

Allegra X-15R and X-12 Series



Brilliance
at every turn.





Allegra X-15R

The best swinging bucket in the industry.

MAXIMUM PERFORMANCE.

If you operate a high-throughput lab, get to know the Allegra X-15R. Able to process up to 148 tubes per cycle, this centrifuge delivers maximum performance and cost-effective versatility through a wide variety of available adapters and accessories.

With a powerful, built-in refrigeration system, the X-15R cools more quickly and maintains temperature even at maximum speed. The ARIES Smart Balance Rotor System automatically detects and corrects up to 50 grams of imbalance, which not only reduces mid-run interruptions to save you time, but also helps to minimize instrument wear and tear. BioCertified[†] rotors, covers, and canisters are tested to ensure potentially harmful aerosols are contained, maintaining safety in the lab.



Allegra X-12 Series

Streamlined, simplified separations.

STREAMLINE AND SIMPLIFY.

You'll also find the ARIES Smart Balance Rotor System, as well as BioCertified[†] rotors, covers, and canisters, in our Allegra X-12 Series. Offering maximum capacity, these centrifuges are ideal for streamlining general separations. Whether harvesting cells or preparing blood samples, the X-12 will simplify your processes and save you time.

SPECIFICATIONS

	ALLEGRA X-15R	ALLEGRA X-12R	ALLEGRA X-12
Swinging Bucket Rotor	4,750 rpm; 5,250 x g	3,750 rpm; 3,270 x g	3,750 rpm; 3,270 x g
Fixed-Angle Rotor	10,200 rpm; 11,400 x g	10,200 rpm; 11,400 x g	10,200 rpm; 11,400 x g
Refrigeration	CFC-Free	CFC-Free	Constant Controlled Temp.
Maximum Capacity	3 Liters	3 Liters	3 Liters
Run Time	Timed (Up to 99 hrs, 59 min) Hold, Pulse (Short Run)	Timed (Up to 99 hrs, 59 min) Hold, Pulse (Short Run)	Timed (Up to 99 hrs, 59 min) Hold, Pulse (Short Run)
Drive	Brushless Induction	Brushless Induction	Brushless Induction
Programs	10 user-defined	10 user-defined	10 user-defined
Accel/Decel Profiles	10/11	10/11	10/11
Temperature Range	-10° C to +40° C (selectable)	-10° C to +40° C (selectable)	Preset to +20° C*
Dimensions (cm), Lid Closed	76.2W x 62.2D x 34.3H	76.2W x 62.2D x 34.3H	76.2W x 62.2D x 34.3H
Dimensions (in), Lid Closed	30W x 24.5D x 13.5H	30W x 24.5D x 13.5H	30W x 24.5D x 13.5H
Weight Without Rotor	128 kg (283 lb)	121 kg (267 lb)	121 kg (267 lb)
Electrical	208V, 60 Hz; 230V, 50 Hz; 200V, 50/60 Hz	208V, 60 Hz; 230V, 50 Hz; 200V, 50/60 Hz	208V, 60 Hz; 230V, 50 Hz; 200V, 50/60 Hz

* The constant-temperature Allegra X-12 utilizes a refrigeration system pre-set to +20° C.

ORDERING INFORMATION

BENCHTOP CENTRIFUGE MODELS						
Part No.	Description	Max Speed (rpm)		Max RCF, x g		
		Swinging Bucket/Fixed Angle		Swinging Bucket/Fixed Angle		
392932	Allegra X-15R Centrifuge, 208V, 60 Hz	4,750 / 10,200		5,250 / 11,400		
392934	Allegra X-15R Centrifuge, 230V, 50 Hz	4,750 / 10,200		5,250 / 11,400		
392933	Allegra X-15R Centrifuge, 200V, 50/60 Hz	4,750 / 10,200		5,250 / 11,400		
392302	Allegra X-12R Centrifuge, 208V, 60 Hz	3,750 / 10,200		3,270 / 11,400		
392304	Allegra X-12R Centrifuge, 230V, 50 Hz	3,750 / 10,200		3,270 / 11,400		
392303	Allegra X-12R Centrifuge, 200V, 50/60 Hz	3,750 / 10,200		3,270 / 11,400		
392472	Allegra X-12 Centrifuge, 208V, 60 Hz	3,750 / 10,200		3,270 / 11,400		
392474	Allegra X-12 Centrifuge, 230V, 50 Hz	3,750 / 10,200		3,270 / 11,400		
392473	Allegra X-12 Centrifuge, 200V, 50/60 Hz	3,750 / 10,200		3,270 / 11,400		
ROTORS						
Part No.	Description	Max Speed (rpm)		Max RCF, x g		
Swinging Bucket Rotors		X-15R/X-12 Series		X-15R/X-12 Series		
369702	4x750 mL Swinging Bucket Rotor Assembly w/o Bucket Covers (SX4750)	4,750 / 3,750		5,250 / 3,270		
369704	4x750 mL ARIES Self-Balancing Swinging Bucket Rotor Assembly w/o Bucket Covers (SX4750A)	4,750 / 3,750		5,250 / 3,270		
392806	Microplate Carriers for SX4750 & SX4750A Rotor (SX4750 _μ) 16 Standard Microplates, 4 Deep-Well/Square-Well Plates, 4 of 96-Tube Racks (Set of 2), Includes Carriages	4,450 / 3,750		4,060 / 2,885		
Fixed-Angle Rotor						
369735	BioCertified [®] 6x100 mL Fixed-Angle Rotor (FX6100)	10,200		11,400		
Rotor Accessories						
392805	BioCertified [®] Covers for SX4750 & SX4750A Round Buckets (Set of 2)	—		—		
393070	BioCertified [®] Covers for SX4750 _μ Microplate Carriers (Set of 2)	—		—		
392873	Replacement Carriages for SX4750 _μ Microplate Carriers, Includes Rubber Pads (Set of 2)	—		—		
TUBE AND BOTTLE ADAPTERS						
Part No.	Tube Volume (mL)	Tube Diameter (mm)	Color	Bucket Capacity	Rotor Capacity	Pack of
Adapters for Direct Centrifugation of Cell Culture Flasks (SX4750 & SX4750A)						
369292	Corning T-75 cm ² Canted-Neck Cell Culture Flasks**	—	Orange	1	4	2
369295	Corning T-25 cm ² Canted-Neck Cell Culture Flasks**	—	Green	2	8	2
Modular Disk Adapters for Tubes (SX4750 & SX4750A)						
359469 & 359148	3 mL & 5 mL	10 & 12	Blue	37	148	2 & 4
343108	Tube Decanter for P/N 359469 & 359148	—	—	—	—	Each
354511	1.5 mL Adapter Plate for P/N 359469 & 359148	—	—	—	—	Each
359478 & 359157	3 mL & 5 mL	13	Tan	30	120	2 & 4
359470 & 359149	7 mL & 10 mL	14	Orange	24	96	2 & 4
359471 & 359150	12 mL	16	Purple	19	76	2 & 4
359472 & 359151	15 mL (Conical)	18	Green	14	56	2 & 4
359473 & 359152	15 mL & 20 mL (Round)	18	Green	14	56	2 & 4
359475 & 359154	30 mL (Round) & 50 mL (Conical)	30	Light Green	4	16	2 & 4
393266 & 393267	50 mL (Conical)	29	Grey	7	28	2 & 4
359474 & 359153	50 mL (Round)	29	Yellow	7	28	2 & 4
359476 & 359155	50 mL (Round)	35	Dark Blue	4	16	2 & 4
359477 & 359156	100 mL	44	Brown	2	8	2 & 4

(Tube and bottle adapters continued on back)

** Corning is a trademark of Corning Incorporated.

† Validation of microbiological containment was done at an independent, third-party facility (CAMR, Porton Down, UK or USAMRIID, Ft. Detrick, MD, USA). Improper use or maintenance may affect seal integrity and, thus, containment.

BIOC 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.

TUBE AND BOTTLE ADAPTERS (continued)

Part No.	Tube Volume (mL)	Tube Diameter (mm)	Color	Bucket Capacity	Rotor Capacity	Pack of
Adapters for Bottles (SX4750 & SX4750A)						
356983	230 mL (Conical)-Insert for Use with 349946	62	Orange	1	4	Each
349946	250 mL Flat Bottom	62	Yellow	1	4	Each
349849	250 mL (Corning Conical)	62	Warm Red	1	4	Each
349945	500 mL Flat Bottom	70	Tan	1	4	Each
349846	750 mL Flat Bottom (Can also be used with 500 mL Corning Conical Tubes with Corning Cushion)	96	Blue	1	4	Each
BioCertified[†] Aerosolve Canisters and Tube Racks (SX4750 & SX4750A)						
359481 & 359232	BioCertified [†] Aerosolve Canisters	—	—	—	—	2 & 4
354495	1.5 mL	11	White	24	96	Each
359482 & 359160	3 mL & 5 mL	12	Blue	24	96	2 & 4
359489 & 358993	5 mL	12	Tan	24	96	2 & 4
359483 & 359161	10 mL	14	Orange	18	72	2 & 4
359484 & 359162	12 mL, 3 mL & 5 mL	16, 12	Purple	12, 6	48, 24	2 & 4
344517	15 mL	14	White	10	40	Each
359485 & 359163	15 mL & 20 mL (Round), 3 mL & 5 mL	18, 12	Green	12, 6	48, 24	2 & 4
359487 & 358991	15 mL (Conical), 3 mL & 5 mL (Round)	17, 12	Light Green	6, 6	24, 24	2 & 4
359486 & 359164	50 mL, 3 mL & 5 mL	29, 12	Yellow	4, 4	16, 16	2 & 4
359488 & 358992	50 mL (Conical), 3 mL & 5 mL (Round)	30, 12	Lime Green	4, 4	16, 16	2 & 4
356985	230 mL (Conical)	62	Orange	1	4	2
Adapters for Tubes (FX6100)						
344497	1.5 mL/2.0 mL (must be inserted into P/N 392830)	—	White	—	36	Each
392824	10 mL	16	White	—	6	6
392823	15 mL	18	White	—	6	6
392270	15 mL (Conical)	17	Beige	—	6	6
392822	30 mL	25.7/25	White	—	6	6
392821	35 mL	29	White	—	6	6
392830	50 mL	29	Blue	—	6	6
392268	50 mL (Conical)	30	Beige	—	6	6

Accommodates 100 mL, 94 mL, 85 mL, 80 mL and 70 mL tubes without adapters, directly in rotor cavity.

With over three-quarters of a century in developing scientific instruments, Beckman Coulter has earned the sobriquet of gold standard from our customers around the globe. We've built a reputation of innovation by challenging convention and redefining standards of excellence. The Allegra suite of instruments is no exception. These general purpose, benchtop mainstays perform with endurance under the most exacting standards—their sheer longevity is a form of built-in sustainability. And each instrument is a precision-made beauty worthy of the Beckman Coulter name.

[†] Validation of microbiological containment was done at an independent, third-party facility (CAMR, Porton Down, UK or USAMRIID, Ft. Detrick, MD, USA). Improper use or maintenance may affect seal integrity and, thus, containment.

BIO **Certified** 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 logo, Aerosolve and Allegra 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

B2011-12485-5K-LC

© 2011 Beckman Coulter, Inc.

PRINTED IN U.S.A.

