



AQUIOS CL Load & Go Flow Cytometer and Reagent Kits

Empowering Life Changing Decisions



ACCELERATING
answers

 **BECKMAN
COULTER**
Life Sciences

Enabling 24/7 flow cytometry service

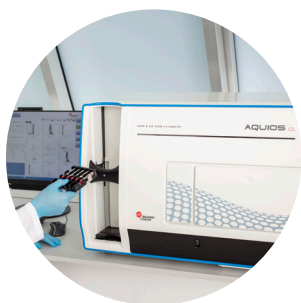


Compact size

with dimensions W: 32" (82cm),
D: 22" (56cm), H: 22" (56cm)
to easily fit your lab space.

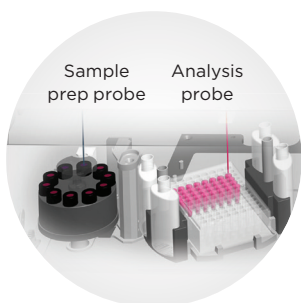
When patients are waiting for test results, few things are more important than your lab's efficiency.

To help optimize lab efficiency and productivity—while maintaining the highest standards of accuracy—consider the AQUIOS CL flow cytometry system, which provides:



“Load & Go” Cytometry

- Makes sample loading as easy as using a hematology analyzer
- Simply load primary samples and receive results



Automated sample preparation

- Eliminates the need for manual sample prep
- Enables simultaneous sample prep and data acquisition



More time for other crucial tasks

- Lets your team focus on data review and validation of results

Boost your efficiency with true walk-away automation

The AQUIOS CL Load & Go Flow Cytometer helps streamline your assays so your laboratory can perform at its highest potential. This closed, fully automated system enables:

Random-access sample loading

Cassette autoloader lets you load 8 cassettes at one time (with up to 5 sample tubes each) for continuous loading/unloading—**no batching required**

Separate single-tube loader for STAT samples, open vial samples or irregular tube sizes



On-board sample preparation

Automated sample preparation matches samples to test requests, prepares them in 96-well microplates with no manual pipetting, and launches protocols

Pipeline sample processing enabled by technology that uses 2 multitasking probes.

Barcoded reagents for tracking/monitoring reagent consumption as samples are processed

Cap piercing to help prevent sample contamination

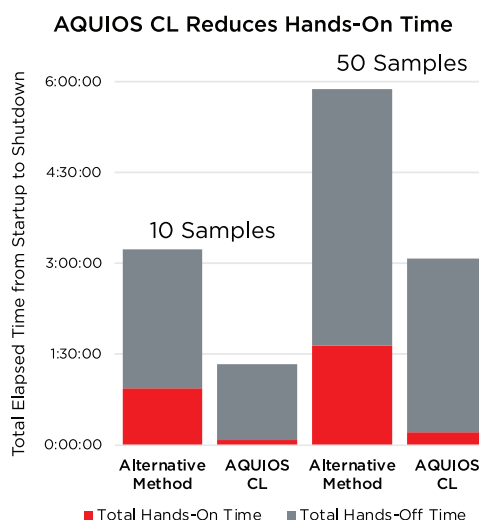
Error minimization

Digital data transfer via barcode scanner and bidirectional LIS-link

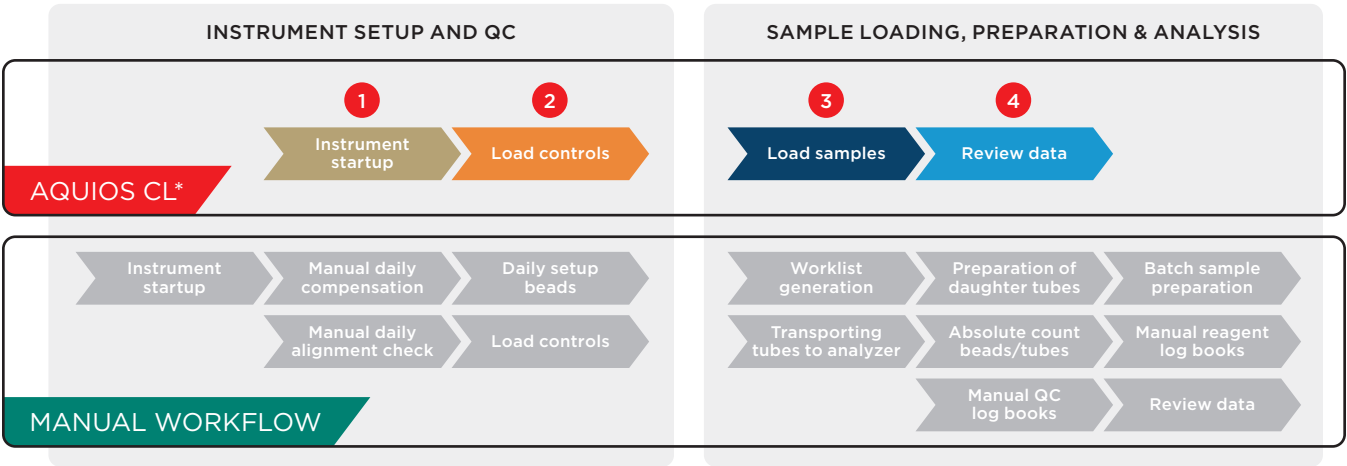
Comprehensive quality control requiring no manual QC while generating a full, easily reviewable audit trail

More time for techs to:

- **Analyze data and review results**
- **Perform value-adding tasks**
- **Develop/hone new skills**



When you have routine, low-complexity assays to analyze, you can increase laboratory efficiency by automating your processes.



*Only steps that require manual intervention are shown

"The machine is able to analyze leucocytes in a sample of whole blood in only 20 minutes. In addition, the data obtained can be exported directly to a hospital file system. This allows cellular diagnostics in a 24/7 context with high reproducibility."

Leo Koenderman, Ph.D.
University Medical Center, Utrecht

"Since it is so simple and intuitive, training my techs was a breeze. You can train them in less than an hour. So, we were able to cross-train most of the techs in the area in very short time. That makes everything efficient in the laboratory."

Alexandra Amador, MT (ASCP), CHT (ABHI)
Manager Histocompatibility Lab
University of Miami

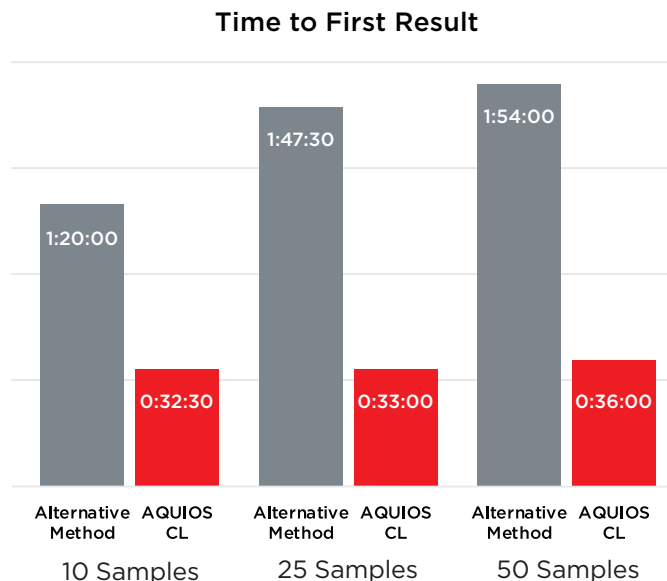
Deliver critical results faster

The AQUIOS CL system expedites results by providing:

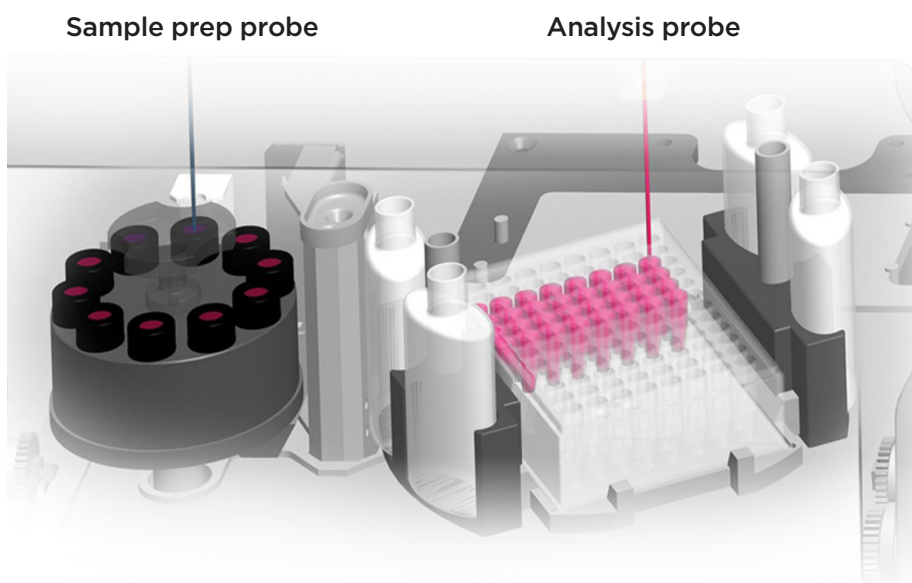
- Capabilities for 24/7 flow cytometry service
- Full bidirectional LIS connectivity
- Smart track reagent monitoring
- Analysis as soon as samples are ready
- Digital data transfer

Time to first result* was analyzed for lymphocyte subset samples on AQUIOS CL vs. an alternative method (BD FACSCalibur plus SPA II system)

* Sum of cytometer startup & cleaning, QC & sample prep and time for analysis



Two probes enable faster processing



A system that supports your desire for multitasking

- Continuous loading (no batching required)
- Simultaneous sample prep & acquisition
- Analyze samples as soon as they're ready

Minimize human error

The AQUIOS CL system provides:

- Real-time sample monitoring
- Single Platform Technology (SPT) for lymphocyte subset analysis/enumeration with no need for counting beads
- Reviewable audit trail
- User alerts via email/text in case of a QC fail
- Intelligent analysis algorithm that sets all gates and regions automatically
- Daily instrument quality control
 - Measure 2 levels of cellular process controls
 - No setup beads, alignment beads or compensation reagents
- Reagent tracking via barcodes

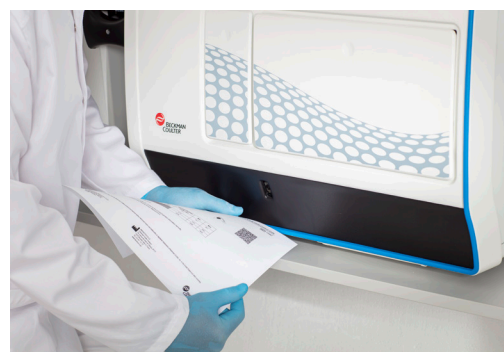


Reduce Error-Prone Steps Up to 87.5%

Real-time specimen quality control

Automatically checks every patient sample for:

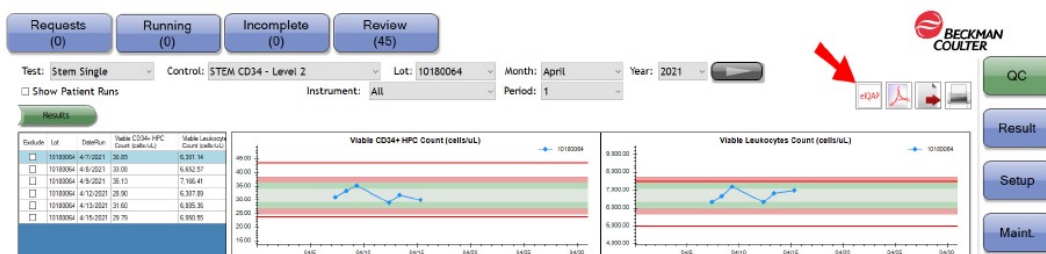
- Stable event rate vs. time
- Stable mean channel of key cell populations
- FL compensation
- Separation quotient
- Correct location/appearance of all cell populations
- Appearance of increased debris due to under-lysis of RBCs



Transfer target values from the assay sheet with the barcode scanner

Interlaboratory Quality Assurance Program (IQAP)

This unique program lets you further reinforce accuracy by comparing results from control runs with those from peer groups.

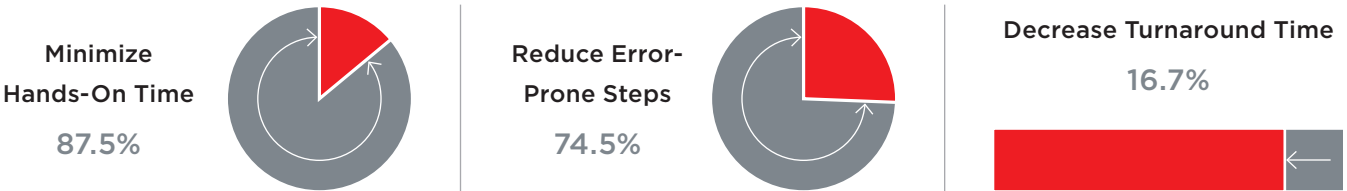


Ensure consistent performance with AQUIOS reagent kits

AQUIOS TETRA KIT

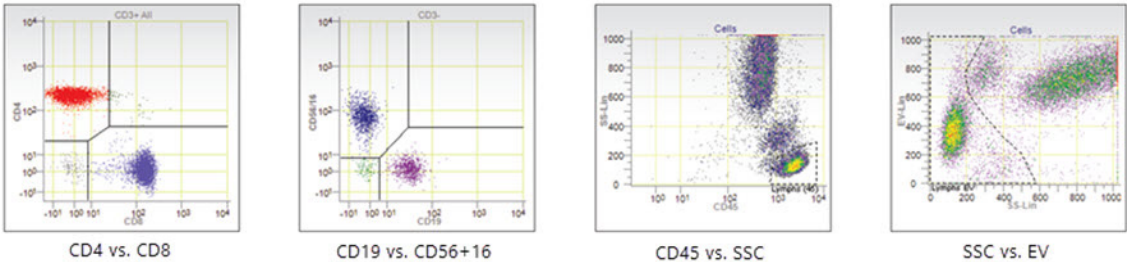
Intuitive lymphocyte subset analysis

- Analyze T, B and NK lymphocytes in whole blood
- Use STAT feature for urgent requests
- Enumerate without a need for counting beads
- Get peer group comparison reports for your laboratory's QC data using our Interlaboratory Quality Assurance Program



AQUIOS CL system (■) running AQUIOS Tetra-1 was compared to the FC 500 and TQ Prep systems (■), running CYTO-STAT tetraCHROME CD45-FITC/CD4-RD1/CD8-ECD/CD3-PC5 for 50 test samples.

	FITC	PE	ECD	PC5
AQUIOS Tetra-1	CD45	CD4	CD8	CD3
AQUIOS Tetra-2+	CD45	CD56+16	CD19	CD3



AQUIOS STEM KIT



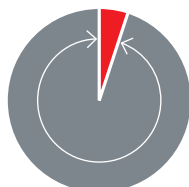
Evolution of the “Gold Standard”

- Fully automates ISHAGE-compliant CD34+ progenitor cell analysis
- STAT capabilities for urgent requests
- AQUIOS STEM software includes choice of 3 panels:
 - Duplicate + negative control
(CD45-FITC/CD34-PE/7-AAD, CD45-FITC/CD34-PE/7-AAD, CD45-FITC/CD34-CTRL/7-AAD)
 - Duplicate W/O negative control
(CD45-FITC/CD34-PE/7-AAD, CD45-FITC/CD34-PE/7-AAD)
 - Single test
(CD45-FITC/CD34-PE/7-AAD)



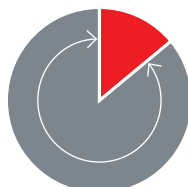
Minimize
Hands-on Time

95%



Reduce Error-
Prone Steps

87.5%

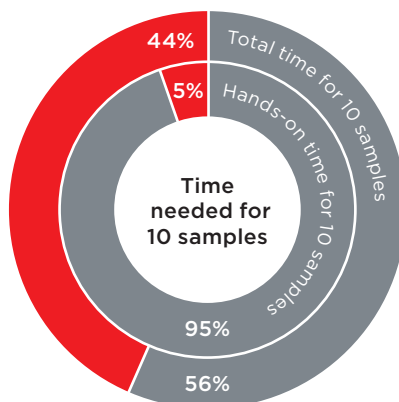
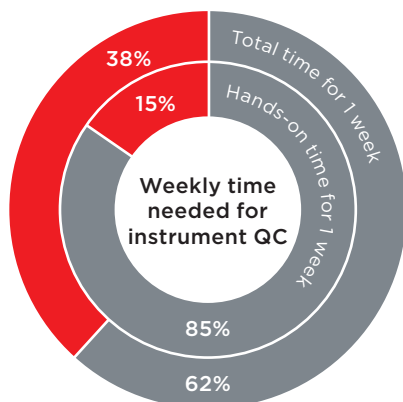


Decrease Turnaround Time

23%



AQUIOS STEM system (■) was compared directly to the FC 500 and Stem-Kit (■) to run 10 test samples.



Decrease time for weekly
instrument QC and sample runs

While time-to-result is largely determined by incubation and lysing times, hands-on time could be reduced by up to 95% compared to an alternative method.

■ AQUIOS CL ■ Alternate method

Add even more flexibility with AQUIOS Designer Software (ADS)

Using AQUIOS Designer Software with your AQUIOS CL System, enables you to:

- Create custom applications that integrate sample preparation and analysis
- Tailor your specific assay requirements
- Utilize user-defined reagent combinations

AQUIOS
Designer Software
LOAD & GO FLOW CYTOMETRY



"It's truly a Load & Go System. For us to be able to just load the system and have the samples upfront for routine analysis which for our uses was T cells, B cells and NK cells, was extremely useful. My overall level of satisfaction is that this is an excellent system."

Philip Ruiz, MD, Ph.D., Director
Transplantation Laboratories and Immunopathology, University of Miami



Designed for Sustainability

At Beckman Coulter Life Sciences, we are dedicated to pioneering sustainable solutions that empower a greener future for generations to come. We are integrating sustainability into the development of new products and packaging.



Service and support—by flow cytometrists FOR flow cytometrists

Our service/support team members include:

- Experts holding MS, MD & Ph.D. degrees
- Immunologists
- Cell biologists
- Virologists...and more

And they're a great source of information we use to improve our current portfolio and innovate new instruments.

BeckmanConnect

- Remote desktop sharing (view/control)
- File transfer (to and from instrument controller PC)
- Downloadable and customer installable
- Customer must approve all remote activity



Scan here—or contact your local Beckman Coulter Life Sciences sales representative



Ordering information

Systems

Part Number	Description
B39101	AQUIOS CL Flow Cytometry System with Uninterruptible Power Manager (100-120V)
B39102	AQUIOS CL Flow Cytometry System with Uninterruptible Power Manager (220-240V)

Reagents and beads

Part Number	Description
B25697	AQUIOS Sheath Solution, 10L
B25698	AQUIOS Cleaning Agent, 500 mL
C48097	AQUIOS Cleaning Agent, 500 ml (REACH)
B23536	AQUIOS Sodium Hypochlorite Solution, (4 X 50 mL)
B23538	AQUIOS Lysing Reagent Kit, 100 tests
B23502	AQUIOS Deep Well Plate (case of 50)
6605359	Flow-Check Beads, 3x10mL

AQUIOS STEM kits and reagents

Part Number	Description
B77691	AQUIOS STEM Kit, 50 tests
C43667	AQUIOS STEM CD34+ Control Cells, 2 levels, 15 tests per level
C89793	AQUIOS STEM Upgrade Kit

AQUIOS TETRA kits and reagents

Part Number	Description
B23533	AQUIOS Tetra-1 Panel, 50 Tests
B23534	AQUIOS Tetra-2+ Panel, 50 Tests
B23535	AQUIOS IMMUNO-TROL Cells (2x3.0 mL)
B25700	AQUIOS IMMUNO-TROL Low Cells (2x3.0 mL)

NOTE: AQUIOS and reagents are CE-IVD. The AQUIOS CL Flow cytometer in combination with the AQUIOS Tetra-1 Panel has been accepted for the WHO list of prequalified IVD products.

Choose Beckman Coulter Life Sciences

for Benchmark Expertise
and Innovation

Since 1935, Beckman Coulter has driven innovation. We remain committed to shaping flow cytometry technology that fits seamlessly into your lab's workflow and provides an optimal user experience. When you choose a Beckman Coulter solution you receive a high level of expertise, innovation, and quality assurance.

Contact your local Beckman Coulter sales representative.



Scan for more information on
AQUIOS CL Load & Go Flow
Cytometer and Reagent Kits



CE 2797

© 2025 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo and the Beckman Coulter product and service marks mentioned herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries. All other trademarks are the property of their respective owners.

For Beckman Coulter's worldwide office locations and phone numbers, please visit Contact Us at beckman.com
2024-GBL-EN-107086-V1

