



CERTIFICATE



This is to certify that



Sartorius Lab Instruments GmbH & Co. KG

Otto Brenner Straße 20
37079 Göttingen
Germany

has implemented and maintains a
Quality and Environmental Management System.

Scope:

Development, production, sales and service of mechatronic products and services for weighing technology, lab water systems and process automation, service of pipettes, calibration laboratories

Through an audit, documented in a report, it was verified that the management system fulfills the requirements of the following standards:

ISO 9001 : 2015

ISO 14001 : 2015

Certificate registration no.	000440 QM15 UM15
Valid from	2018-08-15
Valid until	2021-08-14
Date of certification	2018-08-15



DQS GmbH

Stefan Heinloth
Managing Director