Smart2Pure Consumables Replacement Information

External Pretreatment

Pretreatment is important for the life of the RO membrane. One or more of the following pretreatment cartridges may be used depending on feed water quality. Check the boxes below for the external pretreatment installed before this system.

Note: Frequency of replacement may vary based on feed water supply.
□ 06.5452 Hardness stabilizer: Replace 6-12 months

□ 06.5201 5 µm filter with carbon cartridge 3-4 months

□ 06.5101 10" 1 micron prefilter 4-6 months

□ 06.5204 10" 5 μm filter with hardness stabilizer cartridge 6-12 months

	Consumable	Part #	Estimated Life Time *	How You Will Know To Change
System Consumables	Smart2Pure 3 LPH Pretreatment Cartridge (RO Membrane)	09.2003	Up to 12 months	Ultrapure cartridge exhausting quickly
	Smart2Pure 6 LPH Pretreatment Cartridge (RO Membrane)	09.2006	Up to 12 months	Ultrapure cartridge exhausting quickly
	Ultrapure Cartridge	09.1020	Up to 12 months	Purity drops below acceptable limits "Limit value Cond." alarm and LED on
	0.2 µm Filter For System And Hand Remote Dispenser	09.1003	Up to 12 months	 Dispensing water flow is reduced Ultrapure cartridge is replaced Bacteria is present in product water
	Ultrafilter	50133981	When applicable – up to 2 years	Dispensing water flow is reducedReplace every 2 years
	UV Lamp For Type I System	09.1002	When applicable – up to 2 years	UV lamp is not lit Replace every 2 years
Reservoir Consumable	Reservoir Vent Filter	22.0091	Up to 12 months	Replace once a year
Cleaning	Cleaning Solution	CMX25	n/a	Ultrapure cartridge is replaced Minimum once a year Bacteria is present in system

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.

Smart2Pure System



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

Barnstead Smart2Pure 3LPH & 6LPH System Quick Reference Guide





Display Functionality Keys and Display Readout



Note:

Avoid draining the tank completely to prevent

introducing air to the system.

Displays real time conductivity/resistivity of product water or temperature.

Display Line: LED illuminates next to the current measurement mode in μ S/cm, Mohm-cm and C.

Limit Value Line: LED illuminates next to the error condition when a fault has occurred.

- Cond.: Conductivity level is out of range.
- Temp.: Temperature of pure water exceeded limit.
- Status Sensor: Error of conductivity and temperature measuring cell.

Up/Quit button increases adjustable values, silences acoustic alarms.

Enter button confirms changes, adjusts the cursor position.

Down button decreases adjustable values.

Menu button: Access to system display or limit value parameters. Press **Menu** to go to setting, ↑ ↓ to change setting. Press **Enter** to save changes.

Note: Parameters scroll up or down, not in continuous loop.

- Press Menu 1X: Set display purity units and temperature, recommended setting: MΩxcm or μS/cm.
- Press Menu 2X: Set purity limit value in conductivity or OFF.
- Press Menu 3X: Set temperature limit value, set to "OFF" in units with UV.

Operation

Start Up

 On/Off Power: To power up/down the system, simply unplug power supply.

Daily Operation

- When Water Is Being Dispensed: System is in run mode.
 Water is recirculated through system and tank continuously for optimal water quality.
- When Water Is NOT Being Dispensed After 20 Minutes:
 System automatically switches to standby mode, system and tank re-circulates ~10 minutes every 30 minutes.
- Nights and Weekends: Keep system power on and allow it to automatically switch to standby mode.

Key Components

- Boost Pump: 3 minute delay upon start up. Turns on when storage tank is low. Water will flow from concentrate line when tank is filling.
- External Pretreatment: Protects reverse membrane (RO)/ pretreatment cartridge.
- Units With UV Light: UV light is on during recirculation mode, water temperature can increase with recirculation. Suggested setting for temperature limit = OFF.

Dispensing Options

- Type I Water: Dispense Type 1 water by manually turning the flow control knob counter-clockwise to open to the desired flow.
 Note: Do not over tighten dispense knob to close.
- Water can be dispensed from right side of system using supplied valve or optional remote hand dispenser.

Maintenance

- Water in storage tank should be turned over every 1-2 days.
- Replace external prefilters, pretreatment cartridge (RO membrane), UV light, UF, and tank vent filter on regular set schedule.
- Clean system before replacing ultrapure cartridge and final filter when conductivity limit value is reached or minimum of once a year.
- When cleaning Smart2Pure, use disinfection adapter supplied with unit.
- When pretreatment cartridge (RO membrane) is replaced, discard first tank of water.

Common Alerts

- Audible alarm and red LED when limit value is reached for purity, temperature or sensor.
- Press Up/Quit button to silence alarm.
- Purity Alarm: Change ultrapure cartridge.
- Temperature Alarm: Drain water from system to reduce water temperature. If alarm continues, contact service.
- Sensor Error: Contact service.

