



Sort Smarter

The S3e Cell Sorter provides exceptional sort purity, recovery, and yield without compromising performance and sensitivity. The S3e Cell Sorter is equipped with one or two lasers and up to four-color detection, making it ideal for sorting cells expressing fluorescent proteins, such as green fluorescent protein, or cells labeled with fluorescent markers.

The S3e Cell Sorter offers researchers:

- **Simplified instrument setup** — ProDrop™ technology automates drop delay calculation and droplet break-off monitoring, enabling precise 1–4 color sorting with minimal training
- **Hands-free alignment** — AutoGimbal™ System features fully automated nozzle tip alignment that is driven by software
- **Compact design** — only 70 x 65 x 65 cm with onboard complete fluidics and Peltier solid-state temperature control
- **Intuitive interface** — user-friendly software for effortless instrument control and sort logic definition

Visit bio-rad.com/web/S3eSortSmarter for more information.



BIO-RAD

Safety – Minimizing Risks

While the S3e Cell Sorter has an interlocked sort chamber door to protect from aerosols formed by the sort stream in the collection area, safety is always a key concern. Adhering to National Institutes of Health (NIH) biosafety standards, a biosafety cabinet is the primary containment system recommended. The S3e Cell Sorter fits seamlessly in the custom-designed S3e Biosafety System Class I. Fully integrated with the ProSort™ Software, the system airflow is controlled and the HEPA filter is monitored in real time. The S3e Biosafety System Class I is uniquely designed with four magnetically attached vinyl walls, which allows easy access to the instrument inside for service or routine maintenance. This option is available to be compliant with NIH standards. Feel confident in protecting both the user and the surrounding environment when performing sort experiments in an S3e Biosafety System Class I.

Support – Technical and Service Care

Rest assured that Bio-Rad's recognized technical and product service support is focused on customer needs. Experienced service engineers will provide either on-call or on-site product and instrument support for instruments under warranty or service contract. It is Bio-Rad's long-term commitment to science and the scientific community that stands behind each Bio-Rad-labeled product.

Instrument support includes:

- Instrument installation
- Installation qualification and operational qualification
- Service contracts and preventive maintenance plans
- Technical application support
- Time and material billable services

Technical support scientists can be reached via our support hotline at **1-800-424-6723**.

Specifications

Droplet frequency	37–43 kHz
Sorting type	True jet-in-air for high-performance sorting
Sorting rate	No hardware limitations for sort rate, limited only by droplet frequency and application
Sorting purity	>99% pure
Nozzle size	100 µm
Sorting direction	2-way sorting
Sorting collection	Up to 5 x 5 ml sample tubes each direction Up to 5 x 1.5 ml tubes each direction Microscope slides 8-well strip each direction
Lasers	Single laser: 488 nm 100 mW Single-laser upgrade: 488 nm 100 mW plus 561 or 640 nm 100 mW Dual laser: 488 and 561 nm 100 mW; 488 and 640 nm 100 mW
Detection	Forward scatter (FSC) with photomultiplier tube (PMT) Side scatter (SSC) with PMT Up to 4 fluorescence detectors with PMT Minimum resolution: 0.5 µm
Sensitivity	<125 MESF for fluorescein isothiocyanate and phycoerythrin
Temperature range	Sample and collection temperature control system: 4–37°C Peltier solid-state system
Fluidics	Onboard fluidics and dilution of ProFlow™ Sort Grade 8x Sheath Fluid with deionized water
Data format	FCS 3.1
Dimensions (W x D x H)	70 x 65 x 65 cm (2.3 x 2.1 x 2.1 ft)

Ordering Information

Catalog #	Description
145-1005	S3e Cell Sorter , 488 nm 100 mW laser, includes 2 fluorescence detectors with filters, AutoGimbal System, 2 fluidic containers with connectors and tubing (sheath, water, waste), power cord, ProSort Software
145-1006	S3e Cell Sorter , 488 and 561 nm 100 mW lasers, includes 4 fluorescence detectors with filters, AutoGimbal System, 2 fluidic containers with connectors and tubing (sheath, water, waste), power cord, ProSort Software
145-1008	S3e Cell Sorter , 488 and 640 nm 100 mW lasers, includes 4 fluorescence detectors with filters, AutoGimbal System, 2 fluidic containers with connectors and tubing (sheath, water, waste), power cord, ProSort Software
145-1078	S3e Biosafety System Class I
145-1029	S3e Cell Sorter (488 nm) with S3e Biosafety System Class I
145-1030	S3e Cell Sorter (488/561 nm) with S3e Biosafety System Class I
145-1032	S3e Cell Sorter (488/640 nm) with S3e Biosafety System Class I
145-1065	S3e Accessory Kit , includes 100 µm nozzle tip, 2 nozzle O-rings, 2 nozzle alignment disks, 1 ml syringe, 2 neutral density filters (1.0), 2 mm hex driver, spanner wrench
145-1082	ProFlow Sort Grade 8x Sheath Fluid , 5 x 4 L containers, preservative free
145-1083	ProFlow Sort Grade Water , 5 x 4 L containers, sterile-filtered, endotoxin-tested water
145-1084	S3e Fluidic Container , 3 x 4 L containers, gamma-irradiated, sterile
145-1085	ProLine™ Rainbow Beads , 1 x 5 ml bottle, mixture of 3.0–3.4 µm beads dyed with 8 different fluorescence intensities
145-1086	ProLine Universal Calibration Beads , 3 x 5 ml bottles



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 43 1 877 89 01 **Belgium** 03 710 53 00 **Brazil** 55 11 3065 7550
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 49 89 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 (0) 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026674 55 05
Taiwan 886 2 2578 7189 **Thailand** 1800 88 22 88 **United Kingdom** 020 8328 2000