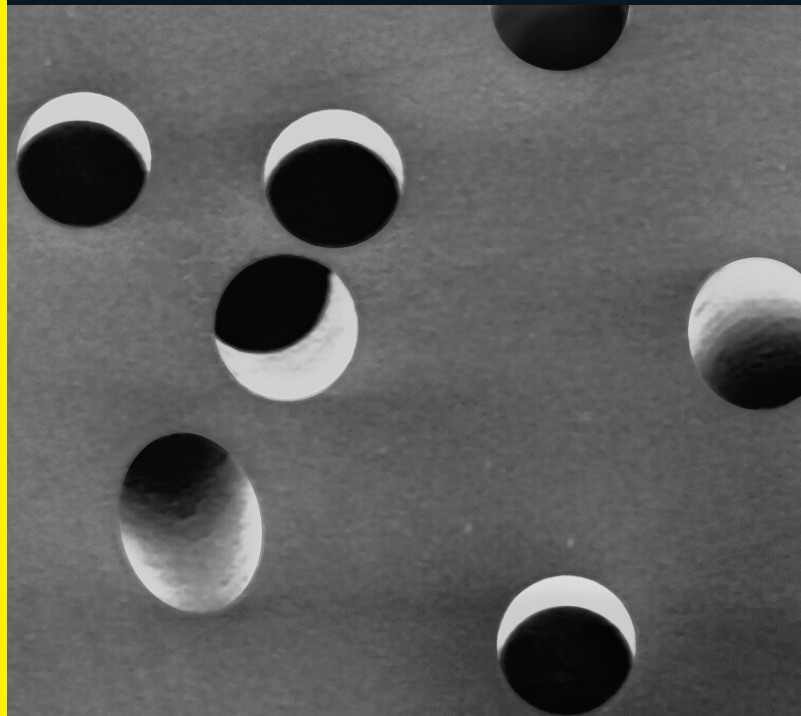


Polycarbonate Track-Etched Filters



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
Autoclavable (30 min at 121 °C)	Yes
Temperature resistance (max. continuous operating temp.)	140 °C
pH range	1 - 13

Typical Results

Specifications Grade	23058	23007	23006	23004	23A42	23015
Pore size (µm)	0.1	0.2	0.4	0.8	5	15
Thickness (µm)	25	25	25	25	11	37
Pore density (cm ⁻²)	6.0E+08	5.0E+08	1.50E+08	4.0E+07	4.0E+05	5.0E+04
Water flow rate (mL/min/cm ² at 0.7 bar)	≥ 0.5	≥ 10	≥ 30	≥ 40	≥ 900	≥ 1200
Air flow rate (L/min/cm ² at 0.7 bar)	≥ 0.2	≥ 1	≥ 3	≥ 7.5	≥ 50	≥ 100
Water bubble point (bar)	≥ 7.0	3.5	2.0	0.6	N/A	N/A
Burst strength (bar)*	≥ 0.7	≥ 0.7	≥ 0.7	≥ 0.7	N/A	N/A

* Pressure required to cause 1 cm² of non-supported membrane to rupture

Ordering Information

Grade	Pore Size (μm)	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
23006	0.4	76	50	23006--76-----K
23004	0.8	25	100	23004--25-----N
23A42	5	47	100	23A42--47-----N
23015	15	47	100	23015--47-----N

Other pore sizes and diameters are available on request.

Germany

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

USA

Sartorius Corporation
3874 Research Park Dr.
Ann Arbor, MI 48108
Phone +1 734 769 1600

 For additional information,
visit www.sartorius.com