

# MET ONE HHPC+ Series Handheld Particle Counters

---



**Simple**

**Fast**

**MET ONE**

# MET ONE HHPC+ series handheld particle counter.

Counting 00:00:14 Infinite	
Location 1	Counts/ft <sup>3</sup>
Size	Cumulative
0.3 $\mu$ m	155.8
0.5 $\mu$ m	77.9
1.0 $\mu$ m	13.0
2.0 $\mu$ m	13.0
5.0 $\mu$ m	0.0
10.0 $\mu$ m	0.0
2.170 L	

Stopped Infinite	
Location 1	Counts/L
Size	Cumulative
0.5 $\mu$ m	0.0
5.0 $\mu$ m	0.0



## Routine monitoring

The easy-to-read high resolution display is fully configurable, presenting only the data you want in a clear large font.

## Light weight and comfortable for those hard to reach HEPA filter leaks

At only 1.5 lbs the MET ONE HHPC+ slim design allows single handed operation. Perfect for effortless troubleshooting of filter leaks in mini-environments and workstations.

## Cleanroom classification

Quickly check or validate ISO Class 5 (FED STD Class 100) or higher cleanrooms and controlled environments.

## Visual particle trending

Save analysis time with on-the-spot process profiling to pinpoint particle events.



# Easy data download

## USB Cable • Memory Stick • Ethernet

### Easy network connectivity

Download to Excel® via Ethernet using industry standard browsers (Internet Explorer®, Safari®, Firefox®, etc.) CFR 21 Part 11 compliant.

### Use a memory stick – grab the data and go!

Plug a USB memory stick in the MET ONE HHPc+ to transfer your particle data direct to a PC.

### USB cable direct to PC

Plug it in like a camera, your data appears in Excel®.

### Leave your MET ONE HHPc+ in the cleanroom on the charge cradle

Your handheld is fully charged, and ready for use at all times. The unique cradle provides PC access to particle counting data via Ethernet or USB cable.



Sample Volume	Sample Units	Count Mode	Concentration	Counts normalized	
2.83 L	Cum.	Counts/L	0.3µm	0.5µm	
2.83 L	Cum.	Counts/L	2	1	
2.83 L	Cum.	Counts/L	0	0	
2.83 L	Cum.	Counts/L	0	0	
2.83 L	Cum.	Counts/L	0	0	
2.83 L	Cum.	Counts/L	0	0	
2.83 L	Cum.	Counts/L	0	0	
2.83 L	Cum.	Counts/L	0	0	
0:01:00	2.83 L	Cum.	Counts/L	0	0
0:01:00	2.83 L	Cum.	Counts/L	0	0
0:01:00	2.83 L	Cum.	Counts/L	2	1
0:01:00	2.83 L	Cum.	Counts/L	0	0
0:01:00	2.83 L	Cum.	Counts/L	1	1
0:01:00	2.83 L	Cum.	Counts/L	0	0
0:01:00	2.83 L	Cum.	Counts/L	0	0
0:01:00	2.83 L	Cum.	Counts/L	0	0
0:01:00	2.83 L	Cum.	Counts/L	1	1



Transfer cleanroom particle data by easy-to-use USB Memory Stick and Ethernet electronic data transfer, you'll be spending less time in the cleanroom and more time on the results!

## Specifications

<b>Particle Size Range</b>	Size channels per ISO 14644-1 (FS 209E) industry standard. HHPC 6+ and HHPC 3+ (0.3µm version) - 0.3µm - 10µm HHPC 2+ and HHPC 3+ (0.5µm version) - 0.5µm - 5µm
<b>Number of Channels</b>	HHPC 6+ Channel 1 fixed 0.3µm. Up to 5 additional size channels: 0.5µm, 1.0µm, 2.0µm, 5.0µm, 10.0µm. HHPC 3+ (0.3µm) Channel 1 fixed 0.3µm. Up to 2 additional size channels: 0.5µm, 1.0µm, 2.0µm, 5.0µm, 10.0µm. HHPC 3+ (0.5µm) Channel 1 fixed 0.5µm. Up to 2 additional size channels: 1.0µm, 2.0µm, 5.0µm. HHPC 2+ Channel 1 fixed 0.5µm. 1 additional size channel: 1.0µm, 2.0µm, 5.0µm.
<b>Weight</b>	Ultra-lightweight at 1.5 lbs for comfortable handheld use.
<b>Flow Rate</b>	2.83 lpm. Take ISO 14644-1 class 5 compliant 0.3µm samples in a minute
<b>Communications</b>	Access particle counting data direct to PC via Ethernet and USB cable, or use USB memory stick. No device drivers or software required.
<b>Display</b>	Large 3.5" high resolution 320 x 240 color for super clear viewing with intuitive icons for easy navigation.
<b>Display Configuration</b>	User selectable size channels with super-large font options.
<b>Display Data Modes</b>	Traditional tabular particle counts or trend graph for easy particle profiling.
<b>Enclosure</b>	High impact polycarbonate ABS for exceptional durability.
<b>Dimensions</b>	10.7" x 3.9" x 2.1" Contoured for convenient single handed operation.
<b>Battery Duration</b>	Over 10 hours with typical use model. Minimum 5.5 hours continuous sampling.
<b>Battery Charging</b>	3.5 hours
<b>Data Storage</b>	10,000 records for the most demanding applications. (5,000 records when secure CFR 21 Part 11 compliant files enabled)
<b>Counting Efficiency</b>	HHPC 3+ (0.3µm version) and HHPC 6+ 50% @ 0.3µm and 100% for particles > 0.45µm (per ISO 21501) HHPC 2+ and HHPC 3+ (0.5µm version); 50% @ 0.5µm and 100% for particles > 0.75µm (per ISO 21501)
<b>Zero Count Level</b>	1 count / 5 minute (per JIS B9921)
<b>Concentration Limits</b>	4,000,000 / ft <sup>3</sup> @ 10% coincidence loss HHPC 6+, HHPC 3+ (0.3µm), 2,000,000 / ft <sup>3</sup> @ 10% coincidence loss HHPC 2+, HHPC 3+ (0.5µm)
<b>Count Modes</b>	Raw counts, N/ft <sup>3</sup> , N/m <sup>3</sup> , N/L in Cumulative or Differential mode
<b>Security</b>	Administrator password controlled (optional) CFR 21 Part 11 compliant Excel compatible files.
<b>Alarms</b>	User selected particle channels and limits
<b>Operating Environment</b>	50°F to 104°F (10°C to 40°C) / < 95% non-condensing
<b>Storage Environment</b>	14°F to 122°F (-10°C to 50°C) / Up to 98% non-condensing
<b>Accessories Included</b>	Power supply, Universal 110-240V Zero count inlet-filter and zero count filter adaptor 6' USB cable Cradle for ease of charging and convenient USB and Ethernet communications (HHPC 6+ only) Carry case (HHPC6+ and HHPC3+ models only)

For more information on the MET ONE HHPC+ handheld particle counter, please visit us at [www.particle.com](http://www.particle.com).

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

