

Balance Test Report

Testing against
Manufacturer, Customer or
OIML R76 Tolerances



Benefits

- Possibility to test against different tolerance settings
- Proof of whether the balance is working within the tolerance settings
- Multiple possibilities to adapt the test to customer needs
- Easy-to-read certificates

Product Information

Test equipment must work within the specified tolerance ranges to achieve successful quality assurance. For this purpose, equipment must be tested at suitable time intervals. Since there are no general normative specifications for non-automatic balances that specify a tolerance, for example depending on the resolution of the device, the manufacturer's tolerances or the user's tolerance specifications are needed when evaluating balances. With a balance test report, balances can be tested against different tolerance requirements. The measurement results are documented in an easily readable certificate.

Specification

- Calibration against manufacturer, customer or OIML R76 tolerance settings
- Examination of key performance parameters of the balance
- Depending of the test report the repeatability, eccentric loading, adjustment accuracy, accuracy and linearity is assessed
- When testing against manufacturer tolerances environmental conditions at the installation place are taken into account
- When testing against customer tolerances the possibility is given to test against warning and action | intervention limits
- The possibility is given also to test balances without type approval against tolerances based on the OIML R76
- Clear assessment of whether the balance is working inside or outside the tolerance limits
- Easy-to-read display of all results in the report

Customer Prerequisites

- When creating a balance test report against customer tolerances the tolerances must be specified in advance
- Device is at the maintenance site and has been sufficiently acclimatized
- The device is freely accessible

Optional Services

- Calibration certificate with ISO | IEC 17025 accreditation
- USP <41> certificate
- USP <1251> test report
- Minimum weight certificate according Eur. Pharmacopoeia (2.1.7.)

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