



## Cell Sorting Made Simple

The S3 cell sorter is expertly engineered to be affordable without compromising high performance and sensitivity. The S3 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 S3 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
- **Reduced carryover** — dual-position loading stage ensures the sample line is clean between sort runs
- **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

For more information, visit us on the Web at [www.bio-rad.com/cellsorter](http://www.bio-rad.com/cellsorter).



**BIO-RAD**

## Specifications

Acquisition rate	Up to 100,000 events/sec
Sorting rate	Up to 30,000 events/sec
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	Primary: 488 nm 100 mW Optional: 561 nm 100 mW
Detection	Forward scatter (FSC) with PMT Side scatter (SSC) with PMT Up to 4 fluorescence detectors with PMT Minimum resolution: 0.5 µm
Sensitivity	<125 MESF for FITC and PE
Temperature range	Sample and collection temperature control system: 4–37°C Peltier solid state system
Fluidics	Onboard fluidics and dilution of ProFlow™ 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
145-1001	<b>S3 Cell Sorter</b> , 488 nm 100 mW laser, includes 2 fluorescence detectors with filters, 2 fluidic containers with connectors and tubing (sheath, water, waste), power cord, ProSort™ software
145-1002	<b>S3 Cell Sorter</b> , 488 and 561 nm 100 mW lasers, includes 4 fluorescence detectors with filters, 2 fluidic containers with connectors and tubing (sheath, water, waste), power cord, ProSort software
145-1065	<b>S3 Cell Sorter Accessory Kit</b> , 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
145-1081	<b>ProLine™ Calibration Beads</b> , 3 x 5 ml bottles
145-1082	<b>ProFlow 8x Sheath Fluid</b> , 5 x 4 L containers, preservative free
145-1083	<b>ProFlow Sort Grade Water</b> , 5 x 4 L containers, sterile-filtered, endotoxin-tested water
145-1084	<b>S3 Fluidic Containers</b> , 3 x 4 L containers, gamma-irradiated, sterile
145-1085	<b>ProLine Rainbow Beads</b> , 1 x 5 ml bottle, mixture of 3.0–3.4 µm beads dyed with 8 different fluorescence intensities



**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699  
**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** 089 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** 03 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** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 026 674 55 05  
**Taiwan** 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000