

Customized Picus® Bundled with a Molecular Diagnostic System

The Company has been at the forefront of the industry for decades, significantly contributing to public health improvements and fostering a healthier society through its diagnostic solutions. These solutions are crucial for healthcare professionals and industry leaders, enabling them to make decisions that positively impact patient outcomes and consumer safety. Key technologies: immunoassays, molecular biology and microbiology.

Testing Challenge

- Supply preanalytical tools for preparing samples in a Diagnostic system for pathogen detection
- Reduce user-based variance and support on reproducible results, over multiple pipetting steps
- Interference-free filter tips to ensure optimal testing performance
- Global service capability that could be accessed by all end user

The Solution

- Electronic pipette based on the highly versatile Picus® unit with customized protocols including mixing, multi-dispensing and electronic tip ejection
- Optimally fitting DNase-, RNase-, human DNA-free Safetyspace® filter tips
- Global service for maintenance and calibration

The Value

- Sartorius' team of experts allows the Company to focus on its core business, which is the delivery of specific, sensitive and reliable diagnostics to patients
- Steady supply of instruments and consumables secured by regular follow-ups and updated forecasts
- Continued maintenance, calibration and repair of the pipettes



For more information,
visit www.sartorius.com

SARTORIUS