

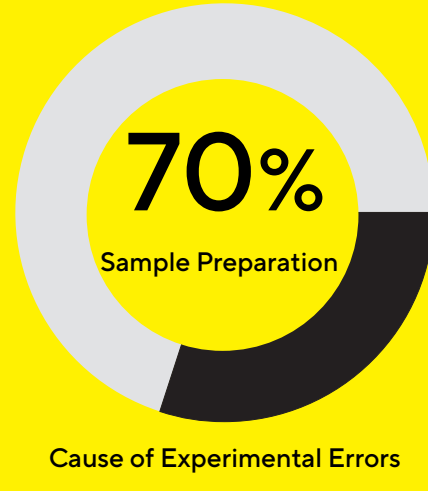
It's Simple: Good Sample Prep Delivers Good Analytical Data



Why Sample Prep Matters

Sample preparation accounts for **up to 70%** of overall errors.

Without high-quality sample preparation workflows, even the most sophisticated instruments can't produce reliable data.



Sensitive Techniques Need Quality Sample Prep

In chromatography, mass spectrometry, spectroscopy, and elemental analysis, even **small errors in prep** can distort the results, impacting reporting and release decisions.

What Is at Stake:

- Calibration curves
- Quantitative accuracy
- Detection limits
- Elemental ratios
- Peak area and retention time

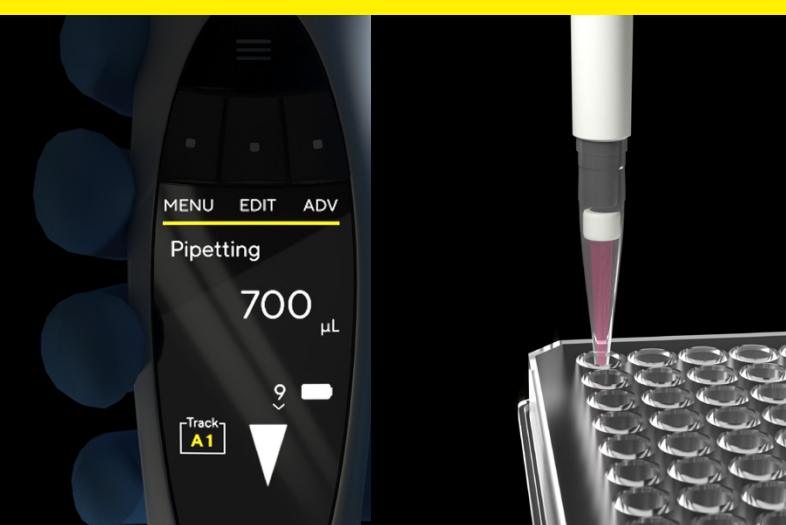


Reduce Risk, Boost Confidence

Manual Pipetting

Tacta® Pipettes and Optifit Tips

Ergonomically designed for precise pipetting



Sample Dilution

Cubis® II Balances and Picus® 2 Pipettes

Equipped with workflow apps to ensure complete traceability

Sample Filtration

Minisart® Syringe Filters

Reliable removal of particulates without analyte loss



Sample Purification

Sartobind® Chromatography Units

More efficient purification of antibodies, viral vectors, nucleic acids, and more



Sample Concentration

Vivaspin® Centrifugal Filters

Faster sample concentration with high target recoveries and easier retentate collection



Pipette Calibration

SpeedCal Mobile

ISO-compliant pipette calibration for multichannel units



A clean prep means I spend less time troubleshooting and more time generating results I trust!

Sartorius tools are built to simplify, standardize, and safeguard your workflow.

- ✓ Reliable
- ✓ Reproducible
- ✓ Regulatory-Ready

