

Calibration Certificate with ISO | IEC 17025 Accreditation

Electronical Weighing Instrument



Benefits

- Reassurance during quality audits
- Calibrations according calibration guideline EURAMET cg-18
- Traceability certificate for your measurement results
- Confidence in the accuracy and quality of your measurement results
- International acceptance and recognition

Product Information

The demands on the accuracy and reliability of laboratory data are constantly increasing. Laboratory scales must be calibrated regularly in order to achieve precise results and meet regulatory requirements. Our Sartorius Service has the right calibration for your needs. Calibration certificates with ISO | IEC 17025 accreditation are internationally recognized and provide the necessary security for your testing materials monitoring. The calibration certificates ensure traceability to national standards and thus confidence in the accuracy of the measurement results. The uncertainty of measurement is also displayed in detail.

Specification

- Accreditation of the calibration laboratory by an independent accreditation body (e.g. DAkkS, UKAS, ANAB)
- Calibration in accordance with the Calibration Guideline EURAMET cg-18
- The repeatability, eccentric loading and error of indication is measured
- Specially adapted measurement of multi range and multi interval balances
- Calculation and indication of the measurement uncertainty
- Indication of the uncertainty of measurement in use
- Easy-to-read display of all measured values in the calibration
- Affixation of a calibration mark
- Guaranteed traceability certificate to the national standard, as required by many quality management standards (e.g. ISO 9001)
- Wide international recognition and acceptance of the calibration certificate

Customer Prerequisites

- If a certificate of conformity is required, the customer tolerances must be defined in advance
- Device is at the maintenance site and has been sufficiently acclimatized
- The device is freely accessible

Optional Services

- As found (initial) calibration to document the balance data before the maintenance
- Multiple possibilities to customise the measurements of repeatability and the test - "error of indication"
- Electronic signature for added security for electronic storage of calibration certificate
- Conformity certificate with confirmation that customer-specific process tolerances and process requirements have been fulfilled
- Minimum sample weight certificate, showing which minimum sample quantities can be measured for different process accuracies and safety factors
- USP <41> certificate
- USP <1251> test report
- Balance test report

You are interested?


Get in touch with us right away. We will be happy to provide you with a non-binding service offer. You can also find further information on our website: www.sartorius.com/en/services

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
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 contacts, visit
www.sartorius.com