

Certificate of Conformity

Compliance Certificate to DIN ISO 3696, Grade 2, System Series: Arium® Advance EDI

Parameter	Requested Specification	Arium® Advance EDI
pH-value at 25 °C	not applicable	not applicable
Electrical conductivity, in $\mu\text{S}/\text{cm}$ at 25 °C, max.	1.0	< 1.0
Oxidizable matter, Oxygen content in mg/l, max	0.08	< 0.08
Absorbance at 254 nm and 1 cm optical path length, absorbance units, max.	0.01	< 0.01
Residue after evaporation on heating at 110 °C in mg/kg, max.	1	< 1
Silica (SiO_2) content in mg/l, max.	0.02	< 0.02

Execution and Analysis Procedure

The water analysis was executed by SYNLAB Analytics & Services Germany GmbH, an independent company for analytical analysis, based on the DIN ISO 3696. The samples were taken directly from the product water outlet. The tests were performed with the Arium® Advance EDI (H2O-EDI-1-T), without final filter, fed with municipal water (04416 Markkleeberg).



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