

MET ONE 3411

Portable Airborne Particle Counter



CHARACTERIZED
by ingenuity.

 **BECKMAN
COULTER**
Life Sciences

MET ONE 3411

PORTABLE AIRBORNE PARTICLE COUNTER

Do you trust
the results?

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

Does your particle
counter contaminate
your cleanroom?



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

Ensured 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 ensures the repeatability of particle measurements between multiple instruments; critical for process benchmarking and troubleshooting activities.

Specifications

Number of Size Channels	6	
Particle Sizes	0.1 μ, 0.2 μ, 0.3 μ, 0.5 μ, 1.0 μ, 5.0 μ	
Flow Rate	1.0 CFM (28.3 LPM)	
Zero Count	ISO 21501-4 and JIS B 9921: 1 count or less in 5 minutes, 95% UCL	
Coincidence Loss	5% at 50,000 /ft3 per ISO 21501-4 method	
Counting Efficiency	50% at 0.1 μm; 100% for particles > 0.15 μm per ISO 21501-4 and JIS B 9921	
Light Source	Helium Neon Laser, 5 mW Max Power at 632.8 nm	
Pump Type	Patented multi-lobe closed loop controlled rated for continuous use	
Display	¼ VGA Color TFT touch screen	
Printer	High speed thermal	
Language	English, French, German, Italian, Spanish, Korean*, Japanese*, Chinese (Simplified and Traditional)*	
Maximum Count Displayed	9,999,999 displayed	
Delay Time	15 seconds to 23 hours 59 minutes 59 seconds	
Sample/Hold Times	1 second to 23 hours 59 minutes 59 seconds	
Count Alarms	1 to 9,999,999 counts	
Data Storage	5,000 samples, scrollable on Historical Data review screen (FIFO or Overflow)	
Locations	ID: 0 to 999; NAME: Alphanumeric, appears on printout	
Outputs	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)	
Communication Protocol	Modbus TCP, Modbus RTU, Serial FX	
Inputs	Air Velocity Probe RH/Temperature Probe	
Auto CDA Purge	Purge solenoid activated by connection to CDA	
Battery Type	Lithium ion smart battery; rechargeable, ejectable, and hot-swappable	
Battery Quantity Included	2	
Battery Operating Time	3 hours minimum	
Battery Recharge Time	6.75 hours minimum, 10 hours maximum	
Power	24 VDC with 100-240 VAC, 50/60 Hz adapter	
Enclosure Material	Easy to clean passivated Stainless Steel	
Size	33 W x 55.9 D x 22.9 H cm (13 x 22 x 9 inches) including protrusions, handles, feet, etc.	
Weight	Without battery: 15.9 kg (35 lbs), Battery: 0.66 kg (1.45 lbs)	
Environment	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
Optional Accessories	-40° to 50°C (-40° to 122°F); 10 to 90% relative humidity, non-condensing	

For more information, please contact:



© 2020 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks used herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com

PART-6601SB03.20