

Arium® Ultrafilter

Ultrapure water, free from endotoxins, DNases and RNases

- High flow rates
- Integrity tested
- Long service lives
- Certified quality

Description

The hollow-fiber ultrafilter utilizes crossflow technology to reliably remove bacterial endotoxins, micro-organisms and particles, as well as DNases and RNases from the ultrapure water.

The filters are developed and produced in accordance with a DIN EN ISO 9001|DIN EN 46 001 certified quality assurance system which fulfills the prerequisites of the Quality System Regulation (Regulation on Quality Systems of the FDA) 21 CFR Part 820. During the manufacturing process, every unit is integrity tested to meet the highest quality standards and safety regulations.



Technical Specifications | Ordering Information

Materials	
Membrane	Polysulfone
Composite material	Polyurethane (PUR)
Housing, caps	Polycarbonate (PC)
Plugs	Polypropylene (PP)
Efficient membrane surface	2.1 m ²
Max. operating pressure	3 bar at room temperature
Sterilization	200 ppm sodium hypochlorite, 45 min, max. 1 x/week
Tank volume	
Lumen	152 ml
Filtrate side	306 ml

Retention of bacteria and endotoxins	
<i>Brev. diminuta</i>	LRV 7 – 10
<i>E. coli</i> O55:B5 endotoxin	LRV > 3,5
Natural endotoxins	LRV > 3,0
Dimensions of the fibers	
Interior diameter	215 µm
Wall thickness	50 µm
Molecular weight cut-off (MWCO)	5,000 (= 5 kD)
Order number	Description
611CDU5	Arium® Ultrafilter, 1 pc

Intended Use

Device type:

- Arium® Pro VF and Pro UF