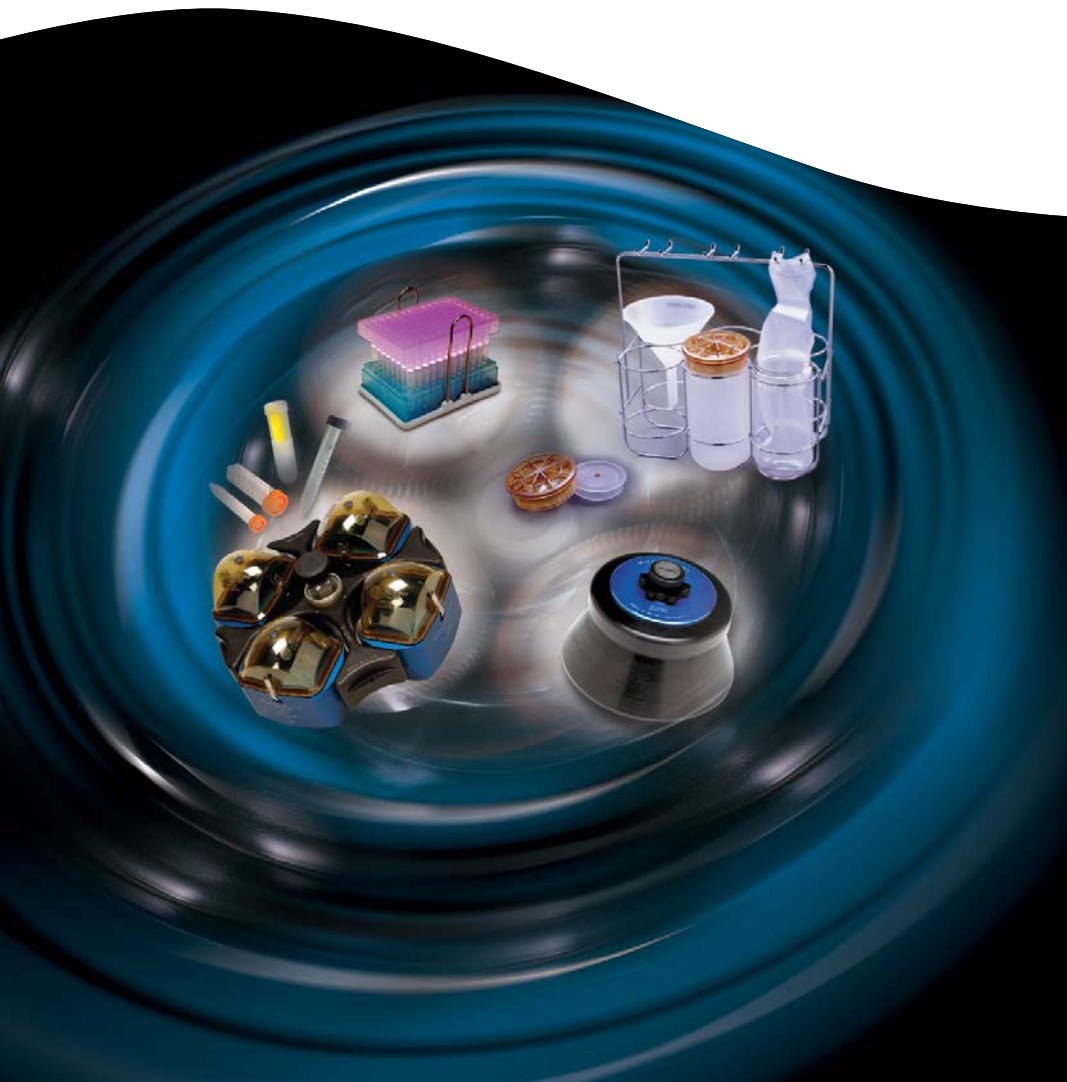




*A name you can trust for
dependability and value*

AVANTI® J SERIES
CENTRIFUGE SYSTEMS



A name you can trust for

THE AVANTI® ADVANTAGE

The Avanti series is designed, developed, produced and supported by Beckman Coulter, a company with more than 50 years of continuous excellence and leadership in centrifugation.

Building on this great history, the Avanti delivers dependability and value as the most productive and versatile high-performance centrifuge system with market-leading innovations that include:

- BioSafe systems
- High-torque drive design
- Lightweight rotors
- Ergonomic functionality
- Energy efficiency

Avanti —

A name you can trust for dependability and value.

APPLICATION VERSATILITY

The Avanti Series High-Performance Centrifuge system is an incredibly powerful and versatile tool for a wide range of separations. Whether you need high-throughput bioprocessing of up to 9 liters in a single run, fast microplate-based sample preparation, or high g-force purification of subcellular particles, the Avanti is the best solution. In fact, with its versatility, the Avanti can replace two or more older centrifuges.



SHORTER RUN TIMES — WITH HIGH-TORQUE SWITCHED RELUCTANCE (SR)[†] DRIVE

- Less waiting with fast accel and decel rates
- Imbalance tolerant drive allows eye-balancing of tubes
- Zero maintenance, brushless, SR drive technology

COMPLETE BIOSAFE[®] SYSTEMS FOR LABORATORY SAFETY

- Full line of BioSafe labware for any application
- Certified BioSafe rotors ensure containment of liquids and aerosols
- BioSafe centrifuges with HEPA filtration prevent infectious materials from being expelled into the lab environment



Biosafety

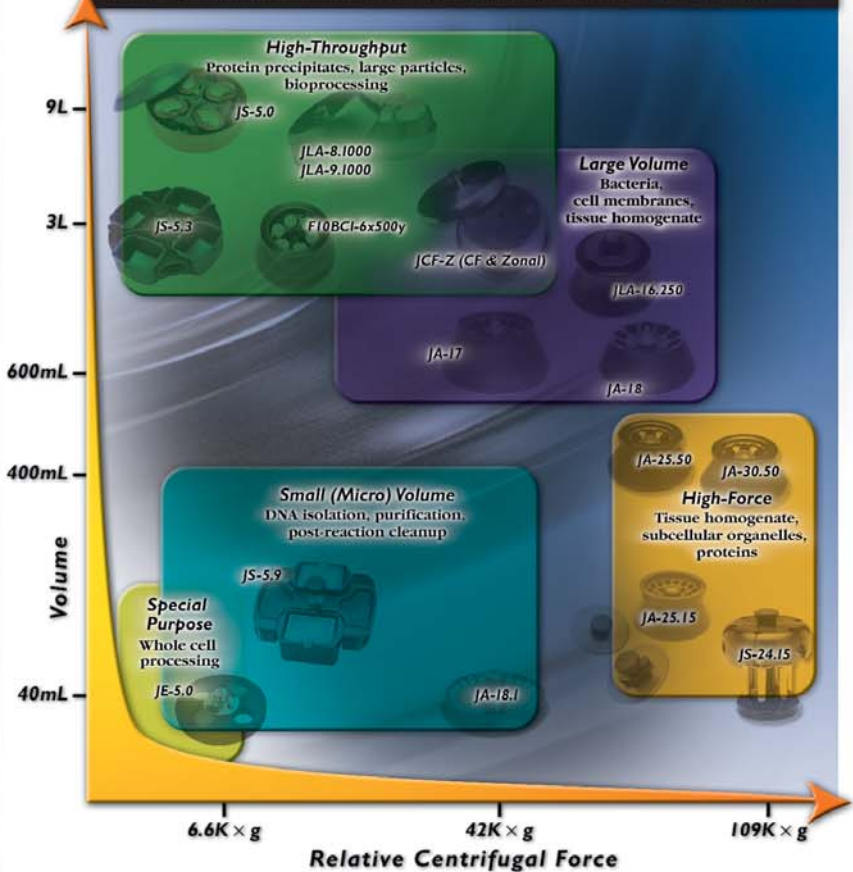
*Validation of microbiological containment was done at an independent, third-party testing facility, Health Protection Agency, Porton Down. Improper use or maintenance may affect seal integrity and, thus, containment.

LARGER VOLUMES & GREATER SEPARATION FORCES WITH "SMART FRS"[®] (FRICTION REDUCTION SYSTEM)

- Zero-maintenance partial vacuum system reduces demand on refrigeration system and drive for quieter, more economical operation
- FRS delivers higher rotor speeds, faster acceleration, and more precise temperature control
- FRS reduces laboratory air conditioning requirements, due to reduced heat output

dependability and value

Beckman Coulter's High-Performance Rotor Systems Break Through the Traditional Limits of High-Speed Separations!





AVANTI® J-301

Ultracentrifuge power from a High-Performance system



AVANTI® J-26 XP/XPi

The only High-Performance Centrifuge your lab may ever need



AVANTI® J-E

The economical performer that fits right in your lab



AVANTI® J-25/J25i

Higher productivity than any other conventional High-Speed Centrifuge

The Avanti® J Series is an important part of a broad spectrum of centrifuges — including High Performance, Ultracentrifuge and Benchtop — for a complete range of applications, including biosafety. As the global leader in centrifugation, Beckman Coulter designs and manufactures centrifuge systems offering rotors, adapters, bottles, microplate carriers, tubes, accessories and software that increase your productivity.

* BioSafe is a term intended to describe the enhanced biosafety features of Beckman Coulter products. Validation of microbiological containment was done at an independent, third-party testing facility, Health Protection Agency, Porton Down. Improper use or maintenance may affect seal integrity and, thus, containment.
† Online technical support available in North America
‡ SR Drives are manufactured by Beckman Coulter with technology licensed from Switched Reluctance Drives Limited

For additional information, please contact your local Beckman Coulter representative or visit our web site at:

www.beckmancoulter.com/centrifugefirst



Developing innovative solutions in Systems Biology.

Innovate SIMPLIFY Automate

Beckman Coulter, Inc. • 4300 N. Harbor Boulevard, Box 3100 • Fullerton, California 92834-3100
Sales & Service: 1-800-742-2345 • Telex: 678413 • Fax: 1-800-643-4366 • www.beckmancoulter.com

Worldwide Biomedical Research Division Offices:

Australia (61) 2 9844-6000 Canada (905) 819-1234 China (86) 10 6515 6028 Eastern Europe, Middle East, North Africa (41) 22 994 07 07
France 01 49 90 90 00 Germany 49 21 513335 Hong Kong (852) 2814 7431/2814 0481 Italy 02-953921 Japan 03-5404-8359
Mexico (55) 560-57770 Netherlands 0297-230630 Singapore (65) 6339 3633 South Africa, Sub-Saharan Africa (27) 11-805-2014/5 Spain 91 3836080
Sweden 08-564 85 900 Switzerland 0800 850 810 Taiwan (886) 2 2378 3456 Turkey 90 216 309 1900 U.K. 01494 441181 U.S.A. 1-800-742-2345