



THE INGENIX SUITE

Bolster data integrity, lab efficiency, and compliance with data and fleet management software

Tracking tools and equipment can consume substantial time and resources in the lab. Regulatory compliance and collaboratively used equipment add an additional level of complexity. Many labs face the problem of mounting paperwork and data management to ensure they meet data integrity, and compliance requirements necessary for continued operations. In addition, when equipment is shared between multiple laboratory groups, centralized reports are needed.

Together, this places considerable work on laboratory staff. Ingenix Suite by Sartorius is designed to address this problem by providing a central, digital database for storage of compliance and routine calibration data and measurement readouts for easy access and organization. The Ingenix Suite Fleet Manager is a platform now enhanced by the newly introduced Ingenix Suite Advanced Pipette Calibration Module.

UTILIZING THE ADVANCED PIPETTE CALIBRATION MODULE

The Advanced Pipette Calibration Module, which is ISO 8655:2022 compliant, offers improved digital calibration management for greater efficiency in the lab. The module can be used with pipettes from any other brand in addition to Sartorius. The Advanced Pipette Calibration Module can manage your pipette calibration data automatically, organizing your readouts and reports, keeping the lab compliant with current regulations, and scheduling calibration reminders ahead of time, preventing a lapse in calibration assessments. Ingenix Suite can also be used in conjunction with Laboratory Information Management System (LIMS) or as a standalone platform.

KEEPING PIPETTE WORK COMPLIANT WITH INGENIX SUITE

Two modules complete the Ingenix Suite for pipettes. The Pipette Check Module enables a convenient means to check the performance of your pipettes during the daily routine, before sensitive applications. The GxP Compliance Module can be used to track overall system- and instrument-level events around all weighing measurements and balance or pipette calibrations. It includes the full audit trail of all audit relevant events and provides the ability to review and approve digital reports, according to the 21 CFR part 11 and EMA Annex 11 needs

WHY CHOOSE INGENIX SUITE

Using the Ingenix Suite modules together offers maximum benefits. Streamlined digital management expedites categorization and facilitation of calibration and compliance documents and identifies equipment problems. Ingenix Software also makes inter-lab sharing of equipment and data easier by creating one localized file repository for documents associated with a piece of equipment. Increase the efficiency, performance, and reliability of your data by utilizing Ingenix Suite to track calibration and compliance documents for pipettes and balances.

SARTORIUS

To learn more, visit:
[www.sartorius.com/en/products/
lab-data-fleet-management-software](http://www.sartorius.com/en/products/lab-data-fleet-management-software)