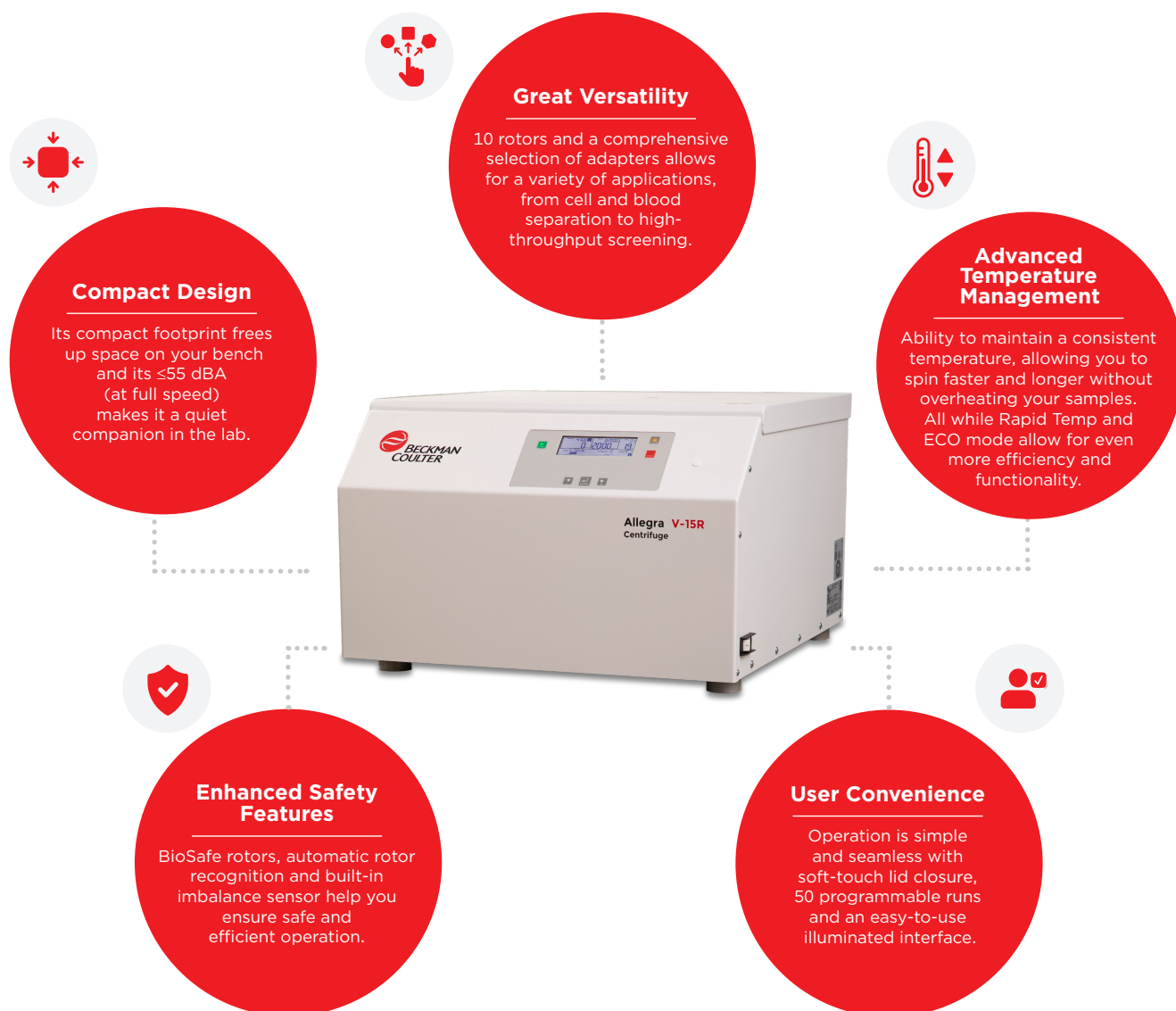




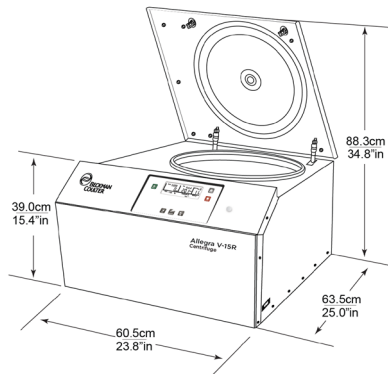
Allegra V-15R Benchtop Centrifuge

The Allegra V-15R benchtop centrifuge brings more versatility to your bench. This compact, refrigerated centrifuge is ideal for evolving labs looking to:

- Free up precious benchtop space
- Support multiple workflows with just one instrument



SPECIFICATIONS



Feature	Allegra V-15R Benchtop Centrifuge	
Max RCF	20,412 x g	
Fixed Angle	20,412 x g	
Swinging Bucket (120 V / 240 V)	4,264 x g / 4,651 x g	
Plates (120 V / 240 V)	4,932 x g / 5,495 x g	
Max Speed	13,500 rpm	
Capacity	4 x 750 mL / 60 x 15 mL / 28 x 50 mL / 4 x 4 plates	
Acceleration / Deceleration	10 / 10	
Number of Programs	50	
Set Temperature Range	-10 to 40° C	
Refrigerant	CFC-Free R452A	
Noise Level	≤55 dBA at max speed	
Dimensions	39 x 60.5 x 63.5 cm / 15.4 x 23.8 x 25 in (H x W x D)	
Ordering Information	General Use ¹	IVD ²
220 - 240V 50 Hz	C63126	C63127
200V 50/60Hz, 208V 60Hz	C63128	C63129
120V 60Hz	C63124	C63125
220 - 240V 50Hz with CCC Plug	C63161	-

Mobile Cart for Allegra V-15R

Cart Dimensions	58.1 x 60.5 x 63.5 cm / 22.9 x 23.8 x 25 in (H x W x D)
Cart with Centrifuge Dimensions	97.1 x 60.5 x 63.5 cm / 38.2 x 23.8 x 25 in (H x W x D)
Ordering Information	Description
Order No.	C63177



¹ General Use. Not intended or validated for use in the diagnosis of disease or other conditions.

² In Vitro Diagnostics Use.

APPLICATION PACKAGES

Cell Culture Package

Includes: Instrument, VS 4.750 swinging-bucket rotor, 15 mL and 50 mL conical tube adapters (qty of 4 of each)

Wash, clarify, concentrate, harvest and separate cells and tissues from cultures. From mammalian to bacterial cells and everything in between, our Cell Culture Application Package is designed to deliver the results you need. Run 15 mL or 50 mL tubes. Ideal for processing cell cultures in large-, medium- and mini-scales. Increase laboratory safety with available BioCertified bucket covers (not included).



	General Purpose ¹	IVD ²
120V, 60 Hz	C63438	C63439
220-240V, 50 Hz	C63436	C63437
200V, 50/60 Hz, 208V, 60 Hz	C63440	C63441

Blood Sample Preparation Package

Includes: Instrument, VS 4.750 swinging-bucket rotor, BioSafe bucket covers, and 12.8 mm and 13.8 mm and 17.5 mm adapters (qty of 4 each)

Processing up to 148, 12.8 mm diameter blood tubes, 92, 13.8 mm diameter blood tubes or 76, 17.5 mm diameter blood tubes per run. Our Blood Sample Preparation packages deliver the speed and flexibility your busy laboratory needs.



	General Purpose ¹	IVD ²
120V, 60 Hz	C63420	C63421
220-240V, 50 Hz	C63418	C63419
200V, 50/60 Hz, 208V, 60 Hz	C63422	C63423

¹ General Use. Not intended or validated for use in the diagnosis of disease or other conditions.
² In Vitro Diagnostics Use.

High Capacity Blood Sample Package

Includes: Instrument VS 4.750-Hex swinging-bucket rotor and adapters 13.5 mm, and 17.5 mm round bottom blood tubes (qty of 4 each)

Maximize throughput and capacity—processing up to 144, 13.5 mm diameter blood tubes or 100, 17.5 mm diameter blood tubes per run.

	General Purpose ¹	IVD ²
120V, 60 Hz	C63426	C63427
220-240V, 50 Hz	C63424	C63425
200V, 50/60 Hz, 208V, 60 Hz	C63428	C63429



Swinging-Bucket Package

Includes: Instrument and VS 4.750 swinging-bucket rotor

Achieve great versatility all in one rotor with the ability to swing 4x750 mL bottles or other common tubes and bottles with our selection of adapters.

	General Purpose ¹	IVD ²
120V, 60 Hz	C63386	C63387
220-240V, 50 Hz	C63384	C63385
200V, 50/60 Hz, 208V, 60 Hz	C63388	C63389



BioSafe Swinging-Bucket Package

Includes: Instrument, VS 4.750 swinging-bucket rotor and BioSafe bucket covers (qty 4)

Minimize contamination from broken or leaking tubes with BioSafe bucket covers, contributing to laboratory safety. Each cover seals with an O-ring and is autoclavable up to 121° C for 20 minutes.

	General Purpose ¹	IVD ²
120V, 60 Hz	C63444	C63445
220-240V, 50 Hz	C63442	C63443
200V, 50/60 Hz, 208V, 60 Hz	C63446	C63447



¹ General Use. Not intended or validated for use in the diagnosis of disease or other conditions.

² In Vitro Diagnostics Use.

Microplate Swinging-Bucket Package

Includes: Instrument and VS 4.750-96 swinging-bucket rotor

Ideal for high-throughput microplate applications, enabling speedy preparation of your samples for downstream applications including UV-VIS absorbance, fluorescence, luminescence, PCR and real-time PCR.

	General Purpose ¹	IVD ²
120V, 60 Hz	C63450	C63375
220-240V, 50 Hz	C63448	C63449
200V, 50/60 Hz, 208V, 60 Hz	C63376	C63377



Deep Well Plate Swinging-Bucket Package

Includes: Instrument and VS 2.5-96 swinging-bucket rotor with BioSafe bucket covers

Ideal for high-throughput and high g-force microplate applications, as well as recovering protein precipitates, soluble proteins and elements in supernatant released from cells or tissues cultured in plates. Our Deep Well Package supports laboratory safety with the option to use BioSafe bucket covers. Each cover seals with an O-ring and is autoclavable up to 121° C for 20 minutes.

	General Purpose ¹	IVD ²
120V, 60 Hz	C63380	C63381
220-240V, 50 Hz	C63378	C63379
200V, 50/60 Hz, 208V, 60 Hz	C63382	C63383



Fixed-Angle Rotor Package

Includes: Instrument and VFC 8.50 rotor

High-speed rotor that can spin 8 x 50 mL conical tubes for a variety of applications. Adapters allow for the spinning of 50 mL round-bottom tubes, 25-30 mL skirted tubes and 15 mL conical tubes.

BioSafe lid minimize contamination from broken or leaking tubes and helps ensure laboratory safety.

	General Purpose ¹	IVD ²
120V, 60 Hz	C63432	C63433
220 - 240V, 50 Hz	C63430	C63431
200V, 50/60 Hz, 208V, 60 Hz	C63434	C63435



¹ General Use. Not intended or validated for use in the diagnosis of disease or other conditions.
² In Vitro Diagnostics Use.

ROTOR AND ADAPTER SPECIFICATIONS AND ORDER INFORMATION

VS 4.750 Swinging-Bucket Rotor

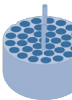
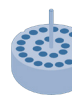


Item No.	C63142
Rotor Capacity	4 x 750 mL
Rotor Angle	90°
Max. RPM (120V / 240V)	4,500 / 4,700
Max. RCF (120V / 240V)	4,264 / 4,651
Permissible Net Weight	4 x 1,000 g
Max. Speed for 4° C Operation	4,700

Tube Specifications							
Tube Type	Flat	Flat	Conical	Flat	Conical	Conical	Round
Tube Volume	750 mL	500 mL	500 mL	250 mL	200 /230/ 250 mL	200 / 250 mL	100 mL
Tubes per Rotor	1 x 4	1 x 4	1 x 4	1 x 4	1 x 4	1 x 4	2 x 4
Max. Tube Diameter (mm)	100	85	95.5	62	61 / 61 / 62	61 / 61	44
Max. Tube Height (mm)	127	135	147	120	145 / 130 / 139.5	118 / 137	100
Required Adapter Specifications							
Item No.	Not Required	C63170	C63171	C63401	C63399	C63169	C63398
Max. RPM	4,700	4,700	4,700	4,700	4,700	4,700	4,700
Max. RCF	4,651	4,577	4,651	4,577	4,651	4,651	4,577
Quantity per Pack	—	2	2	2	2	2	2

*Not compatible with BioSafe Lids (C63153)

Tube Specifications								
Tube Type	Round	Round	Conical	Conical	Skirted	Skirted	Conical	Conical
Tube Volume	50 mL	50 mL	50 mL	50 mL	50 mL	25 - 30 mL	15 mL	15 mL
Tubes per Rotor	4 x 4	7 x 4	5 x 4	7 x 4	5 x 4	7 x 4	12 x 4	15 x 4
Max. Tube Diameter (mm)	35	29	29	29	29	25	16.5	16.5
Max. Tube Height (mm)	100	107	117	117	117	90 - 107	120	120
Required Adapter Specifications								
Item No.	C63397	C63166	C63167	C63168*	C63373	C63372	C63164	C63165*
Max. RPM	4,700	4,700	4,700	4,700	4,700	4,700	4,700	4,700
Max. RCF	4,577	4,577	4,602	4,602	4,503	4,330	4,267	4,267
Quantity per Pack	2	2	2	2	2	2	2	2

Tube Specifications					
Tube Type	Blood Collection	Blood Collection	Blood Collection	Blood Collection	Micro Tube
Tube Volume	10 mL	7 mL	5 mL	10 mL	2 mL
Tubes per Rotor	20 x 4	23 x 4	37 x 4	19 x 4	26 x 4
Max. Tube Diameter (mm)	16	13	12	16	11
Max. Tube Height (mm)	118	107	108	108 - 131	42
Required Adapter Specifications					
Item No.	C63174	C63173	C63172	C63188	C63163
Max. RPM	4,700	4,700	4,700	4,700	4,700
Max. RCF	4,528	4,503	4,453	4,528	4,478
Quantity per Pack	2	2	2	2	2

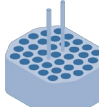
*Not compatible with BioSafe Lids (C63153)

Accessory Specifications			
Item No.	C63152	C63153	961930
Accessory Type	Bucket	BioSafe Lid	O-ring
Max. RPM (120V / 240V)	4,500 / 4,700	-	-
Max. RCF (120V / 240V)	4,264 / 4,651	-	-
Quantity per Pack	2	2	2

VS 4.750-Hex Swinging-Bucket Rotor



Item No.	C63143
Rotor Capacity	4 x 25 x 10 mL
Rotor Angle	90°
Max. RPM and flip 120V / 240V	4,300 / 4,700
Max. RCF (120V / 240V)	3,748 / 4,478
Permissible Net Weight	4 x 900 g
Max. Speed for 4° C Operation	4,700

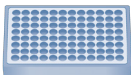
Tube Specifications				
Tube Type	Blood Collection	Beckman Coulter AutoMate 2500	Blood Collection	Beckman Coulter AutoMate 2500
Tube Volume	10 mL	10 mL	7 mL	5 - 7 mL
Tubes per Rotor	25 x 4	16 x 4	36 x 4	25 x 4
Max. Tube Diameter (mm)	16	16	7	13
Max. Tube Height (mm)	91 - 131	57	13.5	57
Required Adapter Specifications				
Item No.	B30603	B32767	C63176	B32766
Max. RPM	4,700	4,700	4,700	4,700
Max. RCF	4,379	4,354	4,379	4,404
Quantity per Pack	2	2	2	2

Accessory Specifications	
Item No.	C63154
Accessory Type	Bucket
Max. RPM (120V / 240V)	4,300 / 4,700
Max. RCF (120V / 240V)	3,748 / 4478
Quantity per Pack	2

VS 4.750-96 Swinging-Bucket Rotor



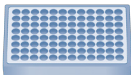
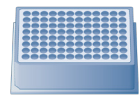
Item No.	C63144
Rotor Capacity	4 x 4 x 96 mL
Rotor Angle	90°
Max. RPM (120V / 240V)	4,500 / 4,700
Max. RCF (120V / 240V)	3,561 / 3,884
Permissible Net Weight	4 x 500 g
Max. Speed for 4° C Operation	4,700 RPM

Tube Specifications		
Tube Type	Microtiter Plate	Deep Well Plates
Tube Volume	MTP	DWP
Plates per Rotor	4 x 4	1 x 4
Max. Plate Dimensions (mm)	86 x 128	86 x 128
Accessory Specifications		
Item No.	C63156	C63157
Accessory Type	Bucket Assembly (includes carrier)	Carrier
Max. RPM (120V / 240V)	4,500 / 4,700	—
Max. RCF (120V / 240V)	3,561 / 3,884	—
Quantity per Pack	2	2

VS 2.5.96 Swinging-Bucket Rotor



Item No.	C63145
Rotor Capacity	2 x 5 x 96 mL
Rotor Angle	90°
Max. RPM (120V / 240V)	5,400 / 5,700
Max. RCF (120V / 240V)	4,932 / 5,495
Permissible Net Weight	4 x 520 g
Max. Speed for 4° C Operation	5,600 RPM

Tube Specifications				
Tube Type	Microtiter Plate		Deep Well Plates	
Tube Volume	MTP		DWP	
Plates per Rotor	5 x 2		1 x 2	
Max. Plate Dimensions	86 x 128		86 x 128	

Accessory Specifications				
Item No.	C63160	C63158	C63159	C63181
Accessory Type	Bucket Assembly (includes carriage)	Carriage	Biosafe Lids	Gasket
Max. RPM (120V / 240V)	5,400 / 5,700	--	--	--
Max. RCF (120V / 240V)	4,932 / 5,495	--	--	--
Quantity per Pack	2	2	2	2

VF 48.2 Fixed-Angle Rotor **BIOC**



Item No.	C63136
Rotor Capacity	48 x 2 mL
Rotor Angle	25° (Outer Row) / 53° (Inner Row)
Max. RPM	13,500
Max. RCF	20,412
Max. Speed for 4° C Operation	13,000

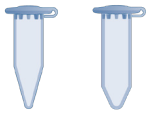
Tube Specifications				
Tube Type	Microtube	Microtube	Microtube	Microtube
Tube Volume	1.5 - 2 mL	0.5 mL	0.4 mL	0.2 mL
Tubes per Rotor	1 x 48	1 x 48	1 x 48	1 x 48
Max. Tube Diameter (mm)	10.7	7.9	5.8	5.85
Max. Tube Height (mm)	42	28	47	20
Required Adapter Specifications				
Item No.	Not Required	C63416	C63415	C63414
Max. RPM	13,500	13,500	13,500	13,500
Max. RCF	20,412	20,004	20,004	19,596
Quantity per Pack	—	12	12	12

Accessory Specifications			
Item No.	C63146	A57892	C63179
Accessory Type	BioSafe Lid	Large O-ring	Small O-ring

VF 100.2 Fixed-Angle Rotor



Item No.	C63137
Rotor Capacity	100 x 2 mL
Rotor Angle	45°
Max. RPM	6,500
Max. RCF	7,713 (Outer Row) 7,145 (Inner Row)
Max. Speed for 4° C Operation	6,500

Tube Specifications				
Tube Type	Microtube	Microtube	Microtube	Microtube
Tube Volume	1.5 - 2 mL	0.5 mL	0.4 mL	0.2 mL
Tubes per Rotor	1 x 100	1 x 100	1 x 100	1 x 100
Max. Tube Diameter (mm)	10.7	7.9	5.8	5.85
Max. Tube Height (mm)	42	28	47	20
Required Adapter Specifications				
Item No.	Not Required	C63416	C63415	C63414
Max. RPM	6,500	6,500	6,500	6,500
Max. RCF	7,713	7,619	7,619	7,524
Quantity per Pack	—	12	12	12

Accessory Specifications		
Item No.	C63147	A57892
Accessory Type	Lid	Large O-ring

VFC 24.15 Fixed-Angle Rotor



Item No.	C63137
Rotor Capacity	24 x 15 mL
Rotor Angle	30° (Outer Row) 40° (Inner Row)
Max. RPM	9,000
Max. RCF	11,431
Max. Speed for 4° C Operation	9,000

Tube Specifications	
Tube Type	Conical
Tube Volume	15 mL
Tubes per Rotor	1 x 24
Max. Tube Diameter (mm)	16.5
Max. Tube Height (mm)	120
Required Adapter Specifications	
Item No.	Not Required
Max. RPM	9,000
Max. RCF	11,431
Quantity per Pack	—

Accessory Specifications			
Item No.	C63148	A57892	368300
Accessory Type	Lid	Large O-ring	Small O-ring

VFC 8.50 Fixed-Angle Rotor



Item No.	C63139
Rotor Capacity	8 x 50 mL
Rotor Angle	25°
Max. RPM	11,360
Max. RCF	15,032
Max. Speed for 4° C Operation	11,360

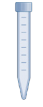
Tube Specifications				
Tube Type	Conical	Round	Skirted	Conical
Tube Volume	50 mL	50 mL	25 - 30 mL	15 mL
Tubes per Rotor	1 x 8	1 x 8	1 x 8	1 x 8
Max. Tube Diameter (mm)	29.5	29	25	16.5
Max. Tube Height (mm)	117	104	90 - 107	120
Required Adapter Specifications				
Item No.	Not Required	C63371	C63369	C63417
Max. RPM	11,360	11,360	11,360	11,360
Max. RCF	15,032	13,586	13,586	14,887
Quantity per Pack	—	2	2	2

Accessory Specifications			
Item No.	C63149	C69755	C63179
Accessory Type	BioSafe Lid	Large O-ring	Small O-ring

VF 6.94 Fixed-Angle Rotor



Item No.	C63140
Rotor Capacity	6 x 94 mL
Rotor Angle	25°
Max. RPM	10,000
Max. RCF	11,872
Max. Speed for 4° C Operation	10,000

Tube Specifications									
Tube Type	Round	Conical	Round	Round	Round	Round	Conical	Round	Microtube
Tube Volume	94 mL	50 mL	50 mL	30 mL	30 mL	15 mL	15 mL	10 mL	1.5 mL
Tubes per Rotor	1 x 6	1 x 6	1 x 6	1 x 6	1 x 6	1 x 6	1 x 6	2 x 6	6 x 6
Max. Tube Diameter (mm)	38	29.5	29	25.3	23.6	17.5	16.5	16.1	11
Max. Tube Height (mm)	102	117	104	89	106	106	120	81.1	40
Required Adapter Specifications									
Item No.	Not Required	C63412	392830	C63410	C63411	C63408	C63409	C63407	344497 in 392830
Max. RPM	10,000	10,000	10,000	10,000	10,000	10,000	10,000	10,000	10,000
Max. RCF	11,872	11,872	11,536	9,968	11,088	10,976	11,760	8,624	8,669
Quantity per Pack	—	2	6	2	2	2	2	2	6

Accessory Specifications			
Item No.	C63150	C69755	C63179
Accessory Type	BioSafe Lid	Large O-ring	Small O-ring

VF 6.250 Fixed-Angle Rotor **BIOC**



Item No.	C63141
Rotor Capacity	6 x 250 mL
Rotor Angle	30°
Max. RPM	5,450
Max. RCF	4,824
Max. Speed for 4° C Operation	5,450

Tube Specifications					
Tube Type	Flat	Round	Conical	Conical	Microtube
Tube Volume	250 mL	50 mL	50 mL	15 mL	1.5 mL
Tubes per Rotor	1 x 6	1 x 6	1 x 6	4 x 6	9 x 6
Max. Tube Diameter (mm)	62	29	29	16.5	10.7
Max. Tube Height (mm)	120	104	117	120	42
Adapter Specifications					
Item No.	Not Required	C63403	C63405	C63404	C63406
Max. RPM	5,450	5,450	5,450	5,450	5,450
Max. RCF	4,824	4,558	4,691	4,790	4,258
Quantity per Pack	—	2	2	2	2

Accessory Specifications			
Item No.	C63151	C69757	C63179
Accessory Type	BioSafe Lid	Large O-ring	Small O-ring

LABWARE SPECIFICATION TABLE

Bottles

Item No.	Type	Package Quantity	Closure	Material	Nominal	Diameter	Length
C63392	Flat-Bottom, Wide-Mouth	2	Screw-cap	Polypropylene	750 mL	100 mm	127 mm
C63162	Flat-Bottom, Wide-Mouth	6	Screw-cap	Polypropylene	500 mL	85 mm	135 mm
356011	Flat-Bottom, Wide-Mouth	6	Cap Assembly	Polypropylene	250 mL	62 mm	120 mm
356013	Flat-Bottom, Wide-Mouth	6	Cap Assembly	Polycarbonate	250 mL	62 mm	120 mm
358275	Flat-Bottom, Wide-Mouth	6	Open-top	Polycarbonate	250 mL	62 mm	120 mm
358326	Flat-Bottom, Wide-Mouth	6	Open-top	Polypropylene	250 mL	62 mm	120 mm

Tubes

Item No.	Type	Package Quantity	Closure	Material	Nominal	Diameter	Length
355643	Round-Bottom Tube	25	Open-top	Polypropylene	94 mL	38 mm	102 mm
363081	Round-Bottom Tube	6	Cap Assembly	Polycarbonate	85 mL	38 mm	104 mm
363082	Round-Bottom Tube	6	Screw-cap	Polypropylene	80 mL	38 mm	104 mm
361693	Round-Bottom Tube	24	Cap Assembly	Polycarbonate	50 mL	29 mm	104 mm
361694	Round-Bottom Tube	24	Cap Assembly	Polypropylene	50 mL	29 mm	104 mm
357002	Round-Bottom Tube	25	Screw-Cap	Polycarbonate	50 mL	29 mm	104 mm
357003	Round-Bottom Tube	25	Screw-Cap	Polypropylene	50 mL	29 mm	104 mm
363647	Round-Bottom Tube	25	Open-top	Polycarbonate	50 mL	29 mm	104 mm
357007	Round-Bottom Tube	25	Open-top	Polypropylene	50 mL	29 mm	104 mm
363664	Round-Bottom Tube	25	Snap-On Cap	Polycarbonate	50 mL	29 mm	104 mm
357000	Round-Bottom	6	Cap Assembly	Polycarbonate	50 mL	29 mm	104 mm
357001	Round-Bottom	6	Cap Assembly	Polycarbonate	50 mL	29 mm	104 mm
357005	Round-Bottom Tube	25	Snap-On Cap	Polypropylene	50 mL	29 mm	104 mm
363075	Round-Bottom, Graduated Tube	8	Open-top	Polycarbonate	50 mL	29 mm	104 mm
340196	Microtube	500	Snap-On Cap	Polyethylene	1.8 mL	11 mm	45 mm
357448	Microtube	500	Snap-On Cap	Polypropylene	1.5 mL	11 mm	38 mm
343169	Microtube	500	Separate Cap	Polypropylene	1.5 mL	11 mm	40 mm



BioSafe and BioSafety are terms intended to describe the enhanced biocontainment features of our products. 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 (TÜV NORD CERT GmbH, Große Bahnstraße 31, D-22525 Hamburg, Germany, or Public Health England, Porton Down, UK). Improper use or maintenance may affect seal integrity and, thus, containment.

© 2022 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](https://www.beckman.com)
22.01.3762.CENT