



MD8 Airscan®

Active sampling of airborne viruses, bacteria, yeasts and fungi

Product Information

Monitoring ambient air in cleanrooms and controlled areas in the (c)GMP environment for viable microorganisms is a routine task in the pharmaceutical | biotechnology industry. As a rule, collection of data and subsequent analysis of the results permit trends to be recognized early and problems (OOS*) with bioburden levels exceeding the limits to be prevented.

* Out of specification



Active Air Monitoring in Cleanrooms and Controlled Areas

Active monitoring of viable airborne microorganisms is an important instrument for ensuring product quality and safety. Sampling of at least 1 m³ of air is required by regulation standards. In this procedure, care must be taken that turbulence is not caused by, for example, devices such as active air samplers and that sampling is performed isokinetically.

The following areas need to be monitored using active sampling of airborne microorganisms:

In the pharmaceutical | biotechnology industry:

- Cleanrooms (grades A – D; ISO classes 5 – 9)
- Sterility testing isolators
- Filling lines
- Controlled areas, e.g. quality control (QC) laboratories
- Warehouses | raw materials depots
- Personnel airlocks and aseptic transfer chambers for materials

In the food and beverage industry:

- Manufacturing areas and production-related areas
- Quality control (QC) laboratories
- Warehouses | raw materials depots

Diagnostics | Medicine:

- Cleanroom or controlled area
- Filling lines
- Manufacturing areas and production-related areas
- Quality control (QC) laboratories
- Warehouses | raw materials depots

Quality Control

As Sartorius is a recognized supplier of products and services for the pharmaceutical industry, product quality and safety are its number one priorities. To meet stringent quality requirements, 100% final testing of the MD8 Airscan® is performed in manufacturing among the quality control checks carried out. In addition, every stainless steel sampling head is calibrated and adjusted after assembly.

Technical Specifications

Types of MD8 Airscan® Stainless Steel Sampling Heads

Stainless steel sampling head with bayonet connector

Specially designed as a "stand-alone" version"

- Type of connector: Bayonet
- Flow capacity: Maximum flow rate capacity: up to 7 m³/h; adjustable in increments of 0.1 m³/h
- Saves the calibration data and actively regulates air flow
- Order number for sampling head: 16746-01--SHB
- Order number for bayonet package version: 16746SHBCOM

Contents of package version:

- Command unit with 3 m connecting cable
- Stainless steel sampling head with bayonet flange connector
- Aluminum filter holder*



Stainless steel sampling head with Tri-Clamp connector

Specially designed for integrating into lines

- Type of connector: Tri-Clamp, 50.5 (TC 50.5)
- Maximum flow rate capacity: up to 7 m³/h; adjustable in increments of 0.1 m³/h
- Saves the calibration data and actively regulates air flow
- Order number for sampling head: 16746-01--SHT
- Order number for Tri-Clamp package version: 16746SHTCOM

Contents of package version:

- Command unit with 3 m connecting cable
- Stainless steel sampling head with Tri-Clamp connector
- Stainless steel adapter, for connection of TC 50.5 to filter holder
- Aluminum filter holder*



MD8 Airscan® Command Unit

- Plastic material PA6 with 30% glass fiber
- RS-232 interface port for connecting a printer
- 2 status LEDs for displaying operating status and errors
- Stores the test results
- Order number of the command unit: 16746-01--COM



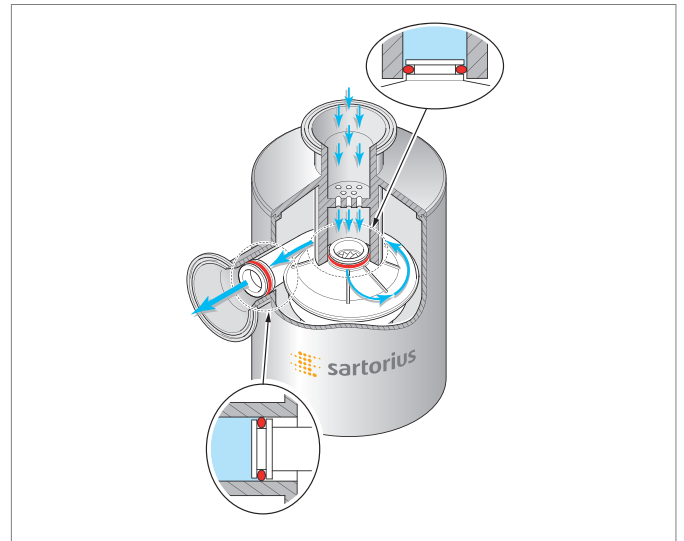
*The filter holder is also available in an AISI 316L stainless steel version (German equivalent standard 1.4404).

Disinfection | Sterilization

- Stainless steel sampling heads (bayonet and TC versions)
- IPA (70%)
- Ethanol (70%)
- Peracetic acid (1%)
- Dismozon® (1%)
- Sodium hypochlorite (5%)
- Formaldehyde (4%)
- Flowpath can be disinfected or sterilized | inline with vaporized hydrogen peroxide

Command unit

- IPA (70%)
- Ethanol (70%)
- Peracetic acid (1%)
- Dismozon® (1%)
- Sodium hypochlorite (5%)
- Formaldehyde (4%)



Specifications of the Stainless Steel Sampling Heads, Bayonet and TC Versions

- Electropolished stainless steel: AISI 316L (German equivalent standard) 1.4404 or better

General Information

	Unit	Sampling Head with Bayonet Connector	Sampling Head with Tri-Clamp Connector (TC 50.5)
Dimensions (L × W × H)	mm inches	128 × 112 × 137 ~5 × 4.4 × 5.4	128 × 112 × 154 ~5 × 4.4 × 6
Weight	kg lb	2.6 ~6	2.6 ~6
Power supply			
Only through MD8 Airscan® command unit			
Material			
Gaskets		Silicone	Silicone
Surfaces		Stainless steel	Stainless steel

Ambient Conditions

	Unit	Specification
Installation site		
Usual laboratory spaces and cleanrooms as defined by ISO 5 to ISO 9; use in controlled areas: Cleanroom classification A–D	m above sea level ft. above sea level	<3,000 <9,843
Required clearance for air exhaust outlet	mm inches	> 100 4
Ambient temperature		
Operation	°C	4 – 40
Storage and transportation	°C	4 – 45

Specifications of the Stainless Steel Sampling Heads, Bayonet and TC Versions

	Unit	Specification
Max. temperature of motor in the sampling head	°C	65
Relative humidity		
Operation	%	50–80
Storage	%	20–90
Electromagnetic compatibility according to DIN EN 61326-1: Electrical equipment for measurement, control and laboratory use – EMC requirements – Part 1		
Defined immunity to interference: Suitable for use in industrial areas		
Limitation of interference	Class	B

Command Unit Specifications

Technical Specifications

Dimensions and Weight

	Unit	Specification
Dimensions (L×W×H)	mm inches	282×102×107 11×4×4.2
Weight	kg lbs	1.4 ~3.1

Ambient Conditions

	Unit	Specification
Installation site		
Usual laboratory spaces and cleanrooms as defined by ISO 5 to ISO 9; use in controlled areas: Cleanroom classification grades A–D	m above sea level ft. above sea level	< 3,000 < 9,843
Required clearance for air exhaust outlet	mm inches	> 100 4
Ambient temperature		
Operation	°C	4–40
Storage and transportation	°C	4–45
Max. temperature of motor in the sampling head	°C	65
Relative humidity		
Operation	%	50–80
Storage	%	20–90
Electromagnetic compatibility according to DIN EN 61326-1: Electrical equipment for measurement, control and laboratory use – EMC requirements – Part 1		
Defined immunity to interference: Suitable for use in industrial areas		
Limitation of interference	Class	B

Measurement Parameters

	Unit	Specification
Adjustable volume: adjustable in increments of 0.01 m ³		
For gelatin membrane filters	m ³	0.01–9.99
For cellulose nitrate membrane filters	m ³	0.01–1.10
Adjustable air flow: adjustable in increments of 0.1 m ³		
With gelatin membrane filters	m ³ /h	1.8–7
With cellulose nitrate membrane filters	m ³ /h	1.8–3

Interfaces

RS-232 interface port for connecting a printer

Suitable printer: Sartorius YDP30 (baud: 9600; parity: odd; data bit: 8; stop bit: 1)

Power Supply

Power Supply Unit	Unit	Rating Specification
Type: Sartorius power supply, model 1000018304		
Primary		
Voltage	V_{AC}	100 – 240 \pm 10%
Frequency	Hz	50 – 60
Power consumption, max.	A	2.0
Secondary		
Voltage	V_{DC}	24 \pm 5%
Power, max.	A	6.25
Short-circuit protection: Electronic		
Protection class according to DIN EN/IEC 60950-1		I
Installation site according to DIN EN/IEC 60529-1		m above sea level ft. above sea level
		< 3,000 < 9,843
Installation category (overvoltage) according to IEC 60950-1		II
Contamination level according to IEC 60950-1		2
Connecting cable for power supply		Class
		B
Connecting cable according to DIN EN/IEC 60320-1/C14: country-specific, 3-pin; can be plugged in at both ends		
Socket according to DIN EN/IEC 60320-1/C14: 3-pin		
Additional data: See label on the power supply		




Command Unit and Sampling Head	Unit	Rating Specification
Power supply for control unit: Only via Sartorius power supply, model 1000018304		
Voltage	V_{DC}	24 \pm 5%
Power consumption	A	5.0
Power supply for sampling head: Only through command unit		

Ordering Information

Recommended Accessories and Consumables








Accessories

This table shows a sampling of the range of accessories that can be ordered. Please contact Sartorius for information on additional products.

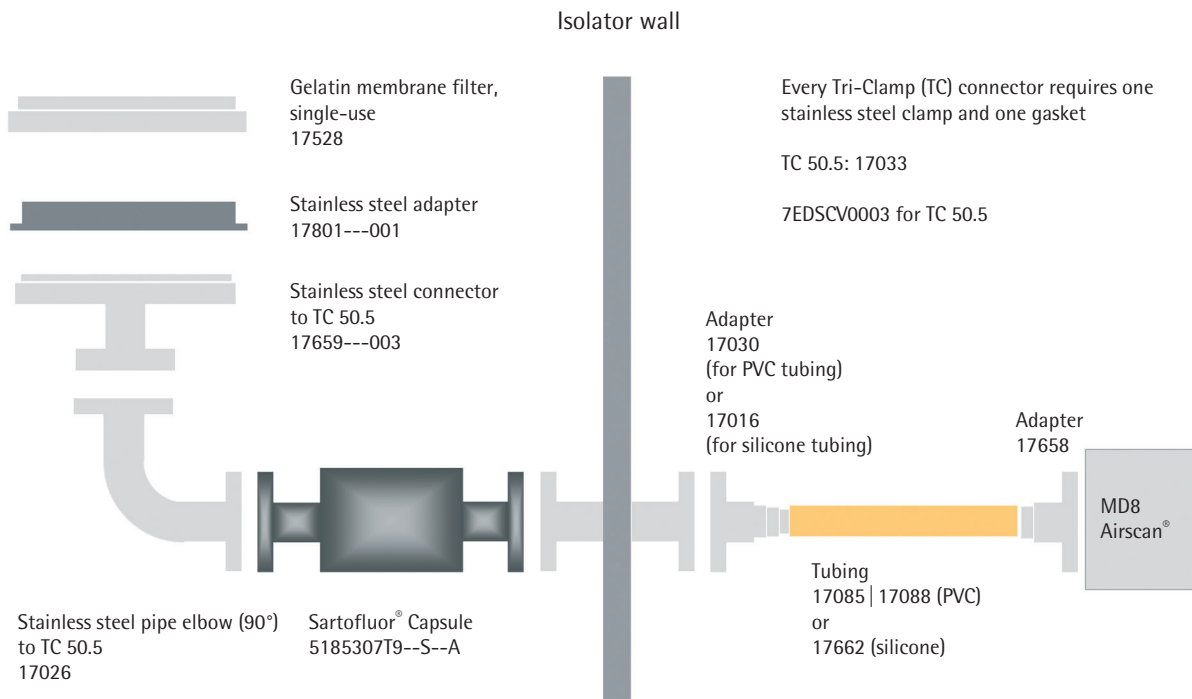
	Product	Quantity	Order Number
	Adapter for gelatin membrane filter	1	17801
	Adapter, Tri-Clamp (TC 50.5) to bayonet, stainless steel	1	17659---003
	Adapter, bayonet to Tri-Clamp (TC 50.5), stainless steel	1	1ZAD-0025
	Adapter, TC 50.5 to single-step DN 25 hose barbs, stainless steel. Can be used on the air outlet for conducting exhaust air outside the area.	1	17016
	Clamp for 1" to 1½" Sanitary flange	1	17033

Consumables

This table shows a sampling of the range of consumables that can be ordered. Please contact Sartorius for information on additional products.

	Product	Quantity	Order Number
	Gasket for Tri-Clamp TC 50.5 connector, silicone	2	7EDSCV0003
	Sartofluor® capsule, sterilizing-grade air filter, 0.2 µm, 0.2 m² filter area	4	5181307T9--SS--A
	PVC tubing, 2 m (~6.5 ft.)	1	17085
	PVC hose, 5 m (~16 ft.)	1	17088
	Silicone tubing, 1 m (~3.3 ft.)	1	17662
	Quick-connect cable	1	69898530
	Cable for YDP30 printer	1	YCC01-0041M3

Installation of the MD8 Airscan® | Isolator



Sartorius Service

We ensure the quality of your results

From equipment qualification IQ | OQ to regular maintenance: We offer you just the right choice of services to ensure that you comply with the high regulatory requirements right from the beginning, especially in GMP environments. Not only that: we will take care that your air sampler operates reliably over the long term.

Get more information at
www.sartorius.com/service

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen

Phone +49.551.308.0
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan

Phone +33.1.70.62.50.00
Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles

Phone +32.2.756.06.90
Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki

Phone +358.9.755.951
Fax +358.9.755.95.200

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi

Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11

Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Via Torino 3/5
20814 Varedo (MB)

Phone +39.0362.5557.11
Fax +39.0362.5557.99

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001
Fax +31.30.60.52.917

info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn

Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg

Phone +7.812.327.53.27
Fax +7.812.327.53.23

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)

Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800
Fax Spain +34.913.589.623
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)

Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ

Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716

Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. AVALOS 4251
B1605ECS Munro
Buenos Aires

Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900
Fax +55.11.4362.8901

Canada

Sartorius Canada Inc
1173 North Service Road West, D4
Oakville, ON L6M 2V9

Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605

Phone +52.55.5562.1102
Fax +52.55.5562.2942

leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima

Phone +51.1.441 0158
Fax +51.1.422 6100

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175

Phone +61.3.8762.1800
Fax +61.3.8762.1828

China

Sartorius (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China

Phone +86.21.6878.2300
Fax +86.21.6878.2882

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong

Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India

Phone +91.80.4350.5250
Fax +91.80.4350.5253

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408
Fax +81.3.3740.5406

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684

Phone +65.6872.3966
Fax +65.6778.2494

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310

Phone +66.2643.8361-6
Fax +66.2643.8367



◀ www.sartorius.com