

# 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.

# Technical Specifications

Material	Diameter	Weight (g/m <sup>2</sup> )	Wall Thickness (mm)	Air Flow at 2 bar (L/m <sup>2</sup> / sec)
Cellulose	19×90	3	1.3	15
	22×80	2.5	1.3	18
	22×100	2.5	1.3	18
	25×60	2.5	1.3	18
	25×70	3	1.3	20
	25×80	3.5	1.3	20
	25×100	5	1.4	30.5
	26×60	2.5	1.3	15
	28×60	3	1.3	15
	28×80	3.7	1.3	20
	28×100	4.5	1.3	25
	30×80	3.8	1.3	21
	30×100	5	1.3	25
	33×60	3.2	1.3	15
	33×80	4.3	1.3	23
	33×90	4.6	1.3	30
	33×94	5	1.3	30
	33×100	5.5	1.3	32
	33×118	6.3	1.3	35
	33×130	7	1.3	37
33×205	12	1.5	60	
35×150	9	1.3	43	
40×100	7.2	1.7	40	
40×123	10	1.7	45	
43×123	13	1.8	50	
Glass fiber	19×90	1.6	1.2	22
	22×80	2	1.6	21
	25×100	2.7	1.6	25
	26×60	1.8	1.5	18
	30×100	3	1.5	33
	33×94	3.5	1.5	29
	43×123	6.5	1.7	50

# Ordering Information

Order No.	Description (Material, Grade, Inner Diameter×Length)
FT-1201-019090	Cellulose, C300, 19×90 mm
FT-1201-022080	Cellulose, C300, 22×80 mm
FT-1201-022100	Cellulose, C300, 22×100 mm
FT-1201-025060	Cellulose, C300, 25×60 mm
FT-1201-025070	Cellulose, C300, 25×70 mm
FT-1201-025080	Cellulose, C300, 25×80 mm
FT-1201-025100	Cellulose, C300, 25×100 mm
FT-1201-028060	Cellulose, C300, 28×60 mm
FT-1201-028080	Cellulose, C300, 28×80 mm
FT-1201-028100	Cellulose, C300, 28×100 mm
FT-1201-030080	Cellulose, C300, 30×80 mm
FT-1201-030100	Cellulose, C300, 30×100 mm
FT-1201-033060	Cellulose, C300, 33×60 mm
FT-1201-033080	Cellulose, C300, 33×80 mm
FT-1201-033090	Cellulose, C300, 33×90 mm
FT-1201-033094	Cellulose, C300, 33×94 mm
FT-1201-033100	Cellulose, C300, 33×100 mm
FT-1201-033118	Cellulose, C300, 33×118 mm
FT-1201-033130	Cellulose, C300, 33×130 mm
FT-1201-033205	Cellulose, C300, 33×205 mm
FT-1201-035150	Cellulose, C300, 35×150 mm
FT-1201-040100	Cellulose, C300, 40×100 mm
FT-1201-040123	Cellulose, C300, 40×123 mm
FT-1201-040150	Cellulose, C300, 40×150 mm
FT-1201-043123	Cellulose, C300, 43×123 mm
FT-1201-026060	Cellulose, C300, 26×60 mm

Order No.	Description (Material, Grade, Inner Diameter×Length)
FT-1204-019090	Glass Fiber, G400, 19×90 mm
FT-1204-022080	Glass Fiber, G400, 22×80 mm
FT-1204-025100	Glass Fiber, G400, 25×100 mm
FT-1204-026060	Glass Fiber, G400, 26×60 mm
FT-1204-030100	Glass Fiber, G400, 30×100 mm
FT-1204-033094	Glass Fiber, G400, 33×94 mm
FT-1204-043123	Glass Fiber, G400, 43×123 mm



Should the size you require not be listed, we invite you to contact us. Our team is eager to provide assistance and tailor our products to your exact specifications.

## Germany

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

## USA

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

 For more information, visit  
[www.sartorius.com](http://www.sartorius.com)

Specifications subject to change without notice.

© 2024 Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Status: 02 | 2024