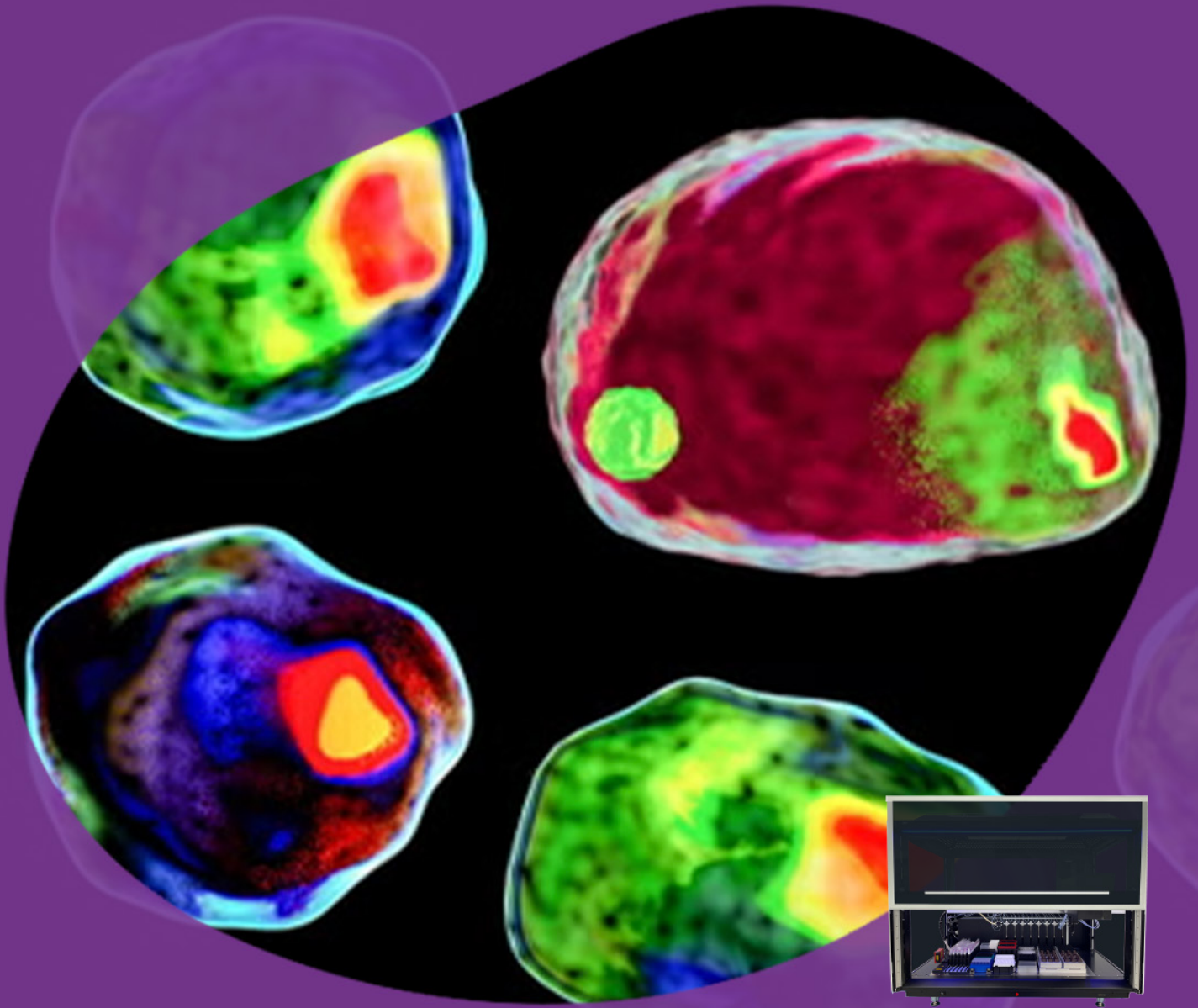


Scientist-driven flow
cytometry solutions
supporting data quality
and reliability.



revvity

Flow cytometry reagents,
software, and automation

Advancing your research with automated cocktail and sample preparation, combined with a wide-ranging antibody catalog.

Our knowledge and expertise with fluorophores, chemistries, and antibody conjugation have helped us become a preeminent expert on flow cytometry. Our scientists are crafting dyes that ignite discoveries by emitting in the far red, filling previously unoccupied spectral spaces, and providing enhanced stability and/or brightness.

Antibodies, fluorophores, and buffers

Explore our fluorophore options below to see which one best fits your panel.

Fluorophores

Fluorophore	Excitation Max (nm)	Emission Max (nm)	Filter	Brightness
Ultraviolet laser (355 nm)				
Spark PLUS UV395™	355	385	379/28	4
Spark UV™ 387	355	388		2
StarBright™ UltraViolet 575	340	569	586/15	4
StarBright™ UltraViolet 605	340	609	610/20	5
StarBright™ UltraViolet 665	340	669	670/30	5
StarBright™ UltraViolet 740	343	737	740/35	5
StarBright™ UltraViolet 795	340	792	820/60	5
Violet laser (405 nm)				
Spark Violet™ 423	405	423	450/50	3
BV421™	405	421	450/50	5
Pacific Blue™	410	455	450/50	1
Spark PLUS V450™	405	452	450/50	4
Spark PLUS V475™	404	475	450/50 or 460/22	4
Spark Violet™ 500	393	500	510/50	1
BV510™	405	510	510/50	2
Spark Violet™ 538	399	538		1
BV570™	405	570	585/42	2
BV605™	405	603	610/20	4
BV650™	405	645	660/20	4
BV711™	405	711	710/50	5
BV750™	405	750	780/60	3
BV785™	405	785	780/60	4

StarBright™ Dyes are powered by Bio-Rad.

Fluorophore	Excitation Max (nm)	Emission Max (nm)	Filter	Brightness
Blue laser (488 nm)				
Spark PLUS B488™	488	520	530/30	5
KIRAVIA Blue 520™	488	520	530/30	4
Alexa Fluor® 488	495	519	530/30	3
FITC	493	525	530/30	3
Spark Blue™ 515	506	511	530/30	2
Spark PLUS B550™	516	540	530/30	3
Spark Blue™ 550	516	540	530/30	1
Spark PLUS B574™	506	574	570/40	4
Spark Blue™ 574	506	574	570/40	1
PerCP	482	675	695/40	1
PerCP/Cyanine5.5	482	690	695/40	2
PerCP/Fire™ 780	482	774	780/60	4
PerCP/Fire™ 806	482	802	780/60	4
Yellow/Green laser (532, 561 nm)				
PE	565	575	585/20	5
Spark PLUS YG581™	565	581	575/26 or 585/15	4
Spark YG™ 581	565	581	575/26 or 585/15	2
Spark YG™ 593	565	593		3
PE/Dazzle™ 594	565	610	610/20	5
PE/Fire™ 640	565	639		4
PE/Cyanine5	565	670	660/20	5
PE/Fire™ 700	565	695	710/50	5
PE/Fire™ 744	565	743	780/60	5
PE/Cyanine7	565	774	780/60	4
PE/Fire™ 810	565	806	780/60	4

Fluorophore	Excitation Max (nm)	Emission Max (nm)	Filter	Brightness
Red laser (633 nm)				
APC	650	660	660/20	4
Alexa Fluor® 647	650	668	660/20	4
Spark NIR™ 685	660	685		3
Spark Red™ 718	697	711	720/45	4

Fluorophore	Excitation Max (nm)	Emission Max (nm)	Filter	Brightness
Alexa Fluor® 700	696	719	720/45	2
APC/Cyanine7	650	774	780/60	2
APC/Fire™ 750	650	774	780/60	2
APC/Fire™ 810	787	807		3

Antibodies

Antibodies and targets at glance

Our catalog features over 22,000 antibodies and kits for flow cytometry applications. We supply Research Use Only (RUO) products, Analyte Specific Reagents (ASR) products, and products manufactured according to GMP guidelines. Adherence to quality is supported by our ISO 13485:2016 and MDSAP certifications.

Surface markers are crucial for phenotyping in flow cytometry, as combinations of these markers identify cell populations. For example, mature T cells are identified by TCR, CD3, and either CD4 or CD8. For other cell types, other important surface markers include TRBC1, kappa and lambda light chains, and OX-40.

Intracellular targets include proteins like cytokines, chemokines, and growth factors, some of which require protein transport inhibitors for detection. Key intracellular markers are MPO and Ki-67, which are vital for signaling pathways and protein regulation.

Nuclear targets reveal cell fate information through transcription factors and genetic content. T helper cells, for instance, are characterized by transcription factors like FOXP3, T-bet, and GATA3. Histones are also assessed for epigenetic modifications.

Reagents

Sample preparation solutions	
Buffer	Function
True-Stain™ Multi-Fluor Buffer	Prevents interactions between fluorescent reagents when staining cells with multiple BV or BUV antibody conjugates without reducing their signal.
TruStain FcX™	Fc receptor blocking solution that prevents non-specific binding of antibodies to cells.
Lymphopure™	Density gradient media for mononuclear cell (PBMC) isolation.
RBC Lysis Buffer (10x)	Concentrated buffer for red blood cell lysis. Contains ammonium chloride, potassium carbonate, and EDTA.
RBC Lysis/Fixation Solution (10x)	Concentrated buffer for red blood cell lysis and fixation of remaining leukocytes.
Fixation buffers	
Buffer	Function
Fixation Buffer	Paraformaldehyde-based buffer used to fix cells.
FluoroFix™ Buffer	Gentler fixation buffer with a lower percentage of paraformaldehyde specialized for use with tandem dyes.
Storage buffers	
Buffer	Function
Cyto-Last™ Buffer	Specially formulated buffer for prolonged storage of cytokine producing cells.



[Learn more](#)

Discover our flow cytometry reagents.

Enhance flow cytometry reliability with automated cocktail preparation

In the realm of flow cytometry, the reliable preparation of reagents is crucial for obtaining precise and consistent outcomes. Flow cytometry is a critical tool in research, facilitating in-depth analysis of cell populations by detecting surface and intracellular markers. Preparing the antibody cocktail reliably is crucial for determining the quality of the collected data.

Fontus liquid handler

Fontus™ workstation streamlines and enhances the antibody cocktail preparation and staining process by focusing on these core aspects. This leads to more reliable and reproducible results, better resource utilization, and improved operational efficiency in research.

- Barcode reader for sample traceability
- Cooling blocks for antibody vial temperature control
- Flexible deck size based on customer's throughput and workflow
- Tinted Plexiglas for light blocking
- Automated handling of antibody reagents and cocktail reagent preparation
- Sub 5% CVs and 1 μ L to 5 mL sample working volumes
- Minimal dead volume with liquid level detection
- High throughput with dual pipetting arms

Why automate?

- Provides standardization of flow cytometry sample preparation
- Improved operator safety through minimal sample interaction
- Reduced operator variability and potential risk of errors
- Open system allows for flexible laboratory testing protocols
- Versatile for routine and complex flow protocols
- Compatible with multiple manufacturers' antibody vials
- Enables highly skilled staff to concentrate on more complex operations



Precision

- Reliable dispensing technologies that allow for precise volumes
- Minimize dead volume
- Reduce human error and variability in measurement

Standardization

- Consistency in preparation process
- Enhance result reproducibility

Enhance productivity

- Increase walkway time
- Focus on value-added activities
- Higher throughput

Traceability

- Automatic audit trail

Streamlined data capture

Flexibility

- Software and deck customizable
- Deck QC identifies discrepancies pre-run
- Optional HEPA filter

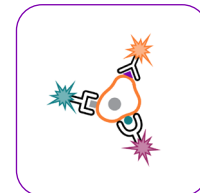


[Learn more](#)

Fontus automated
workstation

BioLegend from Revvity's flow cytometry data analysis solution

Experience our cloud-based, innovative, and easy-to-use flow cytometry data analysis software. Move beyond single-computer access, plug-ins, and installation headaches – use the cloud to access, analyze, and manage your data from anywhere, at any time. Experience our fully featured and highly interactive user interface equipped with all the standard features you'd expect, including built-in high-dimensional analysis methods and AI-powered tools.



Why CytoScribe?



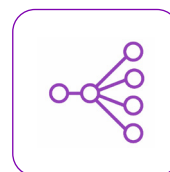
Fully cloud-based and accessible anywhere

- No installation required, runs directly in a web browser
- Accessible from any standard computer
- No hardware upgrades and software maintenance needed
- Enables continuous innovation



Interactive and intuitive user experience

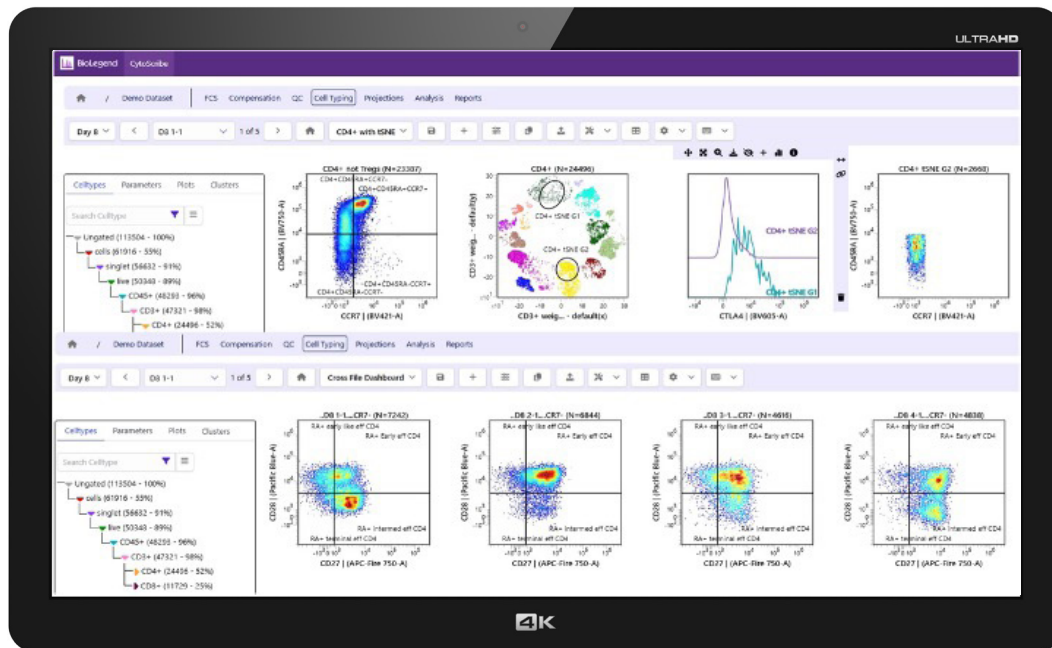
- Live, interactive single-cell data visualization
- Easy-to-use interface for beginners and experts
- Save and reuse gating strategies and layouts
- Synchronize gating across multiple samples



Advanced analytics made accessible

- High-dimensional analysis tools:
 - Clustering (PARC, Leiden, FlowSOM)
 - Dimensionality reduction (UMAP, t-SNE, weighted t-SNE)
- GPU-powered computation for performance
- Simplifies complex data exploration for bench scientists

Interactive cell typing interface



Researcher collaboration and data sharing made easy

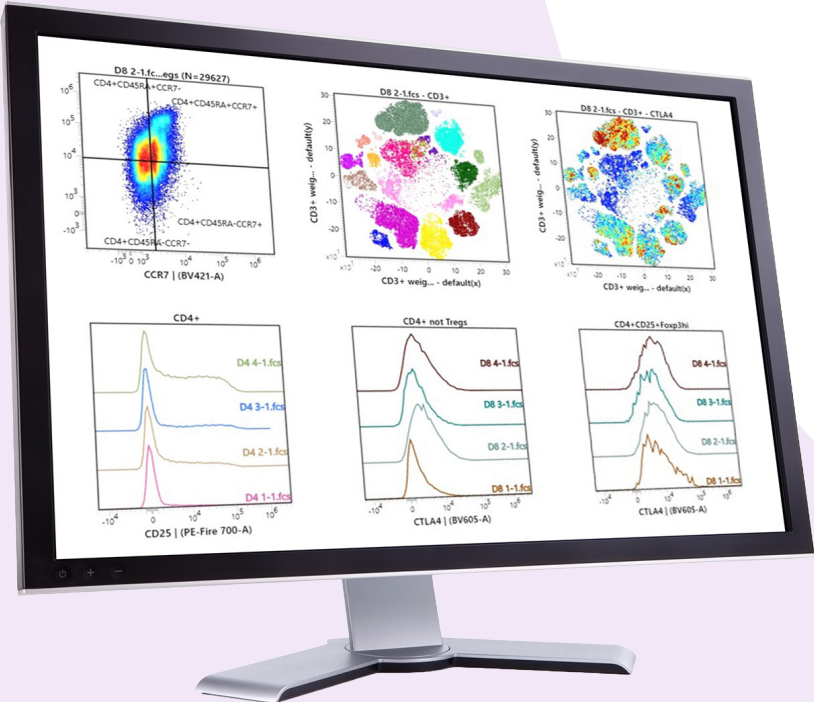
End-to-end flow cytometry workflow support

- Supports the entire analysis workflow, from QC to reporting
- Built-in tools for:
 - Data QC (e.g., EMDcut, PeacoQC)
 - Compensation & spectral unmixing
 - Gating hierarchies and statistics
 - Dimensionality reductions
- Export-ready outputs (PowerPoint images, Excel tables)

CytoScribe is scalable to your research needs

Options for site-wide licenses and custom storage configurations make it perfect for labs of all sizes.

Feature	Free	Premium	Enterprise
Standard cell analysis tool	✓	✓	✓
Dimensionality reductions	✓	✓	✓
Share with support	✓	✓	✓
Collaboration (team sharing)		✓	✓
Experiment design tool		✓	✓
CytolQ: AI gating		✓	✓
Custom AI models			✓
Dedicated server			✓
Cloud storage	2 GB	50 GB (scalable)	Customized



Learn more

Get your free account

Precision where it matters: supporting your flow cytometry workflow with Revvity solutions

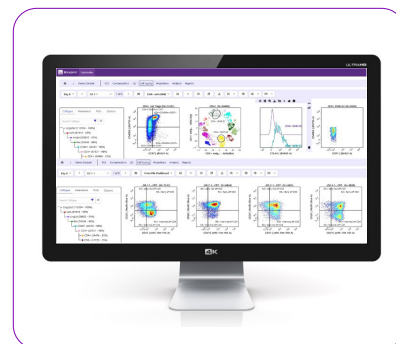
Explore Revvity's fully integrated approach, combining automation, reagents, and data solutions to support your end-to-end flow cytometry workflow.



Fontus™ liquid handler



BioLegend reagents



BioLegend CytoScribe

Precision and standardization: achieve reproducible results in your research workflows with our automated liquid handlers like the Fontus systems for reliable and precise reagent preparation while minimizing variability and errors.

Enhanced productivity: Free up your valuable time for critical analysis by automating cocktail prep and data workflows, allowing efficient high-throughput research operations.

Comprehensive reagent support: Take advantage of BioLegend's comprehensive reagent catalog, including antibodies, fluorophores, and buffers tailored for advanced flow cytometry applications.

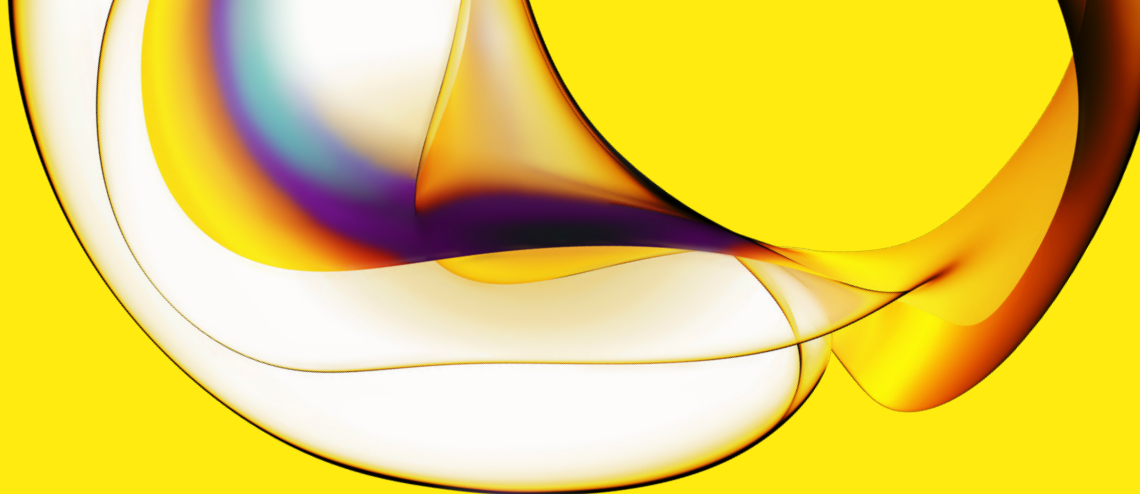
Intuitive data analysis with CytoScribe: A cloud-based, user-friendly platform offering powerful AI-driven tools to simplify high-dimensional data visualization, collaboration, and reporting.

Flexibility for every workflow: Open systems and scalable solutions designed to support the specific needs of your lab, from small-scale studies to enterprise-level operations.



Learn more

Visit the
BioLegend website



www.revivity.com

revvity

Copyright ©2026, Revvity, Inc. All rights reserved.
Revvity is a trademark of Revvity, Inc. All other trademarks are the property of their respective owners.

1843541

For research use only.
Not for use in diagnostic procedures.