



Clinical flow cytometry solutions

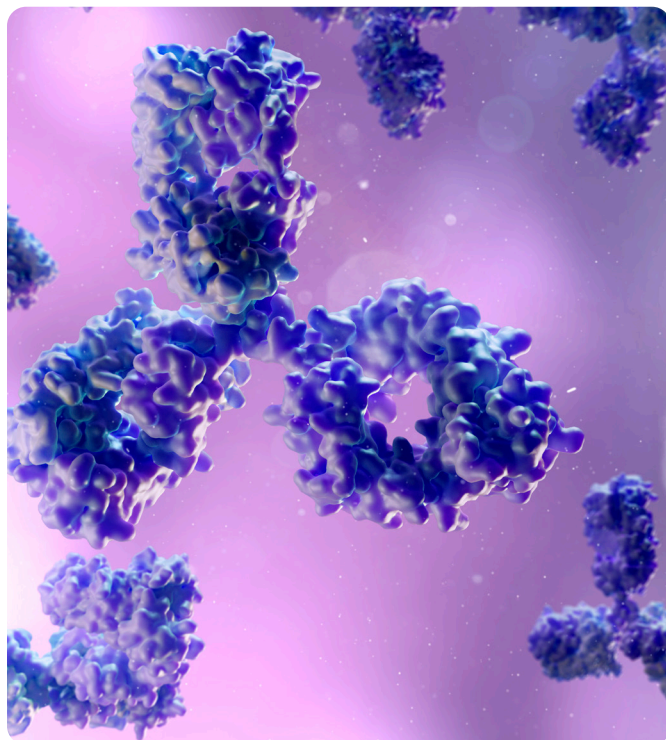
Combined power of automated cocktail preparation and clinical reagents.

Single color antibodies for clinical applications - Analyte specific reagents* (ASR)

Wide range of high-quality ASR antibodies:

- 300+ antibodies for flow cytometry applications
- Premium antibodies ensuring dependable results and precise reporting

BioLegend ASR products are manufactured in accordance with US 21 CFR Part 820 in an ISO 13485:2016 and MDSAP-certified facility following GMP compliant procedures to ensure the highest quality standards.



* Analytical and performance characteristics of these reagents are not established.

Single color ASR antibodies for clinical applications

	405				488							633	
	Pacific Blue®	Spark Violet™ 423	Spark Violet™ 500	Spark Violet™ 538	FITC	PE	PE/Cyanine5	PE/Cyanine7	PE/Dazzle™ 594	PerCP	PerCP/Cyanine5.5	APC	APC/Fire™ 750
CD1a						•						•	
CD2	•				•	•		•			•	•	•
CD3	•		•		•	•		•	•	•	•	•	•
CD4	•	•			•	•		•	•	•	•	•	•
CD5	•				•	•		•			•	•	•
CD7	•				•	•		•	•			•	
CD8	•				•	•		•		•	•	•	•
CD9					•						•	•	•
CD10					•	•		•	•		•	•	•
CD11b	•				•	•		•			•	•	•
CD11c	•					•		•			•	•	
CD13						•		•	•		•	•	
CD14	•				•	•		•	•		•	•	•
CD15	•				•	•		•					
CD16	•				•	•		•	•		•	•	•
CD19	•	•			•	•	•	•			•	•	•
CD20	•				•	•		•			•	•	•
CD21													•
CD22											•		
CD23					•	•		•			•	•	
CD24						•					•		•
CD25					•	•		•			•	•	•
CD26					•	•		•					
CD27	•				•	•		•			•	•	•
CD28	•					•		•			•	•	
CD30					•	•						•	
CD31					•								
CD33					•	•	•	•			•	•	
CD36					•								•
CD38	•				•	•		•	•		•	•	•
CD41					•	•						•	
CD42b						•							
CD45	•			•	•	•		•	•	•	•	•	•
CD45RA					•			•	•			•	•
CD45RO					•	•						•	
CD49d													•
CD56 (NCAM)	•				•	•		•	•		•	•	•
CD58						•							
CD59					•								
CD61					•	•							
CD62L	•				•	•			•			•	
CD62P					•	•							
CD64	•				•	•		•	•		•		•
CD66b					•								
CD69						•						•	
CD71					•			•				•	•
CD79a					•	•						•	
CD79b						•						•	•
CD81	•				•								
CD94					•							•	
CD95						•							
CD99					•	•							
CD103					•	•						•	
CD117						•		•	•		•	•	•
CD123	•					•		•		•	•	•	
CD127						•		•			•		•
CD134 (OX40)						•							
CD138					•								
CD159a								•					
CD159c						•							
CD154								•					
CD183						•							
CD185						•						•	
CD197(CCR7)						•							
CD200						•		•				•	
CD203c						•		•					
CD223												•	
CD279(PD-1)						•							
CD303												•	
CD304	•					•							
CD309												•	
CD335 (NKp46)								•					
Cytochrome c					•								
GranzymeA						•							
HLA-DR	•	•			•	•		•	•	•	•	•	•
IFN-gamma						•						•	
IgD						•							
IgG1-IsoCtr	•				•	•		•	•		•	•	•
IgG2a-IsoCtr	•				•	•		•	•		•	•	•
IgG2b-IsoCtr	•				•			•			•	•	
IgLC-Kappa	•				•	•						•	
IgLC-Lambda					•	•						•	•
IgM-μ					•	•						•	
IgM-IsoCtr					•								
IL-4						•							
TCRalpha-beta					•	•						•	
TNF-a						•							

Precision at every step

Automated flow cytometry reagent cocktail preparation

Streamline your laboratory's workflow with cutting-edge technology

In the realm of flow cytometry, the accurate preparation of reagents and cocktails is crucial for obtaining precise and consistent outcomes, determining the quality of the collected data.

Precision

- Reliable dispensing technologies that enable 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

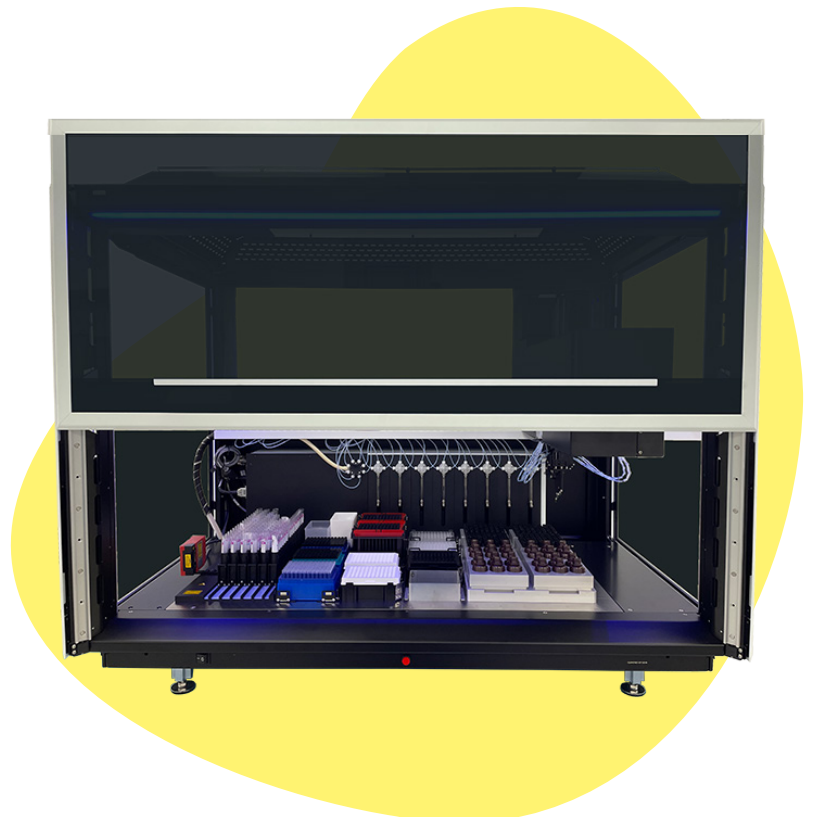
Traceability

- Automatic audit trail
- Streamlined data capture

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 and clinical laboratories.

- 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 µL to 5 mL sample working volumes
- Minimal dead volume with liquid level detection
- High throughput with dual pipetting arms

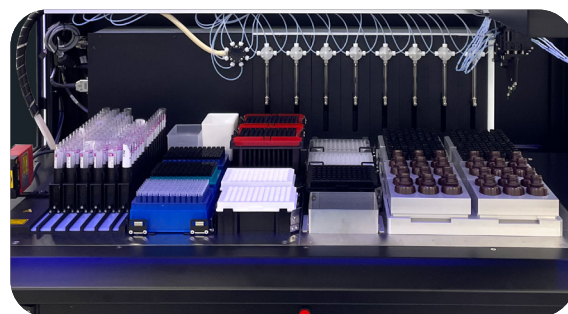


Revvity's solution: precision in automated liquid handling meets high reagent quality

Combine the power of Fontus automation with BioLegend's antibody reagents to streamline your workflow, increase throughput, and achieve consistency in your flow cytometry assay.

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

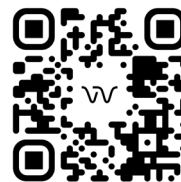


Why BioLegend flow cytometry reagents?

- Wide range of high-quality antibodies
- ASR antibodies are manufactured in an ISO 13485:2016 and MDSAP-certified facility
- Custom Services:
 - Custom conjugations offered
 - Custom antibody cocktails
- Extensive selection of fluorophores to choose from.



For more information, contact
your local representative or
visit our website:



Fontus cytometry
cocktail prep workstation



BioLegend ASR
antibodies



revvity