



The Versatile Rotor Solution.

AllSpin JS-5.3 Rotor

FOR AVANTI® J-E AND J-26XP CENTRIFUGES



The highly versatile AllSpin rotor is ideally suited for Cellular Analysis, Genetic Analysis, and Protein Analysis providing you with a wide assortment of labware for your application requirements.

The 4-place design provides easy sample handling and ready access for time savings. Productivity is further enhanced with fast acceleration and deceleration from powerful Avanti SR* drive technology. And Beckman Coulter's Friction Reduction System delivers precise temperature control and energy saving results at up to 6,130 $\times g$ in both the Avanti J-26XP and Avanti J-E High Performance Centrifuges.¹

CELLULAR ANALYSIS: Up to 2L (4 x 500 mL) conical tube capability for efficient separation of cells, cell membranes, tissue homogenates, and blood cell components.


GENETIC ANALYSIS: Up to 24 (4 x 6 plates) 96-well plate capability for improved productivity in DNA isolation, purification, and post-reaction cleanup.

PROTEIN ANALYSIS: Up to 72 (4 x 18 place) 15 mL and 28 (4 x 7 place) 50 mL conical tube capability for fast preparation of protein/virus precipitates, and centrifugal filtration.

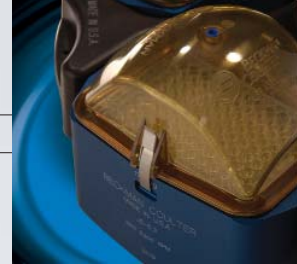
BIO SAFETY: Biocontainment from 5 mL tubes to 250 mL bottles and from PCR*** plates to deep well 96-well plates in a single rotor.



AllSpin JS-5.3 Rotor Parts and Accessories

Part No.	Description	Max. RPM
368690	Swinging-Bucket Rotor (Rotor Assembly)	5,300 rpm
969314	Swinging-Bucket Rotor Package (Rotor Assembly plus Microplate Carrier Kit - part no. 368914)	5,300 rpm
368417	Aeroseal™ Covers for JS-5.3 AllSpin rotor (set of 2)* 	5,300 rpm
368706	Bucket Upgrade Kit (set of 4 blue buckets)*	5,300 rpm

*rotors with black buckets require Bucket Upgrade Kit 368706



Pelleting Bacteria/Mammalian Cells

Part No.	Description	Max. RCF, tubes/bottles
392077	250 mL round-bottom bottle; 230 mL conical* bottle adapter	6,670 x g (5,300 rpm)
392078	500 mL conical bottle adapter (not compatible with Aeroseal Covers)	6,870 x g (5,300 rpm)
392079	250 mL conical bottle adapter; 60 x 172 mm	6,870 x g (5,300 rpm)

*requires a cushion (part no. 356983—set of 4) to be placed at bottom of adapter

Clearing Media — Cell Culture

Part No.	Description	Max. RCF, tubes/bottles
392073	15 mL round-bottom adapter (20 places), 17 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392074	50 mL round-bottom tube adapter (8 places), 29 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392075	15 mL conical tube adapter (18 places), 17 mm tubes	4,000 x g (4,200 rpm)
392076	50 mL conical tube adapter (7 places), 30 x 115 mm	4,000 x g (4,200 rpm)
392077	250 mL round-bottom bottle; 230 mL conical* bottle adapter	6,670 x g (5,300 rpm)
392078	500 mL conical bottle adapter (not compatible with Aeroseal Covers)	6,870 x g (5,300 rpm)
392079	250 mL conical bottle adapter; 60 x 172 mm	6,870 x g (5,300 rpm)

*requires a cushion (part no. 356983—set of 4) to be placed at bottom of adapter

Cell Membranes — Binding Studies

Part No.	Description	Max. RCF, tubes/bottles
392071	5 mL round-bottom tube adapter (33 places), 13 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392072	10 mL round-bottom adapter (24 places), 16 mm tubes	Use manufacturer's recommendation for max g-force
392073	15 mL round-bottom adapter (20 places), 17 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392074	50 mL round-bottom tube adapter (8 places), 29 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392075	15 mL conical tube adapter (18 places), 17 mm tubes	4,000 x g (4,200 rpm)
392076	50 mL conical tube adapter (7 places), 30 x 115 mm	4,000 x g (4,200 rpm)

Lysate/Tissue Homogenates — Animal Liver, Human Biopsy

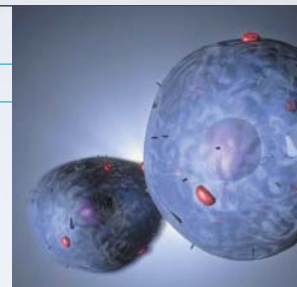
Part No.	Description	Max. RCF, tubes/bottles
392071	5 mL round-bottom tube adapter (33 places), 13 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392077	250 mL round-bottom bottle; 230 mL conical* bottle adapter	6,670 x g (5,300 rpm)
392078	500 mL conical bottle adapter (not compatible with Aeroseal Covers)	6,870 x g (5,300 rpm)
392079	250 mL conical bottle adapter; 60 x 172 mm	6,870 x g (5,300 rpm)

*requires a cushion (part no. 356983—set of 4) to be placed at bottom of adapter

Blood Cell Components — Blood Component Separation

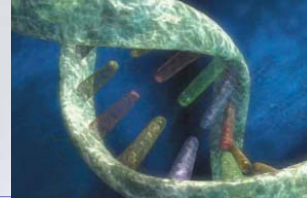
Ficoll-Paque[†] • AccuSpin System[†] • Histopaque[†] • I-Step 1.077/295[†] • I-Step monocytes • I-Step polymorphs

Part No.	Description	Max. RCF, tubes/bottles
392073	15 mL round-bottom adapter (20 places), 17 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392074	50 mL round-bottom tube adapter (8 places), 29 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392075	15 mL conical tube adapter (18 places), 17 mm tubes	4,000 x g (4,200 rpm)
392076	50 mL conical tube adapter (7 places), 30 x 115 mm	4,000 x g (4,200 rpm)



DNA Sample Preparation

Isolation & Purification — Microplate Format



Part No.	Description	Max. RCF, microplates
368914	Microplate Carrier Kit (required for all microplate applications)	
609844	24 polystyrene 96-well plates 300µL/well (nonsterile) (4 x 6 plates)	6,130 x g (5,300 rpm)
140504	4 Square-well polypropylene plates 2 mL/well (4 x 1 plates)	6,130 x g (5,300 rpm)
**	4 PCR plates and (4 x 1 plates)	6,130 x g (5,300 rpm)

Isolation & Purification — 8-well Strip

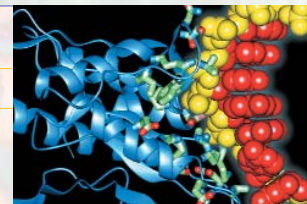
Part No.	Description	Max. RCF, microplates
368914	Microplate Carrier Kit (required for all microplate applications)	
267001	4 deep-well polystyrene 96-well plates 1 mL/well (nonsterile—nonstacking)	6,130 x g (5,300 rpm)
267004	4 deep-well polystyrene 96-well plates 1 mL/well (sterile—nonstacking)	6,130 x g (5,300 rpm)
267006	8 deep-well polypropylene 96-well plates 1 mL/well (nonsterile) (4 x 2 plates)	6,130 x g (5,300 rpm)
267007	8 deep-well polypropylene 96-well plates 1 mL/well (sterile) (4 x 2 plates)	6,130 x g (5,300 rpm)

Post-reaction cleanup — Filter Assembly

Part No.	Description	Max. RCF, microplates
368914	Microplate Carrier Kit (required for all microplate applications)	
**	4 PCR plates (4 x 1 plates)	6,130 x g (5,300 rpm)
**	4 DNA Kit 96-well filtration plates and (4 x 1 plates)	6,130 x g (5,300 rpm)

Virus PEG Precipitation

Part No.	Description	Max. RCF, tubes/bottles
392071	5 mL round-bottom tube adapter (33 places), 13 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392072	10 mL round-bottom adapter (24 places), 16 mm tubes	Use manufacturer's recommendation for max g-force



Proteins — Centrifugal Filtration

Part No.	Description	Max. RCF, tubes/bottles
392071	5 mL round-bottom tube adapter (33 places), 13 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392072	10 mL round-bottom adapter (24 places), 16 mm tubes	Use manufacturer's recommendation for max g-force
392073	15 mL round-bottom adapter (20 places), 17 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392074	50 mL round-bottom tube adapter (8 places), 29 mm tubes, variable lengths	Use manufacturer's recommendation for max g-force
392075	15 mL conical tube adapter (18 places), 17 mm tubes	4,000 x g (4,200 rpm)
392076	50 mL conical tube adapter (7 places), 30 x 115 mm	4,000 x g (4,200 rpm)



ORDERING INFORMATION:



	PART No.	DESCRIPTION
Avanti J-E Compact Centrifuge	369001	50/60 Hz, 208/240 V
	369003	50 Hz, 230 V
	369005	50/60 Hz, 200 V
BioSafe[†] Avanti J-E	969352	50/60 Hz, 208/240 V
	A20698	50 Hz, 230 V
	A20699	50/60 Hz, 200 V
Avanti J-26XP	393124	50/60 Hz, 200/208/240 V
	393125	50 Hz, 230 V
	393126	50 Hz, 220 V, 3-phase
Avanti J-26XPI	393127	50/60 Hz, 200/208/240 V
	393128	50 Hz, 230 V
	393129	50 Hz, 220 V, 3-phase
BioSafe[†] Avanti J-26XP	A25577	50/60 Hz, 200/208/240 V
	A25578	50 Hz, 230 V
	A25579	50 Hz, 220 V, 3-phase
BioSafe[†] Avanti J-26XPI	A25574	50/60 Hz, 200/208/240 V
	A25575	50 Hz, 230 V
	A25576	50 Hz, 220 V, 3-phase

¹The AllSpin JS-5.3 Rotor is also fully compatible with the J-20XP and J-20 Series High Performance Centrifuges.

References:

* 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.

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

** Plates available from outside manufacturers
All trademarks are the property of their respective owners

*** PCR is licensed under US patent numbers 4,683,202, 4,683,195, 4,965,188 and 5,075,216 or their foreign counterparts owned by Hoffman-La Roche and F. Hoffman-La Roche Ltd.

† Ficoll-Paque is a registered trademark of Pharmacia Fine Chemicals, Inc. AccuSpin and Histopaque are registered trademarks of Sigma Chemical Company. 1-Step is a registered trademark of Accurate Chemical and Scientific Corporation

†† 50 mL conical adapter (part no. 356966)
When used with part no. 392077, Max RCF is 6,670 x g

††† 15 mL conical adapter (part no. 356964)
When used with part no. 392077, Max RCF is 6,670 x g



Developing innovative solutions in Systems Biology.

Innovate Automate
SIMPLIFY

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