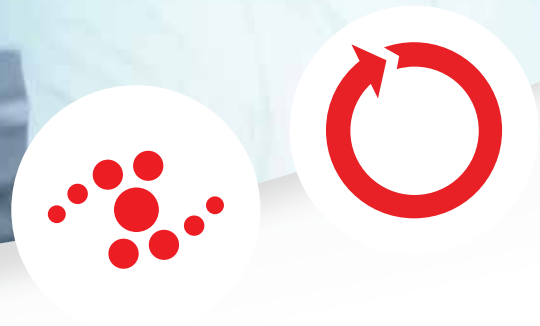


BRILLIANCE
at every turn.

BECKMAN COULTER THE POWER BEHIND YOUR RESEARCH



When you acquire a centrifuge or particle characterization device, we believe it's just the beginning of your relationship with us. That's why behind every Beckman-built instrument, we've mobilized a world-class Service team with the knowledge and experience to optimize instrument performance and maximize critical system uptime to keep your research going.

We make it easy to protect your investment during and beyond the warranty period. With flexible service plan options and features, you choose the level of support to best meet your needs and budget.



Find the right plan for you

Plan Benefits	Beckman Coulter Service Plans				
	WARRANTY PERIOD		POST-WARRANTY PERIOD		
	Warranty	Start-Up Care Premium Services	Prevention Plus	Protective	Comprehensive
Certified parts, labor	●	●	20% Off	●	●
Travel expenses ¹	●	●	20% Off	●	●
Annual preventive maintenance		1 PM	1 PM	1 PM	1 PM
Annual health check ²		●		●	●
Onsite response time guarantee ³	●	2 DAYS		3 DAYS	2 DAYS
Remote technical support	●	●	●	●	●
Software & Engineering updates ⁴		●	20% Off	●	●
Annual basic operator training ⁵		●	20% Off	●	●
Operational qualification (OQ)		AVAILABLE		AVAILABLE	AVAILABLE
Application Support ⁶	AVAILABLE	AVAILABLE	AVAILABLE	AVAILABLE	●
PROService Remote Connectivity ⁷	●	●	●	●	●
Relocation support		●			●



Need help choosing the right protection plan?

We're here to help:

(800) 742-2345, Option 3 | protection@beckman.com

1. Within 100 miles of a Beckman Coulter Service hub.
2. Proactive service check performed by a Beckman Coulter field service engineer to address any system or service performance issues.
3. Priority response guaranteed. Contact your local service representative for guaranteed response times based on your location.
4. Reliability updates include hardware, software and instrument modifications to recommended levels.
5. Services performed at your facility by a Beckman Coulter field service engineer.
6. Beckman Coulter application scientist will provide basic onsite support, at customer request. Advanced application support packages available. Contact your Beckman Coulter sales representative for more information.
7. PROService-enabled instruments include the Navios, Navios EX, Gallios, AQUIOS and Biomek i-Series liquid handling systems.



© 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

MULTI-3467FLY02.18