

The **ideal choice** for cell culture
and general purpose separations.

Allegra 6 Series Benchtop Centrifuges

Blood Banking
Capillary Electrophoresis
Cell Analysis
Centrifugation
Genomics
Lab Automation
Lab Tools
Particle Characterization



The Allegra 6 Series Benchtop Centrifuges provide versatile, high-capacity performance for a wide variety of cell culture, plasma and general purpose separations.



4 x 750 mL
GH-3.8
Swinging-
bucket Rotor

GH-3.8 Swinging-bucket Rotor

- 4 x 750 mL Maximum Capacity
- 11 Different Adapters — Spin From 1.5 mL Tubes to 750 mL Bottles
- Bucket Covers Certified* for Biocontainment
- Up to 3,210 x g

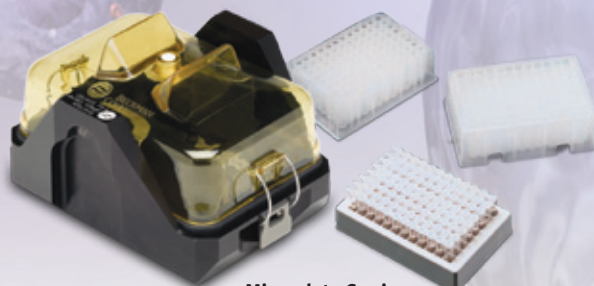
Aerosolve Canisters and Lids for Maximum Protection— Certified Biocontainment

- Clear Plastic Body and Special Adapters Allow Tube Inspection While Canister Remains Sealed
- Autoclavable up to 121°C
- Available adapters from 1.5 mL Tube to 230 mL Conical Bottle



Microplate Carriers — Perfect for Cell Adhesion

- For Use with GH-3.8 Swinging-bucket Rotor
- Holds 16 Standard Microplates, 4 Deep-well (1.8 mL) or 4 Square-well (2.0 mL) Plates
- Up to 2,980 x g



Microplate Carriers
for GH-3.8 Rotor



Modular Disk
Adapters for
GH-3.8 Rotor

Specifications

	Allegra 6R	Allegra 6	Allegra 6KR
Refrigeration	Yes, CFC-Free	No	Yes, CFC-Free
Maximum Speed	6,100 rpm (Fixed-angle) 3,750 rpm (Swinging-bucket)	6,100 rpm (Fixed-angle) 3,750 rpm (Swinging-bucket)	6,100 rpm (Fixed-angle) 3,750 rpm (Swinging-bucket)
Maximum <i>g</i> -Force	5,710 x <i>g</i> (Fixed-angle) 3,210 x <i>g</i> (Swinging-bucket)	5,710 x <i>g</i> (Fixed-angle) 3,210 x <i>g</i> (Swinging-bucket)	5,710 x <i>g</i> (Fixed-angle) 3,210 x <i>g</i> (Swinging-bucket)
Maximum Capacity	4 x 750 mL	4 x 750 mL	4 x 750 mL
Run Time	Timed (0-30 min), Hold	Timed (0-30 min), Hold	Timed (0-30 min), Hold
Drive	Direct, Brush Type	Direct, Brush Type	Direct, Brush Type
Noise at Max Speed	≤ 65 dBa	≤ 65 dBa	≤ 65 dBa
Operating Temperature	-5°C to +25°C	Ambient	-5°C to +25°C
Set Temperature Range	-10°C to +40°C	N/A	-10°C to +40°C
Dimensions (cm) Lid Closed	76W x 66D x 39.4H	54.6W x 66D x 40.6H	54.6W x 66D x 67.3H
Dimensions (in) Lid Closed	30W x 26D x 15.5H	21.5W x 26D x 16H	21.5W x 26D x 26.5H
Weight without Rotor	75.8 kg (167 lbs)	51.7 kg (114 lbs)	90.3 kg (199 lbs)

Ordering Information

Part No.	Description
366816	Allegra 6R Centrifuge, 120V/60Hz
366802	Allegra 6 Centrifuge, 120V/60Hz
366830	Allegra 6KR Centrifuge, 120V/60Hz

Rotors and Accessories

Part No.	Description	Max Speed (rpm)	Max RCF, <i>g</i>
360581	4 x 750 mL Swinging-bucket Rotor (GH-3.8) -11 Different Adapters Available	3,750	3,210
360585	Bucket Covers for GH-3.8 only (Set of 2)		
359232	Aerosolve Canisters for GH-3.8 only (Set of 4) -10 Different Adapters Available		
392806	Microplate Carriers (Set of 2) -Each Carrier Accommodates up to 4 Standard Microplates or 1 Deep-well Plate	3,750	2,890
367054	6 x 250 mL Fixed-angle Rotor (GA-6) -5 Different Adapters Available	6,100	5,710

For a complete listing of adapters, visit www.beckmancoulter.com/centrifuge



* **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 (Health Protection Agency, Porton Down, UK or USAMRIID, Ft. Detrick, MD, USA). Improper use or maintenance may affect seal integrity and, thus, containment.

Beckman Coulter, the stylized symbol, Allegra and Aerosolve are trademarks of Beckman Coulter, Inc. and are registered in the USPTO.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at www.beckmancoulter.com

B2010-11848

© 2011 Beckman Coulter, Inc.

PRINTED IN U.S.A.