



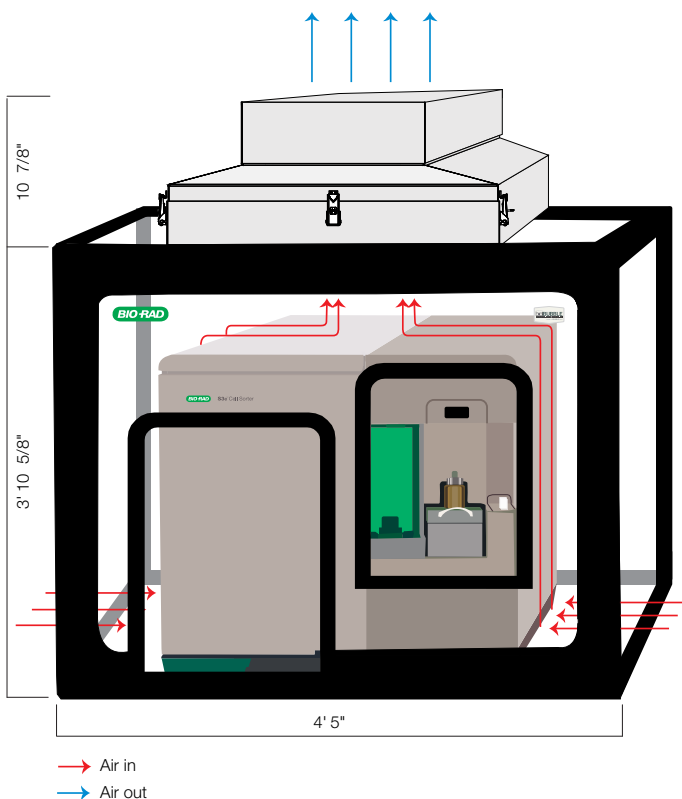
- **Unique design** provides an easy method to access all sides of the instrument for cell sorting, cleaning, and service
- **Designed to reduce vibration** impact to the cell sorter inside
- **Affordable** alternative to traditional large and expensive biosafety hoods offered for cell sorters

Introducing the S3e™ Biosafety System Class I



The S3e Biosafety System is a Class I aerosol containment hood that works in conjunction with ProSort™ Software, providing users maximum protection from aerosols created during the cell sorting process.

The S3e Biosafety System Class I



Airflow is drawn from all four sides and through the access doors when opened, allowing maximum air circulation. The air within the system is exchanged approximately six to eight times a minute. The HEPA filter used in the system provides 99.9997% efficiency at 0.3 μ m. If the airflow through the HEPA filter falls below required limits for containment, the software will alert the user and stop the sort.

Benchtop Design

The S3e Biosafety System's small footprint complements the compact S3™ and S3e Cell Sorters and allows quick access and easy interaction with the instrument. The system is easily assembled and light enough to move if necessary.

ProSort Software Controlled

The S3e Biosafety System is fully integrated and monitored by the ProSort Software. To maintain containment, the fan speed is regulated based on the cell sorter operation mode and the airflow rate through the HEPA filter.

Fully Accessible

Vinyl walls with magnetic sealing are easily opened on all four sides to access any part of the S3 or S3e Cell Sorter for cleaning and maintenance.

Quiet and Energy Efficient

The S3e Biosafety System generates low vibrations and sound pressure levels. Specialized fans and low backpressure from the HEPA filter reduce energy usage.

S3e Biosafety System Class I

Serial #: 779BR0002

Firmware: 000A

Status: OK

On

Exhaust

Filter Life : 100.00%

Temperature : 29.30°C



ProSort Software includes an icon that allows users to view the status of the S3e Biosafety System. The software monitors HEPA filter life and the temperature of the system while it's running.