



arium® Cartridge Sets

for ultrapure water systems



Specifications

Materials of Construction

Housing:	Virgin polypropylene
Connecting screws:	Stainless steel
Purification media:	Spherical catalytical, activated carbons. Semiconductor-grade ultrapure mixed bed ion exchange resin.

Description

The new arium cartridge sets, for arium ultrapure water purification systems, are designed to produce the highest purity water and exceed all specifications of the ASTM Type 1 standard.

Each cartridge set, consisting of one pretreatment and one polishing cartridge, is specifically formulated for each particular arium system. The exceptional quality of the purification media ensures the best product water specs in the industry.

The cartridges are optimized to remove organic and inorganic contaminants. Highly porous activated carbon combined with efficient ion exchange resins enables long cartridge life with unparalleled performance, producing consistent, reproducible results.

In addition, Sartorius Stedim Biotech purification media has been specifically developed to produce water with the lowest total organic carbon (TOC) content for arium systems utilizing ultraviolet (UV) radiation technology.

All cartridges are engineered so that water flows from top-to-bottom through the ion exchange medium, which prevents separation of the mixed bed resin and allows full use of each cartridge, with no wasted media.

Order Number	Product Name	Description	Capacity to 18.2 MΩ × cm ultrapure water as CaCO ₃ in Grain	Capacity to 18.2 MΩ × cm ultrapure water in eq
H2O-A-PACK	Analytical Kit	Cartridge Set for arium® pro/611 VF or UV ultrapure water systems	965	1.25
H2O-B-PACK	Biological Kit	Cartridge Set for arium® pro/611 UF ultrapure water systems	1141	1.48
H2O-E-PACK	Elemental Kit	Cartridge Set for arium® pro/basic/611 DI ultrapure water systems	1268	1.64
H2O-U-PACK	Universal Kit*	Cartridge Set for arium® ultrapure water systems with tap water feed	965	1.25

* Tap water cartridges should be used only after the feed water specifications have been checked and following consultation with the responsible Sartorius application specialist.

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Japan +81.3.3740.5407

Specifications subject to change
without notice. Printed and copyrighted
by Sartorius Stedim Biotech GmbH. | W
Publication No.: SLG2034-e11084
Order No.: 85030-518-89
Ver. 08 | 2011