



# J-Lite JLA-12.500 Rotor

## The Next Revolution

Built for the Avanti JXN-26 and JXN-30 centrifuges, the JLA-12.500 Rotor improves RCF, increases bottle volume and boasts a biosafe design, making it the ideal choice for lysate clarification, protein precipitation and other biopharma workflows.

### KEY BENEFITS

#### 1 Best-in-Class Separation Performance

- Operates up to 26,890 x *g*
- 30° rotor angle provides a smaller pellet surface area, potentially reducing pellet disturbance during handling and decanting
- Aluminum construction provides sample temperature control

#### 2 BioC (BioCertified)\* **BIOC**

- Dual-locking lid assembly provides sample containment without the use of individual canisters
- Fluid containment annulus prohibits liquid from escaping into the centrifuge chamber

#### 3 Supports Sterile Workflows

- Aluminum construction allows the rotor to be autoclaved at 121°C for 60 minutes
- Capable of being autoclaved repeatedly and at a longer duration than composite rotors

#### 4 Provides Workflow Flexibility

- Supports numerous applications
- Operating volume (50-400 mL) supports sample density up to 1.2 g/mL
- Compatible with existing J-Lite adapters and numerous bottles and tubes



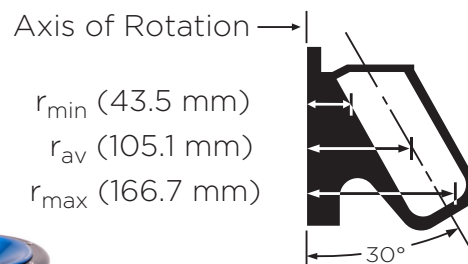
**BRILLIANCE**  
*at every turn.*

\*Validation of microbiological containment was done at an independent third-party testing facility (CAMR, Porton Down, UK, or USAMRIID, Ft. Detrick, MD, USA, or TÜV NORD CERT GmbH Große Bahnstraße 31 Hamburg, Germany).

## Specifications

Part Number	C55767
Max Speed (rpm)	12,000
Max Force (x g)	26,890
k-Factor	2,363
Fluid k-Factor*	1,277
No. of Tubes x Volume (mL)	6 x 500
Max Tube Size (mm)	69 x 160
Rotor Capacity (mL)	3,000
Instrument Compatibility	Avanti JXN-26, Avanti JXN-30

\*k-factor of maximum sample volume at maximum speed in PP bottle (361691).



## Tubes and Bottles

Nominal Dimensions and Volume	Part Number	Tube Description	Max Fill Volume (mL)	Required Adapter P/N	Max Speed/ RCF
69 x 160 mm 500 mL	361691	Polypropylene bottle with cap	400	None	12,000 rpm
					26,890 x g
69 x 160 mm 500 mL	361690	Polycarbonate bottle with cap	400	None	10,010 rpm
					18,711 x g
62 x 120 mm 250 mL	356011	Polypropylene wide mouth bottle with cap	250	362750	10,020 rpm
					18,263 x g
62 x 120 mm 250 mL	356013	Polycarbonate wide mouth bottle with cap	250	362750	10,020 rpm
					18,263 x g
29 x 104 mm 50 mL	357000	Polycarbonate round bottom tube with cap and plug	42	356996	10,040 rpm
					16,011 x g
29 x 104 mm 50 mL	357001	Polypropylene round bottom tube with cap and plug	42	356996	10,040 rpm
					16,011 x g
29 x 104 mm 50 mL	357002	Polycarbonate round bottom tube with cap	34	356996	10,040 rpm
					16,011 x g
29 x 104 mm 50 mL	357003	Polypropylene round bottom tube with cap	34	356996	10,040 rpm
					16,011 x g

Maximum speeds listed are guidelines only. These speeds have been achieved in testing at Beckman Coulter, but, because of variations in materials and chemicals that the product touches, no guarantee of performance or fit is expressed or implied.

When running less than max fill volume, speed must be reduced to 10,000 rpm maximum. When running at max speed, the minimum operating set temperature is 4°C. The useful service life of the 361691 Bottle Assembly has been tested to 100 hours (1 hour cycles) when running at the full rated speed/RCF at 25°C but will largely depend on handling and actual usage. Usage includes but is not limited to run duration, speed settings, temperature settings, cleaning and sterilization methods, and chemicals used. Bottle assembly components that show signs of visible wear, damage, stress, or deformation should be replaced. Do not use damaged components.

## Adapters to Support Tubes From Other Manufacturers

Type and Volume	Tube Description	Required Adapter*
Conical, 50 mL	Polypropylene conical tube with cap	356965
Conical, 15 mL	Polypropylene conical tube with cap	356960

\*Use manufacturers' recommendations for tube maximum fill volumes and speed/RCF limits

## Rotor Accessories

Description	Part Number
Rotor Lid Assembly	C72536
Lid O-ring Kit	C70015
Rotor Cleaning Kit	339558
Silicone Vacuum Grease (1 oz)	335148
Spinkote Lubricant (2 oz)	306812



Software version 2.3.232 and firmware version 4.17.0 or above are required for operation. Contact your local sales representative for information if operating an earlier version.

©2021 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](http://beckman.com)  
 21.04.1356.CENT