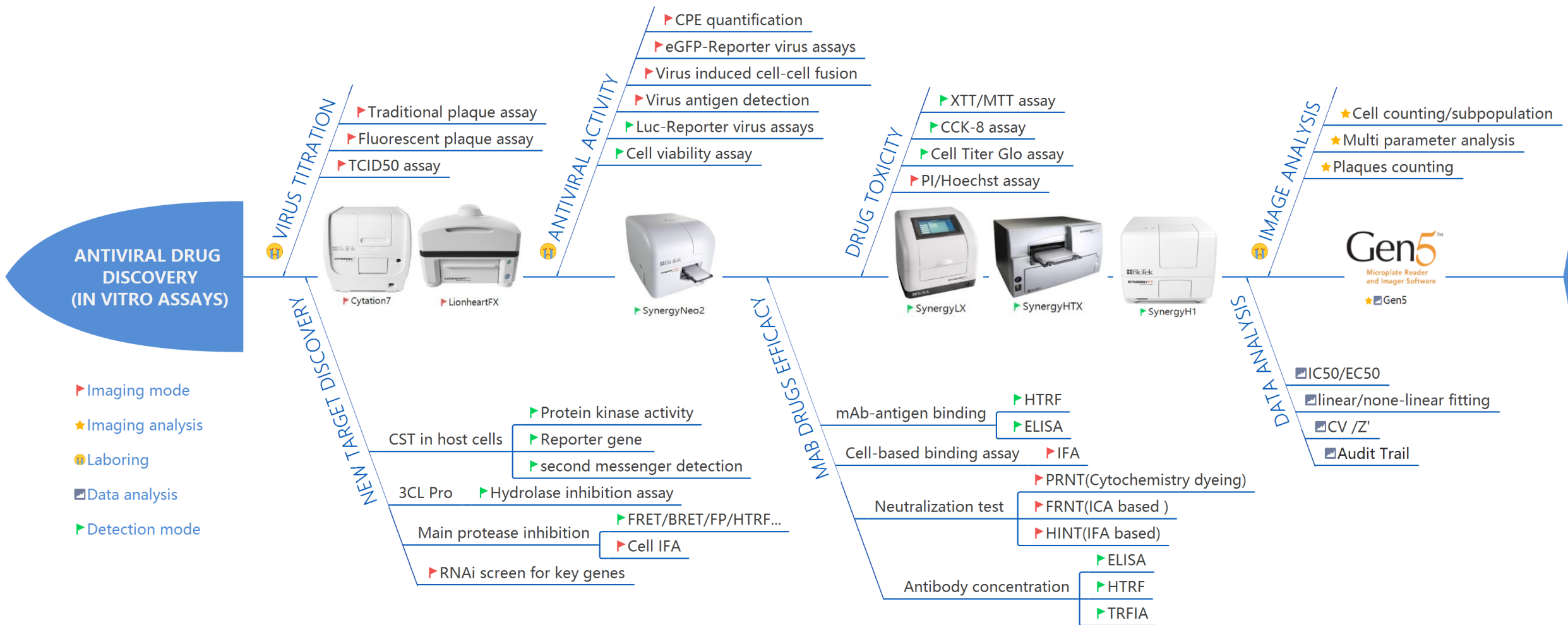
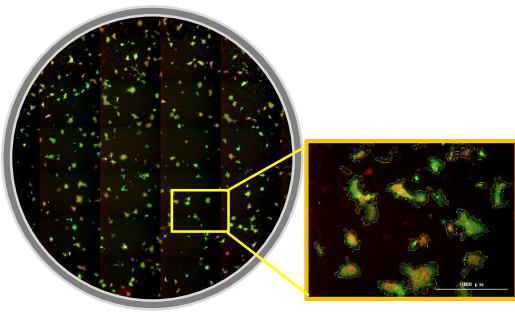


In Vitro Assays for Antiviral Drug Development: Solutions from BioTek

During an infectious virus outbreak, global medical workers are anxious to find effective therapeutic drugs to suppress the virus. Antiviral drugs can include vaccines, small molecular compounds, peptide inhibitors, nucleoside analogues, natural products and monoclonal antibodies. No matter what the drug development strategy is, in vitro assays in microplates are essential. The following shows the main types of microplate-based assays and detection methods.



Cytation™ 7 cell imaging multi-mode reader with WFOV can easily capture the viral plaques in a whole well. Gen5 software is used to count the plaques.



High through-put imaging with 6- to 1536-well plates.

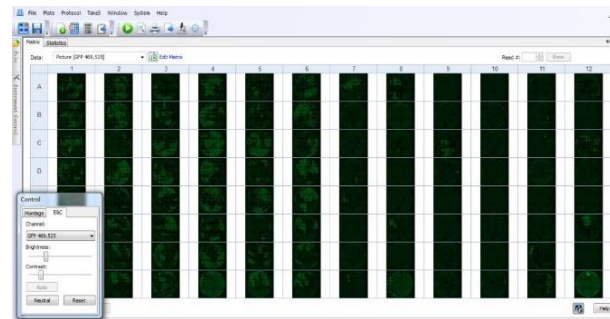
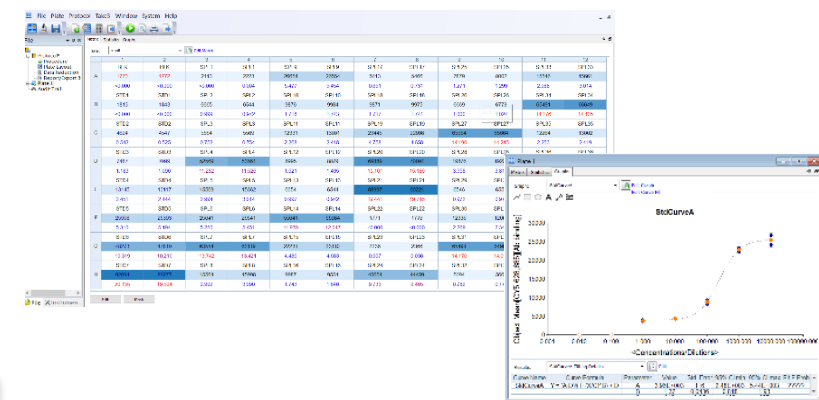


Image & data analysis with Gen5 software. Linear and non-linear curve fitting, EC₅₀/IC₅₀ calculation, statistics and data management.



Liquid Handling Solutions

There are many repetitive procedures in the anti-coronavirus drug in vitro assays, including cell plating, fluid handling, high-throughput reading/imaging requirements, etc. BioTek's washers, dispensers and automation products are versatile and can be combined into different automatic solutions to help researchers by reducing the repetitive manual work, while ensuring stable and reliable data.

Automation Solutions

50™ TS



50 TS is a compact microplate washing system with easy touch operation for washing full or partial plates.

405™ TS



405 TS is a high throughput washer with many patented technologies.

MultiFlo™ FX



MultiFlo FX automates rapid microplate dispensing and washing, as well as gentle media exchange and dispensing for suspension cells and 3D cultures.

BioStack™



BioStack is a compact and versatile microplate stacker compatible with BioTek's washers, dispensers, readers and imaging systems.

BioSpa™ Automation



BioSpa Live Cell Analysis System automates many applications in multiple plates for continuous live cell imaging and analysis. The system includes BioSpa 8 Automated Incubator, Cytation imaging readers and BioTek liquid handling instruments. Together, they automate kinetic live cell workflows in multiple plates for hours, days or weeks.