

Consumables

Arium® Pro Cartridge Sets

Pretreatment and post-treatment cartridge using top-down technology

- High performance capacity thanks to efficient ion exchange resins
- Effective adsorption of impurities through high-grade activated carbon
- Optimized flow pattern, prevents separation of the mixed resin bed
- Patented connection process simplifies the replacement of consumables



Description

The cartridge sets are optimized for the removal of both organic and inorganic constituents. Every set has been designed specifically to match the unit and delivers ultrapure water that even exceeds the ASTM type 1 quality standard. This consistent level of high-quality water ensures optimal reproducibility of your results.

Optimized purification materials such as highly effective activated carbon coupled with a efficient ion-exchange resins

deliver long-lasting performance and thereby ensure long maintenance intervals.

The top-down flow technology produces ideal purification kinetics and prevents any mixing of cleaning media. The cartridge was designed with the applicable standards for flow rate in the cross section and contact time with the medium in mind.

Technical Specifications | Ordering Information

Materials	
Housing	High-purity polypropylene
Fixing screws	Stainless steel
Cleaning media	Spherical, catalytic activated carbons Ultrapure mixed bed ion exchange resin
Feed water requirements	see "Technical Specifications" page 2

Exchange capacity for 18.2 MΩ × cm ultrapure water relative to CaCO₃

	[Grain]	[Equivalent]
Analytical Kit	965	1.25
Biological Kit	1,14	1.48
Elemental Kit	1,268	1.64
Universal Kit	965	1.25

* With the Universal Kit, Arium® Pro can be directly fed with untreated drinking water to produce ultrapure water. The appropriate Sartorius application specialists should be consulted to check the feed water specifications.

Order number	Description
H2O-A-PACK	Analytical Kit, Arium® Pro cartridge set for biological, chemical-analytical and standard ultrapure water applications, 1 pc
H2O-B-PACK	Biological Kit, Arium® Pro cartridge set for biological ultrapure water applications, 1 pc
H2O-E-PACK	Elemental Kit, Arium® Pro cartridge set for standard ultrapure water applications, 1 pc
H2O-U-PACK*	Universal Kit, Arium® Pro cartridge set for untreated feedwater*, 1 pc

Intended Use

Device type:

- H2O-A-PACK - Arium® Pro VF and Pro UV
- H2O-B-PACK - Arium® Pro UF
- H2O-E-PACK - Arium® Pro and Pro DI
- H2O-U-PACK* - Arium® Pro, Pro DI, Pro UF, Pro UV and Pro VF