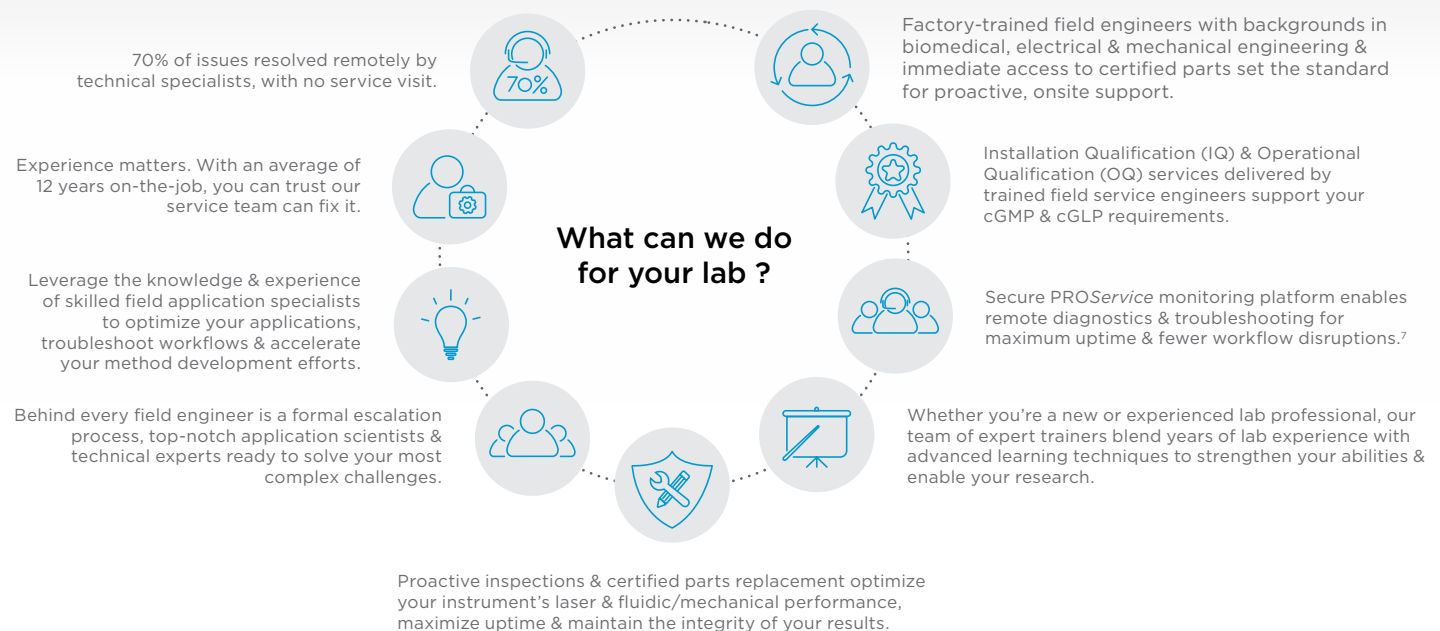


EVERY  
*event matters.*

# BECKMAN COULTER THE POWER BEHIND YOUR WORKFLOWS



With every Beckman Coulter service plan, you can count on original service parts and expert support to ensure your flow cytometer is operating within specification. Exchange downtime and workflow disruption risks for optimized system performance and quality results with a range of protection plans. With flexible 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 <sup>1</sup>	●	●	20% Off	●	●
Annual preventive maintenance		1 PM	1 PM	1 PM	1 PM
Annual health check <sup>2</sup>		●		●	●
Onsite response time guarantee <sup>3</sup>		●		●	●
Remote technical support	●	●	●	●	●
Software & Engineering updates <sup>4</sup>		●	20% Off	●	●
Annual basic operator training <sup>5</sup>		●	20% Off	●	●
Operational qualification (OQ)	AVAILABLE	AVAILABLE	AVAILABLE	AVAILABLE	AVAILABLE
Application Support <sup>6</sup>	●	●	AVAILABLE	●	●
PROService Remote Connectivity <sup>7</sup>	●	●	●	●	●
Relocation support		●			●



Need help choosing the right protection plan?

We're here to help:

(800) 742-2345, Option 3 | [protection@beckman.com](mailto: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](http://beckman.com)

FLOW-3433FLY01.18