

MET ONE HHPC+ Series Handheld Particle Counters

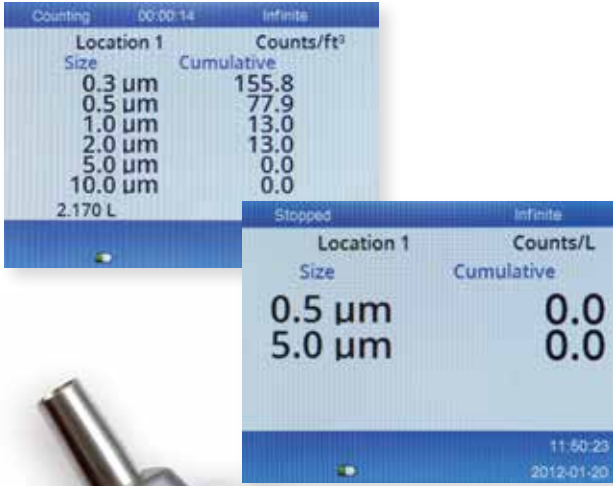


Simple

Fast

MET ONE

MET ONE HHPC+ series handheld particle counter.



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

CFR 21 Part II secure 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 HHPG+ 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 HHPG+ 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.



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

| | |
|------------------------------|---|
| Particle Size Range | 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 |
| Number of Channels | 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. |
| Weight | Ultra-lightweight at 1.5 lbs for comfortable handheld use. |
| Flow Rate | 2.83 lpm. Take ISO 14644-1 class 5 compliant 0.3 µm samples in a minute |
| Communications | Access particle counting data direct to PC via Ethernet and USB cable, or use USB memory stick. No device drivers or software required. |
| Display | Large 3.5" high resolution 320 x 240 colour for super clear viewing with intuitive icons for easy navigation. |
| Display Configuration | User selectable size channels with super-large font options. |
| Display Data Modes | Traditional tabular particle counts or trend graph for easy particle profiling. |
| Enclosure | High impact polycarbonate ABS for exceptional durability. |
| Dimensions | 10.7" x 3.9" x 2.1" Contoured for convenient single handed operation. |
| Battery Duration | Over 10 hours with typical use model. Minimum 5.5 hours continuous sampling. |
| Battery Charging | 3.5 hours |
| Data Storage | 10,000 records for the most demanding applications. (5,000 records when secure CFR 21 Part 11 compliant files enabled) |
| Counting Efficiency | 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) |
| Zero Count Level | 1 count / 5 minute (per JIS B9921) |
| Concentration Limits | 4,000,000 / ft ³ @ 10% coincidence loss (HHPC 6+ and 3+), 2,000,000 / ft ³ @ 10% coincidence loss (HHPC 2+) |
| Count Modes | Raw counts, N/ft ³ , N/m ³ , N/L in Cumulative or Differential mode |
| Security | Administrator password controlled (optional) |
| Alarms | User selected particle channels and limits |
| Operating Environment | 50°F to 104°F (10°C to 40°C) / < 95% non-condensing |
| Storage Environment | 14°F to 122°F (-10°C to 50°C) / Up to 98% non-condensing |
| Accessories Included | 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.



CLASS 1 LASER PRODUCT
Contact manufacturer for complete compliance details



Contact us to place an order, request information or receive technical support.



Service packages and extended warranty up to 5 years.



On site technical support.



Seminars and workshops: Practical and hands on training.



Regular customer information by post and email.



www.particle.com up to date and secure, with downloads, information and e-shop.

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