

Improved Recovery of Airborne Microorganisms

■ In clean areas using a portable air sampler



Abstract

A new portable air monitoring system, recently developed at Millipore Corporation, provides improved recovery of airborne microorganisms. The M Air T™ system is based on the Anderson principle and uses a sieve with about 1,000 micro-perforations, which reduces the potential for overlapping colonies and minimizes the desiccation of the medium.

The system incorporates a unique cassette design that employs a larger agar volume and depth along with a flat surface, which, in combination with a stepwise increase in air velocity, allows consistent particle size impaction. Up to 1,000 L (1 m³) of air can be sampled within seven minutes. The cassette surface rapidly rehydrates after impaction thereby renewing the water activity and reducing further effects of drying on organisms present.

The M Air T System was compared to the United States Pharmacopeia's recommended method, Slit-to-Agar (STA) sampling, and two other portable systems in two test areas. Recoveries were equivalent to the Slit-to-Agar impactor and significantly higher than the other portable systems in the test area with a significant microbial population. Recoveries were equivalent

for all samplers in the M 3.5 (Class 100) clean room. The M Air T System provides an alternative to other sampling devices for monitoring airborne microorganisms where both portability and sensitivity are needed.

Introduction

Microbiological monitoring of air is a critical component of the environmental monitoring program of any pharmaceutical, biotherapeutic, or medical device manufacturer. Air monitoring is also a Good Manufacturing Practice (GMP) requirement for both the United States and Europe.

Air monitoring provides information on the quality of the processing environment during manufacturing and enables the study of microbiological air quality trends. A critical component of air monitoring is the selection and validation of the microbiological air sampler.

The selection of a sampler typically entails comparison to an "industry standard." Such a comparison is difficult as the mechanisms of particle capture and the particle capture cutoff-size may vary among air samplers. Further, to obtain accurate results the devices must be oriented to minimize positional bias and air turbulence due to sampler exhaust.

The purpose of this study was to compare the performance of a new microbial air sampler (M Air T), to the U.S. industry standard (Slit-to-Agar or STA) and two other widely used air samplers. The samplers were tested in areas with two different levels of cleanliness.

MILLIPORE

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 m³ (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 m³) and in a Class 100 clean room (room volume 70 m³). All samplers were placed at least 1 m from the ceiling and 1 m from each other (except between positions 2 and 3 in the clean room, which were 0.76 m apart). 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. Samples from the clean room were enumerated only at the completion of the two-step incubation.

Statistical Evaluation

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 (α 0.05). An orthogonal pairwise analysis was performed on: M Air T versus STA, RCS High Flow versus RCS PLUS, and grouped M Air T/STA versus grouped RCS High Flow/RCS Plus. The results were also confirmed by a non-parametric analysis using the ranks of the square root of the counts (rank of the count). In addition, a prevalence analysis was performed on the clean room data. In this case, all zero cfu results were coded as "0" and all positive cfu results were coded as "1." All data were analyzed using Minitab release 12.2 (Minitab, Inc., State College, Pennsylvania).

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 ³)	6.5 min (140 L/min 1 st 500 L, 180 L/min 2 nd 500 L)	20 min (50 L/min)	10 min (100 L/min)	20 min (50 L/min)
Water loss after sampling 1 m ³ *	4.2%	4.3%	10%	13%
Particle diameter cutoff size (d ₅₀)	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 at both incubation enumeration points and a significant difference among test positions. The P-values for these analyses are shown in Tables 2 and 3.

Overall results for the clean room showed a significant difference in test position but no significant difference in recovery among air samplers. Because of the high number of zero cfu results, the rank of the square root of the counts (rank of the counts) was used in the clean room analysis. The orthogonal analysis was performed on the final time-point counts for the clean room. As expected, the clean room results indicated no significant difference among pairs of samplers (Table 4 and Figure 1).

The overall analysis for the unclassified room suggested a significant difference in recovery for certain pairs of samplers. An orthogonal 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 2).

Table 2.

P-Values for DOE Overall Analysis

	Unclassified Room 48 hr at 35 °C SQRT* of count (n=479)	Unclassified Room 72 hr at 25 °C SQRT of count (n=478)	Clean Room Rank of count (n=480)
P-value tester	0.000	0.000	0.628
P-value position	0.032	0.035	0.000
MSE	1.385	1.499	10778

*SQRT denotes square root.

Table 3.

P-Values for Orthogonal Analysis

	M Air T vs. STA	High Flow vs. PLUS	M Air T/STA vs. High Flow/PLUS
Clean Room (rank of count)			
Sample size	240	240	480
P-value tester	0.243	0.598	0.714
P-value position	0.033	0.004	0.000
MSE	2675	2586	11049
Unclassified room, 48 hr at 35 °C (square root of count)			
Sample size	239	240	479
P-value tester	0.772	0.675	0.000
P-value position	0.078	0.250	0.032
MSE	1.495	1.359	1.388
Unclassified Room, 72 hr at 25 °C (square root of count)			
Sample size	239	239	478
P-value tester	0.166	0.552	0.000
P-value position	0.043	0.282	0.035
MSE	1.848	1.190	1.507

Table 4.

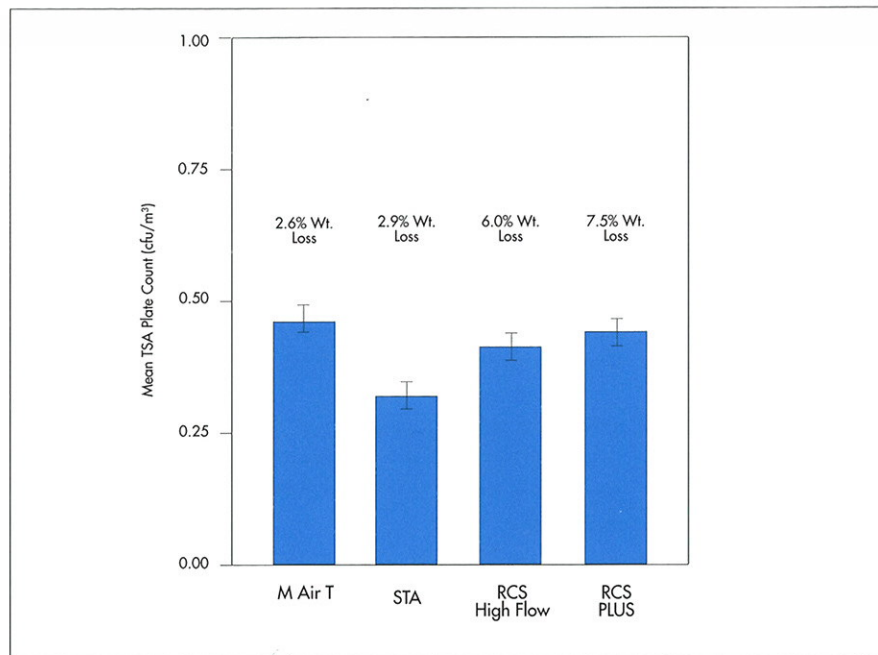
Prevalence of Positive Results for Three Microbial Air Samplers Against the Industry Standard Slit-to-Agar in a Clean Room Environment (Class M3.5)*

	STA +	STA -	Total
M Air T +	10	26	36
M Air T -	19	65	84
Total	29	91	
RCS PLUS +	13	23	36
RCS PLUS -	16	68	84
Total	29	91	
RCS High Flow +	5	27	32
RCS High Flow -	24	64	88
Total	29	91	

*Sampler recovery agar with colony forming units was scored as a positive. Sampler recovery agar without growth was scored as a negative.

Figure 1.

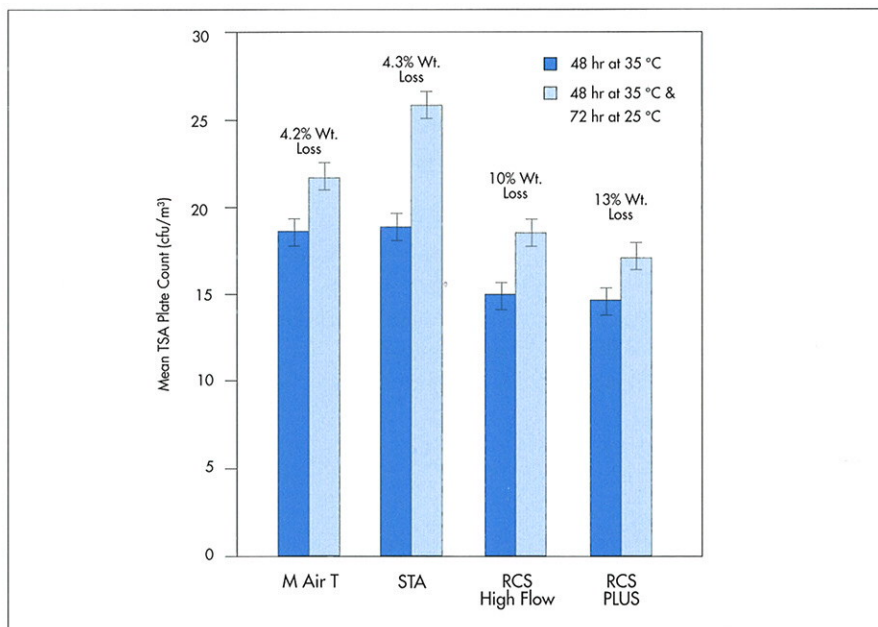
Mean Recovery and Standard Error for Four Different Microbial Air Samplers Operated Simultaneously in a Clean Room (M 3.5).



Conditions: 53% relative humidity; 48 hr incubation at 35 °C followed by 72 hr incubation at 25 °C; 120 samples per sampler.

Figure 2.

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.

Discussion

A significant increase in plate counts was observed during the two-stage incubation of 48 hour at 35 °C followed by 72 hour at 25 °C for the unclassified area. Since there were so few colonies recovered in the clean room, the plates were only counted at the end of the two-stage incubation.

The percent weight loss 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. However, this relationship was not established in the clean room environment. While the RCS PLUS and the RCS High Flow had higher percent weight loss, the four samplers were not significantly different for recovery.

The lack of relationship between percent weight loss and recovery in the clean room is due to the low incidence of contamination. The mean untransformed plate count for the samplers in the clean room ranged from 0.32 to 0.45 cfu (n=120 per sampler). The mean final incubation plate count (untransformed) in the unclassified room ranged from 18 to 26 cfu. The clean room trials resulted in a significant amount of zero cfu results. The low counts, coupled with the zero cfu results, may not enable the prediction of differences in recovery among samplers or establish a relationship between recovery and percent weight loss in the clean room environment.

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.
- The low incidence of microbial colonies in the clean room study contributed to no significant difference among samplers in this environment.
- 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

Presented 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)

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

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 69 60
Fax 08-628 64 57

SWITZERLAND

Volketswil:
Tel. (01) 908-30-60
Fax (01) 908-30-80
Lausanne:
Tel. (021) 648 51 18
Fax (021) 648 51 37

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 or an affiliated company
Lit. No. TB002EN00 Printed in U.S.A. 8/99
99-218

Copyright 1999 Millipore Corporation or an affiliated company, Bedford, MA. All rights reserved.

♻️ Printed with soy inks on recycled paper.