rinse Time Press the ▲ /▼ keys to highlight Rinse Time and press the *f2 (select)* key. Use the numeric keypad to enter the time (in seconds) that the electrodes will be rinsed in each rinse beaker and press the *f2 (accept)* key. The rinse time can be set from 1 to 59 seconds.
- Stirrer Speed Press the ▲ /▼ keys to highlight Stirrer Speed and press the *f2* (select) key. Press the ▲ /▼ keys to highlight Off, 1, 2, 3, 4 or 5 as the stirrer speed of the stirrer probe and press the *f2* (accept) key.
- Samples Press the ▲ / ▼ keys to highlight Samples and press the *f*2 (*select*) key.
  - To enter the number of samples that will be run on the autosampler, press the ▲ /▼
     keys to highlight Number of Samples and press the *f2 (accept)* key. Use the numeric keypad to enter the number of samples and press the *f2 (accept)* key.
  - To set the sample ID, press the ▲ /▼ keys to highlight Sample ID and press the f2 (accept) key. Press the ▲ /▼ keys to highlight Off, Manual or Auto Incremental and press the f2 (accept) key. If Auto Incremental is selected, use the alphanumeric keypad to enter the initial sample ID and then press the f2 (accept) key. Refer to the meter pullout guide for the location of the letter keys on the numeric keypad. The meter will save the entered sample ID and add a sequential number after it.
  - Once all of the sample settings are entered, press the f1 (back) key.
- Once all of the autosampler settings are entered, press the f1 (back) key.

## Running the Autosampler

- 1. Set all of the autosampler parameters in the autosampler setup menu. Refer to the Autosampler Setup Menu section for detailed instructions.
- 2. Set the meter display so only the channel or channels that will be run on the autosampler are displayed. Press the channel key to select a dual channel display for running both channels on the autosampler, a single channel display for running channel 1 on the autosampler or a single channel display for running channel 2 on the autosampler.

**Note:** The meter will log the data points for both channels during the autosampler run if the dual channel display is selected. The meter will log the data points for only one channel if the single channel display is selected.

- Ensure that all of the electrodes are connected to the meter inputs that correspond with the channel(s) that will be run on the autosampler. If the stirrer probe, Cat. No. 096019, will be used with the autosampler, connect the stirrer probe to the stirrer input on the back of the autosampler.
- 4. Set the method(s) for the channel(s) that will be run on the autosampler. Methods must be assigned to the channel(s) that will be run on the autosampler before the autosampler can be run. Make sure to assign the measurement mode and electrode ID for the channel(s) that will be run on the autosampler.

The operator may select Current Settings as the method and the meter will use the current meter settings to run on the autosampler. If Current Settings is selected as the method, calibrations cannot be run on the autosampler and the samples will be measured using the last saved calibration.

**Note:** A method can be selected when the autosampler setup parameters are entered, but any method editing must be done from the method list. Calibration values from the last calibration performed in the method are stored to the method, so calibrations can be performed on the autosampler.

- 5. To begin an autosampler run, press the *auto* key, verify that all of the autosampler setup parameters are correct and press the *f3 (start)* key.
  - a. During the autosampler run, press the *f1 (pause)* key to pause the run. Press the *f1 (resume)* key to continue the autosampler run at the point where it was paused, press the *f2 (auto)* key to access the autosampler setup parameters or press the *f3 (end)* key to cancel and end the autosampler run.
  - b. Press the *f*3 (*end*) key to abort the autosampler run. The autosampler arm will return to home position and the meter will return to the measurement mode.
- 6. Once the autosampler run is complete, another autosampler run can be started or the meter can be used as normal. The method that was selected for the previous autosampler run will be used for the next run, unless the operator changes the method in the autosampler setup menu.

Error Code	Description
E036	Bad Value Entered – The value that was entered into the meter was outside of the allowable range of values. Verify the value range and reenter the value.
E200 E201	Communication Error – There is an error in communication between the meter and autosampler. Check cable connections, communication protocols and meter settings and then retry running the autosampler. The baud rate of the meter must be set to 1200 to run the autosampler.
E202	Autosampler Jam – The autosampler cannot properly move the arm assembly. Press the $f1$ (pause) key and clear any obstructions from the autosampler. If the jam does not clear, turn the autosampler off and wait a minute before turning it back on. The autosampler should return to the home position.
E203	Unstable Readings – The measurements taken while using the autosampler were unstable. Check the electrode for proper function and ensure that the electrode cables are properly connected.

#### Autosampler Error Messages

For assistance with Orion products, contact Technical Support by email at <u>wai.techservbev@thermofisher.com</u> or by phone – within the United States call 1-800-225-1480 and outside the United States call +1-978-232-6000 or fax +1-978-232-6031.

For additional product information, contact your local authorized dealer, local Thermo Scientific Orion technical sales representative or contact us using the Water and Laboratory Products (WLP) information on the page back of this manual.



# **SECTION 9** Customer Services

### **Troubleshooting Guide**

#### **Electrode Stability Test**

- In the measurement mode, press the *setup* key to enter the setup menu. Press the ▲ /▼ keys to highlight Instrument Parameters and press the *f2* (*select*) key. Press the ▲ /▼ keys to highlight Troubleshooting and press the *f2* (*select*) key.
- Press the ▲ /▼ keys to highlight Stability Test and press the *f*2 (accept) key. Press the ▲ /▼ keys to highlight Channel 1 or Channel 2 and press the *f*2 (accept) key.
- 3. Make sure that the electrode or probe to be tested is connected to the channel that was selected in the previous step and press the *f*2 (*yes*) key.
- 4. Place the electrode or probe in an appropriate solution with a stable temperature and press the *f*2 (*yes*) key.
  - A pH 7 buffer is recommended for a pH electrode.
  - A calibration standard with a moderate concentration of the ion of interest is recommended for an ion selective electrode (ISE).
  - An ORP standard is recommended for an ORP electrode.
- 5. The meter will perform the stability test and show the mV per minute drift and noise on the display.
- 6. Press the f1 (escape) key to return to the measurement mode.
### **Meter Self Test and Checkout**

A meter self-diagnostic test is performed automatically when the meter is powered on and can be manually activated from the setup menu. The self test that is performed during the meter power up does not require any operator action. If the self test is activated from the setup menu, the operator must press the meter keys when prompted during the test.

#### **Meter Self Test**

- In the measurement mode, press the *setup* key to enter the setup menu. Press the ▲ /▼ keys to highlight Instrument Parameters and press the *f2* (*select*) key. Press the ▲ /▼ keys to highlight Troubleshooting and press the *f2* (*select*) key.
- 2. Press the ▲ / ▼ keys to highlight Self Test and press the *f*2 (accept) key.
- 3. Disconnect all of the electrodes and probes from the meter, attach both BNC shorting caps to the two BNC inputs and press the *f*2 (*yes*) key.
- 4. Press the f2 (yes) key to start the self test.
- 5. The meter will perform an accuracy test. When the meter displays Accuracy Test Passed, press the *f*2 (*next*) key.
- 6. The meter will perform an EPROM test. When the meter displays E2Prom Test Passed, press the *f2* (*next*) key.
- 7. The meter will perform a keypad test. Press each of the blinking keys shown on the display in the order that they are shown. When the meter displays Keypad Test Passed, press the *f*2 (*next*) key.
- 8. The meter will display Self Test Passed. Press the *f1 (escape)* key to return to the measurement mode.

Note: If any problems are found during the self test, refer to the Meter Error Codes section.

### Meter Checkout

- After the self test is complete, the meter will return to the measurement mode. Keep all of the electrodes disconnected from the meter and keep the BNC shorting caps attached to the BNC inputs on the meter.
- Set the measurement mode for both channels to pH. The meter should read a steady 7.000 pH ± 0.002 pH on both channels. If the meter does not read a steady 7.000 ± 0.002, perform a one point pH calibration with the BNC shorting caps attached to the BNC inputs. Set the pH value to 7.000 and the slope value to 100.0. Refer to the <u>pH Calibration</u> <u>Procedure</u> section for detailed instructions.
- In the measurement mode, the meter should read a steady 0.0 mV ± 0.2 mV. If the meter does not read a steady 0.0 mV ± 0.2 mV with the BNC shorting caps attached to the BNC inputs, contact Technical Support.

### **User Reset**

If the user reset is performed, all of the meter settings will be deleted and reset to the default settings. The data log, calibration log and methods will not be deleted with the user reset.

- In the measurement mode, press the *setup* key to enter the setup menu. Press the ▲ / ▼ keys to highlight Instrument Parameters and press the *f*2 (*select*) key. Press the ▲ / ▼ keys to highlight Troubleshooting and press the *f*2 (*select*) key.
- 2. Press the ▲ / ▼ keys to highlight User Reset and press the *f*2 (accept) key.
- 3. Press the f2 (yes) key to perform the reset.
- 4. After the user reset is done, the meter will enter the EZ Startup menu and then proceed to the measurement mode.

### **Factory Configuration**

If the factory configuration is performed, all of the current meter settings will be deleted and reset to the default settings and the data log, calibration log and methods will be deleted.

- In the measurement mode, press the setup key to enter the setup menu. Press the ▲ /▼ keys to highlight Instrument Parameters and press the f2 (select) key. Press the ▲ /▼ keys to highlight Troubleshooting and press the f2 (select) key.
- 2. Press the ▲ /▼ keys to highlight Factory Config and press the *f*2 (accept) key.
- 3. Press the f2 (yes) key to perform the reset.
- 4. Use the alphanumeric keypad to enter the PIN number and press the *f2 (accept)* key to perform the reset. The default PIN number is 11111111.
- 5. Press the *f1 (measure)* key, the meter will enter the EZ Startup menu and then proceed to the measurement mode.

### **About Meter**

- In the measurement mode, press the *setup* key to enter the setup menu. Press the ▲ /▼ keys to highlight Instrument Parameters and press the *f2* (*select*) key. Press the ▲ /▼ keys to highlight Troubleshooting and press the *f2* (*select*) key.
- 2. Press the ▲ /▼ keys to highlight About Meter and press the *f*2 (accept) key.
- 3. The meter serial number and software version will be displayed.
- Press the *f1 (back)* key to return to the troubleshooting settings, press the *f1 (back)* key to return to the instrument parameter settings and press the *f1 (back)* key to return to the measurement mode.

### **Electrode Condition Icon**

The electrode condition icon indicates the performance of the pH electrode or ion selective electrode (ISE), based on the last saved calibration and electrode measurement stability.

lcon	Definition of Icon – pH Electrode	Definition of Icon – ISE Electrode
1 U	Electrode condition is good and the electrode slope is 90 to 110 %.	Electrode condition is good and the electrode slope is 0 to 199 mV/decade (cation) or 0 to -199 mV/decade (anion).
ſ	Electrode condition is fair and the electrode slope is 80 to 120 %.	N/A
۵.	Electrode condition is bad and the electrode slope is less than 80 % or greater than 120 %.	Electrode condition is bad and the electrode slope has the wrong sign or the value is wrong. Verify that the correct electrode ID was selected in the
۲	Consult the electrode user guide for instructions on how to clean, condition and troubleshoot the electrode.	setup menu. Consult the electrode user guide for instructions on how to clean, condition and troubleshoot the electrode.

## Assistance

For assistance with Orion products, contact Technical Support by email at <u>wai.techservbev@thermofisher.com</u> or by phone – within the United States call 1-800-225-1480 and outside the United States call +1-978-232-6000 or fax +1-978-232-6031.

For additional product information, contact your local authorized dealer, local Thermo Scientific Orion technical sales representative or contact us using the Water and Laboratory Products (WLP) information on the page back of this manual.

# Warranty

For the most current warranty information, visit www.thermoscientific.com/water.

## Meter Error Codes

### **Over Range Measurements**

If the measurement is flashing 9999 and over range, the measurement value is outside of the allowable measurement range. Make sure that the electrode is connected to the appropriate meter input and the electrode parameters are entered correctly in the setup menu. Clean the electrode according to the electrode user guide and re-calibrate the electrode with new buffers or standards. If the error persists, perform the meter self test.

Error Code	Description	
E001	Display Error – If this error occurs, contact Technical Support.	
E002	Hardware Error – If this error occurs, contact Technical Support.	
E003	<b>Channel 1 Input Error</b> – Disconnect all of the electrodes from the meter, connect the two BNC shorting caps to the channel 1 and channel 2 BNC inputs and repeat the self test.	
E004	<b>Channel 2 Input Error</b> – Disconnect all of the electrodes from the meter, connect the two BNC shorting caps to the channel 1 and channel 2 BNC inputs and repeat the self test.	
E005	<b>Reference Input Error</b> – Make sure that all of the electrodes are disconnected from the meter (including the reference input), connect the two BNC shorting caps to the channel 1 and channel 2 BNC inputs and repeat the self test.	
E006	<b>Ground Input Error</b> – Make sure that all of the electrodes are disconnected from the meter (including the ground input), connect the two BNC shorting caps to the channel 1 and channel 2 BNC inputs and repeat the self test.	
E007	<b>Keypad Failure</b> – Repeat the self test. When the meter prompts, press each of the keys one at a time, including the power key, within five seconds of the meter prompts. If the error persists, contact Technical Support.	
E008	Hardware Error – If this error occurs, contact Technical Support.	
E009	Hardware Error – If this error occurs, contact Technical Support.	
E020	<b>Measurement Out of Range</b> – Perform the meter self test, clean the electrode according to the electrode user guide and re-calibrate the electrode with new buffers or standards.	
E027	<b>Data Transfer Error</b> – A communication error between the meter and computer occurred. Check the cable connections and communication protocols. Make sure that the meter is set to the correct baud rate and the computer function is enabled in the setup menu.	
E028	Incremental Techniques Error – Make sure that the correct incremental technique was used and all of the parameters were entered correctly. The measured millivolt difference may not be great enough to permit the calculation of the unknown concentration. Increase the concentration of the standard or dilute the sample. The concentration of the unknown may have been calculated to be less than 0.0001. If the unknown was diluted, use a fresh aliquot of undiluted or less diluted unknown. The electrode ID may be incorrect or the electrode may not have been prepared correctly, filled with the proper filling solution or connected to the appropriate meter input. If the error persists, measure a standard with a known concentration and the same background as the sample. Adjust the standard concentration and volume until accurate results are obtained.	
E032	<b>Printer Communication Error</b> – A communication error between the meter and printer occurred. Check the cable connections and communication protocols. Make sure that the meter is set to the correct baud rate for the printer and the printer function is enabled in the setup menu. The printer may be out of paper, the printer paper may be jammed or there may be another printer error causing the printer not to be ready.	

Error Code	Description	
E035	<b>Incorrect Password Entered</b> – The password that was entered into the meter was incorrect. Verify the password and reenter it.	
E036	<b>Bad Value Entered</b> – The value that was entered into the meter was outside of the allowable range of values. Verify the value range and reenter the value.	
E038	<b>Calibration Log Full</b> – The calibration log is full and must be cleared before any additional calibration data can be saved.	
E039	<b>Calibration Log Empty</b> – There is no calibration saved in the calibration log and a calibration needs to be performed before any calibration data can be viewed.	
E040	Data Log Full – The data log is full and must be cleared before any data points can be saved.	
E041	<b>Data Log Empty</b> – There are no data points saved in the data log. The data log needs to be enabled in the setup menu and a measurement needs to be logged before any data points can be viewed.	
E105	<b>pH Auto Buffer Recognition Error</b> – The millivolts measured for a pH buffer during calibration are outside of the required range for automatic buffer recognition. Review the pH calibration procedure and check that the correct buffer set was selected in the setup menu. Clean the electrode according to the electrode user guide. Re-calibrate the electrode with fresh buffers.	
E107	<b>pH Calibration Standard Error</b> – The millivolts measured during calibration are the same for two buffers or the millivolts for the two buffers are too close together. There must be at least a 5 mV difference between the two buffers. Review the pH calibration procedure and verify that the electrode was placed in the proper buffers at the appropriate times. Clean the electrode according to the electrode user guide. Re-calibrate the electrode with fresh buffers.	
E109	<b>Bad pH Slope or Calibration Offset</b> – Clean the electrode according to the electrode user guide. Re-calibrate the electrode with fresh buffers.	
E306	<b>ISE Automatic Blank Error</b> – Review the ISE calibration procedure and verify that the electrode was placed in the proper standards at the appropriate times. Clean the electrode according to the electrode user guide. Re-calibrate the electrode with fresh standards. If the error persists, disable the autoblank feature in the setup menu and re-calibrate the meter without a zero standard. See the Automatic Blank Correction Feature section for more information.	
E307	<b>ISE Calibration Standard Error</b> – The millivolts measured during calibration are the same for two standards or the millivolts for the two standards are too close together. There must be at least a 20% difference between the entered standard values. Review the ISE calibration procedure and verify that the electrode was placed in the proper standards at the appropriate times. Clean the electrode according to the electrode user guide. Re-calibrate the electrode with fresh standards.	
E309	<b>Bad ISE Slope</b> – Verify that the correct electrode ID was selected in the setup menu. Clean the electrode per the electrode user guide. Re-calibrate the electrode with new standards.	
E511	<b>ORP Calibration Standard Error</b> – The millivolts measured for the ORP standard during calibration are outside of the required range for an ORP calibration. Review the ORP calibration procedure and verify that the electrode was placed in the standard at the appropriate time and the Thermo Scientific ORP standard was used. Clean the electrode according to the electrode user guide. Re-calibrate the electrode with fresh ORP standard.	

## Notice of Compliance

This equipment generates, uses, and can radiate radio frequency energy and if not installed and used in accordance with the instruction manual, may cause interference to radio communications. It has been tested and found to comply with the limits for a Class A computing device pursuant to Subpart J of Part 15 of FCC Rules, which are designed to provide reasonable protection against such interference when operated in a commercial environment. Operation of this equipment in a residential area is likely to cause interference, in which case the user, at his own expense, will be required to take whatever measures may be required to correct the interference.

"This digital apparatus does not exceed the (Class A) limits for radio noise emissions from digital apparatus set out in the Radio Interference Regulations of the Canadian Department of Communications."

"Le présent appareil numérique n' émet pas de bruits radioélectriques dépassant les limites applicables aux appareils numériques (de la class A) prescrites dans le Règlement sur le brouillage radioélectrique édicté par le ministère des Communications du Canada."

## WEEE Compliance

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the symbol shown here.



Further information on compliance with these directives, the recyclers in your country, and information on Thermo Scientific Orion products that may assist the detection of substances subject to the RoHS Directive are available by contact us using the Water and Laboratory Products (WLP) information on the back page of this manual.

## **Declaration of Conformity**

Manufacturer: Thermo Fisher Scientific Inc Address: Ayer Rajah Crescent Blk 55 #04-16/24 Singapore 139949 Singapore

#### Hereby declares that the following product:

Orion Dual Star pH/ISE Benchtop Meter is rated 100 to 240 VAC, 50/60 Hz, 0.5A

**Equipment Class:** 

Measurement, control and laboratory Orion Dual Star meters are EMC Class A

#### Conforms with the following directives and standards:

EN61326-1:2013	Electromagnetic Compatibility (EMC Directive)
	Electrical equipment for measurement, control and laboratory use -
	EMC requirements
EN61010-1:2010	Safety Standards
UL61010-1:2012	Safety requirements for electrical equipment for measurement,
CAN/CSA C22.2 No. 61010-1:2012	control and laboratory use - general requirements.



Cheow Kwang Chan QA/Regulatory Manager Place and Date of Issue: 14 May 2015 Singapore

## Real Time Clock Battery Maintenance

The Orion Dual Star meter uses an internal BR-1225 type, 3 V lithium coin cell battery for real time clock operation. If the real time clock function is lost, the battery needs to be replaced. Since the battery is inside of the meter housing, only qualified technicians should attempt to replace the battery.

### **Replacing and Disposing of the Battery**

- 1. Disconnect the universal power adapter from the meter and the wall outlet.
- 2. Orientate the meter so the display is face down on a soft, nonmetallic, non-scratching surface. The bottom of the meter (with the pullout guide) should be closest to you and the top of the meter (with the electrode and cable inputs) should be farthest from you.
- Use a medium size Phillips head screwdriver to unscrew the four screws on the back of the meter. Remove the back cover of the meter. Be careful not to spill any liquids, dust or other foreign matter onto the exposed circuit board.
- 4. Locate the battery holder. It is a round, small, coin size battery holder that is located in the left bottom corner of the circuit board and has a reference designator BAT41 marked in white on the circuit board.
- 5. Use a small flat screwdriver to pry out the battery from the holder. Slide the new battery, from the side, into the battery holder and be careful not to break the metal tab that holds the battery in place.
- 6. Place the back cover on the meter and fasten the screws.
- 7. Dispose of the used battery in safe and environmentally responsible way, following all applicable federal, state and local regulations.

**Important Note:** Replace the battery only with a new BR-1225 type, 3 V lithium coin cell battery. Do not substitute any other battery type.

**Battery Recommendations**: Use caution, lithium can cause fire, explosion and burns. Follow all applicable regulations for disposal of products containing lithium. Do not tamper with the batteries or try to open them. Do not crush, recharge, incinerate or heat the batteries above 212 °F (100 °C). Do not expose the batteries to heat sources. Burning the batteries will result in corrosive and toxic lithium hydroxide gas release.

# Ordering Information

Cat. No.	Description	
2115000	Orion Dual Star pH/ISE benchtop meter with 927007MD Orion stainless steel ATC probe, electrode stand, USB cable, universal power supply, literature & test certificate	
2115001	Orion Dual Star pH/ISE benchtop meter kit with 8102BNUWP ROSS Ultra glass-body pH electrode, 927007MD Orion stainless steel ATC probe, 910104 Orion pH 4.01 buffer (475 mL), 910107 Orion pH 7.00 buffer (475 mL), 910110 Orion pH 10.01 buffer (475 mL), 810001 ROSS storage solution (475 mL), electrode stand, USB cable, universal power supply, literature & test certificate	
2115101	Orion Dual Star pH/ISE benchtop meter & electrode set with 8102BNUWP ROSS Ultra glass-body pH electrode, 927007MD Orion stainless steel ATC probe, electrode stand, USB cable, universal power supply, literature & test certificate	
2115102	Orion Dual Star pH/ISE benchtop meter & Triode set with 8157BNUMD ROSS Ultra Triode epoxy-body pH/ATC electrode, electrode stand, USB cable, universal power supply, literature & test certificate	
2115201	Orion Dual Star pH/ISE benchtop meter pH & ammonia kit with 9512BNWP Orion ammonia ion selective electrode, 8172BNWP ROSS Sure-Flow glass-body pH electrode, 927007MD Orion stainless steel ATC probe, 096019 Orion Star series stirrer probe, 951207 Orion 100 ppm ammonia standard (475 mL), 951211 Orion ammonia ISA ionic strength adjuster (475 mL), 810199 ROSS pH buffer and solution kit, electrode stand, USB cable, universal power supply, literature & test certificate	
2115202	Orion Dual Star pH/ISE benchtop meter pH & chloride kit with 9617BNWP Orion chloride ion selective electrode, 8172BNWP ROSS Sure-Flow glass-body pH electrode, 927007MD Orion stainless steel ATC probe, 096019 Orion Star series stirrer probe, 941708 Orion 1000 ppm chloride standard (475 mL), 940011 Orion chloride ISA ionic strength adjuster (475 mL), 810199 ROSS pH buffer and solution kit, electrode stand, USB cable, universal power supply, literature & test certificate	
2115203	Orion Dual Star pH/ISE benchtop meter pH & fluoride kit with 9609BNWP Orion fluoride ion selective electrode, 8172BNWP ROSS Sure-Flow glass-body pH electrode, 927007MD Orion stainless steel ATC probe, 096019 Orion Star series stirrer probe, 040908 Orion 10 ppm fluoride standard premixed with TISAB II (475 mL), 040906 Orion 1 ppm fluoride standard premixed with TISAB II (475 mL), 040906 Orion 1 ppm fluoride standard premixed with TISAB II (475 mL), 940909 Orion TISAB II total ionic strength adjustment buffer (3.8 L), 810199 ROSS pH buffer and solution kit, electrode stand, USB cable, universal power supply, literature & test certificate	
2115204	Orion Dual Star pH/ISE benchtop meter pH & nitrate kit with 9707BNWP Orion nitrate ion selective electrode, 8172BNWP ROSS Sure-Flow glass-body pH electrode, 927007MD Orion stainless steel ATC probe, 096019 Orion Star series stirrer probe, 930707 Orion 100 ppm nitrate standard (475 mL), 930711 Orion nitrate ISA ionic strength adjuster (475 mL), 810199 ROSS pH buffer and solution kit, electrode stand, USB cable, universal power supply, literature & test certificate	
2115205	Orion Dual Star pH/ISE benchtop meter pH & sodium kit with 8611BNWP ROSS sodium ion selective electrode, 8172BNWP ROSS Sure-Flow glass-body pH electrode, 927007MD Orion stainless steel ATC probe, 096019 Orion Star series stirrer probe, 841108 Orion 1000 ppm sodium standard (475 mL), 841109 Orion KAP sodium standard & ISA solution (475 mL), 841111 Orion sodium ISA ionic strength adjuster (475 mL), 841113 Orion sodium electrode conditioning solution (475 mL), 841101 Orion sodium electrode storage solution (475 mL), 810199 ROSS pH buffer and solution kit, electrode stand, USB cable, universal power supply, literature & test certificate	

Cat. No.	Description	
STARA-BEA	Meter-attachable electrode stand, includes electrode arm, holder & meter bracket	
STARA-HB	Freestanding weighted base for use with meter-attachable electrode stand	
090043	Swing arm electrode stand, includes electrode arm, holder & weighted base	
810017 Storage sleeve and base for 12mm diameter electrodes		
1010003	Universal power adapter for Star A series meters	
1010004	RS232 computer interface cable and printer interface cable set	
1010053	Star series RS232 computer cable	
090045	BNC shorting cap	
1010006	Orion Star series inkjet printer with RS232 printer interface cable	
1010008	Ink cartridge for the Orion Star series inkjet printer	
1010009	Replacement paper for the Orion Star series inkjet printer	
096019	Orion Star stirrer probe, pin tip connector	
810199	Orion ROSS All-in-One pH buffer kit, includes 475 mL bottle each of pH 4.01, 7.00, 10.01 buffers; ROSS pH electrode storage solution (810001); 30 mL bottle of pH electrode cleaning solution D; and pH electrode storage bottle	
910199	All-in-One pH buffer kit, includes 475 mL bottle each of pH 4.01, 7.00, 10.01 buffers; pH electrode storage solution (910001); and pH electrode storage bottle	
927007MD	Orion ATC temperature probe with stainless steel body, MiniDIN connector	
927005MD	Orion ATC temperature probe with epoxy body, MiniDIN connector	
928007MD	Orion micro ATC temperature probe with stainless steel tip, MiniDIN connector	
8102BNUWP	ROSS Ultra glass-body refillable pH electrode, BNC connector	
8156BNUWP	ROSS Ultra epoxy-body refillable pH electrode, BNC connector	
8172BNWP	ROSS Sure-Flow glass-body refillable pH electrode, BNC connector	
8165BNWP	ROSS Sure-Flow epoxy-body refillable pH electrode, BNC connector	
8302BNUMD	ROSS Ultra Triode glass-body refillable pH/ATC electrode, BNC & MiniDIN connectors	
8157BNUMD	ROSS Ultra Triode epoxy-body refillable pH/ATC electrode, BNC & MiniDIN connectors	
8107BNUMD	ROSS Ultra Triode epoxy-body gel-filled pH/ATC electrode, BNC & MiniDIN connectors	
8135BNUWP	ROSS Ultra epoxy-body refillable pH electrode with flat surface bulb, BNC connector	
8163BNWP	ROSS spear tip glass-body refillable pH electrode, BNC connector	
8103BNUWP	ROSS Ultra semi-micro glass-body refillable pH electrode, BNC connector	
8115BNUWP	ROSS Ultra semi-micro epoxy-body refillable pH electrode, BNC connector	
8220BNWP	ROSS micro glass-body refillable pH electrode, BNC connector	
810199	ROSS All-in-One pH buffer and storage solution kit, includes pH 4, 7, 10 buffers; ROSS storage solution; pH electrode cleaning solution; pH electrode storage bottle	
810001	ROSS pH electrode storage solution, 475 mL	
910001	Standard pH electrode storage solution, 475 mL	
910168	Orion pH 1.68 buffer, 475 mL	

Cat. No.	Description
910104	Orion pH 4.01 buffer, 475 mL
910105	Orion pH 5.00 buffer, 475 mL
910686 Orion pH 6.86 buffer, 475 mL	
910107 Orion pH 7.00 buffer, 475 mL	
910918	Orion pH 9.18 buffer, 475 mL
910110	Orion pH 10.01 buffer, 475 mL
910112	Orion pH 12.46 buffer, 475 mL
9678BNWP	Orion Sure-Flow ORP epoxy-body refillable electrode, BNC connector
9180BNMD	Orion Triode ORP/ATC epoxy-body refillable electrode, BNC & MiniDIN connectors
967901	Orion ORP standard solution, 475 mL
967961	Orion ORP standard solution, 5 x 60 mL
9512HPBNWP	Orion high-performance ammonia ion selective electrode, BNC connector
9512BNWP Orion standard ammonia ion selective electrode, BNC connector	
951007 Orion 1000 ppm ammonia standard, 475 mL	
951211	Orion ammonia ionic strength adjuster (ISA) with pH-indicating blue dye, 475 mL
951210	Orion low level ammonia ISA with pH-indicating blue dye, 475 mL
9609BNWP	Orion fluoride ion selective electrode, BNC connector
940907	Orion 100 ppm fluoride standard, 475 mL
940909	Orion TISAB II total ionic strength adjustment buffer for fluoride analysis, 3.8 L
040906	Orion 1 ppm fluoride standard premixed with TISAB II, 475 mL
040907	Orion 2 ppm fluoride standard premixed with TISAB II, 475 mL
040908	Orion 10 ppm fluoride standard premixed with TISAB II, 475 mL
9707BNWP	Orion nitrate ion selective electrode, BNC connector
920707	Orion 1000 ppm nitrate standard, 475 mL
930711	Orion nitrate ionic strength adjuster (ISA), 475 mL
930710	Orion nitrate interference suppressor solution, 475 mL
8611BNWP	ROSS sodium ion selective electrode, BNC connector
841108	Orion 1000 ppm sodium standard, 475 mL
841111	Orion sodium ionic strength adjuster (ISA), 475 mL

Visit <u>www.thermoscientific.com/water</u> for a complete listing of all available Thermo Scientific Orion meters, electrodes, solutions and accessories.

# Meter Specifications

Orion Dual Star pH/ISE Benchtop Meter Operating Conditions		
Operating Ambient Temperature	5 to 45 °C	
Operating Relative Humidity	5 to 85%, non-condensing	
Storage Temperature	-20 to 60 °C	
Storage Relative Humidity	5 to 85%, non-condensing	
Pollution	Degree 2	
Overvoltage	Category II	
Meter Weight	1.15 kg (2.5 lb)	
Meter Dimensions (H x W x D)	10 cm x 18 cm x 23 cm (3.9" x 7.1" x 9.1")	
Regulatory and Safety	CE, TUV 3-1, FCC Class A	
Power Rating	DC input: 9VDC, 1A	
Shock and Vibratian	Shock: drop test in packaging per ISTA #1A	
	Vibration: shipping/handling per ISTA #1A	
Enclosure (designed to meet)	IP54	
Warranty	3 years	

Universal Power Adapter Operating Conditions	
Operating Ambient Temperature	0 to 50 °C
Operating Relative Humidity	0 to 90%, non-condensing
Storage Temperature	-20 to 75 °C
Storage Relative Humidity	0 to 90%, non-condensing
Pollution	Degree 2
Overvoltage	Category II

Orion Dual Star pH/ISE Benchtop Meter Specifications	
Measurement Channels	2 – pH, mV, RmV, ORP or ion concentration with temperature
Display	Graphic LCD with adjustable backlight (5 levels)
View Options	View 1 or 2 measurement channels simultaneously
Measurement Modes	Auto-Read, On Ready, At Time Intervals, Continuous, Value Change In Measurement and Timed Reading
Data Log Memory	1000 data points with time and date stamp
Data Log Functions	Automatic data logging with Auto-Read, On Ready, At Time Intervals, Value Change In Measurement and Timed Reading; manual data logging with Continuous measurement mode
Methods	20 (10 per channel with calibration data, password protection)
Calibration Log	10 calibrations per channel with time and date stamp

Computer Software		Orion Dual Star Com data transfer software
	Range	- 2.000 to 19.999
	Resolution	0.1 / 0.01 / 0.001
	Relative Accuracy	±0.002
	Calibration Points	1 to 6
рН	Auto-Buffer Recognition	US/NIST (pH 1.68, 4.01, 7.00, 10.01, 12.46 buffers) EURO/DIN (pH 1.68, 4.01, 6.86, 9.18 buffers)
	Calibration Editing	Yes
	Calibration Options	Calibration graph with point to point slopes and average slope
	Input Impedance	> 10^12 ohms
	Electrode Status	Onscreen indicator for good, fair or bad electrode status
	Range	0.0001 to 19900
	Resolution	0.0001 minimum, 1 to 3 significant digits (user selectable)
	Relative Accuracy	$\pm 0.2$ mV or $\pm 0.05\%$ of reading, whichever is greater
	Units	ppm, mg/L, %, M or unitless
ISE (Ion Concentration)	Calibration Points	2 to 6
,	Calibration Editing	Yes
	Incremental Techniques	Single known addition & subtraction, double known addition & subtraction
	Advanced Features	Segmented (point-to-point) slope, non-linear selectable auto- blank, low concentration range stability
	Range	±1999.9 mV
	Resolution	0.1 mV
mV / RmV	Relative Accuracy	$\pm 0.2$ mV or $\pm 0.05\%$ of reading, whichever is greater
	Relative mV Mode	Yes
	E <sub>H</sub> ORP Mode	Yes
	Range	-5 to 105 °C, 22 to 221 °F
Temperature	Resolution	0.1 °C, 0.1 °F
	Relative Accuracy	±0.1 °C
	Offset Calibration	1 point
	Source Options	Manual or automatic with ATC probe
Sancar Innuts	BNC (2)	pH electrode, ORP electrode or ion selective electrode (ISE)
	Pin Tip (2)	Reference electrode
Sensor inputs	8 pin MiniDIN (2)	ATC temperature probe
	Stirrer Jack	Orion Star series stirrer probe
Outputs	Ports	USB, RS232



## **SECTION 10** Advanced Meter Features

## Multipoint Calibration Curve Feature

The Orion Dual Star meter utilizes a point-to-point scheme for multipoint calibrations in both pH and concentration modes. In this technique, straight line segments are drawn between each successive pair of calibration points. Separate slope and Eo values are calculated using the adjacent pair of data points. For example, in a five point calibration, four slopes and five Eo values are calculated to correspond to the four line segments that connect the data. Concentration or pH values for samples are calculated according to the segment that their potential values fall into. Those above or below the range of calibration are calculated by extrapolation from the top or bottom most segment.

The displayed slope is an average slope of the separate segments. The slope may be outside the range normally expected for an electrode. For example, if the slope of one segment is low, then the average will be lower than theoretical, even though the response may be Nernstian over most of the range.

The following figure is an example of a typical calibration curve using this method.



### Automatic Blank Correction Feature

The Orion Dual Star meter has a feature known as automatic blank correction. In this technique, the meter decides whether blank correction is the best measurement strategy by analyzing the electrode response during a multipoint calibration. A separate blank does not have to be run. Graphically, automatic blank correction is equivalent to drawing a smooth curve through the lowest three points of the multipoint calibration and extrapolating to zero concentration based on the assumption of Nernstian electrode behavior.

Multipoint calibrations at the lower limit of detection are desirable when the response of an ion selective electrode is non-linear and cannot be characterized with a one or two point calibration. This is usually seen as a low electrode slope. Generally, the electrode is behaving in a Nernstian manner but the effect of a blank is being observed. See the figure below.





This blank may be a true reagent blank, traces of analyte ion in the reagents or it may be the "mud" value of the electrode. It could also be an interference in the reagents that becomes apparent at low levels of analyte ion, or it could be any combination of these effects.

The expanded version of the Nernst equation traditionally used for blank correction is as follows:

 $E = E_0 + S \times \log (C + b)$  where b is the blank

In a multi-point calibration, a set of equations is generated and the relationship between them evaluated. For example the equations generated in a three point calibration would be as follows:

 $E_{1} = E_{0} + S \times \log (C_{1} + b)$  $E_{2} = E_{0} + S \times \log (C_{2} + b)$  $E_{3} = E_{0} + S \times \log (C_{3} + b)$ 

The meter evaluates the relationships between the three potentials  $E_1$ ,  $E_2$  and  $E_3$  and the three concentrations  $C_1$ ,  $C_2$  and  $C_3$ . If the relationships indicate that blank correction is desired, a blank will automatically be calculated and the non-linearity will be corrected for in a Nernstian manner. If the appropriate conditions are not met, the blank is set at zero and each segment of the multipoint calibration is treated independently.

When all three of the following conditions are met, blank correction is invoked.

- 1. The concentration of the first standard is zero, or the slope of the electrode between the first and second standards is less than the slope between the second and third.
- 2. Potential differences between points are significant. For example,
- 3.  $E_3 E_1 > 10 \text{ mV}$
- 4. The blank correction algorithm converges at reasonable blank and slope values. Conditions in steps 1 and 2 prevent failure to converge in most situations. However, the slope will attain any value necessary for convergence to a calculated blank value of 3 x C<sub>3</sub>.

When blank correction is implemented, the slope value for the electrode may be outside the range of values normally considered acceptable during an ordinary calibration. For best results, calibration standards should be close in range to the expected sample concentrations and should bracket the expected sample concentration.

When conditions 1, 2 and 3 are not met, the calibration data is handled by the multipoint calibration method described earlier.

**Note:** When a calibration standard is defined as zero and the meter cannot calculate a blank, an error code E306 will be displayed.

In calibrations with more than three points, a combination of methods are utilized. Automatic blank correction is used if the lowest three points satisfy the criteria and multipoint calibration is used for the other points. The slope displayed after calibration is an average of the slope calculated in the automatic blank correction algorithm and the slope(s) for each additional segment.

#### **Error Code E306**

#### With the Autoblank Function On

This error occurs during a three or more point calibration when one solution is defined as zero concentration and the actual concentration of the blank is substantially greater than the standard with the third highest concentration. Since the blank cannot be calculated, the meter cannot use the autoblank function and cannot use the zero concentration standard in the two point calibration segment. To clear the error, calibrate the meter without using a zero concentration standard.

#### With the Autoblank Function Off

This error occurs during a one or two point calibration if one standard is defined as zero concentration. To clear the error, calibrate the meter without using a zero concentration standard.

## **Isopotential Point Feature**

To take advantage of the temperature compensation feature for concentration measurements in the Orion Dual Star meter, it is necessary to experimentally determine the isopotential point of the electrode in use. The slope of all electrodes changes with temperature. This is due in fact to the slope term, S, in the Nernst equation. The slope is actually 2.3 RT/nF, where R and F are constants, n is the charge on the measured species and T is the temperature in degrees Kelvin. Examples of calibration curves at varying temperatures are shown in the figure below.

These curves intersect at the isopotential point, which is the concentration at which the potential of the electrode does not vary with temperature. If the isopotential point is known or can be measured experimentally, and if the meter has the means of adjusting the isopotential point, temperature compensation for an ion selective electrode is possible. Temperature compensation can be performed for one and two point calibrations.



### **Determining an Isopotential Point**

- 1. Prepare several standards with concentration ranges over the measuring range of the electrode.
- 2. Measure the millivolt value of the standards at room temperature, about 20 to 25 °C.
- 3. Measure the millivolt value of each standard at 75 °C.
- 4. Measure the millivolt value of each standard at 10 °C.
- On semi-logarithmic graph paper, plot the concentration values on the log axis versus the millivolt values on the linear axis, as any calibration curve would be drawn (see the figure on the previous page).
- 6. The lines will intersect at the isopotential point. Read the concentration off the graph for this point from the log axis. This value is the isopotential point.
- 7. See the *Setup Menu* section for instructions on how to adjust the isopotential point on the Orion Dual Star meter.

## Automatic pH Buffer Recognition Feature

The Orion Dual Star meter is capable of automatically recognizing pH 1.68, 4.01, 7.00, 10.01 and 12.46 buffers or pH 1.68, 4.01, 6.86, and 9.18 buffers during a pH calibration, depending on the pH buffer set that is selected in the setup menu. During a calibration, the meter uses the selected buffer set and the raw mV reading of the pH electrode in the buffer to recognize and display the buffer value at the measured temperature. The raw mV reading of the pH electrode in the buffer in order for the meter to automatically recognize the buffer.

### USA pH Buffer Set

Buffer	mV Range
1.68	+285 to +345 mV
4.01	+207 to +147 mV
7.00	- 30 to + 30 mV
10.01	-207 to -147 mV
12.46	-293 to -353 mV

### EURO pH Buffer Set

Buffer	mV Range
1.68	+285 to +345 mV
4.01	+207 to +147 mV
6.86	+38 to -22 mV
9.18	-99 to -159 mV

### Testing a pH Electrode for Automatic Buffer Recognition

To verify that the raw mV reading of the pH electrode in use is  $\pm$  30 mV from the theoretical mV reading of the buffer, and therefore verify that the pH electrode in use is capable of performing automatic buffer recognition, perform the following procedure.

- 1. Set the measurement mode to mV in the setup menu. Prepare the pH electrode according to the instructions in the pH electrode user guide. Ensure that the pH electrode is connected to the meter input that corresponds with the channel that will be measured.
- 2. Rinse the pH electrode with distilled water, blot it dry with a lint-free tissue and place the pH electrode into a pH 4.01 buffer at 25 °C.
- 3. Wait for the measurement to stabilize and record the mV value of the pH 4.01 buffer when the measurement is stable. The read type selected in the EZ Startup menu or setup menu will determine how the mV measurements are displayed by the meter.
- 4. Remove the pH electrode from the pH 4.01 buffer, rinse the pH electrode with distilled water, blot it dry with a lint-free tissue and place it into a pH 7.00 buffer at 25 °C.
- 5. Wait for the measurement to stabilize and record the mV value of the pH 7.00 buffer when the measurement is stable.
- 6. The mV reading of the pH electrode in pH 4.01 buffer should be +207 to +147 mV and the mV reading of the pH electrode in pH 7.00 buffer should be -30 to +30 mV. If the mV readings are in the correct ranges, the pH electrode is capable of performing automatic buffer recognition. If the mV readings of the pH electrode are not in the correct ranges, the values of the pH buffers will need to be manually entered during a pH calibration.

#### thermoscientific.com/water

© 2015 Thermo Fisher Scientific Inc. All rights reserved. Microsoft and Excel are either registered trademarks or trademarks of Microsoft Corporation. All other trademarks are the property of Thermo Fisher Scientific and its subsidiaries.

Water and Lab Products

North America Toll Free: 1-800-225-1480 Tel: 1-978-232-6000 Info.water@thermofisher.com

Germany Tel: (49) 6184-90-6000 info.water.uk@thermofisher.com

China Tel: (86) 21-68654588 wai.asia@thermofisher.com India Tel: (91) 22-4157-8800 wai.asia@thermofisher.com

Singapore Tel: (65) 6778-6876 wai.asia@thermofisher.com Japan Tel: (81) 045-453-9175 wai.asia@thermofisher.com

Australia Tel: (613) 9757-4300 In Australia (1300) 735-295 InfoWaterAU@thermofisher.com

