

MicroPure Consumables Replacement Information

Consumable	Part #	Estimated Life Time *	How You Will Know To Change
Ultrapure Cartridge	09.1006	Up to 12 months depending on feed water quality	<ul style="list-style-type: none"> Purity drops below acceptable purity level "Lim. val. pure w." displayed Change cartridge message displayed – cartridge exceeded maximum time in the system
0.2 µm Final Filter	09.1003	Up to 12 months	<ul style="list-style-type: none"> Dispensing water flow is reduced 09.1006 ultrapure cartridge is replaced
Ultrafilter	50133981	When applicable – up to 2 years	<ul style="list-style-type: none"> Low flow Replace at least every 2 years
UV Lamp	09.1002	When applicable – up to 2 years	<ul style="list-style-type: none"> "UV duration" alert displayed Replace at least every 2 years
Vent Filter – Micropure ST Units Only	22.0091	When applicable – up to 12 months	<ul style="list-style-type: none"> Replace at least once a year
Cleaning Solution	CMX25	n/a	<ul style="list-style-type: none"> Bacteria present in water When 09.1006 ultrapure cartridge is replaced Requires reusable cleaning cartridge

Note: The actual life time of the consumables depend on the quality of the feed water and the amount of water used by the system.

MicroPure System



Quick Reference Guide

Thermo Scientific

Barnstead MicroPure/MicroPure ST
Ultrapure Water System Quick Reference Guide



Please refer to operation manual for complete instructions

thermoscientific.com/purewater

© 2016 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. Specifications, terms and pricing are subject to change. Not all products are available in all countries. Please consult your local sales representative for details.



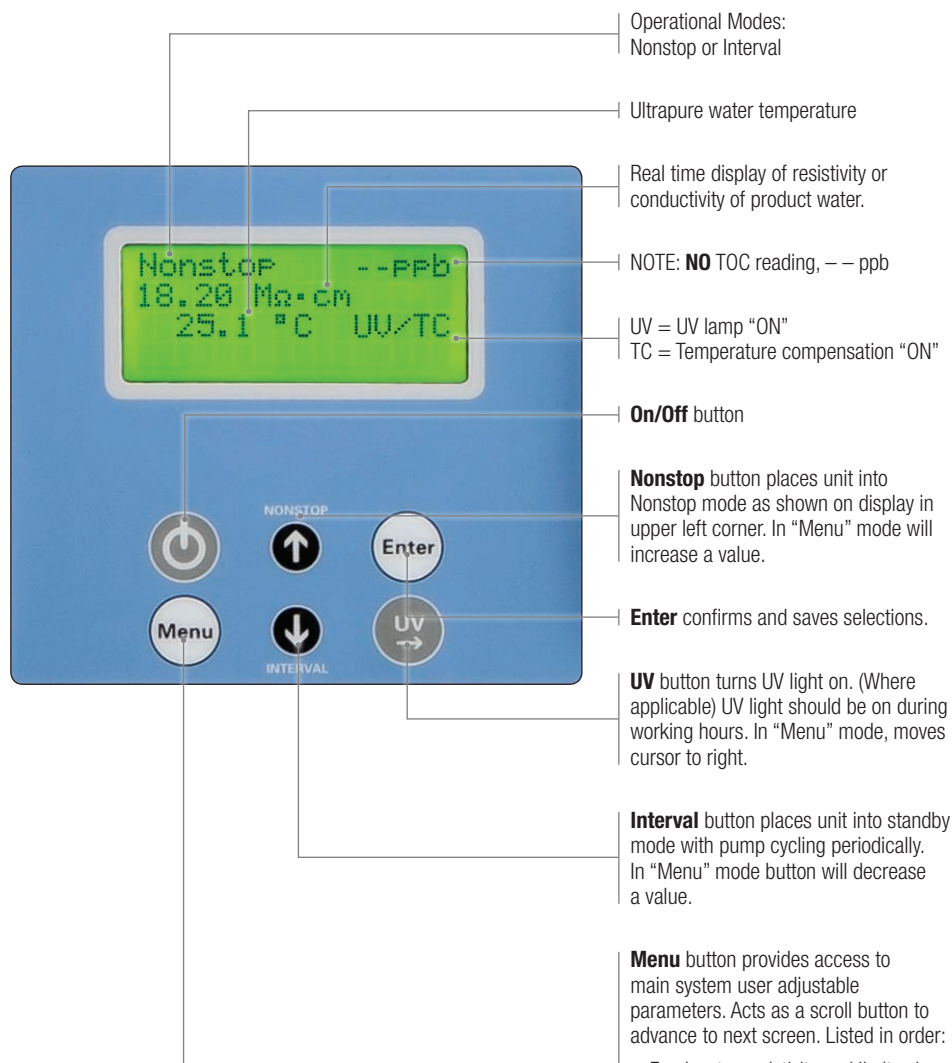
Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand

BRWPQGMICROPURE 0416

Thermo
SCIENTIFIC

Display Functionality Keys and Display Readout



Note:

MicroPure ST units only. Manually add pretreated water to storage tank in back of system. Avoid draining the 5.5 L tank completely to prevent introducing air to the system.

Operation

- **Nonstop:** Pump runs continuously. Use this mode during working hours to ensure highest purity water. To put system in Nonstop mode from Interval mode, press the Nonstop button.
- **Dispensing** in Nonstop mode only. Dispense by turning flow control knob counter clockwise to open to the desired flow.
- **Interval:** Standby mode. Unit will automatically switch into interval after 2 hours of non-use.
 - In interval mode, pump cycles for 5 min (adjustable) out of every 30 min.
 - **Nights and Weekends:** Keep system power on and allow it to automatically switch to interval mode.
 - Displayed purity may change when system is not recirculating.
 - Manually put system in Interval mode by pressing On/Off button twice.
- **Rinse:** Flush of unit – manually activated for 30 seconds (adjustable) in user menu. Rinse system if unit has been off or unused for a long weekend or holiday.
 - Activate rinse by pressing menu button until "Rinse" is displayed and press "Enter".
 - **Automatic Rinses:** Occur when system switches from Interval to Nonstop and when pump recirculates during Interval mode.

Maintenance

Refer to operation manual for complete maintenance instructions.

- Regular maintenance is important for consistent purity.
- Change ultrapure cartridge and 0.2 μm final filter when purity has dropped to limit value or a minimum of once a year.
- Disinfect system when changing the ultrapure cartridge. Use reusable disinfection cartridge, 09.1102 with the cleaning solution.
- Change UV light and/or ultrafilter on a regular set schedule.

Common Alerts

Check operation manual for full list of alerts.

- **Lim. Val. Pure W.:** Displayed when purity has reached ultrapure water limit value. Replace ultrapure cartridge.
- **Limit Value Feed:** Displayed when feed water purity limit is out of range. Check feed water supply.
- **Change Cartridge:** Cartridge exceeded maximum time in the system. Disinfect and replace ultrapure cartridge.
- **UV Duration:** UV light has exceeded maximum time in the system. Replace UV light.

