Data Log

VERSA STAR meters offer a 2000 point data log. Each point includes measurements from one to four channels with the date and time. When the data log function is turned on in the Instrument Settings setup menu, the read type determines how the displayed measurements are saved to the data log.

- 1. In the measurement mode, press the setup key
- 2. Press the *◄* or *▶* key to highlight Log View and press the f3 (Select) key.
- 3. Press the \triangleleft or \triangleright key to highlight Data Log.
- 4. Press the \blacktriangle or \checkmark key to highlight View and press the f3 (Select) key.
- 5. Press the \blacktriangle or \checkmark key to highlight a point and press the enter key to view detailed data for that point. Press the ✓ or ► key to scroll through additional data log points.
- 6. Press the measure (esc) key to return to the measurement mode.

thermoscientific.com/water

© 2014 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries

Tel: (91) 22-4157-8800

Tel: (86) 21-68654588

Tel: (81) 045-453-9175

wai asia@thermofisher.com

wai asia@thermofisher.com

wai.asia@thermofisher.com

Australia

Tel: (613) 9757-4300

in Australia (1300) 735-295

InfoWaterAU@thermofisher.com

India

China

Japan

Water Analysis Instruments

North America

Toll Free: 1-800-225-1480 Tel: 1-978-232-6000 info.water@thermo.com

Netherlands

Tel: (31) 020-4936270 info water uk@thermo.com

Singapore Tel: (65) 6778-6876

wai asia@thermofisher.com

Calibration Log

The calibration log contains up to the ten most recent calibrations per pH. RmV. ORP. ISE, incremental technique, conductivity, DO and RDO parameter.

- 1. In the measurement mode, press the setup key.
- Press the ◄ or ► key to highlight Log View and press the f3 (Select) key.
- 3. Press the \triangleleft or \blacktriangleright key to highlight Cal Log.
- 4. Press the \blacktriangle or \checkmark key to highlight the desired calibration parameter and press the f3 (Select) key.
- 5. Press the \blacktriangle or \checkmark key to view each calibration for the selected parameter.

SCIENTIFIC

A Thermo Fisher Scientific Brand

6. Press the measure (esc) key to return to the measurement mode.



Thermo Scientific Orion

VERSA STAR Conductivity/TDS/Salinity/Resistivity Temperature Module

This literature provides basic instructions on operating the Thermo Scientific[™] Orion[™] VERSA STAR[™] meter when the VERSA STAR conductivity measurement module is installed. For comprehensive information on meter and module system setup, operation and advanced features, please refer to the VERSA STAR user manual available on the VERSA STAR literature CD or www.thermoscientific.com/water.

Preparation

- 1. Prepare the universal power adapter, install the meter-attached electrode stand and verify that the conductivity module is connected 8 pin to the meter. If it is not connected, insert the module into an available MiniDIN channel on the back of the meter.
- Prepare the electrodes as instructed in the electrode user manuals. For improved movement and control, place the electrodes into the electrode stand.
- 3. Connect the conductivity probe cable to the 8 pin MiniDIN input on the module.





IS-68X591203-B 0514

- 4. If using the Thermo Scientific[™] Orion[™] Star[™] stirrer probe, attach the stirrer probe cable to the STIRRER 1 input (default input for channels 3 & 4) or STIRRER 2 input (default input for channels 1 & 2) on the meter.
- 5. Set up the work area with calibration standards, rinse water, samples and other supplies.
- 6. If applicable, connect the meter to an external device using the appropriate cable. A USB cable is included with the meter for computer interfacing.
- 7. Power on the meter by connecting the universal power adapter to the meter and power outlet.

Conductivity Mode Setup Menu

Within the Channel 1, 2, 3 & 4 setup menus are Method, Mode and Temperature submenus, which can be used to customize measurement settings and parameters for the selected channel.

- 1. In the measurement mode, press the setup key.
- Press the ◄ or ► key to highlight the appropriate <u>Channel</u> setup menu and press the f3 (Select) key.
- 3. Press the \blacktriangleleft or \blacktriangleright key to highlight <u>Mode</u> and press the *f3 (Select)* key.
- Press the ▲ or ▼ key to highlight <u>Conductivity</u> and press the *f3* (*Select*) key.
- 5. View and update the displayed menu options.

Conductivity Setup Menu

Default values are in bold. The mV, Relative TDS, Salinity and Resistivity menus are similar with options relevant to the mode.

Read Type	AutoRead, Timed, Single-Shot, Continuous
Cell K	0.475
Temp. Compensation	Linear, nLFn, nLFu, EP, Off
Ref. Temp	5°C, 10°C, 15°C, 20°C, 25°C
Temperature Coeff.	2.10
Stability	Smart, Fast, Medium, Slow

Averaging	Off, Automatic Smart
Alarm	Limit (Off), CalDue (Off), Set Point (Off)
Cell Type	Standard, USP
Electrode SI No	(no value)
Sample ID	Off, Manual, Auto Increment

Conductivity Calibration

- The Orion 100 μ S/cm, 1413 μ S/cm and 12.9 mS/cm conductivity standards can be automatically recognized during calibration when the nominal cell constant (Cell K) of the conductivity probe is entered in the Conductivity Mode setup menu.
- 1. In the measurement mode, press the *f1 (Cal)* key to start the calibration.
 - a. When measuring more than one channel, press the

 ▲ or ▼ key to highlight the channel to be calibrated
 and press the *f3 (Select)* key.
- Rinse the conductivity probe and any other electrodes in use with deionized water, blot dry with a lint-free tissue and place into the standard.
- 3. When the conductivity probe and standard are ready, press the *f3 (Start)* key.
 - The stirrer probe will start stirring and continue until the reading stabilizes.
- 4. Wait for the conductivity value on the meter to stabilize and perform one of the following actions:
- a. Press the *f2 (Accept)* key to accept the displayed value, or
- b. Press the *f3 (Edit)* key, press the *f3 (Clear)* key and use the numeric keypad to enter the conductivity value of the standard at the measured temperature.
 Press the *f2 (Accept)* key to confirm the entered value.
- Press the *f2 (Next)* key to proceed to the next standard and repeat steps 2 through 4 or press the *f3 (Cal Done)* key to save and end the calibration.

6. The meter will display the calibration summary and export the data to the calibration log. Press the *measure (esc)* key to proceed to the measurement mode.

Measurement

The read type selected for each channel will determine how measurements are displayed and saved (data log must be on in the Instrument Settings setup menu).

Auto-Read – Press the *measure (esc)* key to start a measurement. When the measurement is stable, the AR icon will stop flashing and the measurement will be locked on the display and saved to the data log. Press the *measure (esc)* key to take a new measurement.

Timed – Measurements are continuously updated on the display and saved to the data log at the pre-set time interval until the measurement mode is exited.

Single-Shot – Press the *measure (esc)* key to start a measurement. When the pre-set wait time is reached, the measurement is locked on the display and saved to the data log. Press the *measure (esc)* key to take a new measurement.

Continuous – Measurements are continuously updated on the display. Press the *log/print* key to save the measurement to the data log.

Press the *channel* key until the desired combination of measurement channels are shown.

- Rinse the electrodes with deionized water or appropriate solution, blot dry with a lint-free tissue and place into the sample.
- 2. Start the measurement and wait for the reading to stabilize or reach the set time.
 - a. When using the Auto-Read or Single-Shot read type, press the *measure (esc)* key to start the measurement and stirrer probe.
 - b. When using the Timed or Continuous read type, measurements will start immediately; press the *stirrer* key to start and stop the stirrer probe.
- 3. Once the measurement is stable or reaches the set time, record all applicable parameters.

- 4. Remove the electrodes from the sample, rinse, dry and place into the next sample.
- 5. Repeat steps 2 through 4 for all samples. When all samples have been measured, store the electrodes as instructed in the electrode user manuals.

Example Conductivity Measurement – Single Channel Display



Customizing the Display

Select the information shown in the measurement mode.

- 1. In the measurement mode, press the *setup* key.
- 2. Press the ► key until <u>Instrument Settings</u> is highlighted and press the *f3 (Select)* key.
- <u>Display</u> will be highlighted in the left column. Press the
 key to also highlight <u>Display View</u> and press the *f3* (*Select*) key.
- Press the ▲ or ▼ key to highlight an item and press the *f3 (Select)* key to check (display) or uncheck (hide).
- 5. Press the *f1 (Done)* key. Press the *measure (esc)* key to return to the measurement mode.