



SWC Safety Weighing Cabinet

Personnel Safety and
Reliable Weighing Results
From a Single Source

Simplifying Progress

SARTORIUS

Greater Protection and Reliability in Weighing Hazardous Powdered Compounds

For researchers and laboratory technicians, the risks involved with handling toxic and high-potency compounds are considerable. Weighing even the smallest sample quantities can expose them to potential hazards.

Appropriate protective measures must be taken to protect the health of all laboratory staff when they weigh hazardous compounds in powdered form.

The Sartorius SWC Safety Weighing Cabinet offers important advantages in design and construction over traditional laboratory hoods: The SWC creates a contained area around the laboratory balance which prevents any air or finely powdered particulates from escaping into the operating personnel's work area. At the same time, the constant inlet air velocity and the special cabinet construction both keep the air current practically turbulence-free, and, as a result, ensure consistent and repeatable weighing results.

The balance and weighing cabinet are perfectly matched to each other. With its SWC Safety Weighing Cabinet, Sartorius has succeeded in uniting two otherwise contradictory requirements: maximum personnel protection and reliable weighing results.

The Safety Weighing Cabinets are available in four different dimensions for special applications, such as using a second laboratory balance in the cabinet or for unusually high structures:

SWC900	W 890 × D 750 × H 510 mm
SWC1200	W 1,230 × D 750 × H 510 mm
SWC900T	W 890 × D 750 × H 770 mm
SWC1200T	W 1,230 × D 750 × H 770 mm

Each of the four basic models consists of:

Safety Weighing Cabinet with a separate HEPA filter unit, data-logging alarm, lighting unit, waste disposal system (on one side), airflow smoke test kit and antistatic cleaning wipes.

Sartorius SWC Safety Weighing Cabinets comply with the requirements of EN14175.

The clear acrylic glass construction affords the analyst a perfect view of the balance and the sample to be weighed. This feature is particularly important for working with extremely small sample quantities. All surfaces are easy to clean and compatible with conventional cleaning agents.





The extremely stable base made of black granite absorbs vibrations that can affect the balance. The analyst can instantly see even the slightest spillage. Contaminated materials can be disposed of safely in a double-bag system.



The calibrated airflow alarm system with audit trail function sounds an alarm whenever the airflow velocity drops below a defined level.



The filter unit with an HEPA H14 filter is free-standing to effectively prevent vibrations. The "bag in, bag out" filter system allows the filter to be changed quickly and contamination-free.

Safety Weighing Cabinets

Model with filter unit	Model without filter unit	Dimensions in mm (width × depth × height)
SWC900	SWC900NF	890 × 750 × 510
SWC1200	SWC1200NF	1,230 × 750 × 510
SWC900T	SWC900TNF	890 × 750 × 770
SWC1200T	SWC1200TNF	1,230 × 750 × 770

Accessories

YWCF02	Carbon filter for solvent vapors
YWCF03	Box for carbon filters; fixes to the fan filter box
YWCG01	Disposal chute for waste; fits to side of cabinet
YWCG02	Disposable chute bags, qty. of 100
YWCG03	Silencer unit option; attaches to exhaust of fan
YWCG04	Airflow smoke test kit
YWCG07	Anti-static decontamination wipes

YWCG16	Printer table for mounting on the Safety Weighing Cabinet
YWT10	Laboratory bench: fits SWC900, SWC900T and SWC900NF
YWT11	Laboratory bench: fits SWC1200, SWC1200T and SWC1200NF

Other accessories for our Safety Weighing Cabinets are available on request.




🌐 Find out more
For more information, please visit
www.sartorius.com

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further contacts, visit
www.sartorius.com