



BACTair™

A New Method for Microbiological Air Monitoring



Description

As regulations become more stringent, the microbial monitoring of ambient air is gaining increasing importance in today's world. For example, microorganisms can have a major impact on product quality and the production process, and can even pose health hazards.

Sartorius has developed a new system for sampling airborne microorganisms that allows impactation onto culture media plates, where the plates function directly as collection heads.

This new method uses the AirPort MD8 air sampler to pump the air stream, as this sampler has proved its efficacy for years in the gelatin membrane filter method.

Technical Specifications

BACTair™ culture media plates

Material	Polystyrene
Dimensions	116 × 24 mm
Number of impactation holes	400 holes, each 0.47 mm Ø
High retention of particles	> 0.65 µm

AirPort MD8 Air Sampler for BACTair™ culture media plates

Flow rate control	By integrated vane anemometer
Flow rate	125 l/min
Predefined sampling volume	25, 50, 100, 250, 500, 750 and 1000 liters
User-definable sampling volumes	From 10 to 2000 l in 5 l-increments
Total sampling time with one battery charge	Approx. 4 hours at 125 l/min
Power consumption of fan and electronics	Approx. 10 watt
Noise level	< 48 dB (A)
Weight	Approx. 2.5 kg
Dimensions	300 × 135 × 165 mm
CE approved	yes

Power Supply of Air Sampler

Rechargeable battery pack	NiMH 16.8 V/3800 mA
Charger	Input 100-240 V/ 47-63 Hz/600 mA Output 24 V/ 1000 mA
Charging time	Approx. 4.5 h when the battery is fully discharged

The air sampler can also be operated while charging.

Ordering Information

BACTair™ culture media plates

Description	Order Number
BACTair™ culture media plates with TSA agar, 110 mm, sterile and individually packaged, 10 units	14320-110----ACD
BACTair™ culture media plates with Sabouraud agar, 110 mm, sterile and individually packaged, 10 units	14321-110----ACD

Air Sampler

Description	Order Number
AirPort MD8 Air Sampler for BACTair™ culture media plates incl. charger	16757

Accessories for the AirPort MD8 Air Sampler BACTair™ culture media plates

Description	Order Number
Adapter for BACTair™ culture media plates on the AirPort MD8 air sampler	17803
Covers for BACTair™ culture media plates, 10 × 2 units individually, sterile packaged	1ZPX-D0002
Charger	6989525

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany
Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius.com
USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Russian Federation +7.812.327.5.327
Japan +81.3.4331.4300

Specifications subject to change
without notice. Printed and copyrighted
by Sartorius Stedim Biotech GmbH. | W
Publication No.: SL-2047-e140103
Order No.: 85032-533-89
Ver. 01 | 2014