

**Medfuture Variable Volume Single Channel
Pipette *1-5 ml***

Catalog number: L55010

Unit size: Each

Applications

This single-channel pipette is easy to use. The ergonomic design of low plunger pressure avoids repetitive strain injuries, light weight, easy to control with full hand, reducing hand fatigue. It is calibrated for precision and accuracy. Easily rotate the piston button to select dispense volume.