

Certificate of Conformity

Compliance Certificate to DIN ISO 3696, Grade 1 Water

Parameter	Requested Specification	arium® pro ¹ & comfort ²
pH-value at 25 °C	not applicable	not applicable
Electrical conductivity, in µS/cm at 25 °C, max.	0.1	0.055
Oxidizable matter, Oxygen content in mg/l, max	not applicable	not applicable
Absorbance at 254 nm and 1 cm optical path length, absorbance units, max.	0.001	< 0.001
Residue after evaporation on heating at 110°C in mg/kg, max.	not applicable	not applicable
Silica (SiO ₂) content in mg/l, max.	0.01	< 0.01

Execution and Analysis Procedure


The water analysis was executed by RH GmbH, an independent company for analytical analysis, based on the DIN ISO 3696/DIN 38404-C3 and DIN ISO 3696. The samples were taken directly from the water dispense outlet.

¹ Determined with DI water (Goettingen)

² Determined with municipal water (Goettingen)

19.05.2015

Date



Dr. Susan Gerlachhausen
Director of Quality

Sartorius Lab Instruments GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen, Germany

www.sartorius.com