

Simplify Virus Particle Quantification with the Virus Counter[®] Plus Platform

The Virus Counter[®] Plus platform is purpose built to rapidly and precisely quantify total virus particles in process development and manufacturing. The Virus Counter[®] Plus platform featuring the new Virus Counter[®] Plus high-throughput instrument, highly specific reagents and integrated software, make virus particle quantification rapid, precise and simple.

- Accelerate development times with real-time virus particle titer insights for applications in vaccines and viral therapies
- Improve viral vector product yield by optimizing development and manufacturing processes
- Increase product quality by in-depth sample analysis by quantification of total intact particles



Rapid, Direct, Biologically Relevant Virus Quantitation



Rapid Results

Get virus titer insights rapidly to accelerate your time to result and process development timelines



Simple Workflows

Virotag® assay kits offer ready-to-use, specific and flexible quantification solutions with simple, no-wash assays. Labeling reagents allow the direct enumeration of virus particles through fluorescence



High Precision

High instrument reproducibility for confidence in your data. Performance verification solutions ensure optimal instrument performance



Integrated Data Analysis

Integrated software directs instrument function, automated analysis in a 96-well format, and data analysis. New advanced signal detection improves sensitivity and precision



Robust Design

Easy to clean surfaces and redesign of critical components including the optical deck support high reliability and availability

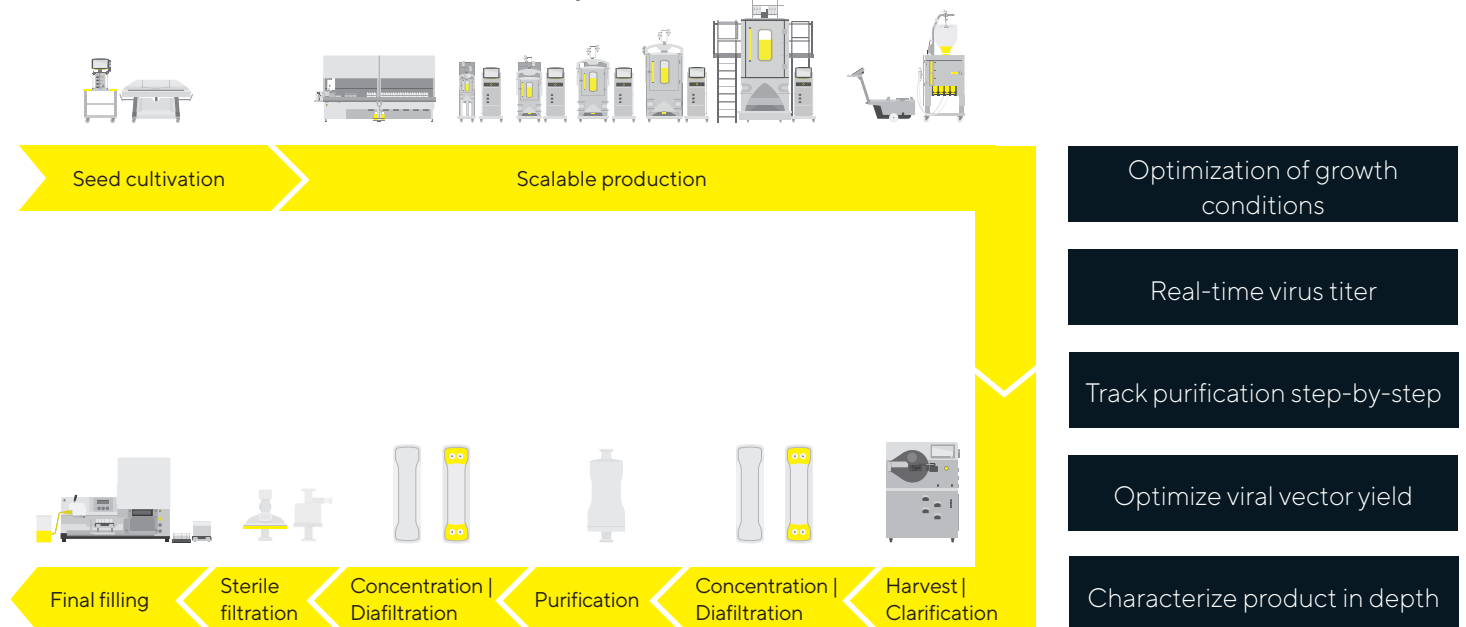


Continuous Development

The platform addresses existing and evolving customer needs by continuously expanding the Virotag® reagent portfolio and adding new features to the system

Results You Can Count On Throughout Your Process

Precise results in minutes instead of hours or days



Simple, No Wash Sample Staining Allows Rapid Quantification



Pipette Virotag® reagent to viral vector samples



Incubate 30 min - 1 h in the dark



Multiple sample (96-well plate) processing in Virus Counter® Plus instrument with walk-away automation



Analyze viral vector titer in simple steps directly in the Virus Counter® Plus software

Virus Counter[®] Reagent Kits

Virotag[®] DY Reagents

Universal Dual Stain Labeling of Viruses

- Broad dye-based quantification approach
- Identifies and detects a wide range of viruses using simultaneous staining of proteins and nucleic acids
- Optimized for enveloped viruses
- Compatible with single- and double-stranded DNA and RNA viruses

Virotag[®] AB Reagents

Antibody Based Labeling of Viruses

- Highly specific detection reagent based on antibody recognition of epitopes on viral vectors
- Detection of total viral particles independent of nucleic acid content
- Quantification of virus-like particles (VLP)
- Enumeration of both enveloped and non-enveloped viruses

Virotag [®] DY Reagents	Compatible Virus viral vector
Virotag [®] DY ENV	All enveloped viruses (e.g., Lentivirus, Influenza, Retrovirus, Baculovirus)

Virotag [®] AB Reagents	Compatible Virus viral vector
Virotag [®] BCVB	Baculovirus
Virotag [®] INVA	Influenza virus A
Virotag [®] INVB	Influenza virus B
Virotag [®] VSVG	Lentivirus (VSV-G pseudotyped)
Virotag [®] XXX	Different Virus types in development

Additional Virotag[®] reagents are currently in development and will be released soon.

Customer Testimonials

“The Virus Counter[®] has proven to be a vital tool in obtaining accurate, quantifiable, and timely results in our efforts of continual process improvement and intensification. Viral titers using the traditional plaque assay typically would take a few days, and many labor hours, but with the VC assay, our sample-to-result time has dramatically decreased to less than one hour.” Novavax

“Virus Counter[®] results significantly correlated with both plaque assay and qRT-PCR. These results demonstrated that the VC is an easy, fast, and consistent method to quantify filoviruses in stock preparations.”
USAMRIID

For Further Information Contact Your Sartorius Sales Representative


The Virus Counter[®] Plus platform is for research use or further manufacturing use only – not for use in therapeutic or diagnostic procedures. They are not for in vitro diagnostic use nor are they medical devices. Drug manufacturers and clinicians are responsible for obtaining the appropriate IND/BLA/NDA approvals for clinical applications. Class 1 Laser Product.

Germany

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

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

 For further contacts, visit
www.sartorius.com