

Tacta® & Picus® Pipettes

Reliable and Ergonomic Pipetting

Sartorius simplifies pipetting and offers repeatable results time after time. The pipette and tips are designed to fit together and let you produce repeatable results at the level you need them to be.

Features

- Intuitive and simple user experience
- Low pipetting and tip ejection forces that reduce the risk of repetitive strain injury (RSI)
- Uniquely lightweight and compact
- Advanced safety features: Optilock, Safe-cone filters, password protection, calibration reminders

Benefits

- Achieve reliable results
- Be kind to your hand, by choosing Tacta® and Picus® pipettes
- Speed up your work with electronic Picus® & Picus® Nxt



Discover more about
Tacta® and Picus® Pipettes



Please scan!