

QC Micro Webinar Series

**Lonza**

---

## **Lonza's Paperless EM Success & Acquisition of MODA**

Rapid Testing Solutions / Lonza / 20 July 2010

# Today's Presenters

---

- **Douglas Danne**

Head, Rapid Testing Solutions, Lonza Bioscience

- **Bob Toal**

Segment Manager , Informatics, Lonza Wayne (formerly MODA)

- **Michael Goetter**

Director, Informatics Strategy, Lonza Wayne (formerly MODA)

- **Susan Harrison**

Senior Manager, Microbiology, Lonza Walkersville

## 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

---

- Lonza's acquisition of MODA Technology Partners
- MODA™ solution overview
- Paperless EM case study: Lonza's Walkersville facility
  - Update: Lean Six Sigma Initiative using MODA™ product line
  - ROI: Financial, Compliance, User Adoption

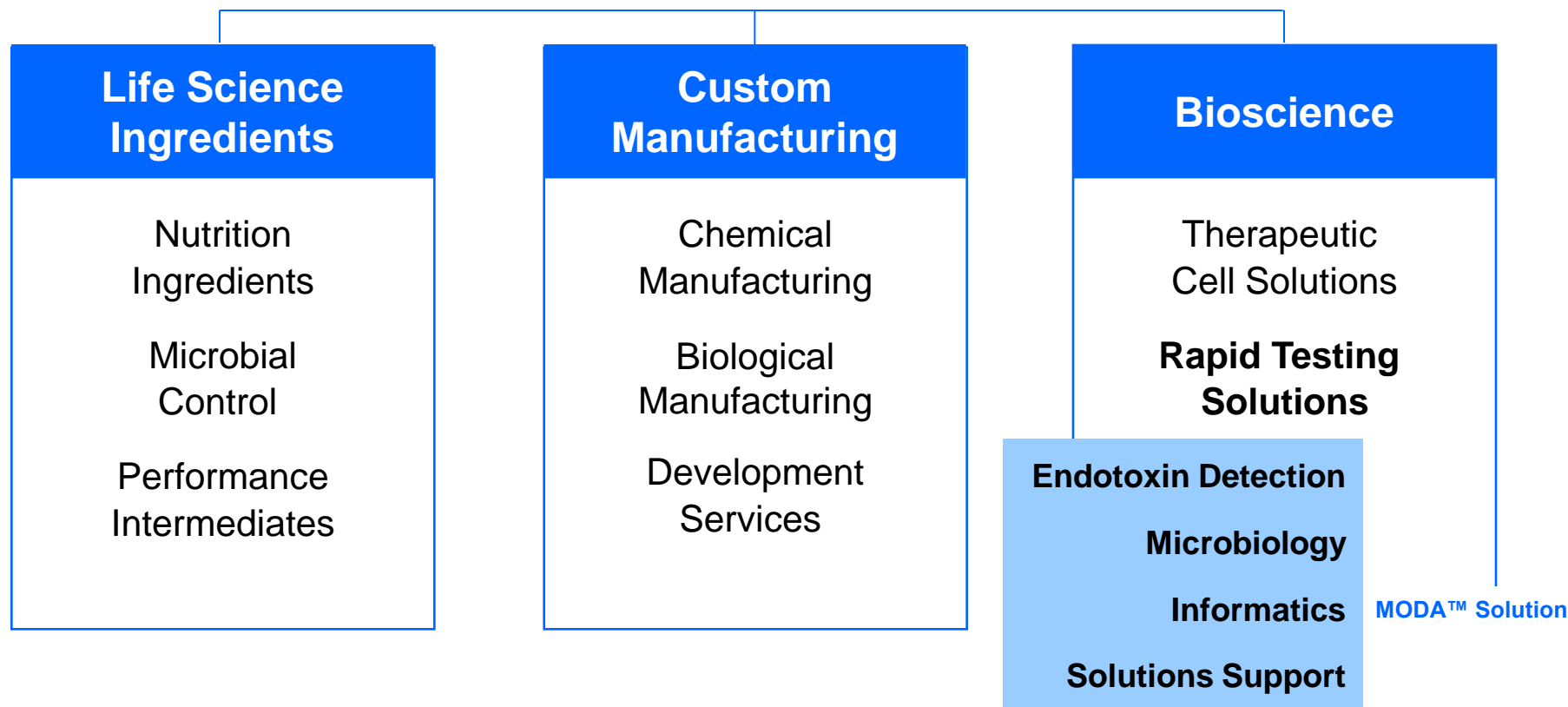
# Lonza

## Lonza Acquisition of MODA

# Lonza's Life-Science Platform

---

## Lonza



# Rapid Testing Solutions

---

- **Endotoxin Detection** - LAL-based assays, sustainable recombinant Factor C technology, plate readers and robotic liquid handling systems
- **Microbiology** - Includes the microCompass™ system, a qPCR-based method that detects bacteria, yeast and mold in less than six hours with applications for sterility and bioburden testing
- **Informatics** - Integrated and flexible software solutions, including the WinKQCL® system and now the MODA™ solution providing results on demand, helping customers make intelligent decisions
- **Solutions Support** - Service programs, equipment and software validation, LIMS support, instrument training and lab testing services

# Walkersville, MD (USA)

---



**Lonza Walkersville, Inc.**

- Area 2.8 ha
- Employees: Over 500
- Bioscience Headquarters
- Plants/Products
  - Liquid and powder media preparation and filling
  - Cell therapy and cell discovery
  - Rapid testing service
  - LAL processing
  - Contract services
- Products
  - Powdered media
  - Liquid media
  - Primary cells
  - Cell culture reagent
  - Endotoxin detection kits
  - Microbiology testing kits



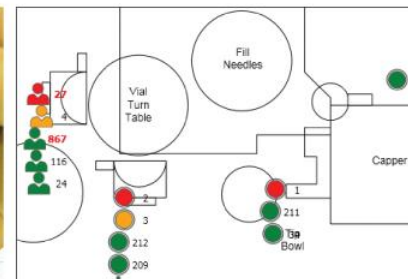
# QC automation just got greener!

MODA is now a Lonza company. More science. Less paper.™

Lonza's Rapid Testing Solutions now features mobile and wireless technologies by MODA—a paperless QC Micro solution. What started as a company-wide initiative to eliminate waste, reduce lead and labor time, and improve overall compliance, inspired a robust informatics platform for the life sciences industry. Automate your data collection. Gain improved insight into your manufacturing operation. Capture and upload test sample data for EM, utilities, and products at point of sample. Track and trend your data for immediate analytical reporting. Even see results through our visualization tools.

**Rapid Testing Solutions. Go paperless today!**

Explore our offerings at [www.lonza.com/rts](http://www.lonza.com/rts)



# Lonza

---

## MODA™ Solution

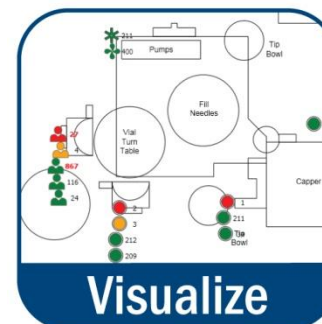
# MODA™ Solution Value Proposition

## ■ More science. Less paper.™

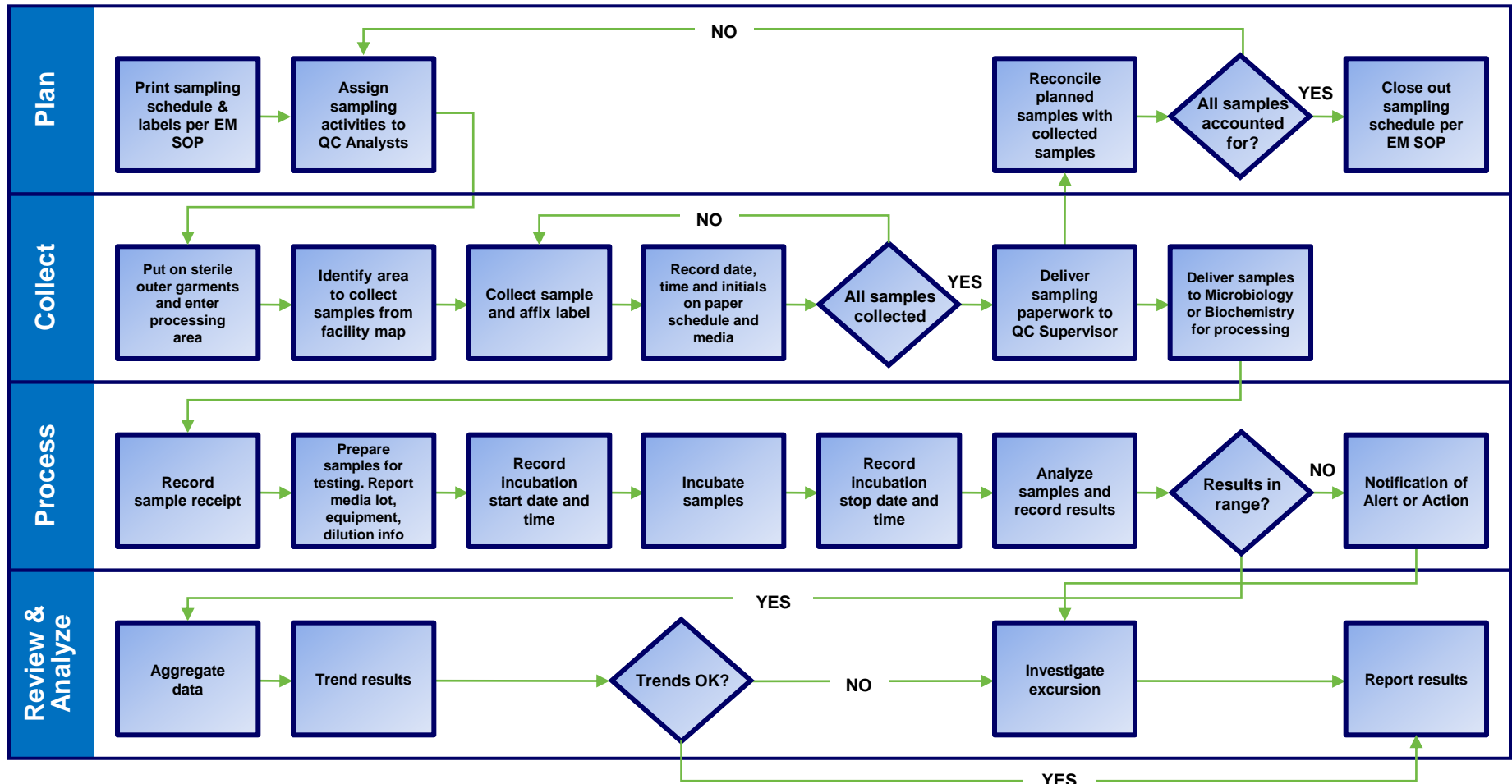
- Quickly move from paper-intensive QC Monitoring & Analysis

## ■ Increase operational efficiency, improve quality, reduce costs

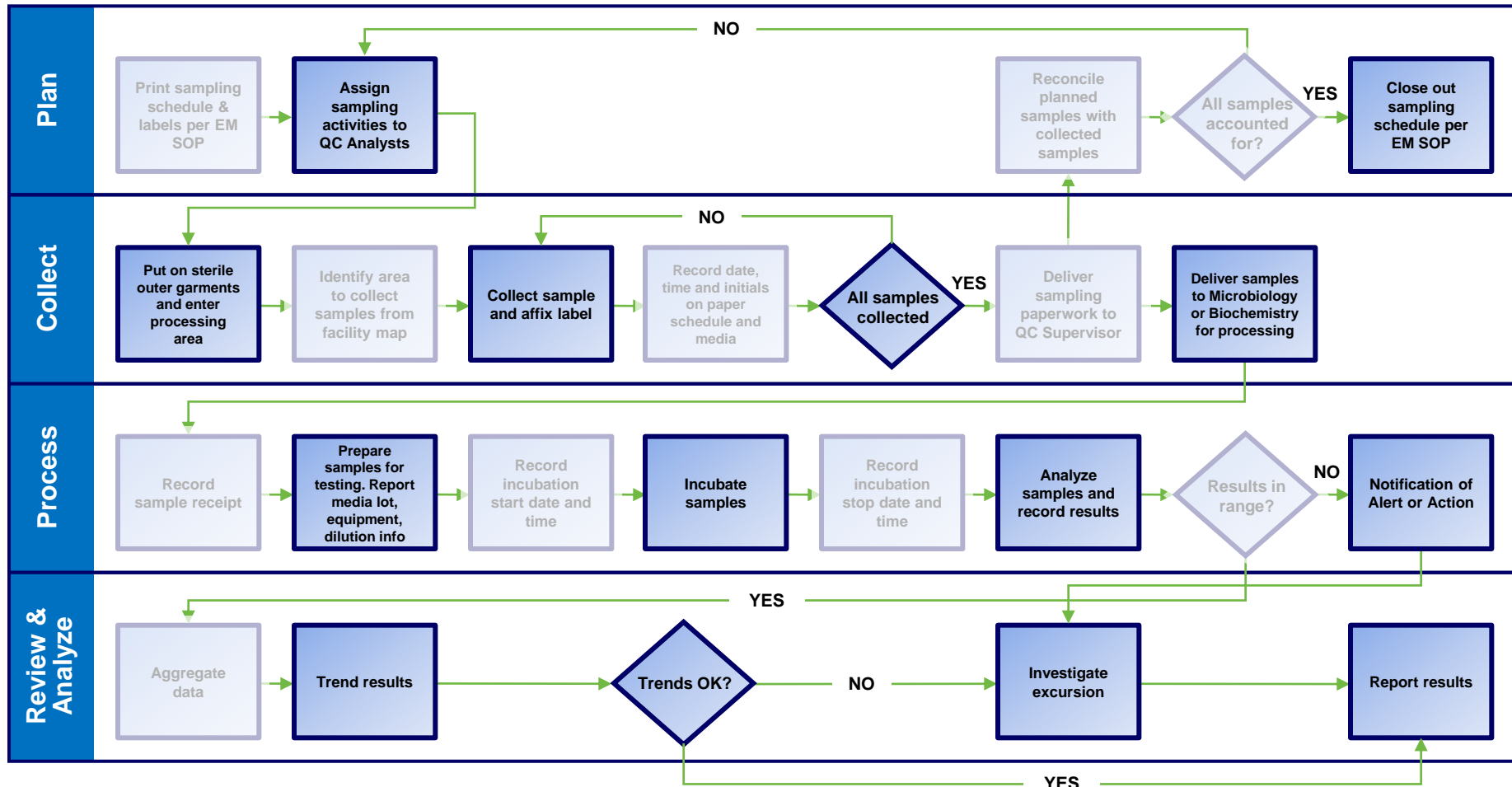
- MODA-EM™ platform offers mobile computing technology & advanced visualization tools






# The Paper-based QC Process



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



# Paperless Efficiency Example

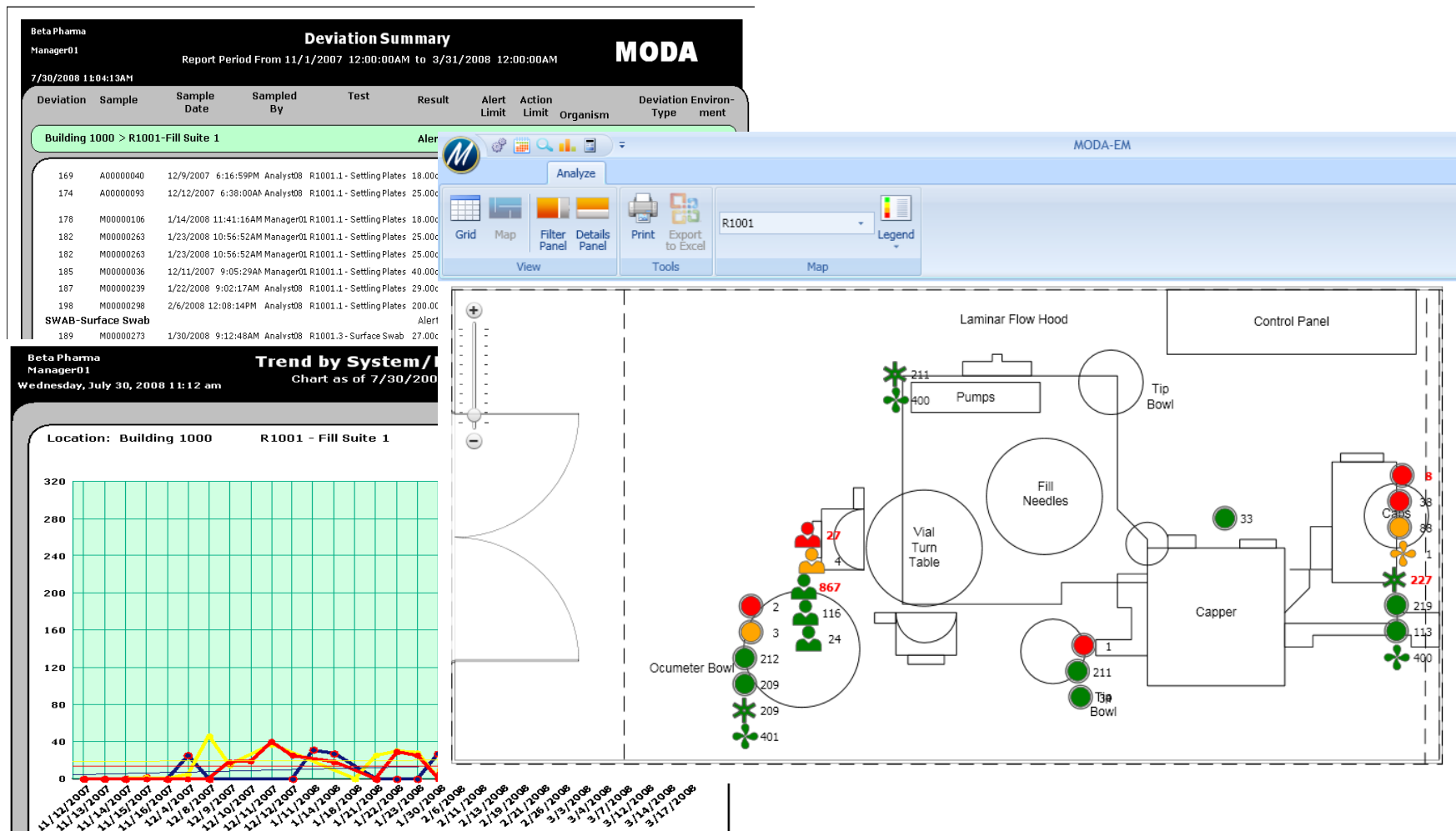
Record on paper template	Record on plate	Reconcile media and template	Enter data into Excel/LIMS
1: BB50, SS43, Water, IE5 2: BB50, SS43, RODAC, IP72/IP82 3: BB50, SS44, Water, IE5/IE7 4: BB50, SS44, Water, IE9/SP2 5: BB50, SS65, RODAC, IP74			
6 hours		1 hour	1 hour

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

	Paper	8 hours
–	Paperless	4 Hours
=	<b>Savings</b>	<b>4 Hours</b>

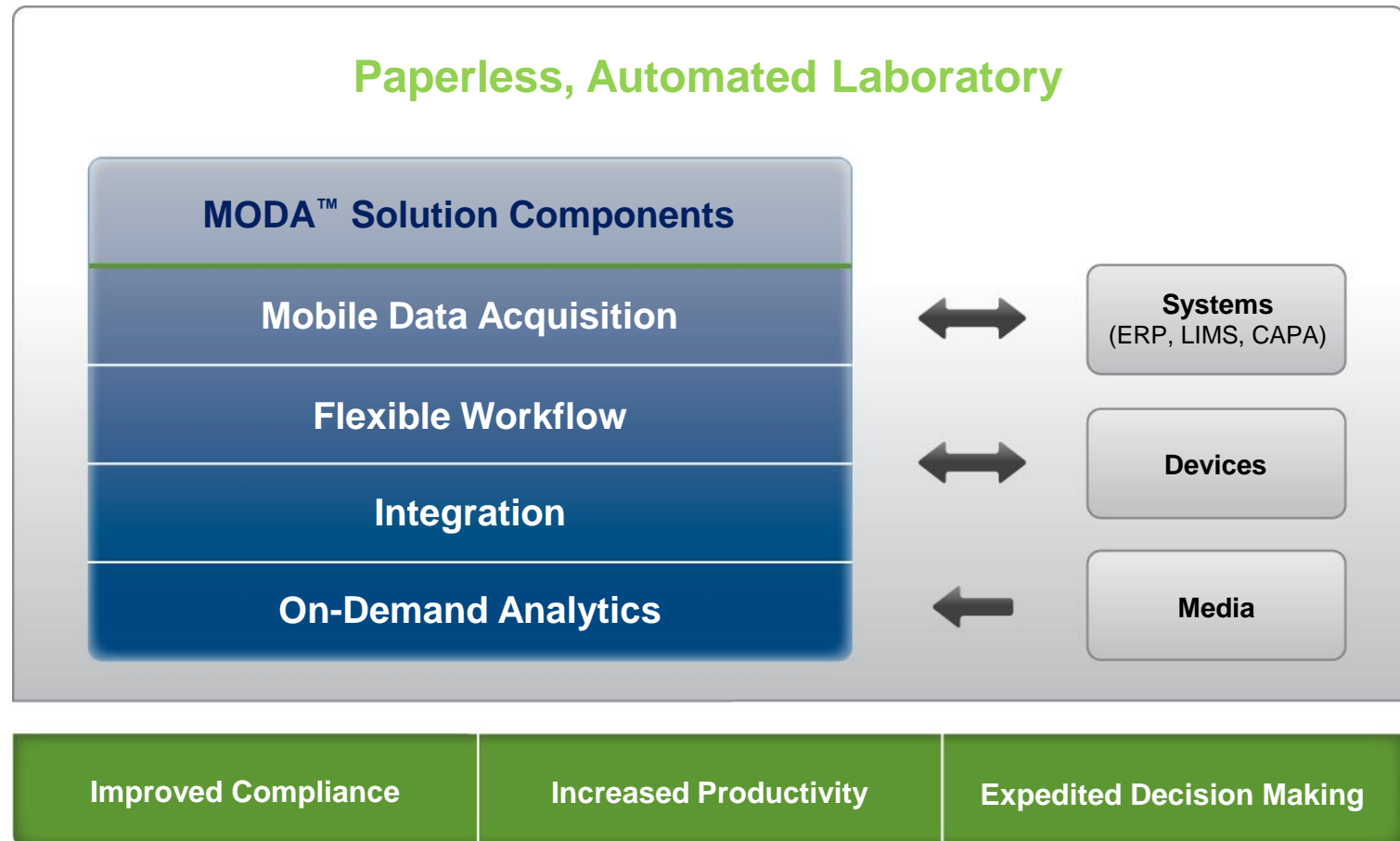


# Automated Reports & Analytics



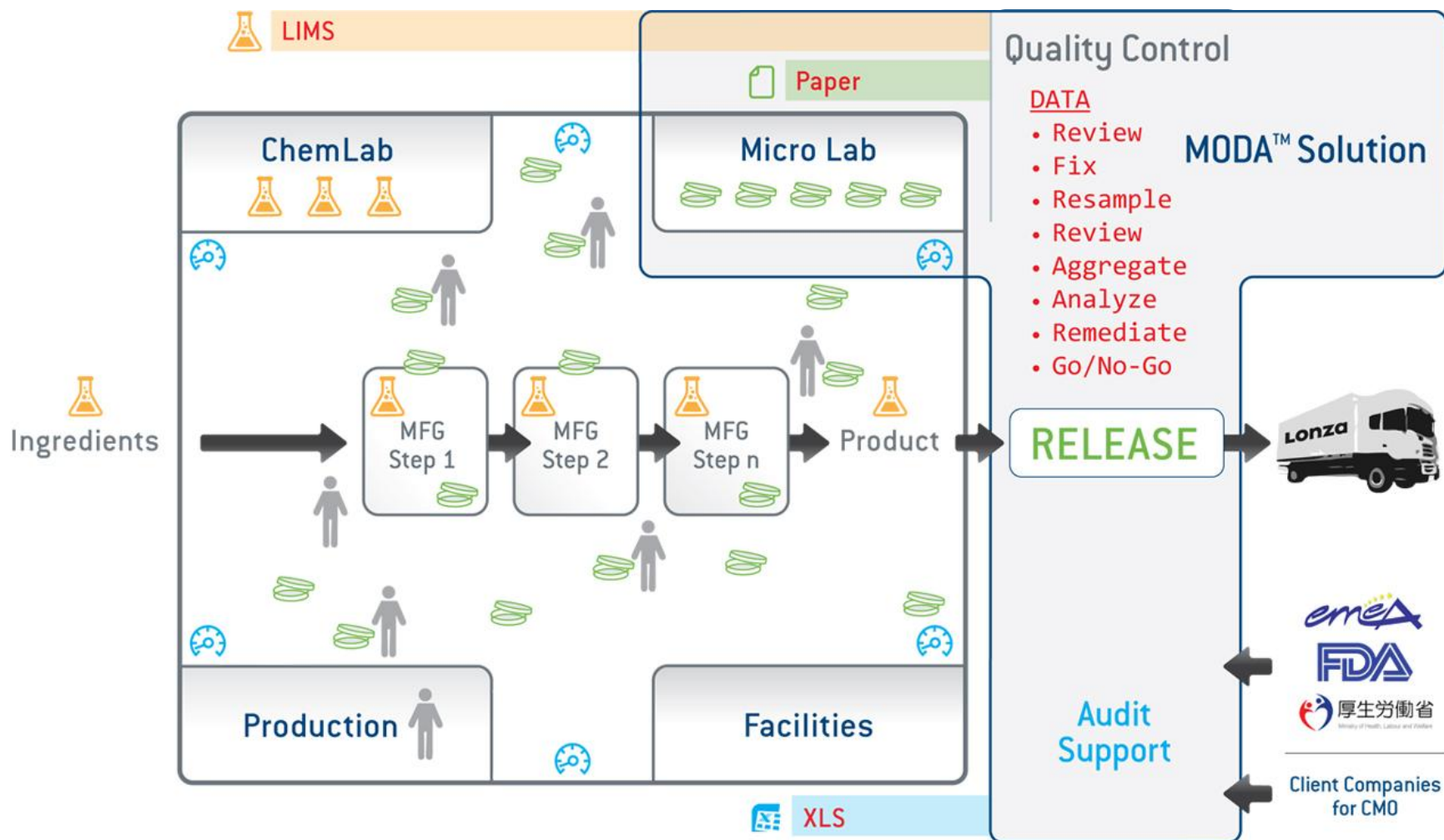
# Now: MODA™ Solution Concept

---



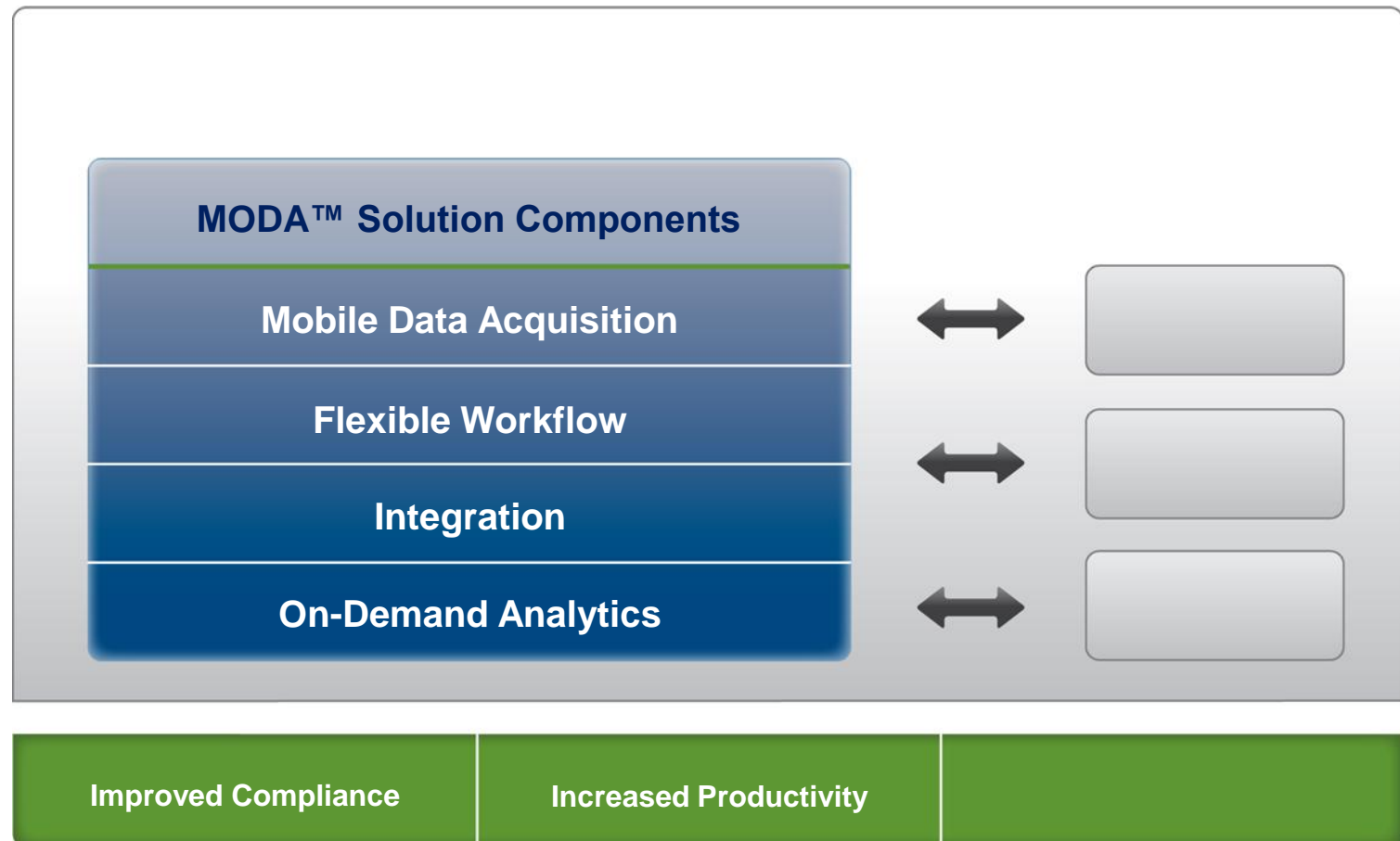


# Now: MODA™ Solution Concept

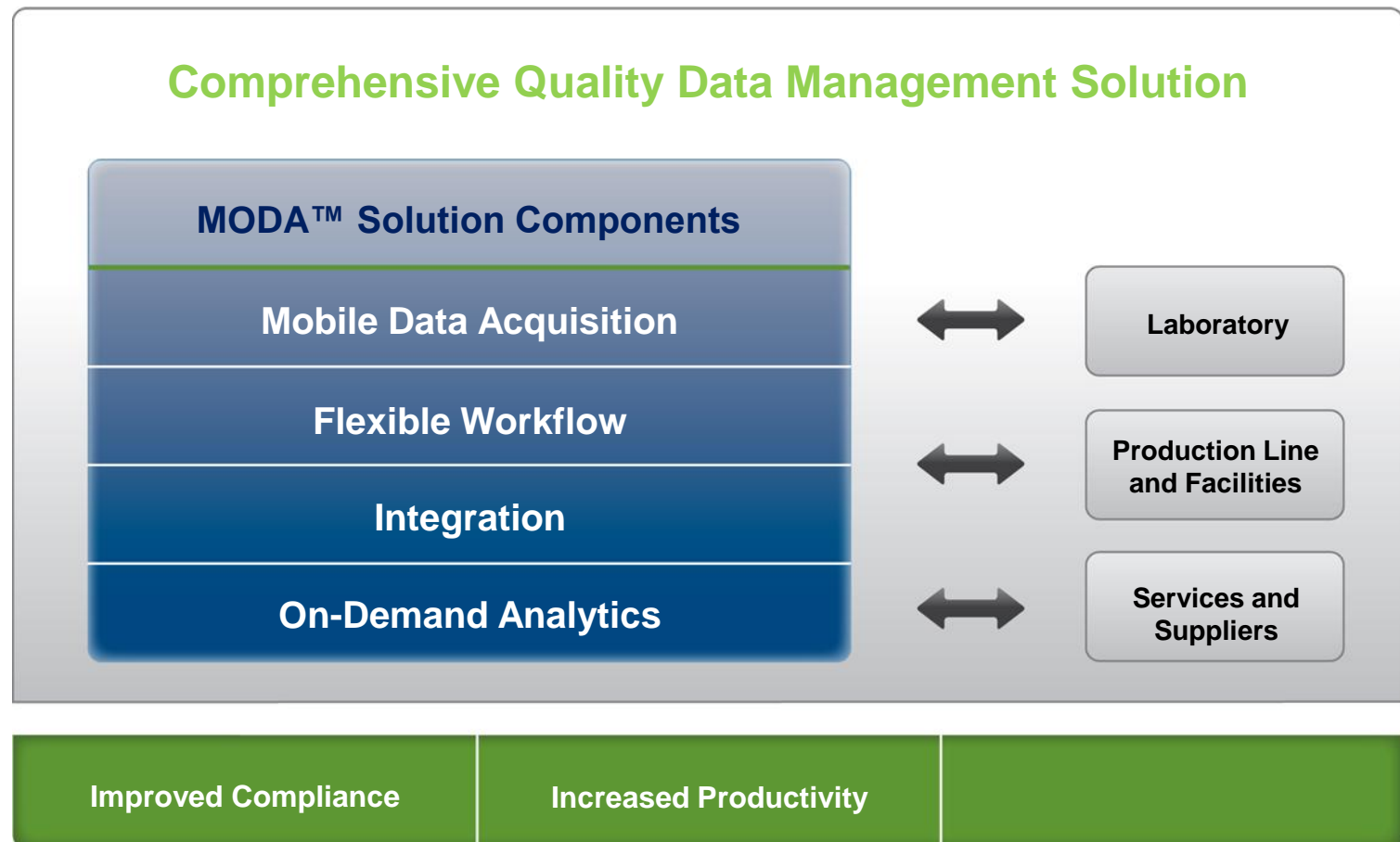


# Looking Forward: MODA™ Solution Vision

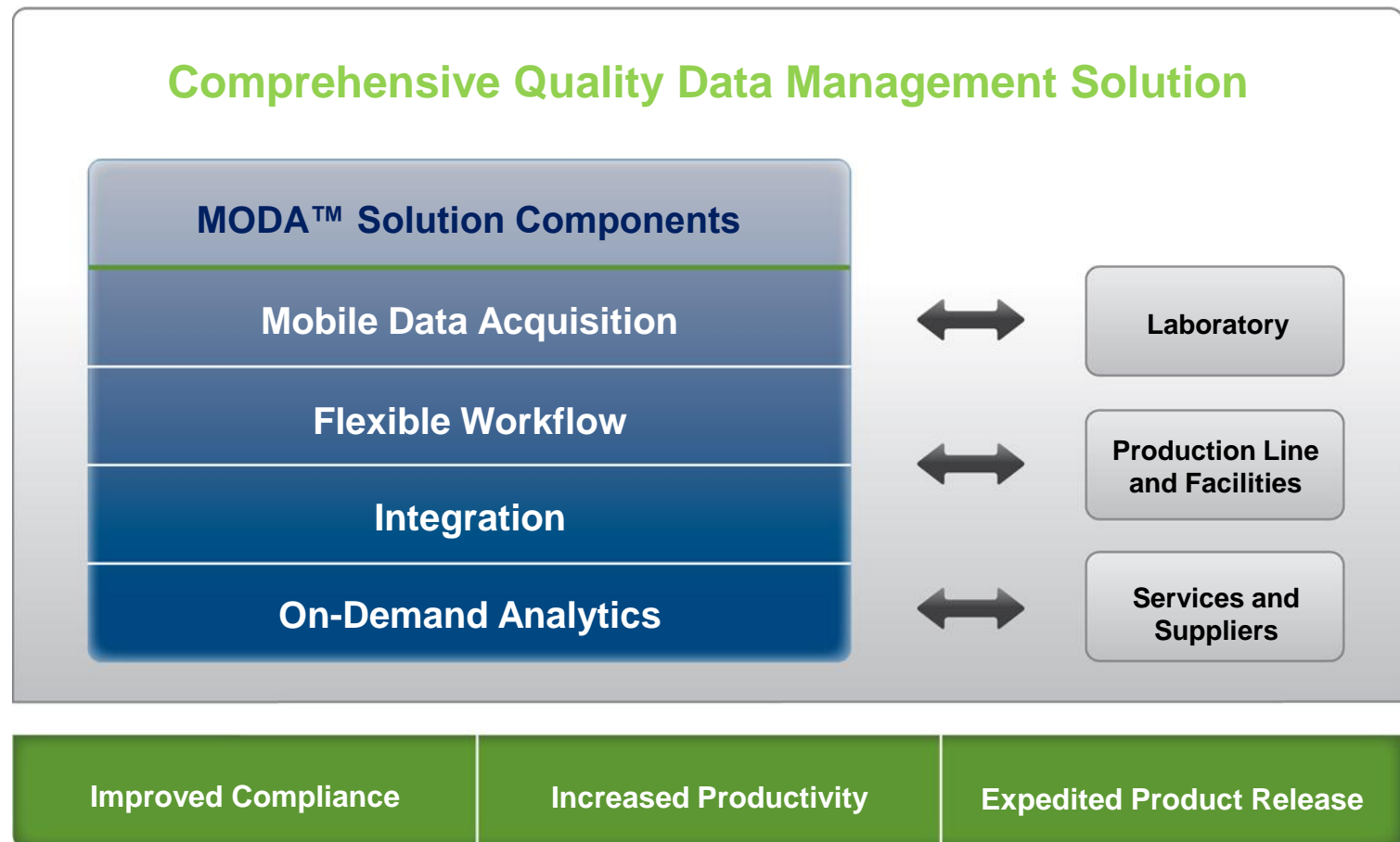
---



# Looking Forward: MODA™ Solution Vision



# Looking Forward: MODA™ Solution Vision





# Walkersville Case Study - MODA-EM™ Platform

# Walkersville MODA™ System Implementation

---

- 12-2007: Beta-Site implementation
  - 3 cleanrooms using Environmental Monitoring
  - Previous MODA software version
  
- 04-2009: Site-wide Environmental Monitoring implementation
  - 13 cleanroom suites
  - Current MODA software version
  
- 04-2010: Utility Monitoring implementation
  - 5 current water loops
  - 3 new loops added in Q2

## Before MODA™ Solution – Recurring Issues

---

- Paper based system (labor intensive)
- Duplication of effort
  - Completing paperwork
  - Data entry
- Timely data review (volume)
- Paper-Based MS Access database
  - Limited selection of reports
  - Excessive time spent generating trending reports
  - Not validated or CFR21 Part 11 compliant
    - Requires data verification

## Paper-Based System vs. MODA™ Solution

---

### ■ Paper-Based System

Particle-generating paper  
autoclaved into clean room

Data capture onto paper

Redundant Entry of data  
information in access/excel

Numerous Corrections for  
Omissions and Errors

Trending is manual, time-  
consuming and limited in scope

### ■ MODA™ Solution

✓ **Clean room compatible:**  
***“No paper”***

✓ **Real-time electronic data  
capture**

✓ **One-step data capture of  
information**

✓ **Reduces corrections through  
mandatory fields/dropdowns**

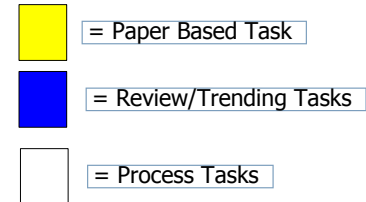
✓ **System generates a variety of  
automatic trend reports**



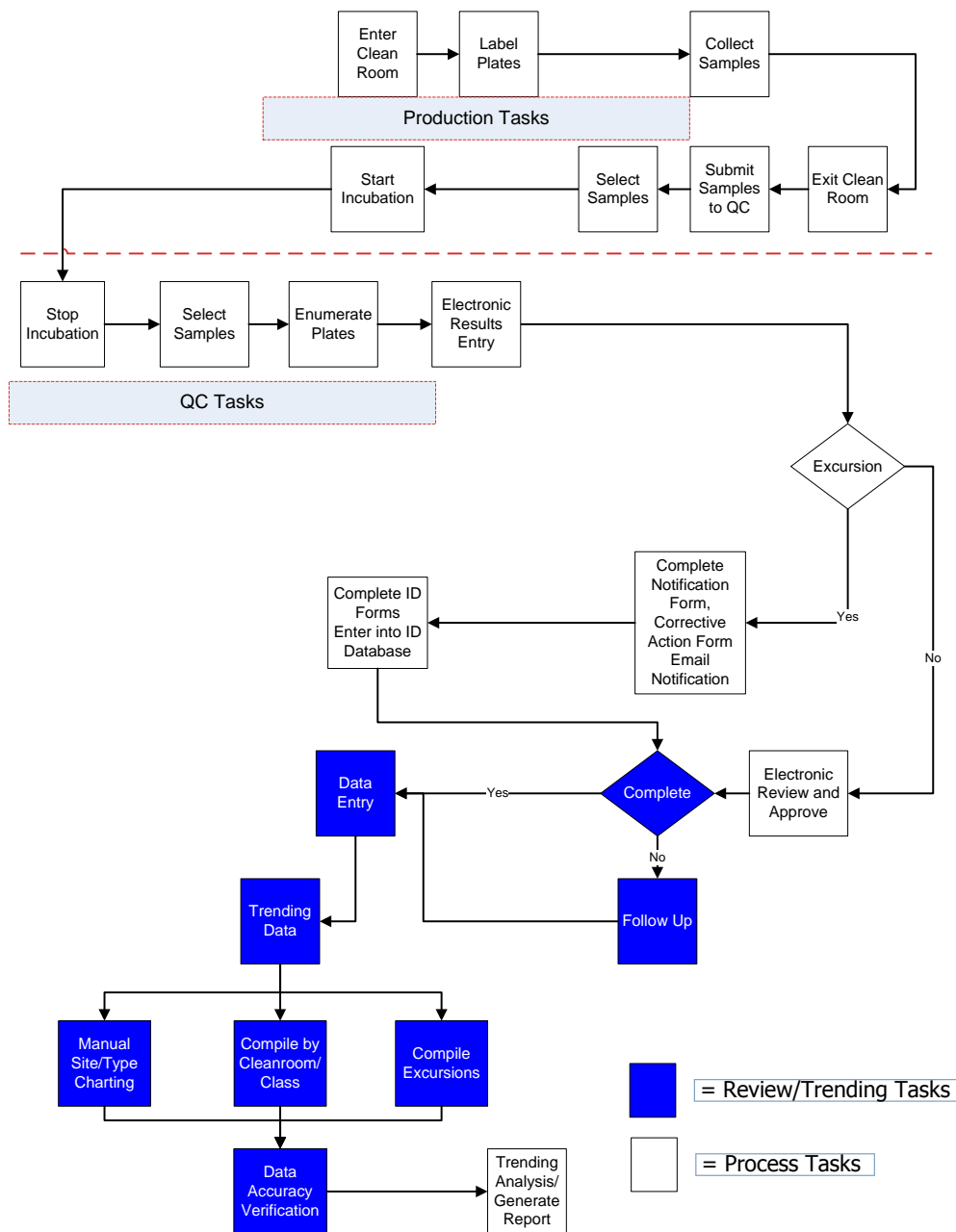
## **MODA™ Solution Improves EM Processes**

---

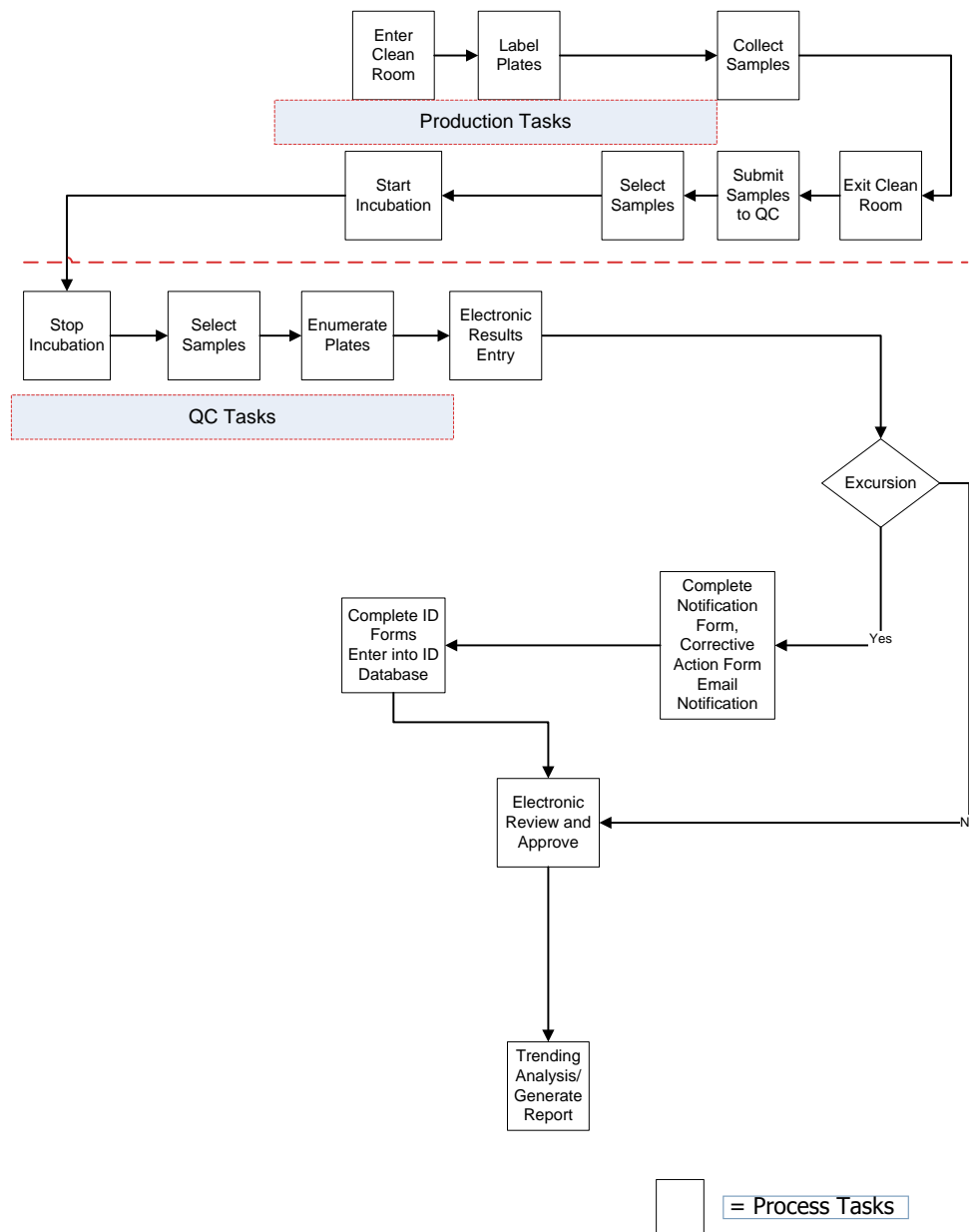
- Eliminates several steps from the paper-based EM Process
- Reduces Task and Process Time
  - Task Time – Hands-on time to complete each task
  - Process Time – Turn Around Time
- Data readily available
- Improves overall compliance



# Process Flow without Paper Tasks



# Process Flow with MODA™ Solution



# Financial Benefits

---

## ■ Estimated

- 1.3 hours/day x 3 beta sites = 3.9 hours/day
- 1.3 hours/day x 11 clean room suites = 14.3 hours/day
- 18.2 hours/day ÷ 8 hours day/analyst = 2.25 FTEs

## ■ Realized

- Staff Reduction from 11 to 8 FTE's (~25%)
- Miscellaneous savings (binders, paper, ink, particle-counter tape, autoclave costs, archiving and storage of data)
- Shift of resource to other value-added activities

## ■ Favorable ROI

## cGMP Compliance

---

### ■ Paper-Based System

Significant number of documentation errors

Long turn around time for processing data

Paper presents particulate and microbial contamination risk

GMP Compliance Risk

### ■ MODA™ Solution

✓ **Eliminates documentation errors**

✓ **Short-turn around time for processing data**

✓ **Clean room compatible equipment reduces contamination risk**

✓ **Sustainable GMP compliance**

## 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**



## Benefits of MODA™ Solution

---

- ✓ 21 CFR Part 11 compliant
- ✓ Instant data entry eliminating paper transpositions
- ✓ Generation of timely trending reports
- ✓ Equipment Interface (MetOne Particle counter)
- ✓ Eliminates duplication of efforts in data entry and trending
- ✓ Storage of paper documentation eliminated
- ✓ Clean room compatible equipment
- ✓ Provide Superior Quality System to our clients (potential revenue generation)
- ✓ Severability: Continues sampling process without wireless internet connection
- ✓ Added bonus: DMS access through PC tablets
  - ✓ i.e. Documentum, NovaManage

# Now: 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



# Now: Automated Lab Processing



# Today's 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](#)

## Today's 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](#)

**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

---

## Week of Nov 8

- 2-day QC Automation Training Workshop, Walkersville, MD

## See Lonza and MODA™ solution in action

- Sep 21-24: ILMAC, Basel
- Sep 28-30: ScanLab, Copenhagen
- Oct 5-8: Biotechnica , Hannover, Germany
- Oct 18: Quality in BioPharma (QBC), Raleigh, NC
- Oct 25-28: PDA Microbiology, Washington DC
- Nov 4: Paperless Lab, Bonn, Germany
- Nov 14-19: AAPS, New Orleans
- Nov 15-17: IVT ACE, Europe

## Sep 2010

- Feature Article in Pharma Technology

# Upcoming Webinars

---

## Sept 2010

- IT Perspective on Automated Quality Control

## Late Fall (topics under consideration)

- Advancements in Automated Water Testing
- Non-sterile Manufacturing Quality Control
- Client Case Studies



# Wrap-up

---

## Personal Consultation

Bob Toal - Segment Manager, Informatics

Phone: +1 484 253 1000 x133

Email: [robert.toal@lonza.com](mailto:robert.toal@lonza.com)

## Learn More

Web: [www.lonza.com/rts](http://www.lonza.com/rts)

# **Lonza**

---

**Thank You**