

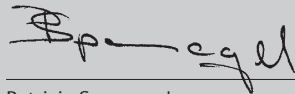
# Certificate of Conformity

## ASTM Compliance Certificate to ASTM D1193-06 Type 1 Water

### arium<sup>®</sup> Laboratory Water Systems Models: pro DI, pro UV, pro UF, pro VF, pro, comfort I, comfort II, mini & mini plus

We hereby certify that the above mentioned products are capable of producing Ultrapure Water that meets or exceeds the current published ASTM D1193-06 Standard Specification for Reagent Water Type 1.

Sartorius further certifies that the above mentioned products are in compliance with ASTM D1193-06 regarding the mixed bed ion exchange resin and the 0.2 µm membrane filter as installed.



Patricia Spannagel  
Quality Management



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

[www.sartorius.com](http://www.sartorius.com)