

S3e™ Cell Sorter (488 nm)

Laser	Filter	Fluorophore or Dye	Excitation Maximum, nm	Emission Maximum, nm
488	FL1 525/30	EGFP	488	507
		ReadiLink 492/516 (1351002)*	492	516
		CFDA-SE (1351201)*	492	517
		DyLight 488	493	518
		FITC	495	519
		Alexa Fluor 488	499	519
		SYTOX Green	504	523
		CytoTrack™ Green 511/525 (1351203)*	511	525
		EYFP	514	527
		488	FL2 540/LP	PE
DsRed	554			586
PE-CF594	496			612
PI	536			617
ReadiDrop™ PI (1351101)*	538			617
PE-Texas Red	496			619
PE-Alexa Fluor 610	496			626
ReadiDrop 7-AAD (1351102)*	548			648
7-AAD	548			648
PE-Alexa Fluor 647	496			669
PE-Cy5	496			670
PE-Cy5.5	490			693
PerCP-Cy5.5	482			694
PE-Cy7	496			767

* 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 488 nm S3e Cell Sorter.

7-AAD, 7-aminoactinomycin D; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; EGFP, enhanced green fluorescent protein; EYFP, enhanced yellow fluorescent protein; FITC, fluorescein isothiocyanate; PE, phycoerythrin; PerCP, peridinin chlorophyll protein; PI, propidium iodide.

Alexa Fluor, SYTOX, and Texas Red are trademarks of Life Technologies Corporation. Cy is a trademark of GE Healthcare. DyLight is a trademark of Thermo Fisher Scientific Inc. ReadLink is a trademark of AAT Bioquest, Inc.

S3e™ Cell Sorter (488/561 nm)

Laser	Filter	Fluorophore or Dye	Excitation Maximum, nm	Emission Maximum, nm
488	FL1 525/30	EGFP	488	507
		ReadiLink 492/516 (1351002)*	492	516
		CFDA-SE (1351201)*	492	517
		DyLight 488	493	518
		FITC	495	519
		Alexa Fluor 488	499	519
		SYTOX Green	504	523
		CytoTrack™ Green 511/525 (1351203)*	511	525
		EYFP	514	527
488	FL2 586/25	PE	496, 565	578
		RFP	554	584
488	FL3 615/25	PE-CF594	496, 565	612
		PI	536	617
		ReadiDrop™ PI (1351101)*	538	617
		PE-Texas Red	496, 565	619
488	FL4 655/LP	ReadiDrop 7-AAD (1351102)*	548	648
		7-AAD	548	648
		PE-Cy5	496, 565	670
		PE-Cy5.5	490, 565	693
		PerCP-Cy5.5	490	695
		PE-Cy7	496, 565	785
561	FL2 586/25	SYTOX Orange	540	570
		ReadiLink 555/570 (1351003)*	555	570
		PE	496, 565	578
		tdTomato	554	581
		RFP	554	584
		DsRed	554	586
561	FL3 615/25	mCherry	587	610
		ReadiLink 594/610 (1351004)*	594	610
		PE-CF594	496, 565	612
		PE-Texas Red	496, 565	613
		PI	536	617
		ReadiDrop PI (1351101)*	538	617
		DyLight 594	593	618
		PE-Texas Red	496, 565	619
561	FL4 655/LP	ReadiDrop 7-AAD (1351102)*	548	648
		7-AAD	548	648
		PE-Alexa Fluor 647	496, 565	669
		PE-Cy5	496, 565	670
		PE-Cy5.5	490, 565	693
		PE-Cy7	496, 565	785

* 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 488/561 nm S3e Cell Sorter.

7-AAD, 7-aminoactinomycin D; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; EGFP, enhanced green fluorescent protein; EYFP, enhanced yellow fluorescent protein; FITC, fluorescein isothiocyanate; PE, phycoerythrin; PerCP, peridinin chlorophyll protein; PI, propidium iodide; RFP, red fluorescent protein.

Alexa Fluor, SYTOX, and Texas Red are trademarks of Life Technologies Corporation. Cy is a trademark of GE Healthcare. DyLight is a trademark of Thermo Fisher Scientific Inc. ReadiLink is a trademark of AAT Bioquest, Inc.

S3e™ Cell Sorter (488/640 nm)

Laser	Filter	Fluorophore or Dye	Excitation Maximum, nm	Emission Maximum, nm
488	FL1 526/48	EGFP	488	507
		ReadiLink 492/516 (1351002)*	492	516
		CFDA-SE (1351201)*	492	517
		DyLight 488	493	518
		FITC	495	519
		Alexa Fluor 488	499	519
		SYTOX Green	504	523
		CytoTrack™ Green 511/525 (1351203)*	511	525
		EYFP	514	527
488	FL2 593/40	CytoTrack Yellow 542/556 (1351204)*	542	556
		PE	496	578
		RFP	554	584
		DsRed	554	586
		PE-CF594	496	612
		PI	536	617
		ReadiDrop™ PI (1351101)*	538	617
		PE-Texas Red	496	619
		PE-Alexa Fluor 610	496	626
488	FL3 670/30	ReadiDrop 7-AAD (1351102)*	548	648
		7-AAD	548	648
		PE-Alexa Fluor 647	496, 653	669
		PE-Cy5	496, 650	670
		PE-Cy5.5	490	693
		PerCP-Cy5.5	490	695
488	FL4 700/LP	PE-Cy7	496	785
640	FL3 670/30	CytoTrack Red 628/643 (1351205)*	628	643
		APC	650	660
		ReadiLink 633/655 (1351005)*	633	655
		DyLight 633	638	658
		SYTOX Red	640	658
		Alexa Fluor 647	650	668
ReadiLink 647/674 (1351006)*	647	674		
640	FL4 700/LP	APC-Cy5.5	650	694
		ReadiLink 680/701 (1351007)*	680	701
		Alexa Fluor 680	679	702
		DyLight 680	692	712
		ReadiLink 700/713 (1351008)*	700	713
		Alexa Fluor 700	696	719
APC-Cy7	650	774		

* 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 488/640 nm S3e Cell Sorter.

7-AAD, 7-aminoactinomycin D; APC, allophycocyanin; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; EGFP, enhanced green fluorescent protein; EYFP, enhanced yellow fluorescent protein; FITC, fluorescein isothiocyanate; PE, phycoerythrin; PerCP, peridinin chlorophyll protein; PI, propidium iodide; RFP, red fluorescent protein.

Alexa Fluor, SYTOX, and Texas Red are trademarks of Life Technologies Corporation. Cy is a trademark of GE Healthcare. DyLight is a trademark of Thermo Fisher Scientific Inc. Readilink is a trademark of AAT Bioquest, Inc.