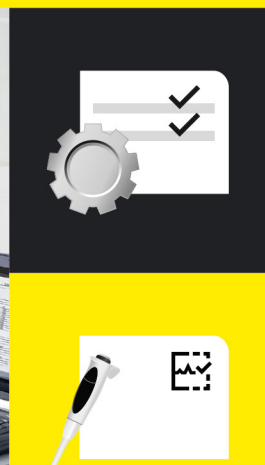
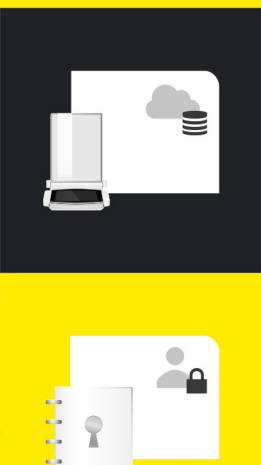


Ingenix Suite - Advanced Pipette Calibration Module

Pipette calibration software for an easy and efficient up-to-date process management

Highlights

- ISO 8655:2022 and CFR 21 Pt. 11 compliant
- Secure SQL database
- Large and growing pipette database
- Languages: English, French, German, Italian, Mandarin



Product Information

The Advanced Pipette Calibration Module is an integral part of the Ingenix Suite software family when it comes to simplifying workflow for your pipette calibration.

The convenient interface offers a guided workflow for your pipette calibration tasks. It allows easy scheduling of calibration intervals and management of the respective pipette fleet.

As every other module, it can be installed and run as a stand-alone software solution for your pipette calibration tasks, or in combination within the Ingenix Suite family of modules.

It is completely compliant with ISO8655:2022, ISO17025 and 21 CFR Pt. 11. Network capabilities with automatic transfer to a SQL database provide central data management and the highest level of security standards.

Technical Specifications

Advanced Pipette Calibration Module

License Period	1 or 5 years
Balance Connections	1 (included) – additional balance connections optional
Supported Sartorius Balances	Speedcal Mobile, Cubis® MPS, Cubis® I, Cubis® II
ISO 8655 Compliance	Yes
21 CFR Pt. 11 Compliance	Yes
SQL Data Base with MS SQL Database	Yes
Languages	English, French, German, Italian and Mandarin
User Management	Yes (LDAP and/or standalone)
Role Management	Yes
GxP Compliance (Audit trail)	Yes
e-Signatures	Yes
Comprehensive Pipette Database	Yes
Data Import of Pipette Database(s)	Yes (CSV import)
Import of Environmental Data	Yes, supported for Sartorius climate modules.
Balances from Multiple Vendors	Supported
Number of Connected Balances	Unlimited
Users per License	Unlimited
Pipette Tip Management	Yes
Report Management	Yes (variable filter functions e.g. date, pipette etc.)
Minimum System Requirements	Laptop with following specifications required: <ul style="list-style-type: none">▪ Processor: 1 gigahertz (GHz) or faster processor or SoC▪ RAM: 2 GB for 64-bit▪ Hard disk space: 12 GB▪ Graphics card: DirectX 9 or later▪ Display resolution: 1024 x 768▪ RS232 ports: For serial connection (optional)

Ordering Information

Description	Order No.
Ingenix Suite Advanced Pipette Calibration	
1 year license	DSX1PC01LWGADC01
5 year license	DSX5PC01LWGADC01
Ingenix 1 Balance licence (optional)	
1 year license	DSX1PC--LWGSBL--
5 year license	DSX5PC--LWGSBL--
Optional Add-ons	
Ingenix Suite Pipette Advanced Calibration Validation Document	DSX-PC01LWGSVM03

How can I get a demo?



Option A: Fill out a request form on the Ingenix Suite Website



Option B: Email us at DigitalSolutionsLPS@Sartorius.com

How can I request a quote?



Option A: Visit our Sartorius eShop for Ingenix Suite



Option B: Contact your local Sales Representative

Discover More



Lab Balance Calibration Solutions



Pipetting and Pipette Tips




Ingenix Suite

Germany

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

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com