

CHARACTERIZED  
*by ingenuity.*

# CERTIFIED CALIBRATION

Ensure count accuracy and reproducibility



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

When it comes to an **ISO 11171 or 21501-3** calibration for your HIAC instrument, no one knows the technology better than our GMP-trained service engineers.

For added protection, HIAC instrument are equipped with built-in safeguards and security features designed to protect the integrity of your calibration routines and system performance. With controlled access, our service engineers are fully equipped to properly calibrate your system—including making adjustments to correct any system deficiencies, for an audit-quality calibration.

**If your service engineer isn't factory-trained and using CoreCAL software, HIAC-embedded security passcodes & NIST-traceable media, your HIAC liquid particle counter may not be calibrated to mandatory test parameters.<sup>1</sup>**



- HIAC PODS+ & 8011+ liquid particle counters are designed & manufactured to be calibrated in full accordance with the ISO 11171 calibration standard.
- Oil particle counter calibrations are performed to provide reporting capability to ISO4406, SAE AS4059, NAS1638, GOST, NAVAIR, DEF STAN & ASTM reporting standards.
- Fully accredited ISO 11171 calibrations are performed for customers requiring calibration in accordance to all clauses & annexes pertaining to the ISO 11171.
- HIAC 9703+ instruments are designed & manufactured to comply with ISO 21501-3 & USP requirements for pharmaceutical manufacturing, microelectronics, precision machines & medical operations.
- For counting accuracy, pharmaceutical particle counter calibrations are performed using particle count standard reference validation (USP) using NIST-traceable latex spheres/standards.
- Detailed calibration certificate & supporting data package included.

## Audit-quality calibration services for your HIAC particle counting equipment.

- Certified calibration certificate and supporting data package
- Service continuity & peace of mind with six-month, annual or customized service options—performed in accordance with applicable ISO standards<sup>2</sup> & requirements using NIST-traceable equipment & calibration media
- Factory-trained service engineers knowledgeable in quality guidelines enforced by the FDA & ISO
- Engineer technical training records available on request
- Onsite or return-to-depot calibration service options
- Five depots across North America
- Discounted parts & labor<sup>3</sup>

### 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 & validation of newly installed, moved or modified equipment
- Flexible service & repair programs to avoid the high cost of instrument failure
- Customized service plans designed to unique customer requirements
- Remote phone support
- Instructor-led & online training options
- Applications & workflow support

## Repeatability, advanced sensitivity & resolution top the list of the most important factors for today's particle system operators.

What do particle size analyzer users have to say? Lab Manager, September 2017.

<sup>1</sup> CoreCAL software is routinely monitored and updated for guaranteed accuracy and alignment with ISO 11171 and 21501 requirements. If working with an authorized third-party, please ensure CoreCAL version 4.5 is being used to calibrate your system (December 2017).

<sup>2</sup> ISO 17025 calibrations are performed at our ISO 17025-accredited facility.

<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 HIAC particle counter.



© 2018 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-3276FLY02.18