

Certificate of Conformity

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

Parameter	Requested Specification	Arium® Pro
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 Sartorius 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® Pro VF (H2OPRO-VF-T-TOC), without final filter, fed with DI water.



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: 08 | 2022