



Global Service Center – Services & Pricing

A service request must be submitted via the online form. An RA number and shipping/decontamination form will be emailed once your online request is received and processed. Instruments received without an RA number & signed decontamination form will be returned without service.

MET ONE AIR CALIBRATION

USD \$

A-CAL - Models 2100 • 2200 • 237H • 3411	1,975.00
B-CAL - Models 3413 • 3415 • 3423 • 3425 • 3445	1,550.00
C-CAL - Models: 237 A/B • 2400/2408 • 33XX • HHPC (For HHPC+ series - see C1-CAL option)	1,410.00
C1-CAL - Models: HHPC+ • 6000/7000 series	1,410.00
D-CAL - Models: 227 A/B • 228 • 229 • R4700 • R4800 • R4900	1,300.00
E-CAL - Models: 3413+, 3415+, 3423+, 3425+, 3445+	3,295.00
ADD ISO 21501-4 CAL - Includes Counting Efficiency & Resolution <i>Available For Models: 3413, 3415, 3423, 3425, 3445, 6000 & 7000 series</i>	640.00
ADD ISO 17025 CAL - <i>Available For Models: 3400 series (except 3411), HHPC+ series, 6000 & 7000 series</i>	640.00
Relative Humidity & Temperature Probe calibration	765.00

HIAC LIQUID CALIBRATION

PSL WATER CALIBRATION HRLD SENSORS (Includes Data)	2,075.00
ADD ISO 17025 CAL - HRLD sensor PSL ONLY	640.00
PSL WATER MICROCOUNT CALIBRATION MC-05, MC-100 (Includes Data)	2,120.00
MTD OIL CALIBRATION HRLD models: 100 • 150 • 600	2,365.00
ULTRAPURE UP-100	3,130.00
ROC VERIFICATION (Pass/Fail Verification only)	750.00
PODS OR PODS+	2,365.00
GLYCOUNT PORTABLE	2,620.00

ANATEL TOC CALIBRATION

CALIBRATION (includes calibration, battery, lamp, filter) - Models: A1000, PAT700, A643, TOC600	5,240.00
CALIBRATION only - Models: A1000, PAT700, A643, TOC600	4,500.00
SYSTEM SUITABILITY - Models: A1000, PAT700, A643, TOC600	1,785.00
AS RECEIVED DATA - Models: A1000, PAT700, A643, TOC600	615.00
CALIBRATION RESISTOR	360.00

REPAIR - ALL PRODUCTS

EVALUATION FEE	630.00
----------------	--------

SERVICE PRIORITY SELECTION

PRIORITY - Next in line for bench calibration or evaluation of repair. Overnight return shipping for customers within the United States. Standard return shipping for International customers.	1,250.00
--	----------



© 2024 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
2024-AMER-EN-105236-v1

