SARTORIUS

Tacta® Mechanical Pipettes

Superior Ergonomics

- Easy to clean, with only three parts to disassemble
- Optiload feature in both single and multichannel models for easy and light tip loading
- Optilock's dual-function volume locking prevents accidental volume changes
- Easy-to-read, four-digit volume display helps to set exact volumes
- Color-coding of volumes to ease the selection of corresponding pipette tips
- Safe-Cone Filters for models above 10 μL with convenient built-in filter ejection mechanism

Simplifying Progress

