

Sample
preparation
made easy.



revvity

CF-350 Centrifuge

Universal sample preparation made easy.

The CF-350 is a universal laboratory centrifuge designed for use in medical, industrial, and scientific laboratories for separating substances with different relative densities by centrifugal force. The maximum rotational speed of 15,000 rpm gives the centrifugal force of 21,630 rcf.

A wide usage of programming options and electronic operation control allow for trouble-free use. With a special drive, quiet, maintenance-free operation with no carbon dust pollution is guaranteed.

The centrifuge has user-friendly options that make operation and standard settings easier. Built-in error-detecting functions keep the user from entering incorrect values and monitor the whole operation.

The centrifuge has the possibility to save programs. In the program memory, you can save up to 100 different programs. The centrifuge always keeps the last run program in its memory for an unlimited amount of time, allowing the program to be restarted at any time, even if the centrifuge was turned off during operation. All important operation parameters can be seen at a glance.

The settings are entered with knobs and keys on the control panel. The centrifuge's interior is also easy to clean.

User manual, rotor key, and power cord included with instrument. All rotors, buckets, and inserts are sold separately.

Part number:
CF-350-120/220/220UK



Specifications

Power rating	CF-350-120 - 120V, 50/60 Hz
	CF-350-220 - 220V, 50/60 Hz
	CF-350-220UK - 220V, 50/60 Hz
Max. power consumption	500 Watts
Max. rotation speed	200 to 15,000 rpm, in steps of 10 rpm
Maximum G force	21,630 rcf
Maximum capacity	6 x 50 mL (Angle Rotor) 4 x 200 mL (Swing out Rotor)
Rotors available	13
Time	10 s to 99 min and continuous, step 10 s
Programs	100
Acceleration levels	10
Deceleration levels	10 (Level 0 without braking)
Display	LED
Noise level (at max. speed)	≤ 66 dBA
Fast spin	Yes
Lid locking mechanism	Motorized
Lid Lock	2 Latches
Motor type	Induction brushless motor
Permitted ambient conditions	2 to 35 °C, Humidity: 75 %, non-condensing
Dimensions	W: 15.9" (40.5 cm), D: 19.7" (50.0 cm), H: 12.8" (32.5 cm) (H: lid open 23.9" (60.6 cm))
Weight	63.9 lbs. (29 kg)
Warranty	One year
Certification	CE

Rotors



**24 x 2 mL
Microcentrifuge Tube
Rotor with Lid**

Part number: 228442

- Rotor type: RA 24/2 angle rotor with lid
- Maximum capacity: 24 x 1.5/2.0 mL microcentrifuge tubes - optional 0.2 mL or 0.5 mL adapters for PCR tubes (283451 or 283349)
- Maximum tube dimensions: 11 x 40 mm
- Maximum speed: 15,000 rpm
- Maximum centrifugal force: 21,630 rcf
- Radius: 86 mm
- Angle: 45°
- Minimum acceleration time: 22 s
- Minimum braking time: 23 s
- Rotor and lid autoclavable: Yes



30 x 12 mL Tube Rotor
Part number: 454655

- Rotor type: RA 30/12 angle rotor
- Maximum capacity: 30 x 12 mL blood collection tubes - optional A5 adapters for 75 to 90 mm tubes
- Maximum tube dimensions: 17 x 120 mm
- Maximum speed: 4,400 rpm
- Maximum centrifugal force: 3,200 rcf
- Radius: 148 mm
- Angle: 45°
- Minimum acceleration time: 21 s
- Minimum braking time: 21 s
- Rotor and shields autoclavable: yes



**24 x 2 mL
Microcentrifuge Tube
Rotor with Aerosol Lid**

Part number: 375944

- Rotor type: RA 24/2 angle rotor with aerosol tight lid
- Maximum capacity: 24 x 1.5/2.0 mL microcentrifuge tubes - optional 0.2 mL or 0.5 mL adapters for PCR tubes (283451 or 283349)
- Maximum tube dimensions: 11 x 40 mm
- Maximum speed: 15,000 rpm
- Maximum centrifugal force: 21,630 rcf
- Radius: 86 mm
- Angle: 45°
- Minimum acceleration time: 21 s
- Minimum braking time: 21 s
- Rotor and lid autoclavable: yes



6 x 50 mL Tube Rotor
Part number: 448914

- Rotor type: RA 6/50 angle rotor with buckets
- Maximum capacity: 6 x 50 mL conical centrifuge tubes or 6 x 15 mL conical centrifuge tubes with adapters A15 (462092)
- Maximum tube dimensions: 29 x 120 mm
- Maximum speed: 4,400 rpm
- Maximum centrifugal force: 3,200 rcf
- Radius: 148 mm
- Angle: 45°
- Minimum acceleration time: 20 s
- Minimum braking time: 21 s
- Rotor and buckets autoclavable: yes



**30 x 2 mL
Microcentrifuge Tube
Rotor with Lid**

Part number: 228451

- Rotor type: RA 30/2 angle rotor with lid
- Maximum capacity: 30 x 1.5/2.0 mL microcentrifuge tubes - optional 0.2 mL or 0.5 mL adapters for PCR tubes (284351 or 283349)
- Maximum tube dimensions: 11 x 40 mm
- Maximum speed: 13,000 rpm
- Maximum centrifugal force: 18,520 rcf
- Radius: 98 mm
- Angle: 45°
- Minimum acceleration time: 21 s
- Minimum braking time: 21 s
- Rotor and lid autoclavable: yes



**24 x Glass Hematocrit
Rotor**
Part number: 246701

- Rotor type: RH 24 hematocrit rotor with reading device
- Maximum capacity: 24 x glass hematocrit capillary tubes
- Maximum tube dimensions: 1.6 X 75 mm
- Maximum speed: 13,000 rpm
- Maximum centrifugal force: 18,520rcf
- Radius: 98 mm
- Angle: 90°
- Minimum acceleration time: 20 s
- Minimum braking time: 21 s
- Rotor and lid autoclavable: yes



30 x 15 mL Tube Rotor

Part number: 454656

- Rotor type: RA 30/15 angle rotor with shields
- Maximum capacity: 30 x 15 mL open conical tubes or 21 x 15 mL closed conical tubes
- Maximum tube dimensions: 17 x 120 mm
- Maximum speed: 4,400 rpm
- Maximum centrifugal force: 3,200 rcf
- Radius: 148 mm
- Angle: 45°
- Minimum acceleration time: 21 s
- Minimum braking time: 21 s
- Rotor and shields autoclavable: yes



4 x 8 PCR Strip Rotor

Part number: 273335

- Rotor type: RA 6/PCR angle PCR strip rotor
- Maximum capacity: 6 x 8 PCR strips
- Maximum tube dimensions: 9 x 22 mm
- Maximum speed: 13,000 rpm
- Maximum centrifugal force: 18,520 rcf
- Radius: 98 mm
- Angle: 48°
- Minimum acceleration time: 22 s
- Minimum braking time: 21 s
- Rotor and lid autoclavable: yes



16 x 5 mL Microcentrifuge Tube Rotor with Aerosol Lid

Part number: 403055

- Rotor type: RA 16/5 angle rotor with aerosol tight lid
- Maximum capacity: 16 x 5.0 mL microcentrifuge tubes
- Maximum tube dimensions: 16.5 X 75 mm
- Maximum speed: 13,000 rpm
- Maximum centrifugal force: 18,520 rcf
- Radius: 98 mm
- Angle: 45°
- Minimum acceleration time: 22 s
- Minimum braking time: 22 s
- Rotor and lid autoclavable: yes



2 Place Microtitration Plate Rotor

Part number: 270848

- Rotor type: RS 4/100 four place swing out rotor
- Maximum capacity: 4 x 100 mL centrifuge tubes
- Maximum tube dimensions: 44 x 100 mm
- Maximum speed: 4,400 rpm
- Maximum centrifugal force: 3,200 rcf
- Radius: 91 mm
- Angle: 90°
- Minimum acceleration time: 21 s
- Minimum braking time: 22 s
- Rotor and lid autoclavable: yes
- Adapters: buckets B100 (168721) lids for buckets BL100 (168781) and B100XL (322022)



4 Place 100 mL Swing Out Rotor

Part number: 450653

- Rotor type: RS 4/200 four place swing out rotor
- Maximum capacity: 4 x 200 mL centrifuge tubes
- Maximum tube dimensions: 57 x 105 mm
- Maximum speed: 4,400 rpm
- Maximum centrifugal force: 3,200 rcf
- Radius: 91.5 Mm
- Angle: 90°
- Minimum acceleration time: 21 s
- Minimum braking time: 22 s
- Rotor and lid autoclavable: yes
- Adapters: buckets B200 (424173) lids for buckets BL200 (412627)



4 Place 200 mL Swing Out Rotor

Part number: 383672

- Rotor type: RA 24/2 angle rotor with lid
- Maximum capacity: 24 x 1.5/2.0 mL microcentrifuge tubes - optional 0.2 mL or 0.5 mL adapters for PCR tubes (283451 or 283349)
- Maximum tube dimensions: 11 x 40 mm
- Maximum speed: 15,000 rpm
- Maximum centrifugal force: 21,630 rcf
- Radius: 86 mm
- Angle: 45°
- Minimum acceleration time: 22 s
- Minimum braking time: 23 s
- Rotor and lid autoclavable: yes



Cytology Rotor

Part number: 349771

- Rotor type: RC 12 cytology rotor
- Maximum capacity: 12 x 6 mL cytology funnels
- Maximum funnel volume: 6 mL
- Maximum speed: 2,500 rpm
- Maximum centrifugal force: 1,030 rcf
- Radius: 123 mm
- Angle: 90°
- Minimum acceleration time: 19 s
- Minimum braking time: 21 s
- Rotor autoclavable: yes

Rotor tube compatibility

Rotor	RA 24/2 & 24/2 AERO			RA 30/2			RA 8/15			RA 16/5	RM 2/3	
Rotor order no.	228442 or 375944			228451			448807			403055	270848	
Tube type	PCR	PCR	Micro	PCR	PCR	Micro	Blood Collection		Conical	Micro	Standard MTP	
Tube image												
Nominal vol. (mL)	0.2/0.4	0.5	1.5 to 2	0.2/0.4	0.5	1.5 to 2	5	7-10	10-12	15	5	96 x 350 µL max.
Dimensions (mm) (D x L)	6 x 18	8 x 30	11 x 38	6 x 18	8 x 30	11 x 38	13 x 75	13 x 100	16 x 100	17 x 120	16.5 x 75	128 x 86 x 15 128 x 86 x 45
Adapter image			-			-		-	-	-	-	-
Adapter	A0.2	A0.5	-	A0.2	A0.5	-	A5	-	-	-	-	-
No. tube/rotor	24	24	24	30	30	30	8	8	8	8	16	6 or 2
No. tube/adapter	24	24	0	30	30	0	8	0	0	0	-	-
Order no.	238451	284349	-	238451	284349	-	450281	-	-	-	-	168790

Rotor	RA 30/15				RA 30/12				RA 6/50		RH 24	RA 6/PCR
Rotor order no.	454656				454655				448914		246701	273335
Tube type	Blood Collection		Conical		Blood Collection		Conical	Conical	Conical	Capillaries	8-Well PCR Strip	
Tube image												
Nominal vol. (mL)	5	7 - 10	10 - 12	15	5	7 - 10	10 - 12	15	15	50	0.075	0.2
Dimensions (mm) (D x L)	13 x 75	13 x 100	16 x 100	17 x 120	13 x 75	13 x 100	16 x 100	17 x 120	17 x 120	30 x 115	1.6 x 75	-
Adapter image		-	-	-		-	-	-		-	-	-
Adapter	A5	-	-	-	A5	-	-	-	A15	-	-	-
No. Tube/rotor	30	30	30	21	30	30	30	15	6	6	24	48
No. Tube/adapter	30	0	0	0	30	0	0	0	6	0	-	-
Order no.	450283				450283	-	-		462092	-	-	

Rotor	RS 4/100												
Rotor order no.	450653												
Tube type	PCR	PCR	Micro	Open Glass	Blood Collection	Open Glass	Conical	Open Glass	Open Glass	Blood Collection	Conical	Blood Collection	
Tube image													
Nominal vol. (mL)	0.2	0.5	1.5 to 2	5	5	25	50	50	100	10 - 12	15	5 - 7	
Dimensions (mm) (D x L)	6 x 18	8 x 30	11 x 38	12 x 75	12 x 75 13 x 100	24 x 100	30 x 115	36 x 100	44 x 100	16 x 100	17 x 120	13 x 75 13 x 100	
Insert image													
Inserts (4/pk)	I100 12/5	I100 12/5	I100 12/5	I100 12/5	I100 7/5-7	I100 2/25	I100 1/50	I100 1/50	I100 1/100	I100 6/10-12	I100 2/15	I100 12/5-7XL	
No. Tubes/insert	12	12	12	12	7	2	1	1	1	6	2	12	
No. Tubes/rotor	48	48	48	48	28	8	4	4	4	24	8		
Insert order no.	168722	168722	168722	168722	168737	168724	168725	168725	168726	168723	168785		
Autoclavable	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes		
Adapter/reducer	A0.2-0.4	A0.5	A1.5				A50						
Adapter/reducer image													
Radius (mm)	162	162	162	162	163	161	161	161	161	161	163	151	
G force (rcf)	3505	3505	3505	3505	3527	3484	3484	3484	3484	3484	3527	3267	
Adapter order no.	168782	168783	168784				168738					322015	

Rotor	RC 12									
Rotor order no.	349771									
Cytology slide	Rotoclip JC390	Rotofunnel Reusable JC406	Rotofunnel Disposable JC320-500	Filtercards JC407	Cytoslides JC311	ECOfunnel Reusable JC313	ECOfilter JC345	ECOSlides JC314	ECOseal JC315	
Nominal vol. (mL)	-	0.5	0.5	-	-	5	-	-	-	-
Sedimentation fields	-	1 Field round 6 mm	1 Field round 6 mm	-	-	1 Field 22 x 15 mm	-	-	-	-
Positions/rotor	12									
Quantity	4 (3 sets required)	12	500	200	100	12	100	100	100	100
Order no.	312491	496090	403270	484130	484136	503420	503417	503419	503418	











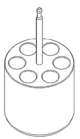
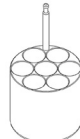
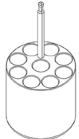
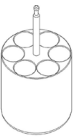
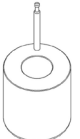

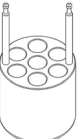

Rotor	RS 4/200								
Rotor order no.	383672								
Tube type	Glass	Blood Collection	Blood Collection BD	Blood Collection Thermo	Blood Collection	Skirted	Conical	Blood Collection BD	Conical
Tube image									
Nominal vol. (mL)	5	5-7	5-7	5-7	10-12	30	50	10-12	15
Dimensions (mm) (D x L)	13 x 75	13 x 75 13 x 100	13 x 75 13 x 100	13 x 75 13 x 100	12 x 75 16 x 100	24 x 100 25 x 110	30 x 115	16 x 100	17 x 120
Insert image									
Inserts (4/pk)	I200 12/5	I200 6/5-7	I200 7/5-7	I200 9/5-7	I200 6/10	I200 1/30	I200 1/50	I200 7/10	I200 3/15
No. Tubes/insert	12	6	7	9	6	1	1	7	3
No. Tubes/rotor	48	24	28	36	24	4	4	28	12
Insert order no.	424176	424177	424178	424179	424180	424182	424175	424181	424174
Autoclavable	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes	Yes
Radius (mm)	156.2	157.2	157.2	155.2	156.2	156.2	157.2	156.2	159.2
G force (rcf)	3380	3402	3402	3358	3380	3380	3402	3380	3445


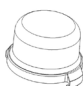
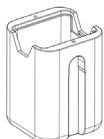


Image		
Code	B200	BL100
Description	Buckets B200	Lids for B200
Quantity	4	4
Autoclavable	Yes	No
Order No.	424173	412627

Image			
Code	B100	BL100	B100 XL
Description	Buckets B100	Lids for B100	Buckets B100 XL
Quantity	4	4	4
Autoclavable	Yes	No	Yes
Order No.	168721	168781	322022



Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA
www.revivity.com

For research use only. Not for use in diagnostic procedures.

For a complete listing of our global offices, visit www.revivity.com
Copyright ©2024, Revvity, Inc. All rights reserved.

www.revivity.com

1414106