

How Safe is the Air Indoors?
Stay on Top of Airborne COVID 19
and Keep Safe

Simplifying Progress

SARTORIUS

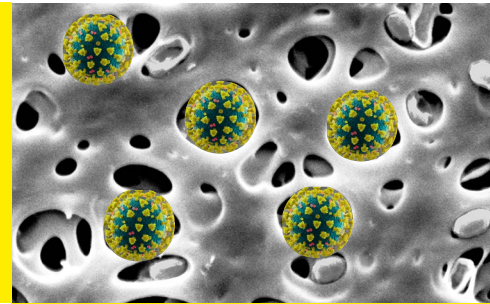
Worried about Virus Recovery From the Air?

We (Dis)solve Your Problem!

The soluble air filter for 99.94%* recovery of airborne viruses

The need to monitor the airborne viral load in critical areas is more important than ever and can be utilized as an early warning system of the pandemic spread.

The unique Sartorius Gelatine Airfilters combined with the patented membrane-dissolving technology enables you to detect every single virus that was retained on your membrane.



* for T3 coli phage at 80% rel. humidity by Gelatine Membrane Filters

Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

I. Air Sampling

For more information please click on the [+](#)



47 mm Gelatine Filter [+](#)



Portable Air Sampler:
MD8 Airport [+](#)



MD8 Airport Adapter [+](#)



Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

I. Air Sampling


For more informa

47 mm Gelatine Filter

Gelatine filters paired with the MD8 Air Samplers are used for collection of airborne microbes and viruses. Gelatine filters are pre-sterilized by gamma irradiation and ensure near absolute retention of microbes and viruses (99.9995% for *B. subtilis niger*; 99.94% for T3 coli phages).

Since the filters are made of 100% gelatine, they are fully soluble which provides easy access to the retained microbes and viruses for subsequent analysis.



47 mm Gelatine Filter 

Portable Air Sampler: 
MD8 Airport

MD8 Airport Adapter 

Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

I. Air Sampling

For more information

Portable Air Sampler: MD8 AirPort
The MD8 air sampler is designed to sample the smallest viruses and microorganism in the air by using the unique Gelatine Membrane Filters (GMF). With its digital interface, all needed parameters, such as flow rate and sampling volume, can easily be adjusted.



47 mm Gelatine Filter +

Portable Air Sampler: MD8 AirPort +

MD8 Airport Adapter +

Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

I. Air Sampling

For more informa

MD8 Airport Adapter

The geometry of this stainless steel adaptor is specifically designed to support the 47 mm Gelatin Membranes and optimized to deliver appropriate flow rates.

After sampling, the membrane is easily transferred touch-free to the dissolving chamber (Microsart®@solve) via an inner lining of glue.



47 mm Gelatine Filter +

Portable Air Sampler: MD8 Airport +

MD8 Airport Adapter +

Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

II. Dissolving the Membrane

For more information please click on the [+](#)



Microsart®@solve [+](#)



Dissolve the Membrane [+](#)



Centrifuge Adapter [+](#)

Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

II. Dissolving t For more informa

Microsart®@solve

Within this closed chamber your gelatine membrane can securely be transported from the sampling location to the lab and is ready to be dissolved afterwards.



Microsart®@solve



Dissolve the Membrane



Centrifuge Adapter



Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

II. Dissolving the Membrane

For more information

Dissolve the Membrane

Add 1.7 ml solvent into the chamber and incubate for 10 min at 37°C to fully dissolve your Gelatine Membrane filter. If the goal is to perform virus infectivity studies through transfection, simply dissolve the membrane in the appropriate culture medium.



Microsart®@solve



Dissolve the Membrane



Centrifuge Adapter



Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

II. Dissolving the Membrane

For more information, see the Sartorius website.

Centrifuge Adapter

After incubation, centrifuge your dissolved membrane into a standard 2 ml reaction tube. Sartorius provides you with the appropriate centrifuge adapter for your application.



Microsart®@solve



Dissolve the Membrane



Centrifuge Adapter

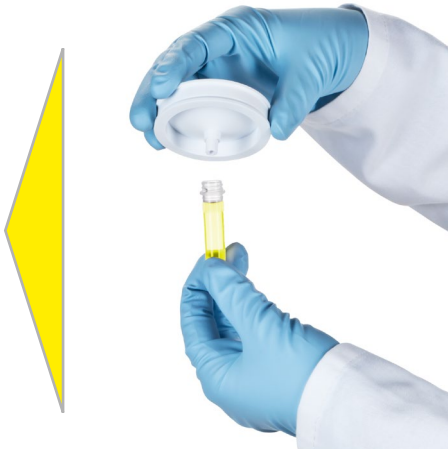


Workflow -

Collection of Airborne Viruses Using Gelatine Membrane Filtration

With the help of the portable air sampler MD8 AirPort, the air of all high-contamination risk areas can be sampled for coronavirus, as well as any other virus. After sampling, the membrane filter can simply be dissolved in minimal volumes of water or buffer. This aids RNA sample preparation for rapid testing, using methods such as PCR.

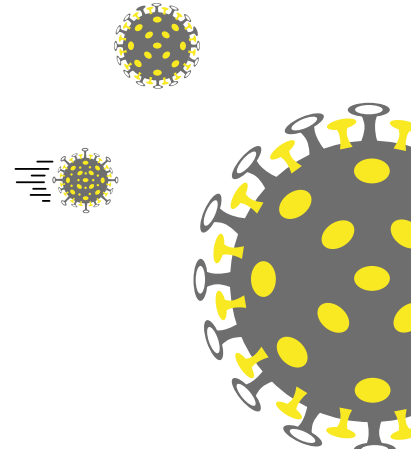
III. Virus Detection



Your Entire Sample in
Just One Tube



Ready-to-use for
Subsequent Analysis.



Virus Detection via PCR

Learn More

Find application data for virus detection, as well as relevant scientific publications for airborne COVID 19 detection using Sartorius Gelatine Membrane Filtration on our website.

Contact us for a quote, call, or demo:

Visit Our Website:



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

Germany

Sartorius GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius North America Inc.
5 Orville Drive, Suite 200
Bohemia, NY 11716
Toll-Free +1 800 368 7178