



sartorius

Sartorius Membranes and Papers for Blotting Applications

Blotting Membranes – Features & Benefits

- High membrane surface area for high binding capacity & no sample loss
- Exceptionally low background allowing longer exposure times & better results
- High membrane stability for easy handling



Blotting Papers – Features & Benefits

- Made of high-purity cotton linters for uniform buffer flow and resulting blots
- No additives to avoid any interference during the transfer
- Supplied in sheets, rolls as well as in customized sizes to save time and avoid any waste

Blotting Membranes



Sartorius nitrocellulose membranes are available in two pore sizes, 0.22 μm and 0.45 μm . Both versions combine the advantages of high protein binding capacity with low background and high membrane stability, which ensures easy handling. Due to its large surface area, the 0.22 μm membrane version is recommended for small proteins.

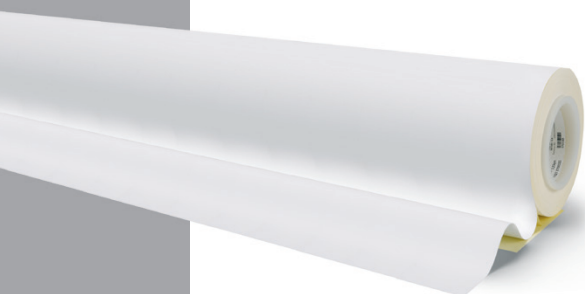
Sartorius blotting membranes are ideal for western blotting, DNA blotting as well as dot or slot blots. They have been optimized for all protein blotting systems, such as electrotransfer, semi-dry or simple capillary blotting.

Specifications Blotting Membranes

Description	11327	11306
Material	Cellulose Nitrate	Cellulose Nitrate
Pore size (μm)	0.22	0.45
Thickness (μm)	130	130
Water flow rate (mL/[min.cm ² bar])	27	70
Bubble point with water (bar)	4.4	2.4
Extractables in water (%)	< 1	< 1
Burst pressure (bar)	0.8	0.2
Binding capacity for IgG ($\mu\text{g}/\text{cm}^2$)	200	200

Ordering Information

Grade	Roll Size	Order Number
11327	30 cm \times 3 m	11327-----41BL
11306	30 cm \times 3 m	11306-----41BL



Blotting Papers

Sartorius blotting papers are made of cotton linters without any additives in order to completely prevent potential interference with transfer processes. They are available in a choice of different weights and thicknesses as well as in different formats to cover the majority of blotting applications. An overview of standard sizes for the most commonly used paper grades is provided in the following. Customized formats and other blotting grades are available on request.



Specifications Blotting Papers

Description	FN 100	BF 3	BF 4
Weight (g/m ²)	195	330	550
Thickness (mm)	0.35	0.76	1.30
Capillary rise (mm/30 min)	115	30	-
Capillary rise (mm/10 min)	70	130	160

Ordering Information

Size (mm)	Qty./Box	Order Number
Grade FN 100		
200 \times 200 mm	100	FT-2-527-200200N
260 \times 410 mm	100	FT-2-527-260410N
460 \times 570 mm	100	FT-2-527-460570N
580 \times 600 mm	50	FT-2-527-580600K
580 \times 680 mm	50	FT-2-527-580680K
Grade BF 3		
135 \times 155 mm	100	FT-2-520-135155N
200 \times 200 mm	100	FT-2-520-200200N
460 \times 570 mm	50	FT-2-520-460570K
580 \times 600 mm	50	FT-2-520-580600K
Grade BF 4		
110 \times 170 mm	25	FT-2-521-110170G
150 \times 150 mm	25	FT-2-521-150150G
580 \times 580 mm	25	FT-2-521-580580G
580 \times 600 mm	25	FT-2-521-580600G

Sartorius Lab Instruments
GmbH & Co. KG
Weender Landstrasse 94-108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

Specifications subject to change without notice.
Printed and copyrighted by
Sartorius Stedim Biotech GmbH. | W
Publication No.: SL-1112-e150702
Order No.: 85037-545-62
Ver. 07 | 2015

www.sartorius.com

