

Essen BioScience Pipeline™ and Peripette™ Specifications

Dispense Speeds	4, 6 and 8 mL/s approx.
Programmable Volumes	
0.1 mL – 99.9 mL	0.1 mL increments
100 mL – 9999 mL	1.0 mL increments
Memory	8 user presets
Remote Control	RS-232 Serial Foot Switch
Precision	
1 mL Dispense	CV < 2.5% *
> 2 mL Dispense	CV < 2% *
Accuracy - Typical	± (2.0% + 0.12 mL)
Source Bottle	
Maximum Base Diameter	4.75" (120mm)
Maximum Height	10.5" (266mm)
Power	90-250 VAC 50-60 Hz 1.5 Amp at 120V max
Dimensions (H x W x D)	14" x 11.25" x 7" (355mm x 285mm x 178mm)
Weight	24.5 lbs. (9.15 Kg)

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

Ordering Information

Pipeline™	
Left-Side Configuration	Catalog # 4354
Peripette™	
Shelf pack (qty 50)	Catalog # 4351
Case (qty 600)	Catalog # 4352