

Certificate of Conformity

Compliance Certificate to DIN ISO 3696, Grade 1, System Series: Arium® Pro

Parameter	Requested Specification	Arium® Pro
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. The tests were performed with the Arium® Pro VF (H2Opro-VF-T), without final filter, fed with DI water (Goettingen).



Mansoor Mobarak
Plant Manager Quality

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0
www.sartorius.com

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG.
Status: 02 | 2022