

Certificate of Conformity

Compliance Certificate to to ASTM D1193-06, Type II, System Series:
System Series: Arium® Comfort II, Type II Water

Parameter	Requested Specification	Arium® Comfort II
Production Process	Distillation or alternatively Ion exchange, reverse osmosis and organic absorption	Ion exchange, reverse osmosis and organic absorption
Electrical conductivity, in $\mu\text{S}/\text{cm}$ at 25 °C, max.	1.0	1.0
pH	not applicable	not applicable
TOC $\mu\text{g}/\text{L}$ (max.)	50	< 50
Sodium $\mu\text{g}/\text{L}$ (max.)	5	< 5
Chloride $\mu\text{g}/\text{L}$ (max.)	5	< 5
HBC cfu/mL (max.)	not applicable	not applicable
Endotoxin, EU/mL (max.)	not applicable	not applicable
Total Silica $\mu\text{g}/\text{L}$ (max.)	3	< 3

Execution and Analysis Procedure

The analysis were performed according to the ASTM D1193-06. The conductivity and TOC tests were executed by Sartotius quality lab with calibrated equipment. Sodium, chloride and silica testing was performed by SYNLAB Analytics & Services Germany GmbH, an independent company for analytical analysis. The samples were taken directly from the product water outlet. The tests were performed with the Arium® Comfort II (H2O-II-1-TOC-T), without final filter, fed with tap water Goettingen, TOC 1,000 ppb and conductivity 280 $\mu\text{S}/\text{cm}$.



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