

Explore new dimensions in flow cytometry analysis

Kaluza analysis software

Genomics

Cell Analysis

Particle Characterization

Capillary Electrophoresis

Lab Automation

Centrifugation

Bioseparation

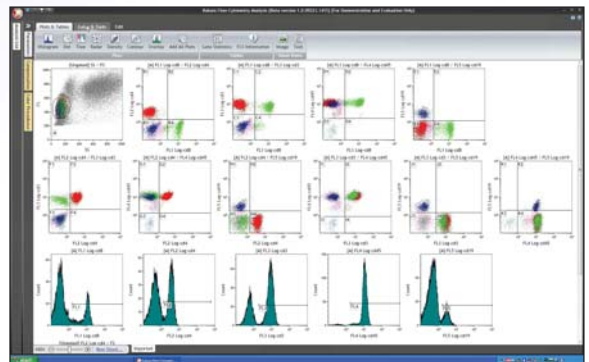
Lab Tools

Today, more colors require more plots, events, protocol complexity and ultimately increased analysis time. Kaluza software challenges today's solutions with a new vision of what flow cytometry analysis should be. Designed to simply, efficiently and quickly analyze multicolor, multiparametric data, Kaluza proficiently manages your analysis from the cytometers of today and tomorrow.

Intelligent Workspace Design

Kaluza's innovative dynamic workspace allows flexibility and freedom of organizing displays with minimal effort providing an intuitive environment for data exploration.

- Analyze directly from report or plot sheet
- Drag & drop plots into a report page
- Multiple report layouts per sample



Data replay is made quick and easy with automatic loading of the original plots, gates, and colors from files with embedded analysis protocols.

Radial Menu

Spend less time "hunting" for options and more time exploring your data. Kaluza's novel use of a radial menu provides one-click access to all plot and region editing eliminating the need to search for features.

On-plot access to

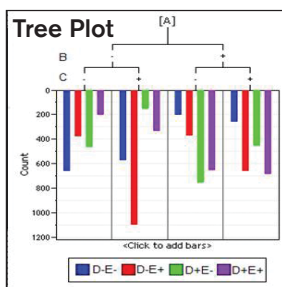
- Creating gates
- Formatting
- Annotating
- Statistics



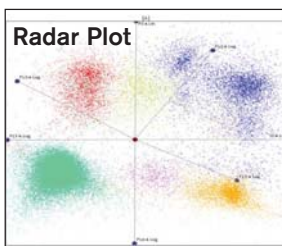
Multiparametric Plots

Kaluza's innovative multiparametric plots revolutionize the way multicolor data is analyzed.

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.



The radar plot provides multidimensional view of data without drawing a single gate. Clusters easily appear revealing relevant and rare population typically not seen using traditional bivariate plots and gating strategies

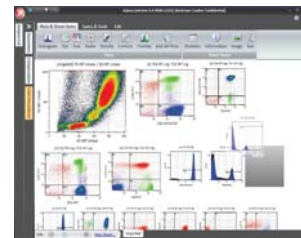


- Simplifies multiparametric data analysis into interactive display
- Enables quick pattern recognition analysis
- Analyze all events in a single view

Plot Management

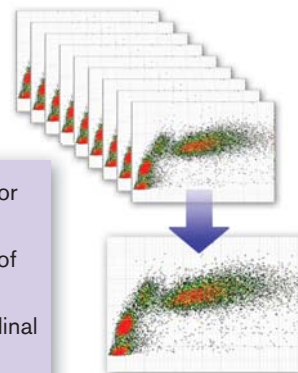
Kaluza's unique Auto-Layout design intelligently reconfigures your workspace for you. Minimizes the clicks required to create and maintain your workspace to save you time and frustration.

- Custom plot sizing
- Easy creating of plots
- Zoom in and out
- Guided drag & drop



Data Merge

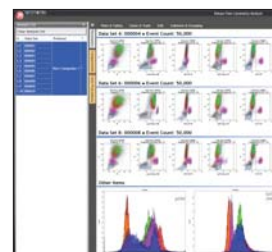
The data merge feature enables better measurement of rare populations and facilitates analysis of longitudinal and time course studies



- Easily combine multiple files for quality analysis
- Enables better measurement of rare populations
- Facilitates analysis of longitudinal and time course studies

Multi-File Analysis

Analysis of multiple datasets is as simple as a drag & drop into the intuitive analysis list. Composite analysis allows you to quickly and easily change compensation and create overlays with multiple files.



Ordering Information

Kit description	Part Number
Kaluza Software Kit Single User	A82959
Kaluza Software Kit 3 User Bundle	A85788
Kaluza Software Kit 5 User Bundle	A82960
Kaluza Software Kit 5 User Network	A84175
Kaluza Software Kit 10 User Network	A82961
Kaluza Software Kit 1 year Single User renewable license	Call for info

Space for
FSC Logo

Printed with soy ink

Kaluza is for Research Use Only. Not for use in diagnostic procedures.

Beckman Coulter and the stylized logo are registered trademarks of Beckman Coulter, Inc. Kaluza is a trademark of Beckman Coulter, Inc.

Australia, Gladesville (61) 2 9844-6000 Austria, Vienna (43) 1 291 01 241 Canada, Mississauga (1) 905 819 1234
 China, Shanghai (86) 21 3865 1000 Croatia, Zagreb (38) 51 489 9003 Czech Republic, Prague (420) 272 01 73 32
 Eastern Europe, Middle East, North Africa, South West Asia: Switzerland, Nyon (41) 22 365 3707 France, Villepinte (33) 1 49 90 90 00
 Germany, Krefeld (49) 2151 33 35 Hong Kong (852) 2814 7431 India, Mumbai (91) 22 3080 5000 Italy, Cassina de' Pecchi, Milan (39) 02 953921
 Japan, Tokyo (81) 3 5530 8500 Korea, Seoul (82) 2 404 2146 Latin America (1) (305) 380 4709 Malaysia, Kuala Lumpur (60) 3 5621 4793
 Mexico, Mexico City (001) 52 55 9183 2800 Netherlands, Woerden (31) 348 462462 Poland, Warszawa (48) 22 366 0180 Puerto Rico (787) 747 3335
 Singapore (65) 6339 3633 South Africa/Sub-Saharan Africa, Johannesburg (27) 11 564 3203 Spain, Madrid (34) 91 3836080
 Sweden, Bromma (46) 8 564 85 900 Switzerland, Nyon (41) 0800 850 810 Taiwan, Taipei (886) 2 2730 2500 Turkey, Istanbul (90) 216 570 17 17
 UK, High Wycombe (44) 01494 441181 USA, Brea, CA (1) 800 742 2345