

# Full Care Lease-to-Own Program

Combines the Power to Purchase and Service at a Fixed Monthly Rate



## Increase Your Lab Efficiency, Not Your Lab Spend

Full Care is a flexible leasing option that combines the latest technology from Beckman Coulter Life Sciences with comprehensive service support including parts, on-site support and maintenance all with zero upfront investment and one fixed monthly payment.

## Eligible for a wide variety of Beckman Coulter instruments:

All products, with the exception of hand-held Particle Counters, Microfuges, and Clinical Flow Analyzers.



## Benefits of Full Care Leasing Program:

- **Direct Lease from Beckman Coulter**  
Full Care Program is a direct financing option from Beckman Coulter to you. No need to work with a 3rd party leasing company.
- **Flexible Financial Solution**  
Cost-effective lease that makes critical instrumentation more affordable. Structured to fit your budget requirements.
- **Increased Efficiency**  
Improved workflow effectiveness and uptime with the latest technology from Beckman Coulter combined with comprehensive service and ongoing maintenance.

## Ownership Following Full Care Lease Term:

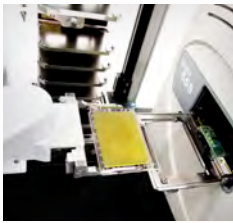
- **Purchase the instrument** for \$1 buyout price to acquire the title. A service agreement would need to be purchased for continued ongoing maintenance and coverage.



# Beckman Coulter Care

## Full Care Lease-to-Own Program

With one low monthly fixed cost, your Full Care Package includes:



### Equipment

A brand new, state-of-the-art instrument from Beckman Coulter



### Service

Locked in service rate for five years that includes on-site support, preventive maintenance and repairs including labor and travel



### Parts

Repair or replacement of certified and tested factory parts

### Full Care Lease-to-Own Program

available with one of the following service options:

	Protective Combine a preventive plan with repair service	Comprehensive Full proactive care for optimized performance
Equipment Lease	●	●
On-site Repair	●	●
Preventive Maintenance	●	●
Software & Engineering Updates	●	●
Phone Support	●	●
Calibration Service <small>(For Particle Counting models only)</small>	1 Visit	2 Visits
On-site Response Time Guarantee	3 Days	2 Days
Annual Basic Operator On-site Training	2 Hours	4 Hours
Rotor Inspection		●

● Included

Instrument coverage and response time vary by region.

To Learn More Please Contact:



Only monthly payments are accepted. Early termination requires payment of all past due invoices and the remaining balance on the contract. MOD Category 1 & 2 are included as updates.

© 2020 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service names mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. Echo is a trademark or registered trademark of Labcyte 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](http://beckman.com)

MULTI-7952SB09.20