# Paperless QC Micro Webinar Series



# Introducing Hach's MET ONE Compliance Workflow System™ for Paperless Air Monitoring Powered by MODA

Presented by

Scott Hjelmervik

Systems Director Hach Company

**Bob Toal** 

Vice President
MODA Technology Partners

# Agenda



- Introduction to Hach Company and MODA Technology Partners
- Air Particle Monitoring in the Pharmaceutical Industry
- Paper-based vs. Paperless Air Particle Monitoring Processes
- MET ONE Compliance Workflow System Demo
- Summary of Features and Benefits
- Questions and Answers

A copy of this presentation and recording to this live session will be available for download at <a href="https://www.modatp.com">www.modatp.com</a> and <a href="https://www.modatp.com">www.modatp.c

#### MET ONE Air Particle Counters



 The Hach MET ONE brand is the world's leading series of airborne particle counters for regulatory compliance







MET ONE HHPC
Handheld particle counter
Used for troubleshooting

MET ONE 3400 Series Cleanroom classification Environmental monitoring Short-term online sampling

MET ONE 6000 & 7000 Series for production monitoring Grade A & B areas

# MODA Technology Partners



- Provide mobile data acquisition solutions that automate regulated manufacturing processes
  - Pharmaceuticals to consumer products
- Office locations
  - Wayne, PA Headquarters (suburban Philadelphia)
  - London, England
- World-wide sales in North America, Europe, and Asia
- Strategic partners

# MODA: Value Proposition

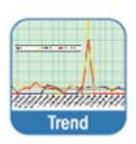


- More science. Less paper.
  - Quickly move from paper-intensive
     QC Monitoring & Analysis
- Increase operational efficiency, improve quality, reduce costs
  - MODA-EM™ offers mobile computing technology and advanced visualization tools











# MET ONE and MODA Partnership



- Environmental monitoring is the most labor intensive aspect of many particle monitoring programs and it generates the most paper
- Since 2007, MET ONE and MODA have worked together to develop an automated solution for environmental monitoring for our customers



















# MET ONE & MODA Partnership

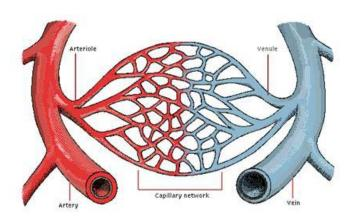


- MET ONE Compliance Workflow System (MET ONE CWS)
  - Affordable and low-impact entry point for paperless quality control monitoring and analysis of non-viable particle counts
- Leverages MODA-EM specific features:
  - Scheduling & Workflow
  - Data Acquisition and Storage
  - Alert and Action Notifications
  - Reporting and Analytical Trending Tools
- Now available for all existing and future MET ONE 3400 series air particle counters from Hach Company

#### Why do we care about Particle Counts?



- Particles as "contamination" negatively affect a process or product
  - Chemical composition
  - Stability
  - Purity
  - Safety
  - Reliability



- Pharmaceutical, Biotechnology
  - Particles in injection could cause occlusion of blood vessels

Red Blood cells are about 5 µm

Capillary (5 to 10 µm)

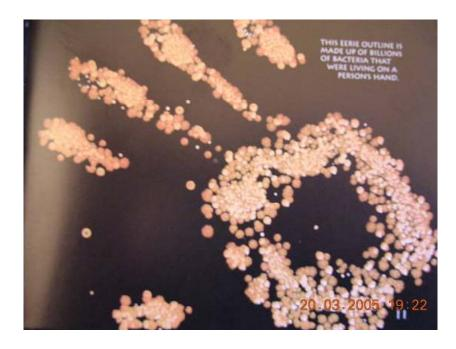
Large veins (10 to 50 µm)

- Viable particles in injection can trigger infection
- Possibility of reaction to foreign substances (RES/allergic reaction)

# Why monitor for particles?

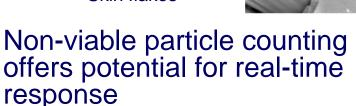


Greatest concern is for viable microorganisms



Non-viable particle count monitoring is a critical component of the total Environmental Monitoring program

- Technology is not available today to measure viable counts in real time
  - Requires incubation time
- Total "non-viable" particle counts used as a surrogate
- Non-viable counts
  - Includes all types of airborne material
    - Solid particles
    - Fibers
    - Microorganisms
    - Skin flakes

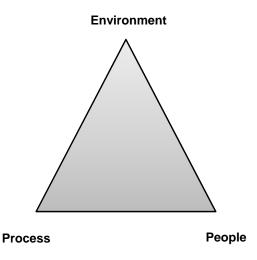


#### Non-Viable Particle Count Monitoring



- Classification of Cleanrooms formal process to ensure the environment meets class standard
  - Rigid standards apply (ISO14644)
- Aseptic Process Monitoring

   – risk-based
   process to ensure that the people, process
   and the environment remain in control
   during the desired task
  - cGMP/GMP Guidance from FDA/EMEA
- Environmental Monitoring risk-based process to verify that the environment is ready to carry out the desired task
  - Inferred guidance, not rigid by regulation



#### Cleanroom and Clean Zone Classification



INTERNATIONAL ISO 14644-1 STANDARD Cleanrooms and associated controlled environments — Part 1: Classification of air cleanliness Salles propres et environnements maîtrisés apparentés — Partie 1: Classification de la propreté de l'air

- Classification is the formal process of qualifying the environment by the number of particles using a standard method (ISO 14644)
  - Performed on a regular basis but <u>not</u> <u>frequently</u>
    - Grade A areas: Six months, "ISO 4.8"
    - Grade B areas: Six months, ISO 5
    - Grade C, D areas: Annually, ISO 7/8
  - Standards define minimum number of points
    - Based on area of cleanroom or clean zone
  - Standards define minimum amount of air to be sampled
    - Minimum volumes for statistically valid samples (typically 1 minute at 1 cfm)
    - Grade A requires minimum 1 m3 (Annex 1)

Classification is a rigid protocol

### ISO 14644 Classification Limits



Class	Number of Particles per Cubic Meter by Micrometer Size					
	0.1 μm	<b>0.2</b> μm	<b>0.3</b> μm	<b>0.5</b> μm	1 μm	5 μm
ISO 1	10	2				
ISO 2	100	24	10	4		
ISO 3	1,000	237	102	35	8	
ISO 4	10,000	2,370	1,020	352	83	
ISO 5	100,000	23,700	10,200	3,520	832	29
ISO 6	1,000,000	237,000	102,000	35,200	8,320	293
ISO 7				352,000	83,200	2,930
ISO 8				3,520,000	832,000	29,300
ISO 9				35,200,000	8,320,000	293,000



Typically performed with portable counters

# Aseptic Process Monitoring



- Governed by EMEA and FDA Guidance
  - For critical (Grade A) areas and Grade B
  - Recommends continuous monitoring
- Number of sample points based on risk assessment
  - Minimum locations is well defined by common practice
- Location of points based on risk assessment
  - Some guidance on sampling location
- Sample volumes defined by need for event detection

Published guidance, industry practice

May be manual or automated remote process



# Aseptic Process Monitoring



- Pharmaceutical Industry Compliance Guidance
  - FDA cGMP Guidance for Industry
    - Sterile Drug Products
       Produced by Aseptic
       Processing
  - EU GMP Annex I (EMEA)
    - Manufacture of Sterile Medicinal Products

**EU GMP Annex 1 is more rigid than FDA cGMP Guidance** 



EUROPEAN COMMISSION ENTERPRISE AND INDUSTRY DIRECTORATE-GENERAL

Consumer goods

Brussels, 14 February 2008

EudraLex
The Rules Governing Medicinal Products in the European Union

Volume 4
EU Guidelines to
Good Manufacturing Practice
Medicinal Products for Human and Veterinary Use

#### Annex 1 Manufacture of Sterile Medicinal Products

Document History		
Previous version dated 30 May 2003, in operation since	September 2003	
Revision to align classification table of clean rooms, to include guidance on media simultations, bioburden monitoring and capping of freeze-dried vials	November 2005 to December 2007	
Date for coming into operation and superseding	01 March 2009 <sup>1</sup>	

Note: Provisions on capping of freeze-dried vials should be implemented by 01 March 2010.

Commission Européenne, B-1049 Bruxelles / Europese Commissie, B-1049 Brussel - Belgium. Telephone: (32-2) 299 11 11

# **Environmental Monitoring**



- Determine readiness of room to carry out designated task
- Performed whenever relevant activity will occur daily work
- Most labor intensive

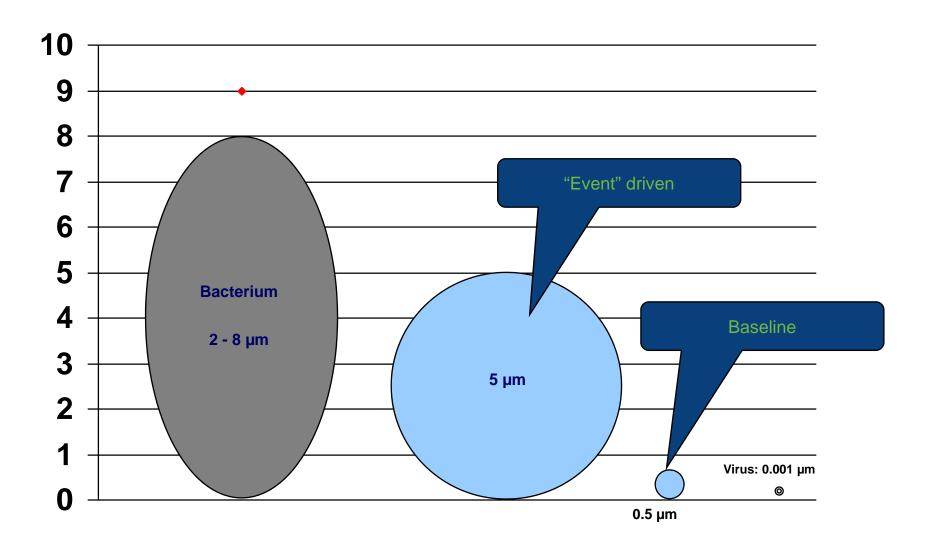


- Number of sample points defined by risk assessment
  - Activity to be performed
  - Risk to product
- Frequency and volume of sample points defined by risk assessment
  - Frequent enough to show control
  - Frequent enough to manage risk of product contamination

Inferred guidance, not rigid by regulation

# Particle Monitoring





#### Manual Monitoring: A Paper Based Process





- Technicians manually record the time, placement and calibration check of devices on paper
- Data is transcribed and reconciled manually from paper printouts to log books and spreadsheets
- Analysts spend days or weeks formatting data in spreadsheets to provide basic analysis and trends, especially when data needs correlation with viable monitoring activity

#### Key Issues with Current EM Methodologies



- The majority of non-viable particle monitoring performed is a time-consuming manual process governed by company specific written SOPs
- This process:
  - Is paper-based and error-prone
  - Is labor intensive
  - Requires high level expertise to understand and follow protocols
  - Results in slow corrective actions (and few pre-emptive actions)

There is a better way ... MET ONE CWS

# MET ONE Compliance Workflow System





#### MET ONE CWS Consists of:

MET ONE 3400 Portable Particle Counter

MET ONE CWS Software (Powered by MODA)

Laptop or tablet PC for online data collection

Server PC for Database

#### MET ONE CWS Software Offers



- Automated scheduling to drive workflow
- Paperless capture of portable particle count data (wired or wireless)
- Immediate notification of non-compliant events/trends
- Comprehensive trending, mapping and reporting tools
- Comprehensive data visualization tools
- Ability to export data to other systems (LIMS, SPC Tools)
- 21 CFR Part 11 compliant





# **Demonstration**

MET ONE Compliance Workflow System

#### MET ONE CWS Software



- Automated scheduling to drive workflow
  - Frequency based for EM routines
  - On-demand
- Paperless capture of portable particle count data (wired or wireless)
  - Run MET ONE 3400 directly from PC tablet or laptop
  - Supports disconnected operation
  - No paper tape and associated manual data entry and validation
- Immediate notification of non-compliant events/trends
  - Automatic email notification when alert or action limits are exceeded
  - Configure for below-limit excursions or frequency-based events

#### MET ONE CWS Software



- Comprehensive trending, mapping and reporting tools
  - Quickly generate required reports
  - Trend charts and statistical summaries
  - Ad-hoc analysis for determining root cause
- Comprehensive data visualization tools
  - Particle count data overlaid on room diagrams
  - Workflow-driven ISO 14644 sampling: more than calculations
    - ISO sampling pattern overlaid on facility blueprint as SOP
    - Samples automatically controlled for Grade A/B/C/D areas
- Automatic compliance checking for equipment calibration
- Data export to other systems (LIMS, SPC Tools)
- Allows you to be 21 CFR Part 11 compliant
  - Including E-signatures, data security, audit trail

# MET ONE Compliance Workflow System Requirements



#### Minimum

- Met One 3400 particle counters
- One laptop or tablet PC for each counter
- One server computer
- Wired or wireless Ethernet network

#### Optional

Bar code reader for each tablet PC

All MET ONE 3400 Particle Counters are designed to comply with ISO 21501



#### MET ONE CWS: Products and Services



#### MET CWS software license

 Support for up to six MET ONE 3400 counters

or

Support for more than six MET
 ONE 3400 counters

#### **AND**

- 1st year software support agreement
- Optional accessories
  - Bar code readers

#### Optional MET ONE services

- Validation documentation package
- Pre-installation consultation services
- On-site installation, commissioning and training
- On-site validation assistance
- Data integration with in-house LIMS system
  - Real-time updates or historical data

## **Upgrade Path to MODA-EM**



- Provides all MET ONE CWS features
- Automated, paperless monitoring
  - Viables testing
  - Utilities (water, compressed gases)
  - Product testing
- Additional reports for full QC Micro coverage
  - Trend particulate count data versus other EM and Utility monitoring results
  - Organism reports, personnel monitoring reports
  - Visualization of entire spectrum of QC Micro tests on facility maps
  - Additional device integration available
- Upgrade accomplished via configuration updates
  - No need for new software installation
  - Minimizes validation impact

## Like what you see?



#### Choose the MET ONE CWS Package:

- Automation of the non-viable particle count data collection only
- Lower cost to implement and validate
- Incremental "low-risk" approach to proving benefits of full EM automation
- Available from Hach Company

#### With an upgrade path ...

- Choose the full MODA-EM Package:
  - Automation of the entire environmental monitoring program
  - All elements of QC Microbiology (viable, non-viable, utilities, etc)
  - Available from MODA Technology Partners

# Summary



- MET ONE Compliance Workflow System
  - Automated scheduling to drive workflow
  - Paperless capture of portable particle count data (wired or wireless)
  - Immediate notification of non-compliant events/trends
  - Comprehensive trending, mapping and reporting tools
  - Comprehensive data visualization tools
  - Data export to other systems (LIMS, SPC Tools)
  - Complete support services
- Now available for all existing and future MET ONE 3400 series air particle counters from Hach Company



Bob Toal MODA btoal@modatp.com (484) 253-1000 x133

# **Questions & Answers**

→ 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.

A copy of this presentation and recording to this live session will be available for download at <a href="https://www.modatp.com">www.modatp.com</a> and <a href="https://www.modatp.com">www.modatp.c