



# Global Service Center – Services & Pricing

A service request must be submitted via the online form. A return authorization (RA) number and shipping/decontamination form must be assigned before shipping to Beckman Coulter Life Sciences. Instruments received without an RA number and signed decontamination form will be returned without service.

## MET ONE AIR CALIBRATION

A-CAL - Models 2100 • 2200 • 237H • 3411	\$ 1,818
B-CAL - Models 3413 • 3415 • 3423 • 3425 • 3445	\$ 1,489
ADD ISO 21501-4 CAL (UNDER B-CAL OPTION)	\$ 588
C-CAL - Models 237 A/B • 2400/2408 • 33XX • HHPC/HHPC+ • 6000/7000	\$ 1,348
D-CAL - Models 227 A/B • 228 • 229 • R4700 • R4800 • R4900	\$ 1,030
E-CAL - Models 3413+ • 3415+ • 3423+ • 3425+ • 3445+ (Includes ISO-21501-4)	\$ 3,235
Relative Humidity & Temperature Probe calibration	\$ 765

## HIAC LIQUID CALIBRATION

PSL WATER CALIBRATION HRLD SENSORS (Includes Data)	\$ 2,059
PSL WATER MICROCOUNT CALIBRATION MC-05, MC-100 (Includes Data)	\$ 2,117
MTD OIL CALIBRATION HRLD 100 • 150 • 600	\$ 2,177
ULTRAPURE UP-100	\$ 2,882
ROC VERIFICATION (Pass/Fail Verification only)	\$ 384
PODS OR PODS+	\$ 2,177
GLYCOUNT PORTABLE	\$ 2,412

## ANATEL TOC CALIBRATION

CALIBRATION A1000 S10, S20, S20P	\$ 2,305
CALIBRATION A1000 XP & XPS	\$ 2,470
CALIBRATION A643-P & A643-S	\$ 2,006
CALIBRATION TOC600	\$ 1,835
CALIBRATION PAT700	\$ 3,152
SYS-SUIT A1000 S10, S20, S20P, XP & XPS	\$ 3,270
SYS-SUIT A643	\$ 1,495
SYS-SUIT PAT700 & TOC 600	\$ 1,365
AS RECEIVED DATA A1000 S10, S20, S20P, XP, XPS, A643-P, A643-S	\$ 1,160
AS RECEIVED DATA PAT700 & TOC 600	\$ 1,160
UV LAMP A1000 S10, S20, S20P	\$ 1,070
UV LAMP A1000 XP & XPS	\$ 1,656
UV LAMP A643 & TOC600	\$ 909
UV LAMP PAT700	\$ 1,058
CALIBRATION RESISTOR	\$ 330
C80 REPAIR SERVICE	\$ 1,717

## REPAIR - ALL PRODUCTS

FLOW/PUMP ISSUE (Air units only)	\$ 5,818
HRLD SENSOR ISSUE	\$ 4,411
INSTRUMENT WILL NOT TURN ON	\$ 4,017
DISPLAY SCREEN REPLACEMENT	\$ 1,891
CALIBRATION ALARM (Calibration not included)	\$ 6,845
EVAL FEE / REPAIR OTHER - (Anatel/TOC only select Evaluation Option)	\$ 598

## SERVICE PRIORITY SELECTIONS

<b>RUSH</b> - Next on bench service with overnight return shipping (Note: available for calibration service or evaluation of repair only) <b>Excludes: Anatel, UP100, CCPHA, 3400 STD, MC-100/200 &amp; INT'L units</b>	\$ 1,000
---	----------



©2023 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](https://beckman.com)  
2023-AMER-EN-100728-v1