

Go beyond cell counting to detailed analysis.

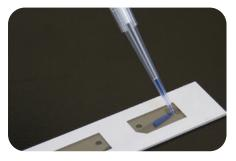
Cellometer Spectrum Image Cytometry System

revvity

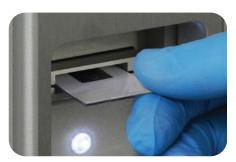
The Cellometer® Spectrum is designed for 20 µL cell-based assays

- All-in-one system
 Basic cell counting, primary cell viability, and cell-based assays.
- Dual-fluorescence for accurate primary cell viability
 No interference from red blood cells. Analyze bone marrow, peripheral blood, and cord blood without lysing.
- Unique algorithms for advanced cell analysis
 Determine concentration and viability of hepatocytes, adipocytes, and other sophisticated cell types.
- Fast results
 Obtain cell images, counts, size measurements and viability calculations in < 30 seconds per sample.

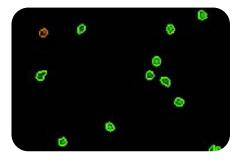




1. Pipette 20 µL



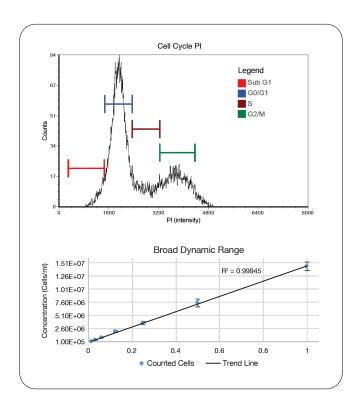
2. Insert slide & count



3. Get images & data

Simple 20 µL cell-based assays

- Accurate Cell Counting
- Versatile Analysis and Applications
- User-changeable Fluorescent Modules
- Pre-qualified Reagents
- Pre-defined Instrument Settings
- Assay-Specific data templated
- Export to FCS Express* (Flow like data analysis)
 - GFP
 - Apoptosis
 - Cell cycle
 - Cell mediated cytotoxicity
 - Cell counting & viability



User-changeable fluorescence optics modules**

Cellometer assays use specific optic modules for maximum performance and discrimination between fluorescence channels. Example filter modules are listed below. Additional filter modules available upon request.

Optics module	Fluorophores	Nucleic acid stains
S1-534-470 Ex: 470nm (452-498nm) Em: 534nm (510-560nm)	GFP Calcein FITC AlexaFluor® 488	AO (acridine orange) SYTO®9, SYTO®13 SYTOX®Green SYTO®BC
S1-605-527 Ex: 527nm (498-550nm) Em: 605nm (570-638nm)	AlexaFluor® 546 AlexaFluor® 555 Cy3® PE (R-phycoerythrin) RFP	PI (propidium iodide) EB (ethidium bromide) SYTOX® Orange
S1-655-527 Ex: 527nm (525-558nm) Em: 655nm (642-665nm)	RFP 7-AAD Nile Red	PI (propidium iodide) EB (ethidium bromide)

^{**} This table is a partial list of compatible fluorophores, nucleic acid stains, and fluorescent proteins. Please contact technical support regarding compatibility of other reagents.

^{*} FCS Express Flow Cytometry software is a product of De Novo Software and is included with the Cellometer Spectrum





 $^{^{\}ast}$ Sytox, Alexa Fluor, and Cy are trademarks of Life Technologies.