



# Back to the lab

## Laboratory equipment checklist

Here are a few practical reminders to keep in mind for maintaining your equipment

---

Check the qualified state of your equipment- schedule if overdue

---

Schedule your next preventative maintenance, calibrations, or certifications

---

Consult equipment manuals for start-up recommendations- available on [www.thermofisher.com](http://www.thermofisher.com)

---

Check stock and order replacement supplies for equipment like tubing, bottles, tubes, filters and reagents

---

Check expiration dates on reagents for replacement

---

Label reagent/sample vials (including open date, expiration date)

---

Replace bench top protectors and contamination control mats

---

Calibrate pipettors or send them to a service

---

Wipe down the handles and exterior of the equipment

---

Remember to clean corners of the lab and top of and under equipment where dust can collect

---

Activate the sterilization cycles on your CO<sub>2</sub> Incubators before start-up per manual

---

Order cleaning supplies for the lab- 70% ethanol and quaternary ammonium disinfectants

---

Check gas supply levels to your equipment and schedule replacement

---

Clean the interior the equipment - BSCs, incubators, shakers, centrifuges, and baths - with recommended disinfectant (remember no bleach)

---

Wipe bench top surfaces, pipettors, pipet tip boxes with ethanol/DNase/RNase

---

Autoclave water for your sterile water needs like baths and incubators

Clean gaskets by removing ice, frost, or condensation

Inspect gaskets for damage, replace if needed

Eliminate cardboard storage around refrigerators and freezers because can get wet and breed fungi.

Test the operation on your tabletop equipment prior to use

Order recommended replacements parts and consumables for annual maintenance.

Replace HEPA filters in your CO<sub>2</sub> incubators

Drain and refill your pH and ion selective electrodes. Change the storage solution

Clean or replace the front grill filter on your ULT

If capable, download data and verify last two months of ULT performance

Check LN<sub>2</sub> levels in all manual dewars per manual or auto-fill dewars

Verify tight connections (water, electrical, and CO<sub>2</sub>/LN<sub>2</sub> back-up) to your equipment

Clean down your SpeedVac system, change out the CryoCool from the cold trap and change the oil from your oil pump

Clean condenser fins with compressed air in heated/refrigerated bathes per manual

Inspect your centrifuge rotors for any damage or pitting

Record the maintenance of your equipment in your equipment log

Validate the calibration of your setting with secondary probes or monitors



For more information, visit  
[thermofisher.com/backtothelab](https://www.thermofisher.com/backtothelab)

**ThermoFisher**  
SCIENTIFIC