

# S3e™ Cell Sorter (405/488/561 nm lasers)

Fluorophore or Dye	Laser	Ex Max, nm	Em Max, nm	Fluorescence Emission Color			
				FL1 417–477	FL2 510–540	FL3 573–598	FL4 600LP
DyLight 405	405	400	420	•			
Alexa Fluor 405	405	401	421	•			
BV421	405	407	421	•			
Horizon V450	405	404	448	•			
VivaFix™ 410/450 (1351112)*	405	410	450	•			
Pacific Blue	405	401	452	•			
CytoTrack™ Blue 403/454 (1351202)**	405	403	454	•			
ReadiLink 405/454 (1351011)***	405	403	454	•			
DAPI	405	359	461	•			
ECFP	405	434	477	•			
EGFP	488	488	507		•		
VivaFix 408/512 (1351113)*	405	408	512		•		
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		•		
VivaFix 398/550 (1351114)*	405	398	550			•	
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 (405/488/561 nm)	12007058	FL1: 447/60, FL2: 525/30, FL3: 586/25, FL4: 600LP					

**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 405/488/561 nm S3e Cell Sorter.

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

\*\* Catalog number for cell proliferation assay.

\*\*\* Catalog number for antibody labeling kit.

7-AAD, 7-aminoactinomycin D; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; DAPI, 4',6'-diamidino-2-phenylindole dihydrochloride; ECFP, enhanced cyan fluorescent protein; 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, Pacific Blue, and Texas Red are trademarks of Thermo Fisher Scientific Inc. BD Horizon and BD Horizon Brilliant are trademarks of Becton, Dickinson and Company (BD Horizon V450 and BD Horizon Brilliant Violet 421 are referred to as Horizon V450 and BV421, respectively). Cy is a trademark of GE Healthcare. ReadiLink is a trademark of AAT Bioquest, Inc.

