

# Beckman Coulter MET ONE FMS – Continuous Viable and Non-Viable Particle Monitoring Systems



Beckman Coulter offers scalable solutions for viable and non-viable continuous particle monitoring systems to

- Enhance data integrity
- Simplify 21 CFR Part 11 Compliance
- Reduce risk of human error with automation
- Reduce likelihood of audit findings

## MET ONE Non-Viable Remotes

- MET ONE 6000 for central vacuum system
- MET ONE 6000P – self pumped for minimal impact to infrastructure. Ideal for BSCs and Isolators.
- MET ONE 7000 – specifically designed for aseptic monitoring with NEMA enclosure and local vacuum control



## Viable Active Air Remotes

- iVAS - viable self-pumped active air samplers
- 50 or 100 LPM air flow, continuous or interval sampling to your 90 mm sample plates



## MET ONE Non-Viable Portable

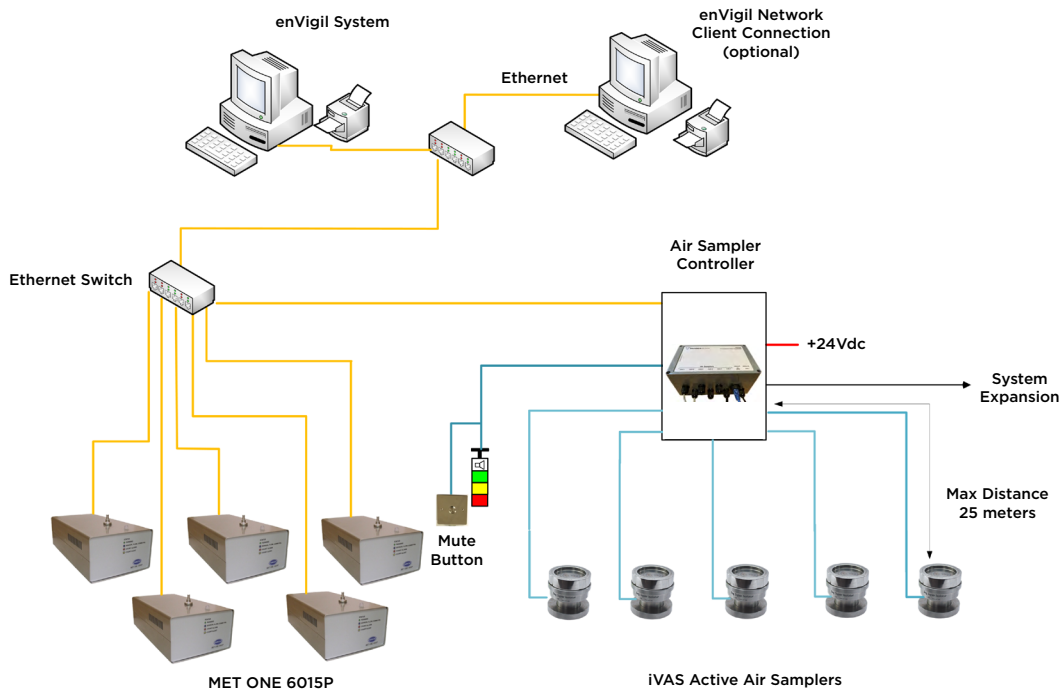
- Cleanroom Classification and Routine Environmental Monitoring for non-viable particles
- With LDAP and internal audit trail
- Remote control and configuration via web-browser
- Also used for continuous monitoring in BSCs and hoods with local on-screen control



**CHARACTERIZED**  
*by ingenuity.*

 **BECKMAN  
COULTER**  
Life Sciences

## enVigil Particle Monitoring System with MET ONE 6015P Particle Counters and Active Air Samplers



**Figure 1.** This is an example of a non-viable and viable self pumped system with enVigil PnP. Systems are configurable and expandable to meet your needs, including central vacuum sensors and larger systems running on virtual servers.

### Scalable Software for Small to Large systems-

#### enVigil PnP Software

- Pre-packaged software that is easy to set up and configure
- Ideal for smaller systems, such as isolators
- Non-viable and viable capability
- Temperature/ relative humidity and differential pressure sensor capability
- Validation Documents included
- Up to 20 air particle counters

#### enVigil FMS Software

- Ideal for customized and very large systems
- Non-viable and viable capability
- Drivers for multiple environmental sensors
- Customizable Software with built in workflow
- Batch Processing
- Multiple reporting and trending options
- Validation documents available
- Industrial Server Interface design

Contact your local Beckman Particle rep for more information:



© 2021 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](http://beckman.com)  
21.04.1467.PCC