



Hollow Fiber Ultrafilter 611CDU5



Specifications

Materials

Membrane	Polysulfone
Potting material	Polyurethane (PUR)
Housing, caps	Polycarbonate (PC)
Plugs	Polypropylene (PP)

Effective membrane area
2.1 m²

Connectors Inlet | Outlet
Acc. to EN 1283

Maximum Transmembrane Pressure (TMP)
0.8 bar at 0°C.
Maximum operating pressure up to 3 bar (45 psi) at room temperature.

Sanitization

200 ppm sodium hypochlorite, 45 min.
Frequency: not more than once a week.

Priming volume

Lumen	152 ml
Filtrate side	306 ml

Fiber dimensions

Inner diameter	215 µm
Wall thickness	50 µm
Molecular Weight Cut off (MWCO)	5.000

Order information

Order code	611CDU5
Weight	0.24 kg
Quantity per package	1 unit

Description

Disposable, self-contained, hollow fiber ultrafilter designed for arium® 611 Laboratory Water Systems. High flow rate, low extractable ultrafilter suitable for laboratory use.

Applications

Removal of bacterial endotoxins, bacteria and particles from ultrapure water with a resistivity of up to 18.2 megohms × cm at 25°C.

Quality

The filters are designed and manufactured in a quality management system in compliance with DIN EN ISO 9001/DIN EN 46 001 and the FDA Quality System Regulation, 21 CFR Part 820.

100% integrity tested prior to release.

Bacterial and Endotoxin Retention

Bacteria:
Brevundimonas diminuta
ATCC 19146 19146 in saline lactose broth.
Cell diameter approx. 0.3 mm 7-10 LRV*

Endotoxin:
E. coli O55:B5 endotoxin (Whittaker, USA) > 3.5 LRV*

Natural endotoxins:
Endotoxins in water from various sources > 3.0 LRV*

* LRV = Logarithmic reduction value;

$$LRV = \log_{10} \frac{\text{Number of organisms in challenge suspension}}{\text{Number of organisms in filtration}}$$

Performance

Filtration flow Q _F (ml/min)	Pressure drop Δp (mmHg)	
	Solution 21°C	Solution 37°C
300	43 – 64	29 – 43
500	71 – 107	48 – 71
700	100 – 150	67 – 100

Determined with 0.9 % sodium chloride solution. Pressure drop Δp is measured between fluid inlet and filtrate outlet.

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/arium

USA +1.800.3687178
UK +44.1372.737100
France +33.1.69192100
Italy +39.055.634041
Spain +34.91.3586102
Japan +81.3.37405407

Specifications subject to change
without notice. Printed and copyrighted
by Sartorius Stedim Biotech GmbH
W/sart-000 · G
Publication No.: SLG2028-e07086
Order No.: 85030-515-94