



sartorius

Filter Papers for the Laboratory and Industry



turning science **into solutions**



Table of Contents

4	Filter papers – An Introduction
5	Product Overview
6	Ash-free Filter Papers for Quantitative and Gravimetric Analyses
8	Wet-strengthened Filter Papers for Qualitative Analyses
10	High-purity Filter Papers for Qualitative Analyses
12	Smooth Filter Papers for Qualitative & Technical Analyses
14	Crêped Filter Papers for Qualitative & Technical Analyses
16	Paper Boards for the Filtration and Absorption of Liquids
18	Seed Testing Papers
20	Filter Papers for the Sugar Industry
21	Diatomaceous Earth Filter Paper
22	Phase Separating Paper
23	Nonwovens
24	Weighing Paper
24	Lens Cleaning Paper
25	Surface Protection Paper
26	Blotting Papers
27	Blotting Membranes
28	Chromatography Papers
30	Glass Microfiber Filters With Binder
32	Glass Microfiber Filters Without Binder
34	Pre-heated Quartz Microfiber Filters
35	Quality Control Test Methods
36	Index of Grades



Filter Papers – An Introduction

High-grade filter papers are indispensable for routine work in laboratory and industrial applications. Sartorius supplies you with a broad range of filter papers for myriad filtration tasks and supports you in solving all your filtration challenges.

With this catalog, we invite you to familiarize yourself with our broad product range. Here, you will find typical examples intended to help you quickly select the filter paper that is right for your application.

Our Product range covers:

- Quantitative, qualitative filter papers
- Technical papers and boards
- Blotting and chromatography papers
- Glass and quartz microfiber filters
- And many other paper grades for special applications

Quality Assurance and Quality Control

Sartorius pays particular attention to continuous in-process quality control; additionally, regular checks and exact analyses of raw material and of each individual finished product assure constant high quality and product uniformity.

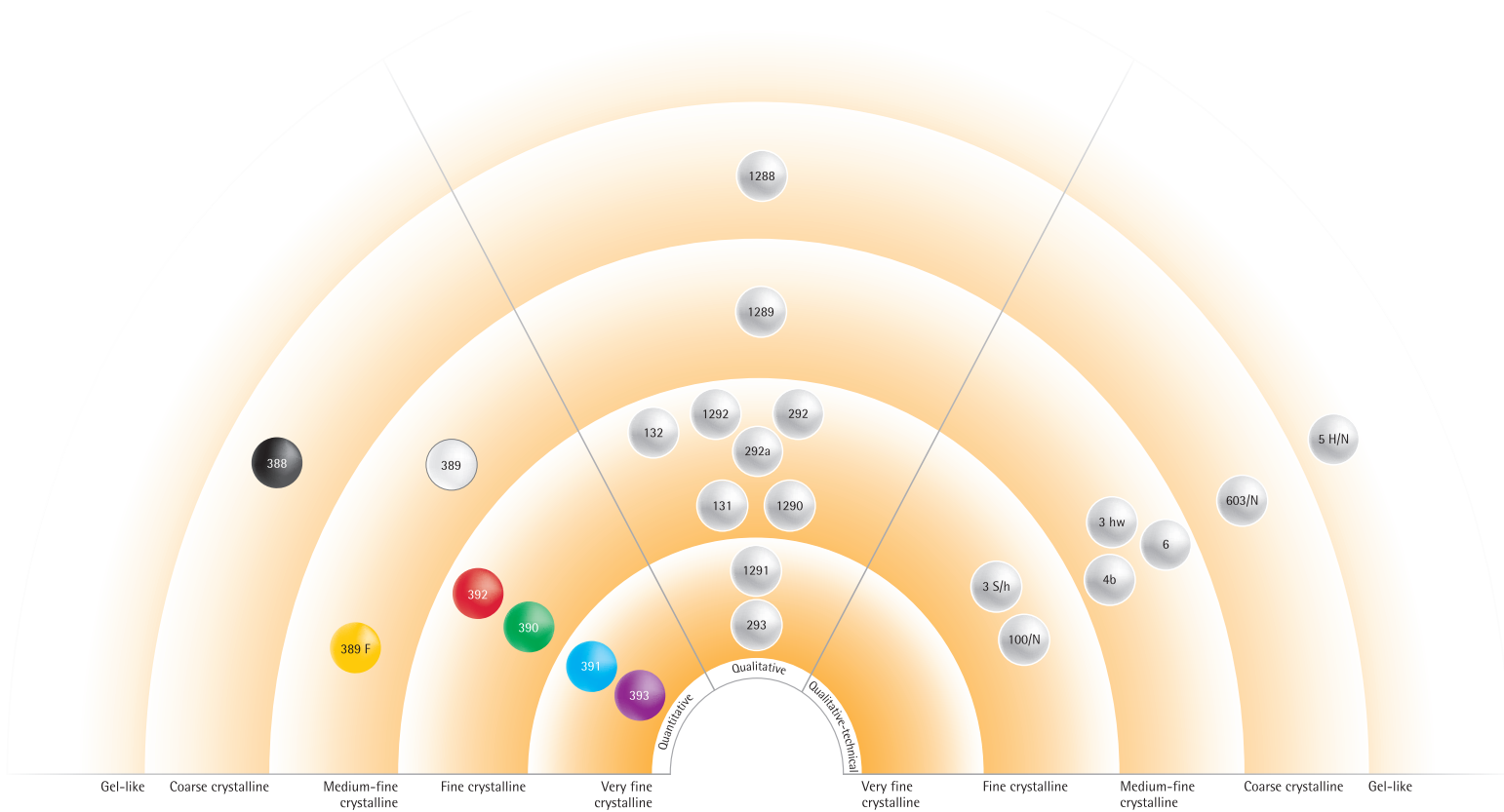
The paper mill meets the requirements set by the ISO 9001 quality management system and the ISO 14001 environmental management system.

How Do Filter Papers Work?

Filter papers are actually depth filters. Various parameters influence their effectiveness: Mechanical particulate retention, absorption, pH, surface properties, thickness and strength of the filter paper as well as the shape, density and quantity of particles to be retained. The precipitates deposited on the filter form a "cake layer", which – depending on its density – increasingly affects the progress of a filtration run and decisively affects the retention capability. For this reason, it is essential to select the right filter paper to ensure effective filtration. This choice also depends on the filtration method to be used, among other factors. In addition, the amount and properties of the medium to be filtered, the size of the particulate solids to be removed and the required degree of clarification are all decisive in making the right choice.



Product Overview



Quantitative Filter Papers



Black dot

Fast filtering, wide-pore, loose structure, ash-free, wet-strengthened



White dot

Medium fast filtering, medium- to wide-pore, ash-free, wet-strengthened



Yellow dot

Medium fast filtering, medium- to wide-pore, low-fat content, ash-free, wet-strengthened



Red dot

Medium fast filtering, medium density ash-free, wet-strengthened



Green dot

Slow filtering, narrow-pore, dense, ash-free, wet-strengthened



Blue dot

Very slow filtering, fine-pore, very dense, ash-free, wet-strengthened



Purple dot

Very slow filtering, very fine-pore, very dense, ash-free, wet-strengthened

Ash-free Filter Papers for Quantitative and Gravimetric Analyses

These filter papers are used for quantitative and gravimetric analyses as well as for pressure or vacuum filtration. They are made out of 100 % cotton linters with an α -cellulose content of > 98 % and are acid-washed to make the papers ashless and achieve high purity.

In gravimetric applications, the cake layer is calcined and the residue quantified. For quantitative analysis of the filtrate, the filter paper must not give off any foreign substances. This guarantees that no test results are falsified. That is why it is important that the filters are ash-free.

For some quantitative analyses, the cake layer has to be mechanically removed from the filter (for example, with a water jet or a spatula). The filter must be wet-strengthened so that it doesn't break when the cake layer is removed.










Application Examples

Application	Grade
Determination of ash content	388
Gravimetric analysis of metals	388
Analysis of alkaline earth carbonates	389
Determining the fate content in natural raw materials	389 F
Gravimetric analyses in power plants	392
Filtration of fine precipitates	390
Filtration of fine-grained precipitates	391, 393
Blaine Test for cement (EN 196-6:2010)	392, 391, 390, 389

- Made of 100 % cotton linters
- Ash-free (Ash content $\leq 0,01$ % according to DIN 54370)
- Wet-strengthened
- Color-coded box for easy selection
- Supplied in rolls, sheets, discs and folded filters

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Particle retention (µm)	Filtration (s)*	Precipitates	Properties
 388	84	0.21	12–15	10	Coarse crystalline	Wide-pore, loose structure, fast filtering
 389	84	0.19	8–12	20	Medium-fine crystalline	Medium- to wide-pore, medium fast filtering
 389 F	84	0.19	8–12	20	Medium-fine crystalline	Medium- to wide-pore, medium fast filtering
 392	84	0.17	5–8	50	Fine crystalline	Medium dense, medium fast filtering
 390	84	0.16	3–5	100	Fine crystalline	Narrow-pore, dense, slow filtering
 391	84	0.15	2–3	180	Very fine crystalline	Fine-pore, dense, very slow filtering
 393	100	0.17	1	300	Very fine crystalline	Very fine-pore, very dense, very slow filtering

* See test methods, page 35

Ordering Information



Filter Discs, 100 pieces

Ø in mm	Grade 388	Grade 389	Grade 389 F	Grade 390	Grade 391	Grade 392	Grade 393
55	FT-3-101-055	FT-3-102-055	FT-3-112-055	FT-3-103-055	FT-3-104-055	FT-3-105-055	FT-3-127-055
70	FT-3-101-070	FT-3-102-070		FT-3-103-070	FT-3-104-070	FT-3-105-070	FT-3-127-070
90	FT-3-101-090	FT-3-102-090	FT-3-112-090	FT-3-103-090	FT-3-104-090	FT-3-105-090	FT-3-127-090
110	FT-3-101-110	FT-3-102-110	FT-3-112-110	FT-3-103-110	FT-3-104-110	FT-3-105-110	FT-3-127-110
125	FT-3-101-125	FT-3-102-125	FT-3-112-125	FT-3-103-125	FT-3-104-125	FT-3-105-125	FT-3-127-125
150	FT-3-101-150	FT-3-102-150	FT-3-112-150	FT-3-103-150	FT-3-104-150	FT-3-105-150	FT-3-127-150
185	FT-3-101-185	FT-3-102-185	FT-3-112-185	FT-3-103-185	FT-3-104-185	FT-3-105-185	FT-3-127-185
240	FT-3-101-240	FT-3-102-240		FT-3-103-240	FT-3-104-240	FT-3-105-240	FT-3-127-240



Folded Filters, 100 pieces

Ø in mm	Grade 388	Grade 389	Grade 389 F	Grade 390	Grade 391	Grade 392
110	FT-4-101-110	FT-4-102-110		FT-4-103-110	FT-4-104-110	FT-4-105-110
125	FT-4-101-125	FT-4-102-125		FT-4-103-125	FT-4-104-125	FT-4-105-125
150	FT-4-101-150	FT-4-102-150		FT-4-103-150	FT-4-104-150	FT-4-105-150
185	FT-4-101-185	FT-4-102-185	FT-4-112-185	FT-4-103-185	FT-4-104-185	FT-4-105-185
240	FT-4-101-240	FT-4-102-240			FT-4-104-240	



Sheets in 580 × 580 mm, 100 pieces

Grade 388	Grade 389	Grade 390	Grade 391	Grade 392	Grade 393
FT-2-101-580580	FT-2-102-580580	FT-2-103-580580	FT-2-104-580580	FT-2-105-580580	FT-2-127-580580

Other dimensions are available on request

Wet-strengthened Filter Papers for Qualitative Analyses

These qualitative filter papers are essentially used for analytical purposes and routine analyses, whenever no gravimetric analyses are required. They are wet-strengthened and can be used for pressure and vacuum filtration. They are made of refined pulp and linters with an > 95% α -cellulose content, are very pure with an ash content $\leq 0.1\%$.

Application Examples

Application	Grade
Must analysis	1288
Routine filtration for malt analysis	1289
Rapid filtration of fine precipitates	1292
Analysis of coffee extracts	1290
Tannin solutions	1291
Clarification of wine	293



- Made of refined pulp and cotton linters with an > 95% α -cellulose content,
- Ash content $\leq 0,1\%$ according to DIN 54370
- Wet-strengthened
- Supplied in rolls, sheets, discs and folded filters

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Particle retention (μ m)	Filtration (s)*	Precipitates	Properties
1288	84	0.21	12 – 15	10	Coarse crystalline	Wide-pore, loose structure, fast filtering
1289	84	0.21	8 – 12	20	Medium-fine crystalline	Medium- to wide-pore, medium fast filtering
1292	84	0.17	5 – 8	50	Fine crystalline	Medium dense, medium fast filtering
1290	84	0.15	3 – 5	100	Fine crystalline	Narrow-pore, dense, slow filtering
1291	84	0.15	2 – 3	180	Very fine crystalline	Fine-pore, dense very slow filtering
293	80	0.15	1 – 2	300	Very fine crystalline	Very fine-pore, very dense very slow filtering

* See test methods, page 35

Ordering Information



Filter Discs, 100 pieces

Ø in mm	Grade 1288	Grade 1289	Grade 1290	Grade 1291	Grade 1292	Grade 293
55	FT-3-206-055	FT-3-207-055	FT-3-208-055	FT-3-209-055	FT-3-210-055	FT-3-211-055
70	FT-3-206-070	FT-3-207-070	FT-3-208-070	FT-3-209-070	FT-3-210-070	FT-3-211-070
90	FT-3-206-090	FT-3-207-090	FT-3-208-090	FT-3-209-090	FT-3-210-090	FT-3-211-090
110	FT-3-206-110	FT-3-207-110	FT-3-208-110	FT-3-209-110	FT-3-210-110	FT-3-211-110
125	FT-3-206-125	FT-3-207-125	FT-3-208-125	FT-3-209-125	FT-3-210-125	FT-3-211-125
150	FT-3-206-150	FT-3-207-150	FT-3-208-150	FT-3-209-150	FT-3-210-150	FT-3-211-150
185	FT-3-206-185	FT-3-207-185	FT-3-208-185	FT-3-209-185	FT-3-210-185	FT-3-211-185
240	FT-3-206-240	FT-3-207-240	FT-3-208-240	FT-3-209-240	FT-3-210-240	



Folded Filters, 100 pieces

Ø in mm	Grade 1288	Grade 1289	Grade 1290	Grade 1291	Grade 1292	Grade 293
110	FT-4-206-110	FT-4-207-110	FT-4-208-110	FT-4-209-110	FT-4-210-110	
125	FT-4-206-125	FT-4-207-125	FT-4-208-125	FT-4-209-125	FT-4-210-125	FT-4-211-125
150	FT-4-206-150	FT-4-207-150	FT-4-208-150	FT-4-209-150	FT-4-210-150	FT-4-211-150
185	FT-4-206-185	FT-4-207-185	FT-4-208-185	FT-4-209-185	FT-4-210-185	FT-4-211-185
240	FT-4-206-240	FT-4-207-240	FT-4-208-240	FT-4-209-240	FT-4-210-240	FT-4-211-240



Sheets in 580 × 580 mm, 100 pieces

Grade 1288	Grade 1289	Grade 1290	Grade 1291	Grade 1292	Grade 293
FT-2-206-580580	FT-2-207-580580	FT-2-208-580580	FT-2-209-580580	FT-2-210-580580	FT-2-211-580580

Other dimensions are available on request

High-Purity Filter Papers for Qualitative Analyses

These paper grades are used for analytical purposes that require a low ash content. Grades 292 and 292a are especially suitable for soil analyses because they are low in nitrogen. For phosphate or sodium determination, we recommend grades 131 and 132.



Application Examples

Application	Grade
Malt filtration according to EBC standards	292
Determination of nitrogen content in soils	292, 292a
Determination of phosphate and sodium content in soils	131, 132

- Pure cotton linters or cotton linters with refined pulp
- No additives, such as wet-strengthening agents
- Supplied in rolls, sheets, discs and folded filters

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Particle retention (µm)	Filtration (s)*	Material
292	87	0.18	5–8	45	Cotton linters, low-nitrogen and nitrates, ash content ≤ 0.06% according to DIN 54370
292a	97	0.19	4–7	60	Cotton linters, low-nitrogen and nitrates, ash content ≤ 0.06% according to DIN 54370
132	80	0.17	5–7	55	Cotton linters and refined pulp, low-phosphate and low-potassium, ash content < 0.02% according to DIN 54370
131	80	0.16	3–5	100	Cotton linters and refined pulp, low-phosphate and low-potassium, ash content < 0.02% according to DIN 54370

* See test methods, page 35

Ordering Information



Filter Discs, 100 pieces

Ø in mm	Grade 131	Grade 132	Grade 292	Grade 292a
55		FT-3-329-055	FT-3-205-055	FT-3-215-055
70		FT-3-329-070	FT-3-205-070	FT-3-215-070
90		FT-3-329-090	FT-3-205-090	FT-3-215-090
110		FT-3-329-110	FT-3-205-110	FT-3-215-110
125	FT-3-351-125	FT-3-329-125	FT-3-205-125	FT-3-215-125
150		FT-3-329-150	FT-3-205-150	FT-3-215-150
185		FT-3-329-185	FT-3-205-185	FT-3-215-185
240		FT-3-329-240	FT-3-205-240	FT-3-215-240



Folded Filters, 100 pieces

Ø in mm	Grade 131	Grade 132	Grade 292	Grade 292a
110	FT-4-351-110	FT-4-329-110	FT-4-205-110	FT-4-215-110
125	FT-4-351-125	FT-4-329-125	FT-4-205-125	FT-4-215-125
150	FT-4-351-150	FT-4-329-150	FT-4-205-150	FT-4-215-150
185	FT-4-351-185	FT-4-329-185	FT-4-205-185	FT-4-215-185
240		FT-4-329-240	FT-4-205-240	FT-4-215-240



Sheets in 580 × 580 mm, 100 pieces

Grade 292	Grade 292a
FT-2-205-580580	FT-2-215-580580

Other dimensions are available on request

Smooth Filter Papers for Qualitative & Technical Analyses

These filter papers are used for routine analyses like clarification, determination of substances, but also as discs with a center hole for technical applications. Grades with a wet burst resistance > 30 kPa are referred to as wet-strengthened and are therefore suitable for pressure or vacuum filtration. White and bright particles can be easily detected with the black paper grade 918, due to the color contrast for example for the detection of fluorine or silicon in water or the detection of mycelium in mildews.

Application Examples

Application	Grade
Routine work in the lab	3 hw
Degassing beer before analysis	6
Determination of the sugar content	100/N
Clarification of clear or dyed liquids	3 m/N
Water Absorption test for mortar according to EN 1015-18	3 S/h
Durum wheat flour and semolina – Determination of yellow pigment content (ISO 11052:1994)	918
Clarification and brightening of dull and dark urines	69 k
Polarimetric sugar determination	69 k



- Made of refined pulp and cotton linters with an > 95% α -cellulose content
- Ash content between 0.1 – 0.15% (grade 100/N < 0.1%)
- Wet-strengthened
- Supplied in rolls, sheets, discs and folded filters as well as customer-specific cuts

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Filtration (s)*	Particle Retention (μ m)	Wet burst resistance (kPa)*	Properties
10	120	0.33	8	11 – 16	30	Fast filtering, wide-pore
6	80	0.17	15	10 – 13	30	Fast filtering
3 w	65	0.14	15	9 – 13	15	Medium fast filtering
3 hw	65	0.14	20	8 – 12	40	Medium fast filtering
C 140	140	0.30	20	7 – 11	> 50	Medium fast filtering
4 b	75	0.15	22	8 – 12	30	Medium fast filtering
3 m/N	65	0.14	30	7 – 10	30	Medium fast filtering
50 S	120	0.22	30	6 – 10	20	Medium fast filtering
100/N	85	0.18	30	6 – 8	80	Medium fast filtering, low ammonium, potassium & sodium content
3 h	65	0.13	35	6 – 9	15	Medium fast to slow filtering
918	85	0.17	45	8 – 10		Medium fast to slow filtering, black paper
460/N	90	0.17	50	3 – 4	30	Medium fast to slow filtering
3 S/h	200	0.36	55	5 – 7	15	Medium fast to slow filtering, narrow-pore
69 k	155	0.36	65	7 – 9	20	Slow filtering, narrow-pore, black activated charcoal paper

* See test methods, page 35

Ordering Information



Filter Discs

Ø in mm	Grade 10 (50 pieces)	Grade 100/N (100 pieces)	Grade 3 h (100 pieces)	Grade 3 hw (100 pieces)	Grade 3 m/N (100 pieces)	Grade 3 S/h (50 pieces)
55	FT-3-352-055	FT-3-328-055	FT-3-302-055	FT-3-303-055	FT-3-305-055	FT-3-307-055
70	FT-3-352-070	FT-3-328-070	FT-3-302-070	FT-3-303-070	FT-3-305-070	FT-3-307-070
90	FT-3-352-090	FT-3-328-090	FT-3-302-090	FT-3-303-090	FT-3-305-090	FT-3-307-090
110	FT-3-352-110	FT-3-328-110	FT-3-302-110	FT-3-303-110	FT-3-305-110	FT-3-307-110
125	FT-3-352-125	FT-3-328-125	FT-3-302-125	FT-3-303-125	FT-3-305-125	FT-3-307-125
150	FT-3-352-150	FT-3-328-150	FT-3-302-150	FT-3-303-150	FT-3-305-150	FT-3-307-150
185	FT-3-352-185	FT-3-328-185	FT-3-302-185	FT-3-303-185	FT-3-305-185	FT-3-307-185
240	FT-3-352-240	FT-3-328-240	FT-3-302-240	FT-3-303-240	FT-3-305-240	FT-3-307-240

Ø in mm	Grade 3 w (100 pieces)	Grade 4 b (100 pieces)	Grade 6 (100 pieces)	Grade 69 k (100 pieces)	Grade 918 (100 pieces)	Grade C 140 (50 pieces)
55	FT-3-308-055	FT-3-309-055	FT-3-312-055		FT-3-607-055	
70	FT-3-308-070	FT-3-309-070	FT-3-312-070	FT-3-326-070		
90	FT-3-308-090	FT-3-309-090	FT-3-312-090	FT-3-326-090	FT-3-607-090	FT-3-356-090
110	FT-3-308-110	FT-3-309-110	FT-3-312-110	FT-3-326-110		
125	FT-3-308-125	FT-3-309-125	FT-3-312-125	FT-3-326-125		
150	FT-3-308-150	FT-3-309-150	FT-3-312-150	FT-3-326-150		
185	FT-3-308-185	FT-3-309-185	FT-3-312-185	FT-3-326-185		FT-3-356-185
240	FT-3-308-240	FT-3-309-240	FT-3-312-240			



Folded Filters, 100 pieces

Ø in mm	Grade 10	Grade 100/N	Grade 3 h	Grade 3 hw	Grade 3 m/N
110	FT-4-352-110		FT-4-302-110	FT-4-303-110	FT-4-305-110
125	FT-4-352-125		FT-4-302-125	FT-4-303-125	FT-4-305-125
150	FT-4-352-150	FT-4-328-150	FT-4-302-150	FT-4-303-150	FT-4-305-150
185	FT-4-352-185		FT-4-302-185	FT-4-303-185	FT-4-305-185
240	FT-4-352-240	FT-4-328-240	FT-4-302-240	FT-4-303-240	FT-4-305-240
270	FT-4-352-270	FT-4-328-270	FT-4-302-270	FT-4-303-270	FT-4-305-270
320	FT-4-352-320	FT-4-328-320	FT-4-302-320	FT-4-303-320	FT-4-305-320
385	FT-4-352-385		FT-4-302-385	FT-4-303-385	FT-4-305-385

Ø in mm	Grade 3 S/h	Grade 3 w	Grade 4 b	Grade 6	Grade C 140
110		FT-4-308-110	FT-4-309-110	FT-4-312-110	FT-4-356-110
125		FT-4-308-125	FT-4-309-125	FT-4-312-125	FT-4-356-125
150		FT-4-308-150	FT-4-309-150	FT-4-312-150	FT-4-356-150
185		FT-4-308-185	FT-4-309-185	FT-4-312-185	FT-4-356-185
240	FT-4-307-240	FT-4-308-240	FT-4-309-240	FT-4-312-240	FT-4-356-240
270	FT-4-307-270	FT-4-308-270	FT-4-309-270	FT-4-312-270	FT-4-356-270
320	FT-4-307-320	FT-4-308-320	FT-4-309-320	FT-4-312-320	FT-4-356-320
385		FT-4-308-385	FT-4-309-385	FT-4-312-385	



Sheets in 580 × 580 mm, 100 pieces

Grade 10	Grade 100/N	Grade 3 h	Grade 3 hw	Grade 3 m/N
FT-2-352-580580	FT-2-328-580580	FT-2-302-580580	FT-2-303-580580	FT-2-305-580580

Grade 3 S/h	Grade 3 w	Grade 4 b	Grade 460/N	Grade 6
FT-2-307-580580	FT-2-308-580580	FT-2-309-580580	FT-2-332-580580	FT-2-312-580580

Other dimensions are available on request

Crêped Filter Papers for Qualitative & Technical Analyses

Crêped filter papers are mostly used for the rapid filtration of relatively coarse precipitates; because of their crêped structure they provide a larger filtration area than smooth filter papers. Grades with a wet burst resistance > 30 kPa are referred to as wet-strengthened and are therefore suitable for pressure or vacuum filtration. Below you will find an overview of the most commonly used grades.



Application Examples

Application	Grade
Cooking oils	39/N
Vinegar filtration	39/N
Fine filtration of tar	17/N
Galvanic baths	34/N
Prefilters for transformer oil	6 S/N
Filtration of sugar solutions	603/N
Filtration of essential oils	5 H/N
Juice filtration	67/N

- Made of refined pulp and cotton linters with an > 95% α -cellulose content
- Ash content between 0.1 – 0.15%
- Wet-strengthened
- Supplied in rolls, sheets, discs and folded filters as well as customer-specific cuts

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Filtration (s)*	Wet burst resistance (kPa)*	Air resistance (mbar)*	Properties
5 H/N	85	0.28	3	≥ 40		Very fast filtering, wide-pore
34/N	60	0.20	4	40	1.5	Very fast filtering
37/N	135	0.50	4	≥ 70	1.9	Very fast filtering, wide-pore
FT 55	55	0.15	5	20		Very fast filtering
1602/N	70	0.23	5	≥ 30		Very fast filtering
39/N	180	0.65	5	90	2.5	Very fast filtering, wide-pore
39/N	300	0.95	5	120	2.5	Very fast filtering, wide-pore
603/N	75	0.25	8	≥ 50		Fast filtering
6 S/N	145	0.55	12	≥ 90		Medium fast filtering
601/N	65	0.19	13	≥ 30		Medium fast filtering
67/N	160	0.65	13	≥ 70	5.5	Medium fast filtering
17/N	90	0.30	20	≥ 30		Medium filtering

* See test methods, page 35

Ordering Information



Filter Discs

Ø in mm	Grade 5 H/N (100 pieces)	Grade 6 S/N (50 pieces)	Grade 601/N (100 pieces)	Grade 603/N (100 pieces)	Grade 37/N (50 pieces)	Grade 39/N, 180 g/m ² (50 pieces)
47	FT-3-423-047				FT-3-480-047	
70		FT-3-314-070			FT-3-480-070	
90	FT-3-423-090	FT-3-314-090		FT-3-335-090	FT-3-480-090	
110	FT-3-423-110	FT-3-314-110	FT-3-354-110	FT-3-335-110	FT-3-480-110	FT-3-483-110
125	FT-3-423-125	FT-3-314-125	FT-3-354-125	FT-3-335-125	FT-3-480-125	
150	FT-3-423-150	FT-3-314-150	FT-3-354-150	FT-3-335-150	FT-3-480-150	
185	FT-3-423-185	FT-3-314-185	FT-3-354-185	FT-3-335-185	FT-3-480-185	FT-3-483-185
240	FT-3-423-240	FT-3-314-240	FT-3-354-240	FT-3-335-240	FT-3-480-240	
320			FT-3-354-320	FT-3-335-320		



Folded Filters, 100 pieces

Ø in mm	Grade 5 H/N	Grade 6 S/N	Grade 603/N	Grade 34/N	Grade 37/N	Grade 39/N, 180 g/m ²
125	FT-4-423-125	FT-4-314-125	FT-4-335-125	FT-4-478-125	FT-4-480-125	
150	FT-4-423-150	FT-4-314-150	FT-4-335-150		FT-4-480-150	FT-4-483-150
185	FT-4-423-185	FT-4-314-185	FT-4-335-185		FT-4-480-185	FT-4-483-185
240	FT-4-423-240	FT-4-314-240	FT-4-335-240		FT-4-480-240	FT-4-483-240
270	FT-4-423-270	FT-4-314-270	FT-4-335-270			
320	FT-4-423-320	FT-4-314-320	FT-4-335-320	FT-4-478-320	FT-4-480-320	
385	FT-4-423-385					FT-4-483-385
500	FT-4-423-500	FT-4-314-500			FT-4-480-500	FT-4-483-500



Sheets in 580 × 580 mm, 100 pieces

Grade 5 H/N	Grade 6 S/N	Grade 601/N	Grade 603/N
FT-2-423-580580	FT-2-314-580580	FT-2-354-580580	FT-2-335-580580
Grade 17/N	Grade 37/N	Grade 39/N, 180 g/m ²	Grade FT 55
FT-2-321-580580	FT-2-480-580580	FT-2-483-580580	FT-2-348-580580

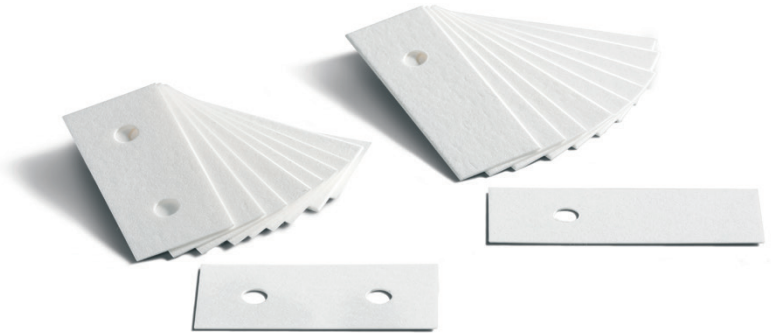
Other dimensions are available on request

Paper Boards for the Filtration and Absorption of Liquids

Among other applications, these boards are used for the filtration of cooking and transformer oils, galvanic baths and as base paper for further impregnation with certain reagents. Grades with a wet burst resistance > 30 kPa are referred to as wet-strengthened and are therefore suitable for pressure or vacuum filtration.



- Made of refined pulp or cotton linters
- Smooth
- Supplied in sheets, discs and as well as customer-specific cuts



Application Examples

Application	Grade
Cytocards	151
Transformer oils	C 250
Clarifications of galvanic baths	C 350
Blotting paper for determination of water absorptiveness according to COBB (EN 20535)	A 250
Cooking oil	C 251
Paper air fresheners	157
Fragrance test cards	C 160

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Filtration (s)*	Air resistance (mbar)*	Capillary rise (mm/10 min)*	Dry burst resistance (kPa)*	Wet burst resistance (kPa)*	Water capacity (%)
C 160	160	0.30	40	25	80		≥ 50	
S 165	165	0.34		35	≥ 40			
C 250	250	0.48	40	25	100		≥ 80	
C 251	250	> 0.55	8	3.5	180	≥ 200	≥ 80	
A 250	250	0.45			70			180
C 300	300	0.55	40	25	100		≥ 80	
1339	315	0.63		42	60	≥ 500	≥ 230	
C 350	350	0.63	40	25	110		≥ 80	
C 350L	360	0.75		30	80		≥ 200	
C 450	450	0.95	40	25	110		≥ 50	
151	460	1.00		19	120	≥ 400		
1220	475	1.10	200		120			
K12	520	1.60		2.5	200	≥ 250	≥ 50	
157	700	1.80		8	150			
SEK 770	770	1.00						400

* See test methods, page 35

Ordering Information



Sheets in 580 × 580 mm, 100 pieces

Grade C 160	Grade C 250	Grade C 251	Grade C 300
FT-2-343-580580	FT-2-344-580580	FT-2-355-580580	FT-2-345-580580
Grade C 350	Grade C 450	Grade S 165	
FT-2-346-580580	FT-2-347-580580	FT-2-368-580580	

Other dimensions are available on request

Seed Testing Papers

These papers satisfy the requirements for the determination of germination capability according to ISTA (International Seed Testing Association) and are ideal for ensuring optimal moisture content for the most diverse types of seeds and germination forms. Their pH ranges between 6.0 and 7.5, they are wet-strengthened and their special structure prevents fine seed roots from growing through the paper. The colored papers are produced with dyes that do not influence the growth of roots. These papers are mainly used to count more easily very fine and white roots.



Application Examples

This method is mainly applied with corn, sugar beets, wheat, barley and various grasses, but can also be used for all other seed types.

PP ("Pleated Paper") Method

The pleated paper is placed in a box; the seeds are distributed among the folds of the pleated paper and covered with a wrapping strip to keep the seeds moist. The pleated papers have 50 double folds that are 20 mm in depth; usually, 2 seeds are placed in each fold. Both white and grey papers are available. Colored paper makes it easier to count white seed species.

Technical Specifications & Ordering Information

Grade	Properties	Weight (g/m ²)*	Thickness (mm)*	Size (mm)	Qty per box	Order No.
20	Pleated strips, white	110	0.22	2,000 × 110	1,008**	FT-2003532000110
20, grey	Pleated strips, grey	110	0.22	2,000 × 110	1,008**	FT-2003662000110
4 b	Wrapping strips	75	0.15	110 × 580	100	FT-2-309-110580
6	Wrapping strips	80	0.17	110 × 580	100	FT-2-312-110580

* See test methods, page 35

** 112 rods à 9 pleated strips

BP ("Between Paper") Method

One wetted paper sheet is laid on top of a second, the seeds are placed on the double sheet which is then rolled up.

Application Examples

The method is used for peas and oats, among others.

Technical Specifications & Ordering Information

Grade	Properties	Weight (g/m ²)*	Thickness (mm)*	Size (mm)	Qty per box	Order No.
1750	Sheets, white	90	0.20	220 × 400	500	FT-210607-220400
1755	PE-coated sheets	82	0.19	190 × 400	500	FT-210608-190400
39/N	Crêped white paper	180	0.65	580 × 580	100	FT-2-483-580580

* See test methods, page 35



TP ("Top of Paper") Method

The seeds are placed on the paper (discs or sheets) and then transferred either to petri dishes or plastic boxes. By supplying the filter with water, wick papers are used for constant moistening during the Jacobsen method. They are also supplied as blue and yellow papers to make it easier to count white seed species.

Application Examples

The method is applied to small seeds like clover species, for example.

Technical Specifications & Ordering Information

Grade	Properties	Weight (g/m ²)*	Thickness (mm)*	Size (mm)	Qty per box	Order No.
94/N	Smooth, white wick paper	100	0.20	30 × 60	100	FT-210374-030160
50 S	Smooth, white filter discs	120	0.22	95	50	FT-3-353-095
50 S	Smooth white paper sheets	120	0.22	225 × 225	100	FT-2-353-225225
C 140	Smooth white paper	140	0.30	240 × 400	100	FT-2-356-240400
6 S/N	Crêped white paper	145	0.55	150 × 580	100	FT-2-314-150580
193	Smooth, yellow paper sheets	160	0.32	120 × 300	100	FT-2-381-120300
193	Smooth, yellow paper sheets	160	0.32	110 × 170	1000	FT-2-381-110170
190	Smooth, blue wick paper	300	0.65	25 × 133	100	FT-2-378-025133
190	Smooth, blue paper sheets	300	0.65	90 × 133	100	FT-2-378-090133
194	Smooth, dark blue paper	430	0.68	200 × 350	100	FT-3-383-200350
191	Smooth, blue paper	700	1.35	140 × 200	100	FT-2-379-140200
192	Smooth, dark blue paper	720	1.45			

Other paper grades and dimensions are available on request

* See test methods, page 35

Filter Papers for the Sugar Industry

In the sugar industry, filter papers are used in laboratories to assay sugar beet or cane sugar. The sugar beets are mashed and further analyzed according to the aluminum sulfate method. Potassium, nitrogen, sodium and saccharose content are measured using a spectrophotometer or the likes. These papers are wet-strengthened and either smooth or crêped. They are made of cellulose or a mixture of cellulose and diatomaceous earth.

Grade 100/N is not only supplied as discs or folded filters, but also on rolls for VENEMA systems.

Technical Specifications

Grade	Properties	Weight (g/m ²)*	Thickness (mm)*	Filtration (s)*	Wet burst resistance (kPa)*	Order No.
603/N	Crêped paper, very fast filtering	75	0.25	8	≥ 50	See page 15
6 S/N	Crêped paper, very fast filtering	145	0.55	12	≥ 90	See page 15
601/N	Crêped paper, fast filtering	65	0.19	13	≥ 30	See page 15
3 hw	Smooth paper, medium fast filtering	65	0.14	20	40	See page 13
470	Diatomaceous earth filter paper, slow filtering	140	0.32	80	30	See page 21
100/N	Smooth paper, medium fast filtering, low phosphate and low sodium	85	0.18	30	80	See below on rolls or page 13 as sheets, discs or folded filters

* See test methods, page 35

Ordering Information

Venema Rolls, Grade 100/N

Width	Length	Qty per box	Order No.
150 mm	1,000 m	1 roll	FT-1-328-1501000
240 mm	1,000 m	1 roll	FT-1-328-2401000



Diatomaceous Earth Filter Paper

Grade 470 papers are made of cellulose and diatomaceous earth and offer a much better separating capability than pure cellulose papers at the same rate of filtration. This grade quickly retains the finest particles at high flow rates.



Application Examples

Clarification of beer, wine, urine during spectrophotometric or refractometric tests

Filtration of the finest, semi-colloidal precipitates, e.g. those consisting of proteins, clay or cold-precipitated barium

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Filtration (s)*
470	140	0.32	80

* See test methods, page 35

Ordering Information



Filter Discs, 100 pieces

Ø in mm	Order No.
90	FT-3-606-090
110	FT-3-606-110
125	FT-3-606-125
150	FT-3-606-150
185	FT-3-606-185



Folded Filters, 100 pieces

Ø in mm	Order No.
125	FT-4-606-125
150	FT-4-606-150
185	FT-4-606-185
240	FT-4-606-240
320	FT-4-606-320

Phase Separating Paper

Grade 480 is impregnated with stabilized silicon, thus rendering it hydrophobic: It retains water, but allows solvents to flow through. The flow stops automatically when the entire solvent has passed through. In many applications, this phase separator paper eliminates the need to use separating funnels.

Application Examples

Filtration of organic solvents that have been contaminated with water to separate the aqueous from the organic phase

Filtration of extracting solvents in clinical or medical labs

Separation of emulsions that are formed during the extraction of aqueous plant or drug solutions



Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*
480	85	0.19

* See test methods, page 35

Ordering Information



Filter Discs, 100 pieces

Ø in mm	Order No.
70	FT-3-602-070
90	FT-3-602-090
110	FT-3-602-110
125	FT-3-602-125
150	FT-3-602-150
185	FT-3-602-185



Folded Filters, 100 pieces

Ø in mm	Order No.
90	FT-4-602-090
125	FT-4-602-125
150	FT-4-602-150
185	FT-4-602-185
270	FT-4-602-270

Nonwovens

These nonwoven grades are made of rayon or polyester and are available in different weights. They can be used for the filtration or prefiltration of viscous solutions containing particles visible with the naked eye.



Application Examples

Testing of mortars containing mineral binders; determination of water retentivity of freshly mixed mortar by the filter plate method according to DIN 18555-7 : FT-3-01308-185

Food-processing industry: Detection of contaminants in dairy products, baby food : FT-3-01324-032, FT-3-01324-047

Petrochemical industry: Filtration of viscous solutions to retain the biggest particles before further analysis with a viscosimetry: FT-3-01304-025

Metal-processing industry: Non-woven belt filters (rolls) for the removal of particles from liquids.

Technical Specifications

Grade	Material
2601	Rayon (viscose) supplied in pore sizes from 75 to 115 μm
2701	Polyester supplied in pore sizes from 60 to 140 μm
2602	Rayon (viscose) Polyester

Ordering Information

Grade, Weight	Thickness (mm)*	Mean pore size (μm)	Max. pore size (μm)	Format	Size	Qty per box	Order No.
2601, 20 g/m ²	0.19	75	200	Filter discs	185 mm	100	FT-3-01308-185
2601, 60 g/m ²	0.46	50	95	Filter discs	25 mm	100	FT-3-01304-025
2601, 60 g/m ²	0.46	50	95	Filter discs	47 mm	100	FT-3-01304-047
2701, 60 g/m ²	0.20	60	80	Roll	380 mm x 40 m	1	FT-1-01323-38040
2602, 150 g/m ²	0.80	-	-	Filter discs	25 mm	1,000	FT-3-01324-025
2602, 150 g/m ²	0.80	-	-	Filter discs	32 mm	1,000	FT-3-01324-032
2602, 150 g/m ²	0.80	-	-	Filter discs	47 mm	1,000	FT-3-01324-047

Weighing Paper

Grade 605 weighing paper is made of transparent smooth parchment that is ideal for the weighing of viscous, semi-crystalline or solid substances.



- Keeps the balance pan clean
- Allows easy transfer of weighed substances
- Nitrogen-free, smooth, transparent and soluble
- Supplied in sheets, discs as well as weighing boats

Technical Specifications & Ordering Information

Grade	Material	Weight (g/m ²)*	Thickness (mm)*	Format	Size	Qty per box	Order No.
605	Nitrogen-free parchment paper	20	0.02	Sheets	100 × 100 mm	1,000	FT-2-605-100100R
605	Nitrogen-free parchment paper	20	0.02	Boats	58 × 10 × 10 mm	100	FT-2-605-581010

* See test methods, page 35

Lens Cleaning Paper

Grade 2113 lens cleaning paper is a thin, non-linting silk tissue paper used for cleaning very sensitive surfaces, such as optical glasses or lenses without scratching them.



- Prevents scratching of surfaces
- Supplied in sheets as well as customer-specific cuts

Technical Specifications & Ordering Information

Grade	Material	Weight (g/m ²)*	Format	Size	Qty per box	Order No.
2113	Non-linting silk paper	13	Sheets	120 × 120 mm	500	FT-2-1111-120120

* See test methods, page 35

Surface Protection Paper

LabSorb is a highly absorptive grade of paper coated on one side with polyethylene. Used with the cellulose side up, the paper absorbs liquids, which are stopped by the polyethylene layer and thus prevented from soaking through. Used with the polyethylene side up, the paper is highly useful for recovery of valuable or toxic liquids.



Technical Specifications

- Weight: 140 g/m²
- Water capacity: 140 %

Application Examples

Preventing radioactive contamination of work surfaces in radiochemical laboratories

Recovering spilled solutions containing expensive reagents

Protecting laboratory bench surfaces from spillage or splashes of liquids by preventing absorption and seepage of these liquids into work surfaces

Lining animal cages for protection and hygiene

Reducing the risk of objects breaking after falling on hard surfaces

Ordering Information

Grade	Format	Size	Qty per box	Order No.
LabSorb	Roll	400 mm × 50 m	1	FT-1-601-400050
LabSorb	Roll	400 mm × 100 m	1	FT-1-601-400100
LabSorb	Roll	460 mm × 50 m	1	FT-1-601-460050
LabSorb	Roll	600 mm × 50 m	1	FT-1-601-600050
LabSorb	Roll	600 mm × 100 m	1	FT-1-601-600100
LabSorb	Sheets	460 × 570 mm	50	FT-2-601-460570K
LabSorb	Sheets	480 × 600 mm	50	FT-2-601-480600K

Blotting Papers

These blotting papers are made from the purest raw materials with the maximum degree of absorptiveness and cellulose content. They are available in a choice of different weights and thicknesses as well as in different formats to cover the majority of blotting applications. Furthermore, they are the ideal complement to the Sartorius nitrocellulose blotting membranes available in two pore sizes, 0.22 μm and 0.45 μm .



Application Examples

Application	Grade
To cover the gel membrane sandwich in the buffer tank	BF 1
For gel wicking and drying, capillary blotting using Western, Southern or semidry techniques	BF 2
To increase and maintain the transport of liquid from the buffer and as buffer reservoir in capillary and semidry blotting methods	BF 3
To transfer DNA or RNA according to the Southern technique or semidry blotting of proteins	BF 4

- 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

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Capillary rise (mm/10 min)*	Capillary rise (mm/30 min)*
BF 1	90	0.16	80	140
BF 2	195	0.35	70	115
BF 3	330	0.76	130	
BF 4	550	1.30	160	

* See test methods, page 35

Ordering Information



Sheets

Grade	Size (in mm)	Qty per box	Order No.
BF 1	580 × 600	100	FT-2-518-560600N
BF 2	80 × 90	100	FT-2-519-080090N
BF 2	130 × 210	100	FT-2-519-130210N
BF 2	200 × 200	100	FT-2-519-200200N
BF 2	460 × 570	100	FT-2-519-460570N
BF 2	580 × 600	100	FT-2-519-580600N
BF 3	135 × 155	100	FT-2-520-135155N
BF 3	200 × 200	100	FT-2-520-200200N
BF 3	460 × 570	50	FT-2-520-460570K
BF 3	580 × 600	50	FT-2-520-580600K
BF 4	110 × 170	25	FT-2-521-110170G
BF 4	150 × 150	25	FT-2-521-150150G
BF 4	580 × 580	25	FT-2-521-580580G
BF 4	580 × 600	25	FT-2-521-580600G

Other dimensions are available on request

Blotting Membranes

Sartorius blotting membranes are ideal as a complement to the blotting papers 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.



- 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

Technical Specifications

Description	11327	11306
Material	Cellulose Nitrate	Cellulose Nitrate
Pore size (µm)	0.22	0.45
Thickness (µm)	120	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 (µg/cm ²)	200	200

Ordering Information



Rolls

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

Chromatography Papers

Chromatography papers are made of 100% cotton linters. These highly pure papers are not only ideal for blotting & chromatography, but also for a wide range of absorption applications like those common in the life sciences and diagnostics.



Application Examples

Application	Grade
The most commonly used chromatography paper	FN 100
Analytical paper for routine and repetitive separations	FN 1
Routine analysis of proteins in serum (e.g. human albumin)	FN 3
Blotting paper in analysis sets	FN 8
Antibiotic test strips	FN 30
Testing of mortars containing mineral binders; determination of water retentivity of freshly mixed mortar by the filter plate method according to DIN 18555-7: FT-210501-190190	27 CH

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Capillary rise (mm/30 min)*
FN 1	90	0.19	145
FN 2	125	0.24	145
FN 3	90	0.19	95
FN 4	125	0.24	95
FN 5	90	0.18	60
FN 7	150	0.32	145
FN 7a	200	0.41	145
FN 8	280	0.55	170
FN 30	320	0.90	240
FN 100	195	0.35	115
27 CH	700	1.3	170

* See test methods, page 35

Ordering Information



Sheets

Grade	Size (in mm)	Qty per box	Order No.
FN 1	400 × 400	100	FT-2-501-400400N
FN 1	470 × 580	100	FT-2-501-470580N
FN 1	580 × 600	100	FT-2-501-580600N
FN 2	580 × 600	100	FT-2-502-580600N
FN 3	300 × 580	100	FT-2-503-300580N
FN 3	460 × 570	100	FT-2-503-460570N
FN 3	580 × 600	100	FT-2-503-580600N
FN 4	580 × 600	100	FT-2-504-580600N
FN 5	200 × 200	100	FT-2-505-200200N
FN 5	580 × 600	100	FT-2-505-580600N
FN 7	460 × 570	50	FT-2-507-460570K
FN 7	580 × 600	50	FT-2-507-580600K
FN 7a	460 × 570	50	FT-2-508-460570K
FN 7a	580 × 600	50	FT-2-508-580600K
FN 8	470 × 580	50	FT-2-509-460580K
FN 8	580 × 600	50	FT-2-509-580600K
FN 30	254 × 305	100	FT-2-526-254305N
FN 30	580 × 600	25	FT-2-526-580600G
FN 100	76 × 102	100	FT-2-527-076102N
FN 100	200 × 200	100	FT-2-527-200200N
FN 100	260 × 410	100	FT-2-527-260410N
FN 100	460 × 570	50	FT-2-527-460570K
FN 100	460 × 570	100	FT-2-520-460570N
FN 100	580 × 600	50	FT-2-527-580600K
FN 100	580 × 680	50	FT-2-527-580680K

Other dimensions are available on request

Glass Microfiber Filters With Binder

These filters are mostly used either for monitoring air and gas or as prefilter. They are manufactured with synthetic binding agents to ensure that the filter has a defined strength. They are mechanically and chemically stable, have a temperature resistance up to 180°C and – depending on the binding agent used – are either hydrophobic or hydrophilic.



Application Examples

Application	Grade
Prefiltration	13400, MG 1387/1
Smoke number measurement	MG 227
Gas monitoring	MG 1387/1
Air pump protection	MG 227/1/60

- Mechanically and chemically stable
- Temperature resistant up to 180°C
- Supplied as discs or sheets

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Penetration 0.3 µm (%)	Pressure drop 5.3 cm/s (Pa)	Binding agent
MG 227/1/60	60	0.32	< 0.5	260	Hydrophobic
13430	220	1.25	0.02	360	Hydrophilic
13400	73	0.39	0.015	363	Hydrophilic
MG 227	75	0.40	< 0.01	350	Hydrophobic
MG 400	80	0.38	< 0.001	400	Hydrophilic
MG 1387/1	90	0.37	≤ 0.003	400	Hydrophilic

* See test methods, page 35

Ordering Information



Filter Discs

Ø in mm	MG 227/1/60 (100 pieces)	13430**	13400**	MG 227 (100 pieces)	MG 1387/1 (50 pieces)
13			13400--13-----S		
16			13400--16-----S		
20			13400--20-----S		
25			13400--25-----Q		
42			13400--42-----Q		
44			13400--44-----Q		
45			13400--45-----Q		FT-3-01125-045
47		13430--47-----S	13400--47-----Q	FT-3-01120-047	FT-3-01125-047
50			13400--50-----Q		FT-3-01125-050
55				FT-3-01120-055	FT-3-01125-055
80			13400--80-----N		
100		13430-100-----K	13400-100-----K		
110			13400-110-----K	FT-3-01120-110	FT-3-01125-110
120			13400-120-----K		
124			13400-124-----K		
125					FT-3-01125-125
127		13430-127-----K	13400-127-----K		
130		13430-130-----K	13400-130-----K		FT-3-01125-130
150	FT-3-01124-150				

Other dimensions are available on request

** K= 50 pieces,
N= 100 pieces,
Q = 500 pieces,
S= 200 pieces



Glass Microfiber Filters Without Binder

Binder-free glass microfiber filters are recommended for analytical and gravimetric analyses and also as prefilters. These filters combine fast flow rates with high load capacity and the retention of very fine particles; they are biologically inert, are resistant to most chemicals and withstand temperatures up to 500 °C (grade 550-HA & MG 169 up to 550 °C).



- Manufactured from 100% borosilicate glass
- 100% binder free
- pH stable
- Withstand temperatures up to 500 °C (Grade MG 550-HA & MG 169 up to 550 °C)
- Supplied as discs or sheets

Application Examples

Application	Grade
Prefiltration	13440, MGB, MGD
Analysis of suspended solids in wastewater according to EN 872	MGC
Analysis of suspended solids in wastewater according to 2540D	MG 550-HA, MG 169
Clarification of buffer & reagent solutions	MGA
Clarification of protein solutions	MGF
Air Monitoring, PM10	MG 160

Technical Specifications

Grade	Weight (g/m ²)*	Thickness (mm)*	Penetration 0.3 µm (%)**	Particle retention in liquids (µm)	Filtration speed (mL/min)*	Fulfills the requirements in EN 872:2005 (weigh loss)
MGA	54	0.23	< 0.001	1.6	510	yes
MGB	143	0.70	< 0.001	1.0	210	
MGC	52	0.26	< 0.001	1.2	335	yes
MGD	120	0.53	< 0.1	2.7	920	
MGF	75	0.38	< 0.001	0.7	110	
MGG	64	0.28	≤ 0.001	1.5	600	
13440	88	0.44		0.7	120	yes
MG 160	75	0.35	0.002	1.2	400	
MG 550-HA	65	0.27		1.5	400	

* See test methods, page 35

** Measurement according to EN 143 (0.3 µm, 5.3 cm/s, paraffin oil)

Ordering Information



Filter Discs

Ø in mm	MGA (100 pieces)	MG 160 (50 pieces)	MGB (50 pieces)	MGC (100 pieces)	MGD (50 pieces)
13					FT-3-1104-013
20	FT-3-1101-020				
21			FT-3-1102-021	FT-3-1103-021	
25	FT-3-1101-025		FT-3-1102-025	FT-3-1103-025	FT-3-1104-025
37	FT-3-1101-037	FT-3-01110-037			
47	FT-3-1101-047	FT-3-01110-047	FT-3-1102-047	FT-3-1103-047	FT-3-1104-047
50	FT-3-1101-050	FT-3-01110-050	FT-3-1102-050	FT-3-1103-050	FT-3-1104-050
55	FT-3-1101-055		FT-3-1102-055	FT-3-1103-055	
70	FT-3-1101-070	FT-3-01110-070	FT-3-1102-070	FT-3-1103-070	FT-3-1104-070
80	FT-3-1101-080				
90	FT-3-1101-090	FT-3-01110-090	FT-3-1102-090	FT-3-1103-090	FT-3-1104-090
100	FT-3-1101-100	FT-3-01110-100	FT-3-1102-100	FT-3-1103-100	FT-3-1104-100
110	FT-3-1101-110	FT-3-01110-110	FT-3-1102-110	FT-3-1103-110	FT-3-1104-110
125	FT-3-1101-125		FT-3-1102-125	FT-3-1103-125	FT-3-1104-125
150	FT-3-1101-150		FT-3-1102-150	FT-3-1103-150	FT-3-1104-150
185	FT-3-1101-185			FT-3-1103-185	
240	FT-3-1101-240		FT-3-1102-240		FT-3-1104-240
293					FT-3-1104-293

Ø in mm	MGF (100 pieces)	MGG (100 pieces)	MG 550-HA (100 pieces)	13440***
20		FT-3-1106-020		
24			FT-3-01147-024	
25	FT-3-1105-025	FT-3-1106-025		
37		FT-3-1106-037		
42				13440--42-----Q
44				13440--44-----Q
47	FT-3-1105-047	FT-3-1106-047	FT-3-01147-047	13440--47-----Q
50	FT-3-1105-050	FT-3-1106-050	FT-3-01147-050	13440--50-----Q
55	FT-3-1105-055	FT-3-1106-055	FT-3-01147-055	
60		FT-3-1106-060		
70	FT-3-1105-070	FT-3-1106-070	FT-3-01147-070	
90	FT-3-1105-090	FT-3-1106-090	FT-3-01147-090	
100				13440-100-----K
110	FT-3-1105-110	FT-3-1106-110	FT-3-01147-110	
125	FT-3-1105-125	FT-3-1106-125	FT-3-01147-125	
130				13440-130-----K
142	FT-3-1105-142			
150	FT-3-1105-150	FT-3-1106-150		
240	FT-3-1105-240			

Other dimensions are available on request

*** Q = 500 pieces
K = 50 pieces

Pre-heated Quartz Microfiber Filters

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.

Application Examples

Analysis of dust levels according to
EN 13284 1:2001

Emission monitoring at high temperatures
(air pollution)

Analysis of hot and acidic gases

Trace element analysis

Analytical and gravimetric analyses

- High-purity filters with the lowest trace metal values
- Extreme temperature resistance up to 900°C
- Exceptional chemical resistance
- Excellent weight and dimensional stability
- Biologically inert
- Certificate on trace elements available for every batch

Technical Specifications

Grade	Material	Weight (g/m ²)*	Thickness (mm)*	Penetration 0.3 µm (%)**	Tensile strength, dry, MD (N/m)	Pressure drop 5.3 cm/s (Pa)
Q3400	100% Quartz microfiber silicium dioxide (SiO ₂)	85	0.43	< 0.002	200	450

* See test methods, page 35

** according to EN 143 (0.3 µm, 15 cm/s, paraffin oil)

Ordering Information



Filter Discs

Size (in mm)	Qty per box	Order No.
20	25	Q3400--20-----G
25	25	Q3400--25-----G
30	25	Q3400--30-----G
37	25	Q3400--37-----G
45	25	Q3400--45-----G
47	25	Q3400--47-----G
50	25	Q3400--50-----G
82	100	Q3400--82-----N
90	100	Q3400--90-----N
142	50	Q3400-142-----K
150	50	Q3400-150-----K

Other dimensions are available on request



Quality Control

Test Methods

Basis Weight According to DIN EN ISO 536

The basis weight is determined by weighing a paper sheet that is between 500 cm² and 1000 cm² in size on a calibrated paper scale showing an accuracy of +/- 0.5%. The basis weight is expressed in grams per square meter (g/m²).

Thickness According to DIN EN ISO 20534

The thickness is measured using a thickness meter or gauge readings and is expressed in millimeters.

Filtration Speed (s)

The time required to filter 10 mL of distilled water at 20 °C through a free-hanging, fully-wetted filter disc with a diameter of 110 mm folded in quarters. The filtration rate is expressed in seconds.

Filtration Speed (ml/min – Herzberg)

The time required to filter distilled water at 20 °C through a filter surface of 10 cm² and at a constant pressure of 5 cm water column. The filtration rate is expressed in ml/min.

Ash Content According to DIN 54370

The ash content is the residue determined after ignition of 10 g of filter paper at 800 °C in a platinum crucible. The ash content is expressed in percent.

Tensile Strength According to DIN EN ISO 1924-2

A continually increasing load is applied vertically to a paper strip measuring 15 mm in width and 180 mm in length. The tensile strength is defined as the stretching force necessary to break the piece and measured lengthwise and in the transverse direction. The tensile strength is expressed in N/15 mm.

Dry Bursting Strength According to DIN ISO 2758

A paper with a surface area of 10 cm² is clamped over and subjected to increasing pressure from a rubber diaphragm. The bursting strength is the pressure reading at the time of rupture. The bursting strength is expressed in kilopascal (kPa).

Wet Burst Resistance According to DIN ISO 3689

A paper with surface area of 10 cm² is immersed in water and then clamped over a rubber diaphragm. The paper is subjected to evenly increasing pressure from the rubber diaphragm. The bursting strength is the pressure reading at the time of rupture. The wet bursting strength is expressed in kilopascal (kPa).

Air Resistance

Air resistance is the pressure drop that occurs after filtration of a defined air stream (270 L/h and | or 75 cm/s at 10 cm²) through a filter paper. The air resistance is expressed in mbar.

Capillary Rise According to DIN ISO 8787 (Klemm Method)

The capillary rise is defined as the height to which a paper strip measuring 15 mm in width and 250 mm in length, whose narrow side is immersed in prefiltered distilled water (20 °C), is wetted after 10 or 30 min. After this test period of 10 and 30 min., the wetted part of the strip is measured in mm. The capillary rise is expressed in mm per 10 min and | or 30 min.

Index of Grades

Grade	Middler number	Description	Page	Grade	Middler number	Description	Page
10	352	Qualitative-technical paper, smooth	12–13	● 388	101	Quantitative filter paper	7
100/N	328	Qualitative-technical paper, smooth	12–13, 20	○ 389	102	Quantitative filter paper	7
1120, 475 g/m ²	10389	Paper board	16–17	● 389 F	112	Quantitative filter paper	7
1288	206	Qualitative filter paper	8–9	● 390	103	Quantitative filter paper	7
1289	207	Qualitative filter paper	8–9	● 391	104	Quantitative filter paper	7
1290	208	Qualitative filter paper	8–9	● 392	105	Quantitative filter paper	7
1291	209	Qualitative filter paper	8–9	● 393	127	Quantitative filter paper	7
1292	210	Qualitative filter paper	8–9	3 h	302	Qualitative-technical paper, smooth	12–13
131	351	Qualitative filter paper	10–11	3 hw	303	Qualitative-technical paper, smooth	12–13, 20
132	329	Qualitative filter paper	10–11	3 m/N	305	Qualitative-technical paper, smooth	12–13
1339	441	Paper board	17	3 S/h	307	Qualitative-technical paper, smooth	12–13
13400	n.a.	Glass microfiber filter, with binder	30–31	3 w	308	Qualitative-technical paper, smooth	12–13
13430	n.a.	Glass microfiber filter, with binder	30–31	460/N	332	Qualitative-technical paper, smooth	12–13
13440	n.a.	Glass microfiber filter, binder-free	32–33	470	606	Diatomaceous earth filter	20–21
151	449	Paper board	16–17	480	602	Phase separating paper	22
157	437	Paper board	16–17	4 b	309	Qualitative-technical paper, smooth	12–13, 18
1602/N	342	Qualitative-technical paper, créped	14	50 S	353	Germination test paper Qualitative-technical paper, smooth	12, 19
1750	10607	Germination test paper	19	5 H/N	423	Qualitative-technical paper, creped	14–15
1755	10608	Germination test paper	19	6	312	Qualitative-technical paper, smooth	12–13, 18
17/N	321	Qualitative-technical paper, créped	14–15	6 S/N	314	Qualitative-technical paper, creped	14–15, 19
190, light blue	378	Germination test paper	19	601/N	354	Qualitative-technical paper, creped	14–15, 20
191, blue	379	Germination test paper	19	603/N	335	Qualitative-technical paper, creped	14–15, 20
192, dark blue	380	Germination test paper	19	605	605	Parchment paper	24
193, yellow	381	Germination test paper	19	67/N, 160 g/m ²	477	Qualitative-technical paper, creped	14–15
194, dark blue	383	Germination test paper	19	69 K	326	Qualitative-technical paper, smooth	12–13
20	00353	Germination test paper	18	918	607	Qualitative-technical paper, smooth, black	12–13
20 grey	00366	Germination test paper	18	94/N	10374	Germination test paper Qualitative-technical paper, smooth	19
2113	1111	Lens cleaning paper	24	A 250	412	Paper board	16–17
2601, 20 g/m ²	01308	Nonwovens, viscose	23	BF 1	518	Blotting paper	26–27
2601, 60 g/m ²	01304	Nonwovens, viscose	23	BF 2	519	Blotting paper	26–27
2602, 150 g/m ²	01324	Nonwovens, viscose, polyester	23	BF 3	520	Blotting paper	26–27
2701, 60 g/m ²	01323	Nonwovens, polyester	23	BF 4	521	Blotting paper	26–27
27 CH	10501	Chromatography paper	28	C 140	356	Qualitative-technical paper, smooth	12–13, 19
292	205	Qualitative filter paper	10–11	C 160	343	Filter paper for liquid absorption	16–17
292a	215	Qualitative filter paper	10–11				
293	211	Qualitative filter paper	8–9				
34/N, 60 g/m ²	478	Qualitative-technical paper, créped	14–15				
37/N, 135 g/m ²	480	Qualitative-technical paper, créped	14–15				
39/N, 180 g/m ²	483	Qualitative-technical paper, créped	14–15, 19				
39/N, 300 g/m ²	487	Qualitative-technical paper, créped	14				

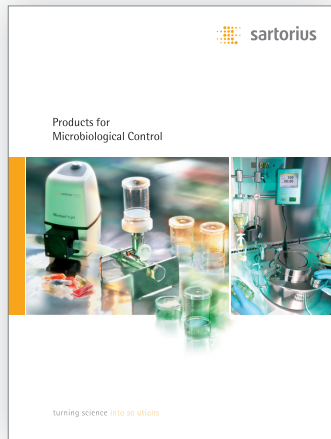
Grade	Middler number	Description	Page
C 250	344	Paper board	16 – 17
C 251	355	Paper board	16 – 17
C 300	345	Paper board	17
C 350	346	Paper board	16 – 17
C 350L	340	Paper board	17
C 450	347	Paper board	16 – 17
FN 1	501	Chromatography paper	28 – 29
FN 100	527	Chromatography paper	28 – 29
FN 2	502	Chromatography paper	28 – 29
FN 3	503	Chromatography paper	28 – 29
FN 30	526	Chromatography paper	28 – 29
FN 4	504	Chromatography paper	28 – 29
FN 5	505	Chromatography paper	28 – 29
FN 7	507	Chromatography paper	28 – 29
FN 7a	508	Chromatography paper	28 – 29
FN 8	509	Chromatography paper	28 – 29
FT 55	348	Qualitative-technical paper, creped	14 – 15
K 12	438	Paper board	17
LabSorb	601	Surface protection paper	25
MG 1387/1	01125	Glass microfiber filter, with binder	30 – 31
MG 160	01110	Glass microfiber filter, binder-free	32 – 33
MG 227	01120	Glass microfiber filter, with binder	30 – 31
MG 227/1/60	01124	Glass microfiber filter, with binder	30 – 31
MG 400	01123	Glass microfiber filter, with binder	30
MG 550-HA	01147	Glass microfiber filter, binder-free	32 – 33
MGA	1101	Glass microfiber filter, binder-free	32 – 33
MGB	1102	Glass microfiber filter, binder-free	32 – 33
MGC	1103	Glass microfiber filter, binder-free	32 – 33
MGD	1104	Glass microfiber filter, binder-free	32 – 33
MGF	1105	Glass microfiber filter, binder-free	32 – 33
MGG	1106	Glass microfiber filter, binder-free	32 – 33
Q3400	n.a.	Quartz microfiber filter	34
S 165	368	Filter paper for liquid absorption	17
SEK 770	419	Paper board	17

■ Did You know ...

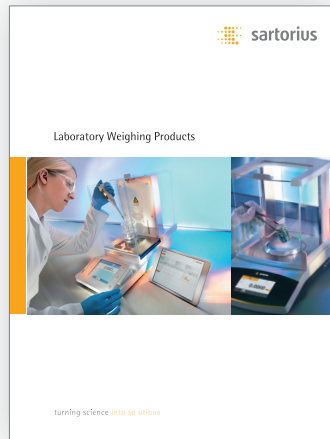
... that you can order more lab catalogs?

Get an overview of our high-quality laboratory instruments, high-grade consumables and excellent services.

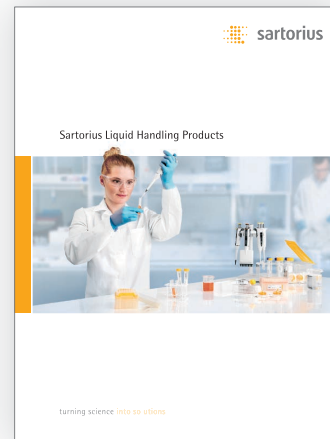
Our product portfolio focuses on laboratory instruments, such as lab balances, pipettes and laboratory water purification systems. Moreover, we offer the widest range of consumables like laboratory filters and pipette tips.



Products for Microbiological Control
Publication No.: SL-1530-e
Order No.: 85034-538-81



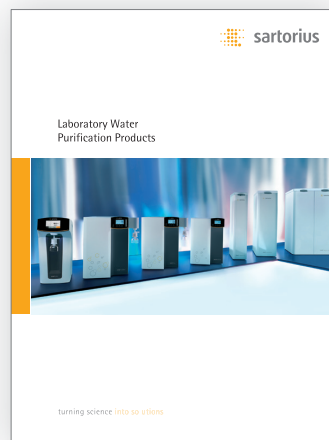
Laboratory Weighing Products
Publication No.: WL-0007-e
Order No.: 98649-018-13



Sartorius Liquid Handling Products
Publication No.: SUL0002-e
Order No.: 85032-542-25



Laboratory Filtration Products
Publication No.: SLU0006-e
Order No.: 85032-543-83



Laboratory Water Purification Products
Publication No.: SL-0002-e
Order No.: 85037-548-54



Download our lab catalogs from our website:
www.sartorius.com/lab-catalogs

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49.551.308.0

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan
Phone +33.1.70.62.50.00

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965760.0

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.90

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358.9.755.951

Hungary

Sartorius Hungária Kft.
Kagyló u. 5.
2092 Budakeszi
Phone +3623.457.227

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11
Phone +353.1.8089050

Italy

Sartorius Italy S.r.l.
Via Torino 3/5
20814 Varedo (MB)
Phone +39.0362.5557.11

Netherlands

Sartorius Netherlands B.V.
Phone +31.30.60.53.001
info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.6473830

Russian Federation

LLC "Sartorius RUS"
Vasilyevsky Island
5th line 70, Lit. A
199178 St. Petersburg
Phone +7.812.327.53.27

Spain & Portugal

Sartorius Spain, S.A.
Avda. de la Industria, 32
Edificio PAYMA
28108 Alcobendas (Madrid)
Phone Spain +34.913.586.095
Phone Portugal +351.800.855.800

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)
Phone +41.44.746.50.00

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ
Phone +44.1372.737159

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1.631.254.4249
Toll-free +1.800.635.2906

Argentina

Sartorius Argentina S.A.
Int. A. Ávalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900

Canada

Sartorius Canada Inc
1173 North Service Road West, D4
Oakville, ON L6M 2V9
Phone +1.905.569.7977
Toll-Free +1.800.668.4234

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepotzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepotzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Avenue Alberto del Campo 411
Floor 12 - The Office
15076 - San Isidro, Lima
Phone +51.1.441 0158

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800

China

Sartorius (Shanghai) Trading Co., Ltd.
3rd Floor, North Wing, Tower 1
No. 4560 Jinke Road
Zhangjiang Hi-Tech Park
Pudong District
Shanghai 201210, P.R. China
Phone +86.21.6878.2300

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong
Phone +852.2774.2678

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250

Japan

Sartorius Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan
Phone +81.3.3740.5408

Malaysia

Sartorius Malaysia Sdn. Bhd
Lot L3-E-3B, Enterprise 4
Technology Park Malaysia
Bukit Jalil
57000 Kuala Lumpur, Malaysia
Phone +60.3.8996.0622

Singapore

Sartorius Singapore Pte. Ltd
10 Science Park Rd
The Alpha #02-13/14
Singapore Science Park II
Singapore 117684
Phone +65.6872.3966

South Korea

Sartorius Korea Ltd.
8th Floor, Solid Space B/D,
PanGyoYeok-Ro 220, Bundang-Gu
SeongNam-Si, GyeongGi-Do, 463-400
Phone +82.31.622.5700

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310
Phone +66.2643.8361-6



◀ www.sartorius.com