

## 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  $\mu\text{m}$  and 0.45  $\mu\text{m}$ . 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  $\mu\text{m}$  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.

## Typical values

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

## Ordering Information

	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 <sup>2</sup> )	195	330	550
Thickness (mm)	0.35	0.76	1.30
Capillary rise (mm/30 min)	115	-	-
Capillary rise (mm/10 min)	70	130	160

## Ordering Information

Size (mm)	Qty./Box	Order Number
<b>Grade FN 100</b>		
200 × 200 mm	100	FT-2-527-200200N
260 × 410 mm	100	FT-2-527-260410N
460 × 570 mm	100	FT-2-527-460570N
580 × 600 mm	50	FT-2-527-580600K
580 × 680 mm	50	FT-2-527-580680K
<b>Grade BF 3</b>		
135 × 155 mm	100	FT-2-520-135155N
200 × 200 mm	100	FT-2-520-200200N
224 × 224 mm	100	FT-2-520-224224N
460 × 570 mm	50	FT-2-520-460570K
580 × 600 mm	50	FT-2-520-580600K
<b>Grade BF 4</b>		
110 × 170 mm	25	FT-2-521-110170G
150 × 150 mm	25	FT-2-521-150150G
580 × 580 mm	25	FT-2-521-580580G
580 × 600 mm	25	FT-2-521-580600G




**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 contacts, visit  
[www.sartorius.com](http://www.sartorius.com)