

# Ingenix® Suite

Elevate Accuracy, Efficiency,  
and Compliance  
With Various Equipment  
Brands in Your Lab



## Highlights

- Open On-Premises Software Solution with Modular design giving maximum flexibility to purchase only what is needed
- Optimizes lab weighing and pipette calibration workflows
- Pipette Calibration Process meets the requirements of ISO 8655:2022 and supports ISO 17025
- Supports compliance with 21 CFR Part 11 and EMA Annex 11
- IQOQ Documentation, Installation and Qualification Services available

## Customizable Modularity for Maximum Flexibility

- Balance Device Manager Module: Manages digital data of Cubis® II MCA lab balances .
- External Balance Calibration Module: Digitally manages Cubis® II MCA lab balance calibrations using external weight sets.
- Pipette Check Module: Quick check of all pipettes for accuracy and precision.
- Advanced Pipette Calibration Module: Simplifies pipette calibration process per latest ISO standards.
- GxP Compliance Module: Compliance with 21 CFR Part 11 and EMA Annex 11 requirements



Learn More:




| Feature                                               | Description                                                                                                                                                                                                                                        |
|-------------------------------------------------------|----------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Centralized Device and Device Management</b>       | Simplifies access to all your equipment and data in one place and through the automated data transfer between device and computer it eliminates manual data entry and reduces error rates.                                                         |
| <b>Automated Job Manager</b>                          | Schedule regular reoccurred Jobs. With specific Sartorius Balances it is possible to send Jobs automatically to the assigned balances.                                                                                                             |
| <b>Compliance Support</b>                             | Helps ensure adherence to 21 CFR Pt. 11 and EMA Annex 11 standards.                                                                                                                                                                                |
| <b>Full Audit Trail</b>                               | The GxP Compliance Module ensures a detailed audit trail for individual or all devices. Advanced filtering options allow for the efficient identification of relevant events based on user, timestamp, completed actions, and additional criteria. |
| <b>Standalone Software and ELNs/ LIMSs Connection</b> | Can be used as stand-alone software but also enables connection to existing laboratory systems.                                                                                                                                                    |
| <b>Brand Flexibility</b>                              | Compatible with all pipette brands and models, whether electronic or mechanical, single or multi-channel, for seamless management of all pipettes across different labs.                                                                           |
| <b>E-Mail Notification</b>                            | Notifications directly via email whenever a maintenance or calibration job is scheduled, streamlining the process, and ensuring timely upkeep.                                                                                                     |
| <b>Role Management</b>                                | Enables the customization of user roles with specific access levels e.g. system configuration, access to groups, and data sets, to manage different responsibilities within the lab.                                                               |
| <b>User Management</b>                                | Supports LDAP or local user management exclusively, depending on your IT infrastructure.                                                                                                                                                           |
| <b>Customizable Report Template</b>                   | By selecting specific sections from report templates, the final report can be refined and streamlined to ensure it meets specific needs and requirements. Upload multiple logos and add standard statements.                                       |
| <b>Flexible File Export Formats</b>                   | File export/saving in pdf, xlsx and xml to enable archiving and ELN/LIMS integration.                                                                                                                                                              |
| <b>Online and Offline Activation</b>                  | Enables license activation online or offline through MySartorius.com. For detailed description look at the manual page 26.                                                                                                                         |
| <b>Scalability</b>                                    | Provides the capability to connect an unlimited number of pipettes and balances.                                                                                                                                                                   |
| <b>Create and send diagnostic files</b>               | Create and send diagnostic files to Sartorius for efficient service process.                                                                                                                                                                       |

## Germany

Sartorius Lab Instruments GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen  
Phone +49 551 308 0

## USA

Sartorius Corporation  
3874 Research Park Dr.  
Ann Arbor, MI 48108  
Phone +1 734 769 1600

 For further information, visit  
[www.sartorius.com](http://www.sartorius.com)