

## Certificate of Conformity

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

Parameter	Requested Specification	Arium® Comfort I
Production Process	Purify to 20 µS/cm by dist. or equiv., followed by mixed bed DI, 0.2 µm filtration.	Purify to 20 µS/cm by dist. or equiv., followed by mixed bed DI, 0.2 µm filtration.
Electrical conductivity, in µS/cm at 25 °C, max.	0.0555	0.0555
pH	not applicable	not applicable
TOC µg/L (max.)	50	< 50
Sodium µg/L (max.)	1	< 1
Chloride µg/L (max.)	1	< 1
HBC cfu/mL (max.)	not applicable	not applicable
Endotoxin, EU/mL (max.)	not applicable	not applicable
Total Silica µg/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 I (H2O-I-2-TOC-T), without final filter, fed with tap water Goettingen, TOC 1,000 ppb and conductivity 280 µS/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