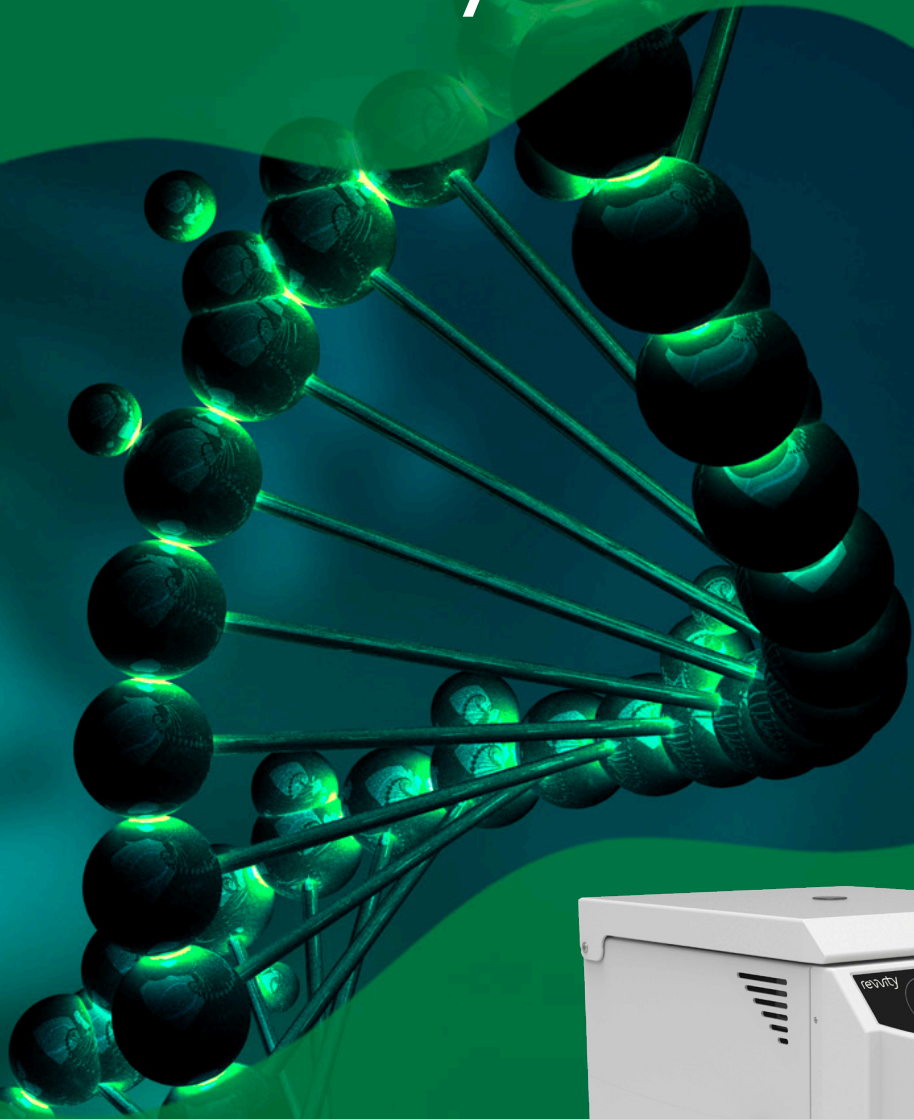


Sample
preparation
made easy.



revvity

CF-220 Centrifuge

Simple and versatile

The CF-220 is an ergonomic and user-friendly clinical centrifuge for separating substances with different relative densities by using centrifugal force. The maximum rotational speed of 15,500 rpm gives the centrifugal force of 23,100 rcf.

With seven different rotors available, it suits various applications. Easy to operate, it streamlines processes while maintaining precision.

The large, illuminated LCD interface is modern and offers great visibility.



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

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

Specifications

Power rating	CF-220-120 - 120V, 50/60 Hz
	CF-220-220 - 220V, 50/60 Hz
	CF-220-220UK - 220V, 50/60 Hz
Max. power consumption	450 watts
Max. rotation speed	200 to 15,500 rpm, in steps of 10 rpm
Maximum G force	23,100 rcf
Maximum capacity	6 x 50 mL
Rotors available	7
Time	10 s to 99 min and continuous, step 10 s
Programs	100
Acceleration levels	10
Deceleration levels	10 (Level 0 without braking)
Display	LCD with backlight
Noise level (at max. speed)	≤ 60 dBA
Fast spin	Yes
Lid locking mechanism	Manual closed, motorized open
Lid Lock	2 Latches
Motor type	Induction brushless motor
Permitted ambient conditions	2 to 35 °C, Humidity: 75 %, Non-Condensing
Dimensions	W: 11" (28.0 cm) D: 14.6" (37.0 cm) H: 8.9" (22.6 cm) (H: lid open 19.4" (49.3 cm))
Weight	31.7 lbs. (14.4 kg)
Warranty	One Year
Certification	CE

Rotors



24 x 2 mL Microcentrifuge Tube Rotor

Part number: 437480

- Rotor type: RA 24/2 angle rotor, transparent plastic 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: 15,500 rpm
- Maximum centrifugal force: 23,100 rcf
- Radius: 86 mm
- Angle: 45°
- Minimum acceleration time: 25 s
- Minimum braking time: 24 s
- Rotor and lid autoclavable: Yes



12 x 5 mL Microcentrifuge Tube Rotor

Part number: 449201

- Rotor type: RA 12/5 angle rotor, transparent plastic lid
- Maximum capacity: 12 x 5 mL microcentrifuge tubes
- Maximum tube dimensions: 16.5 x 75 mm
- Maximum speed: 15,500 rpm
- Maximum centrifugal force 23,100 rcf
- Radius: 82 mm
- Angle: 45°
- Minimum acceleration time: 28 s
- Minimum braking time: 25 s
- Rotor and lid autoclavable: Yes



12 x 12 mL Tube Rotor

Part number: 447993

- Rotor type: RA 12/12 angle rotor with shields
- Maximum capacity: 12 x 12 mL centrifuge tubes or 6 x 15 mL centrifuge tubes - optional A5 adapters for tubes with length 75 to 90 mm (450282)
- Maximum tube dimensions: 17 x 120 mm
- Maximum speed: 6,000 rpm
- Maximum centrifugal force: 4,185 rcf
- Radius: 101.6 mm
- Angle: 35°
- Minimum acceleration time: 18 s
- Minimum braking time: 19 s
- Rotor and shields autoclavable: Yes



8 x 15 mL Tube Rotor

Part number: 438526

- Rotor type: RA 8/15 angle rotor with shields
- Maximum capacity: 8 x 15 mL conical centrifuge tubes - optional A5 adapters for 75 to 90 mm tubes (450281)
- Maximum tube dimensions: 17 x 120 mm
- Maximum speed: 6,000 rpm
- Maximum centrifugal force: 4,185 rcf
- Radius: 101.6 mm
- Angle: 35°
- Minimum acceleration time: 20 s
- Minimum braking time: 23 s
- Rotor and shields autoclavable: Yes



6 x 50 mL Tube Rotor

Part number: 438545

- Rotor type: RA 6/50 angle rotor with shields
- Maximum capacity: 6 x 50 mL conical centrifuge tubes or 6 x 15 mL conical centrifuge tubes with A15 adapters (462092)
- Maximum tube dimensions: 29 x 120 mm
- Maximum speed: 6,000 rpm
- Maximum centrifugal force: 4,185 rcf
- Radius: 104 mm
- Angle: 30°
- Minimum acceleration time: 18 s
- Minimum braking time: 20 s
- Rotor and shields autoclavable: Yes



24 x Glass Hematocrit Rotor

Part number: 449568

- 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: 15,500 rpm
- Maximum centrifugal force: 23,100 rcf
- Radius: 92 mm
- Angle: 90°
- Minimum acceleration time: 28 s
- Minimum braking time: 25 s
- Rotor autoclavable: Yes



4 x 8 PCR Strip Rotor

Part number: 449767

- Rotor type: RA 4/PCR angle PCR strip rotor
- Maximum capacity: 4 x 8 PCR Strips
- Maximum tube dimensions: 9 x 22 mm
- Maximum speed: 15,500 rpm
- Maximum centrifugal force: 23,100 rcf
- Radius: 56 mm
- Angle: 45°
- Minimum acceleration time: 24 s
- Minimum braking time: 23 s
- Rotor autoclavable: Yes

Rotor tube compatibility

Rotor	RA 24/2				RA 12/5	RA 12/12			RA 8/15			
Rotor order no.	437480				449201	447993			438526			
Tube type	PCR	PCR	Micro	Micro	Blood Collection			Conical	Blood Collection		Conical	
Tube image												
Nominal vol. (mL)	0.2/0.4	0.5	1.5 to 2	5	5	7 - 10	10 - 12	15	5	7 - 10	10 - 12	15
Dimensions (mm) (D x L)	6 x 18	8 x 30	11 x 38	16 x 75	13 x 75	13 x 100	16 x 100	17 x 120	13 x 75	13 x 100	16 x 100	17 x 120
Adapter image			-	-		-	-	-		-	-	-
Adapter	A0.2	A0.5	-	-	A5	-	-	-	A5	-	-	-
No. tube/rotor	24	24	24	12	12	12	12	6	8	8	8	8
No. tube/adapter	24	24	0	-	12	0	0	0	8	0	0	0
Order no.	238451	284349	-	-	450282	-	-	-	450281	-	-	-

Rotor	RA 6/50		RH 24	RA 4/PCR
Rotor order no.	438545		449568	449767
Tube type	Conical	Conical	Capillaries	8-Well PCR Strip
Tube image				
Nominal vol. (mL)	15	50	0.075	0.2
Dimensions (mm) (D x L)	17 x 120	30 x 115	1.6 x 75	-
Adapter	A15	-	-	-
No. tube/rotor	6	6	24	32
No. tube/adapter	6	0	-	-
Order no.	462092	-	-	-



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

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

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

www.revvity.com