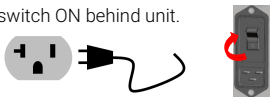


Start Up (Manual Section 1.3.4)

Installation: Do not install or use the pump in an environment exposed to atmospheric agents (rain, snow, ice), dust, aggressive gases, or in explosive environments or those with a high fire risk. During operation, the following environmental conditions must be respected:

- Temperature: from +12 °C to + 45 °C
- Relative humidity: up to 90% (non-condensing)

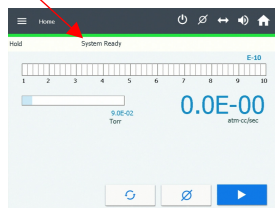
Plug in power cord, flip switch ON behind unit.



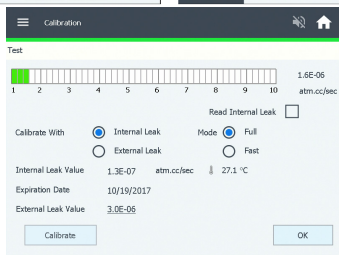
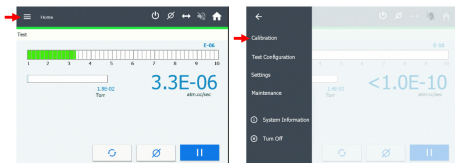
Wait about 7 minutes for "SYSTEM READY."



First time startup will prompt a wizard to help select language, units, date, time, and options. These may be changed any time through the menu screens.



Calibration Set Up (Manual Section 4.1)



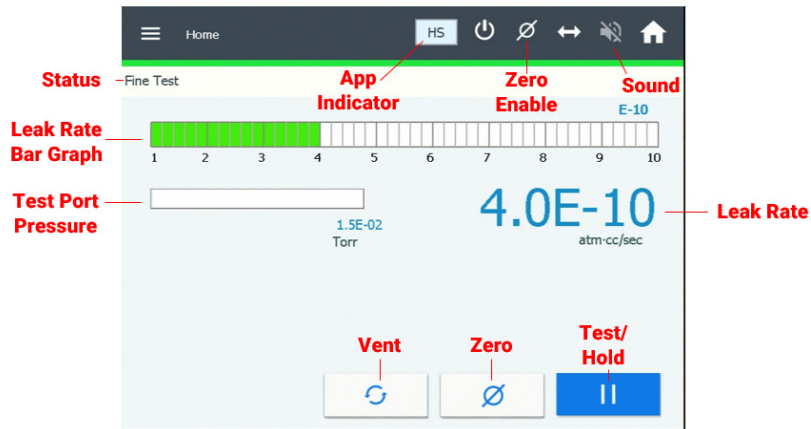
- **Read Internal Leak** - Quick check to see if instrument reads calibrated leak correctly
- **To Calibrate:**
 1. Select Internal Leak and Full.
 2. Tap Calibrate.
 3. Tap OK when calibration is complete.

Menu Navigation (Manual Section 1.5)

Calibration	Calibrate with Internal/External Leak	External leak to be placed on Test Port or vacuum system. Internal leak is provided instrument.
	Calibrate Full/Fast	Full Spectrometer Calibration/ Fast Gain Adjustment
Test Configuration	Set Points	Set Point: Leak Rate, Pressure, Audio
	Advanced Features	High Pressure Test, Fine Test Only, Auto-Zero < 0, Gross Test, Transfer Points (to Test), Stabilization Wait Time
	Application Setup	Sniff Method, Spray Method, High Sensitivity, Auto Sequencer, Split Flow, PPM Sniff Method
Settings	Main Settings	Date and Time, Units, Languages, Touch, and Sound
	Screen Settings	Bar Graph, Range, Analog Output
	Remote Control	Remote Setup with Channel Sync
	Password and Restrictions	Restriction Level, Lock Vent Control, Lock All Control
Maintenance	General	Internal Calibrated Leak Setup, Software Update
	Manual Tuning	Emission, Ion, Gain Settings
	Valve Control	Manually Open and Close Valves
	Pump	Inspect Forepump and Exhaust Pump
	Gauge Calibration	Calibrate Thermocouple Gauge to Vacuum or Atmosphere
	System Defaults	Set to Factory Defaults
System Information	System Information	Software Version, Internal Data
	Current Setup	Set Points and Calibration Type
Turn Off	Turn Off	Turbo will brake. Do NOT move leak detector while turbo is powering down.

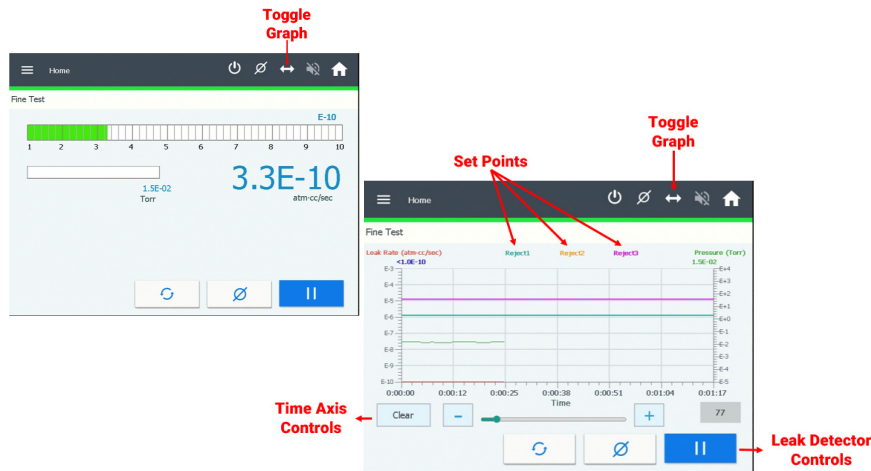


Main Controls (Manual Section 1.5)

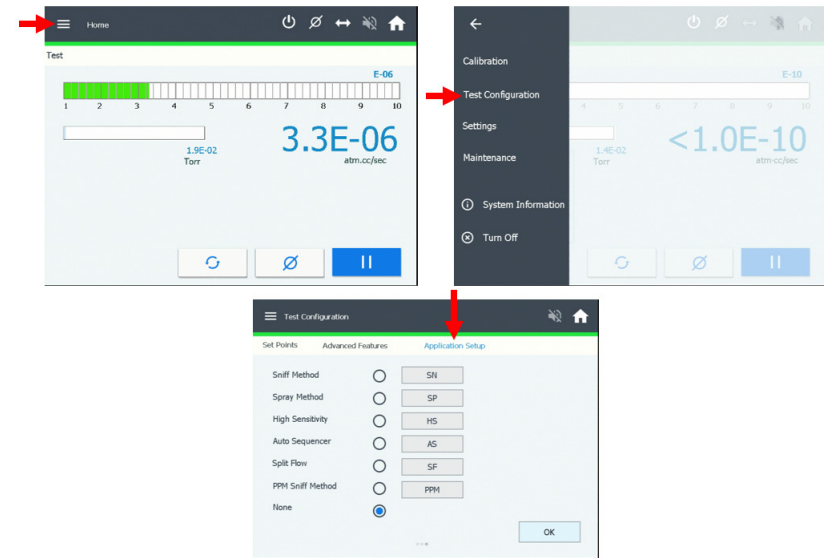


- **Test** - Measure helium leak rate
- **Hold** - Isolate leak detector from connected vacuum system while keeping both under vacuum
- **Zero** - Nullify current leak rate reading and establish a new baseline
- **Vent** - Bring test port to atmospheric pressure

Home Screen Display (Manual Section 1.5 and 3.11)



Application Setup (Manual Section 5.1.3)



Shutdown (Manual Section 1.3.5)

Tap the Power icon and follow the prompts.

