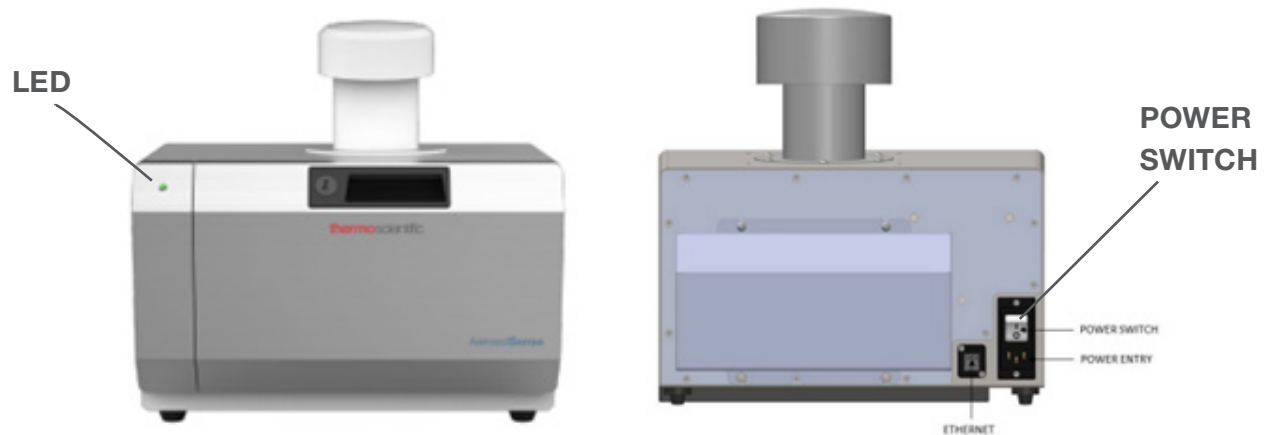


AerosolSense Sampler reset after power down

Procedure for resolving LED and power switch failure

Introduction

The recommended guidance for sample collection with the Thermo Scientific™ AerosolSense™ Sampler 2900 is between 2 and 12 hours and at maximum 24 hours of continuous sample collection. If the sampler is run continuously to sample longer than 24 hours, there is a possibility that the AerosolSense Sampler may power down during sampling and the LED in the top left of the AerosolSense Sampler front panel will not be illuminated despite no power loss to the sampler or toggling of the main power switch. If your AerosolSense Sampler experiences this condition, follow the steps below for resetting your sampler.



Resetting AerosolSense Sampler after power down condition

1. Verify that the main power switch on the back panel of the sampler is in the “ON” position and the power cord is plugged into the sampler.
2. Turn off sampler by toggling main power switch to the “OFF” position.
3. Follow the “Basic Operation” section of the “AerosolSense 2900 Sampler Manual” for removing the sample cartridge.
4. Once removed, the AerosolSense sample cartridge should be placed in a biohazard bag and sent to a lab for testing.

5. After the sampler has been powered off by the main power switch for 5 minutes, toggle the main power switch to “ON”. The sampler is reset and ready for sampling again.
6. Follow the “Basic Operation” section of the “AerosolSense 2900 Sampler Manual” for installing a new cartridge and to start sampling.
7. After performing these steps, if the sampler powers down within a maximum continuous sample time of 24 hours please contact our Technical Support team for further troubleshooting assistance.

Contact Information

Customer Service

1-508-553-1110 and 800-209-8763

orders.apd@thermofisher.com

Technical Support

support.apd@thermofisher.com

USA

27 Forge Parkway
Franklin, MA 02038
support.apd@thermofisher.com
(800) 209-8763

Find out more at [thermofisher.com/aerosolsense](https://www.thermofisher.com/aerosolsense)

ThermoFisher
S C I E N T I F I C