

Preparation of Bioassays and Serial Dilutions in Microplates



How Can QC Labs Achieve Consistent Bioassay Results

Bioassay techniques are essential for quantifying the effects of substances on biological systems, widely used in biotherapeutics, diagnostics, and environmental sciences. Key methods include ELISA for protein and antibody detection and flow cytometry for cell health analysis.



Pipettes and Tips



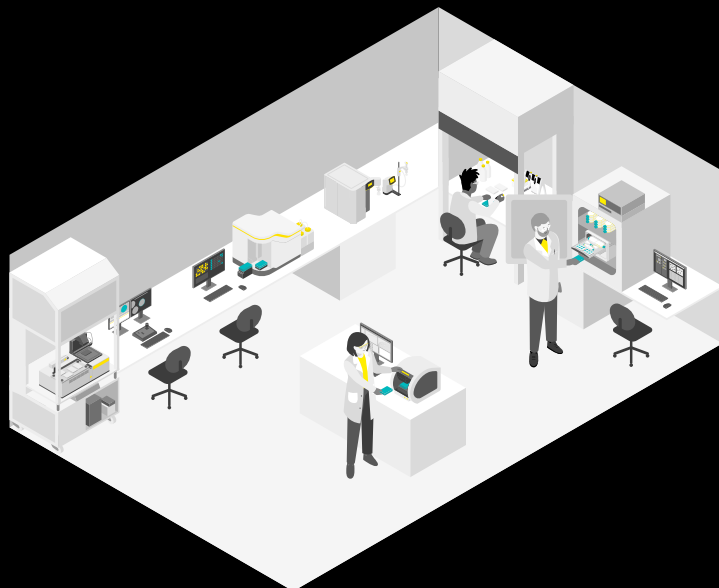
Lab Water Systems



Filtration Devices

Key Needs for Robust Data

- Accurate dispensing across microplate wells
- Contamination-free samples
- Comfortable and easy-to-use lab tools



Picus® 2 Electronic Pipette — Adapt it to Your Needs

Reproducibility

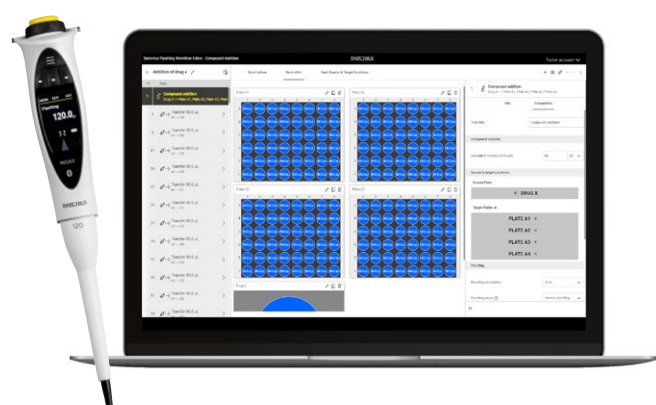
Use predefined, step-by-step protocols to minimize user variance.

No Cross-Contamination

Combine with Safetyspace® Filter tips for contamination-free operation.

Expansive Workflow Library

Access a large selection of protocols with your existing pipette.



Ensure Accuracy with Reliable Pipette Calibration Solutions

Discover More
[ISO 8655 | Sartorius](#)



Minisart® Syringe Filters — Remove Debris and Aggregates



High Recovery

Minimize protein loss with Minisart® NML Cellulose Acetate (CA).

Easier Filtration

Filter samples using less force, thanks to a filtration area of 6.2 cm².

Reduced Waste

Choose bulk packs to reduce waste, with secondary packaging made of cardboard.

Ensure Compliance

Meet ASTM, NCCLS, ISO, and USP standards.

Avoid Contaminants

Keep endotoxins and inorganic salts below detection thresholds.

Fewer Cleaning Cycles

Optimized performance with iJust software: less energy, less feed water

Arium® Pro UF — Ultrapure Water Systems for Cell Cultivation



To learn more, talk to our expert team today



Sample Preparation for QC | Sartorius