

1500 Series A2 Biological Safety Cabinets

Quick-start guide

Initial start-up

Plug the Thermo Scientific™ 1500 Series A2 Class II Biological Safety Cabinet (BSC) into the power outlet, making sure that the supply voltage matches the data plate. When the BSC is powered up, the start-up wizard will provide guidance.

Before using the cabinet

The BSC must be installation tested in accordance with NSF49/ YY0569.

Operating modes

Working mode

Window at operating height, fans energized and airflows stable. Personal and product protections provided.

Standby mode

Window closed, fans energized and sample chamber environment maintained.

Disinfection mode

Window closed, fans and UV lamps energized. Sample chamber disinfected on timed cycle.

Access control

If desired, the administrator can access factory setting and field certification mode following the procedure in the service manual.

Health status



OK



Alarm



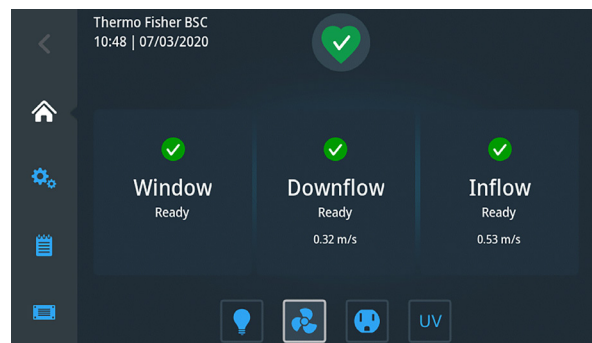
Alert



Snoozed Alarm

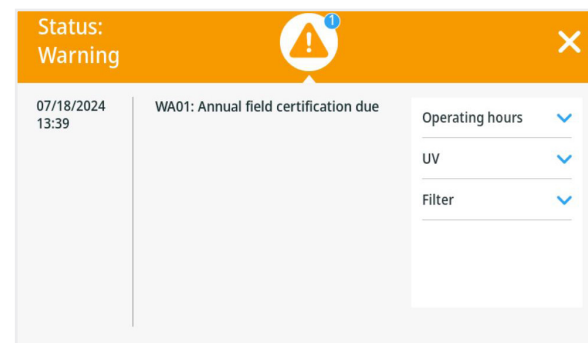
Alarm/alert log

Any alarm or alert that occurs will be recorded into the alarm/alert log. In the log, the maximum number of items is 50. The log can only be displayed on the touchscreen, but cannot be downloaded to a USB drive.



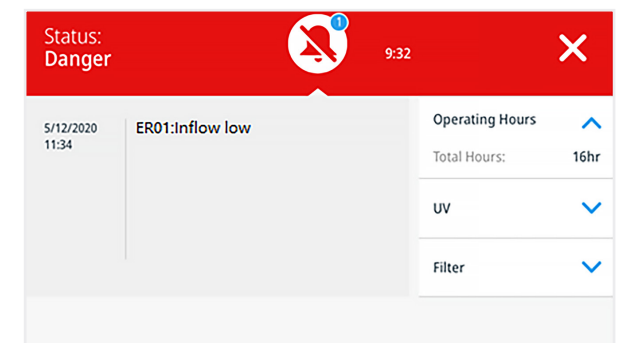
Home screen

- Health status information
- Alarms and alerts
- Settings screen
- Alarm/alert log
- Data plate



Alert screen

- Active past alerts are listed. Tapping the down arrow provides additional information.
- Additional status information is available on the right section of the screen. Tapping the down arrow provides status information by category.



Alarm screen

- There are two main alarms; airflow and window position.
- Active past alarms are listed.
- Tapping the down arrow provides additional information.
- Active alarms will disappear after the condition is cleared.

For technical specifications and safety instructions, including warnings and conditions for safe use, please refer to the user manuals below.
 Made in PRC 1500 A2 BSC Manual: thermofisher.com/1500a2-prc-manual | Made in USA 1500 A2 BSC Manual: thermofisher.com/1500a2-usa-en-manual

Learn more at thermofisher.com/bsc