

# Monitoring Airborne Microorganisms During Food and Beverage Processing

■ Comparison of four microbial air samplers



Air monitoring provides information on the quality of the processing environment during manufacturing and enables the study of microbiological air quality trends (Figure 1).

There are two basic methods of air monitoring: passive and active. Passive monitoring, such as the use of settling plates, can only detect microorganisms that fall onto the surface of the medium. Active air sampling systems draw a volume of air from the test area and, therefore, provide a more representative sample.

*Air monitoring is a GMP requirement in the United States and Europe. It is also an important part of many HACCP compliance programs.*

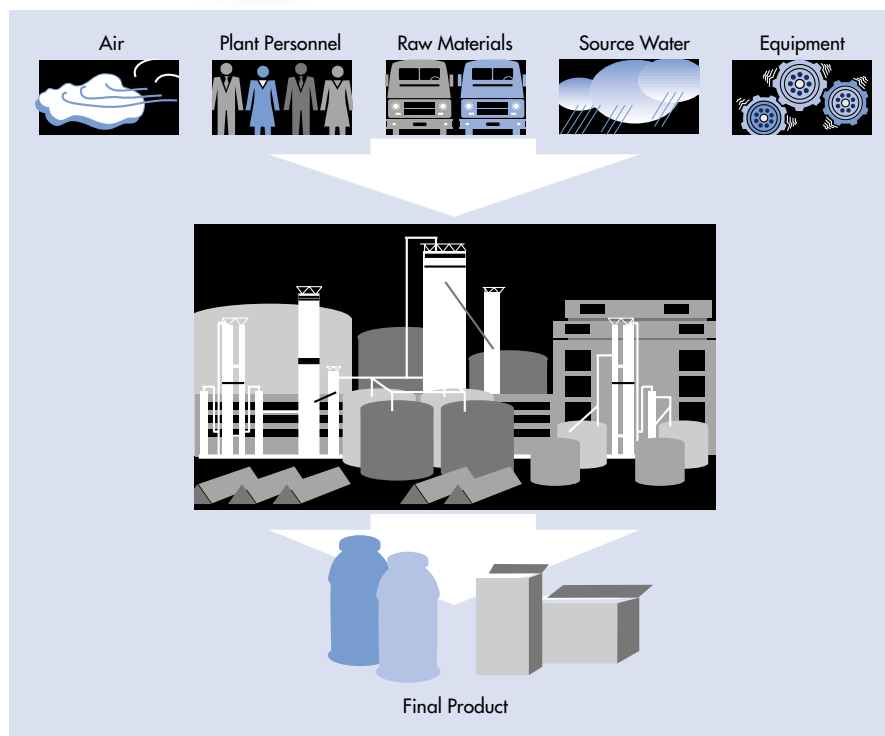


Figure 1. Air is one of the five primary sources of particulate and microbial contamination.

The purpose of this study was to compare the performance of a new, active microbial air sampler (M Air T), to a Slit-to-Agar (STA) sampler and two other widely used air samplers. The M Air T System uses the sieve-impaction sampling method (Figure 2).

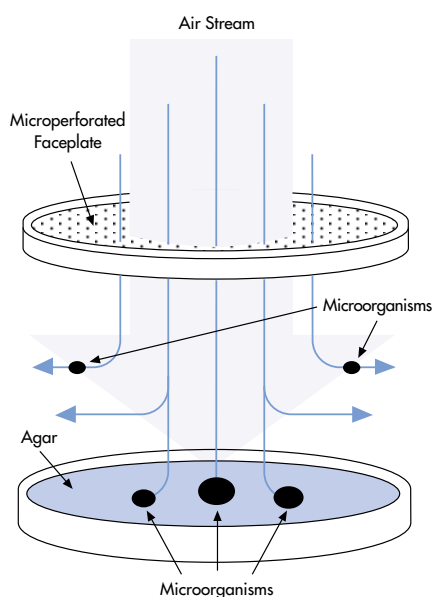


Figure 2. The sieve-impaction sampling method.

## Materials

- Four active microbiological air samplers were selected for evaluation (refer to Table 1 for operating parameters).
  - M Air T Air Tester (Millipore Corporation, Bedford, MA)
  - RCS PLUS Air Sampler (Biotest Diagnostics, Denville, NJ)
  - RCS High Flow Air Sampler (Biotest Diagnostics, Denville, NJ)
  - Slit to Agar Biological Air Sampler, STA 203 (New Brunswick Scientific Co., Edison, NJ)
- Air Current Tubes (Dräger, Sicherheitstechnik, Germany)
- Recovery agar was standard Soybean Casein Digest Agar (also known as Tryptic Soy Agar or TSA). Multiple lots of cassettes, strips, and dishes were used.
  - M Air T TSA Cassettes for M Air T (Millipore Corporation, Bedford, MA)
  - Biotest TSA Test Strips for RCS PLUS and High Flow (Biotest Diagnostics, Denville, NJ)
  - TSA dehydrated powder reconstituted and sterilized according to manufacturer's instructions at 65 mL per 150 mm Petri dish for STA 203 (Difco, Detroit, MI)

## Method

A two-factor, four-level, full-factorial blocked design was used for this study. The factors were Air Tester (M Air T, RCS PLUS, RCS High Flow, STA) and Test Position (P1, P2, P3, P4). The four samplers were tested simultaneously as a group. After all samplers completed sampling 1 m<sup>3</sup> (1,000 L) of air, the samplers were switched for test position. One test block consisted of all air samplers tested in all test positions for a total of 16 tests in four groups. Six test blocks were completed in a single experiment for a total of 96 tests in 16 groups.

Five experiments each were conducted in a controlled-access, but unclassified, testing room (room volume 87 m<sup>3</sup>). All samplers were placed at least 1 m from the ceiling and 1 m from each other. While the samplers operated simultaneously, their orientation was tested for exhaust turbulence and other potential air current disturbances using Dräger air current tubes.

All TSA agar plates, cassettes, and strips were weighed before and after testing. All agar was incubated in humidified environments for 48 hours at 35 °C followed by 72 hours at 25 °C. Samples from the unclassified test room were enumerated at both 48 and 72 hours.

Table 1.

Properties of the Four Microbial Air Samplers

	M Air T (Millipore)	STA-203 (New Brunswick Scientific)	RCS High Flow (Biotest)	RCS PLUS (Biotest)
Particle capture mechanism	Impaction	Impaction	Centrifugal impaction	Centrifugal impaction
Sampling volume (maximum)	1,000 L	3,000 L	1,000 L	1,999 L
Time to sample 1,000 L (1 m <sup>3</sup> )	6.5 min (140 L/min)	20 min (50 L/min)	10 min (100 L/min)	20 min (50 L/min)
	1st 500 L, 180 L/min 2 <sup>nd</sup> 500 L)			
Water loss after sampling 1 m <sup>3</sup> *	4.2%	4.3%	10%	13%
Particle diameter cutoff size (d <sub>50</sub> )	3.5 µm	0.5 µm	2-5 µm	2-5 µm
Aspiration mechanism	Impeller below agar cassette	Vacuum required	Impeller perpendicular to agar strip	Impeller perpendicular to agar strip
Agar volume	34 mL	65 mL	8 mL	8 mL

\*In an unclassified environment with 53% relative humidity. Sample size was 120 per tester.

## Results

Overall results for the unclassified room studies showed a significant difference among samplers and a significant difference among test positions.\*

The overall analysis for the unclassified room suggested a significant difference in recovery for certain pairs of samplers. A pairwise analysis performed on the 48-hour and 72-hour plate counts demonstrated no significant difference between the M Air T and the STA samplers and no significant difference between the RCS PLUS and RCS High Flow samplers. There was a significant difference in recovery between the grouped M Air T/STA and the grouped RCS PLUS/RCS High Flow data (Figure 3).

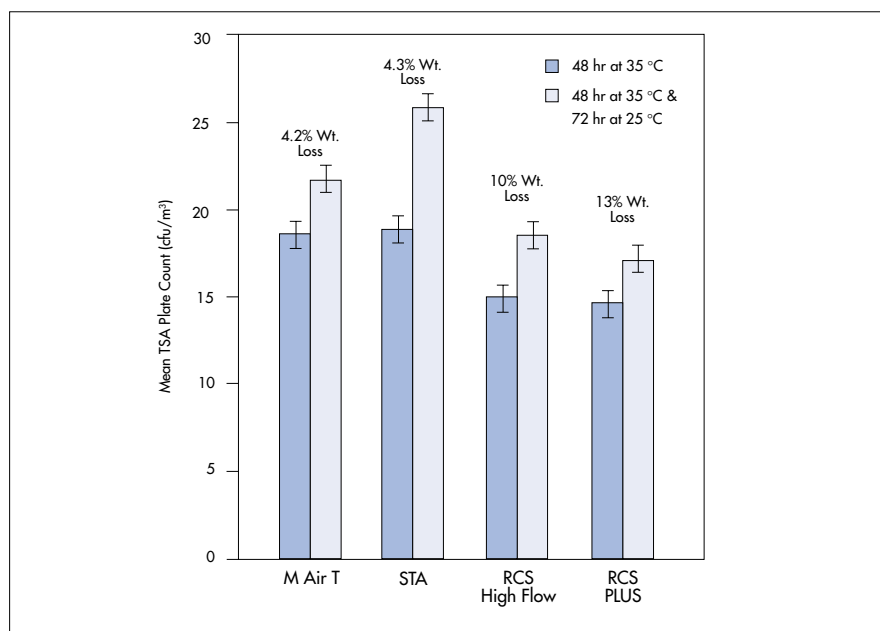
## Discussion

A significant increase in plate counts was observed during the two-stage incubation of 48 hours at 35 °C followed by 72 hours at 25 °C for the unclassified area.

The percent weight loss of the agar in the unclassified room appeared to be related to recovery. M Air T and STA had lower percent water loss and higher microbial recovery as compared to RCS PLUS and RCS High Flow. The mean final incubation plate count ranged from 18 to 26 cfu.

Figure 3.

Mean Recovery and Standard Error for Four Different Microbial Air Samplers Operated Simultaneously in an Unclassified Test Room.



Conditions: 53% relative humidity; 48 hr incubation at 35 °C versus 48 hr incubation at 35 °C followed by 72 hr incubation at 25 °C.

## Conclusions

- Test position was found to have a significant effect upon microbial recovery. By using a proper test design with switching of air samplers among positions, this bias was eliminated from the test results.
- When higher counts are encountered, a statistically significant difference in sampler recovery is observed. In this environment, M Air T is equivalent to STA and more sensitive than RCS PLUS and RCS High Flow.
- Increased sensitivity is important, particularly if the sampler is used in a variety of testing environments or if a test area encounters spikes in microbial counts.

## Authors

Excerpted from a presentation at the 99th American Society for Microbiology General Meeting, May 30-June 3, 1999, Chicago, Illinois. Poster session 153/Q: Aerosols and Air Quality. Poster number Q-230.

Kerry Roche Lentine, Nancy Le,\* Jean Lemonnier,† Agnes Entzmann,‡ and Molly Pickett\* (Millipore Corp., Bedford, Massachusetts, United States,\* St. Quentin-en-Yvelines, France,† and Molsheim, France‡)

## Acknowledgement

The authors would like to acknowledge, with appreciation, the statistical advice of Mark Blanchard (Millipore Corp., Bedford, Massachusetts, United States)

## For Additional Information

For a complete copy of the presentation, request Millipore document TB002EN00.

\* Plate counts were transformed to the square root of the count and analyzed using a general linear ANOVA for a DOE design for all air samplers ( $\alpha$  0.05). All data were analyzed using Minitab release 12.2 (Minitab, Inc., State College, Pennsylvania).

## To Place an Order or Receive Technical Assistance

For additional information call your nearest Millipore office:

In the U.S. and Canada, call toll-free **1-800-MILLIPORE (1-800-645-5476)**

In the U.S., Canada and Puerto Rico, fax orders to **1-800-MILLIFX (1-800-645-5439)**

On the Internet

<http://www.millipore.com>

E-mail: [tech\\_service@millipore.com](mailto:tech_service@millipore.com)

## Millipore Worldwide

### AUSTRALIA

Tel. 1 800 222 111  
or (02) 9888 8999  
Fax (02) 9878 0788

### AUSTRIA, CENTRAL EUROPE, C.I.S., AFRICA, MIDDLE EAST AND GULF

Tel. (43) 1 877-8926  
Fax (43) 1 877-1654

### BALTIC COUNTRIES

Tel. +358 9 804 5110  
Fax +358 9 256 5660

### BELGIUM AND LUXEMBOURG

Tel. (02) 726 88 40  
Fax (02) 726 98 84

### BRAZIL

Tel. (011) 548-7011  
Fax (011) 548-7923

### CANADA

Tel. 1-800-645-5476  
Fax 1-800-645-5439

### CHINA, PEOPLE'S REPUBLIC OF

Beijing:  
Tel. (8610) 6500-8063  
Fax (8610) 6500-7372  
Guangzhou:  
Tel. (8620) 8755-4021  
Fax (8620) 8755-4350  
Hong Kong:  
Tel. (852) 2803-9111  
Fax (852) 2513-0313  
Shanghai:  
Tel. (8621) 5306-9100  
Fax (8621) 5306-0838

### CZECH REPUBLIC

Tel. 02-205 138 41  
02-205 138 42  
Fax 02-205 14294

### DENMARK

Tel. 70 10 00 23  
Fax 70 10 13 14

### FINLAND

Tel. (09) 804 5110  
Fax (09) 256 5660

### FRANCE

Tel. (01) 30.12.70.00  
Fax (01) 30.12.71.80

### GERMANY

Tel. (06196) 494-0  
Fax (06196) 43901

### HUNGARY

Tel. 1-205 9784  
Fax 1-205 9792

### INDIA

Tel. (91) 80-839 46 57  
Fax (91) 80-839 63 45

### ITALY

Vimodrone (Milano):  
Tel. (02) 25.07.81  
Fax (02) 26.50.324  
Roma:  
Tel. (06) 52.03.600  
Fax (06) 52.95.735

### JAPAN

Tel. (03) 5442-9711  
Fax (03) 5442-  
9736 Analytical  
9737 BioProcess  
9734 Lab Water

### KOREA

Tel. (822) 551-0990  
Fax (822) 551-0228

### MALAYSIA

Tel. 603-757 1322  
Fax 603-757 1711

### MEXICO

Tel. (525) 576-9688  
Fax (525) 576-8706

### THE NETHERLANDS

Tel. 076-5022000  
Fax 076-5022436

### NORWAY

Tel. 22 67 82 53  
Fax 22 66 04 60

### POLAND

Tel. 22-669 12 25  
22-663 70 31  
Fax 22-663 70 33

### PUERTO RICO

Tel. (787) 273-8495  
Fax (787) 747-6553

### SINGAPORE

Tel. (65) 842 1822  
Fax (65) 842 4988

### SPAIN AND PORTUGAL

Madrid:  
Tel. 917 283 960  
Fax 917 292 909  
Barcelona:  
Tel. 934 525 530  
Fax 934 516 048

### SWEDEN

Tel. 08-628 6960  
Fax 08-628 6457

### SWITZERLAND

Volketswil:  
Tel. (01) 908-30-60  
Fax (01) 908-30-80  
Lausanne:  
Tel. (021) 641 2550  
Fax (021) 641 2551

### TAIWAN

Taipei:  
Tel. (886-2) 700-1742  
Fax (886-2) 755-3267  
Hsin Chu City:  
Tel. (886-3) 571-0178  
Fax (886-3) 572-9520

### U.K. AND IRELAND

Tel. (01923) 816375  
Fax (01923) 818297

### U.S.A.

Tel. (781) 533-6000  
Fax (781) 533-3110

### IN ALL OTHER COUNTRIES

Millipore Intertech (U.S.A.)  
Tel. +1 (781) 533-8622  
Fax +1 (781) 533-8630

# MILLIPORE

Millipore and M Air T are trademarks of Millipore Corporation.

Lit. No. TB1000EN00 Printed in U.S.A. 8/03 03-255

Copyright 1999, 2003 Millipore Corporation, Billerica, MA. All rights reserved.