S/RICS

Picus® and Picus® NxT

The Most Sophisticated and Ergonomic Electronic Pipettes Ever!

- Ergonomic Design for Reduced Risk of RSI (Repititive Strain Injury)
- Innovative Technology for Reliable Results
- Intuitive User-Interface for Ease of Use
- Comprehensive Range of Pipetting Modes for Speed of Use
- Microwell Plate Tracker for Improved Efficiency and Reliability

Simplifying Progress

