

# Addendum to the Threshold® System User Manual

## Physical Specifications

The following table lists the physical specifications of the Threshold Contaminant Detection System.




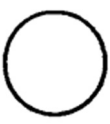


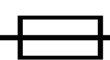
**Table 1-1:** Physical Specifications of the Threshold Contaminant Detection System

Item	Description
Environment	Indoor use only
Power requirements: Threshold	AC Voltage 100/115/220 V, frequency range: 50/60 Hz
Auxiliary Manifold	100 V to 240 V, frequency range: 50/60 Hz
Dimensions	Width 23" Depth 20" Height 8"
Weight	45 Lbs
Ambient operating temperature	15°C to 40°C (59°F to 104°F)
Ambient storage temperature	-5°C to 40°C (23°F to 104°F) continuous; -20°C to 50°C (-4°F to 122°F) transient (up to 10 hours)
Humidity restrictions	0 to 75% (non-condensing)
Altitude restrictions	Up to 2000 m (6562 ft)
Installation category	II
Pollution degree	2
Fuses	250 V, slow-blow UL recognized/CSA/VDE

## Instrument Safety Labels

Safety labels are located on the instrument. Each Safety label consists of a:

- Signal Word panel, which implies a particular level of observation or action (for example, CAUTION or WARNING). If a safety label encompasses multiple hazards, the Signal Word corresponding to the greatest hazard is used.
- Message panel, which explains the hazard and any user action required.
- Safety Alert symbol, which indicates the type of potential personal safety hazard.

Symbol	Indication
	Indicates that product documentation needs to be consulted
	Indicates Alternating Current
	Indicates power on
	Indicates power off
	Indicates the location of the Protective Earth Ground Terminal
	Indicates that you must not discard this electrical/electronic product in domestic household waste
	Indicates fuse

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/productpatents>

Product manufactured by Molecular Devices, LLC.

1311 Orleans Drive, Sunnyvale, California, United States of America 94089.

Molecular Devices, LLC is ISO 9001 registered.

© 2013 Molecular Devices, LLC.

All rights reserved.