SVISCISVS

Our Unique Solutions for the Food and Beverage Industry

Characteristics and Advantages

Filtration of Beverages:

Extensive range of filtration equipment and consumables adapted to your needs

Microbiological Quality Control: Equipment and consumables designed to ensure optimum quality for your products

Expertise:

Recognized experience and knowledge in filtration and microbiological quality control

Compliant To:

Regulations 1935/2004 and 10/2011 of the European Union as well as FDA standards



Simplifying Progress

Our Laboratory Products Range

Laboratory Instruments

- Balances
- Pipettes
- Laboratory water systems
- Moisture analyzers
- Automated filterability index

Microbiological Control

- Filtration manifolds
- Membrane filters
- Single use and ready-to-use filtration units
- Nutrient pad sets and media broths

Consumables

- Syringe filters
- Microbiological test kits
- Pipette tips

Services

- Installation
- Maintenance | Repair
- Qualification (QI | QO)
- Calibration
- Training





Microsart® E-motion automatic membrane dispenser



Filtration manifolds



White, green, or grey membrane filters

Our Solutions for the Filtration of Beverages

Filtration Consumables

- Standard cartridges from 10" to 40"
- Jumbo Star cartridges
- Filtration sheets and modules

Filter Integrity Tester

Automated test device

Filtration Systems

- Manual filtration skids
- Automated industrial filtration skids
- CIP Systems (Clean In Place)
- Housings



the second

Cartridges



Jumbo Star cartridges

Applications

- Wines and sparkling wines
- Water
- Beer
- Spirits
- Fruit juice and soft drinks



Systems



Filtration sheets



Filtration modules



Automated test device



CIP systems



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 further contacts, visit www.sartorius.com