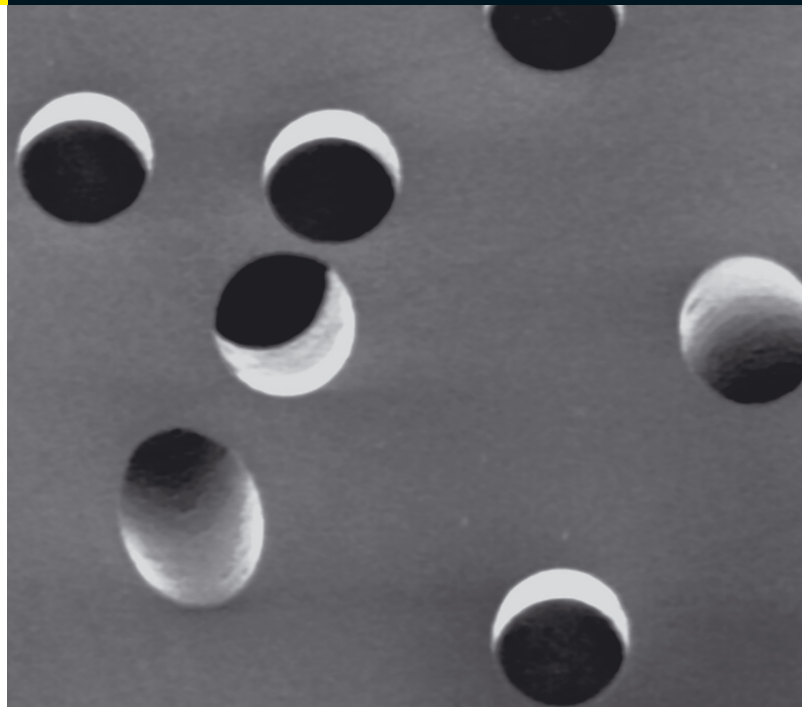


# Polycarbonate Track-Etched Filters



## Benefits

- Sharply defined pore sizes for accurate separation and retention of particles
- Smooth and translucent surface for excellent visibility of the retained particles
- Capture of the particles on the surface of the filter for precise quantification and reading under light microscopes or for SEM

## Product Information

Polycarbonate track-etched filters are an excellent choice when precise surface capture and high sample visibility are required. They are available in a choice of pore sizes and diameters to cover a broad range of applications.

## Description

Polycarbonate track-etched filters are manufactured from high-grade polycarbonate film. This film is irradiated with ions that cross the material by breaking polymer chains. After ion passage, chemical etching dissolves these zones of broken polymer chains to form cylindrical pores of exactly defined sizes and densities.

The results are a well-calibrated pore size, narrow pore size distribution and a uniform structure to retain all particles of interest on the surface of the filter.

## Applications

The major applications for these track-etched filters are particulate analysis, fluid clarification, cytology, cell biology, bioassays, water microbiology and environmental analyses (for example, asbestos).

# Technical Data

## Specifications

Material	Polycarbonate
Color	White, translucent
Structure	Symmetrical
Surface property	Hydrophilic
Binding	Exceptionally low non-specific adsorption
Extractables	Extremely low
Particle release	No
Biologically inert	Yes
Hygroscopic	No
Thickness (µm)	25
Burst strength (bar)*	≥0.7
Autoclavable (30 min at 121°C)	Yes
Temperature resistance (max. continuous operating temp.)	140°C
pH range	1 – 13

\* Pressure required to cause 1 cm<sup>2</sup> of non-supported membrane to rupture

## Typical Results

Specifications   Grade	23058	23007	23006	23004
Pore size (µm)	0.1	0.2	0.4	0.8
Pore density (cm <sup>2</sup> )	6.0E+08	5.0E+08	1.50E+08	4.0E+07
Water flow rate (mL/min/cm <sup>2</sup> at 0.7 bar)	≥0.5	≥10	≥30	≥40
Air flow rate (L/min/cm <sup>2</sup> at 0.7 bar)	≥0.2	≥1	≥3	≥7.5
Water bubble point (bar)	≥7.0	3.5	2.0	0.6

# Ordering Information

Grade	Pore Size ( $\mu\text{m}$ )	Diameter (mm)	Qty./Pkg.	Order Number
23058	0.1	25	100	23058--25-----N
23058	0.1	47	100	23058--47-----N
23007	0.2	25	100	23007--25-----N
23007	0.2	47	100	23007--47-----N
23007	0.2	50	100	23007--50-----N
23006	0.4	25	100	23006--25-----N
23006	0.4	47	100	23006--47-----N
23004	0.8	25	100	23004--25-----N

Other pore sizes and diameters are available on request.

# Sales and Service Contacts

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

## **Germany**

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

## **U.K.**

Sartorius UK Ltd.  
Longmead Business Centre  
Blenheim Road, Epsom  
Surrey KT19 9QQ  
Phone +44 1372 737159