

ZE5 Cell Analyzer Reagent Reference Guide

5-Laser Options	ZE5 Emission Filters			Fluorescent Antibody Conjugates	Antibody Labeling	Viability Dyes (Fixable)	Viability Dyes (Nonfixable)	DNA Content/ Cell Cycle	Cell Proliferation	Apoptosis/Autophagy
	Standard 5 Laser	5 Laser UV Option A	5 Laser UV Option B							
Ultraviolet (355 nm)	387/11	387/11	387/11	SBUV400						
	447/60		460/22	SBUV445		VivaFix 353/442	PureBlu DAPI	PureBlu DAPI PureBlu Hoechst 33342		
	525/50	509/24	509/24	SBUV510			PureBlu DAPI	PureBlu DAPI PureBlu Hoechst 33342		
		577/15	577/15	SBUV575						
		615/24		SBUV605						
	670/30	670/30	670/30	SBUV665			ReadiDrop PI	ReadiDrop PI PureBlu Hoechst 33342		ReadiDrop PI
	700LP	747/33 780LP	747/33 780LP	SBUV740 SBUV795						
Violet (405 nm)	420/10	420/10	420/10	DyLight 405						
	460/22	460/22	460/22	SBV440 SBV475 Pacific Blue	ReadiLink 405/454	VivaFix 410/450	PureBlu DAPI	PureBlu DAPI	CytoTrack Blue 403/454	
	525/50	525/50	525/50	SBV515 Amethyst Orange		VivaFix 408/512				
	615/24	615/24	615/24	SBV570 SBV610						
	670/30	670/30	670/30	SBV670						
	720/60	720/60	720/60	SBV710						
	750LP	750LP	750LP	SBV760 SBV790						
Blue (488 nm)	525/25	525/35	525/35	FITC Alexa Fluor 488 DyLight 488	ReadiLink 492/516 LYNX Rapid Plus DyLight 488 LYNX Rapid Fluorescein	VivaFix 498/521			CytoTrack Green 511/525 CFDA-SE	FAM FLICA Poly Caspase FAM FLICA Caspase-1 FAM FLICA Caspase-2 FAM FLICA Caspase-3/7 FAM FLICA Caspase-6 FAM FLICA Caspase-8 FAM FLICA Caspase-9 FAM FLICA Caspase-10 MitoPT JC-1 (apoptotic) Cathepsin B
	593/52	593/52	593/52	SBB580 SBB615 PE	LYNX Rapid RPE LYNX Rapid Plus Rhodamine LYNX Rapid RPE-Texas Red		ReadiDrop PI	ReadiDrop PI		Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) ReadiDrop PI
	692/80	692/80	692/80	SBB675 SBB700 PE-Cy5 PE-Cy5.5 PE-Alexa Fluor 647	LYNX Rapid PerCP LYNX Rapid PerCP-Cy5.5 LYNX Rapid RPE-Cy5 LYNX Rapid RPE-Cy5.5		ReadiDrop 7-AAD	ReadiDrop 7-AAD		
	750LP	750LP	750LP	SBB765 SBB810 PE-Cy7 PE-Alexa Fluor 750	LYNX Rapid RPE-Cy7					
Yellow/Green (561 nm)	577/15			SBY575 PE DyLight 550	ReadiLink 555/570 LYNX Rapid Plus Cy3 LYNX Rapid Plus DyLight 550 LYNX Rapid RPE LYNX Rapid Plus Rhodamine	VivaFix 547/573				Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) SR FLICA Poly Caspase SR FLICA Caspase-9
	589/15	583/30	583/30	SBY575 PE DyLight 550	ReadiLink 555/570 LYNX Rapid Plus Cy3 LYNX Rapid Plus DyLight 550 LYNX Rapid RPE LYNX Rapid Plus Rhodamine	VivaFix 547/573				Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) SR FLICA Poly Caspase SR FLICA Caspase-9
	615/24	615/24	615/24	SBY605 Texas Red	ReadiLink 594/610 LYNX Rapid RPE-Texas Red	VivaFix 583/603	ReadiDrop PI	ReadiDrop PI		ReadiDrop PI Autophagy Assay Kit Red
	640/20						ReadiDrop 7-AAD	ReadiDrop 7-AAD		
	670/30	670/30	670/30	SBY665 PE-Cy5 PE-Alexa Fluor 647	LYNX Rapid RPE-Cy5		ReadiDrop 7-AAD	ReadiDrop 7-AAD		
	720/60	720/60	720/60	SBY720 PE-Cy5.5	LYNX Rapid RPE-Cy5.5					
	750LP	750LP	750LP	SBY800 PE-Cy7 PE-Alexa Fluor 750	LYNX Rapid RPE-Cy7					
Red (640 nm)	670/30	670/30	670/30	SBR670 APC Alexa Fluor 647 Cy5 DyLight 650	ReadiLink 633/655 ReadiLink 647/674 LYNX Rapid Plus APC LYNX Rapid Plus Cy5 LYNX Rapid DyLight 650	VivaFix 649/660			CytoTrack Red 628/643	Annexin V-APC
	720/60	720/60	720/60	SBR715 Alexa Fluor 700 DyLight 680	ReadiLink 700/713 LYNX Rapid Plus APC-Cy5.5					Pyroptosis 660 Caspase-1 FLICA 660 Caspase-1 FLICA 660 Caspase-3/7 FLICA 660 Poly Caspase
	775/50	775/50	775/50	SBR775 DyLight 755	LYNX Rapid APC-Cy7					
	800LP	800LP	800LP	SBR815						

These reagents are provided as general information. End users must check compatibility with their systems. 7-AAD, 7-aminoactinomycin D; APC, allophycocyanin; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; Cy, cyanine; DAPI, 4',6-diamidino-2-phenylindole dihydrochloride; FITC, fluorescein isothiocyanate; FLICA, fluoro-chrome-labeled inhibitors of caspases; JC-1, tetraethylbenzimidazolylcarbocyanine iodide; LP, long pass; PE, phycoerythrin; PerCP, peridinin chlorophyll; PI, propidium iodide; RPE, R-phycoerythrin; SBB, StarBright™ Blue; SBR, StarBright Red; SBUV, StarBright UltraViolet; SBV, StarBright Violet; SBY, StarBright Yellow; TMRE, tetramethylrhodamine ethyl ester; TMRM, tetramethylrhodamine methyl ester; UV, ultraviolet.

ZE5 Cell Analyzer Reagent Reference Guide

4-Laser Options	ZE5 Emission Filters	Fluorescent Antibody Conjugates	Antibody Labeling	Viability Dyes (Fixable)	Viability Dyes (Nonfixable)	DNA Content/ Cell Cycle	Cell Proliferation	Apoptosis/Autophagy
Violet (405 nm)	420/10	DyLight 405						
	460/22	SBV440 SBV475 Pacific Blue	ReadiLink 405/454	VivaFix 410/450	PureBlu DAPI	PureBlu DAPI	CytoTrack Blue 403/454	
	525/50	SBV515 Amethyst Orange		VivaFix 408/512				
	615/24	SBV570 SBV610						
	670/30	SBV670						
	720/60	SBV710						
	750LP	SBV760 SBV790						
Blue (488 nm)	509/24	FITC Alexa Fluor 488 DyLight 488	ReadiLink 492/516 LYNX Rapid Plus DyLight 488 LYNX Rapid Fluorescein	VivaFix 498/521			CytoTrack Green 511/525 CFDA-SE	FAM FLICA Poly Caspase FAM FLICA Caspase-1 FAM FLICA Caspase-2 FAM FLICA Caspase-3/7 FAM FLICA Caspase-6 FAM FLICA Caspase-8 FAM FLICA Caspase-9 FAM FLICA Caspase-10 MitoPT JC-1 (apoptotic) Cathepsin B
	549/15							
	583/30	SBB580 PE	LYNX Rapid RPE LYNX Rapid Plus Rhodamine		ReadiDrop PI	ReadiDrop PI		Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) ReadiDrop PI
	615/24	SBB615	LYNX Rapid RPE-Texas Red					
	692/80	SBB675 SBB700 PE-Cy5 PE-Cy5.5 PE-Alexa Fluor 647	LYNX Rapid PerCP LYNX Rapid PerCP-Cy5.5 LYNX Rapid RPE-Cy5 LYNX Rapid RPE-Cy5.5		ReadiDrop 7-AAD	ReadiDrop 7-AAD		
750LP	SBB765 SBB810 PE-Cy7 PE-Alexa Fluor 750	LYNX Rapid RPE-Cy7						
Yellow/Green (561 nm)	577/15	SBY575 PE DyLight 550	ReadiLink 555/570 LYNX Rapid Plus Cy3 LYNX Rapid Plus DyLight 550 LYNX Rapid RPE LYNX Rapid Plus Rhodamine	VivaFix 547/573				Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) SR FLICA Poly Caspase SR FLICA Caspase-9
	589/15	SBY575 PE DyLight 550	ReadiLink 555/570 LYNX Rapid Plus Cy3 LYNX Rapid Plus DyLight 550 LYNX Rapid RPE LYNX Rapid Plus Rhodamine	VivaFix 547/573				Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) SR FLICA Poly Caspase SR FLICA Caspase-9
	615/24	SBY605 Texas Red	ReadiLink 594/610 LYNX Rapid RPE-Texas Red	VivaFix 583/603	ReadiDrop PI	ReadiDrop PI		ReadiDrop PI Autophagy Assay Kit Red
	640/20				ReadiDrop 7-AAD	ReadiDrop 7-AAD		
	670/30	SBY665 PE-Cy5 PE-Alexa Fluor 647	LYNX Rapid RPE-Cy5					
	720/60	SBY720 PE-Cy5.5	LYNX Rapid RPE-Cy5.5					
	750LP	SBY800 PE-Cy7 PE-Alexa Fluor 750	LYNX Rapid RPE-Cy7					
Red (640 nm)	670/30	SBR670 APC Alexa Fluor 647 Cy5 DyLight 650	ReadiLink 633/655 ReadiLink 647/674 LYNX Rapid Plus APC LYNX Rapid Plus Cy5 LYNX Rapid DyLight 650	VivaFix 649/660			CytoTrack Red 628/643	Annexin V-APC
	720/60	SBR715 Alexa Fluor 700 DyLight 680	ReadiLink 700/713 LYNX Rapid Plus APC-Cy5.5					Pyroptosis 660 Caspase-1 FLICA 660 Caspase-1 FLICA 660 Caspase-3/7 FLICA 660 Poly Caspase
	775/50	SBR775 DyLight 755	LYNX Rapid APC-Cy7					
	800LP	SBR815						

These reagents are provided as general information. End users must check compatibility with their systems.
 7-AAD, 7-aminoactinomycin D; APC, allophycocyanin; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; Cy, cyanine; DAPI, 4',6-diamidino-2-phenylindole dihydrochloride; FITC, fluorescein isothiocyanate; FLICA, fluorochrome-labeled inhibitors of caspases; JC-1, tetraethylbenzimidazolylcarbocyanine iodide; LP, long pass; PE, phycoerythrin; PerCP, peridinin chlorophyll; PI, propidium iodide; RPE, R-phycoerythrin; SBB, StarBright Blue; SBR, StarBright Red; SBV, StarBright Violet; SBY, StarBright Yellow; TMRE, tetramethylrhodamine ethyl ester; TMRM, tetramethylrhodamine methyl ester.

ZE5 Cell Analyzer Reagent Reference Guide

3-Laser Options	ZE5 Emission Filters			Fluorescent Antibody Conjugates	Antibody Labeling	Viability Dyes (Fixable)	Viability Dyes (Nonfixable)	DNA Content/ Cell Cycle	Cell Proliferation	Apoptosis/Autophagy
	405 nm 488 nm 640 nm	488 nm 561 nm 640 nm	405 nm 488 nm 561 nm							
Violet (405 nm)	420/10		420/10	DyLight 405						
	460/22		460/22	SBV440 SBV475 Pacific Blue	ReadiLink 405/454	VivaFix 410/450	PureBlu DAPI	PureBlu DAPI	CytoTrack Blue 403/454	
	525/50		525/50	SBV515 Amethyst Orange		VivaFix 408/512				
	615/24		615/24	SBV570 SBV610						
	670/30		670/30	SBV670						
	720/60		720/60	SBV710						
	750LP		750LP	SBV760 SBV790						
Blue (488 nm)	509/24	509/24	509/24	FITC Alexa Fluor 488 DyLight 488	ReadiLink 492/516 LYNX Rapid Plus DyLight 488 LYNX Rapid Fluorescein	VivaFix 498/521			CytoTrack Green 511/525 CFDA-SE	FAM FLICA Poly Caspase FAM FLICA Caspase-1 FAM FLICA Caspase-2 FAM FLICA Caspase-3/7 FAM FLICA Caspase-6 FAM FLICA Caspase-8 FAM FLICA Caspase-9 FAM FLICA Caspase-10 MitoPT JC-1 (apoptotic) Cathepsin B
	549/15	549/15	549/15							
	583/30	583/30	583/30	SBB580 PE	LYNX Rapid RPE LYNX Rapid Plus Rhodamine		ReadiDrop PI	ReadiDrop PI		Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) ReadiDrop PI
	615/24	615/24	615/24	SBB615	LYNX Rapid RPE-Texas Red					
	692/80	692/80	692/80	SBB675 SBB700 PE-Cy5 PE-Cy5.5 PE-Alexa Fluor 647	LYNX Rapid PerCP LYNX Rapid PerCP-Cy5.5 LYNX Rapid RPE-Cy5 LYNX Rapid RPE-Cy5.5		ReadiDrop 7-AAD	ReadiDrop 7-AAD		
	750LP	750LP	750LP	SBB765 SBB810 PE-Cy7 PE-Alexa Fluor 750	LYNX Rapid RPE-Cy7					
	Yellow/Green (561 nm)		577/15	577/15	SBY575 PE DyLight 550	ReadiLink 555/570 LYNX Rapid Plus Cy3 LYNX Rapid Plus DyLight 550 LYNX Rapid RPE LYNX Rapid Plus Rhodamine	VivaFix 547/573			
		589/15	589/15	SBY575 PE DyLight 550	ReadiLink 555/570 LYNX Rapid Plus Cy3 LYNX Rapid Plus DyLight 550 LYNX Rapid RPE LYNX Rapid Plus Rhodamine	VivaFix 547/573				Annexin V-RPE MitoPT TMRE MitoPT TMRM MitoPT JC-1 (nonapoptotic) SR FLICA Poly Caspase SR FLICA Caspase-9
		615/24	615/24	SBY605 Texas Red	ReadiLink 594/610 LYNX Rapid RPE-Texas Red	VivaFix 583/603	ReadiDrop PI	ReadiDrop PI		ReadiDrop PI Autophagy Assay Kit Red
		640/20	640/20				ReadiDrop 7-AAD	ReadiDrop 7-AAD		
		670/30	670/30	SBY665 PE-Cy5 PE-Alexa Fluor 647	LYNX Rapid RPE-Cy5					
		720/60	720/60	SBY720 PE-Cy5.5	LYNX Rapid RPE-Cy5.5					
		750/LP	750/LP	SBY800 PE-Cy7 PE-Alexa Fluor 750	LYNX Rapid RPE-Cy7					
Red (640 nm)	670/30	670/30		SBR670 APC Alexa Fluor 647 Cy5 DyLight 650	ReadiLink 633/655 ReadiLink 647/674 LYNX Rapid Plus APC LYNX Rapid Plus Cy5 LYNX Rapid DyLight 650	VivaFix 649/660			CytoTrack Red 628/643	Annexin V-APC
	720/60	720/60		SBR715 Alexa Fluor 700 DyLight 680	ReadiLink 700/713 LYNX Rapid Plus APC-Cy5.5					Pyroptosis 660 Caspase-1 FLICA 660 Caspase-1 FLICA 660 Caspase-3/7 FLICA 660 Poly Caspase
	775/50	775/50		SBR775 DyLight 755	LYNX Rapid APC-Cy7					
	800LP	800LP		SBR815						

These reagents are provided as general information. End users must check compatibility with their systems.
 7-AAD, 7-aminoactinomycin D; APC, allophycocyanin; CFDA-SE, 5(6)-carboxyfluorescein diacetate succinimidyl ester; Cy, cyanine; DAPI, 4',6-diamidino-2-phenylindole dihydrochloride; FITC, fluorescein isothiocyanate; FLICA, fluoro-chrome-labeled inhibitors of caspases; JC-1, tetraethylbenzimidazolylcarbocyanine iodide; LP, long pass; PE, phycoerythrin; PerCP, peridinin chlorophyll; PI, propidium iodide; RPE, R-phycoerythrin; SBB, StarBright Blue; SBR, StarBright Red; SBV, StarBright Violet; SBY, StarBright Yellow; TMRE, tetramethylrhodamine ethyl ester; TMRM, tetramethylrhodamine methyl ester.

BIO-RAD and STARBRIGHT are trademarks of Bio-Rad Laboratories, Inc. in certain jurisdictions. All trademarks used herein are the property of their respective owner. © 2023 Bio-Rad Laboratories, Inc.

Visit bio-rad.com/FlowCytometry and bio-rad-antibodies.com/FlowCytometry for more information.

