

Pre-Heated Quartz Microfiber Filters Grade Q3400



Quartz Microfiber Filters

Properties

- High-purity filters with the lowest trace metal values
- Extreme temperature resistance
- Exceptional chemical resistance
- Excellent weight and dimensional stability
- Biologically inert

Description

The quartz microfiber material of the Sartorius pre-heated filters, grade Q3400, is made of high-purity quartz microfibers without any addition of glass microfibers or binding agents. In addition, the Q3400 filter grade is tempered to remove all chemically combined water and to give the filters excellent weight and dimensional stability. Sartorius filters are especially suitable for emissions monitoring at temperatures of up to 900°C and wherever filters of the highest purity are needed. These quartz microfiber filters are even purer than glass microfiber filters.

Applications

- Emissions monitoring at high temperatures
- Air pollution monitoring (US EPA PM 10*)
- Analysis of hot and acidic gases
- Trace metal element analysis
- Analytical and gravimetric analyses
- Analysis of dust levels according to EN 13284 1: 2001

*US EPA PM 10: United States Environmental Protection Agency – Particulate Matter, Level 10



Specifications

| | |
|---|--|
| Material | 100% high-purity quartz, binder-free, tempered |
| Basis weight | 85 g/m ² |
| Thickness | 0.43 mm |
| Penetration, 0.3 µm 15 cm/s using paraffin oil (EN 143) | <0.002% |
| Efficiency, 0.3 µm 15 cm/s using paraffin oil (EN 143) | 99.998% |
| Temperature resistance | Up to 900°C |

Ordering Information

| Order Number | Ø in mm | Qty. per Box |
|-----------------|---------|--------------|
| Q3400--20-----G | 20 | 25 |
| Q3400--25-----G | 25 | 25 |
| Q3400--30-----G | 30 | 25 |
| Q3400--37-----G | 37 | 25 |
| Q3400--45-----G | 45 | 25 |
| Q3400--47-----G | 47 | 25 |
| Q3400--50-----G | 50 | 25 |
| Q3400--82-----N | 82 | 100 |
| Q3400--90-----N | 90 | 100 |
| Q3400-142-----K | 142 | 50 |
| Q3400-150-----K | 150 | 50 |

Other diameters or formats are available on request.

Germany

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

For further information, visit
www.sartorius.com

USA

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