



MET ONE 6000 Series Remote Airborne Particle Counters

High performing, reliable 0.3 μm sensitivity remote



Be Right™

The MET ONE 6000 series of ISO 21501 compliant counters are the most reliable, top performing, high sensitivity remote particle counters

High performance

0.3 µm sensitivity counter designed for ISO 21501 compliance for data accuracy and reliability

Easy to integrate with any FMS system

Multiple communication options such as Pulse, Serial Modbus, Ethernet, Wireless and Analog

Long Life Laser™ diode

Reduces the need for replacement parts and the overall cost of ownership



Faster trouble shooting and reduced downtime

Instrument can be diagnosed for flow, sensor, communication failures, and count alarm or count alert through built-in or external status light

Easy installation options

Reduces initial installation cost and subsequent reinstallation cost during routine maintenance/calibration work

Built-in flow sensor

Improves operational reliability during continuous monitoring

Port Sizes	Model 6003	0.32 cm (1/8-inch) ID inlet	0.64 cm (1/4-inch) ID outlet
	Model 6005	0.32 cm (1/8-inch) ID inlet	0.64 cm (1/4-inch) ID outlet
	Model 6013	0.64 cm (1/4-inch) ID inlet	0.64 cm (1/4-inch) ID outlet
	Model 6015	0.64 cm (1/4-inch) ID inlet	0.64 cm (1/4-inch) ID outlet
Flow Rate	Model 6003	0.1 CFM (2.83 LPM)	
	Model 6005	0.1 CFM (2.83 LPM)	
	Model 6013	1.0 CFM (28.3 LPM)	
	Model 6015	1.0 CFM (28.3 LPM)	
Sensitivity	Model 6003	0.3 µm at 0.1 CFM (2.83 LPM)	
	Model 6005	0.5 µm at 0.1 CFM (2.83 LPM)	
	Model 6013	0.3 µm at 1.0 CFM (28.3 LPM)	
	Model 6015	0.5 µm at 1.0 CFM (28.3 LPM)	
Range	Model 6003	0.3 µm to 10.0 µm at 0.1 CFM (2.83 LPM)	
	Model 6005	0.5 µm to 10.0 µm at 0.1 CFM (2.83 LPM)	
	Model 6013	0.3 µm to 10.0 µm at 1.0 CFM (28.3 LPM)	
	Model 6015	0.5 µm to 10.0 µm at 1.0 CFM (28.3 LPM)	
Channel Sizes		STANDARD 2-CHANNEL	OPTIONAL 4-CHANNEL*
	Model 6003	0.3 and 0.5 µm	0.3, 0.5, 1.0 and 5.0 µm
	Model 6005	0.5 and 5.0 µm	0.5, 1.0, 5.0 and 10.0 µm
	Model 6013	0.3 and 0.5 µm	0.3, 0.5, 1.0 and 5.0 µm
	Model 6015	0.5 and 5.0 µm	0.5, 1.0, 5.0 and 10.0 µm
Counting Efficiency	Model 6003	50% (± 20%) for 0.3 µm, (100% ± 10% at 1.5 times the minimum sensitivity). Fully complies with ISO21501-4.	
	Model 6005	50% (± 20%) for 0.5 µm, (100% ± 10% at 1.5 times the minimum sensitivity). Fully complies with ISO21501-4.	
	Model 6013	50% (± 20%) for 0.3 µm, (100% ± 10% at 1.5 times the minimum sensitivity). Fully complies with ISO21501-4.	
	Model 6015	50% (± 20%) for 0.5 µm, (100% ± 10% at 1.5 times the minimum sensitivity). Fully complies with ISO21501-4.	

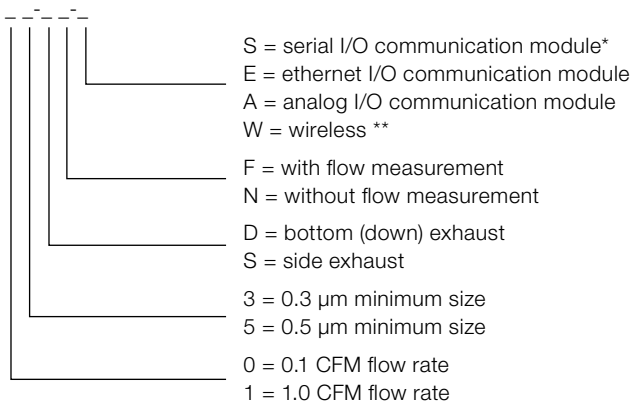
* Pulse models have two channels only

The MET ONE 6000 series remote airborne particle counter offers accurate and reliable continuous particle monitoring. With a sensitivity range of 0.3 or 0.5 to 10 µm, data accuracy and repeatability through ISO 21501 compliant calibration, the MET ONE 6000 series is designed to meet the specific needs of cleanroom operations within the pharmaceutical, semiconductor, hard disk drive and flat panel display industries.

With an array of communication and mechanical installation options, the MET ONE 6000 is easy to integrate with any facility monitoring system. These features also reduce downtime for instrument removal and reinstallation during routine calibration and preventative maintenance cycles. Built with Long Life Laser™ technology, the MET ONE 6000 offers industry leading diagnostic features that reduce troubleshooting time and related downtime costs.

Specifications

Light Source	Long Life Laser™ diode
Weight	0.82 kg (1.8 lbs)
Dimensions	13.56 w x 8.93 d x 12.06 h cm (5.34 x 3.52 x 4.75 inches)
Enclosure	304 stainless steel
Status Indicator	Multi-colored LED for normal status, count alarm, count alert, sensor failure, flow failure or communication failure
Power Requirements	9 to 28 VDC
Power Consumption	Serial and pulse units: 3.3 W; Ethernet unit: 4.3 W; Analog: 3.5 W; Wireless: 7.1 W. Maximum amperage requirement: 1 A
Operating Temperature	10 to 32 °C (50 to 90 °F), 5 to 95% relative humidity, non-condensing
Storage Temperature	-40 to 70 °C (-40 to 158 °F), 5 to 98% relative humidity, non-condensing
Signal Output Options	Pulse Analog 4–20 mA Serial RS-232 Serial RS-485 with Modbus RTU or FXB communication protocol Ethernet or Wireless with TCP/IP protocol
Vacuum Requirements	At least 40.6 cm (16 inches) Hg
Number of Size Channels	Standard 2 or Optional 4 (pulse units have 2 channels only)
Flow Control	Through critical orifice
Coincidence Loss	Model 6003/6005: 5% at 2,000,000 particles/ft ³ (70,600,000 particles/m ³) Model 6013/6015: (except Pulse) 5% at 400, 000 particle/ft ³ (14,000,000 particles/m ³)
False Count Rate	One or less in five minutes
Accessories Included	DIN Rail Mounting Kit, Isokinetic Probe, Male Phoenix Terminal Strip with Cover, Operator Manual
How To Order	20886



Optional Accessories***	Optional 4-Channel Setting (Pulse models have two channels only) Wall Plate Mounting Kit Utility Terminal Box Mounting Kit Purge Filters, 1.0 CFM 1/4 Tube and 0.1 CFM 1/8 Tube Isokinetic Probes, 0.1 CFM and 1.0 CFM Cable, Service Port Connector with RS-232 Converter Relative Humidity/Temperature Probe Brackets, for RH/Temp Probe and LED Indicator Stack/Isokinetic Probe Remote Indicator Light Assembly Cable Assembly, 3 m for Remote Indicator Light Assembly Cable Assembly, RJ-45 Adapter Power Supply Unit, 24 VDC @ 5.0 A, Universal Input for FMS
-------------------------	---

- * If a serial I/O model is ordered, a second part number 2088600-232, 2088600-485, or 2088600-PLS must be ordered that will identify the serial I/O configuration as RS-485, RS-232, or Pulse, respectively.
- ** Wireless models are available in limited number of countries. Contact your Hach sales representative for details.
- *** For a comprehensive parts and accessories list, please refer to the user manual.



Lit. No. 5236
G9 Printed in USA

© Hach Company 2009. All rights reserved.

In the interest of improving and updating its equipment, Hach Company reserves the right to alter specifications to equipment at any time.

At Hach, it's about learning from our customers and providing the right answers.

Keep it pure.

Make it simple.

Be right.

For current price information, technical support, and ordering assistance, contact the Hach office or distributor serving your area.

In the United States, contact:

HACH COMPANY World Headquarters
PO Box 608
Loveland, Colorado 80539 USA
Telephone: 800-866-7889
Fax: 970-461-3914
E-mail: customersupport@hach.com
www.particle.com

U.S. exporters and customers in Canada, Latin America, sub-Saharan Africa, Asia, and Australia/New Zealand, contact:

HACH COMPANY World Headquarters
PO Box 608
Loveland, Colorado 80539 USA
Telephone: 970-663-9760
Fax: 970-461-3914
E-mail: intl@hach.com
www.particle.com

In Europe, the Middle East, and Mediterranean Africa, contact:

HACH LANGE GmbH
Willstätterstraße 11
D-40549 Düsseldorf, GERMANY
Tel: +49 (0) 211 5288-0
Fax: +49 (0) 211 5288-143
E-mail: info@hach-lange.de
www.particle.com



Be Right™