

Calibration Certificate with ISO | IEC 17025 Accreditation

Weights and Weight Sets



Benefits

- Calibration of OIML class E2 weights up to 50 kg and F2 weights up to 100 kg
- Proof of traceability for your test weights
- Confidence in the accuracy and quality of your measurement results
- International acceptance and recognition
- Calibration of test weights from all manufacturers
- Short calibration times

Product Information

When testing your balances you must be able to rely on your test weights. Accurately calibrated test weights are the basis of accurate weighing results. Test weights change over time e.g. due to wear and tear or due to becoming dirty. Test weights must therefore 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

- Calibration in our ISO | IEC 17025 accredited calibration laboratory
- Calibration for weights of the OIML class E2 up to 50 kg and F2 up to 100 kg
- Certification of OIML conformity (compliance with the specified accuracy class)
- Calibration certificate stating the conventional mass and | or the mass with deviation from the nominal value and the measurement uncertainty
- Affixation of a calibration mark on the weights case
- Guaranteed traceability certificate to the national standard
- Wide international recognition and acceptance of the calibration certificate
- Easy-to-read display of all measured values in the calibration

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