

CHARACTERIZED  
*by ingenuity.*

# CERTIFIED CALIBRATION

Ensure count accuracy and reproducibility



## When accuracy is critical, don't trust your calibration to just anyone.

Beckman Coulter is the only service provider with GMP-trained service technicians equipped with the CoreCAL system and test standards required to perform ISO 21501-4 calibrations for your MET ONE equipment.

## Minimize your compliance risk with CoreCAL.

The CoreCAL system from Beckman Coulter is routinely monitored and updated for guaranteed accuracy and alignment with ISO 21501-4 requirements.

Only factory-trained Service Engineers authorized to use a licensed version of CoreCal can identify and correct system deficiencies for audit-ready compliance.

**If your service engineer isn't using CoreCAL and the Beckman Coulter 3400STD transfer standard, it's impossible to complete the following mandatory ISO 21501-4 test parameters:**

- Set / update channel threshold voltages
- Noise test
- Flow rate calibration
- Resolution
- Counting efficiency and accuracy
- @ 50% / @ 100%
- Maximum concentration
- Peak-to-valley
- Sample time error
- False count rate



The CoreCAL system is exclusive to Beckman Coulter service engineers and authorized agents.

## Getting it right matters.

Call or email to learn more: 800-369-0333 | [particlecountingsupport@beckman.com](mailto:particlecountingsupport@beckman.com).

<sup>1</sup>If working with an authorized third-party, ensure CoreCAL version 4.5 is being used to calibrate your system (October 2017).

## Complete peace of mind

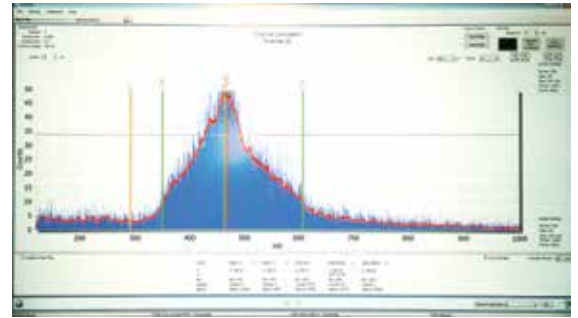
More than 150 engineers across North America are factory trained to calibrate Beckman Coulter instruments to the highest and most exacting standards.

### Audit-quality calibration services for your MET ONE particle counting equipment

- Certified calibration to maintain original system efficiency—only available with the CoreCAL system
- Service continuity and peace of mind with 6-month, annual or customized service options—performed in accordance with ISO 21501-4 and ISO 17025<sup>2</sup> requirements using NIST-traceable equipment and particle media
- ‘As Found’ data available
- Counting resolution and efficiency certificate issued with each instrument
- Expertly-trained service engineers knowledgeable in quality guidelines enforced by the FDA and ISO/IEC 17025
- Engineer technical training records available on request
- Onsite or return-to-depot calibration service options
- Five depots across North America
- Discounted parts and labor<sup>3</sup>

MET ONE 3400, 6000 and 7000 Series systems are designed and manufactured to comply with ISO 21501-4.

If you're older MET ONE equipment, contact us to learn more about upgrading to an ISO 21501-4 compliant counter for improved accuracy and repeatability.



The CoreCAL system is exclusive to Beckman Coulter service engineers and authorized agents.

## Getting it right matters.

Call or email to learn more: 800-369-0333 | [particlecountingsupport@beckman.com](mailto:particlecountingsupport@beckman.com).

### Additional Beckman Coulter services to ensure the integrity of your results and reduce costly downtime:

- IQ/OQ services for expert qualification, compliance and validation of newly installed, moved or modified equipment
- Flexible service and repair programs to avoid the high cost of instrument failure
- Customized service plans designed to unique customer requirements
- Remote phone support
- Instructor-led and online operator training programs
- Application and workflow support

<sup>2</sup>ISO 17025 calibrations performed at our ISO 17025-accredited Loveland, CO depot.

<sup>3</sup>Parts and labor discount applies to services performed during coverage period. For the best performance possible, Beckman Coulter recommends using original replacement parts for the repair and maintenance of your MET ONE or HIAC particle counter.



Only Beckman Coulter certified and trained Field Service Engineers and agents with the Beckman Coulter CoreCAL 3 calibration tool and reference instrument are able to perform fully ISO 21501 compliant calibrations on MET ONE air particle counters.

© 2017 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks mentioned 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](http://beckman.com)

PART-3167FLY11.17