

S3e™ Cell Sorter (405/488/640 nm lasers)

Fluorophore or Dye	Laser	Ex Max, nm	Em Max, nm	Fluorescence Emission Color			
				FL1 417–477	FL2 502–550	FL3 573–613	FL4 650LP
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		•		
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				•
DyLight 650	640	652	672				•
ReadiLink 647/674 (1351006)***	640	647	674				•
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	785				•
PE–Cy7	488	496	785				•

Instrument Filter Configuration

Catalog Number

Filter Set

S3e Cell Sorter (405/488/640 nm)

12007059

FL1: 447/60, FL2: 526/48, FL3: 593/40, FL4: 650LP

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 405/488/640 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; APC, allophycocyanin; 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. ReadLink is a trademark of AAT Bioquest, Inc.

