



a | 5350 Lakeview Parkway South Dr. | Indianapolis, IN 46268  
o | 317.808.4200 or 800.352.4975  
f | 317.808.4395  
w | [beckman.com](http://beckman.com)

March 15, 2022

From: Kelsey Rospos, PCC Product Line Manager  
Re: MET ONE 3400 Obsolescence

Dear MET ONE Customer,

This is an obsolescence letter for the legacy **MET ONE 3400** air particle counter. This is planned in response to obsolescence of the main processor board and the availability of our new MET ONE 3400+ instruments. The MET ONE 3400 product went obsolete on July 1, 2021 in European countries that require RoHS 3 certification, and it went obsolete for the rest of the world on **December 31, 2021**. Legacy MET ONE 3400 instruments are replaced by new MET ONE 3400+ instruments.

We will continue to fully service and support the MET ONE 3400 instrument for a period of five (5) years after the latest obsolescence date up to December 31, 2026. During this time, calibrations, repairs, and reagents required for routine service will be available, however, breakdown repairs cannot be guaranteed if a required service component is made obsolete by our supplier.

Part numbers affected by the MET ONE 3400 obsolescence are listed at the end of this letter.

Please contact your local sales representative to find out more about current MET ONE 3400+ products. For technical questions or details regarding continued service and support, please contact your regional Customer Technical Support (CTS) team.

Thank you,

A handwritten signature in black ink that reads 'Kelsey Rospos'.

Kelsey Rospos  
Product Line Manager, Beckman Coulter Life Sciences  
[KMROSPOS@Beckman.com](mailto:KMROSPOS@Beckman.com)



a | 5350 Lakeview Parkway South Dr. | Indianapolis, IN 46268  
o | 317.808.4200 or 800.352.4975  
f | 317.808.4395  
w | [beckman.com](http://beckman.com)

Part numbers affected by the legacy MET ONE 3400 obsolescence include:

2088900-01	Air Particle Counter MO3413 0.3-10um
2088900-01-W	Air Particle Counter MO3413 0.3-10um WiFi
2088900-06	Air Particle Counter Met One 3413 0.3-10um Gas
2088900-07	Air Particle Counter MO3413 0.3-10um gas WiFi
2088900-05	3413 PARTICLE CNTR, 1CFM W/ASHRAE CAL
2089494-01	MET ONE 3400 PAPERLESS PDF OPTION
2088900-03	Air Particle Counter MO3415 0.5-25um
2088900-02	Air Particle Counter MO3415 0.5-10um
2088900-02-W	Air Particle Counter MO3415 0.5-10um WiFi
2088900-08	Air Particle Counter Met One 3415 0.5-10um Gas
2088900-09	Air Particle Counter MO3415 0.5-10um Gas WiFi
2088900-11	Air Particle Counter MO3423 0.3-10um 50l
2088900-11-W	Air Particle Counter MO3423 0.3-10um 50l WiFi
2088900-12	Air Particle Counter MO3425 0.5-10um 50l
2088900-12-W	Air Particle Counter MO3425 0.5-10um 50l WiFi
2088900-13	Air Particle Counter MO3425 0.5-25um 50l
2083049-11	A3432 MANIFOLD, 115/100VAC w/CONTROLLER
2088900-32	Air Particle Counter MO3445 0.5-10um 100l
2088900-32-W	Air Particle Counter MO3445 0.5-10um 100l WiFi
2088900-03-W	Air Particle Counter MO3415 0.5-25um WiFi
2088900-13-W	Air Particle Counter MO3425 0.5-25um 50l WiFi
2089473-01	MODEL 3413 PORT-AIR PARTICLE COUNTER
2089473-02	Model 3413 Port-Air Particle Counter
2089474-01	Model 3415 Port-Air Particle Counter
2089474-02	Model 3415 Port-Air Particle Counter
2089474-03	3415 W/CHNL SIZES 0.5 to 25 MICRONS
2089475-01	Model 3423 Port-Air Particle Counter
2089475-02	Model 3423 Port-Air Particle Counter
2089476-01	Model 3425 Port-Air Particle Counter
2089476-02	Model 3425 Port-Air Particle Counter
2089477-01	Model 3445 Port-Air Particle Counter
2089477-02	Model 3445 Port-Air Particle Counter
B97964	METONE 3413 AIR PARTICLE COUNTER
2083049-01	A2432 Manifold, 115/100V w/Controller
2083049-02	A2432 Manifold, 220V w/Controller