



# Kaluza C Analysis Software

PROTECTS YOUR DATA AND YOUR REPUTATION.



**VISUALIZE**  
*the possibilities.*



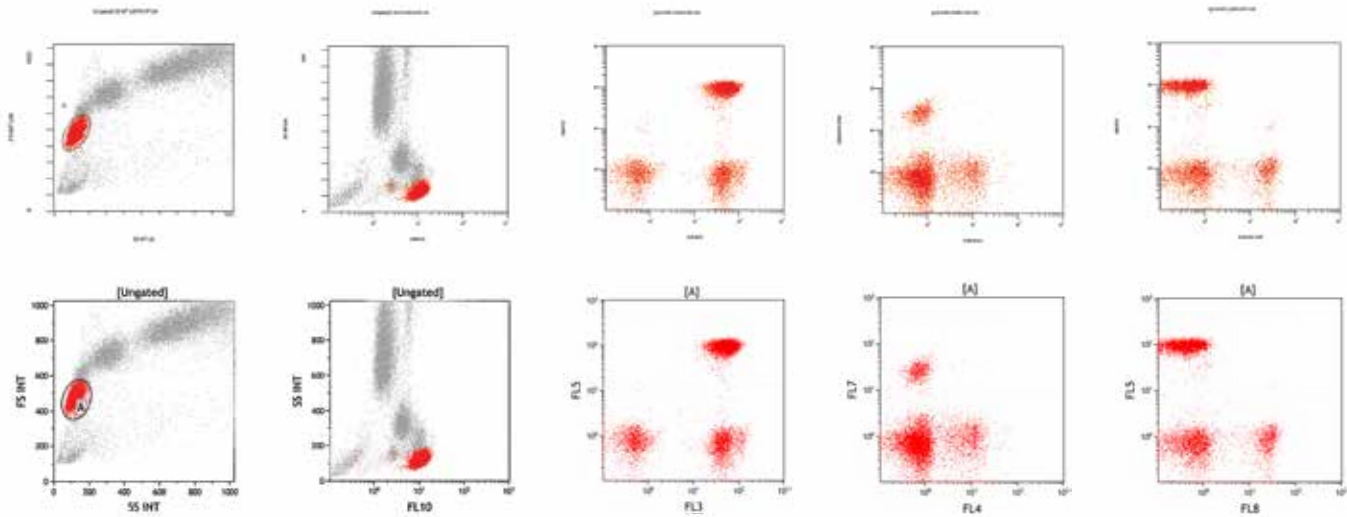
# KALUZA C ANALYSIS SOFTWARE

Kaluza C Flow Cytometry Analysis Software is built upon our successful Kaluza Analysis research use platform. We've turned our attention now to the unique needs and challenges of the clinical flow cytometry laboratory.

Clinical laboratories are under increasing pressure from three sources: reimbursement, regulatory, and resource. Meeting regulatory requirements places demands on both time and expertise. Kaluza C provides features that streamline the clinical QC reporting requirements and address standardization issues in flow cytometry. At the same time it provides the power to process multi-color files of up to 20 million events in real time.

- Listed with the FDA.
- Integration with Laboratory Information Systems (LIS).
- Traceability through user management.
- Specialized QC Report generation functions.
- Loads any listmode file that is compliant with the FCS standard up through version 3.1.

## Analyze Any Flow Cytometry Standard (FCS) Compliant File



Kaluza C uses the FCS \$CYT keyword to identify the flow cytometer used in acquisition. Data scaling preferences for the instrument are set up once upon loading the file into Kaluza. Example shows Navios data using Navios System Software (top row) and after importing into Kaluza Analysis (bottom row). Scatter and fluorescence parameters scaling are identical.

Presented plots and values are for illustrative purposes only and are based on customer reported data. Some information has been modified to simplify the presentation. Individual results may vary.

## Licensing Options

Whether you are looking for an individual timed license or a large site license, we have the solution for you. Special pricing for academic institutions is available. An economical single user 1 year license is available for those that want to try our platform with lower risk. All other licenses are in perpetuity and include future upgrades.

	NON-ACADEMIC PART NUMBER	ACADEMIC PART NUMBER
Kaluza C Analysis Software 1 Year, 1-User (Single)	C10575	C10578
Kaluza C Analysis Software Perpetual, 1-User (Single)	C10574	C10579
Kaluza C Analysis Software, 5-User (Network)	C10577	C10581
Kaluza C Analysis Software, 10-User (Network)	C10576	C10580

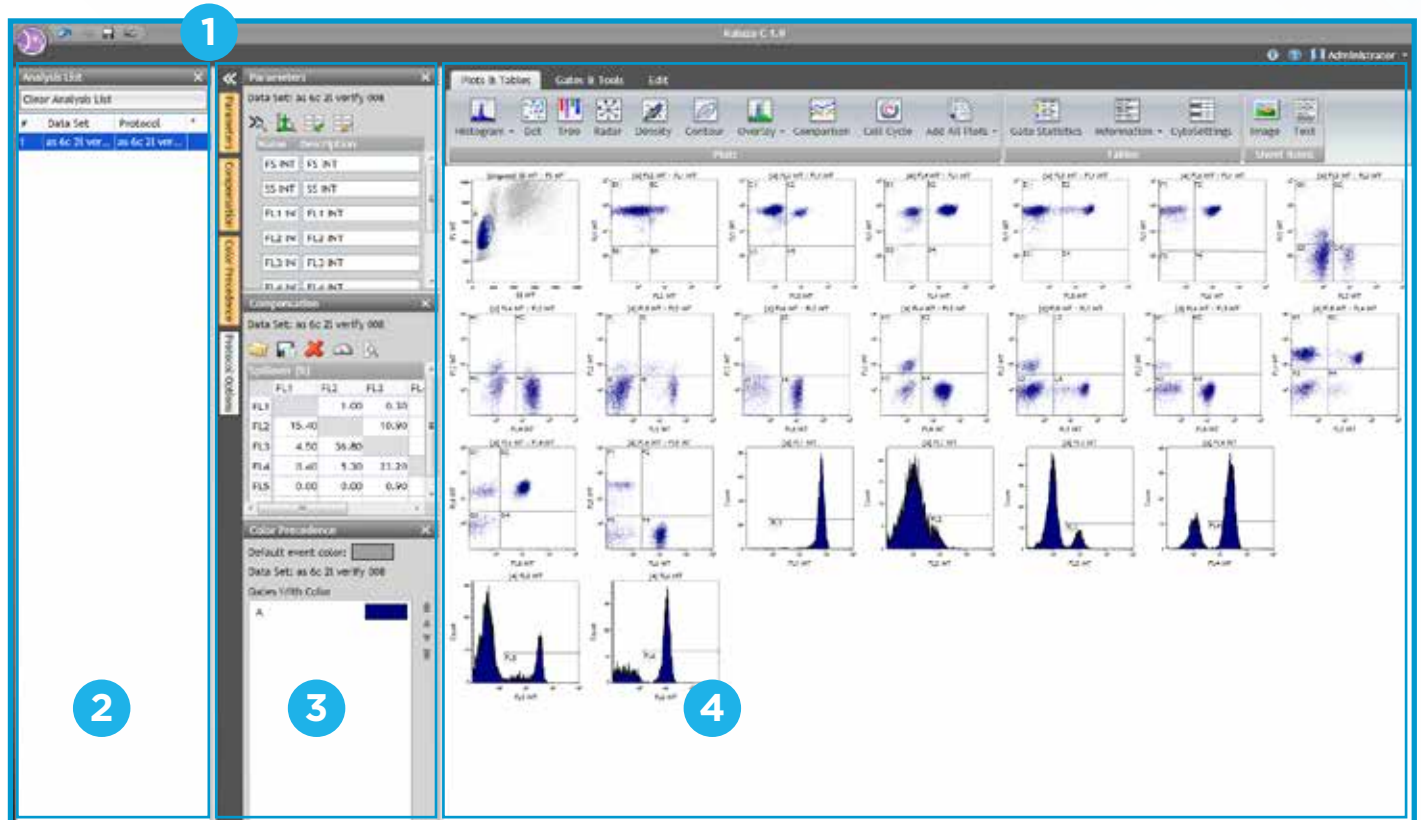
## Minimum Computer Requirements

	KALUZA C ANALYSIS SOFTWARE	NETWORK LICENSE CONTROL CENTER (NLCC)
Required processor	The processor must support the SSE2 Instruction set.	Quad-Core CPU, 3.2 GHz, above.
Required operating system	Windows 7, Enterprise Edition with Service Pack 1 (32 bit or 64 bit) Windows 8.1, Enterprise Edition (32 bit or 64 bit) Windows 10, Enterprise Edition (32 bit or 64 bit)	Windows 7, Enterprise Edition with Service Pack 1 (32 bit or 64 bit) Windows 8.1, Enterprise Edition (32 bit or 64 bit) Windows 10, Enterprise Edition (32 bit or 64 bit)
Required memory	200 MB of disk space to install For 32 bit systems, up to 4 GB of RAM For 64 bit systems, more than 4 GB of RAM	Memory minimum is 4 GB; 8 GB or above is recommended.
Required hard disk space	At least 1 G free space for storing data files.	The following requirements apply to Windows System Drive (c: drive, by default). <ul style="list-style-type: none"> <li>The NTFS format is required for good performance and security.</li> <li>A minimum of 50 GB of available hard disk space of the Windows System Drive.</li> </ul>
Required display	1440 x 900 (Min), Optimized for high-resolution wide screen monitors	No monitor required for the server; standard display for the client.

# SOPHISTICATION MADE SIMPLE

## Efficient Organized Interface

Four simple control panels is all you need to access every aspect of your information!



### 1 Tool Bar

Use the application button to access batch processing, QC report, plugin management, as well as open, save and export. The quick toolbar includes the unlimited undo and redo controls. And the info and help access detailed information including the complete instructions for use.

### 2 Analysis List

Keep track of all of your files, including data files (fcs or lmd), protocols, composite files, or analysis files. Drag and drop to open, reorder, or apply data to protocols. Highlight shows active file and asterisk indicates unsaved work.

### 3 Attributes Pane

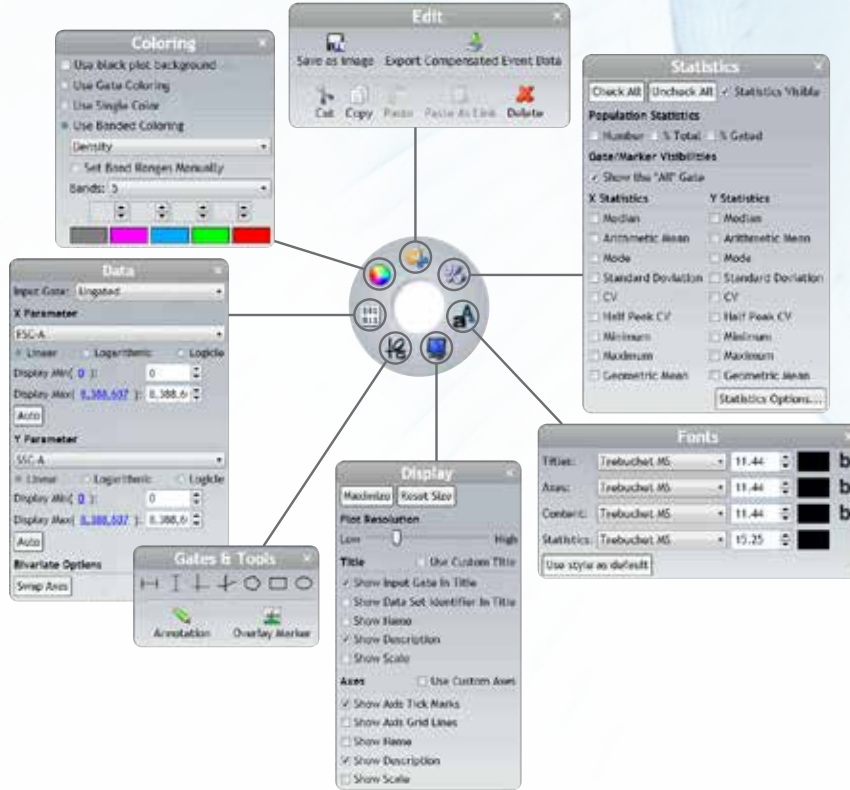
Quickly access to view and edit parameter info, compensation, color precedence, and protocol options. Collapse the panes to maximize analysis space.

### 4 Workspace

Use the ribbon to create plots and tables, add gates, including auto gate, annotations, and adjust compensation. Add additional sheets or reports as you explore your data. Drag and drop plots and use the zoom tool to manage plots viewed in the workspace.

## Context Specific Radial Menu

It took a lot of thought to create an analysis software this simple.  
At the heart of the Kaluza philosophy is the concept that what you need is just a right click away.



**Just Right Click:** Depending on the position of the pointer, different options are available in the radial menu. And the content of the submenu adjusts too.



Right click in the workspace, maybe you want to add an element.



Right click on a plot, maybe you want to add gates.



Right click on a gate, maybe you want to add Boolean logic.



Right click on an information table, maybe you want to adjust text attributes.

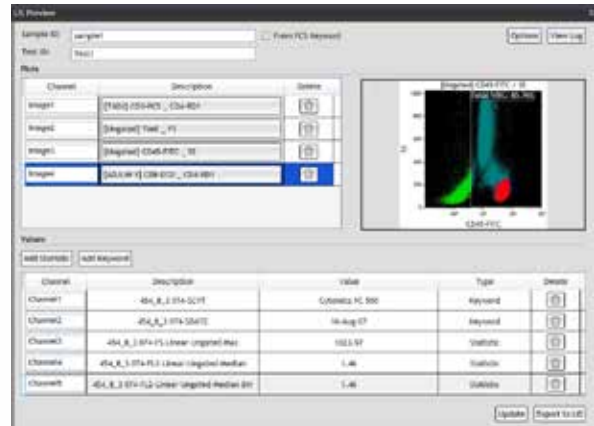
Compare the data table properties sub menu when accessed within the plot or from within a table radial menu. Attributes are different and vary based upon the context.

# PROTECTS YOUR DATA AND YOUR REPUTATION

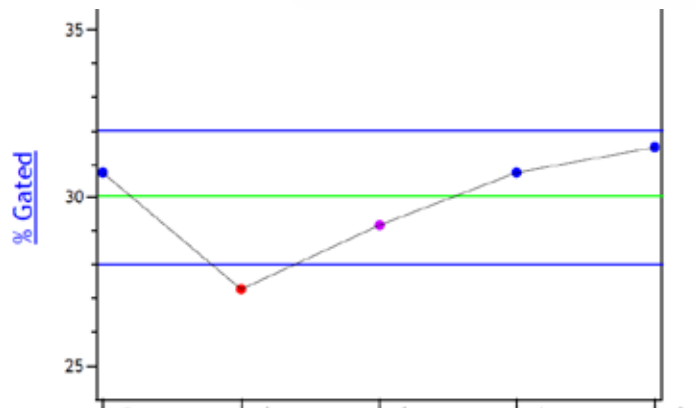
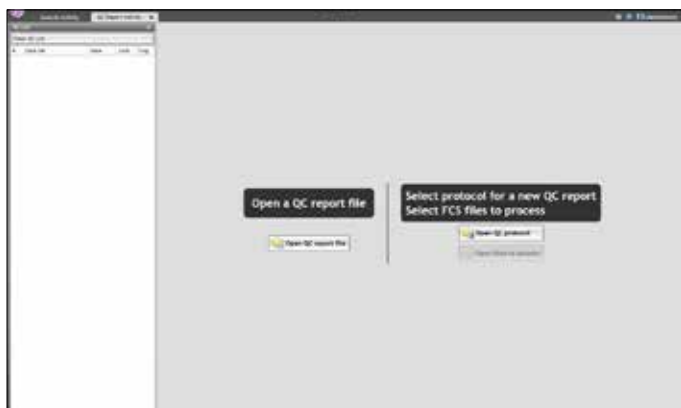
## Reports & LIS Connectivity

Create and edit a Report Sheet, which can include plots, tables, graphics, text, page numbers, and/or a date and time stamp. Quickly arrange plots so that they do not overlap, or align them to the location of your choice. An electronic approval signature can be added to the report sheet. Kaluza C includes an interface that enables the transmission of results, including plots, related statistics, and keywords, to a Laboratory Information System (LIS).



## Quality Control Reports

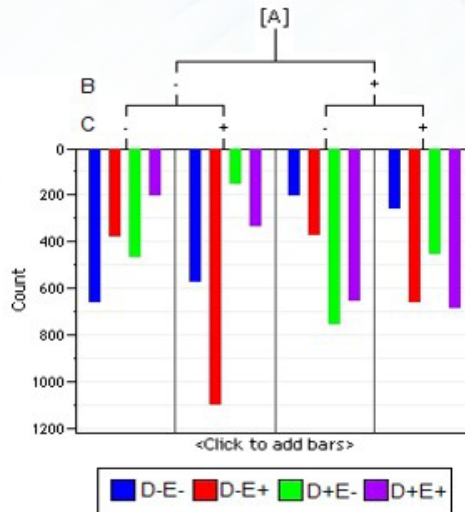
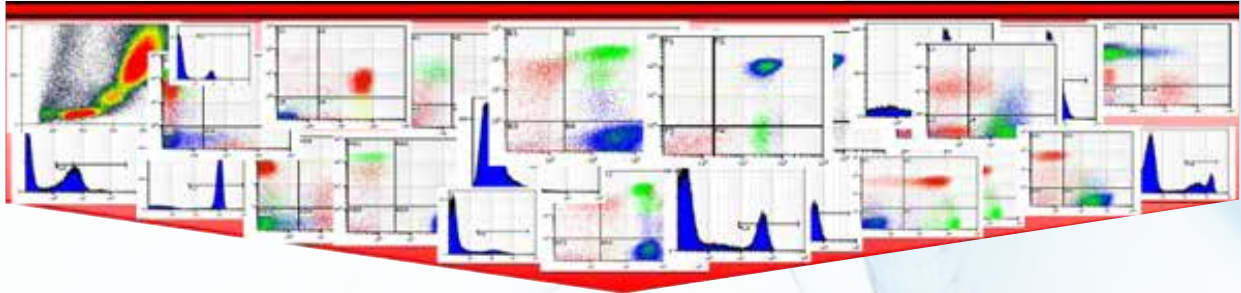
(QC Reports) are designed to help you detect issues that could produce inaccurate results in your data analyses, such as problems with the flow cytometer optics, instrument calibration, reagents, etc. Through the QC Report Activity tab, Kaluza C provides a platform to allow you to create Levey-Jennings plots, which clearly show whether data collected from control samples are either inside or outside of the standard accepted values. A QC Report is made up of one QC Protocol (\*.protocol or \*.qcp) and multiple Data Sets, which can be saved as one QC Report file (\*.qcr).



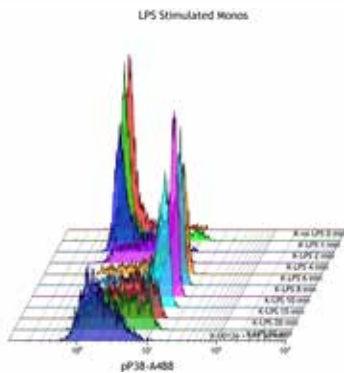
## User Management

User Management controls the level of access of each account set up on the system. There are two types of accounts available in Kaluza C, the Operator and the Administrator accounts. The Administrator account oversees account setup functions and can review the activities for all accounts on the system.

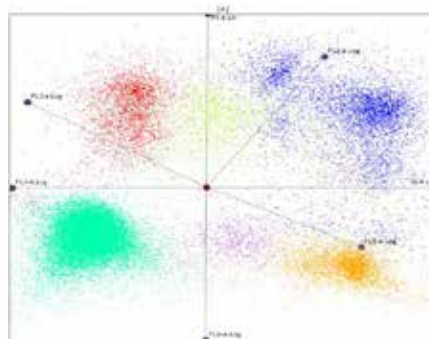
# ANALYSIS TOOLS



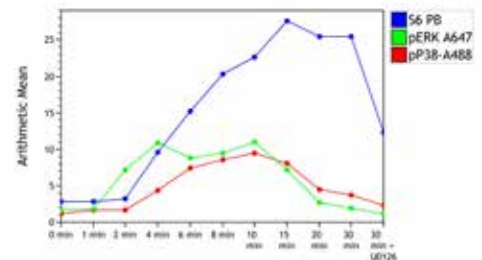
The tree plot can easily represent what typically requires dozens of plots, gates and regions do. Each phenotype can be used as a gate for further data exploration.



Overlay plots show data from multiple plots on a single pane. Density, dot and histograms can all be utilized in an overlay to easily view trends and changes.



The Radar Plot provides a multidimensional view. Clusters appear revealing relationships or rare populations not always apparent using bivariate plots and gating strategies. Animated radar plots can be exported as GIF files.



The comparison plot puts data from multiple plots into a single graph. Use it to view trends over time. Compare samples in a study to controls. Normalize samples or not, the choice is yours.



## Choose Beckman Coulter for Benchmark Expertise and Innovation

For over 80 years Beckman Coulter has driven innovation. We remain committed to shaping flow cytometry technology to fit seamlessly into your lab's workflow and to provide an optimal user experience. When you choose a Beckman Coulter solution you receive the a high level of expertise, innovation, and quality assurance.

Contact your local Beckman Coulter sales representative.



For In Vitro Diagnostic Use. CE marked to the European In Vitro Diagnostic Medical Device Directive 98/79/EC.

© 2017 Beckman Coulter, Inc. All rights reserved. Beckman Coulter, the stylized logo, and the Beckman Coulter product and service marks used herein are trademarks or registered trademarks of Beckman Coulter, Inc. in the United States and other countries.

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

FLOW-3104SB11.17