

AVANTI CENTRIFUGES

DEFINING A NEW ERA OF PERFORMANCE AND POSSIBILITIES.



Brilliance
at every turn.





WORLD-RENOWNED PERFORMANCE
TECHNOLOGICALLY ADVANCED
EASY TO USE HIGH ENDURANCE
CENTRIFUGES WITH MOBILE
CONNECTIVITY* BUILT RIGHT IN
IT'S THE WHOLE PACKAGE

*Mobile connectivity applies to Avanti JXN Series only.



1947 *Model E ultracentrifuge*

1976 *J6—one of the first UL- and CSA-listed centrifuges on the market*

1994 *Avanti J Series of high-performance centrifuges*

2010 *Centrifuge Center of Excellence established in Indianapolis, Indiana*

2011 *Launch of the Optima X Series centrifuges*

2014 *Avanti JXN Series and industry-leading MobileFuge app*

> And to think, it all started in an improvised chemistry lab.

When Arnold O. Beckman's dad let him convert a toolshed into a place to safely perform experiments, he probably had no idea where it would all lead. You know the cliché, necessity is the mother of invention. It's this same resourcefulness that drives Beckman Coulter to develop superior instruments for all ranges of applications and requirements today.

Because of our dedication to the life sciences industry, we created the Avanti Series of centrifuges—specifically for labs performing multiple and varied

applications and who demand versatility from a single instrument. Moreover, the Avanti Series delivers the speed, capacity, and performance you'd expect only from Beckman Coulter.

These high-performance, high-capacity instruments are designed to give you more throughput, faster turns, and greater control over your lab's workflow. And like all of our instruments, each Avanti centrifuge is a precision-made beauty worthy of the Beckman Coulter name.



AVANTI
IT'S BECKMAN BUILT.

Our history is rich with innovations that have benefited the greater good, having found thousands of applications in science, industry, and medicine. It is this heritage that proves our quality stands the test of time. Every piece of our equipment is engineered

Avanti JXN Series

Speed up your workflow.

Ergonomic design

- Low work surface height for easier installation and unloading of rotors
- Foot pedal for hands-free door operation
- Door swings up and out of the way for easy access to the rotor chamber
- Large digital readouts make operation simpler

Gain flexibility with MobileFuge

- Set and monitor run parameters and email diagnostic results
- Control and monitor the centrifuge from your Apple® iOS and Android™ device with MobileFuge
- Get alerts and diagnostics remotely

Intuitive interface

- User-friendly, large LCD screen
- View set and actual run conditions simultaneously
- Select rotor by name, not by arcane number code
- 11 accel/12 decel rates
- Delay start
- Run times to 99:59, Hold, and ω^{2t} integrator for applications that require exquisite reproducibility



with planned durability, so you can trust in reliable performance year after year.

Our equipment isn't the only thing that's exceptional. When you invest in Beckman Coulter solutions, you're backed by an incomparable worldwide

support organization. You can rely on easy Internet or phone access, field service teams, customizable service plans, original Beckman Coulter parts, and more. All to help you maintain your investment and ensure top performance from your lab.

Safe and sustainable

- Friction Reduction System (FRS) allows quick, quiet, and cool runs
- Uses less energy than traditional centrifuges
- BioSafe* options
- Dynamic Rotor Inertia Check
- High-torque, Switched Reluctance (SR)** drive cycle shortens cycle times

Advanced, easy-to-use data management

- Perfect for shared labs or to support GMP environments
- Up to 1,000 user-defined programs
- Set up and store numerous protocols
- Password protection for up to 50 unique user profiles

Versatile across rotors and applications

- Compatible with an extensive library of rotors
- Fixed-angle and swinging bucket options available
- Fully compatible with lightweight J-Lite rotors
- Dynamic Rotor Inertia Check (DRIC) and rotor recognition provide redundant safety checks to assure no rotor can be run above its rated speed
- Temperature control is optimized ensuring sample quality and avoiding unwanted artifacts
- SW and FA rotors capable of greater than 100,000 x g



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

**SR drives are manufactured by Beckman Coulter with technology licensed from Switched Reluctance Drives Limited.

Avanti J-26S XP

One centrifuge. So many possibilities.

The Avanti J-26S XP combines high performance, complete BioSafety* systems and sustainable advantages, making it an incredibly capable choice for any lab environment. This instrument is suitable for elutriation.

Along with high speed, it offers low heat output, imbalance tolerance and low noise. Versatility is delivered by a vast range of separations. Plus, ergonomic design features, such as hands-free door operation, make it easy to use.

For extra peace of mind, the Avanti J-26S XP has an impact-resistant lid, a full containment barrier and automatic rotor identification. Additionally, the resin used to manufacture the centrifuge comes from recycled material, which provides strong impact-, scratch-, and chemical-resistance for the life of the instrument.



J6-MI

It spins volumes.

With six-liter capacity and an innovative range of rotors, carriers, and adapters, this versatile centrifuge can handle multiple high-capacity applications. Ideal for bioprocessing and clinical and blood sample preparation, the J6-MI is the standard for quiet and convenient operation.

Its self-balancing rotor system corrects for imbalances during a run, without shutdown. Its microprocessor-controlled induction motor provides high torque, fast acceleration, and long motor life. And the J6-MI generates speeds and forces up to 6,000 rpm and $6,835 \times g$.



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

Avanti J-HC

Process more, in less time.

The Avanti J-HC is your complete solution for batch bioprocessing. Its capacity enables you to process up to nine liters spun at 5,000 rpm and 7,480 x g or 36 liters per hour when harvesting bacteria. And its selection of large-volume swinging bucket and fixed-angle rotors suits a wide range of processing applications.

Beyond versatility, the J-HC offers impressive levels of BioSafety*. J-Wide labware facilitates rapid harvesting of samples. Exclusive HarvestLine System Liners eliminate time-consuming scraping of samples and allow sealing and storing for future use. And Viewport strobing helps calibrate the instrument and achieves certification for regulatory agency compliance.



Avanti J-E

Compact, yet powerful.

Designed with a smaller footprint so it will fit right in your lab, the Avanti J-E is where you need it, when you need it. It also packs quite a punch.

The Avanti J-E's extensive rotor library maximizes capacity and optimizes g-force for efficient, fast separations and greater variety of applications. Its high-torque, Switched Reluctance (SR) drive technology shortens cycle times, which means less wait time for you. It offers an intuitive, user-friendly interface, as well as lower energy costs.

What's more, the Avanti J-E's low height and contoured front panel help ease the burden of rotor installation and retrieval, for even more convenient operation.





Avanti

As far as your ingenuity can take you.

Beckman Coulter helps create productive and efficient laboratories by providing high-quality centrifuges. Because we believe life-altering research should not be confined by the size of a lab, or the size of its throughput.

With the Avanti Series, we've set an all-new standard for enhanced performance and flexibility. Limits are now a thing of the past. In the footsteps of our founder, Beckman Coulter remains dedicated to your resourcefulness.

Service and support.

When you invest in Beckman Coulter instruments, you're backed by an incomparable support organization. In North America, our 24/7 customer support center directs a team of engineers with extensive product knowledge who have the power to solve your problems quickly and efficiently. And no matter where you are in the world, as a Beckman Coulter customer you have access to experienced, courteous service. Because when your issues are taken care of, you can return your attention to critical tasks.

www.beckmancoultercentrifugation.com



Beckman Coulter, the stylized logo, Avanti and HarvestLine are trademarks of Beckman Coulter, Inc. and are registered with the USPTO. 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 www.beckmancoulter.com

CENT-849SB03.15-A

©2015 Beckman Coulter, Inc.