

## BIOSAFE† CENTRIFUGE SYSTEMS - HEPA FILTERS

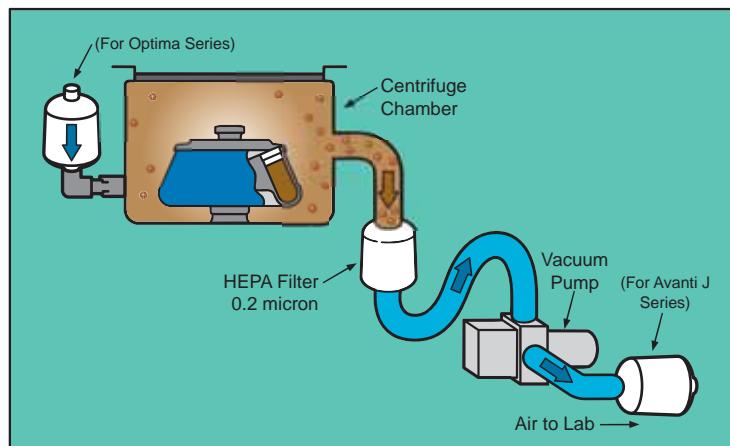
Beckman Coulter introduces a multilevel system of containment for biomedical and microbiological laboratories, Biosafety Levels labs, and vaccine production facilities. Containment of aerosols and spills is achieved in the centrifuge with HEPA (High Efficiency Particle Arresting) filters introduced in the design of the instrument. The HEPA filter provides enhanced aerosol protection from leakage of biohazardous samples such as blood-borne pathogens and other infectious/hazardous material.

HEPA air filtration operates by filtering air passing from the rotor chamber to the vacuum pump, and from the atmosphere to the chamber during venting. HEPA filter kits consist of two disposable filters that have double-layer Emflon® II (Pall) polyvinylidene fluoride (PVDF) membranes for materials compliance in biological tests with a microbial rating greater than 0.2 microns. A solenoid-controlled valve is used to direct the airflow through the filters. Each HEPA filter carries a certificate of test for Pharmaceutical – Grade Sterilizing Filters from the manufacturer.

The Avanti® J Series High Performance Centrifuges have a filter assembly located between the rotor chamber and the vacuum pump. The assembly accommodates two filters: one for evacuation, the other for venting.

The Optima™ Series floor model and tabletop model ultracentrifuges have one filter assembly located between the vacuum pump and diffusion pump for evacuation and a second located on the rotor chamber for venting.

Under normal use, we recommend HEPA filters be replaced once per year by your service rep. Exceptions include clogging, aqueous contact or contamination with sample leakage in the chamber, which requires immediate replacement.



HEPA FILTER FLOW PATH

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

\*All trademarks are the property of their respective owners

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

[www.beckmancoulter.com/biosafe](http://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



