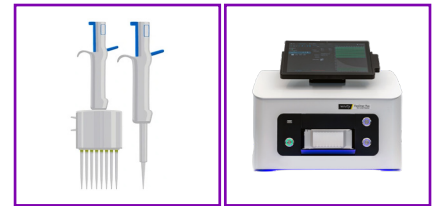




Optimizing antibody dispensing for flow cytometry.



Maximize speed, minimize waste

Introducing FlexDrop Plus

FlexDrop™ Plus is built for precision and speed. Its advanced automation allows for the rapid dispensing of multiple antibodies and reagents, eliminating the bottlenecks and complexity caused by manual pipetting. With its ability to dispense reagents in nanoliter increments, FlexDrop Plus ensures that even the smallest volumes are handled with unmatched accuracy - allowing you to prepare complex cocktails faster and with less waste of expensive antibodies.

How FlexDrop Plus works:

Configure cocktail mix

Select the reagent volumes and types through the simple touchscreen interface or import a CSV file.

Automate dispensing

FlexDrop Plus handles the complex mixing, delivering precise volumes to prepare each cocktail.

Monitor and optimize

Keep track of the entire process with real-time monitoring and post-run report to ensure all dispensing was done correctly using the unique droplet detection of the FlexDrop Plus.

	Manual	FlexDrop Plus
Technology	Pipette	Non-contact nL dispenser
Dispenser	1, 8, 12 fixed channels	1 to 8 flexible channels
Source material	Variable	FlexDrop Source plates
Source capacity	Non-limiting	80 or 500 µL per well
Dead volume	Variable, down to 0 µL	<1 µL
Verification	No	Droplet verification
Precision	100 nL, typical	8 nL
Time*	3 hrs	20 min
Automation compatible	No	Yes

* FlexDrop flow cytometry app note_1568639 APP_v5 - Automated antibody dispensing to improve high-parameter flow cytometry throughput and analysis. (Bosteels et al, 2024).

Optimize your lab efficiency

Save time: Reduce manual pipetting, focus on research.

Save money: Minimize antibody waste, lower costs.

Increase accuracy: Achieve consistent results with precise cocktail preparation.

Key features

Unparalleled accuracy

Allows the dispensing of complex cocktails with multi-reagent transfers as small as nanoliters.

Minimize dead volume

Designed to minimize dead volume of <1 μL , conserving use and cost of your valuable antibodies.

Automated multi-reagent dispensing

Eliminates manual pipetting, enabling faster, more reliable dispensing of up to 8 reagents at a time.

User-friendly interface

Intuitive touchscreen interface allowing quick setup of custom reagent mixes or input CSV files.

Flexible workflow integration

Easily integrated into your existing workflows, as a standalone device or can be integrated with automation.



FlexDrop Plus

- Non-contact dispensing - eliminate carryover and cross-contamination
- Dead volume < 1 μL - conserve valuable reagents and expensive antibodies
- 8 nL or larger dispense volumes and 80 or 500 μL source plate options
- Fast - dispense 10 nL into 96 wells in 10 seconds, into 384 wells in 20 seconds
- Smart - Simultaneously dispense from 1 to 8 wells (a column) to 96, 384 or 1536 well plates
- Built-in droplet verification - verify liquid dispensed

BioLegend reagents

- Wide range of high-quality antibodies
 - Fluorophore conjugated antibodies
 - TotalSeq Antibodies
- Extensive selection of fluorophores to choose from
- Custom conjugations offered
- Custom bottle size option available



Ready to revolutionize your workflow?
Contact us today.

revvity