

Extraction Thimbles

Efficient Solvent-Based
Separation from Solids,
Liquids, and Gases



Benefits

- Composed of high-purity cellulose and glass fiber for uncontaminated results
- Engineered for superior mechanical stability to withstand rigorous use
- Glass fiber thimbles offer remarkable resistance to high temperatures
- Designed for a precise fit across a diverse range of extractors
- Exhibits outstanding chemical resistance, ensuring reliability in various extraction processes.

Product Information

Our extraction thimbles have been designed to be compatible with the Soxhlet apparatus, making them ideal for a variety of applications. These include environmental monitoring tasks such as the separation of particulates like dust and aerosols, as well as the analysis of gas or air streams. They are also indispensable in food quality control processes, where they assist in the extraction of substances such as fats, emulsifiers, and additives. Designed for versatility, our thimbles can be utilized with any extractor, including models like the Tectator, to facilitate efficient and high-yield extraction procedures.

Cellulose and Glass Fiber Material

Both grades are free of binders and offer high purity. This ensures consistent and high overall flow rates. Our glass fiber thimbles are well suited to high temperature applications, whilst cellulose can be used for very sensitive applications. Our thimbles guarantee accuracy for any extraction system with common dimensions. Typical wall thicknesses for our C300 grade are 1.3 mm for inner diameters under 35 mm and 1.7 mm for inner diameters over 35 mm.

Applications

Primarily utilized in applied industrial processes, our thimbles are versatile consumables for various analytical procedures. They are particularly effective for measuring free lipid content in food and beverages, identifying additives in both food products and elastomers, and monitoring soil samples as well as hot air streams. Simply place the thimbles into your preferred extractor and employ the continuous extraction method for reliable and accurate results.

Cellulose Thimbles, Grade C300

Order Number	Inner diameter x Length (mm)	Wall Thickness (mm)	Weight (g/m ²)	Air Permeability at 2 mbar (L/m ² s)	Qty./Pack
FT-1201-019090	19 x 90	1.3	3	15	25
FT-1201-022080	22 x 80	1.3	2.5	18	25
FT-1201-022100	22 x 100	1.3	2.5	18	25
FT-1201-025060	25 x 60	1.3	2.5	18	25
FT-1201-025070	25 x 70	1.3	3	20	25
FT-1201-025080	25 x 80	1.3	3.5	20	25
FT-1201-025100	25 x 100	1.3	3.5	25	25
FT-1201-026060	26 x 60	1.3	2.5	15	25
FT-1201-028060	28 x 60	1.3	3	15	25
FT-1201-028080	28 x 80	1.3	3.7	20	25
FT-1201-028100	28 x 100	1.3	4.5	25	25
FT-1201-030080	30 x 80	1.3	3.8	21	25
FT-1201-030100	30 x 100	1.3	5	25	25
FT-1201-033060	33 x 60	1.3	3.2	15	25
FT-1201-033080	33 x 80	1.3	4.3	23	25
FT-1201-033090	33 x 90	1.3	4.6	30	25
FT-1201-033094	33 x 94	1.3	5	30	25
FT-1201-033100	33 x 100	1.3	5.5	32	25
FT-1201-033118	33 x 118	1.3	6.3	35	25
FT-1201-033130	33 x 130	1.3	7	37	25
FT-1201-033205	33 x 205	1.5	12	60	25
FT-1201-035150	35 x 150	1.3	9	43	25
FT-1201-040100	40 x 100	1.7	7.2	40	25
FT-1201-040123	40 x 123	1.7	10	45	25
FT-1201-040150	40 x 150	1.8	13	50	25
FT-1201-043123	43 x 123	1.8	13	50	25

Glass Microfiber Thimbles, Grade G400

Order Number	Inner diameter x Length (mm)	Wall Thickness (mm)	Weight (g/m ²)	Air Permeability at 2 mbar (L/m ² s)	Qty./Pack
FT-1204-019090	19 x 90	1.2	1.6	22	25
FT-1204-022080	22 x 80	1.6	2	21	25
FT-1204-025100	25 x 100	1.6	2.7	25	25
FT-1204-026060	26 x 60	1.5	1.8	18	25
FT-1204-030100	30 x 100	1.5	3	33	25
FT-1204-033094	33 x 94	1.5	3.5	29	25
FT-1204-043123	43 x 123	1.7	6.5	50	25



If the size you require is not listed, please don't hesitate to reach out to us. Our team is ready to assist you in customizing our products to meet your specific needs.

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

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

 **For further information, visit**
[sartorius.com](https://www.sartorius.com)