

# S3e™ Cell Sorter (488 nm laser)

| Fluorophore or Dye                   | Laser          | Ex Max, nm              | Em Max, nm | Fluorescence Emission Color |              |
|--------------------------------------|----------------|-------------------------|------------|-----------------------------|--------------|
|                                      |                |                         |            | FL1<br>510–540              | FL2<br>540LP |
| 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        | •                           |              |
| PE                                   | 488            | 496                     | 578        |                             | •            |
| DsRed                                | 488            | 554                     | 586        |                             | •            |
| PE–Texas Red                         | 488            | 496                     | 613        |                             | •            |
| ReadiDrop™ PI (1351101)***           | 488            | 538                     | 617        |                             | •            |
| ReadiDrop 7-AAD (1351102)***         | 488            | 548                     | 648        |                             | •            |
| PE–Alexa Fluor 647                   | 488            | 496                     | 668        |                             | •            |
| PE–Cy5                               | 488            | 496                     | 670        |                             | •            |
| PerCP                                | 488            | 482                     | 676        |                             | •            |
| PerCP–Cy5.5                          | 488            | 482                     | 690        |                             | •            |
| PE–Cy5.5                             | 488            | 496                     | 690        |                             | •            |
| PE–Cy7                               | 488            | 496                     | 785        |                             | •            |
| Instrument Filter Configuration      | Catalog Number | Filter Set              |            |                             |              |
| S3e Cell Sorter (488 nm)             | 1451005        | FL1: 525/30, FL2: 540LP |            |                             |              |

**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 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; 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.