

S3e™ Cell Sorter (488/640 nm lasers)

| Fluorophore or Dye | Laser | Ex Max, nm | Em Max, nm | Fluorescence Emission Color | | | |
|--------------------------------------|-------|------------|------------|-----------------------------|----------------|----------------|--------------|
| | | | | FL1 502–550 | FL2 573–613 | FL3 655–685 | FL4 700LP |
| EGFP | 488 | 488 | 507 | • | | | |
| ReadiLink 492/516 (1351002)* | 488 | 492 | 516 | • | | | |
| CFDA-SE (1351201)** | 488 | 492 | 517 | • | | | |
| DyLight 488 | 488 | 493 | 518 | • | | | |
| FITC | 488 | 495 | 519 | • | | | |
| Alexa Fluor 488 | 488 | 499 | 519 | • | | | |
| VivaFix™ 498/521 (1351115)*** | 488 | 498 | 521 | • | | | |
| CytoTrack™ Green 511/525 (1351203)** | 488 | 511 | 525 | • | | | |
| EYFP | 488 | 514 | 527 | • | | | |
| CytoTrack Yellow 542/556 (1351204)** | 561 | 542 | 565 | | • | | |
| PE | 488 | 496 | 578 | | • | | |
| DsRed | 488 | 554 | 586 | | • | | |
| PE–Texas Red | 488 | 496 | 613 | | • | | |
| ReadiDrop™ PI (1351101)*** | 488 | 538 | 617 | | • | | |
| CytoTrack Red 628/643 (1351205)** | 640 | 628 | 643 | | | • | |
| ReadiDrop 7-AAD (1351102)*** | 488 | 548 | 648 | | | • | |
| ReadiLink 633/655 (1351005)* | 640 | 633 | 655 | | | • | |
| VivaFix 649/660 (1351118)*** | 640 | 649 | 660 | | | • | |
| APC | 640 | 650 | 660 | | | • | |
| PE–Alexa Fluor 647 | 488 | 496 | 668 | | | • | |
| Alexa Fluor 647 | 640 | 650 | 668 | | | • | |
| PE–Cy5 | 488 | 496 | 670 | | | • | |
| ReadiLink 647/674 (1351006)* | 640 | 647 | 674 | | | • | |
| DyLight 650 | 640 | 652 | 672 | | | • | |
| PerCP | 488 | 482 | 676 | | | • | |
| PerCP–Cy5.5 | 488 | 482 | 690 | | | • | |
| PE–Cy5.5 | 488 | 496 | 690 | | | • | |
| APC–Cy5.5 | 640 | 650 | 694 | | | • | |
| ReadiLink 680/701 (1351007)* | 640 | 680 | 701 | | | | • |
| Alexa Fluor 680 | 640 | 679 | 702 | | | | • |
| DyLight 680 | 640 | 692 | 712 | | | | • |
| Alexa Fluor 700 | 640 | 696 | 719 | | | | • |
| APC–Cy7 | 640 | 650 | 774 | | | | • |
| PE–Cy7 | 488 | 496 | 785 | | | | • |

Instrument Filter Configuration

Catalog Number

Filter Set

S3e Cell Sorter (488/640 nm)

1451008

FL1: 526/48, FL2: 593/40, FL3: 670/30, FL4: 700LP

Note: This is not a comprehensive list of fluorophores and dyes that can be used with this instrument. Additional options are available for use with the 488/640 nm S3e Cell Sorter.

* Catalog number for antibody labeling kit.

** Catalog number for cell proliferation assay.

*** Catalog number for cell viability assay (live/dead).

7-AAD, 7-aminoactinomycin D; APC, allophycocyanin; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; EGFP, enhanced green fluorescent protein; Em Max, emission maximum; Ex Max, excitation maximum; EYFP, enhanced yellow fluorescent protein; FITC, fluorescein isothiocyanate; PE, phycoerythrin; PerCP, peridinin chlorophyll protein; PI, propidium iodide.

Alexa Fluor, DyLight, and Texas Red are trademarks of Thermo Fisher Scientific Inc. Cy is a trademark of GE Healthcare. ReadiLink is a trademark of AAT Bioquest, Inc.