



## Enhanced Biosafety for Your Lab Users During Centrifugation

Containment of biological agents and infectious substances is an essential element in maintaining a safe environment for biomedical and microbiological laboratories, Biosafety Levels labs, and vaccine production facilities. Our BioCertified\*\* Aerosolve Canisters and Bucket Covers offer:

- A transparent design that enables you to check for broken labware
- Protection from aerosol leakage and spills from rotor buckets
- Biocontainment\* that has been third-party tested
- Bucket covers for 1.5 mL tubes to 750 mL bottles
- Microplate carrier covers for 96-well plate format

Visit [beckman.com](http://beckman.com) to learn more.



\* BioSafe is a term intended to describe the enhanced Biosafety features of our products.



\*\* 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.

© 2020 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service names mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at [beckman.com](http://beckman.com)

CENT-694BADV04.20

