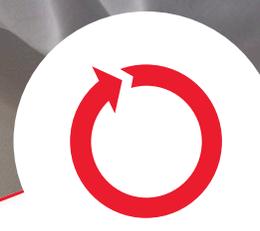


Avanti J-E

COMPACT, YET POWERFUL.



BRILLIANCE
at every turn.

 **BECKMAN
COULTER**
Life Sciences

HIGH-CAPACITY. SMALL FOOTPRINT.

The Avanti J-E is designed to deliver all the power you need, in even the tightest of laboratory spaces.

The Avanti J-E's extensive rotor library maximizes capacity and optimizes g-force for efficient, fast separations and a greater variety of applications. Its high-torque, switched reluctance (SR*) drive technology delivers faster acceleration and deceleration, 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.



Specifications

Maximum Speed/g-Force	21,000 rpm/53,300 x g
Maximum Capacity	4 Liters
Speed Control	±50 rpm of set speed
Set Temperature	-10° C to 40° C in 1° increments
Temperature Control	±2° C of set temperature after equilibration
Ambient Operating Range	15° C to 35° C
Accel/Decel Profiles	2/3
Heat Output	≤6900 BTU/hr (2.0 kW)
Drive Type/Cooling	SR* drive/Air-cooled
Refrigeration	Non-CFC, non-ozone depleting refrigerant
Sound Level	<64 dBa (0.91 m/3 ft from instrument at maximum speed)
Dimensions	63.5W x 80D x 91.4H cm (25W x 31.5D x 36H in) ¹
Weight	267.4 kg (589 lb)



¹Height measured with the door closed.

Note: Only values with tolerances or limits are guaranteed data. Values without tolerances are informative data, without guarantee.

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

ORDERING INFORMATION



Avanti Centrifuge Models

Model	Part No.	Description
Avanti J-E	369001	200/208/240V, 50 Hz
Avanti J-E	369003	230V, 50/60 Hz
Avanti J-E	369005	200V, 50/60 Hz
BioSafe** Avanti J-E	969352	200/208/240V, 50 Hz
BioSafe** Avanti J-E	A20698	230V, 50/60 Hz
BioSafe** Avanti J-E	A20699	200V, 50/60 Hz

Fixed-Angle Rotors

	Rotor Type	Part No.	Max Speed (rpm)	Max RCF, x g	k Factor	No. Tubes x Volume (mL)	Max Tube Size (mm)	Rotor Capacity
BIOC	JA-25.50	363055 Single-Locking Lid [§] 363058 Dual-Locking Lid	21,000	53,300	592	8 x 50	29 x 104	400 mL
BIOC	JA-20.1	342095	20,000	51,500/43,900	371/465	32 x 15	18 x 99	480 mL
	JA-21	334845	21,000	50,400	470	18 x 10	16 x 80	180 mL
BIOC	JA-20	334831	20,000	48,400	769	8 x 50	29 x 104	400 mL
BIOC	JA-17	369691	17,000	39,800	690	14 x 50	29 x 104	700 mL
BIOC	JLA-16.250	363934 Single-Locking Lid [§] 363930 Dual-Locking Lid	16,000	38,400	1,090	6 x 250	62 x 120	1.5 L
BIOC	JA-18	369679	16,000	37,800	716	10 x 100	38 x 102	1 L
BIOC	JA-14.50	B88554 Dual-Locking Lid B88555 Single-Locking Lid [§]	14,000	35,000	787	16 x 50	30 x 115	800 mL
BIOC	JA-14	339247	14,000	30,100	1,764	6 x 250	62 x 120	1.5 L
BIOC	JA-12	360993 Single-Locking Lid [§] 360992 Dual-Locking Lid	12,000	23,200	1,244	12 x 50 Conical	30 x 115	600 mL
BIOC	JLA-10.500	369681	10,000	18,600	2,840	6 x 500	69 x 160	3 L
	JA-10	369687	10,000	17,700	3,610	6 x 500	69 x 160	3 L
BIOC	JLA-9.1000	366754	6,300	8,230	5,192	4 x 1,000	95 x 191	4 L

Swinging Bucket Rotors

	Rotor Type	Part No.	Max Speed (rpm)	Max RCF, x g	k Factor	No. Tubes x Volume (mL)	Max Tube Size (mm)	Rotor Capacity
	JS-13.1	346963	13,000	26,500	1,841	6 x 50	29 x 104	300 mL
BIOC	JS-5.3 AllSpin	368690	5,300	6,870 (bottles) 6,130 (plates)	Various	4 x 500 24 Microplates	Various	2 L

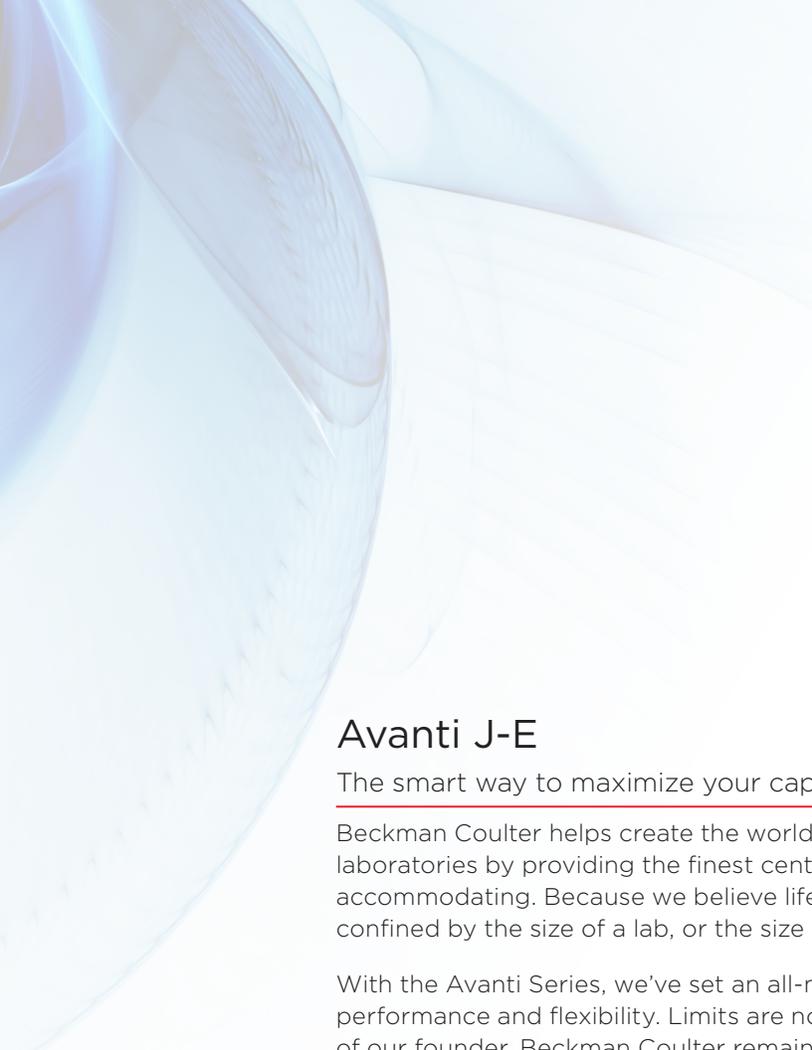
Dual-Locking Lids provide sample containment by enabling the rotor to remain sealed while being transported to a biocontainment hood.

[§]Single-locking lid versions of these rotors are not BioCertified.

****BioSafe and Biosafety** are terms intended to describe the enhanced biocontainment features of our products. Elutriation systems are available. Please contact your sales representative.

BIOC 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 (Health Protection Agency, Porton Down, UK or USAMRIID, Ft. Detrick, MD, USA). Improper use or maintenance may affect seal integrity and, thus, containment.

JS-5.3 is BIOC when using Aeroseal covers.



Avanti J-E

The smart way to maximize your capacity.

Beckman Coulter helps create the world's most productive and efficient laboratories by providing the finest centrifuges available. Finest, and most accommodating. 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.