



arium® 615DI Deionization Cartridges

Water Pre-treatment Systems



Features

- Low water consumption
- Compact design, easy to install
- Reliable procedure
- Conductivity typically < 0.1 $\mu\text{S}/\text{cm}$
- High efficiency

Description

The arium® 615DI deionization cartridges are designed to produce highly deionized water.

These cartridges contain highly effective ion exchange resins to remove cations and anions from water by exchanging H^+ and OH^- . Their uniform and compact mixed bed ensures reliable desalination and optimal utilization of the resin material, reducing conductivity by up to < 0.1 $\mu\text{S}/\text{cm}$.

The arium® line also includes conductivity meters for checking the quality of product water. These convenient options are available in a choice of LED controlled display or digitally controlled display.

The LED controlled conductivity meter enables the user to monitor the conductivity of the product water with a red (> 20 $\mu\text{S}/\text{cm}$) or green (< 20 $\mu\text{S}/\text{cm}$) light. If the water does not meet the user's quality requirements, the LED will light up in red to alert the user.

The digitally controlled conductivity meter lets the user monitor the conductivity by individually defined setpoint limits (0.1–200 $\mu\text{S}/\text{cm}$). The conductivity measurement range is between 0.056 and 200 $\mu\text{S}/\text{cm}$.

Applications

Pre-treatment for:

- Ultrapure water systems
- Autoclaves
- Glassware machinery, etc.

Post-treatment for:

- RO permeate

Specifications

Description of the DI cartridge		615DI004	615DI007	615DI010	615DI014
Material of construction:		Plastic (white)	Plastic (white)	Stainless steel	Stainless steel
Capacity:	at 10° dH	840 l	1,300 l	2,100 l	2,800 l
	at 20° dH	420 l	650 l	1,050 l	1,400 l
Flow:		40 l/h	300 l/h	300 l/h	600 l/h
Dimensions:	Height	40 cm	140 cm	44.5 cm	60 cm
	Diameter	23 cm	23 cm	23.7 cm	23.7 cm
Connectors:	Inlet (male)	3/4"	3/4"	3/4"	3/4"
	Outlet (male)	3/4"	3/4"	3/4"	3/4"
Operating pressure:	Max.	6 bar	6 bar	6 bar	6 bar
	Min.	0.5 bar	0.5 bar	0.5 bar	0.5 bar
Operating temperature: Max.		35°C	35°C	35°C	35°C
Volume of resin:		5 l	12 l	15 l	20 l

Order No.	Description
615DI004	arium® deionization cartridge including set of tube connectors, capacity of up to 420 l at 20°dH
615DI007	arium® deionization cartridge including set of tube connectors, capacity of up to 650 l at 20°dH
615DI010	arium® deionization cartridge including set of tube connectors, capacity of up to 1,050 l at 20°dH
615DI014	arium® deionization cartridge including set of tube connectors, capacity of up to 1,400 l at 20°dH
615ADC1	arium® conductivity meter for deionization cartridges, with digital conductivity display; conductivity measurement range 0.056–200 µS/cm
615ALC1	arium® conductivity meter for deionization cartridges, with LED display (red >20 µS/cm & green <20 µS/cm)
615DIC1	arium® replacement package for deionization cartridges, ion exchange resin mixture in 25 l PE bag

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.: SLG2046-e11022
Order No.: 85032-538-40
Ver. 02 | 2011