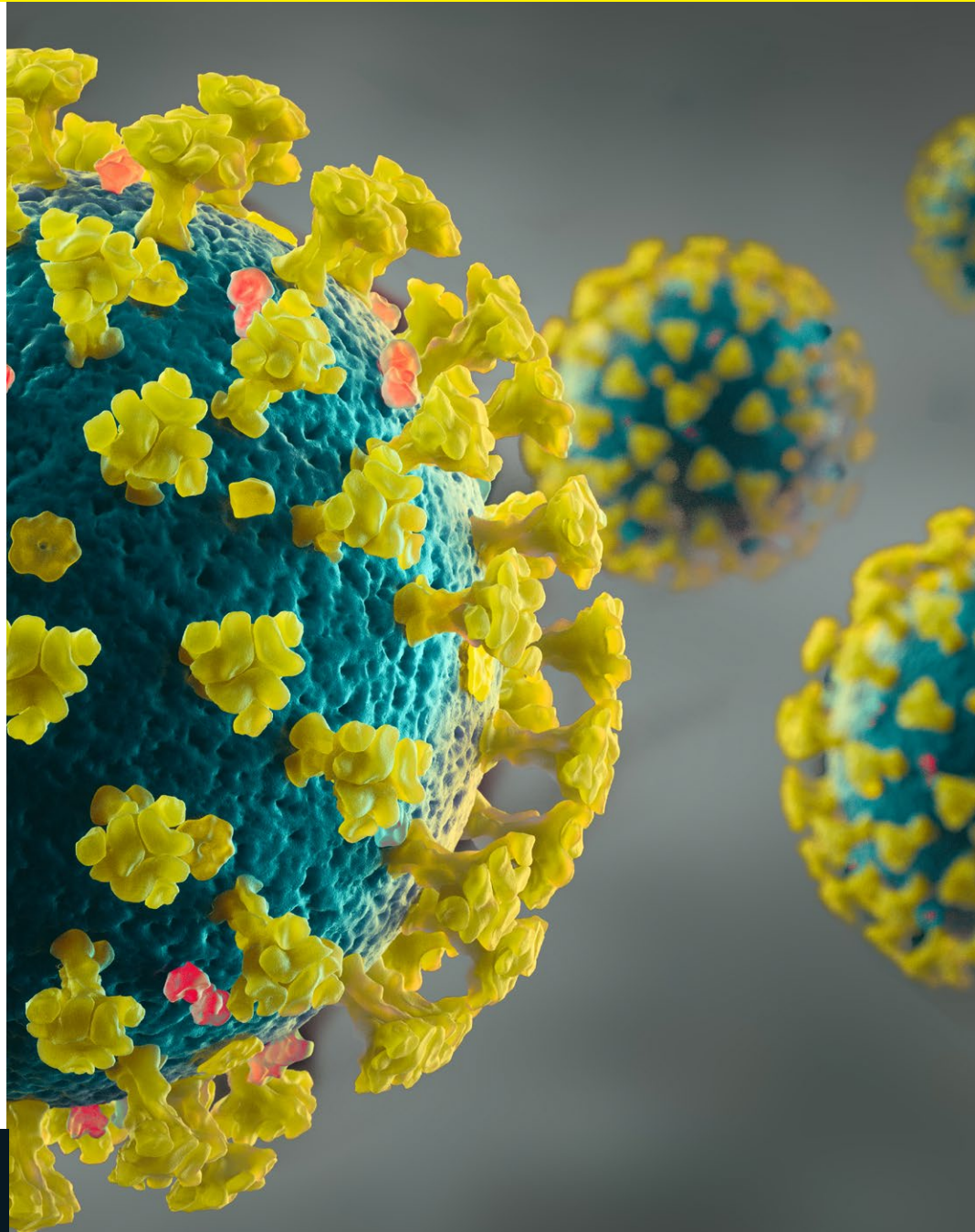


## Sartorius Solutions for the Detection of SARS-CoV-2 Coronavirus in Environmental Samples, Air and Water

To address the need for more insight into COVID-19 in line with CDC and WHO guidelines, we support your efforts with solutions for environmental testing. If you would like to collect viruses from air samples, seek for solutions to concentrate viruses from water samples, or optimize the way you handle your samples, then we offer many solutions to facilitate your work.



### Collection of viruses in Air Samples

Water soluble gelatin membrane filters are the perfect way to quickly capture SARS-CoV-2 in air samples.



Gelatin Membrane sampling heads

With the help of the portable AirPort MD8 air sampler, the air of all high-contamination risk areas can be sampled for coronavirus. After sampling, the membrane filter can be dissolved in minimal volumes of water or buffer, aiding with RNA sample preparation prior to rapid tests methods such as PCR.



MD8 AirPort air sampler

### Concentration of viruses in Water Samples

Ultrafiltration products are ideal to isolate and concentrate viruses or virus like particles from drinking, waste and surface water samples prior to the final detection methods, such as RT-PCR. Sartorius offers ultrafiltration devices for sample volumes from microliters to several liters.

The product range includes several membrane materials and molecular weight cutoffs (MWCO), to cover a wide range of applications.



Vivaspin® Centrifugal Filtration Device from 100 µl to 20 mL



Vivaspin® Centrifugal Filtration Device from 20 mL to 100 mL



Vivaflow® Tangential Flow Filtration devices from 50 mL to several liters

### Pipetting and Sample Transfer

In detecting SARS-CoV-2, RNA preparation and RT-PCR are pipetting intensive protocols, for which precise, accurate and contamination free pipetting is required.



Ergonomically designed Sartorius mechanical Tacta®, electronic pipettes Picus® and Picus® Nxt are Easy-to-clean and resistant making regular decontamination simple during coronavirus testing.



RNase|DNase free pre-sterilized SafetySpace™ Filter tips help protect samples from contamination

### 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](http://www.sartorius.com)