

## Accredited Calibration For the Highest Quality Results



turning science **into solutions**

## Accurate and Reliable Results

The requirements on the accuracy and reliability of your laboratory data are steadily increasing. To ensure precise, dependable and traceable measurement results of the highest quality, your measuring and test equipment needs to be calibrated at regular intervals. We will be more than happy to perform this service for you.

## Your Equipment Is in the Best Hands

Our Sartorius Service organization is accredited for compliance with EN ISO 17025 in many countries and is authorized to issue calibration certificates for balances, weights and pipettes. As a result, we ensure the 100% traceability of your measuring equipment to national standards and confidence in the quality of your measurements and test results.

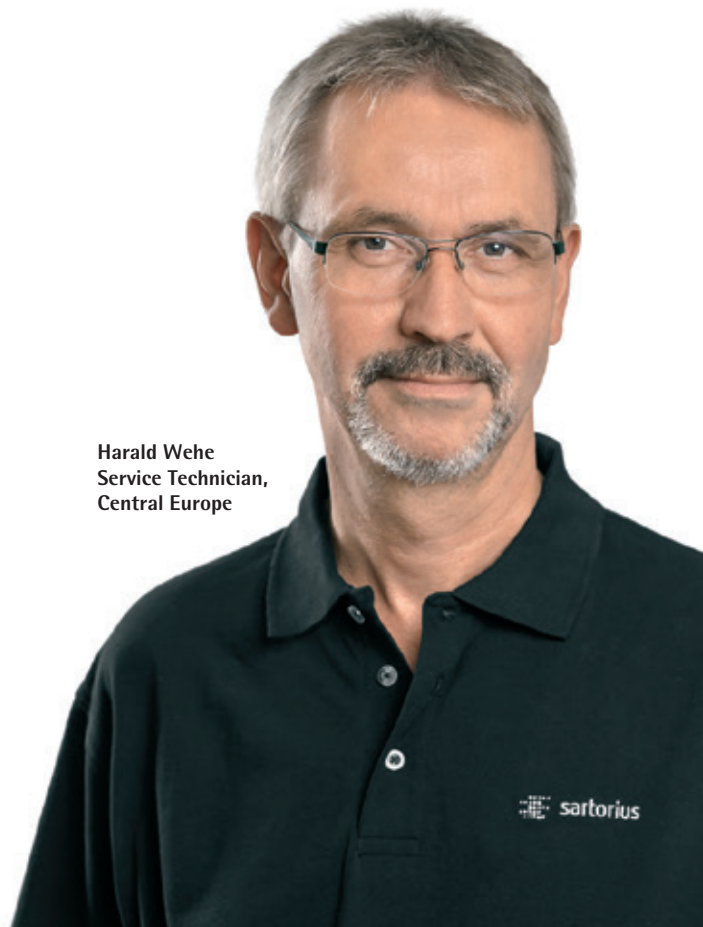
Are you aware that every year we calibrate over 300,000 pipettes and balances worldwide and offer an entire range of services all from a single source? On request, we will also take care of your instruments of different manufacturers' brands.



“ Always be perfectly prepared for the next quality audit.

Our accredited calibration service will eliminate the surprise factor, saving you lots of trouble down the road due to non-conforming equipment.

Harald Wehe  
Service Technician,  
Central Europe



## Accurate and Reliable Results

The requirements on the accuracy and reliability of your laboratory data are steadily increasing. To ensure precise, dependable and traceable measurement results of the highest quality, your measuring and test equipment needs to be calibrated at regular intervals. We will be more than happy to perform this service for you.

## Your Equipment Is in the Best Hands

Our Sartorius Service organization is accredited for compliance with EN ISO 17025 in many countries and is authorized to issue calibration certificates for balances, weights and pipettes. As a result, we ensure the 100% traceability of your measuring equipment to national standards and confidence in the quality of your measurements and test results.

Are you aware that every year we calibrate over 300,000 pipettes and balances worldwide and offer an entire range of services all from a single source? On request, we will also take care of your instruments of different manufacturers' brands.



## Calibration of Laboratory Balances

### Your Advantages:

- Worldwide recognition of your calibration certificates by the International Laboratory Accreditation Cooperation (ILAC)
- Guaranteed traceability to national standards
- Clear presentation of all measured results and of the calculated measurement uncertainty on the calibration certificate
- On-site calibration by specially trained service technicians who use highly accurate calibration weights

## Calibration of Pipettes

### Your Advantages:

- Reliability and compliance with EN ISO 8655
- Guaranteed traceability to national standards
- Precise results: Trust in the accuracy and the best possible performance of your pipette
- The choice is all yours: have us calibrate your pipette either at our laboratory or on-site at your facility
- Added services available on request: Cleaning and maintenance of pipettes

## Calibration of Test Weights

### Your Advantages:

- Calibration of weights of OIML R 111 classes E1, E2, F1 and F2 up to 50 kg
- Guaranteed traceability to national standards
- Inspection and cleaning of your test weights



# Contact

Don't forget to have your balances, pipettes and test weights calibrated at regular intervals.

Just complete the form provided at

[www.sartorius.com/service](http://www.sartorius.com/service)

and we'll get in touch with you right away.

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289  
[www.sartorius.com](http://www.sartorius.com)

Specifications subject to change without notice.  
Copyright Sartorius Lab Instruments GmbH & Co. KG.  
Printed in the EU on paper bleached without chlorine.  
Publication No.: S--1548-e170501  
Order No.: 85037-559-09  
Ver. 05 | 2017