

# S3e™ Cell Sorter Laser Configuration Guide

Fluorophore or Dye	Laser	Ex Max, nm	Em Max, nm	Fluorescence Emission Color						Fluorophore or Dye
				405/488/561	405/488/640	488/561/640	488/561	488/640	488	
DyLight 405	405	400	420	FL1	FL1					DyLight 405
Alexa Fluor 405	405	401	421	FL1	FL1					Alexa Fluor 405
BV421	405	407	421	FL1	FL1					BV421
Horizon V450	405	404	448	FL1	FL1					Horizon V450
Pacific Blue	405	401	452	FL1	FL1					Pacific Blue
DAPI	405	359	461	FL1	FL1					DAPI
ECFP	405	434	477	FL1	FL1					ECFP
EGFP	488	488	507	FL2	FL2	FL1	FL1	FL1	FL1	EGFP
CFDA-SE (1351201)*	488	492	517	FL2	FL2	FL1	FL1	FL1	FL1	CFDA-SE (1351201)*
DyLight 488	488	493	518	FL2	FL2	FL1	FL1	FL1	FL1	DyLight 488
FITC	488	495	519	FL2	FL2	FL1	FL1	FL1	FL1	FITC
Alexa Fluor 488	488	499	519	FL2	FL2	FL1	FL1	FL1	FL1	Alexa Fluor 488
EYFP	488	514	527	FL2	FL2	FL1	FL1	FL1	FL1	EYFP
DyLight 550	561	562	576	FL3		FL2	FL2			DyLight 550
PE	488, 561	496, 565	578	FL3	FL3	FL2	FL2	FL2	FL2	PE
tdTomato	561	554	581	FL3		FL2	FL2			tdTomato
DsRed	488, 561	554	586	FL3	FL3	FL2	FL2	FL2	FL2	DsRed
mCherry	561	587	610	FL4		FL3	FL3			mCherry
PE-Texas Red	488, 561	496, 565	613	FL4	FL3	FL3	FL3	FL2	FL2	PE-Texas Red
PI (1351101)*	488, 561	538	617	FL4	FL3	FL3	FL3	FL2	FL2	PI (1351101)*
7-AAD (1351102)*	488, 561	548	648	FL4	FL4	FL4	FL4	FL3	FL2	7-AAD (1351102)*
APC	640	650	660		FL4	FL4		FL3		APC
PE-Alexa Fluor 647	488, 561	496, 565	668	FL4	FL4	FL4	FL4	FL3	FL2	PE-Alexa Fluor 647
Alexa Fluor 647	640	650	668		FL4	FL4		FL3		Alexa Fluor 647
PE-Cy5	488, 561	496, 565	670	FL4	FL4	FL4	FL4	FL3	FL2	PE-Cy5
DyLight 650	640	652	672		FL4	FL4		FL3		DyLight 650
PerCP	488	482	676	FL4	FL4	FL4	FL4	FL3	FL2	PerCP
PerCP-Cy5.5	488	482	690	FL4	FL4	FL4	FL4	FL3	FL2	PerCP-Cy5.5
PE-Cy5.5	488, 561	496, 565	690	FL4	FL4	FL4	FL4	FL3	FL2	PE-Cy5.5
APC-Cy5.5	640	650	694		FL4	FL4		FL3		APC-Cy5.5
Alexa Fluor 680	640	679	702		FL4	FL4		FL4		Alexa Fluor 680
DyLight 680	640	692	712		FL4	FL4		FL4		DyLight 680
Alexa Fluor 700	640	696	719		FL4	FL4		FL4		Alexa Fluor 700
APC-Cy7	640	650	774		FL4	FL4		FL4		APC-Cy7
PE-Cy7	488, 561	496, 565	785	FL4	FL4	FL4	FL4	FL4	FL2	PE-Cy7

\* Indicates catalog number.

**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 S3e Cell Sorter.



# S3e Cell Sorter Laser Configuration Guide

## Instrument and Reagent Ordering Information

Instrument Filter Configuration	Catalog Number	Filter Set			
		FL1	FL2	FL3	FL4
S3e Cell Sorter (488 nm)	1451005	525/30 (510–540)	540 LP		
S3e Cell Sorter (488/561 nm)	1451006	525/30 (510–540)	586/25 (573–598)	615/25 (602–627)	655LP
S3e Cell Sorter (488/640 nm)	1451008	526/48 (502–550)	593/40 (573–613)	670/30 (655–685)	700LP
S3e Cell Sorter (405/488/561 nm)	12007058	447/60 (417–477)	525/30 (510–540)	586/25 (573–598)	600LP
S3e Cell Sorter (405/488/640 nm)	12007059	447/60 (417–477)	526/48 (502–550)	593/40 (573–613)	650LP
S3e Cell Sorter (488/561/640 nm)	12007052	525/30 (510–540)	586/25 (573–598)	615/25 (602–627)	655LP

Fluorophore or Dye	Catalog Number	Laser	Ex Max, nm	Em Max, nm	Fluorescence Emission Color						Fluorophore or Dye
					405/488/561	405/488/640	488/561/640	488/561	488/640	488	
CFDA-SE	1351201	488	492	517	FL2	FL2	FL1	FL1	FL1	FL1	CFDA-SE
ReadiDrop™ PI	1351101	488, 561	538	617	FL4	FL3	FL3	FL3	FL2	FL2	ReadiDrop™ PI
ReadiDrop 7-AAD	1351102	488, 561	548	648	FL4	FL4	FL4	FL4	FL4	FL4	ReadiDrop 7-AAD
<b>Antibody Labeling Kits</b>											
ReadiLink 405/454	1351011	405	403	454	FL1	FL1					ReadiLink 405/454
ReadiLink 492/516	1351002	488	492	516	FL2	FL2	FL1	FL1	FL1	FL1	ReadiLink 492/516
ReadiLink 555/570	1351003	561	555	570	FL3		FL2	FL2			ReadiLink 555/570
ReadiLink 594/610	1351004	561	594	610	FL4		FL3	FL3			ReadiLink 594/610
ReadiLink 633/655	1351005	640	633	655		FL4	FL4		FL4		ReadiLink 633/655
ReadiLink 647/674	1351006	640	647	674		FL4	FL4		FL4		ReadiLink 647/674
ReadiLink 680/701	1351007	640	680	701		FL4	FL4		FL4		ReadiLink 680/701
<b>Cell Viability Dyes</b>											
VivaFix™ 410/450	1351112	405	410	450	FL1	FL1					VivaFix™ 410/450
VivaFix 408/512	1351113	405	408	512	FL2	FL2					VivaFix 408/512
VivaFix 498/521	1351115	488	498	521	FL2	FL2	FL1	FL1	FL1	FL1	VivaFix 498/521
VivaFix 398/550	1351114	405	398	550	FL3	FL3					VivaFix 398/550
VivaFix 547/573	1351116	561	547	573	FL3		FL2	FL2			VivaFix 547/573
VivaFix 583/603	1351117	561	583	603	FL4		FL3	FL3			VivaFix 583/603
VivaFix 649/660	1351118	640	649	660		FL4	FL4		FL4		VivaFix 649/660
<b>Cell Proliferation Dyes</b>											
CytoTrack™ Blue 403/454	1351202	405	403	454	FL1	FL1					CytoTrack™ Blue 403/454
CytoTrack Green 511/525	1351203	488	511	525	FL2	FL2	FL1	FL1	FL1	FL1	CytoTrack Green 511/525
CytoTrack Red 628/643	1351205	640	628	643		FL4	FL4		FL4		CytoTrack Red 628/643

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 Texas Red 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.

