

**Medfuture Variable Volume Single Channel  
Pipette \*50-200 ul\***Catalog number: L55007  
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.