

Ingenix[®] Suite

Advanced Pipette Calibration Module



Product Information

The Advanced Pipette Calibration Module optimizes calibration for all pipettes, including third-party brands. It centralizes critical information such as pipette ID, status, calibration history, and type, simplifying oversight and maintenance for lab managers. Pipettes can be set up using Ingenix's built-in SQL Lite database or a SQL database setup on a server. The module ensures accurate data recording and management, allowing data review, approval, or invalidation with the integrated GxP Compliance Module.

Automated workflows and scheduled tasks maintain regular calibration intervals, reducing manual intervention and errors. It supports detailed, customizable calibration reports for internal reviews or regulatory audits, ensuring compliance with industry standards. The module supports unlimited number of pipettes and balances, allowing scalable and cost-effective management as your lab grows.



Learn More:




Device/Instrument Compatibility	Balances	<ul style="list-style-type: none">SpeedCal MobileCubis® MPSCubis® I	<ul style="list-style-type: none">Cubis® IIQuintix®Secura®	<ul style="list-style-type: none">MC210SCPA225DME235PSelected models of 3rd party balances
	Pipettes	<ul style="list-style-type: none">Picus®Picus® 2	<ul style="list-style-type: none">Tacta®Mline	<ul style="list-style-type: none">Proline®Proline® PlusAll third-party Pipettes
Features		<ul style="list-style-type: none">Pipetting process according to the requirement of ISO 8655:2022 Part 6, Part7/ 7aExtended Measurement Uncertainty Calculation in accordance with ISO 8655:2022 and ISO/TR 20461:2023 with the option to include lab and operator effectAbility to upload Calibration and Measurement Capability (CMC) TablePipette Calibration scheduler & trackerCustomizable Reports By selecting specific sections from Report TemplatesUser (LDAP or Local) and Role Management with specific access levelsOnline or Offline License Activation Process through MySartorius.comFile Export/Saving Formats (PDF, XML) to enable archiving and ELN/LIMS integration		
System Requirements		<ul style="list-style-type: none">Processor: 2 GHz or faster, 2-coreMemory: 4 GB RAMHard disk: 4 GB of available disk spaceDisplay: 1280 x 768 screen resolution (32-bit requires hardware acceleration for 4K and higher)Graphics: Graphics hardware acceleration requires DirectX 9 or later, with WDDM 2.0 or higher for Windows 10/128 MB graphics memoryOperating system: Windows OS: Windows 10 / 11.NET version: .NET6 and higherServer: MSSQL server 2019 or later		
Product (Item No.)		Mandatory Modules: <ul style="list-style-type: none">Ingenix® Suite Advanced Pipette Calibration Module incl. 1 Balance License 1 Year License (DSX5PC01LWGADC01) or 5 Year License (DSX5PC01LWGADC01) Optional extras: <ul style="list-style-type: none">Ingenix® Suite Balance License for Single Device 1 Year License (DSX1PC-LWGSBL-) or 5 Year License (DSX5PC-LWGSBL-)Ingenix® Suite Software IQ Documents (DSX-PC01LWGSVM01)Ingenix® Suite Advanced Pipette Calibration Module OQ Document (DSX-PC01LWGSVM03)Ingenix® Suite Installation, Qualification & Training Service(DSX-PC01—GXP01)		

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