

Flow Cytometer Buyer's Guide

From Instrument to Configuration



What to Consider When Buying a Flow Cytometer



A flow cytometer is a significant investment for your laboratory and there are many considerations to ensure that you choose an appropriate instrument to meet your users' needs. Things like the number of parameters it can measure, the throughput requirements, and sample handling will all need to be factored into your decision. This Buyer's Guide highlights some key considerations to enable you to narrow down your instrument and configuration options.

Define Your Needs

Having a clear idea of what your users' needs are will pay huge dividends. Not only will it guarantee that you approach your purchase in a practical way, but it will help you justify your budget for new equipment. If you are in a multiuser environment, consider identifying gaps in the current service by surveying your userbase. As a flow cytometer is a substantial investment, make sure you think long-term; anticipating your future needs is almost as important as the here and now.

The Bio-Rad ZE5 Cell Analyzer is designed with flexibility in mind to help adapt to future use. Its integrated plate loader can handle 1.5 and 5 ml tubes as well as plates, and it can be upgraded with the addition of extra lasers and detectors if your needs change. It is automation-ready and ideal for high-throughput screening and routine laboratory applications.

Prioritize Your Applications

Anticipate the different types of applications that your flow cytometer will be routinely used for so you can identify essential features and prioritize laser combinations. For example, if immunophenotyping is a priority, the inclusion of violet (405 nm) and ultraviolet (UV; 355 nm) lasers will help expand the number of compatible fluorophores and enables more flexible panel design.

The ZE5 Cell Analyzer offers a host of features that ensure it is suitable for a wide range of applications, including immunophenotyping, exosome detection, fluorescence resonance energy transfer (FRET), bead arrays, and high-throughput screening. The system can be preconfigured with different laser combinations based on your priority applications (Table 1).

Table 1. Laser priority by application. Larger circles represent higher priority.

Application	Laser					Explanation
	355 nm*	405 nm	488 nm	561 nm	640 nm	
Immunophenotyping	●	●	Included	●	●	UV* and violet lasers provide the highest number of colors.
Small particles**	●	●	Included	●	●	The ZE5 Cell Analyzer can be equipped with a small particle detector triggered off the violet laser. The brightness of phycoerythrin (PE), which is maximally excited by the yellow (561 nm) laser, is also useful for small particle detection.
FRET	●	●	Included	●	●	The yellow and violet lasers give the best choice of compatible FRET pairs.
Novel markers	●	●	Included	●	●	Generally, markers with only a few conjugates will be available in only fluorescein isothiocyanate (FITC), PE, and allophycocyanin (APC). The blue laser (488 nm) covers FITC and PE and the red laser APC.
Bead assays	●	●	Included	●	●	The red laser is often needed for bead arrays and the violet laser gives maximum flexibility for other applications.

* Only available for 4 and 5 laser configurations.

** Also consider the small particle detector.

APC, allophycocyanin; FITC, fluorescein isothiocyanate; FRET, fluorescence resonance energy transfer; PE, phycoerythrin; UV, ultraviolet.

The small particle detector is an additional scatter detector triggered from the violet laser and calibrated to give enhanced resolution of small particles such as bacteria, yeast, and extracellular vesicles. It is included as standard on all five-laser ZE5 Cell Analyzer configurations. Four- and three-laser configurations that have a violet laser are also available with a small particle detector. It is also an option as a field upgrade for four- and three-laser systems equipped with a violet laser.

Count Your Parameters

How many parameters do you need to measure? The number will determine which laser configuration is best for you:

- For 17 or fewer, consider a three-laser configuration
- For 17–24, consider a four-laser configuration
- For 24–27, consider a five-laser configuration that includes a small particle detector

For complex immunophenotyping assays, think about expanding the number of parameters available off the UV laser from five to seven. Note that, in this configuration, yellow laser parameters are reduced from seven to five. Refer to the ZE5 Cell Analyzer Laser and Filter Configuration Guide on pages 6–7 when configuring your ZE5 Cell Analyzer.

Think About Throughput

Ensure that your instrument can handle the volume of samples that you need to process for your current workload, but also factor in a 50% increase in throughput to ensure it can also meet your needs if they evolve. Avoid having to double up to meet throughput demands by choosing a flow cytometer that has additional capabilities, ready for future applications and experiments.

The Bio-Rad ZE5 Cell Analyzer can collect data at a higher rate than any other instrument currently available. The electronics and fluidics make it capable of sampling data at 100,000 events per second without compromising data quality, making rare event detection super-fast. Additionally, in high-throughput mode plates can be processed at speeds previously only available with instruments dedicated to screening: a 96 well plate may take less than 15 minutes and a 384 well plate less than 50 minutes. Combining these two advantages makes it possible to perform complex immunophenotyping panels quickly and maximize productivity.

Consider Automation

The ability to screen more and analyze faster while maintaining high quality, accurate results is key to progressing research. Automating your flow cytometry workflow can dramatically increase productivity and lead to greater progress. As not all instruments are suitable for automation, future proof your workflow by making sure that you opt for an instrument that was designed with automation in mind. Robotic integration and automation features simplify operations, save time, and improve lab performance. Automation providers will be happy to offer further assistance with setting up your workstation.



External Fluidics are especially useful if you are considering automation. The house deionized (DI) connection allows the ZE5 Cell Analyzer to be connected directly to a supply of DI water. The external bulk carboys replace the internal tanks to give up to 24 hr uninterrupted run time.

Automation-ready, the ZE5 Cell Analyzer has been integrated by many providers into a wide range of automation workcells. Large capacity fluidics options ensure uninterrupted 24 hour operation.

Visit bio-rad.com/htscreening for more information on ZE5 Cell Analyzer automation.

Identify Your Installation Site

Laboratory space is often at a premium, especially in more established labs. Think about space below, as well as above, the bench and don't forget to factor in access for servicing. If considering automation, the flexibility of modern robotics allows for workcells to be built in more limited spaces than previously so automation can be possible even in smaller spaces. It's a good idea to contact your local instrument representative as well as your automation provider to seek advice if space is limited.

Unlike many other systems, the ZE5 Cell Analyzer does not need a separate fluidics cart, all the fluidics can be contained within the instrument itself, although larger external tanks are available as well. With a remarkably small footprint, it can be installed on a standard benchtop.

Next Steps

Make a Shortlist and Use Your Network

Once you have a clear idea of what you need, make a shortlist of instruments that match your requirements. Then use your network to gain experience from your peers. Speak to your colleagues as well as the wider userbase on social media, at meetings, and at conferences to get a broad opinion of the instruments on your list.

Talk with a Specialist

Contact your local sales teams and speak with a specialist in flow cytometry within the company. They should ensure that the instrument you have selected meets your requirements and is the right configuration for you. This is your chance to obtain more detailed information, so make sure that you have a list of questions as well as providing as much detail as possible on what you need from the instrument. For a neutral appraisal of the technology, talk to other researchers using the same system you are considering.

Request a Demo

Once you have identified a flow cytometer that meets your requirements, put it through its paces by arranging a demonstration. Involve key stakeholders in the demonstration and perform your most important assays to ensure the machine will deliver what you hope. Plan to have suitable samples ready for your demonstration.

ZE5 Cell Analyzer Laser and Filter Configuration Guide

5-Laser (27 colors)			5-Laser 7 Off UV Option A (27 Colors)			5-Laser 7 Off UV Option B (27 Colors)		
Laser	Filter	Fluorochrome	Laser	Filter	Fluorochrome	Laser	Filter	Fluorochrome
355	387/11	SBUV400, BUV395	355	387/11	SBUV400, BUV395	355	387/11	SBUV400, BUV395
	447/60	SBUV445, AF350, DAPI		509/24	SBUV510, DAPI, Zombie UV		460/22	SBUV445, DAPI, Zombie UV, L/D Blue
	525/50	SBUV510, BUV496		577/15	SBUV575, BUV536		509/24	SBUV510, BUV496
	670/30	SBUV665, BUV661		615/24	SBUV605, BUV615		577/15	SBUV575, BUV563
	700LP	SBUV740, SBUV795, BUV737		670/30	SBUV665, BUV661		670/30	SBUV665, BUV661
405	420/10	BV421	405	747/33	SBUV740, BUV737	405	747/33	SBUV740, BUV737
	460/22	SBV440, Pacific Blue		780 LP	SBUV795, BUV805		780 LP	SBUV795, BUV805
	525/50	SBV515, BV510, CFP		420/10	BV421		420/10	BV421
	615/24	SBV610, BV605		460/22	SBV440, Pacific Blue		460/22	SBV440, Pacific Blue
	670/30	SBV670, BV650		525/50	SBV515, BV510, CFP		525/50	SBV515, BV510, CFP
	720/60	SBV710, BV711		615/24	SBV610, BV605		615/24	SBV610, BV605
	750LP	SBV760, SBV790, BV786		670/30	SBV670, BV650		670/30	SBV670, BV650
488	488/10	Side Scatter	488	720/60	SBV710, BV711	488	720/60	SBV710, BV711
	525/25	FITC, GFP, YFP, AF488, Kiravia520		750LP	SBV760, SBV790, BV786		750LP	SBV760, SBV790, BV786
	593/52	SBB580		488/10	Side Scatter		488/10	Side Scatter
	692/80	SBB675, SBB700, PerCP-Cy5.5		525/35	FITC, GFP, YFP, AF488, Kiravia520		525/35	FITC, GFP, YFP, AF488, Kiravia520
	750LP	SBB765, SBB810		593/52	SBB580		593/52	SBB580
561	577/15	SBY575, PE	561	692/80	SBB675, SBB700, PerCP-Cy5.5	561	692/80	SBB675, SBB700, PerCP-Cy5.5
	589/15	DsRed, tdTomato		750LP	SBB765, SBB810		750LP	SBB765, SBB810
	615/24	SBY605, PE-Dazzle, PE-CF594, mCherry		583/30	SBY575, PE, DsRed, tdTomato, RFP		583/30	SBY575, PE, DsRed, tdTomato, RFP
	640/20	mPlum		615/24	SBY605, PE-Dazzle, PE-CF594, mCherry		615/24	SBY605, PE-Dazzle, PE-CF594, mCherry
	670/30	SBY665, PE-Cy5, PE-AF647		670/30	SBY665, PE-Cy5, PE-AF647		670/30	SBY665, PE-Cy5, PE-AF647
	720/60	SBY720, PE-Cy5.5		720/60	SBY720, PE-Cy5.5		720/60	SBY720, PE-Cy5.5
640	750LP	SBY800, PE-Cy7, PE-AF750	640	750LP	SBY800, PE-Cy7, PE-AF750	640	750LP	SBY800, PE-Cy7, PE-AF750
	670/30	APC, AF647		670/30	APC, AF647		670/30	APC, AF647
	720/60	AF700		720/60	AF700		720/60	AF700
	775/50	APC-Cy7, AF750		775/50	APC-Cy7, AF750		775/50	APC-Cy7, AF750
	800LP	AF790, APC/Fire810		800LP	AF790, APC/Fire810		800LP	AF790, APC/Fire810

Note: This is not a comprehensive list of fluorophores and dyes that can be used with the ZE5 Cell Analyzer.

continues

ZE5 Cell Analyzer Laser and Filter Configuration Guide, continued

4-Laser (24 colors)***			3-Laser (17 Colors)***			3-Laser Option 2 (17 colors)**		
Laser	Filter	Fluorochrome	Laser	Filter	Fluorochrome	Laser	Filter	Fluorochrome
405	420/10	BV421	405	420/10	BV421	488	488/10	Side Scatter
	460/22	SBV440, Pacific Blue		460/22	SBV440, Pacific Blue		509/24	FITC, GFP, AF488, Kiravia520
	525/50	SBV515, BV510, CFP		525/50	SBV515, BV510, CFP		549/15	YFP
	615/24	SBV610, BV605		615/24	SBV610, BV605		583/30	SBB580, PE
	670/30	SBV670, BV650		670/30	SBV670, BV650		615/24	SBB615, PE-CF594, PE-Dazzle
	720/60	SBV710, BV711		720/60	SBV710, BV711		692/80	SBB675, SBB700, PerCP-Cy5.5, PE-Cy5
	750LP	SBV760, SBV790, BV786		750LP	SBV760, SBV790, BV786		750LP	SBB765, SBB810, PE-Cy7, PE-AF750
488	488/10	Side Scatter	488	488/10	Side Scatter	561	577/15	SBY575, PE
	509/24	FITC, GFP, AF488, Kiravia520		509/24	FITC, GFP, AF488, Kiravia520		589/15	DsRed, tdTomato
	549/15	YFP		549/15	YFP		615/24	SBY605, PE-Dazzle, PE-CF594, mCherry
	583/30	SBB580, PE		583/30	SBB580, PE		640/20	mPlum
	615/24	SBB615, PE-CF594, PE-Dazzle		615/24	SBB615, PE-CF594, PE-Dazzle		670/30	SBY665, PE-Cy5, PE-AF647
	692/80	SBB675, SBB700, PerCP-Cy5.5, PE-Cy5		692/80	SBB675, SBB700, PerCP-Cy5.5, PE-Cy5		720/60	SBY720, PE-Cy5.5
	750LP	SBB765, SBB810, PE-Cy7, PE-AF750		750LP	SBB765, SBB810, PE-Cy7, PE-AF750		750LP	SBY800, PE-Cy7, PE-AF750
561	577/15	SBY575, PE	640	670/30	APC, AF647	640	670/30	APC, AF647
	589/15	DsRed, tdTomato		720/60	AF700		720/60	AF700
	615/24	SBY605, PE-Dazzle, PE-CF594, mCherry		775/50	APC-Cy7, AF750		775/50	APC-Cy7, AF750
	640/20	mPlum		800LP	AF790, APC/Fire810		800LP	AF790, APC/Fire810
	670/30	SBY665, PE-Cy5, PE-AF647						
	720/60	SBY720, PE-Cy5.5						
	750LP	SBY800, PE-Cy7, PE-AF750						
640	670/30	APC, AF647						
	720/60	AF700						
	775/50	APC-Cy7, AF750						
	800LP	AF790, APC/Fire810						
3-Laser (20 Colors)***								
Laser	Filter	Fluorochrome						
405	420/10	BV421	405	420/10	BV421	488	488/10	Side Scatter
	460/22	SBV440, Pacific Blue, mTurquoise		460/22	SBV440, Pacific Blue, mTurquoise		509/24	FITC, eGFP, AF488, Kiravia520, mClover
	525/50	SBV515, BV510, CFP, Cascade Yellow		525/50	SBV515, BV510, CFP, Cascade Yellow		549/15	eYFP
	615/24	SBV610, BV605		615/24	SBV610, BV605		583/30	SBB580, PE
	670/30	SBV670, BV650		670/30	SBV670, BV650		615/24	SBB615, PE-CF594, PE-Dazzle
	720/60	SBV710, BV711		720/60	SBV710, BV711		692/80	SBB675, SBB700, PerCP-Cy5.5, PE-Cy5
	750LP	SBV760, SBV790, BV786		750LP	SBV760, SBV790, BV786		750LP	SBB765, SBB810, PE-Cy7, PE-AF750
488	488/10	Side Scatter	488	488/10	Side Scatter	561	577/15	SBY 575, PE
	509/24	FITC, eGFP, AF488, Kiravia520, mClover		509/24	FITC, eGFP, AF488, Kiravia520, mClover		589/15	DsRed, tdTomato, RFP
	549/15	eYFP		549/15	eYFP		615/24	SBY605, PE-Dazzle, PE-CF594, mCherry
	583/30	SBB580, PE		583/30	SBB580, PE		640/20	mPlum
	615/24	SBB615, PE-CF594, PE-Dazzle		615/24	SBB615, PE-CF594, PE-Dazzle		670/30	SBY665, PE-Cy5, PE-AF647
	692/80	SBB675, SBB700, PerCP-Cy5.5, PE-Cy5		692/80	SBB675, SBB700, PerCP-Cy5.5, PE-Cy5		720/60	SBY720, PE-Cy5.5
	750LP	SBB765, SBB810, PE-Cy7, PE-AF750		750LP	SBB765, SBB810, PE-Cy7, PE-AF750		750LP	SBY800, PE-Cy7, PE-AF750
561	577/15	SBY 575, PE	561	577/15	SBY 575, PE	640	670/30	APC, AF647
	589/15	DsRed, tdTomato, RFP		589/15	DsRed, tdTomato, RFP		720/60	AF700
	615/24	SBY605, PE-Dazzle, PE-CF594, mCherry		615/24	SBY605, PE-Dazzle, PE-CF594, mCherry		775/50	APC-Cy7, AF750
	640/20	mPlum		640/20	mPlum		800LP	AF790, APC/Fire810
	670/30	SBY665, PE-Cy5, PE-AF647		670/30	SBY665, PE-Cy5, PE-AF647			
	720/60	SBY720, PE-Cy5.5		720/60	SBY720, PE-Cy5.5			
	750LP	SBY800, PE-Cy7, PE-AF750		750LP	SBY800, PE-Cy7, PE-AF750			

Note: This is not a comprehensive list of fluorophores and dyes that can be used with the ZE5 Cell Analyzer.

* Small particle detector can be included in configuration or omitted from configuration with the option to add later.

** Eligible for an upgrade with a UV laser.

AFxxx, Alexa Fluor; APC, allophycocyanin; BV, Brilliant Violet; BUUV, Brilliant UltraViolet; CFP, cyan fluorescent protein; Cy, cyanine; dsRed, Discosoma Red Fluorescent Protein; FITC, fluorescein isothiocyanate; GFP, green fluorescent protein; eGFP, enhanced green fluorescent protein; eYFP, enhanced yellow fluorescent protein; LP, long pass; PE, phycoerythrin; PerCP, peridinin chlorophyll protein; RFP, red fluorescent protein; SBB, StarBright Blue; SBV, StarBright Violet; SBUV, StarBright UltraViolet; SBY StarBright Yellow; UV, ultraviolet; YFP, yellow fluorescent protein.

Visit [bio-rad.com/cellanalysis](https://www.bio-rad.com/cellanalysis) for a wide range of videos, technical reports, published findings, and other resources describing the Bio-Rad ZE5 Cell Analyzer.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.
© 2022 Bio-Rad Laboratories, Inc.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399
Canada 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23
France 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050
Italy 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 2 3473 4460 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670
The Netherlands 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23
Russian Federation 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23
Switzerland 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

