

Gain Deeper Insight.
Streamline Workflows.
Accelerate Your
Research.



Environmental Testing for Virus Research & Monitoring

Discover Air Sampling Tools for Detection of Viruses in Your Air Indoors.



Airport MD8 Air Sampler



Gelatine Membrane Filters

Efficient Preparation of Wastewater Samples Prior to Virus Detection

Find the optimal ultrafiltration device for virus concentration.



Vivaspin® Centrifugal Concentrators



Vivaspin® Pressure Driven Concentrators



Vivaflow® Crossflow Cassettes

Innovative strategies and technologies are needed to combat global pandemics like COVID-19. Researchers need deeper understanding of the cellular and molecular mechanisms that drive immune responses against infectious diseases. Sartorius provides groundbreaking solutions that simplify complexity and expedite discovery for scientists working on the frontline of viral vaccine research.

- 01 Incucyte® SX5 Live-Cell Analysis System
- 02 iQue®3 Advanced Flow Cytometry Platform
- 03 Octet® Bio-Layer Interferometry (BLI) Platforms



Find out more
For more information, please visit
www.sartorius.com



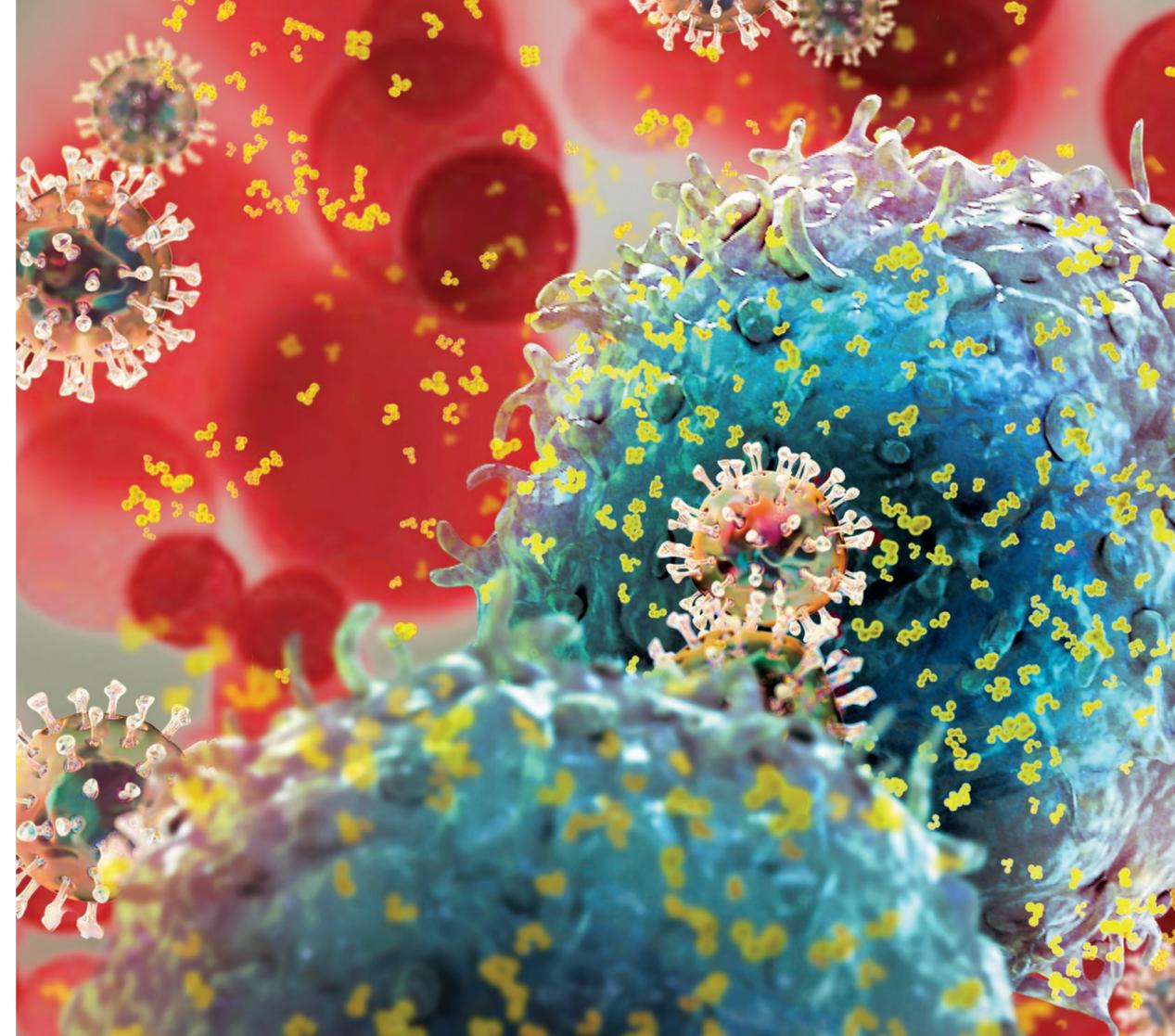
Sales and Service Contacts

For further contacts, visit
www.sartorius.com

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

Specifications subject to change without notice.
Copyright Sartorius Lab Instruments GmbH & Co. KG. All Rights Reserved.
Printed in the EU or US on paper bleached without chlorine.
Version 2021 | 09



Solutions for Viral Vaccine Research

Simplifying Progress

SARTORIUS

Solutions for Viral Vaccine Research

Incucyte® Live-Cell Analysis System



In the race for a vaccine, accelerating biological insight is critical. The Incucyte® Live-Cell Analysis System improves productivity and insight with automated image acquisition and integrated analysis. Accelerate your next discovery with a flexible assay platform that accommodates multiple applications simultaneously with multi-plate throughput.

Viral Replication and Infection Assays

Viral replication and infection assays enable the study of pathogenesis in host cultures and the testing of novel antiviral therapeutics.

Live-Cell Immunocytochemistry

Complement your existing immunocytochemistry protocols and start gaining new insights into the dynamics of surface protein expression.

Cell Health and Viability Assays

Study mechanism of infection through assessing virus modulated cell health, proliferation and viability in real time.

iQue® Advanced Flow Cytometry Platform



The iQue® Advanced Flow Cytometry Platform gets you from samples to actionable results in record time. The iQue platform can be used for viral vaccine research and virology activities such as antibody library screening, immune response assessment, including epitope mapping and antibody neutralization studies.

Antibody Library Screening

Screen antibody libraries against cell surface proteins that play a key role in virus inhibition.

Neutralizing Antibodies

Measure antibody affinity to monitor viral neutralization while concurrently evaluating host cell proliferation.

T Cell Activation and Exhaustion

Identify the extent of T cell activation and markers of T cell exhaustion to provide insight into the immune response following viral infection.



NutriVero™ Flex 10

Next-generation chemically defined serum-free, animal component-free medium.

Virus Counter® for Rapid Virus Quantification



The Virus Counter platform enables rapid, direct and precise quantification of total virus particles. These real-time insights can be used to track virus particles during up- and downstream processes, accelerating development timelines and optimizing product yields.



Microsart® AMP Mycoplasma | Bacteria | Fungi Kit

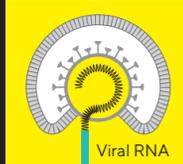
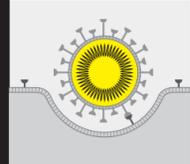
Ensure your cells are free from any microbial contamination with our real-time PCR kits.

Vivaflow® & Vivaspin® Ultrafilters and Vivapure® Membrane Adsorbers

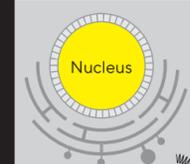


Concentration, purification and diafiltration of proteins, antigens, antibodies and viruses.

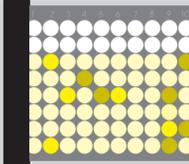
Investigate Host-Virus Interactions and Immunogen Characterization



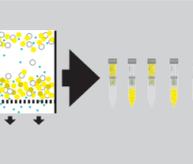
Identify Potential Antigens Profile Immune Responses



Express and Analyze Antigens | Antibodies in Cell Cultures



Harvesting and Concentration



Octet® Bio-Layer Interferometry (BLI) Platforms



Octet® BLI platforms enable fast real-time label-free determination of affinity, kinetics, and antibody | protein concentration analysis. Octet® platforms can be used, for example, to characterize virus-host protein interactions, antibody binding characterization, antibody epitope binning, vaccine titer and potency assays.

- Fluidic-free Dip-and-Read™ measurements help to get results within minutes.
- Crude sample analysis compatibility enables reduced sample preparation times.
- Easily construct your binding and quantitation assays using a range of available biosensor chemistries.

Select Vaccine and Therapeutic Candidates Faster

Measure affinity and binding kinetics of neutralizing antibodies, peptide and small-molecule inhibitors binding to viral proteins to advance best therapeutics candidates.

Assess Epitope Diversity and Coverage

Rapidly categorize and bin antibodies that bind to similar epitope regions using Octet® cross-competition assays.

Measure Antibody and Virus Titers

Octet® assays enable fast at-line testing of antibody and virus titers (e.g., AAV, Influenza, etc.) direct in process development samples.



Ambr® 15 and Ambr® 250

High-throughput cell culture systems with scalable, multiparallel mini-bioreactors and automated functionality accelerate cell line | strain selection and process development.

Sartoclear Dynamics® Lab V kits

Virus harvesting in one step.

