

Ultrafiltration Membranes

Optimize Your Process:
Discs for Sample
Concentration and
Buffer Exchange



Product Information

Our comprehensive range of ultrafiltration membranes offers the flexibility to optimize your concentration and buffer exchange processes for most macromolecules. Available in a variety of diameters, these membranes are ideal for use in stirred cells and custom applications.

Cellulose triacetate (CTA) is a hydrophilic, low-binding membrane that is cast without a substructure that could otherwise trap passing micromolecules. This makes it ideal for analytical applications where the permeate is to be examined, such as sample fractionation, deproteinization, or the removal of other high molecular weight interfering substances.

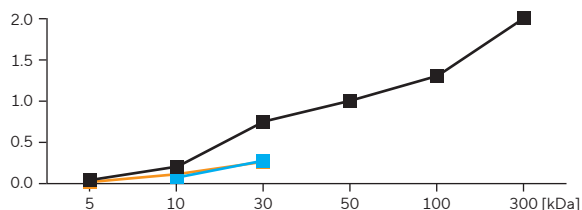
Hydrosart® regenerated cellulose (RC) is a hydrophilic membrane that has been optimized for the biopharmaceutical industry to be non-protein binding and virtually non-fouling, while offering extended chemical resistance. It is often preferred to achieve high recoveries from very dilute or valuable samples, can be sterilized by autoclaving, and is easy to clean.

Polyethersulfone (PES) exhibits no hydrophobic or hydrophilic interactions and delivers excellent performance in research and biotechnology applications. Optimized to provide exceptional flux and high biomolecule recovery, its broad pH and chemical resistance also ensure suitability for a wider range of sample compositions.

Technical Specifications

	CTA	RC	PES
Thickness	120 µm	180 µm	120 µm
Chemical compatibility	pH 4–8	pH 1–13	pH 1–14
Water flux [mL/min/cm ²] ¹	0.11	0.08	0.20
Retention rate ²	90%	99%	95%

Figure 1: Water Flux (mL/min/cm²) for CTA, Orange, RC, Blue, and PES, Black.



Ordering Information

Membrane Material	MWCO	Diameter			
		25 mm	47 mm	63 mm	76 mm
Cellulose triacetate (CTA)	5 kDa	14529--25-----D	14529--47-----D	-	-
	10 kDa	-	14539--47-----D	-	-
	20 kDa	14549--25-----D	14549--47-----D	14549--63-----D	-
Hydrosart® regenerated cellulose (RC)	2 kDa	-	-	14419--63-----D	-
	5 kDa ³	14429--25-----D	14429--47-----D	14429--63-----D	14429--76-----D
	10 kDa	14439--25-----D	14439--47-----D	14439--63-----D	14439--76-----D
	30 kDa	14459--25-----D	14459--47-----D	14459--63-----D	14459--76-----D
	100 kDa	-	14468--47-----D	-	-
Polyethersulfone (PES)	1 kDa	-	14609--47-----D	14609--63-----D	14609--76-----D
	5 kDa	14629--25-----D	14629--47-----D	14629--63-----D	14629--76-----D
	10 kDa ⁴	14639--25-----D	14639--47-----D	14639--63-----D	14639--76-----D
	30 kDa	14659--25-----D	14659--47-----D	14659--63-----D	14659--76-----D
	50 kDa	-	14650--47-----D	-	14650--76-----D
	100 kDa	-	14668--47-----D	14668--63-----D	-
	300 kDa	-	14679--47-----D	-	14679--76-----D

¹ For 10 kDa membranes.

² For cytochrome c (12.4 kDa) on 10 kDa membranes.

³ Also available in 44 mm diameter (order number 14429--44-----D).

⁴ Also available in 43 mm and 150 mm diameters (order numbers 14639--43-----D and 14639-150-----D).

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
sartorius.com