

QC Micro Webinar Series

Lonza

Amylin's Improved Aseptic Gowning and Monitoring Program Using MODA™ Solution

Today's Presenters

- Bob Toal
Segment Manager, Informatics
Lonza Wayne
- Dominick Villani
Manager of QC Microbiology
Amylin Pharmaceuticals

60-Minute Agenda

- Lonza Overview
- 45-min Presentation
- 15-min Interactive Q&A
- Upcoming News/Events
- Wrap-up

A copy of this presentation will be made available through a follow up email.

Webinar Focus Areas

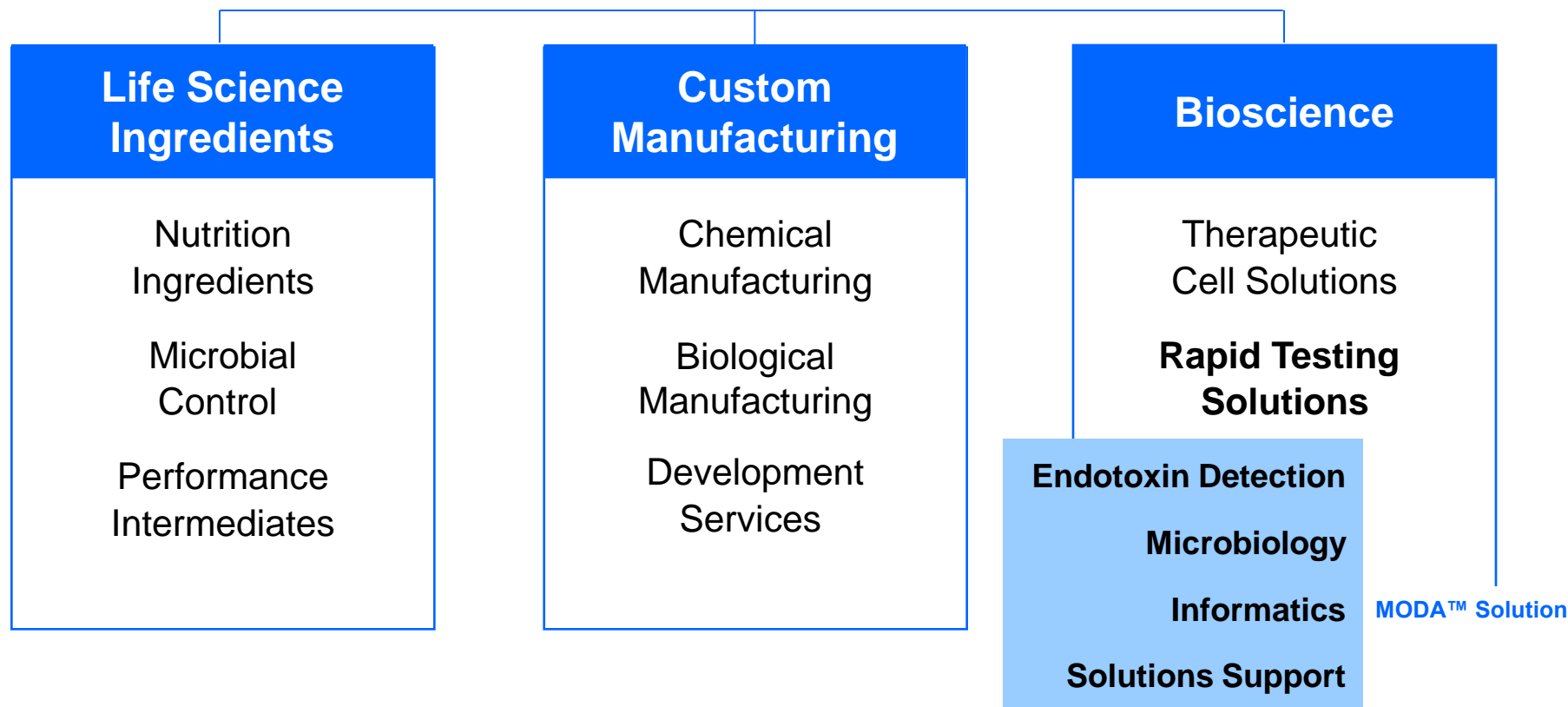
- Industry best practices, strategies, and regulatory guidelines
- Definitions of qualification, re-qualification and routine monitoring
- Use of automation and the MODA™ Solution to meet requirements
- Tracking, trending, and report generation compliance using MODA™ Solution



Lonza and MODA™ Solution Overview

Lonza's Life-Science Platform

Lonza



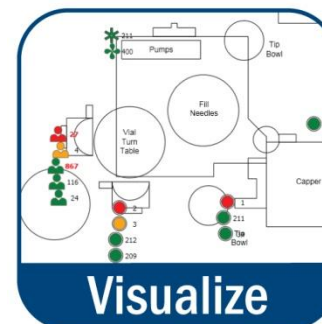
MODA™ Solution Value Proposition

■ More science. Less paper.™

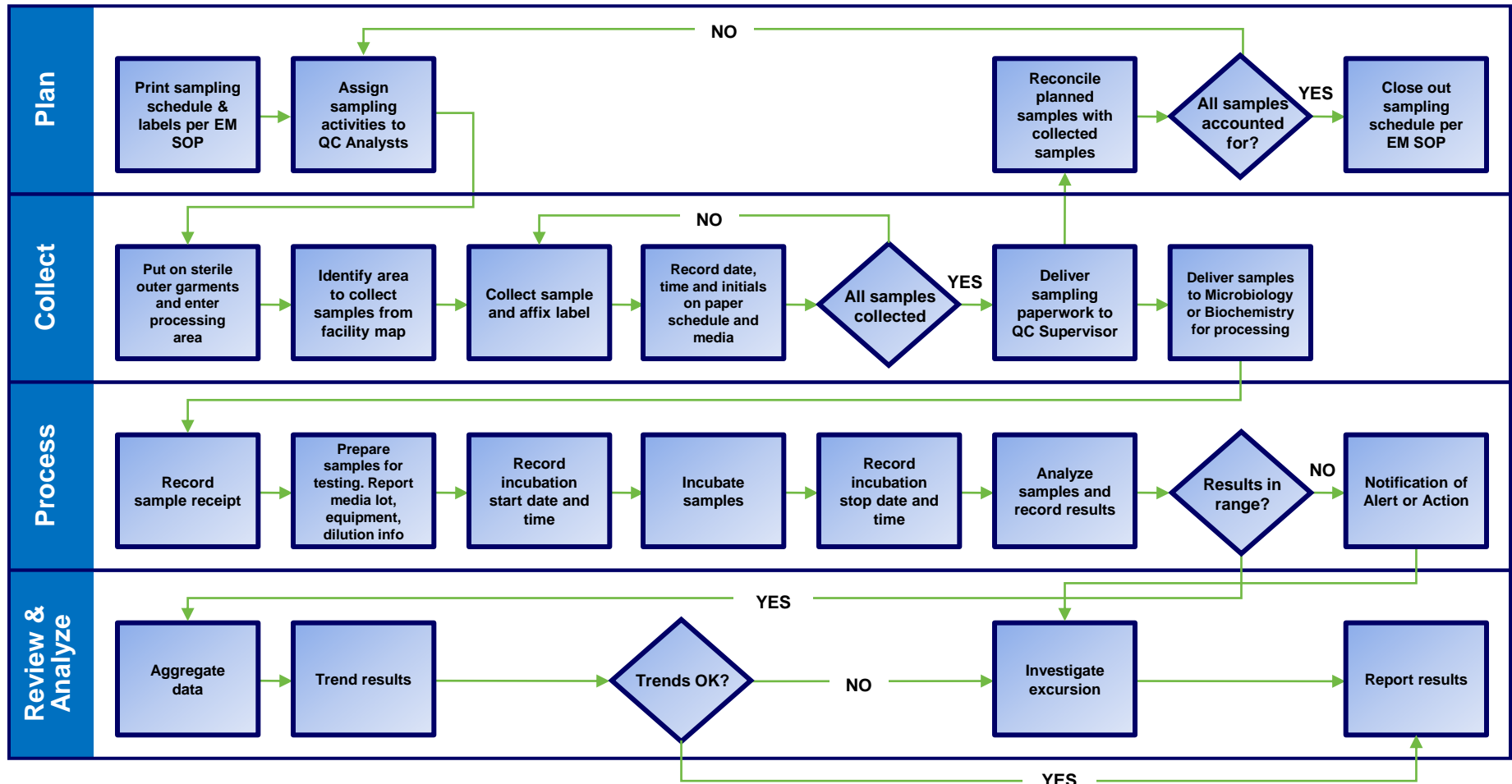
- Quickly move from paper-intensive QC Monitoring & Analysis

■ Increase operational efficiency, improve quality, reduce costs

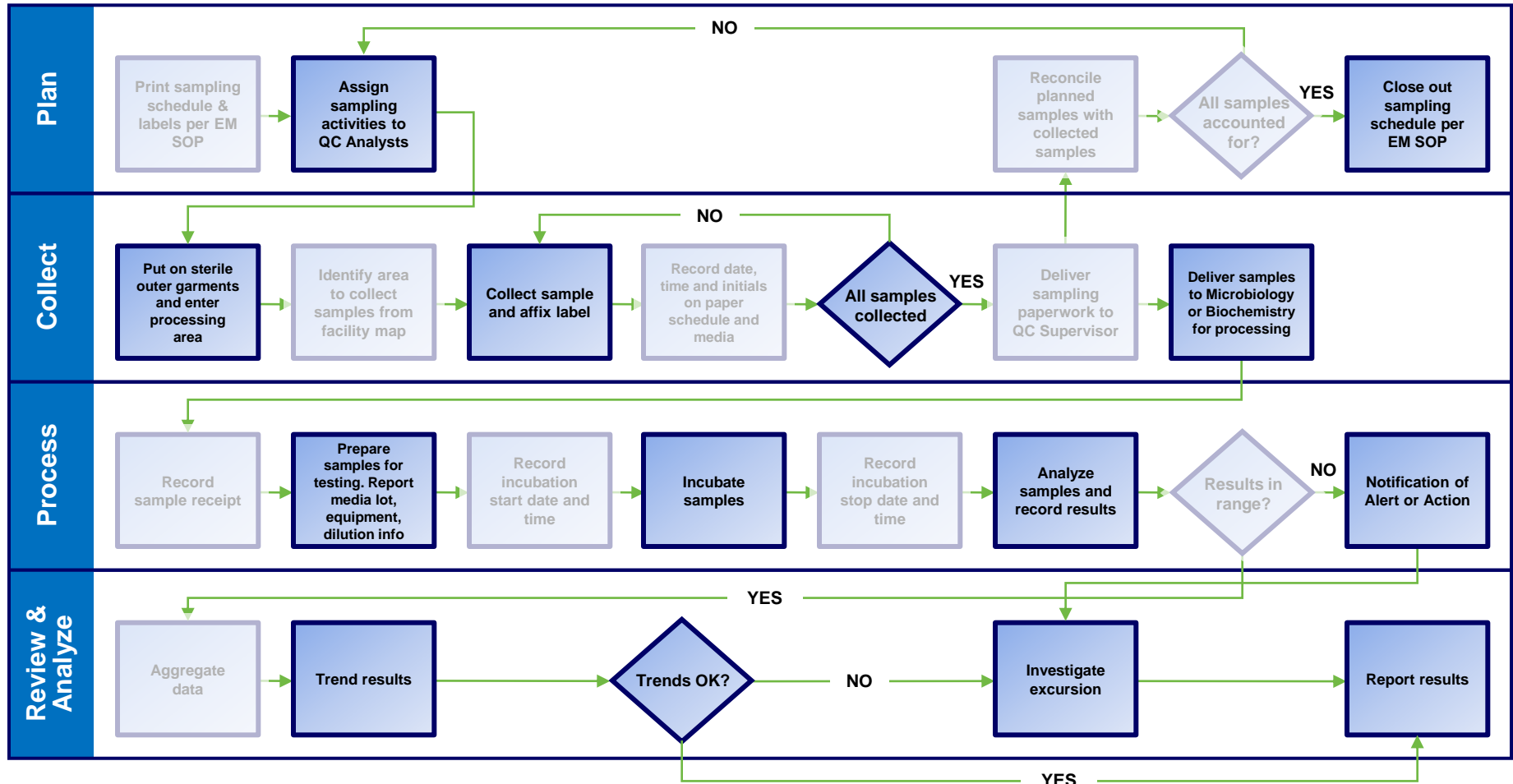
- MODA-EM™ platform offers mobile computing technology & advanced visualization tools






The Paper-based QC Process



The **Paperless** QC Process—11 Steps Removed



Paperless Efficiency Example

Record on paper template	Record on plate	Reconcile media and template	Enter data into Excel/LIMS
1: BB50, SS43, Water, IE5 2: BB50, SS43, RODAC, IP72/IP82 3: BB50, SS44, Water, IE5/IE7 4: BB50, SS44, Water, IE9/SP2 5: BB50, SS65, RODAC, IP74			
6 hours		1 hour	1 hour

Scan room or site	Select sample and print labels	Send to repository
		
4 hours		Finished

	Paper	8 hours
–	Paperless	4 Hours
=	Savings	4 Hours

Automated Field Data Capture



Featured: MODA-FDC™ platform, used for field data capture in clean room areas

- Stainless steel cart
- Ergonomic tablet PC
- Docking station
- Thermal label printer
- Barcode scanner gun
- Proximity reader for RF badges
- Space for equipment
- Space for growth media

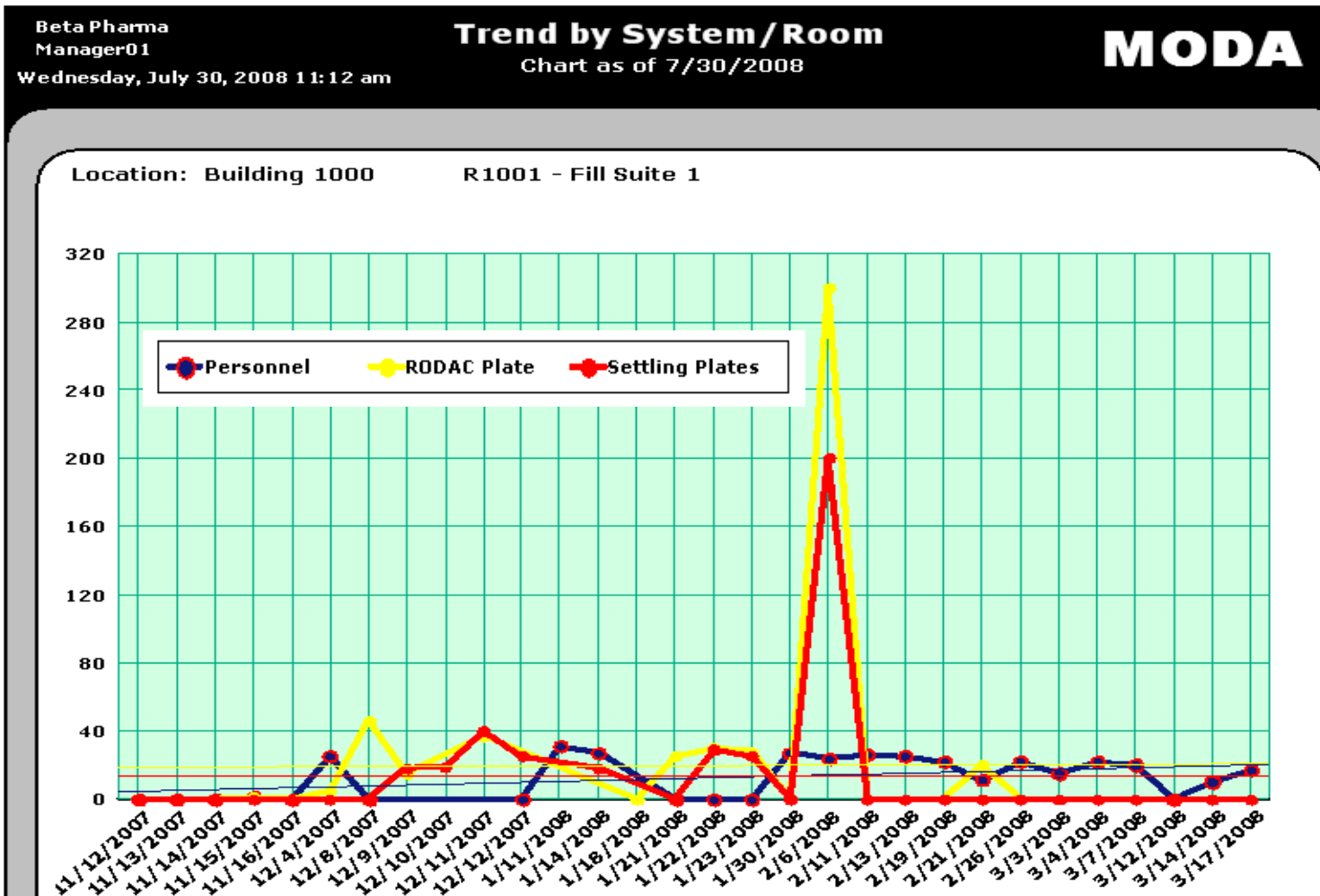




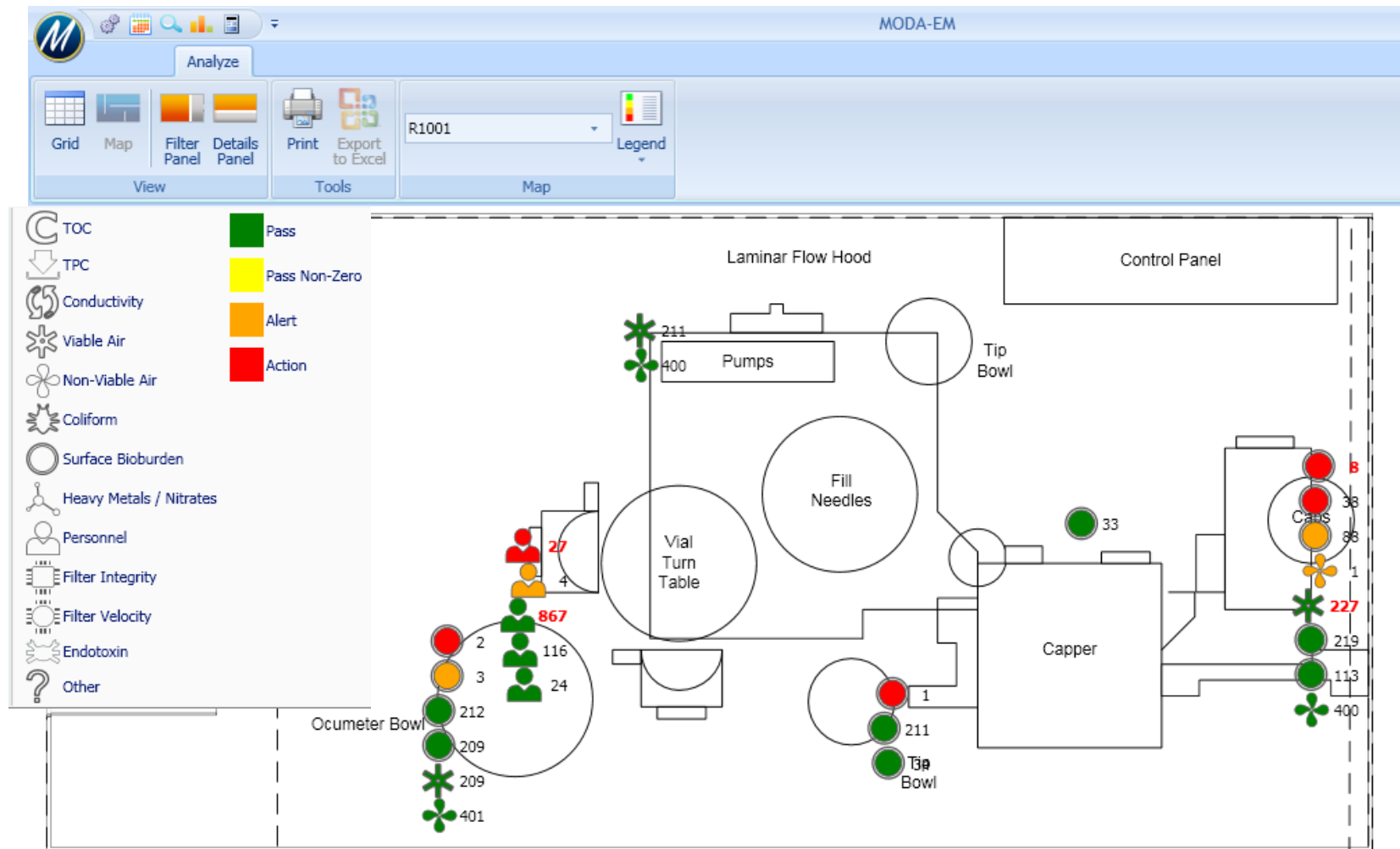
Automated Reporting and Analytics

<div> <div>Beta Pharma</div> <div>Manager01</div> <div>7/30/2008 11:04:13AM</div> </div> <div> <div>Deviation Summary</div> <div>Report Period From 11/1/2007 12:00:00AM to 3/31/2008 12:00:00AM</div> <div>MODA</div> </div>										
Deviation	Sample	Sample Date	Sampled By	Test	Result	Alert Limit	Action Limit	Organism	Deviation Type	Environment
Building 1000 > R1001-Fill Suite 1					Alerts:	6	Actions:	50	Total:	52
169	A00000040	12/9/2007 6:16:59PM	Analyst08	R1001.1 - Settling Plates	18.00cfu	1.00	3.00		>= Action	D
174	A00000093	12/12/2007 6:38:00AM	Analyst08	R1001.1 - Settling Plates	25.00cfu	1.00	3.00	Bacillus licheniformis; Escherichia coli	>= Action	D
178	M00000106	1/14/2008 11:41:16AM	Manager01	R1001.1 - Settling Plates	18.00cfu	1.00	3.00		>= Action	D
182	M00000263	1/23/2008 10:56:52AM	Manager01	R1001.1 - Settling Plates	25.00cfu	1.00	3.00	Escherichia coli	>= Action	D
182	M00000263	1/23/2008 10:56:52AM	Manager01	R1001.1 - Settling Plates	25.00cfu	1.00	3.00	Escherichia coli	>= Action	D
185	M00000036	12/11/2007 9:05:29AM	Manager01	R1001.1 - Settling Plates	40.00cfu	1.00	3.00		>= Action	D
187	M00000239	1/22/2008 9:02:17AM	Analyst08	R1001.1 - Settling Plates	29.00cfu	1.00	3.00		>= Action	D
198	M00000298	2/6/2008 12:08:14PM	Analyst08	R1001.1 - Settling Plates	200.00cfu	1.00	3.00		>= Action	D
SWAB-Surface Swab					Alerts:	0	Actions:	1	Total:	1
189	M00000273	1/30/2008 9:12:48AM	Analyst08	R1001.3 - Surface Swab	27.00cfu	1.00	3.00		>= Action	D
Building 1000 > R1004-Fill Suite 4					Alerts:	0	Actions:	4	Total:	4
PERSON-Personnel					Alerts:	0	Actions:	4	Total:	4
151	R1419009301	11/6/2007 10:32:35PM	Analyst04	R1004.6 - Personnel	6.00cfu	2.00	3.00		>= Action	D
152	R1419009302	11/6/2007 10:32:35PM	Analyst04	R1004.6 - Personnel	6.00cfu	2.00	3.00		>= Action	D
153	R1419009303	11/6/2007 10:32:35PM	Analyst04	R1004.6 - Personnel	6.00cfu	2.00	3.00		>= Action	D
154	R1419009304	11/6/2007 10:32:36PM	Analyst04	R1004.6 - Personnel	6.00cfu	2.00	3.00		>= Action	D

Automated Reporting and Analytics



Automated Reporting and Analytics



Lonza

Our Featured Speaker:
Dom Villani – Amylin Pharmaceuticals

USING MODA™ SOLUTION TO IMPLEMENT AND MAINTAIN A ROBUST AND COMPLIANT ASEPTIC GOWNING QUALIFICATION, RE-QUALIFICATION, AND ROUTINE MONITORING PROGRAM

Presenter: Dominick C. Villani III
07JUN2011

AMYLIN PHARMACEUTICALS

- ◉ Amylin produces sterile products through aseptic processing
- ◉ Amylin implemented and qualified MODA™ in Q1 2009
- ◉ Amylin utilizes MODA™ to track, monitor, and manage all sample plans, sample frequencies, and excursions for classified cleanrooms and critical utilities (i.e. WFI, Clean Steam, Compressed Gases, etc.)
- ◉ Amylin currently uses MODA™ 2.3 and is in the process of upgrading to MODA™ 3.0

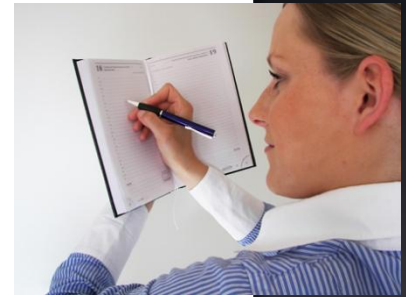
PRESENTER BIOGRAPHY

Dominick Villani

Manager of Quality Control Microbiology at Amylin OH LLC

- ◉ Current responsibilities: Laboratory method development and validation; environmental and utilities monitoring qualifications and routing maintenance programs; and overall Sterility Assurance Programs
- ◉ 13+ years experience with GMP manufacturing operations, QC microbiology activities, and QA oversight of GMP facilities at Amgen, Bioport Corporation, Bristol-Myers Squibb, and Amylin
 - 4 years dedicated to the development and management of EM programs, including aseptic gowning, and sterility assurance programs for sterile products
- ◉ B.S. in Biological Sciences, with an emphasis in Microbiology

AGENDA AND OBJECTIVES



- ◉ Regulatory requirements for a compliant aseptic gowning monitoring program
- ◉ Industry best practices, guidances and standards for a compliant aseptic gowning monitoring program
- ◉ Definitions of qualification, re-qualification, and routine monitoring for aseptic gowning
- ◉ Requirements and guidelines for investigations, immediate actions, and corrective actions
- ◉ MODA™ Solution advantages and solutions for implementing and managing a compliant aseptic gowning monitoring program

ASEPTIC GOWNING INTRODUCTION

- ◎ The quality of a company's product(s) can be directly defined and assessed based on:
 - Robustness of its aseptic gowning program
 - Data produced from that program
- ◎ Good aseptic gowning results signify that:
 - The company and its employees are committed to quality
 - The site and process are in a state of control



ASEPTIC GOWNING INTRODUCTION

- ◎ Aseptic gowning is by far the most important and critical operation for manufacturers of sterile products using aseptic operations.

The program must be:

- Intensive, robust, and rigid, with defined actions and consequences
- Balanced to manage business and compliance expectations
- Evolutionary and current
- Timely



ASEPTIC GOWNING INTRODUCTION

- ◎ The aseptic gowning program must be monitored, managed, and updated in the same way critical utility systems (e.g. WFI) are treated.
 - Sample plans
 - Sample frequencies
 - Sampling methods
 - Timely reporting
 - Data trending



ASEPTIC GOWNING INTRODUCTION

- ◎ Who defines the requirements and components of an aseptic gowning program?
 - Regulatory bodies
 - Industry best practices, guidances, and standards
 - Company management
- ◎ Who defends the requirements and components of an aseptic gowning program?
 - Company management
- ◎ Who is responsible for the requirements, components, and data of program to inspectors?
 - Company management





REGULATORY REQUIREMENTS

◎ Regulatory requirements:

- Legal requirements
- Based on the drug market
- Are minimum requirements



REGULATORY REQUIREMENTS

United States

**21CFR 210 & 211,
“cGMP in
Manufacturing,
Processing, Packing, or
Holding of Drugs and
Finished
Pharmaceuticals”**
•21CFR 211.25 (a) -
Personnel Qualifications
•21CFR 211.28 (a), (b),
and (d) - Personnel
Responsibilities
•21CFR 211.113 (b) -
Control of
Microbiological
Contamination

European Union

**Eudralex Volume 4,
“Good Manufacturing
Practice Medicinal
Products for Human
and Veterinary Use”**
•Chapter 2, “Personnel”
 •2.10, “Training”
 •2.13, 2.15, 2.18,
 “Personnel
 Hygiene”
•Chapter 5,
“Production”
 •5.18 “Prevention of
 Cross-
 Contamination in
 Production”



REGULATORY REQUIREMENTS

Japan

Japanese GMPs, “GMP Guideline for Drugs and Quasi-Drugs (Drug Products) 2005”

- Section 3, “Personnel”
 - 3.10, “Personnel Qualifications”
 - 3.20, 3.22, “Education and Training”
 - 3.30, 3.31, 3.3, “Personnel Hygiene Control”
- Section 7, “Production and In-Process Control”
 - 7.6, “Microbiological Contamination Control”

Harmonization

ICH Harmonised Tripartite Guideline Q7, “Good Manufacturing Practice Guide for Active Pharmaceutical Ingredients”

- Section 3, “Personnel”
 - 3.10, 3.12, “Personnel Qualifications”
 - 3.20, 3.21, 3.24, “Personnel Hygiene”



REGULATORY REQUIREMENTS

◎ Common themes with most observations are:

- Sample frequency
- Sample site rationale
- Sample techniques
- Data tracking and trending
- Employee training



REGULATORY REQUIREMENTS

- ◎ What are the consequences for not implementing and maintaining a robust and compliant aseptic gowning qualification, re-qualification, and routine monitoring program?
 - Observations
 - Warning letters
 - Recalls
 - Consent decrees

INDUSTRY BEST PRACTICES, GUIDANCES AND STANDARDS

- ◉ Industry best practices that are applicable to implementing and managing a compliant aseptic gowning program are:
 - FDA Guidance for Industry section V, “Sterile Drug Products Produced by Aseptic Processing - Current Good Manufacturing Practice”
 - PDA Journal of Pharmaceutical Science and Technology, Technical Report No.13, “Fundamentals of an Environmental Monitoring Program”
 - ISPE, Pharmaceutical Engineering Guides for New and Renovated Facilities, Volume 3, “Sterile Manufacturing Facilities”
 - International Standard (ISO) 13408-1
 - United States Pharmacopeia <1116>



INDUSTRY BEST PRACTICES, GUIDANCES AND STANDARDS

◉ Industry best practices and standards highlight:

- Training
- Personal hygiene and health habits
- Gowning control
- Facilities control
- Aseptic gowning monitoring
- Tracking and trending



ASEPTIC GOWNING DEFINITIONS OF QUALIFICATION, RE-QUALIFICATION, AND ROUTINE MONITORING PROGRAM

◎ Major phases of the program should include:

- Qualification
- Routine monitoring
- Re-qualification



ASEPTIC GOWNING DEFINITIONS OF QUALIFICATION, RE-QUALIFICATION, AND ROUTINE MONITORING PROGRAM

- Routine monitoring activities should include:
 - Routine monitoring
 - Audits



ASEPTIC GOWNING DEFINITIONS OF QUALIFICATION, RE-QUALIFICATION, AND ROUTINE MONITORING PROGRAM

◎ Batch critical interventions

- Sample locations
- Frequency of sampling
- Limits
- Microbiological assessments
- Actions and consequences
- Security



ASEPTIC GOWNING DEFINITIONS OF QUALIFICATION, RE-QUALIFICATION, AND ROUTINE MONITORING PROGRAM

◎ Audits

- Sample locations
- Frequency of sampling
- Limits
- Microbiological assessments
- Actions and consequences
- Security



ASEPTIC GOWNING DEFINITIONS OF QUALIFICATION, RE-QUALIFICATION, AND ROUTINE MONITORING PROGRAM

◎ Re-qualification activities should include:

- Sample locations
- Frequency of sampling
- Limits
- Microbiological assessments
- Actions and consequences
- Security



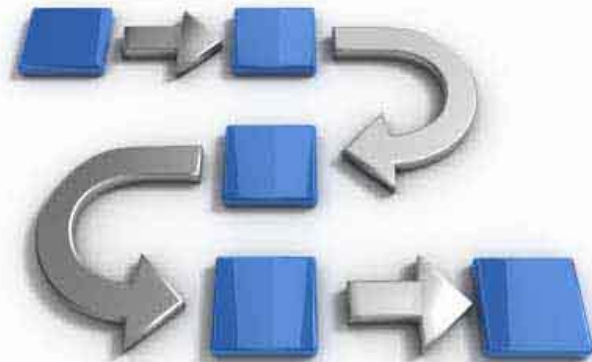
RATIONALE FOR SAMPLE LOCATIONS, SAMPLE FREQUENCY, SAMPLE LIMITS

◎ Definitions of particular importance:

- Alert
- Action
- % Positives

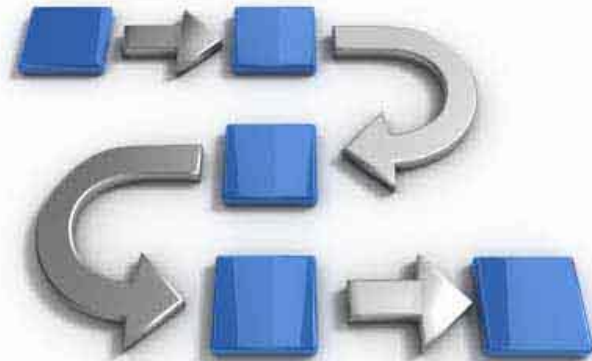
◎ Predecessors to defining locations, frequencies, and limits:

- Sample methods
- Media type
- Equipment type



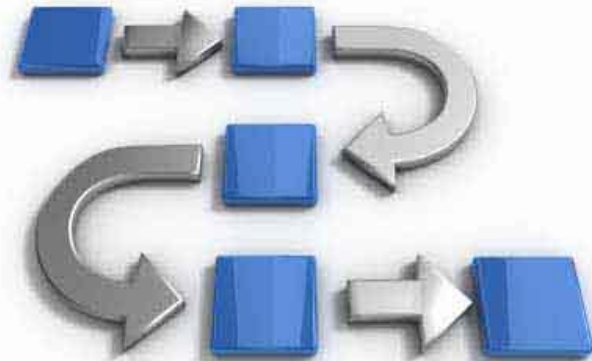
RATIONALE FOR SAMPLE LOCATIONS, SAMPLE FREQUENCY, SAMPLE LIMITS

- ◉ Sample location selection is based on the following factors:
 - Likelihood for contamination
 - Likelihood for contamination transfer
 - Gown attachment points



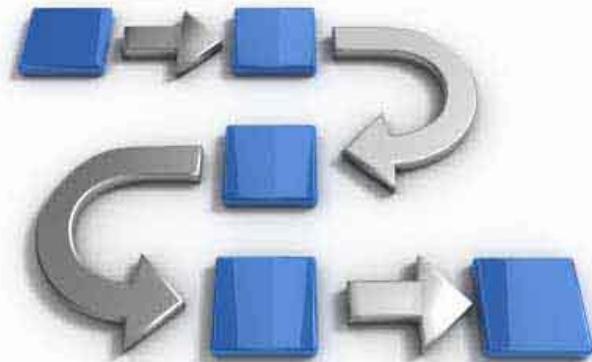
RATIONALE FOR SAMPLE LOCATIONS, SAMPLE FREQUENCY, SAMPLE LIMITS

- ◎ Sample frequencies should be based on:
 - Criticality of tasks
 - Frequency of tasks
 - Trends



RATIONALE FOR SAMPLE LOCATIONS, SAMPLE FREQUENCY, SAMPLE LIMITS

- ◎ Sample limits should be defined and justified:
 - Absence of objectionable microorganisms
 - Alert
 - Action
 - % Positives
 - Total body count



INVESTIGATIONS, IMMEDIATE ACTIONS, AND CORRECTIVE ACTIONS



- ◎ Investigation types include:
 - Alerts
 - Actions
 - % Positives
 - Total body count
- ◎ Investigations into aseptic gowning excursions should include:
 - Immediate response
 - Timely closure
 - Approved by all key stakeholders

INVESTIGATIONS, IMMEDIATE ACTIONS, AND CORRECTIVE ACTIONS



- ◎ Investigations for personnel excursions can be:
 - Difficult and challenging
 - Subjective
 - Adversarial
- ◎ Use the Program
 - Sample locations
 - Frequencies
 - Locations
 - Isolate information

INVESTIGATIONS, IMMEDIATE ACTIONS, AND CORRECTIVE ACTIONS



◎ Immediate actions:

- Actions implemented to prevent nonconforming events from having or continuing to have impact to the process or product

◎ Immediate actions should be:

- Timely
- Diligent
- Be based on a point system

INVESTIGATIONS, IMMEDIATE ACTIONS, AND CORRECTIVE ACTIONS



◎ Corrective actions:

- Actions taken to prevent recurrence of the defined root cause
- A corrective action will always relate to the determined root cause

◎ Preventive Actions:

- Actions taken to prevent the occurrence of an issue

INVESTIGATIONS, IMMEDIATE ACTIONS, AND CORRECTIVE ACTIONS

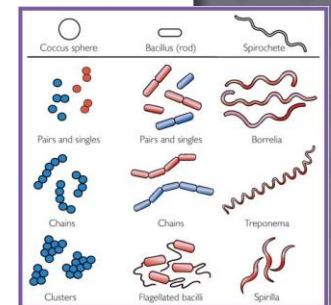


- ◎ Incomplete investigations and root cause determinations may lead to:
 - Implementation of ineffective CAPAs
 - Recurring excursions
 - Compliance vulnerabilities
 - Negative impact to the product and processes

OBJECTIONABLE MICROORGANISMS

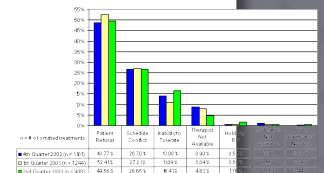
◉ Know your micro flora:

- Frequent microbial identifications
- Define objectionable microorganisms
- When to act and when not to



TRACKING AND TRENDING CRITICAL DATA FOR ASEPTIC GOWNING

- The tracking and trending system must be qualified
- Examples of data that should be tracked and reported include:
 - Excursion rate by area or by person
 - % Positive rate by area or by person
 - Most commonly recovered isolates by area, by person, and by quarter





HOW DOES THE MODA™ SOLUTION
AID IN IMPLEMENTING AND
MAINTAINING A GOOD ASEPTIC
MONITORING PROGRAM?

KEY FUNCTIONALITY FOR AMYLIN

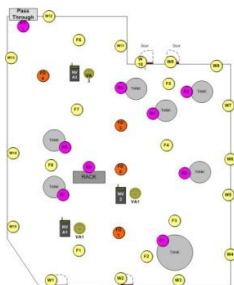


- ◎ **MODA™ Solution provides an automated and compliant solution for:**
 - Defining, Controlling, and Managing Sample Plans
 - Defining, Controlling, and Managing Sample Frequencies
 - Defining and Controlling Security Roles and Rights
 - Forcing and maintaining structured and consistent work flows for samples. Sample Chain of Custody is controlled!
 - Tracking and Controlling Materials, Equipment, Results, and Personnel
 - Assistance in Investigating Aseptic Gowning Excursions
 - Tracking and Trending Critical Data points specific to:
 - Qualifications
 - Batch Monitoring
 - Audits
 - Re-Qualifications

QC PROCESS CHALLENGES

Disparate Data Points

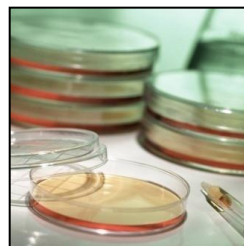
**Air, Surface,
Water**



Personnel



Media



Equipment



“Islands” of Information

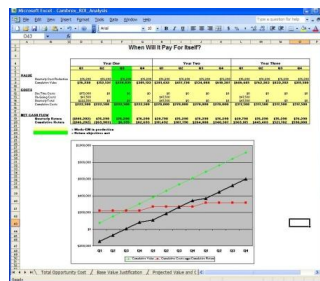
Sample Records



Laboratory



Spreadsheets



Management





WHY IT MATTERS

- ◉ Paper-based, redundant data recording & reconciliation
- ◉ Increased labor and time costs & delays in data analysis/reporting
- ◉ Ineffective data analysis and trending
- ◉ Impacts overall effectiveness of QC program

AUTOMATION IN CLEANROOM



“...cleanroom touch pads or computer terminals that allow for automated data entry IN THE ROOM.”

“...palm-pilot-type of data collection devices... that can directly download to the computer system and allow for direct data transfer without risk of contamination.”

“...real time data for many of the chemistry and microbiology tests that must be performed.”



Source: [Environmental Monitoring A Comprehensive Handbook, Volume 1](#)

AUTOMATION IN QC MICRO LAB



“...analysis and trending of environmental data is essential to aid in the interpretation of process stability and assess overall control performance.”

EM Reports must be “...accurate, traceable, timely, and well-documented.”



Source: [Environmental Monitoring A Comprehensive Handbook, Volume 1](#)



21 CFR PART 11 COMPLIANCE

Paper-Based System

Access database for data storage is not validated

No control of changes in access database

No electronic signature for changes

All users have all privileges

Data stored in binders that are stored on and off site

MODA™ Solution

- ✓ Validated system
- ✓ Audit trail tracks all changes/records
- ✓ Electronic signature for all major steps to enable quick and clear traceability
- ✓ Varying levels of access to system dependent upon job function
- ✓ Data stored on servers that are backed up on a routine basis

TRENDING/REPORT GENERATION COMPLIANCE



Paper-Based System

5 Analysts 8-10 hours per month
after data entry

Limited Scope of Trending
Reports

No Formal way of tracking flora

Difficult to extract client-
specific data

All trending performed by QC
for QA and Manufacturing

MODA™ Solution

- ✓ Quick and efficient trending in real time
- ✓ Wide array of trend report formats
- ✓ Flora can be trended by person, site, room, facility
- ✓ Client-specific reporting for Contract Manufacturing
- ✓ Ease of trending allows others (QA, Manufacturing) to perform their own trending

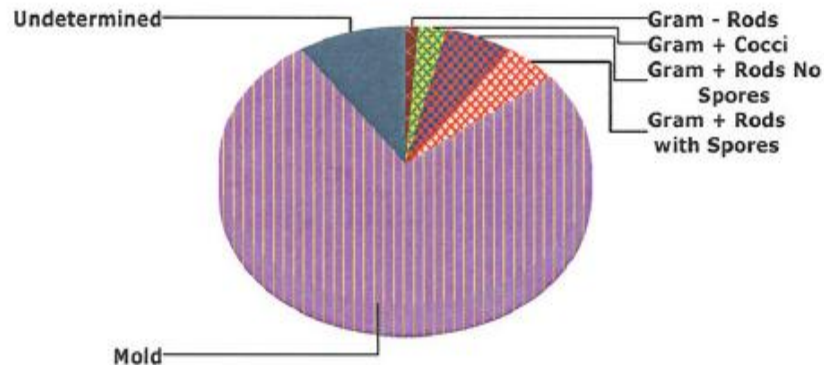
FROM THIS.....



TO THIS.....

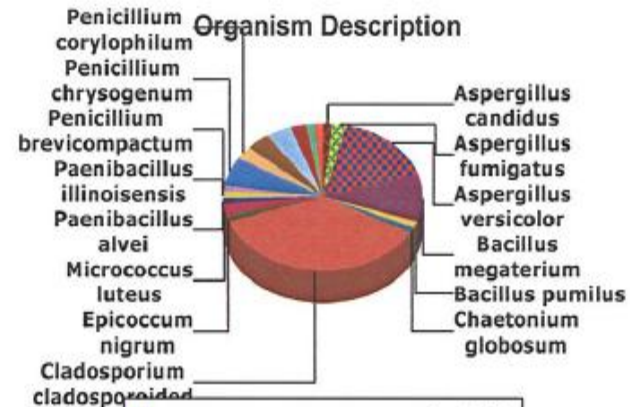


Organism Characterization



+	0	0.0%
Gram - Rods	1	1.2%
Gram + Cocci	2	2.4%
Gram + Rods No Spores	5	5.9%
Gram + Rods with Spores	4	4.7%
Mold	65	76.5%
Undetermined	8	9.4%
Total:	85	100.0%

Organism Description



+	0	0.0%
Aspergillus candidus	1	1.3%
Aspergillus fumigatus	2	2.7%
Aspergillus versicolor	12	16.0%
Bacillus megaterium	8	10.7%
Bacillus pumilus	1	1.3%
Chaetonium globosum	1	1.3%
Cladosporium cladosporioides	27	36.0%
Epicoccum nigrum	1	1.3%
Micrococcus luteus	2	2.7%
Paenibacillus alvei	1	1.3%
Paenibacillus illinoisensis	1	1.3%
Penicillium brevicompactum	1	1.3%
Penicillium chrysogenum	5	6.7%
Penicillium corylophilum	2	2.7%
Penicillium glabrum	3	4.0%
Rhodococcus equi	3	4.0%
Staphylococcus hominis hominis	2	2.7%
Staphylococcus warneri	1	1.3%
Stereum hirsutum	1	1.3%
Total:	75	100.0%



AND THIS



MODA-EM

Review/Approve

Filter Panel Details Panel View

Edit Approve Print Operation

Drag a column header here to group by that column.

	Sampling Location	Sampling Plan	Sampling Site	Sampling Barco	Test Method	Result	Deviation Type	Sampling Date
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	012 - Cellular 124	124-B	A00002502	RODAC - SDA	1	ACTION	9/15/2008 10:55: d
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 115 - RNA Proce	030 - RNA 115 - I	115-A	A00002392	Personnel - Glove	1	ACTION	9/11/2008 12:04: y
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 123 - Cellular Pr	001 - Cellular 123	123-B	B00002154	Personnel - Glove	1	ACTION	9/11/2008 10:23: jk
<input checked="" type="checkbox"/>	Argos Therapeutics -> RCC -> 123 - Cellular Pr	001 - Cellular 123	123-A	A00002316	Viable Air - TSA	1	ACTION	8/29/2008 9:30:0 r
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 123 - Cellular Pr	001 - Cellular 123	123-E	A00002307	RODAC - TSA	6	ACTION	8/29/2008 9:23:0 r
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 123 - Cellular Pr	001 - Cellular 123	123-F	A00002308	RODAC - TSA	11	ACTION	8/29/2008 9:23:0 r
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-A	A00002205	Personnel - Glove	1	ACTION	8/20/2008 10:19: jk
<input checked="" type="checkbox"/>	Argos Therapeutics -> RCC -> 124B - Cellular I	010 - Cellular 124	124B-E	A00002184	RODAC - TSA	22	ACTION	8/20/2008 9:19:0 jk
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-C	A00002174	RODAC - TSA	100	ACTION	8/20/2008 9:12:0 jk
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-E	A00002176	RODAC - TSA	100	ACTION	8/20/2008 9:12:0 jk
<input checked="" type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-F	A00002177	RODAC - TSA	100	ACTION	8/20/2008 9:12:0 jk
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124B - Cellular I	010 - Cellular 124	124B-E	B00001707	RODAC - TSA	6	ACTION	8/19/2008 10:30: r
<input type="checkbox"/>	Argos Therapeutics -> HIV -> 183 - HIV Suite 3	100 - HIV 183 In-	183-B	M00001529	Personnel - Glove	1	ACTION	8/19/2008 10:29: jk
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-C	B00001697R	RODAC - TSA	100	ACTION	8/19/2008 10:00: r
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-E	B00001699R	RODAC - TSA	100	ACTION	8/19/2008 10:00: r
<input checked="" type="checkbox"/>	Argos Therapeutics -> HIV -> 172 - HIV QC La	110 - HIV QC Lab	172-A	A00002165	Personnel - Glove	1	ACTION	8/15/2008 1:41:0 p
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 112 - RNA Proce	020 - RNA 112 - I	112-A	A00002141	Personnel - Glove	3	ACTION	8/14/2008 11:56: jk
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 112 - RNA Proce	020 - RNA 112 - I	112-A	A00002102	Viable Air - TSA	1	ACTION	8/14/2008 8:54:0 jk
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-F	B00001611	Viable Air - TC Co	10	ACTION	8/12/2008 9:08:0 r
<input type="checkbox"/>	Argos Therapeutics -> RCC -> 124 - Cellular Pr	010 - Cellular 124	124-B	B00001548	Personnel - Glove	4	ACTION	8/11/2008 10:01: r

Barcode: A00002316 Sampling Technician: mnorris Sampling Date: 8/29/2008
Location: Argos Therapeutics -> RCC -> 123 - Cellular Processing Room Test Method: Viable Air - TSA

Additional Summary Details:

Negative Control: N
Media/Lot: RM0061 - TSA Rodacs - Release BARCODE : G17206 LOT NUMBER : G17206 LOT EXP DATE : Nov 16 2008 12:00AM VENDOR : Biomerieux
No Test: N
Plan: 001 - Cellular 123 - In-Process
Equipment/Control#: ENV-25 BARCODE : ENV-25 CALIBRATION DATE : Jul 2 2008 12:00AM CALIBRATION VALUE : 0 PERFORMED BY : N/A NEXT CALIBRATION DATE : Jan 31 2009
Site/Class: 123-A - BSC-08
Environment: D
Deviation: ACTION
Person Sampled:
Sampling Date: 8/29/2008
Personnel Site:

Product(s):

Product	Lot
AGS-003-006-002-002	36004-08N01, 08Z01
AGS-003-006-002-002	36004-08Z01, 08N01, 08P01

Approval Date

From: To: Range:

Result Date

From: To: Range:

Sampling Location

Product

Plan

Test Method

NONViable Particulate Count
PERSONG Personnel - Gloves
PERSONS Personnel - Sleeves
RODACSDA RODAC - SDA
RODACTSA RODAC - TSA

Deviation Type

Action
Alert
Pass
Non Zero

Get Results Clear Filters

Ready 119 records were returned.

start MODA-EM 3:59 PM

PAPERLESS QC MICRO AT AMYLIN



- ◎ **MODA™ Solution provides the ability to:**
 - Regulate and control existing and new sample locations
 - Define and set sample frequencies (discrete or standardized)
 - Immediately report nonconforming values
 - Define and control equipment and materials required for sampling and testing activities
 - Track the lifecycle of samples
 - Monitor and track plant flora by rooms, locations, and systems
 - Paperless - barcode scanning

AUTOMATED REPORTING AND TRENDING AT AMYLIN



- ◎ MODA™ Solution makes investigations into aseptic gowning excursions easy. Specifically,
 - Gathers and reports the data (trends overtime, specific areas where microorganisms were recovered, etc.)
 - Removes the human element and reports the facts
 - Data for the classified rooms and systems are housed in the same system

SUMMARY - WHAT DOES SUCCESS LOOK LIKE?

- ◎ A robust and compliant aseptic gowning qualification program is a program that utilizes:
 - A fully qualified 21CFR Part 11 compliant system
 - A system where all data can be tracked and trended together and timely
 - A system that tracks Samples Frequencies, Sample Locations, Sample Plans, Chain of Custody, Materials, and Equipment
 - Standardizes the Requirements and Guidelines for Investigations, Immediate Actions, and Corrective Actions



To submit a question, use the “Q&A” feature of WebEx (bottom right of your screen). If we do not answer a question online, we will be sure to follow up with an e-mail.

Lonza

Questions & Answers

Upcoming News & Events

See Lonza and MODA™ solution in action

- Jun 21-23: IVT ACE, Philadelphia, PA
- Jun 28 – Irish Cleanroom Society, Dublin, Ireland
- Jun 29-Jul 1: Interphex Japan 2011 – Tokyo, Japan
- Aug 1-3 : International Assoc of Food Protection, Milwaukee, WI

Upcoming Webinars (topics under consideration)

- Non-sterile Manufacturing Quality Control
- Advancements in Automated Water Testing
- Process optimization – making the business case
- Client Case Studies

Wrap-up

Personal Consultation

Bob Toal - Segment Manager, Informatics, Lonza Wayne

Phone: +1 484 253 1000 x133

Email: robert.toal@lonza.com

Learn more about MODA Solution: www.lonza.com/moda

Lonza

Thank You