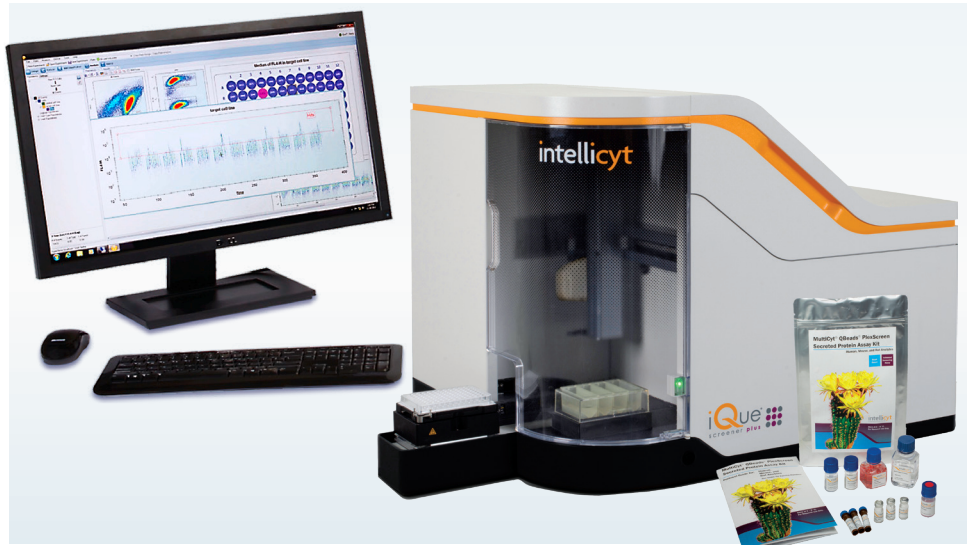


IntelliCyt® iQue Screener PLUS

Fast, multiparameter data analysis from suspension-cell assays



Profile and functionally characterize cells relevant to disease and drug mechanism of action (MOA), and gain the valuable insight needed to make more effective drug development decisions sooner in the process.

The IntelliCyt® iQue Screener PLUS Platform integrates a flow based instrument with enabling and easy to use software, and reagents designed to address key challenges across the suspension-cell screening workflow.

The IntelliCyt Advantage



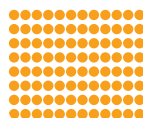
Speed

- Faster plate processing, minutes, not hours.
- Mix and read samples.
- Faster time to results.



Miniaturization

- Consumes less reagents.
- Conserves precious cells.
- Saves money.



Content

- Rich, multiplexed, per-cell content.
- Cell and beads together.
- Secreted protein analysis.



Usability

- Automated workflow.
- Validated reagents.
- Easiest software you will ever love.



Insight

- Link information.
- Run scenarios.
- Create knowledge.
- Make decisions.

IntelliCyt iQue Screener PLUS Platform

The IntelliCyt® iQue Screener PLUS platform is an integrated instrument, software and reagent system that enables rapid, high content, multiplexed analysis of cells and beads in suspension. When speed is a factor or when cells are precious or limited in number this user-friendly system excels. Our unique, software-assisted automation and experiment-based analyses deliver the deep insight needed to answer complex biological questions.

The IntelliCyt iQue Screener PLUS (**Blue-Red** laser configuration) is a phenotypic screening and profiling workhorse that is ideal for applications that require up to 6-color detection, including immunophenotyping, cell health assessment, secreted protein analysis using QBead-based assays and many more applications. Our platform delivers the ForeCyt® Software Workflow Advantage: a single data management workflow from input to output, which means you work faster and work smarter—not harder.

Content is king with the IntelliCyt iQue Screener PLUS (**Violet-Blue-Red** and **Violet-Yellow-Blue** laser configurations). Three-laser configurations offer up to 13-color detection and are ideal for functional and phenotypic applications that demand more choice and flexibility in experimental design. By maximizing the detection and resolution of traditional and/or innovative new reagent dyes, tandem dyes, and fluorescent proteins, IntelliCyt iQue Screener PLUS delivers both high performance multicolor analysis and the ForeCyt Software Workflow Advantage making it hands-down the choice of leaders in immune-based drug discovery, immuno-oncology, and cell therapy applications.

The IntelliCyt iQue Screener PLUS HD (**Blue-Red** configuration) provides ultimate assay miniaturization and is the only high content, per-cell, 1536-well capable suspension screener available.

IntelliCyt iQue Screener PLUS Technical Specifications

	IntelliCyt iQue Screener PLUS Configuration	Blue and Red		Violet, Blue and Red			Violet, Yellow and Blue		
	Lasers	488 nm	640 nm	405 nm	488 nm	640 nm	405 nm	561 nm	488 nm
Detectors	445/45 nm			●			●		
	530/30 nm	●		●	●		●		●
	572/28 nm	●		●	●				
	586/20 nm						●	●	
	615/24 nm			●	●				
	615/20 nm						●	●	●
	660/20 nm						●	●	●
	675/30 nm	●	●	●	●	●			
	780/60 nm	●	●	●	●	●	●	●	
	Forward Light Scatter (relative size)	●			●			●	
	Side Light Scatter (relative granularity)	●			●			●	
Optical	Fluorescence Sensitivity	FITC < 75 MESF; PE < 50 MESF; APC < 20 MESF							
	Minimum Particle Size Detection	0.5 µm							
	Cell Detection Rate	Up to 35,000/second							
	Dynamic Range of Detection*	> 7 Decades							
* This wide dynamic range and a Zoom function permit operation of the system without user adjustments of the voltage or gain of the detectors.									
Sampling	Plate Compatibility	96-well, 384-well or 384-well, 1536-well (iQue Screener PLUS HD BR)							
	Sampling	Continuous Air-gap Delimited							
	Minimum Assay Volume Requirements	10 µL							
	Minimum Sample Aspiration	1 µL							
	Minimum Plate Sampling Time*	< 5 minutes / 96 wells < 20 minutes / 384 wells							
	Carryover	< 2% for typical no-wash assays. Actual amounts are cell and assay dependent and are easily managed by including interwell rinses to reduce carryover to < 0.1%							
	Automated Plate Shaker	Up to 3,000 RPM (Up to 5000 RPM on IntelliCyt iQue Screener PLUS HD BR)							
	Foil-sealed Plate Processing	●							
	Volumetric Cell Counting (< 10% CV)	●							
	* The time required for sampling plates is both sample type and experiment dependent. A range of well-sampling times can be designated from 0.5 seconds – minutes.								
ForeCyt® Software	Auto Compensation	●							
	Real-time Whole-plate Data Analysis	●							
	Dynamic Linked Gating	●							
	Interactive Heat Maps, Profile Maps	●							
	Cross Plate Analysis	●							
	Export Files in FCS, CSV or ForeCyt Formats	●							
	Customizable PDF Data Report	●							
	ForeCyt Enterprise Edition Compatible	●							
Operational	Computer Workstation, Windows Compatible	Xeon Processor, Dual 256 GB SSD (RAID 0), 16GB RAM, 27" Monitor 2560 x 1400							
	Weight (less computer)	205 lbs, 93 kg							
	Dimensions	39" W x 25" D x 26" H • 99 cm W x 63 cm D x 66 cm H							
	Power Requirements	100/115/230 VAC, 50-60Hz							
	Environment Requirements	Temperature: 15–32°C (59–90°F), Relative Humidity: 80% Maximum							
	CE Labeled	●							
	21 CFR Logging Option Compatible	●							
	Robotic Integration Option Compatible	●							
	QMax Refill Module Option Compatible	●							

Learn why IntelliCyt is the choice of leaders in Immuno-Oncology, Antibody Discovery and Immune Target Screening

www.intellicyt.com

For more information or a quote, call us at +1 505-345-9075 or send an email to sales@IntelliCyt.com

©2017 IntelliCyt Corporation. All Rights Reserved. The trademarks used herein are the property of IntelliCyt Corporation or their respective owners.



A SARTORIUS COMPANY

IntelliCyt technology is protected by the following patents and other patents pending: 6,890,487, 6,878,556, 7,368,084, 7,842,244, 8,021,872, 8,268,571, 8,637,261, 8,823,943, 9,012,235, D,722,515