



**sartorius**  
mechatronics

## SWC Safety Weighing Cabinet: Personnel safety and reliable weighing results from a single source



turning science **into solutions**

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.

# 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 1230 × D 750 × H 510 mm
SWC900T	W 890 × D 750 × H 770 mm
SWC1200T	W 1230 × 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 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	1230×750×510
SWC900T	SWC900TNF	890×750×770
SWC1200T	SWC1200TNF	1230×750×770

## Accessories

YWCF02	Carbon filter for solvent vapors	YWCG07	Anti-static decontamination wipes
YWCF03	Box for carbon filters; fixes to the fan filter box	YWCG16	Printer table for mounting on the Safety Weighing Cabinet
YWCG01	Disposal chute for waste; fits to side of cabinet	YWT10	Laboratory bench: fits SWC900, SWC900T and SWC900NF
YWCG02	Disposable chute bags, qty. of 100	YWT11	Laboratory bench: fits SWC1200, SWC1200T and SWC1200NF
YWCG03	Silencer unit option; attaches to exhaust of fan		
YWCG04	Airflow smoke test kit		

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

All of the balances listed below were tested for use in the Safety Weighing Cabinet and achieved their typical repeatability when weighing times are extended accordingly.

Balance series	ME SE	LA Reference	Sartorius CP	Extend ED
Microbalances	ME5 ME36S		CPA2P	
Semi-microbalances	ME235S ME235P		CPA225D	
Analytical balances	ME614S ME414S ME254S	LA310S LA230S LA230P LA120S	CPA324S CPA224S CPA124S CPA64	ED224S ED124S
Precision balances		LA1200S LA620S LA220S LA2000P LA620P LA5200D LA3200D	CPA1003S CPA1003P CPA623S CPA423S CPA323S CPA223S CPA2202S-DS CPA5202S-DS	ED623S ED423S ED323S ED153  All models listed are also available in – CW versions

Sartorius AG  
Weender Landstrasse 94–108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

[www.sartorius-mechatronics.com](http://www.sartorius-mechatronics.com)

Specifications subject to change without notice. Printed in Germany on paper that has been bleached without any use of chlorine.  
W · G  
Publication No.: WL-1008-e08082  
Order No.: 98649-007-78