

# Pipeline™ Dispenser and Peripette™ Consumable



Pipeline™ Dispenser

- Accurate and precise.
- Promotes consistent technique.
- Increases throughput.
- Reduces tiring, repetitive motions.
- Foot switch operation.
- Designed for use in bio-cabinets.

The Pipeline™ dispenser and the Peripette™ consumable provide a fast, accurate and efficient approach to the transfer of sterile fluids. By combining the advantages of a peristaltic pump with the convenience of a sterile consumable, the Pipeline offers an ideal solution for high-throughput research applications.

- Materials meet USP Class VI Gamma-radiation sterilized.
- Non-pyrogenic.
- Compatible with dispensing cells.
- Packaged in easy-to-open pouches.
- Loads quickly into Pipeline.



Peripette™ Consumable

## One-step operation

The Pipeline dispenser eliminates the tedious nature of hand-operated pipetting for large-scale cell culture:

- Three times as fast as manual pipetting.
- Reduces repetitive motion injuries.
- Improved precision and accuracy vs. manual pipetting.

To operate, the user simply snaps in a sterile Peripette (Peristaltic Pipette), lifts the pump head mechanism, inserts the source bottle and lowers the assembly. After setting the desired volume, the user places the destination container under the output spout and initiates the dispenser. Pipeline then performs the fluid transfer in a one-step operation. This saves time and reduces the stirring, repetitive motions (aspirate, move and dispense) as required for hand pipetting. Additionally, the dispenser is ready to repeat the process, as many times as is needed, until the source bottle is empty.

With increased efficiency and high throughput capabilities, Pipeline facilitates repetitive laboratory tasks. For example, Pipeline provides a fast and accurate means for preparing

multiple aliquots of a common reagent. Pipeline is also compatible with dispensing cells and suitable for use in cell culture. In multi-step procedures, one can use an aspirator along with the Pipeline to exchange cell culture media, solutions and reagents without the need to change multiple pipettes. Overall, the Pipeline dispenser enhances laboratory protocols by promoting consistent technique.

## Flexibility

The Pipeline is capable of dispensing different modes:

- A user-defined volume of fluid per switch activation
- Dispensing continuously during switch activation
- Three different dispense speeds tailored to the most demanding cell culture applications

The user can also save and recall up to 8 pre-programmed settings, and the unit can be activated via front-panel switch or foot switch. The accuracy and precision of the dispensed volumes are similar to serological pipettes (see specifications). The Pipeline also provides an RS-232 interface if automation or remote operation is required.

## Convenience

The design of the Pipeline dispenser optimizes the space usage inside the bio-cabinet, requiring a base footprint of less than 7" x 10". When not in use, the pump head can also swivel back out of the way. This head can be configured on either side of the front panel, allowing for convenient positioning on either the left or right wall of the cabinet. In addition, the manual lift operation can accommodate vessels as large as a 2 liter roller bottle.

## Peripette™ Consumable

Peristaltic pump-based systems with flexible tubing have been broadly used in many industries for metering out large volumes of fluid. However, their applications in cell culture have been limited by the complications of maintaining and changing the sterile tubing. These concerns are eliminated by the Peripette, a sterile and easy-to-change consumable. Manufactured from USP Class VI components, the Peripette is gamma-radiation sterilized, non-pyrogenic and completely compatible with standard cell culture reagents and methods, including the transfer of suspended cells.

Ordering information		
Product	Quantity	Cat. No.
Pipeline™ Left-Side Configuration	1 each	4354
Peripette™ Shelf Pack	50	4351
Peripette™ Case	600	4352

## Technical Specifications

DISPENSE SPEEDS	4, 6, and 8 mL/s approximate	
PROGRAMMABLE VOLUMES	0.1 mL – 99.9 mL	<b>0.1 mL increments</b>
	100 ml – 9999 mL	<b>1.0 mL increments</b>
MEMORY	8 user presets	
REMOTE CONTROL	RS-232 Serial Foot Switch	
PRECISION	1 mL Dispense	<b>CV &lt; 2.5%*</b>
	≥2 mL Dispense	<b>CV &lt; 2.0%*</b>
ACCURACY-TYPICAL	± (2.0% + 0.12 mL)	
SOURCE BOTTLE	Max. base diameter	<b>4.75" (120mm)</b>
	Max. height	<b>10.5" (266mm)</b>
POWER	90 x 250 VAC 50-60 Hz 1.5 Amp @ 120V max	
FUSES	5 x 20 mm 1.6 Amp 250V time-delay	
DIMENSIONS (H x W x D)	14" x 11.25" x 8.5" (355 x 285 x 216 mm)	
WEIGHT	24.5 lbs (9.15 Kg)	

\*Each Peripette should be used to dispense no more than 6 liters of fluid.

Learn more at  
[essenbioscience.com/pipeline](https://essenbioscience.com/pipeline)

Or contact us at  
[sales@essenbio.com](mailto:sales@essenbio.com)

Essen BioScience, 300 West Morgan Road, Ann Arbor, Michigan, USA 48108

© 2016 Essen BioScience. All rights reserved. All trademarks are the property of Essen BioScience unless otherwise specified.