

Certificate of Calibration

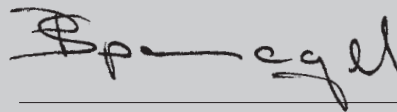
arium® TOC Instrument

This product has passed Sartorius' in-house tests, thus meets Sartorius' stringent quality control standards and the technical specifications in the documentation at the time of shipment from the warehouse.

The measurement standards used are calibrated by comparison to measurement against national standards as defined by the National Institute of Standards and Technology within the limitations of the Institute's Calibration Service, or by accepted values of natural physical constants, or by consensus standards or by other internal measurements of Sartorius.

This product is developed, produced and distributed according to a Quality Management System that is certified for compliance with DIN | ISO 9001.

Please note we recommend a regular calibration verification once a year.



Calibration date

Patricia Spannagel
Quality Manager

Manufactured by
Sartorius Weighing Technology GmbH
Weender Landstrasse 94-108
37075 Goettingen, Germany
www.sartorius.com

