

## 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 impaction 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

<b>BACTair™ culture media plates</b>	
Material	Polystyrene
Dimensions	116 × 24 mm
Number of impaction holes	400 holes, each 0.47 mm Ø
High retention of particles	>0.65 µm
<b>AirPort MD8 Air Sampler for BACTair™ culture media plates</b>	
Flow rate control	By integrated vane anemometer
Flow rate	125 L/min
Predefined sampling volume	25, 50, 100, 250, 500, 750 and 1,000 liters
User-definable sampling volumes	From 10 to 2,000 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
<b>Power Supply of Air Sampler</b>	
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


Description	Order Number
<b>BACTair™ culture media plates</b>	
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
<b>Air Sampler</b>	
AirPort MD8 Air Sampler for BACTair™ culture media plates incl. charger	16757
<b>Accessories for the AirPort MD8 Air Sampler BACTair™ culture media plates</b>	
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	YEPS02-24V0

## 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 information, visit  
[www.sartorius.com](http://www.sartorius.com)