

Thermo Scientific™ Aquanex™ Touch Smart Dispenser

1. Tank level
2. Dispenser ID
3. Dispense volume setting—touch screen
4. System health status indicator
5. Type 1 water temperature
6. Dispensing water purity
7. Dispense and flow speed selector buttons

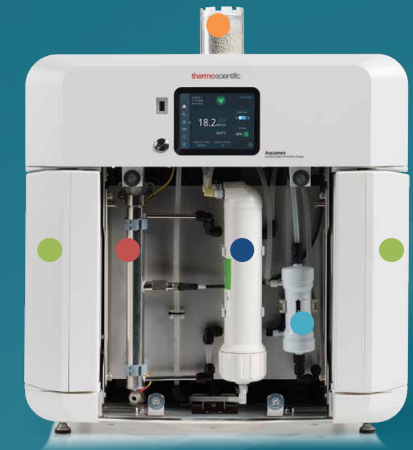


Aquanex system user interface



1. System health/status indicator
2. Unit name, time and date
3. Home screen navigation bar
4. Dispenser volume setting
5. Dispensing water purity
6. Type 1 water temperature
7. Tank level
8. Water type selector
9. Secure mode log in/out

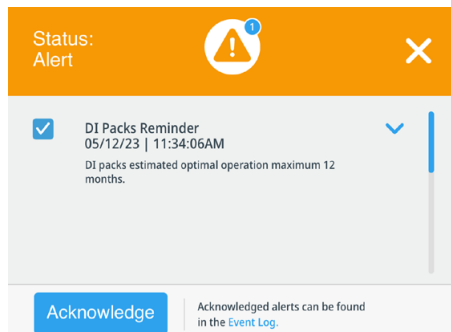
| System consumables | | |
|----------------------------------------------------------------------|---------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Name/catalog number | Estimated lifetime* | When to replace |
| Thermo Scientific™ Reverse Osmosis Cartridge 50157608 | Up to 12 months | <ul style="list-style-type: none"> • “Change RO Assembly” alert on display • DI cartridges exhausting quickly |
| Thermo Scientific™ Aquanex™ DI Cartridge Set 7501050 | Up to 12 months | <ul style="list-style-type: none"> • “Change DI Packs” alert on display • “Water Quality Alert” on display |
| Thermo Scientific™ Barnstead™ Systems Ultrafilters 50133980 | Up to 24 months | <ul style="list-style-type: none"> • “Change Ultrafilter” alert on display • Low flow from dispenser • Pyrogens in product water that don’t improve with disinfection |
| Thermo Scientific™ Barnstead™ Replacement UV Lamps 09.2002 | Up to 24 months | <ul style="list-style-type: none"> • “Change UV” alert on display • Higher TOC levels observed • UV lamp is burned out |
| Thermo Scientific™ CO ₂ Adsorber + Sterile Filter 06.5002 | Up to 12 months | <ul style="list-style-type: none"> • “Change CO₂ Vent Filter” alert on display • Color indicator changed from white to purple |



Dispenser consumables

| Name/catalog number | Estimated lifetime* | When to replace |
|-------------------------------------------------------------------------|---------------------|------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Thermo Scientific™ Barnstead™ Point of Dispense 0.1 μm Filters 50157375 | Up to 6 months | <ul style="list-style-type: none"> • **“Change POU Final Filter” alert on display • Low flow from dispenser • When DI cartridges are replaced • When bacterial growth is suspected in the POU filter |

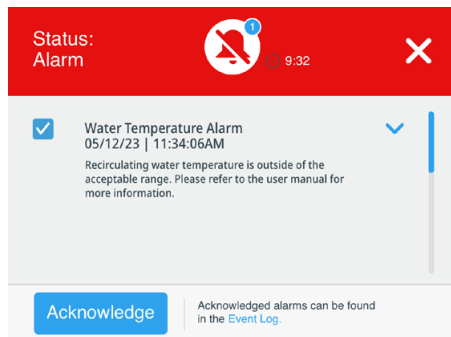
*The actual lifetime of the consumable depends on the quality of the feed water and the amount of water used by the system.



Alerts

A yellow banner message will indicate an alert, but normal operation will not be interrupted. Tap the yellow alert triangle for more information and to acknowledge and clear the alert. Examples include:

- Type 1 water quality
- Feed water pressure low
- Consumable needs replacement



Alarms

A red banner message will indicate an alarm, and the system will automatically enter Safe Mode until acknowledged. Tap the red bell for more information. Examples include:

- Water temperature
- Leak detection
- DI pressure
- RO pressure

Operation

- System will automatically refill tank when level drops below 90%
- System goes into deep sleep after two hours; tap main screen or dispenser screen to wake up
- To change volumetric dispense from Smart Dispenser, press and hold dispenser volume setting
- To change volumetric dispense from main display, tap dispenser volume setting and select desired volume
- When operating in manual dispense mode (∞), use the arrow keys to increase or decrease the flow speed
- Start or stop water flow by pressing the teardrop button on the Aquanex Touch Dispenser
- To dispense Type 2 water, set toggle on main screen to II and use optional Type II hand dispenser

Maintenance

- Consumable status and replacement processes can be accessed within the maintenance menu on the Aquanex home menu
- Detailed instructions on replacement of consumables can be found in the product manual
- Disinfect the Aquanex unit on a regular basis and prior to changing DI cartridges; detailed disinfection instructions are provided in the product manual

Need assistance?

Contact Customer Support

Phone: 866-984-3766

Find out more at thermofisher.com/aquanex

For Laboratory Use. It is the customer's responsibility to ensure that the performance of the product is suitable for customers' specific uses or applications.

© 2019–2023 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

Document 274006-001 REVA EXT6099

Quick reference guide

Thermo Scientific™ Aquanex™ Ultrapure Water Purification System

