



MET ONE 3411 Portable Airborne Particle Counter

MET ONE 3411

PORTABLE AIRBORNE PARTICLE COUNTER

**Does your particle counter
contaminate your cleanroom?**

Do you trust the results?

**How much time do you spend
recording and analyzing the data?**



Introducing the MET ONE 3411 0.1 micron sensitivity portable airborne particle counter. Unlike many particle counters the MET ONE 3411 fan remains off when operating in a standard semiconductor cleanroom environment thus particle generation is virtually eliminated.

Cleanroom friendly monitoring

Reduced particle impact
to cleanroom environment

Unit-to-unit accuracy and reproducibility

Assured through ISO 21501 compliance

Flexible communications

Wireless, Ethernet, Serial and USB

Long continuous operation

Dual hot-swappable batteries

Intuitive touch-screen user interface

Easy area, location, operating
parameter configuration and replication



The MET ONE 3411 is designed to meet the rigorous requirements of ISO-21501-4 to provide the users with unparalleled accuracy. This assures the repeatability of particle measurements between multiple instruments; critical for process benchmarking and troubleshooting activities.

Specifications

<i>Number of Size Channels</i>	6
<i>Particle Sizes</i>	0.1, 0.2, 0.3, 0.5, 1.0, 5.0
<i>Flow Rate</i>	1.0 CFM (28.3 LPM)
<i>Zero Count</i>	ISO 21501-4 and JIS B 9921: 1 count or less in 5 minutes, 95% UCL
<i>Coincidence Loss</i>	5% at 50,000 /ft ³ per ISO 21501-4 method
<i>Counting Efficiency</i>	50% at 0.1 µm; 100% for particles > 0.15 µm per ISO 21501-4 and JIS B 9921
<i>Light Source</i>	Helium Neon Laser, 5 mW Max Power at 632.8 nm
<i>Pump Type</i>	Patented multi-lobe closed loop controlled rated for continuous use
<i>Display</i>	¼ VGA Color TFT touch screen
<i>Printer</i>	High speed thermal
<i>Language</i>	English, French, German, Italian, Spanish, Korean*, Japanese*, Chinese (Simplified and Traditional)*
<i>Maximum Count Displayed</i>	9,999,999 displayed
<i>Delay Time</i>	15 seconds to 23 hours 59 minutes 59 seconds
<i>Sample/Hold Times</i>	1 second to 23 hours 59 minutes 59 seconds
<i>Count Alarms</i>	1 to 9,999,999 counts
<i>Data Storage</i>	5,000 samples, scrollable on Historical Data review screen (FIFO or Overflow)
<i>Locations</i>	ID: 0 to 999; NAME: Alphanumeric, appears on printout
<i>Outputs</i>	USB Client (Version 1.1) USB Host (Version 1.1) RS-485 Ethernet with TCP/IP protocol Wireless with 802.11g protocol (optional) MET ONE 2432 Manifold Auxiliary (alarm and scan probe)
<i>Communication Protocol</i>	Modbus TCP, Modbus RTU, Serial FX
<i>Inputs</i>	Air Velocity Probe RH/Temperature Probe
<i>Auto CDA Purge</i>	Purge solenoid activated by connection to CDA
<i>Battery Type</i>	Lithium ion smart battery; rechargeable, ejectable, and hot-swappable
<i>Battery Quantity Included</i>	2
<i>Battery Operating Time</i>	3 hours minimum
<i>Battery Recharge Time</i>	6.75 hours minimum, 10 hours maximum
<i>Power</i>	24 VDC with 100-240 VAC, 50/60 Hz adapter
<i>Enclosure Material</i>	Easy to clean passivated Stainless Steel
<i>Size</i>	33 W x 55.9 D x 22.9 H cm (13 x 22 x 9 inches) including protrusions, handles, feet, etc.
<i>Weight</i>	Without battery: 15.9 kg (35 lbs), Battery: 0.66 kg (1.45 lbs)
<i>Environment</i>	Operating 10° to 40°C (50° to 104°F); 10 to 90% relative humidity, non-condensing Storage -40° to 50°C (-40° to 122°F); 10 to 90% relative humidity, non-condensing
<i>Optional Accessories</i>	Smart Battery Charger/Calibrator: PN 280-300-5000 Additional Battery: PN 280-120-2024

Beckman Coulter Life Sciences
250 S Kraemer Blvd
Brea, CA 92821 USA
Telephone: 800-866-7889
E-mail: insidesalesgp@beckman.com
www.particle.com