

# Midisart® 2000

## Ready-To-Use Filter for Sterilizing Gases and Venting



### Introduction

Midisart® 2000 is designed for maximum handling ease and safety. The device is lightweight only 20 g – so it will not weigh down or kink the tubing. With a diameter measuring just 64 mm, Midisart® incorporates a filter area of 20 cm<sup>2</sup>, which means that it is “packed” with high flow rate performance power.

Midisart® 2000 filters can be integrity tested (Bubble Point Test, 60% IPA | Water, BPmin. = 1.1 bar for 0.2 µm and BPmin. = 0.9 bar for 0.45 µm). The Midisart® Test Manifold (Order code 1Z-LB-0002) can be used to test 10 filters in parallel which simply leads to a reduction of test time by 90%.

### Product Information

The Midisart® 2000 is an ideal choice for small-scale sterile air | gas filtration applications. Superior flow rates and seamless integration into single-use assemblies provide Midisart® customers with a distinct advantage that directly translates into cost savings. Individual 100% factory integrity testing verifies the quality of every filter prior to use.

## Venting Applications

- Filling and Transfer Vessels
- Fermentation Carboys
- Cell Culture Chambers
- CO<sub>2</sub> incubators (6 to 120 liters)
- Holding and Storage Tanks
- Autoclave venting
- Small Bioreactors

## Technical Data

### Materials

Membrane PTFE – reinforced with polypropylene gauze  
Housing Polypropylene

### Pore Sizes

0.2 µm  
0.45 µm

### Filtration Area

20 cm<sup>2</sup> | 3 square inch

### Max. Differential Pressure for Air

3 bar (in-out)  
2 bar (out-in)

### Housing Diameter

64 mm | 2,5"

### Connectors

Hose Barb, 1/8" NPT, Triclamp, small Hose Barb

## We Prove Your Trust

The single-layer hydrophobic PTFE membrane is self-dewetting and therefore provides high air flow rates even at very low differential pressures over an extended process time. All components were developed and are produced under full control of Sartorius. Furthermore the following indispensable requirements are fulfilled:

- **100% integrity tested prior to release**
- **Clear labeling** with Lot number, individual unit number and "IN" as an indicator for direction of flow. All information lasered on the top part of each filter housing.
- **Retention of ≥ 10<sup>7</sup> Brev. diminuta/cm<sup>2</sup>** in aerosol and liquid

### Sterilization

Max. Temperature 134 °C  
Max. Autoclaving 20 Cycles  
Not suitable for gamma irradiation  
Gamma irradiatable disc filter: Midisart® Sartopore® Air

### Integrity Testing

Min. Bubble Point >1100 mbar | 15.95 psi for 0.2 µm  
Min. Bubble Point >900 mbar | 13.05 psi for 0.45 µm  
Wetting agent: IPA/Water (60 | 40)

### Regulatory Compliance

- Each individual element is tested for integrity in direction of filtration
- 0.2 µm Pore size: Fully validated as sterilizing grade filters according to ASTM current F-838 guidelines
- Designed, developed and manufactured in accordance with an ISO 9001 certified Quality Management System
- USP Plastic Class VI Test
- Non fiber releasing according to 21 CFR

### Technical References

A Validation Guide is available under DIR No. 2729708.

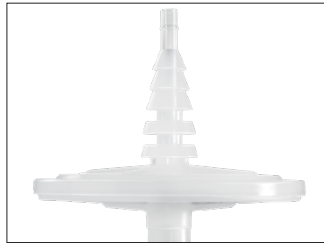
# Ordering Information

Order No.	Pore Size [µm]	Membrane	Connectors	Qty.	Sterile
17804-----E	0.45	PTFE	Hose Barb   Hose Barb	12	Yes
17804-----G	0.45	PTFE	Hose Barb   Hose Barb	25	Yes
17804-----NPE	0.45	PTFE	½"   ½" NPT	12	Yes
17804-----NPG	0.45	PTFE	½"   ½" NPT	25	Yes
17805-----E	0.2	PTFE	Hose Barb   Hose Barb	12	Yes
17805-----G	0.2	PTFE	Hose Barb   Hose Barb	25	Yes
17805-----NPE	0.2	PTFE	½"   ½" NPT	12	Yes
17805-----NPG	0.2	PTFE	½"   ½" NPT	25	Yes
17805-----UPN	0.2	PTFE	Hose Barb   Hose Barb	100	No
17805-----UPQ	0.2	PTFE	Hose Barb   Hose Barb	500	No
17805-----UQN	0.2	PTFE	Hose Barb   Hose Barb	100*	No
17809-----UNN	0.2	PTFE	½"   ½" NPT	100	No
17812-----UNN	0.2	PTFE	½"   Hose Barb	100	No
17805-----TCN	0.2	PTFE	TriClamp   TriClamp	100	No
17877-----UPN	0.2	PTFE	Small Hose Barb   Small Hose Barb	100	No

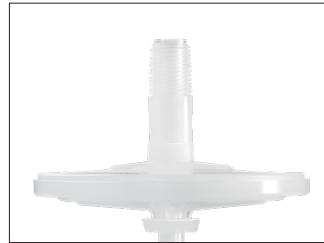
\* = plastic bucket - the bucket with its smooth surface can be easily transferred into clean-room areas without risking particle release from the packaging material.



Standard Hose Barb



Small Hose Barb



½" NPT Thread




TriClamp

**Germany**

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

**USA**

Sartorius Stedim North America Inc.  
565 Johnson Avenue  
Bohemia, NY 11716  
Toll-Free +1 800 368 7178

 For further contacts, visit  
[www.sartorius.com](http://www.sartorius.com)