



GALLIOS

*Flow Cytometry —
Powerful Versatile Performance*

 **BECKMAN
COULTER**
Life Sciences



Gallios Flow Cytometer

Speed, reproducibility, simplicity, reliability...

These are the goals that every flow cytometry lab aims to achieve. Beckman Coulter shows the way with the Gallios Flow Cytometer, a state-of-the-art system designed with your research needs in mind. Gallios delivers consistent and reliable analysis of multi-color assays.



Speed...

With up to four lasers, ten colors and advanced optical design, Gallios delivers high resolution in a short amount of time, with accurate processing of your flow cytometry data.

Reproducibility...

Gallios delivers powerful, consistent performance and data integrity in the most complex analytical environments. Acquire quality data efficiently and accurately, and get stable performance over long periods of time and across a wide range of operating temperatures.

Simplicity...

A multi-sample loading system is designed with busy research labs in mind, and educational support enhances ease and convenience for users.

Reliability...

The optional **PRO Service Remote Diagnostic System** helps keep Gallios operating at peak performance, with minimal downtime. Troubleshoot and resolve any issues promptly and efficiently and, in some cases, identify potential issues that could affect your laboratory productivity.

Enjoy sophisticated technology and flexible options with Gallios Flow Cytometer

Gallios delivers excellent resolution

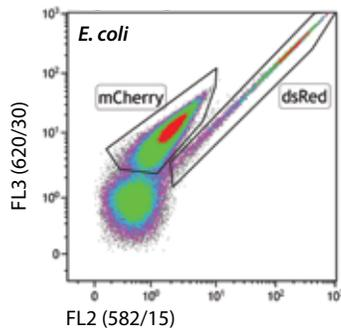
Built on a platform designed with reliability and stability in mind, Gallios samples information at 40MHz and displays data on a 1,048,576 channel scale. These specifications offer resolution of cell/particle characterization without sacrificing analytical speed or data integrity.

An innovative forward-scatter detector enables excellent resolution of submicron particles down to 0.404 μm in diameter. A side-scatter detector incorporates an independently focused, high-performance photodiode with electronic attenuation. Six fluorescence detectors provide simultaneous acquisition of up to six fluorescence signals.

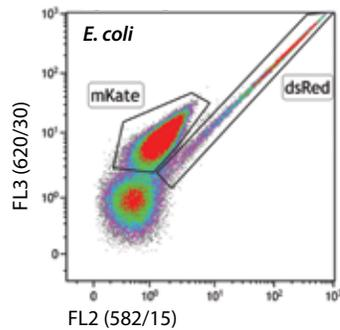
The electronics provide accurate and efficient signal processing at high event rates. A selection of up to 62 parameters can be processed per analysis at acquisition rates of 25,000 events per seconds with high yield. Gallios delivers a stable performance over long periods of time and across a wide range of operating temperatures.

Whether your challenges are dim markers, rare events, or just routine analysis, the Gallios is the cytometer of choice to capture all of your events – collecting four times the information as the cell passes through the interrogation point and displaying the information on a scale with four times the resolution.

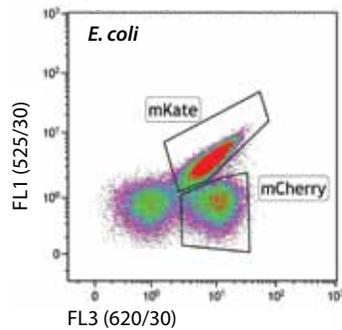
Panel A



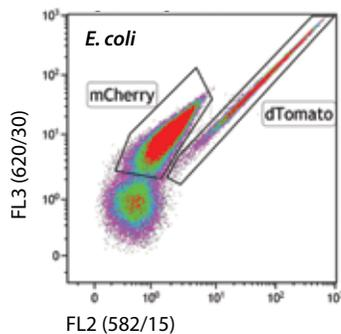
Panel B



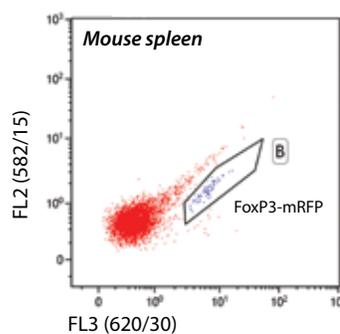
Panel C



Panel D



Panel E



Fluorescent protein detection with 488 and 561 nm excitation. Escherichia coli transfected with mCherry, dsRed, mKate and dTomato could be resolved in FL1, FL2, and FL3 (panels A-D). Mouse spleen cells transfected with FoxP3-mRFP (blue) are shown in panel E.

Gallios delivers flexibility to expand your research

When you incorporate the optional 405nm violet laser and an optional 561nm yellow laser, you get a greater choice of fluorochromes to perform complex multicolor experiments. Easily interchangeable optical filters facilitate detection of a variety of dyes and wavelengths. You can also add up to four additional fluorescence detectors, enabling you to read up to 10 colors concurrently.

Gallios makes it easy to detect fluorescent proteins with the optional 561nm Laser

Gallios is now equipped for an optional 561nm laser to allow for expanded applications and greater choice of fluorochromes for multicolor experiments. With the new 561nm laser system you can analyze multiple fluorescent proteins simultaneously, as well as benefit from a greater ability to detect red fluorescent proteins.

- Analyze red fluorescent proteins such as mCherry and DsRed.
- Take full advantage of improved PE tandem dyes for enhanced detection.
- Versatile for both phenotypic and functional analysis using fluorescent antibodies, dyes and proteins.



State-Of-The Art Optics

The Gallios Flow Cytometer is in a class by itself with its advanced optical design that provides efficient acquisition of superior quality data. The Gallios is built with two highly stable, solid-state lasers in standard red and blue. You can choose the option to add up to four lasers by opting for violet and yellow lasers. Easily interchangeable optical filters facilitate detection of a variety of dyes and wavelengths.

An innovative forward-scatter detector enables superior resolution of submicron particles down to 0.404 μm in diameter. A side-scatter detector incorporates an independently focused, high-performance photodiode with electronic attenuation. Six fluorescence detectors provide simultaneous acquisition of up to six fluorescence signals.

Gallios is available in 2, 3 and 4 laser configurations:

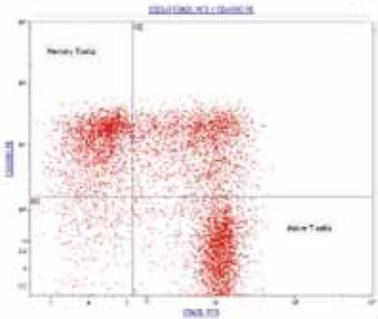
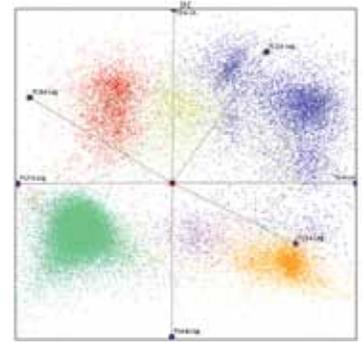
- 6 color, 2 laser (488nm Blue & 638nm Red) (5 + 1)
- 8 color, 2 laser (488nm Blue & 638nm Red) (5 + 3)
- 10 color, 3 laser (488nm Blue, 638nm Red & 405nm Violet) (5 + 3 + 2)
- 10 color, 4 laser (488nm Blue & 561nm Yellow [co-linear], 638nm Red, 405nm Violet) (5 + 3 + 2)

** The optional 561nm yellow laser is co-linear with the blue 488nm laser*



Radar Plot

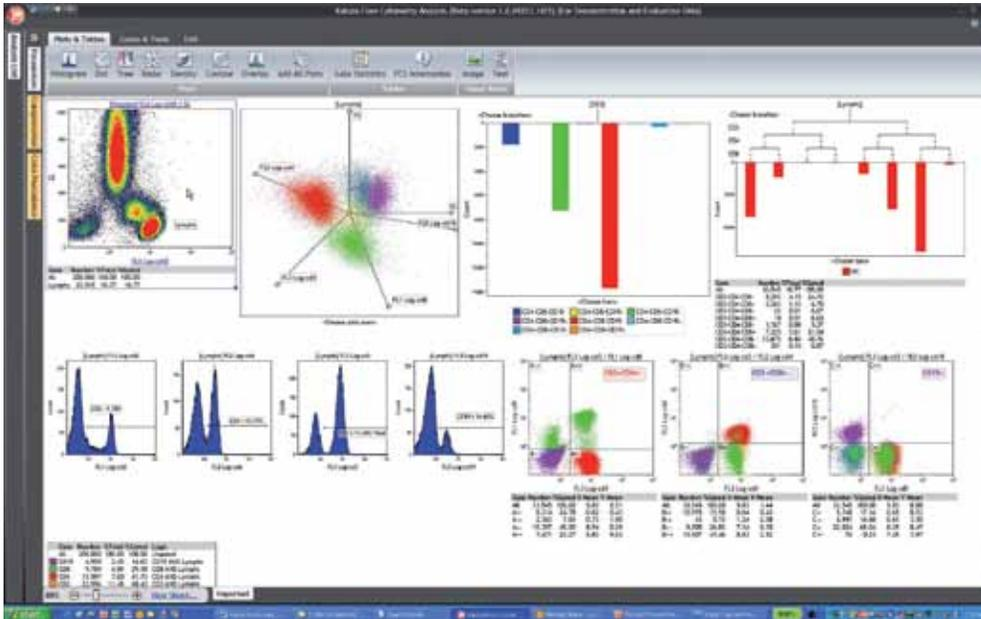
To enable multi-parametric pattern visualization without gating



Logicle Display

Adjustable linear/log transition

Slider bars to adjust compensation



Elevate your data analysis with Kaluza Flow Cytometry Analysis Software

Today, more colors require more plots, events, protocol complexity and ultimately increased analysis time. Kaluza is cutting edge flow cytometry software that performs powerful real-time data analysis of complex multi-color files. Its revolutionary speed (analyzing millions of cells, not thousands) makes analysis of multiple data sets as simple as drag and drop.

Kaluza helps you simplify the management of multiple data sets, enables visualization of high-content data in different spatial dimensions on a single plot, and provides real-time analysis of high-content flow cytometry files. The software features automatic plot organization, a zoom in/out feature to enhance data exploration, and auto-layout, which reconfigures the workspace in order to reduce the process of maintaining an analysis.

Kaluza features a variety of innovative plot types, including tree, 3D radar and overlay histogram as well as standard scatter, histogram and contour plots. Kaluza works with the majority of FCS-compliant files, can operate on either network or stand-alone computers and is supercomputer compatible with the NVIDIA Tesla graphics card.

Download a free 30-day trial version of Kaluza Analysis Software at: www.KaluzaNow.com



Fast, accurate, efficient digital electronics, extraordinary sensitivity and resolution; spatially separated solid state lasers with up to 10 colors give flexibility, accuracy and quality.

Maintain peak performance with **PROService** Remote Diagnostics

Not only do we care about designing and manufacturing quality instrumentation, we also want to ensure that your Gallios Flow Cytometer maintains peak performance. To that end, we created PROService, a remote diagnostic system that enhances the technical support of your flow cytometer.

PROService delivers prompt, efficient troubleshooting and issue resolution, minimizing downtime. In certain circumstances, PROService can preemptively identify potential problems and correct them before they impact your laboratory productivity. Count on Beckman Coulter to provide you with 24-7 support from our expert field consultants.



Gallios Flow Cytometer

Make the most of your flow cytometry capabilities with the powerful versatile performance of Gallios. Contact your Beckman Coulter representative or visit us at: www.GalliosNow.com

Download a free 30-day trial version of Kaluza Analysis Software at: www.KaluzaNow.com

fast track to success.



Gallios and Kaluza are for research use only. Not for use in diagnostics procedures.

Gallios, Kaluza and Beckman Coulter and the stylized logo are trademarks of Beckman Coulter, Inc. and are registered with the USPTO.

NVIDIA is a registered trademark of NVIDIA Corporation.

BR-17146C B2012-13567 © 2013 Beckman Coulter, Inc.

