

Gelatine Filter Disposables

For Quantification and Differentiation
of Airborne Organisms



Product Information

Gelatine filter disposables are individually packaged, presterilized and ready-to-connect units, each consisting of a gelatine membrane filter and a holder. Gelatine filters in conjunction with the MD8 Airscan Air Sampler (gelatine filter method) are used for collection of airborne microbes and viruses.

The gelatin filter method provides the following benefits for airborne organism sampling:

- Absolute retention of organisms since the filters capture all bacteria laden particles in the sample regardless of particle size.
- The filters prevent dessication of the retained organisms due to the filters' residual moisture content.
- Very low bacteria counts can be measured, because the high flow rates allow sampling of large volumes in short time, 1 m³ in < 8 minutes. This makes the method ideal for validating and monitoring the most critical sterile areas in pharmaceutical plants.
- Isokinetic sampling is possible in laminar-flow areas.
- The system is suitable for clean rooms because the disposables may be mounted inside the clean room with the sampling device mounted outside the room.
- The system can be calibrated by an accessory calibration unit.

In combination with the Sartorius MD8 Air samplers, gelatine filter disposables allow a series of air samples to be taken rapidly.

Gelatine filter disposables are easy to use

You only need to remove the disposable from the packaging, place it on the aluminum holder (17801), attach the holder to the MD8 and begin sampling.

Specifications

Technical Specifications of the Disposables	Filter holder: sized to accommodate 80 mm filters Filtration area: 38.5 cm ² Material: recyclable Cyrolite Thermal resistance: 85°C Dimensions: 93 × 16 mm
Technical Specifications of the Gelatine Filters	Material: gelatine Pore size: 3 µm Filter diameter: 80 mm Reaction to water: soluble Thickness: approx. 250 µm Flow rate per cm ² : 2.2 – 3.2 l/min at Δp = 0.05 bar (15 kPa) Thermal resistance: 60°C max.
Maximum Working Conditions for the Gelatine Filters	Temperature: 30°C Humidity: 85%
Sterilization	Gelatine filter disposables are supplied sterile by gamma irradiation.
Retentive Capacity of the Gelatine Filter for Bacteria	99.9995% at an inlet velocity of 0.25 m/s.
Retentive Capacity of the Gelatine Filter for Viruses	99.9% for T1 phages (coli phages) at a relative humidity of 50% and an inlet velocity of 0.3 m/s 99.94% for T3 phages (coli phages) at a relative humidity of 80% and an inlet velocity of 0.3 m/s.



Instruments

Ordering Information

Gelatine Filter Disposables

Order Number	Description	Pore Size	
		[μm]	\varnothing [mm]
17528--80----ACD	Disposable gelatine filter units, sterile, individually packed in 1 PE bag	3	80
17528--80----BZD	Disposable gelatine filter units, sterile, individually packed in 3 PE bags	3	80
17528--80----VPD	Disposable gelatine filter units, sterile, individually packed in 3 PE bags but label on innermost bag	3	80

Aseptic Transfer of Gelatine Membrane Filters (GMF)

Order Number	Description	Pore Size	
		[μm]	\varnothing [mm]
17528-----BFV	Gelatine membrane filter in Biosafe® Bags, sterile, individually packed in 1 PE inside the Biosafe® Bags	3	80



Gelatine Filters in Biosafe® Bags

Gelatine Disc Filter

Order Number	Description	Pore Size	
		[μm]	\varnothing [mm]
12602--37----ALK	Gelatine membrane filter	3	37
12602--47----ALK	Gelatine membrane filter	3	47
12602--47----ALN	Gelatine membrane filter	3	47
12602--50----ALK	Gelatine membrane filter	3	50
12602--50----ALN	Gelatine membrane filter	3	50
12602--80----ALK	Gelatine membrane filter	3	80

Reference Literature

1. Product Information brochure on the MD8 airscan Air Sampler (SLF3001-e).
2. Data Sheet: Calibration Unit (SL-2028-e).
3. Application Notes: Detecting Airborne Influenza Virus A (H3N2)...(SM-4018-e).
4. Application Notes: Collecting Airborne Viruses and Phages Using the Sartorius Gelatine Membrane Filter Method (SLF4028-e).
5. Four translated special reprints from "Bio Tec" on the subject of sampling virus aerosols using the gelatine membrane filter (SM-8030-e, SM-8031-e, SM-8032-e, SM-8034-a).
6. Test Report of the Public Health Laboratory Service, England 1993, Report No. 201/93.



Gelatine Filter Disposables




47 mm Gelatine filter with sampling head

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
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 further information, visit
www.sartorius.com