



LEVEL **3**

Preventive Plus 1

Uptime peace of mind.

Right across the Service value chain, the challenges remain the same. Achieving more from less and for greater value. Beckman Coulter Life Sciences Service understands that proactivity is the key to uptime peace of mind.

Prevent catastrophic failures. Increase workflow productivity.

Sometimes, prevention is the first, the only imperative – the dominant force in managing the risk of major instrument failure and the impact on workflow productivity. But sometimes budget limitations drive that most pressing commercial choice – how can my business protect its most valuable investment and still maintain, or even increase, process throughput in terms of speed, accuracy and sensitivity.

Beckman Coulter Life Sciences Service provides the flexibility to help you make that choice. At the heart of our reputation sits a promise. To achieve maximum uptime without compromising our customers workflows and the productivity challenges they face. And providing the peace of mind every laboratory or facility requires.

Our Plus 1 plan delivers that promise in full.



Preventive Plus 1 Service Plan

The benefits:

- A proactive approach to prevent catastrophic failures and increase your workflow productivity.
- This program includes a single, annual preventive maintenance visit and a discount for other repair calls. Ideal if you have a limited budget that needs to ensure dependable operation and minimal downtime.

Comprehensive

Protective

Preventive Plus 2

Preventive Plus 1

Connect

Start-up Care

Full Care Program





Beckman Coulter Care Packages Overview:

	Year 1		Year 2+				
	Warranty Standard coverage & phone support	LEVEL 1 Start-Up Care Peace of mind when an instrument is new	LEVEL 2 Connect Phone support to assist self servicing of equipment	LEVEL 3 Preventive Plus 1 A proactive approach to preventive care	LEVEL 4 Preventive Plus 2 Greater flexibility and options based on need	LEVEL 5 Protective Combine a preventive plan with repair service	LEVEL 6 Comprehensive Full proactive care for optimized performance
Certified parts, labor and travel ¹	●	●	5% Off	20% Off	20% Off	●	●
Annual Preventive Maintenance ²		1 PM		1 PM	2 PM	1 PM	1 PM
On-site response time guarantee ³		3 Days				3 Days	2 Days
Telephone/email technical support	●	●	●	●	●	●	●
Annual health check visit		●				●	●
Reliability Updates ⁴		●	5% Off	20% Off	20% Off	●	●
Annual basic operator on-site training ⁵		4 Hours	○	○	○	2 Hours	4 Hours
Operational Qualification Level 3		○				○	○
Full Care Leasing Program						○	○

● Included ○ Optional

¹Within 100 miles of a Beckman Coulter service hub. Ask your local service representative about locations.

²For Preventive Plus 2, you can choose 2 PMs, 1 PM and 1 repair call, or 2 repair calls. Labor and travel costs are covered for repair calls. Parts cost will be billed separately.

³Typical for most areas. Contact your local service representatives for guaranteed response times in your location.

⁴Reliability updates include hardware upgrades, software upgrades, and instrument modifications to recommended level.

⁵Training doesn't apply to Flow Cytometry instruments.

Product lines covered by these plans include Centrifuge, Particle Characterization, Automation, and Flow Cytometry. For Centrifuges, Start-Up and Comprehensive plans include rotor inspection and rotor service.

We've designed flexible options to suit the needs of any lab or operation. No matter which package you select, you're making a choice to strengthen your investment.

Need help with the decision or more details? Contact us:

1-800-742-2345 or visit www.beckman.com



© 2016 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