



arium® CellPlus Ultrafilter

For effective removal of endotoxins
in cell culture applications

Advantages

- Effective removal of RNase | DNase
- Reliable removal of endotoxins
- High flow rate performance
- Certified quality
- Sterile-packaged



Product Description

The arium® CellPlus is a point-of-use ultrafilter for efficient removal of endotoxins, RNase, DNase, microorganisms and particles.

Designed for arium® comfort and arium® mini ultrapure water systems, this sterile-packaged ultrafilter provides the highest safety for your critical cell culture applications. A protective bell supplied with the ultrafilter additionally prevents retrograde contamination.

Moreover, the high-grade material selected for arium® CellPlus enables excellent total throughputs and optimal flow rates.

Technical Specifications

Materials	
Membrane	Polysulfone
Composite material	Polyurethane (PUR)
Housing	Acrylonitrile butadiene styrene (ABS)
Protective bell	Polycarbonate (PC)

Typical Specifications	
Flow rate*	Up to 2.0 l/min
Endotoxins	< 0.001 EU/ml
Bacteria	< 1 cfu/100 ml
RNase concentration	< 1 pg/ml
DNase concentration	< 5 pg/ml

* depends on the inlet pressure and type of system

General Specifications	
Dimensions (height × diameter)	169 × 50 mm
Max. operating pressure	6 bar (87 psi)
Max. inlet temperature	50°C
Effective membrane area	0.5 m ²

Ordering Information

Order Number	Quantity per Package
H2O-CUF	1

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius.com

USA Toll-free +1.800.635.2906
UK +44.1372.737159
France +33.1.70.62.50.00
Italy +39.0362.5557.11
Spain +34.913.586.095
Russian Federation +7.812.327.53.27
Japan +81.3.3740.5408

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG.
Printed in the EU on paper bleached without chlorine.
Publication No.: SLG2060-e170203
Order No.: 85037-556-48
Ver. 02 | 2017