

## Kaluza for Gallios

*Easy To Learn & Easy To Use—At All Experience Levels*

Kaluza for Gallios from Beckman Coulter employs cutting-edge technology to make high complexity flow cytometry easy for even the least experienced, while providing powerful tools for the most sophisticated user.

Kaluza for Gallios is built on the strong, innovative foundation of Kaluza Analysis Software.

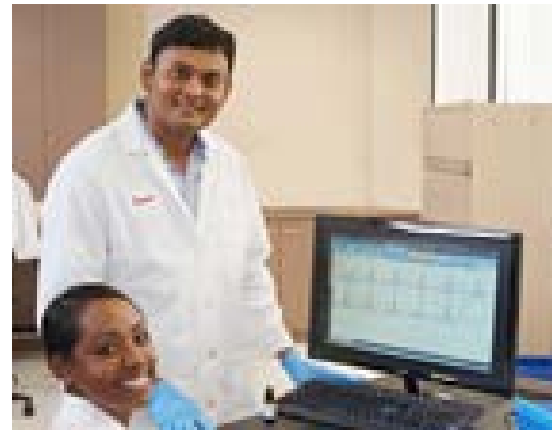
With Kaluza for Gallios, you can power through datasets containing multiple colors, plots and events with speed and precision. The result: your flow cytometry research can find that extra gear to get you to your objective sooner.

With up to four lasers scalable to 10 colors, sophisticated forward and side scatter detectors, state-of-the-art optics, and patented electronics, Gallios delivers powerful, versatile performance. Kaluza for Gallios provides the power to make the most of the data generated.

## Simplify and Speed Up Your Data Acquisition

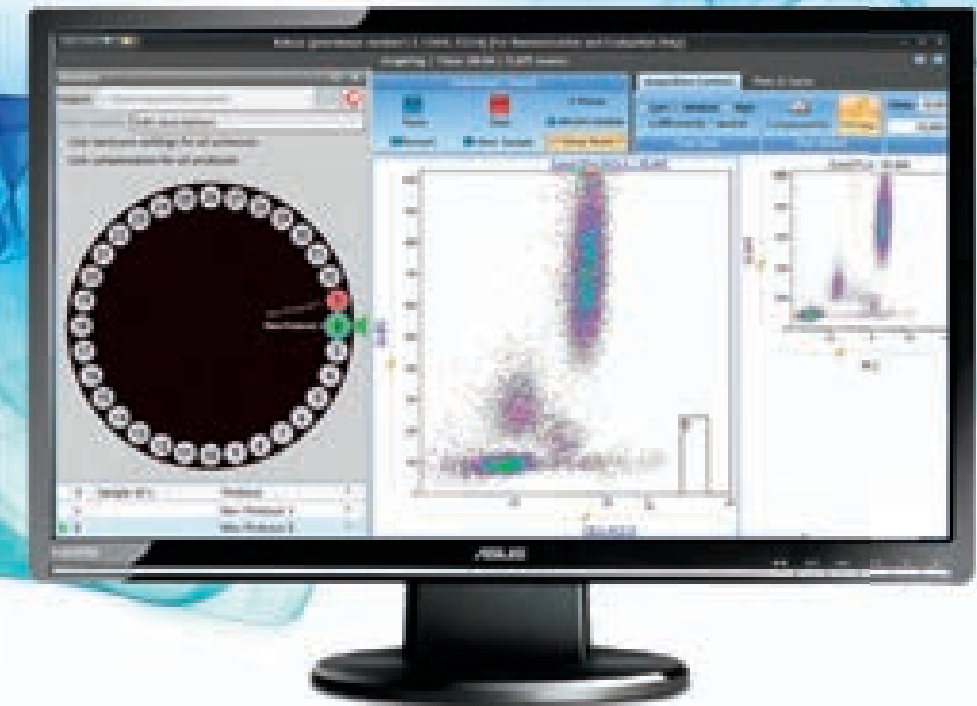
Flow cytometry experiments may be complex, but setting them up doesn't have to be hard. Kaluza for Gallios makes it easy—and fast—to spend more time on research and discovery and less time getting your flow experiments up and running. Learn more about Kaluza for Gallios, and get on the faster road to success.

For more information, go to [www.kaluzasoftware.com](http://www.kaluzasoftware.com)



# Kaluza for Gallios Acquisition Software

Sophisticated Flow Cytometry, Simplified.



Life Sciences

For research use only. Not for use in diagnostic procedures.  
Kaluza, Beckman Coulter and the stylized logo are trademarks of Beckman Coulter, Inc. and are registered with the USPTO. Gallios is a trademark of Beckman Coulter, Inc.

For Beckman Coulter's worldwide office locations and phone numbers, please visit "Contact Us" at [www.beckmancoulter.com](http://www.beckmancoulter.com)

FLOW-57SB04.14-B

© 2014 Beckman Coulter, Inc.

PRINTED IN U.S.A.



**KALUZA**

acquisition software for Gallios



Life Sciences

IT IS EASY TO LEARN, EASY TO USE, AND MAKES SHORT WORK OF SETTING UP AND RUNNING COMPLEX MULTICOLOR FLOW EXPERIMENTS.

## Kaluza for Gallios acquisition software:

### SPEED

- Set up a 10-color experiment quickly.
- Change voltages and compensation quickly, accurately and intuitively.
- Create compensation matrices with ease using compensation wizard.
- Create gates and access editing tools with innovative radial menu that keeps mouse clicks to a minimum.
- Load files into Kaluza for Gallios Analysis Software for more in-depth analysis.
- Configure, load and walk away with flexible carousel control.
- Create a protocol with one click using the “Add All Plots” feature.

### EASE

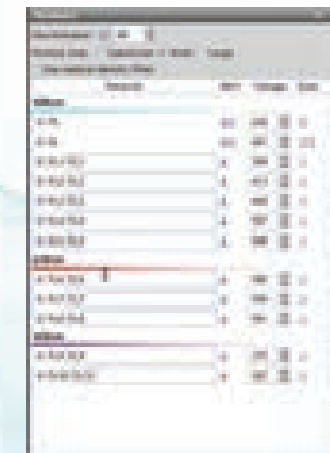
- Save your files to where you specify with no database to impact performance.
- Learn quickly and easily with simple user interface.
- Reconfigure your workspace with auto-layout feature for easy plot creation and sizing, zoom in/out, and guided drag and drop.
- Understand what is happening to your experiment and cytometer with visual management tools.
- Make a mistake? Undo it, with unlimited undo!
- Set voltages and compensation while in Setup Mode, then acquire data.
- Provide as much or as little sample information as you like before running an experiment.

### EFFICIENCY

- Utilizes eight times higher event limit than existing Gallios software.
- Integrate Kaluza for Gallios into your existing lab tools with standard FCS 3.1 files.
- Use offline software to create protocols and worklists to make your instrument time more productive.
- Use Gallios Simulator to train on the use of the instrument so you don't have to run a live experiment every time you train.
- Cleanse on Close means that the cytometer will be ready for action the next morning.
- Have a Gallios? No problem—Kaluza for Gallios will read Gallios protocols.
- Join the Kaluza for Gallios community: User forum to share stories and ask questions.



**Figure 1.** Visual management, drag-and-drop capabilities and linking protocols make configuring a carousel a snap.



**Figure 2.** Change voltage and compensation with easy-to-use sliders.