

GenPure xCAD Plus Consumables Replacement Information

Consumable	Part #	Estimated Life Time *	How You Will Know To Change
Ultrapure Cartridge	09.2005	Up to 12 months	<ul style="list-style-type: none"> Purity drops below acceptable purity level "Lim. val. pure w." displayed "Change cartridge" message displayed – cartridge usage exceeded maximum time in system
0.2 µm Final Filter	09.1003	Up to 12 months	<ul style="list-style-type: none"> Dispensing water flow is reduced 09.2005 ultrapure cartridge is replaced
Ultrafilter	50133980	When applicable – up to 2 years	<ul style="list-style-type: none"> Dispensing water flow is reduced Replace at least every 2 years
UV Lamp	09.2002	When applicable – up to 2 years	<ul style="list-style-type: none"> "UV Intensity" alert (units with TOC only) "UV Duration" alert is displayed
Cleaning Solution	CMX25	n/a	<ul style="list-style-type: none"> Bacteria present in water When 09.2005 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.

GenPure xCAD Plus 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.



Quick Reference Guide

Thermo Scientific

Barnstead GenPure xCAD Plus System
Quick Reference Guide



Please refer to operation manual for complete instructions



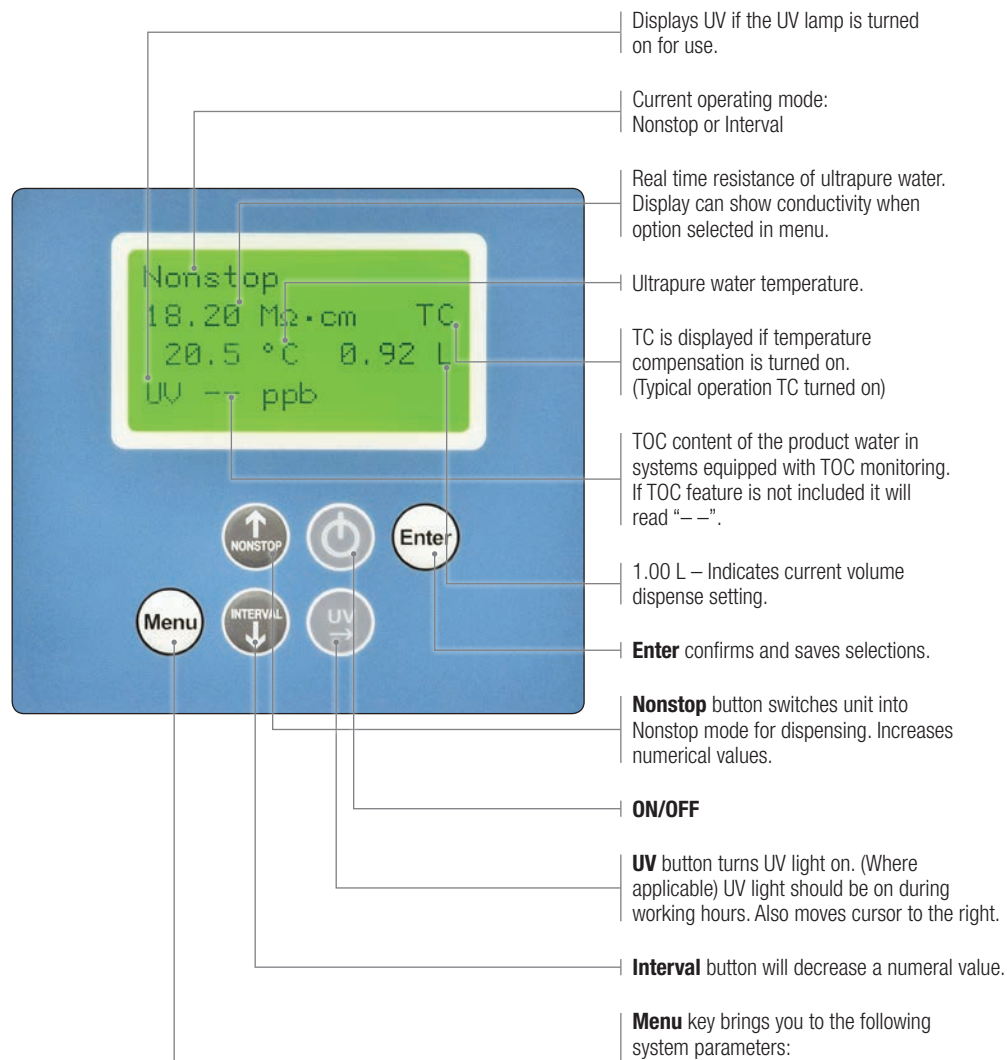
Thermo
SCIENTIFIC

A Thermo Fisher Scientific Brand

BRWPQGGENXCAD 0416

Thermo
SCIENTIFIC

Display Functionality Keys and Display Readout



Maintenance

- Regular maintenance is important for consistent purity. Change ultrapure cartridge and final filter when purity has dropped to limit value or a minimum of once a year.
- Clean system before changing the ultrapure cartridge using reusable, 09.2201 cleaning cartridge.
- Change UV light and/or ultrafilter on a regular set schedule.

Operation

- **Nonstop** mode during working hours for highest purity water. Pump and UV runs continuously.
- **Volumetric Dispense:** In Nonstop mode only. Press Enter or Dispense button, display will show last volume dispensed, press button again to dispense water. While dispensing, pressing Enter or Dispense button will stop water flow.
 - **Change Volume:** Press Enter, then ↓ or ↑ to change volume, UV button to move over to next digit.
- **Manual Dispense:** Press dispense button 2X to start dispensing. Press again to stop. (Change volume on display as necessary to dispense larger volumes of water).
- **Interval:** Standby mode. Unit enters interval upon start up and after period of non-use.
 - In interval mode, pump recirculates water 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:** Manually activate rinse 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. System will rinse for 30 seconds.
 - Interval Rinse Time: System will activate a 0.5 second rinse when system switches from Interval to Nonstop mode and when system initiates recirculation in Interval mode.

Common Alerts

Check operation manual for full list of alerts.

- Screen may turn red indicating an alert condition.
- **Lim. Val. Pure W.:** Displayed when purity has reached ultrapure water limit value. Replace ultrapure polishing cartridge.
- **Limit Value Feed:** Displayed when feed water purity limit is out of range. Check feed water quality.
- **Change Cartridge:** Cartridge exceeds maximum time in the system. Disinfect system and replace ultrapure cartridge. Enter serial number for new cartridge using user code.
- **UV Duration:** UV light has exceeded maximum time in the system. Replace UV lamp.
- **UV Intensity:** Displayed in TOC units only, red screen alert if below 50 % intensity. Replace UV lamp.

