

## Proline<sup>®</sup> Plus Mechanical Pipettes

### Dependable Durability

The Proline<sup>®</sup> Plus mechanical pipette family is designed to offer comfort and quality for your everyday manual pipetting. It combines durable construction with ease and lightness of use, and is therefore the perfect choice for liquid-handling professionals and students alike. In addition, it has the widest pipette range, including fixed volume pipettes for when volumes need to be ready-set to avoid errors.

#### Features and Benefits

- Low pipetting forces that prevent WRULD and improve results in long pipetting series
- Ergonomic finger support minimizes the grip force needed to hold the pipette
- Optiload mechanism in multi-channel models for easy and light tip loading with perfect tip sealing
- Wide variety of adjustable single and multi-channel models as well as fixed single-channel models
- Volume range from 3  $\mu$ L (5  $\mu$ L for fixed) to 10 mL
- Easy volume setting with click stop mechanism
- Big volume display
- Color-coding of different volumes to ease the selection of matching pipette tips
- Safe-Cone Filters available for models > 10  $\mu$ L
- Fully autoclavable without disassembly
- Simple to clean and maintain with only three parts to disassemble



Discover more about Proline<sup>®</sup> Plus  
Mechanical Pipettes



Please scan!