

BENCH SERIES OF BIOSAFE† CENTRIFUGE SYSTEMS

Beckman Coulter introduces a series of BioSafe rotors and labware†† as part of a multilevel approach to containment against biological agents and infectious substances. We confidently stand behind this first line of defense, designed with your applications in mind, to make your job safer.



BIO SAFETY CERTIFIED ROTORS

	PART NO.	ROTOR TYPE	MAX SPEED (rpm)	MAX g FORCE	NO. TUBES x VOLUME	ROTOR CAPACITY
Allegra® X-15R and Allegra X-12 Series	369704	SX4750A ARIES™ Smart Balance Rotor (With Cover P/N 392805)	4,750*	5,250*	4 x 750 mL***	3 L
	369702	SX4750 (With Cover P/N 392805)	4,750*	5,250*	4 x 750 mL***	3 L
	392806	Microplate Carriers for SX4750 Rotor (With Cover P/N 393070)	4,450**	4,060**	4 x 3 std plates 4 x 1 deep well plates	
	369735	FX6100	10,200	11,400	6 x 100 mL	600 mL
Allegra 25R	368308	TS-5.1-500 (With Cover P/N 368472)	5,100	5,530	4 x 500 mL****	2 L
	368298	TA-15-1.5	15,000	25,160	30 x 1.5 mL	45 mL
Allegra X-22 Series	361171	F2402H	15,500	22,065	24 x 1.5 mL	36 mL
Allegra 64R	361171	F2402H	26,000	61,970	24 x 1.5 mL	36 mL
	363000	H6002	12,200	12,400	60 x 1.5 mL	90 mL
Microfuge® 22R	365630	F241.5	14,000	17,500	24 x 1.5 mL	36 mL
	368894	F301.5	14,000	21,920	30 x 1.5 mL	45 mL
	368898	F40.25	14,000	17,500	40 x 0.25 mL	10 mL

* 3,750 RPM & 3,270 x g in Allegra X-12 Series

** 3,750 RPM & 2,890 x g in Allegra X-12 Series

*** Variety of adapters available for tubes 1.5 mL to 750 mL

**** Variety of adapters available for tubes 1.5 mL to 500 mL

LABWARE

	PART NO.	DESCRIPTION
Tube and Bottle Bucket Covers	392805	Covers for 750 mL Round Buckets (set of 2) (for SX4750 and SX4750A Swinging Rotor Buckets)
	393070	Covers for SX4750µ Microplate Carriers (set of 2) (Compatible with SX4750)
	368472	Covers for 500 mL Round Buckets (set of 2) (TS-5.1-500 Rotor)
	359481	Aerosolve™ Cannisters (set of 2) (for SX4750 and SX4750A Rotors)



† BioSafe is a term intended to describe the enhanced Biosafety features of our products

†† Validation of microbiological containment was done at an independent third-party testing facility (CAMR, Porton Down, UK, or USAMRIID, Ft. Detrick, MD, USA)

DECONTAMINATION GUIDELINES

Normal centrifuge operation may involve the use of samples and materials that are infectious or pathogenic. Such biohazardous samples may contain fungi, bacteria, virus, or prions and should not be used in a centrifuge until all necessary safety precautions are taken. It is very important to observe the laboratory safety practices as published by your Laboratory Safety Officer, and/or consult the World Health Organizations Laboratory Biosafety Manual or other appropriate sources. It is sometimes necessary to decontaminate centrifuge system components due to a spill or to comply with general laboratory procedures. Decontamination protocols vary greatly depending on the pathogen type needed to be sterilized, and these must be matched to the chemical and/or heat resistance of the materials used in centrifuges, rotors, and labware. Refer to the individual centrifuge and rotor user manuals and publication IN-175, Chemical Resistances for Beckman Coulter Centrifuge Products, for more information about decontamination procedures.

For additional information, please contact your local Beckman Coulter representative or visit our web site at:

www.beckmancoulter.com/biosafe



Developing innovative solutions in Systems Biology.

Innovate Automate
SIMPLIFY

Beckman Coulter, Inc. • 4300 N. Harbor Boulevard, Box 3100 • Fullerton, California 92834-3100
Sales & Service: 1-800-742-2345 • Telex: 678413 • Fax: 1-800-643-4366 • www.beckmancoulter.com

Worldwide Biomedical Research Division Offices:

Australia (61) 2 9844-6000 **Canada** (905) 819-1234 **China** (86) 10 6515 6028 **Eastern Europe, Middle East, North Africa** (41) 22 994 07 07
France 01 49 90 90 00 **Germany** 49 21 513335 **Hong Kong** (852) 2814 7431/2814 0481 **Italy** 02-953921 **Japan** 03-5404-8359
Mexico (55) 560-57770 **Netherlands** 0297-230630 **Singapore** (65) 6339 3633 **South Africa, Sub-Saharan Africa** (27) 11-805-2014/5 **Spain** 91 3836080
Sweden 08-564 85 900 **Switzerland** 0800 850 810 **Taiwan** (886) 2 2378 3456 **Turkey** 90 216 309 1900 **U.K.** 01494 441181 **U.S.A.** 1-800-742-2345