



ALLEGRA V-15R BENCHTOP CENTRIFUGE

SMALL FOOTPRINT. BIG IMPACT.



BRILLIANCE
at every turn.

 **BECKMAN
COULTER**
Life Sciences

ALLEGRA V-15R FEATURES



Ergonomic Design

Soft-close lid for stress-free closing

Enhanced Safety

Automatic Rotor ID, Rotor Cycle Count and Imbalance Detection

Versatility

Broad rotor and adapter selection to spin commonly used tubes, bottles and plates

Simple and Intuitive

Intuitive, easy-to-read backlit display

Quiet Operation

Quiet operation of ≤ 56 dBA, at max speed, reduces noise pollution

Compact Footprint

Small footprint that can easily fit on the bench

Flexibility of using an optional mobile cart to completely free up the bench

↻ Versatile Performance Meets Simplicity

The Allegra V-15R centrifuge is designed to be a workhorse in the lab. Operation is simple and seamless thanks to an easy-to-use interface and ergonomic design. And with 10 rotor options at your fingertips, you will be ready and able to support multiple workflows with just one instrument.



Enhanced Safety

Automatic rotor recognition and cycle count, built-in imbalance sensor and biosafe rotors and covers protect the centrifuge, your samples and most importantly—you.



User Convenience

Operation is simple and seamless with 50 programmable runs and an easy-to-use interface with illuminated buttons to quickly visualize the status of a run.



Ergonomically Designed

No need to exert unnecessary force with the soft-close lid and low-profile design. Quiet operation of ≤ 56 dBA at max speed makes this centrifuge easy to work next to on the bench.

Greater Temperature Control

With the Allegra V-15R centrifuge, you'll be able to maintain a consistent temperature, allowing you to spin faster and longer without overheating your samples. All while Rapid Temp and Eco Mode allow for even more efficiency and functionality.



Rapid Temp

You no longer need to confirm the speed and time for pre-cooling. With Rapid Temp, the chamber can be pre-cooled to a defined temperature, ensuring optimal sample protection.



ECO Mode

To reduce energy consumption and extend compressor life, ECO Mode deactivates the compressor after a pre-selected duration of up to 8 hours in 30-minute increments. Say goodbye to condensation pooling.



Refrigeration

Uses CFC-free refrigerant and allows you to set a temperature from -10 to 40 °C.

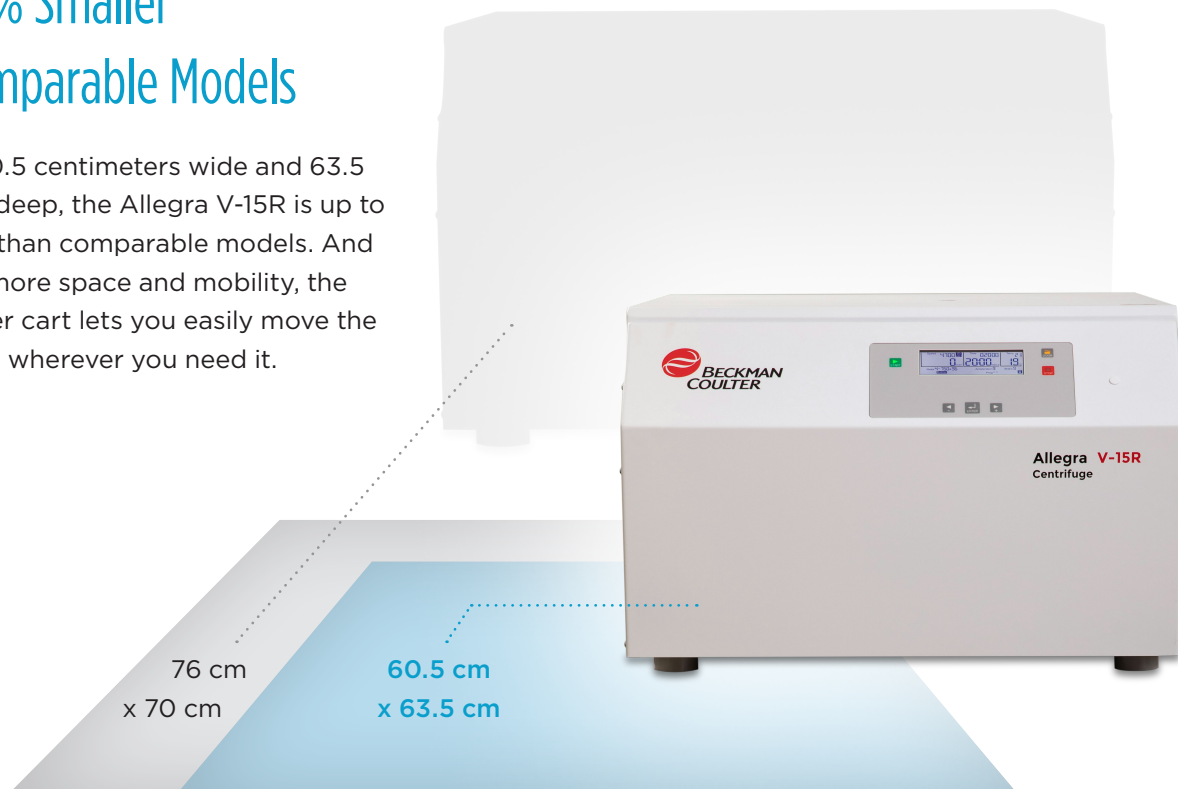


..... *Brilliance At Every Turn*

Free Up Precious Bench Space

Up to 28% Smaller Than Comparable Models

At a mere 60.5 centimeters wide and 63.5 centimeters deep, the Allegra V-15R is up to 28% smaller than comparable models. And if you need more space and mobility, the optional roller cart lets you easily move the centrifuge to wherever you need it.



Roller Cart

Not only does the roller cart give you mobility, it also gives you extra storage for rotors and accessories.

Spin More

With an extensive portfolio of swinging-bucket and fixed-angle rotors, and various adapters that can accommodate a wide range of tubes, bottles and plates, the Allegra V-15R brings more versatility and performance to your bench.

- 10 rotor configurations for a broad spectrum of applications and workflows, from High Throughput Samples all the way to cell and blood separation
- Capacity of up to 4 x 750 mL per run
- Gravitational force of up to 20,412 x g

Swinging-Bucket Rotors

4 x 750 mL Swinging-Bucket Rotor with several configurations.
Easily exchange buckets for microplate carriers or hexagonal carriers.



VS 4.750 **BIOC**

4 x 750 mL round bucket rotor; BioSafe with optional bucket covers
Max speed 240V / 120V:
4,651 x g (4,700 rpm) /
4,264 x g (4,500 rpm)



VS 4.750-96

Swing 4 x 4 microtiter plates or 4 x 1 deep well plates
Max speed 240V / 120V:
3,884 x g (4,700 rpm) /
3,561 x g (4,500 rpm)



VS 4.750-Hex

4 x 25 x 10 mL hexagonal bucket rotor
Max speed 240V / 120V:
4,478 x g (4,700 rpm) /
3,748 x g (4,300 rpm)



VS 2.5-96 **BIOC**

Swing 2 x 5 microtiter plates or 2 x 1 deep well plates; BioSafe when used with included cover
Max speed 240V / 120V:
5,495 x g (5,700 rpm) /
4,932 x g (5,400 rpm)

Fixed-Angle Rotors



VF 48.2 **BIOC**

48 x 1.5/2 mL microfuge tubes with biosafe lid
Max speed: 20,412 x g
(13,500 rpm)



VF 100.2

100 x 1.5/2 mL microfuge tubes
Max speed: 7,713 x g (outer row)
/ 7,145 x g (inner row)
(6,500 rpm)



VFC 24.15

24 x 15 mL conical tubes
Max speed: 11,431 x g
(9,000 rpm)



VFC 8.50 **BIOC**

8 x 50 mL conical tube rotor with BioSafe Lid
Max speed: 15,032 x g
(11,360 rpm)



VF 6.94 **BIOC**

6 x 94 mL tube rotor with Biosafe lid
Max speed: 11,872 x g
(10,000 rpm)



VF 6.250 **BIOC**

6 x 250 mL bottles with Biosafe lid
Max speed: 4,824 x g
(5,450 rpm)

Brilliance At Every Turn

Protecting Your Peak Performance

Service and Preventive Maintenance Agreements

Annual inspection and maintenance of your centrifuge and rotors by one of our factory-trained experts is imperative to ensure your instrument is operating efficiently and safely.

Qualification Services

Our service engineers are extensively trained in Installation Qualification and Operational Qualification (IQ/OQ) procedures and protocols to ensure your instrument is operating to manufacturer's specifications.

Protecting your peak performance is what we do.

Visit beckman.com/service to learn more.



| Beckman Coulter Service Plans | | | | | |
|---------------------------------------------|-----------------|--------------------------------|----------------------|------------|---------------|
| Plan Benefits | 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 | | ● | | ● | ● |
| Remote technical support | ● | ● | ● | ● | ● |
| Software & Engineering updates ⁴ | | ● | 20% Off | ● | ● |
| Annual basic operator training ⁵ | | ● | 20% Off | ● | ● |
| Operational qualification (OQ) | | AVAILABLE | | AVAILABLE | AVAILABLE |
| Relocation support | | ● | | | ● |

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.

Choose Beckman Coulter Life Sciences for Benchmark Expertise and Innovation

Providing 75 years of global leadership in centrifugation, Beckman Coulter Life Sciences designs, manufactures, sells, and services a complete line of centrifuge systems. By offering unique rotors and innovative bottles, tubes and accessories, Beckman Coulter Life Sciences delivers the highest level of expertise, innovation and quality assurance.

Contact your local Beckman Coulter Life Sciences sales representative at [beckman.com](https://www.beckman.com).



BioSafe and BioSafety are terms intended to describe the enhanced biocontainment features of our products.



BioCertified is a term used to describe our products which have been tested and validated to demonstrate containment of microbiological aerosols by an independent, third-party facility (TÜV NORD CERT GmbH, Große Bahnstraße 31, D-22525 Hamburg, Germany, or Public Health England, Porton Down, UK). Improper use or maintenance may affect seal integrity and, thus, containment.



© 2023 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. 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://www.beckman.com)

2023-GBL-EN-103768-v2