



# Thermo Scientific Orion

## Versa Star Pro Meter and USB Printer Interfacing

### General Information

This instruction sheet provides general guidance on interfacing the Thermo Scientific™ Orion™ Versa Star Pro™ meter with a USB printer. Orion Versa Star Pro meters offer expanded USB printer compatibility with new PCL 6 driver hardware for better compatibility with recent, commercially available USB printers. Although we strive to ensure the compatibility of our meters with USB printers, we cannot ensure compatibility with all models.

#### Tested Printer Models:


- Fuji Xerox DocuPrint P225D
- HP LaserJet Pro MFP M225DW

#### Recommended Printer Models:

- Brother HL-L2365DW
- Brother MFC-L2740DW
- Fuji Xerox DocuPrint P255D
- Fuji Xerox DocuPrint P265D
- HP LaserJet M252DW
- HP LaserJet Pro MFP M426FDW

For comprehensive information on meter and module system setup, operation and advanced features, refer to the Orion Versa Star Pro meter user manual available on the Orion Versa Star Pro literature CD and [www.thermofisher.com/water](http://www.thermofisher.com/water).

## Meter to Printer Interface Requirements

| Component     | Requirement  |
|---------------|--|
| Meter         | Thermo Scientific Orion Versa Star Pro meter, serial number V10000 or higher   |
| Printer       | PCL6 driver compatible   |
| Printer Paper | Letter size format   |
| Cable         | Micro AB USB Male (meter connection) to printer-specific USB connection (may vary by printer model)  |
|               | For Tested Printer Models, printer cable requirement is Micro AB USB Male to Standard USB B Type cable:<br> |

**Note:** The printer cable must be purchased separately and the printer connection type may vary by printer model.

## Meter Setup

1. Power on the Orion Versa Star Pro meter.
2. In the measurement mode, press the **setup** key to access the Main Setup menu.
3. Press the **▲**, **▼**, **◀** or **▶** key to highlight the Instrument Settings icon and press the **f3 (Select)** key.
4. Press the **▲** or **▼** key to highlight Communication and press the **◀** or **▶** key to highlight USB PRINTER.
5. Press the **f2 (Page 2)** key, press the **▲** or **▼** key to highlight Printing and press the **◀** or **▶** key to highlight On.
6. Press the **▲** or **▼** key to highlight Print Format and press the **◀** or **▶** key to highlight Printer.
7. Press the **measure (esc)** key to return to the measurement mode.
8. The USB symbol should now be displayed on the top of the screen:

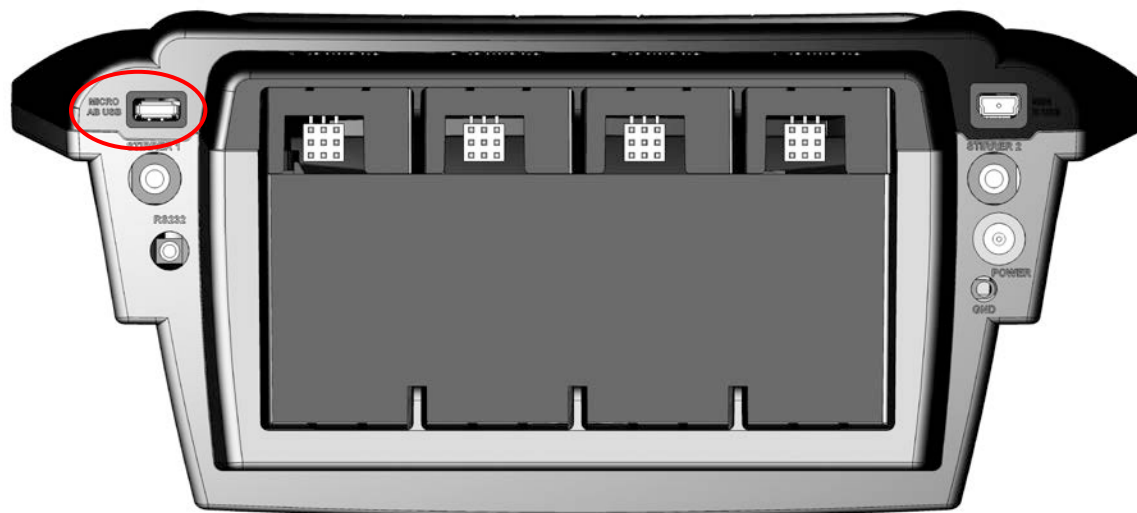


## Printer Setup

1. Unpack the printer and prepare according to the instructions in the printer manual. Add letter size paper to the paper tray.
2. Power on the printer and confirm the printer status is ready.

# Meter to Printer Interfacing

1. Connect the USB cable to the MICRO AB USB connection on the meter and the appropriate USB connection on the printer.



2. Press the **log/print** key on the meter keypad to send the displayed measurement information to the printer. The Printer symbol should be displayed on the top of the screen.



3. The printer should print the measurement information, similar to the following example:

```
-----  
Thermo Scientific (c) 2016  
Orion Versa Star Pro  
Meter S/N          V15454          SW Rev          4.0  
  
12-05-16, 1:26:42 PM  
Channel - 1  
Probe S/n          -----  
User ID            -----  
Mode               pH  
pH                 7.000 pH  
Temperature        25.0 C (MAN)  
Method             100  
CalLog #           0  
Sample ID          -----  
Module S/n #      VA11284  
mV                 0.0 mV  
Average Slope     100.0 t  
Last Cal           -----  
Channel - 2  
Probe S/n          -----  
User ID            -----  
Mode               Conductivity  
Conductivity       0.000 uS/cm  
Cell Constant      0.4750 /cm  
Temperature        25.0 C (MAN)  
Temp. Coefficient  2.10 %/C  
Method             200  
CalLog #           0  
Sample ID          -----  
Module S/n #      VA10941  
Conductance       0.000 uS  
Temp. Reference   25.0 C  
Last Cal           -----  
Operator           _____  
Signature          _____
```

## Ordering Information

| Cat. No.  | Description  |
|-----------|--|
| VSTAR10   | Versa Star Pro pH meter with stand   |
| VSTAR12   | Versa Star Pro pH meter standard kit   |
| VSTAR13   | Versa Star Pro pH meter difficult sample kit                                       |
| VSTAR20   | Versa Star Pro conductivity meter with stand                                       |
| VSTAR22   | Versa Star Pro conductivity meter standard kit                                     |
| VSTAR23   | Versa Star Pro conductivity meter pure water kit                                   |
| VSTAR30   | Versa Star Pro RDO/DO meter with stand   |
| VSTAR32   | Versa Star Pro RDO/DO meter kit  |
| VSTAR40A  | Versa Star Pro pH/ISE meter with stand   |
| VSTAR40A2 | Versa Star Pro pH/ISE meter kit  |
| VSTAR40B  | Versa Star Pro dual channel pH/ISE meter with stand                                |
| VSTAR40B2 | Versa Star Pro dual channel pH/ISE meter kit                                       |
| VSTAR50   | Versa Star Pro pH and conductivity meter with stand                                |
| VSTAR52   | Versa Star Pro pH and conductivity meter kit                                       |
| VSTAR80   | Versa Star Pro pH/LogR meter with stand  |
| VSTAR82   | Versa Star Pro pH/LogR meter difficult sample kit                                  |
| VSTAR83   | Versa Star Pro pH/LogR meter micro sample kit                                      |
| VSTAR90   | Versa Star Pro pH/ISE, conductivity and RDO/DO meter with stand                    |
| VSTAR91   | Versa Star Pro dual channel pH/ISE, conductivity and RDO/DO meter with stand       |
| VSTAR92   | Versa Star Pro pH/ISE, conductivity and RDO/DO meter kit                           |
| VSTAR93   | Versa Star Pro dual channel pH/ISE, conductivity and RDO/DO meter kit              |
| VSTAR94   | Versa Star Pro pH/ISE, conductivity and RDO/DO meter kit with RDO optical DO probe |

## Technical Support

For any questions or if you require assistance, contact our Technical Support Specialists:

- Email [WLP.techsupport@thermofisher.com](mailto:WLP.techsupport@thermofisher.com)
- Within the United States, call **1-800-225-1480**
- Outside the United States, call **+1-978-232-6000**

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 below.

[thermofisher.com/water](http://thermofisher.com/water)

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Water and Lab Products

**North America**  
Toll Free: 1-800-225-1480  
Tel: 1-978-232-6000  
[info.water@thermofisher.com](mailto:info.water@thermofisher.com)

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

**India**  
Tel: (91) 22-4157-8800  
[wai.asia@thermofisher.com](mailto:wai.asia@thermofisher.com)

**Japan**  
Tel: (81) 045-453-9175  
[wai.asia@thermofisher.com](mailto:wai.asia@thermofisher.com)

**China**  
Tel: (86) 21-68654588  
[wai.asia@thermofisher.com](mailto:wai.asia@thermofisher.com)

**Singapore**  
Tel: (65) 6778-6876  
[wai.asia@thermofisher.com](mailto:wai.asia@thermofisher.com)

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

IS-VSPUSBPRT-E RevA 0516

**Thermo**  
SCIENTIFIC  
A Thermo Fisher Scientific Brand