

Success Story

Digital Calibration Certificates (DCCs) Save Time and Improve Your Process

How Webers transfers calibration data from Sartorius DCCs easily and efficiently into its own calibration software

Digital calibration certificates increase process reliability and generate real added value by reducing manual steps and avoiding media discontinuities. Webers shows how this can be achieved using digital calibration certificates (DCCs) for balances.

The Application: Pipette Calibration Service

Webers offers calibration services for pipettes for its customers. They are accredited for this service by the DAkkS in accordance with DIN EN ISO/IEC 17025. Webers uses in-house software that supports both test equipment management and the entire pipette calibration process. Calibration data from laboratory balances used for gravimetric pipette calibration are imported into the software.

The Challenge: Entering laboratory balance calibration data

Whenever a laboratory balance is recalibrated, the relevant data from the calibration certificate must be entered into Webers' internal software. Since this data is critical to quality, it requires particular care. When the data is entered manually, a dual-control (four-eye) principle is essential to identify and prevent transcription errors. As this is a recurring task, it involves a significant amount of manual effort, which Webers sought to reduce without compromising data quality.

The Solution: Using DCC to bridge documents and systems

Sartorius is the first provider to offer calibration certificates for balances as digital calibration certificates (DCCs) in a harmonized, well-documented format. DCCs contain all calibration data in a machine-readable XML file. The key practical advantage is that the DCC is embedded directly in the PDF. As a result, the familiar, human-readable original document is preserved while machine-readable data is simultaneously available for further processing.

Key Benefits at a Glance



Reduced manual data entry: Values are no longer typed in but transferred.



Faster process: Data is immediately available for internal steps and verification.



Greater process reliability: Fewer transmission errors, clear assignment of certificate and device.



Practical hybrid: PDF remains readable and verifiable for users, while the DCC makes the calibration certificate machine-readable and thus automatable.

Here's how it works today: One click instead of columns of numbers

In the solution developed by Webers, the digital calibration certificate is imported directly into internal software. The workflow has been deliberately kept simple, i.e. the user uploads the PDF calibration certificate into the application, and the software automatically extracts the embedded DCC as an XML file and immediately displays the relevant information. A brief plausibility check by the employee remains in place as a safety measure while the time consuming manual data entry is eliminated.

What has improved: More speed, less work

Automated transfer reduces manual routines, shortens throughput times and relieves the workload on specialists. At the same time, the risk of classic transfer errors is reduced, and the documentation remains audit-proof because the original PDF is still available.

Outlook: Developing together

DCC formats continue to evolve, guided by feedback from the application. Webers has already demonstrated how the benefits can be realized today. Sartorius will continue along this path by developing real-world formats that promote interoperability and sharing success stories that help other users get started more easily.

"The DCC is designed to ensure that calibration data flows seamlessly into customer applications, eliminating media discontinuities and unnecessary manual effort."

Julian Haller, Team Lead Applied & Legal Metrology, Sartorius



Ready to simplify your calibration data transfer?

Find out more on our website:

www.sartorius.com/en/services/instrument-service/lab-balances-weighing-instrument-services/digital-calibration-certificate

Germany

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

USA

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



For further information, visit

www.sartorius.com