

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 BSC is powered up, the start up wizard will provide guidance.

Before using 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 procedure in service manual.

Health status



OK



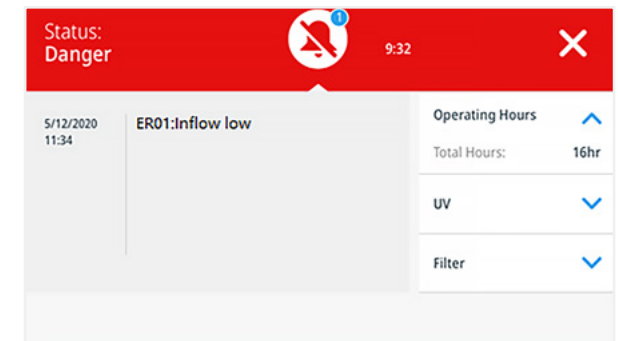
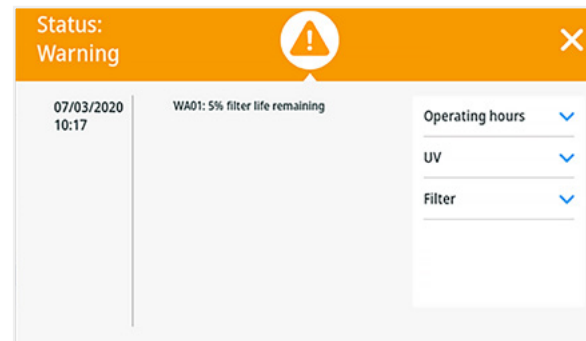
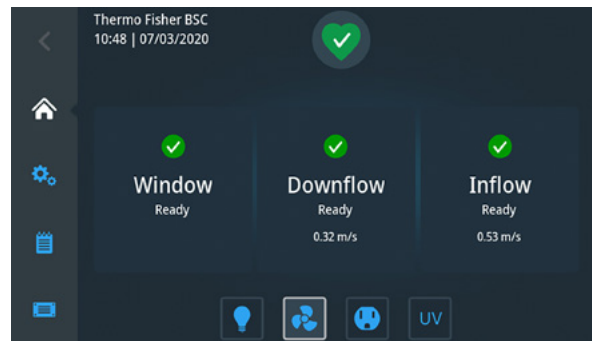
Alarm



Alert



Snoozed Alarm



Home screen

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

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 right selection 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, safety instructions including warnings and conditions for a safe usage, please refer to the user manual:
<https://assets.thermofisher.com/TFS-Assets/LED/manuals/CIC0005677%201500%20BSC%20Op%20Manual%20NSF%20EN%20Rev.%2000.pdf>

Learn more at thermofisher.com/bsc