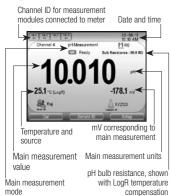
Example pH Measurement – Single Channel Display



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.

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

Water Analysis Instruments

thermoscientific.com/water

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

India

Tel: (91) 22-4157-8800 wai.asia@thermofisher.com

China

Tel: (86) 21-68654588 wai.asia@thermofisher.com

Japan

Tel: (81) 045-453-9175 wai.asia@thermofisher.com

1. In the measurement mode, press the *setup* key.

- Press the ◀ or ▶ key to highlight <u>Log View</u> and press the f3 (Select) key.
- 3. Press the ◀ or ▶ key to highlight Data Log.
- Press the ▲ or ▼ key to highlight <u>View</u> and press the f3 (Select) key.
- Press the ▲ or ▼ 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.

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.
- 2. Press the ◀ or ▶ key to highlight Log View and press the f3 (Select) key.
- 3. Press the ◀ or ▶ key to highlight Cal Log
- 4. Press the ▲ or ▼ key to highlight the desired calibration parameter and press the f3 (Select) key.
- Press the ▲ or ▼ key to view each calibration for the selected parameter.
- 6. Press the *measure (esc)* key to return to the measurement mode.

Australia

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

Thermo scientific

A Thermo Fisher Scientific Brand



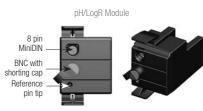
Thermo Scientific Orion

VERSA STAR pH/mV/RmV/ORP and LogR Temperature Module

This literature provides basic instructions on operating the Thermo Scientific™ Orion™ Versa Star™ meter when the Versa Star pH/LogR 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

- Prepare the universal power adapter, install the meter-attached electrode stand and verify that the pH/LogR module is connected to the meter. If it is not connected, insert the module into two available channels 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
- Connect the pH electrode cable to the BNC input on the module and ATC probe cable to the 8 pin MiniDIN input on the module. An ATC probe can be used when calibrating LogR temperature or as the direct temperature source.





- 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 buffers, 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.

pH 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.
- 2. Press the ◀ or ▶ key to highlight the appropriate Channel setup menu and press the f3 (Select) key.
- 3. Press the ◀ or ▶ key to highlight Mode and press the f3 (Select) key.
- 4. Press the ▲ or ▼ key to highlight pH and press the f3 (Select) kev.
- 5. View and update the displayed menu options.

pH Setup Menu

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

Electrode SI No	(no value)
Sample ID	Off, Manual, Auto Increment
Resolution	0.1, 0.01, 0.001
Stability	Smart, Fast, Medium, Slow
Averaging	Off, Automatic Smart
Read Type	AutoRead, Timed, Single-Shot, Continuous
Buffer Group	USA, DIN, Custom

Isopotential Value	Off, On (7.000)
Linear Regression	Off, On
Alarm	Limit (Off), CalDue (Off), Set Point (Off)

Selecting and Calibrating LogR **Temperature**

LogR technology allows most glass-bulb pH electrodes to be used as a temperature source once a LogR temperature calibration is performed with the pH electrode. For the best results, calibrate at two temperature points that bracket the expected sample temperature and are 10 to 20 °C apart. Recalibrate LogR temperature weekly or if the meter loses power.

- 1. In the measurement mode, press the *setup* key.
- 2. Press the ◀ or ▶ key to highlight the appropriate Channel setup menu and press the f3 (Select) key.
- 3. Press the ◀or ▶ key to highlight Temperature and press the f3 (Select) kev.
- 4. Press the ▲ or ▼ key to highlight Temperature Input and press the ▲ or ▼ key to highlight the channel with pH/LoaR module (i.e. Ch2-LoaR).
- 5. Press the ▲ or ▼ key to highlight Temperature Cal and press the f3 (Select) key.
- 6. Place the pH electrode and NIST-traceable thermometers or ATC probe into a calibration solution (pH buffer or tap water) with a known, stable temperature. Moderately stir the solution using the Orion Star stirrer probe or stir plate with stir bar.
- 7. Press the f3 (Start) key and wait for the temperature value on the meter to stabilize.
 - a. The stirrer probe will start stirring and continue until the reading stabilizes.
- 8. Once the temperature value is stable, 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, use the numeric keypad to enter the temperature value and press the f2 (Accept) key to confirm the entered value.
- 9. Press the f2 (Next) key to proceed to the next calibration point and repeat steps 6 through 8 or press the f3 (Cal Done) key to save and end the calibration.

pH Calibration

- 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.
- 2. Rinse the pH electrode and any other electrodes in use with distilled water, blot dry with a lint-free tissue and place into the pH buffer.
- 3. When the electrode and buffer are ready, press the f3 (Start) key.
 - a. The stirrer probe will start stirring and continue until the reading stabilizes.
- 4. Wait for the pH 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 and use the numeric keypad to enter the pH value. Press the f2 (Accept) key to confirm the entered value.
- 5. Press the f2 (Next) key to proceed to the next buffer and repeat steps 2 through 4 or press the f3 (Cal Done) key to save and end the calibration.
 - a. If a one point calibration is performed, press the f3 (Slope Edit) key, press the f3 (Clear) key, use the numeric keypad to enter the new slope value and press the f2 (Accept) key.
- 6. The meter will display the calibration summary including the slope 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.

- 1. Rinse the electrodes with distilled 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.