

Certificate of Conformity

Compliance Certificate to DIN ISO 3696, Grade 3 Water

Parameter	Requested Specification	arium® advance EDI ¹
pH-value at 25 °C	5.0 - 7.5	7.5
Electrical conductivity, in µS/cm at 25 °C, max.	5.0	< 0.2
Oxidizable matter, Oxygen content in mg/l, max	0.4	< 0.4
Absorbance at 254 nm and 1 cm optical path length, absorbance units, max.	not applicable	not applicable
Residue after evaporation on heating at 110°C in mg/kg, max.	2	< 2
Silica (SiO ₂) content in mg/l, max.	not applicable	not applicable

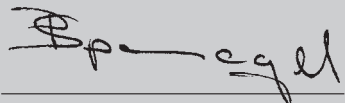
Execution and Analysis Procedure

The water analysis was executed by RH GmbH, an independent company for analytical analysis, based on the DIN ISO 3696. The samples were taken directly from the water dispense outlet and stored in a 1.5 liter glass bottle.

¹ Supplied with municipal water (Göttingen)

06.04.2016

Date



Patricia Spannagel
Plant Manager Quality

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

www.sartorius.com