



Sartorius SPH Safety Powder Hoods



SPH Safety Powder Hoods – Weigh with Confidence



No duct work required. The pre-filter, primary and backup HEPA filters and blower are housed inside the unit. The blower is a brushless, sparkless motor rated for use 24 hours, 7 days a week. Together they ensure that warm, filtered air is returned to the room saving energy costs.

Constructed of polypropylene. The workstation is constructed of tough, chemically-resistant polypropylene. The three sides, top and base are thermally fused to provide a seamless shell - no glue to breakdown and no metal to rust.



Air Flow Manager. For maximum containment, stable and uniform air flow is maintained by the air flow manager. The enclosure's face velocity is constantly monitored and displayed. Any change, often due to laboratory conditions or filter capacity, automatically initiates an adjustment to the blower speed keeping the air flow constant.

Alarms. A low airflow alarm visually and audibly alerts the user to insufficient airflow. The filter alarm monitors the filter life and alerts the user when the HEPA filter is saturated and needs to be changed.



Full protection during filter change. When changing the primary HEPA filter, the backup HEPA filter provides complete containment. Simply remove the pre-filter and primary HEPA filter from inside the enclosure. Bag them and seal the bag safely inside the workstation.

Ergonomically designed. The Sartorius Mechatronics SPH features an ergonomic, energy efficient design with angled face for worker comfort. The clear polycarbonate face can be opened completely to install equipment. The front sash can be raised to a height of 9", while maintaining full containment, allowing for greater accessibility to the workstation.

Perfect fit. The SPH is small enough to fit on your benchtop, but roomy enough to accommodate up to two lab instruments.

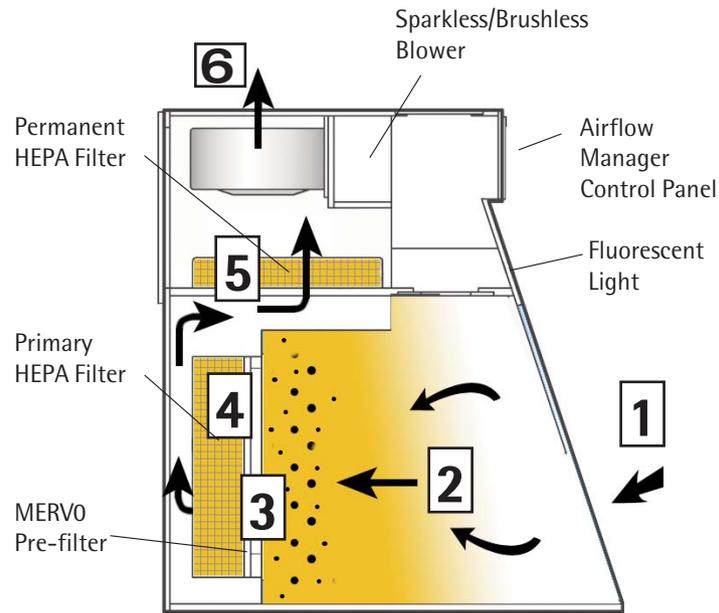
Fully assembled. When you receive the SPH, all you need to do is plug it in and turn it on. All three filters are installed at the factory prior to shipment - where the unit undergoes an extensive quality testing to ensure maximum performance and protection.

Extra features. The florescent light can be turned on for task lighting. Utility ports allow power and communication cords to be connected outside workstation. Lab timer assists with time management.

Disposal Port. This option allows disposal of weighing containers, gloves and other items with traces of powder within the workstation. Simply thread a waste bag through the 4.5" diameter hole and secure inside the workstation. Place contaminated materials inside bag. When full, seal the bag from inside and remove.

Cart provides mobility or extra countertop space. Constructed of 14-gauge epoxy coated steel, these carts are a perfect fit for the SPH. Standard height is under 34" to meet ADA requirements. Choose locking casters to make your workstation portable or stationary feet to expand workspace in the laboratory.

Description of Air Flow



1. Air from the room enters the SPH and is drawn in a laminar flow toward the rear plenum of the enclosure.

2. This air is mixed with the contaminated air which contains powder which is not dense enough to fall to the work surface.

3. Next the air passes through the pre-filter. Here 95% of particles greater than 0.5 microns are trapped. Pre-filters extend the life of the more expensive HEPA filter by capturing large particles that take up precious space in the HEPA filter.

4. The air then enters the primary HEPA filter. Here 99.997% of all particles greater than 0.3 microns are captured. The air that leaves the primary HEPA is filtered and safe to return to the room.

5. The air then passes through the back-up HEPA filter. The back-up HEPA filter is only used during changing of the primary HEPA and comes with a 10-year warranty. Simply stated, if during normal use, the back-up HEPA filter ever needs changing we will send a certified fume hood expert to perform the task at no cost to the customer.

6. Clean, room-temperature air is returned to the room. Compared to traditional fume hoods, energy costs are saved. Instead of taking room air and venting it to the outside, the air is filtered and circulated back into the room.

Features

- Designed specifically for the containment of powder
- Particles are trapped in the pre-filter and HEPA filter
- Save energy by returning filtered, room-temperature air to the room
- Thermally fused Polypropylene to eliminate vibration, provide chemical resistance and wipe-down cleanability.

- Includes HEPA Filter with 99.997% efficiency at 0.3 microns
- Audible and visual alarms warn of any potential problems
- Dynamic air flow volume with real time air velocity feedback
- Continuous HEPA filtration monitoring
- Compatible with MicroBalances
- Modular electronics for easy service

- No installation cost
- Electrical cord and waste disposal ports
- Ductless design enables easy portability
- Available in 32" or 48" widths

Specification	32"	48"
Maximum volume of air filtered	350 CFM	600 CFM
Included Filters	prefilter (1); primary HEPA (1); backup HEPA (1)	prefilter (1); primary HEPA (1); backup HEPA (1)
Electrical Voltage	110V (optional 220V)	110V (optional 220V)
Electrical Current	3.0A (1.5A for 220V)	6.0A (6.0A for 220V)
Electrical Consumption	330W	660W
External Dimensions (L x W x H)	30 x 32 x 32.5" (762 x 813 x 826mm)	30 x 48 x 32.5" (762 x 1219 x 826mm)
Weight	175 lbs. (79.4kg)	250 lbs. (113.4kg)
Minimum Installation Clearance	3 " on right side; 4" on top	3 " on right side; 4" on top
Model No.	SPH32	SPH48
Model No. with Disposal Port	SPH32B	SPH48B
Model No. 220V Option	SPH322	SPH482
Model No. 220V Option with Disposal Port	SPH32B2	SPH48B2

Warranty

1 year

Accessories

	Part No.
Pre-filter for 32" unit (pack of 12)	6009.1
Pre-filter for 48" unit (pack of 2)	6010.1
HEPA for 32" unit	6011.1
HEPA for 48" unit	6012.1
Cart with wheels for 32" unit	6013.1
Cart with wheels for 48" unit	6014.1
Stationary cart for 32" unit	6015.1
Stationary cart for 48" unit	6016.1



Sartorius Mechatronics Corporation

5 Orville Drive

Bohemia, NY 11716

Phone: 631.254.4249

Fax: 631.254.4253

www.sartorius-mechatronics.com

Lit Code SPH