Thermo Fisher

1300 Series A2 Biological Safety Cabinets

Quick start guide

Initial start up

Plug the Thermo Scientific™ 1300 Series Class II, Type A2 Biological Safety Cabinet (BSC) into the power outlet, making sure that the supply voltage matches the data plate.

Before using cabinet

The BSC must be installation tested in accordance with NSF49.

Operating modes

Working mode

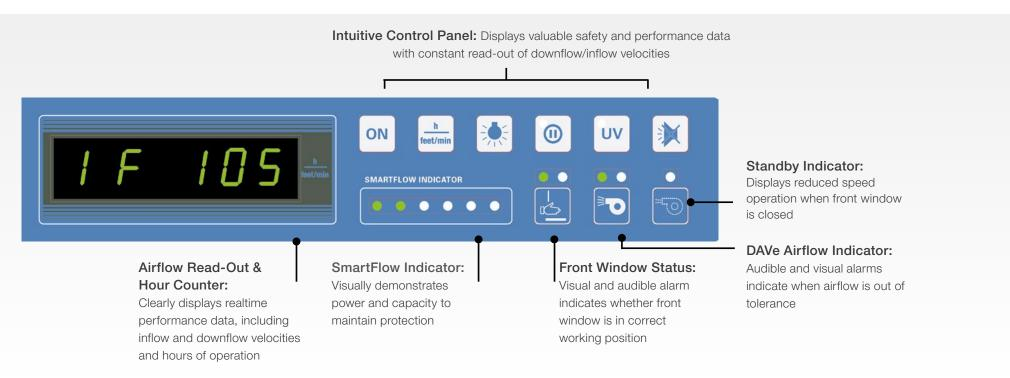
Window at operating height, fans energized and airflows stable. Personal and product protections provided.

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.



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/1300-A2-BSC-User-Manual.pdf



Learn more at thermofisher.com/1300A2

