

Certificate of Conformity

arium® advance, arium® comfort Water Contacting Components

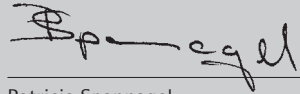
We hereby confirm, that the materials listed below, used in the product water cycle of arium® advance and in the pre-treatment cycle of arium® comfort, are internally tested and released for their compatibility in the production of pure water.

- PP
- Natural PSU (arium® advance EDI & comfort II)
- POM EPDM
- Stainless steel 316
- Stainless steel 1.4401

arium® pro, arium® comfort Water Contacting Components

We hereby confirm, that the materials listed below, used in the ultrapure water cycle of arium® pro and arium® comfort systems, are internally tested and released for their compatibility in the production of ultrapure water.

- PVDF
- EPDM
- POM
- Stainless steel 1.4404
- PEEK



Patricia Spannagel
Quality Management



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

www.sartorius.com