



arium® 611 Cartridge Kits

Disposables for Ultrapure Water



Specifications

Materials of Construction

Housing:	Virgin unfilled polypropylene
Connecting screws:	Stainless steel
Purification media:	20 × 50 mesh granular activated and catalytic activated carbons. Semiconductor-grade ultrapure mixed bed ion exchange resin.

Description

The Sartorius arium® 611 cartridge kits are designed to be installed in the arium® 611 laboratory water purification systems. Each cartridge kit utilizes Sartorius patented technologies to provide ultrapure water that exceeds the ASTM Type 1 Reagent Grade Water Standards. Kits are designed to produce 18.2 MΩ × cm ultrapure water with a low TOC content.

All cartridge kits contain 1 pretreatment pack, 1 polishing pack and 2 final filters (see additional data sheet on the Sartopore 2 150 final filters).

Features and Benefits

The cartridge packs are designed for down-flow operation to promote efficient purification kinetics and to prevent media separation. Inner tube diameters and volumes are engineered to fluid cross-sectional velocity and media bed contact standards.

The packs contain the highest quality granular and catalytic activated carbons. Only new semiconductor (electronics) grade mixed bed ion exchange resin is utilized in the cartridge kits. These kits have some of the highest ion exchange capacity of the industry, resulting in low operating costs.

Application Specific Media Design

Each unique cartridge kit is designed for a specific feed water source and arium® model to provide the highest product water quality available. Special kits are engineered for the different and variable feed water characteristics expected from deionized, reverse osmosis or tap (mains) feed water sources. Purification media recipes are also specifically designed for arium® systems utilizing ultraviolet (UV) radiation technology for the lowest total organic carbon (TOC) content. Ultraviolet systems can produce peroxides and ozone that can damage ion exchange resins. A special catalytic material is used to remove peroxide and ozone downstream of the UV chamber. Kits for low TOC without the UV chamber and high capacity ion exchange are also available.

Laboratory Water Systems

Cartridge Kit for 611DI and 611UF Systems

Order No.	Feed Water Source	Grain Capacity to 1.0 MΩ × cm as CaCO ₃	Grain Capacity to 10 MΩ × cm as CaCO ₃	eq (val)
611CKDI	Deionized EDI	1470	1323	67.4
611CKDO	Deionized EDI	1260	1134	57.8
611CKRI	RO and distilled	1470	1323	67.4
611CKRO	RO and distilled	1260	1134	57.8
611CKHI	RO, distilled, deionized EDI	1630	1467	74.8
611CKTI	Tap mains	1470	1323	67.4
611CKTO	Tap mains	1260	1134	57.8

Cartridge Kit for 611UV and 611VF Systems

Order No.	Feed Water Source	Grain Capacity to 1.0 MΩ × cm as CaCO ₃	Grain Capacity to 10 MΩ × cm as CaCO ₃	eq (val)
611CKDU	Deionized EDI	1140	1026	52.3
611CKRU	RO and distilled	1140	1026	52.3
611CKTU	Tap mains	1140	1026	52.3

Cartridge Kit Order Number Description

6 1 1 C K R U

arium®
product line

Consum-
ables kit

Polishing pack (Pack 2)

I = Low inorganic application

O = Low organic application for 611DI & 611UF systems

U = Low organic application for 611UV & 611VF systems

Pretreatment pack (Pack 1)

R = RO and distilled feed water

D = Deionized | EDI feed water

H = RO, distilled, deionized | EDI feed water,
high capacity ion exchange

T = Tap feed water

Sartorius AG
Weender Landstrasse 94–108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

www.sartorius.com/arium

USA +1.631.2544249
UK +44.1372.737100
France +33.1.69192100
Italy +39.055.634041
Spain +34.91.3586100
Japan +81.3.33295533

Specifications subject to change without notice.
Printed in Germany on paper that has been
bleached without any use of chlorine.
W/sart-105 · G
Publication No.: SLG2034-e03081
Order No.: 85030-518-89