

S3e™ Cell Sorter (488/561 nm lasers)

Fluorophore or Dye	Laser	Ex Max, nm	Em Max, nm	Fluorescence Emission Color			
				FL1 510–540	FL2 573–598	FL3 602–627	FL4 655LP
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	•			
ReadiLink 555/570 (1351003)*	561	555	570		•		
VivaFix 547/573 (1351116)***	561	547	573		•		
DyLight 550	561	562	576		•		
PE	488, 561	496, 565	578		•		
tdTomato	561	554	581		•		
DsRed	488, 561	554	586		•		
VivaFix 583/603 (1351117)***	561	583	603			•	
mCherry	561	587	610			•	
ReadiLink 594/610 (1351004)*	561	594	610			•	
PE–Texas Red	488, 561	496, 565	613			•	
ReadiDrop™ PI (1351101)***	488, 561	538	617			•	
ReadiDrop 7-AAD (1351102)***	488, 561	548	648				•
PE–Alexa Fluor 647	488, 561	496, 565	668				•
PE–Cy5	488, 561	496, 565	670				•
PerCP	488	482	676				•
PerCP–Cy5.5	488	482	690				•
PE–Cy5.5	488, 561	496, 565	690				•
PE–Cy7	488, 561	496, 565	785				•

Instrument Filter Configuration

Catalog Number

Filter Set

S3e Cell Sorter (488/561 nm)

1451006

FL1: 525/30, FL2: 586/25, FL3: 615/25, FL4: 655 LP

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 488/561 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.