

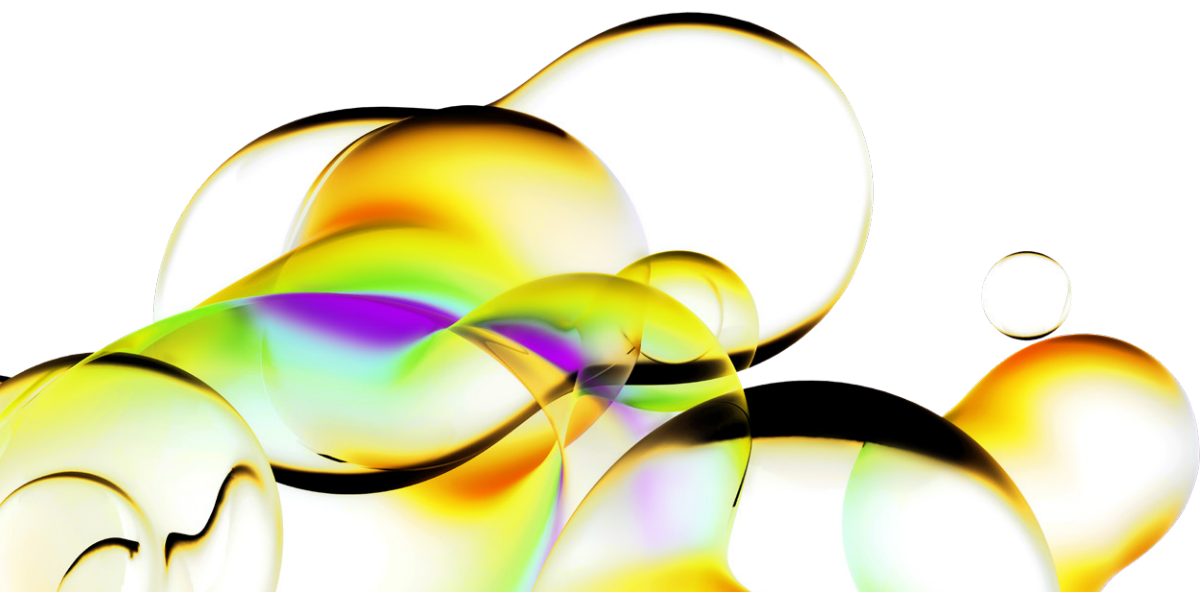
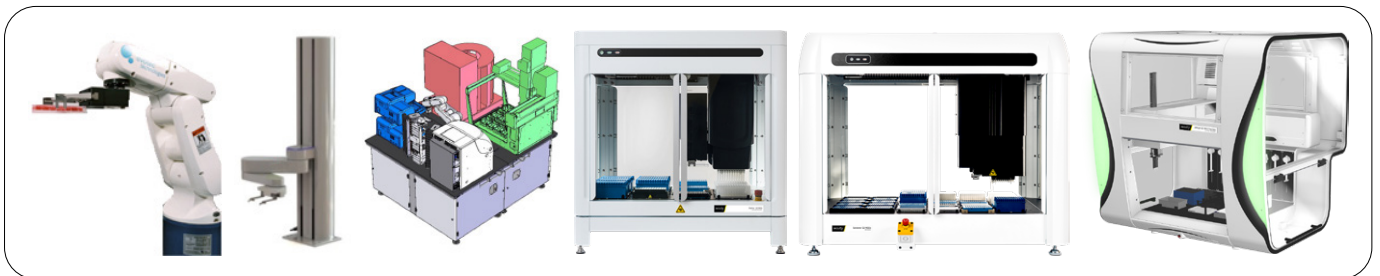
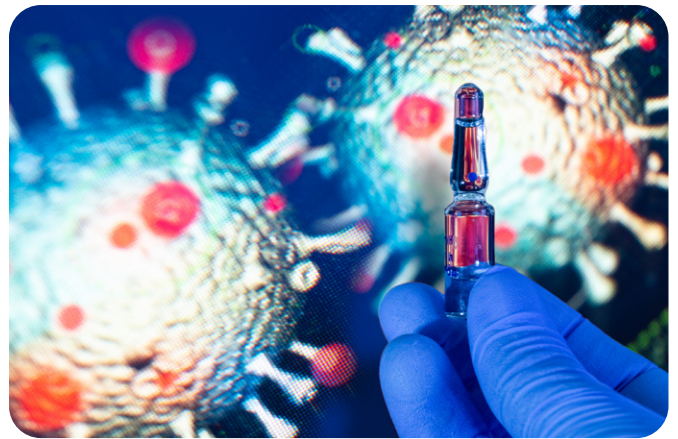


Coming together on a mission.

Revvity offers high-throughput solutions that support many applications in virology and enable researchers to accelerate workflows to meet the needs of rapid vaccine and therapeutics development.

Automation

Whether you are researching virus genetic diversity and tracking mutations or trying to understand differences in human disease severity, our genomics solutions are optimized and automated, supporting you from DNA/RNA extraction up to library quality control.



Virus titer and virus neutralization



EnSight® – Automated virus plaque imaging and analysis using the EnSight multimode plate reader provides a rapid and reliable way for researchers to determine viral titers and perform plaque reduction neutralization assays.

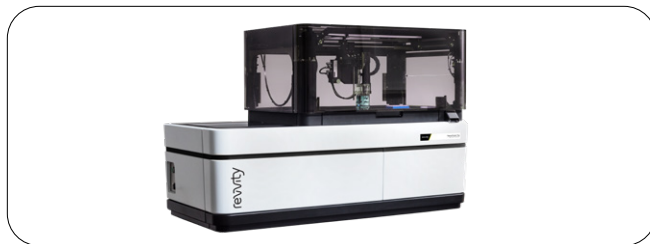


Lites™ – Based on luminescence, Lites Reporter gene assays allow sensitive detection of viral infection or virus neutralization. ATPLite luminescence assays provide simple ways to measure virus induced cell toxicity.

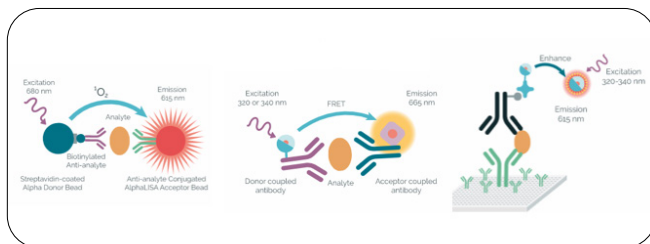
Host-virus interactions



VICTOR® Nivo™, EnVision® – Multimode detection instruments, together with reagents and microplates provide robust and convenient tools to study protein-protein interactions or identify inhibitors of PPIs.



Opera Phenix® Plus – High-content based siRNA screening is a drug target identification method, that when applied to virus research can uncover host cell factors required for virus infection.

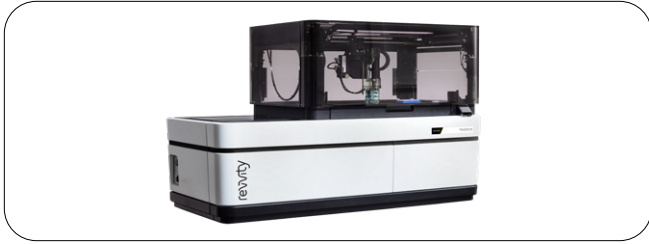


Alpha, DELFIA, Revvity's HTRF® – Revvity's range of reagents allow measuring viral presence, viral attachment, genome integration and virus assembly as well as cytokine responses, inflammatory pathways and mediators, which are critical indicators that need evaluation during vaccine development.

Drug/vaccine discovery and development



EnVision®, and VICTOR® Nivo™ plate readers (with reagent kits) enable detection and characterization of viral proteins and host immune markers.



Opera Phenix Plus and Operetta® CLS high-content screening and analysis platforms allow understanding the pathogenesis of viral infection and the host immune response, and support compound screens for drug discovery or repurposing as well as toxicity studies for drug development.



LabChip® GXII Touch™ – Microfluidic capillary electrophoresis on the LabChip GXII Touch, combined with consumables and reagents provides rapid, reliable and quantitative protein and nucleic acid analysis for a variety of virus related applications: QC of mRNA/DNA/protein vaccines, characterization of antibody therapeutics in between analysis of virus-like particles.



IVIS® – Small animal *in vivo* imaging on the IVIS system allows studying viral infection, biodistribution and dosing *in vivo*. Animals can be studied repeatedly over time, thus limiting the number of animals needed and allowing to understand therapy or vaccination effectively over time

***In vivo* imaging reagents** – *In vivo* imaging reagents enable visualization of important biomarkers *in vivo*, such as those involved in lung inflammation and others. Deep red fluorescent tags can label a variety of molecules, such as drugs or vaccines to study biodistribution *in vivo*.

Analytics Supporting Your Clinical Development Life Cycle

Revvity Clinical Solutions, powered by TIBCO® Spotfire® – provides a single unified platform to support your data analytics – from source to visualization to action. With unrivaled workflow flexibility to support dynamic collaboration, Revvity's Clinical solutions are accelerating the delivery of urgently needed therapeutics to patients.

Reach out to your local sales representative or visit www.revvity.com for more information.

For research use only. Not for use in diagnostic procedures.

The Revvity logo, consisting of the word "revvity" in a lowercase, sans-serif font.