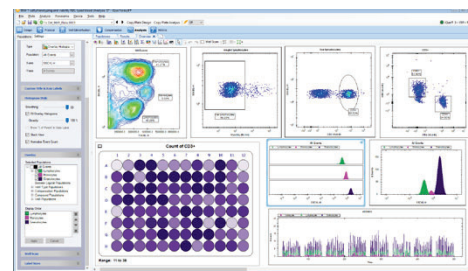


Accelerate your insights with an all-inclusive software package

iQue Forecyt® software provides automated real time analysis with plate level analytics. Unique analysis tools quickly transform even the most complex experimental data sets into actionable results and eliminate the need to perform time consuming analysis of data files on a well-by-well basis.



iQue® Software Editions

iQue Forecyt® Standard Edition – Instrument Control and Analysis



Connected directly to iQue®

Included with every instrument. Used to set up experiments, acquire data, and analyze results on the computer connected to the iQue®.

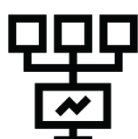
iQue Forecyt® Offline Edition – Data Analysis Without the Lab or Network



Instrument-free analysis powerhouse

Take insights on the go with Offline Edition. Ideal for users who want to analyze data away from the laboratory. No network connection required – import experiments and templates to a locally installed iQue Forecyt® database. This version has no acquisition capabilities, and is for experiment set-up and analysis purposes only.

iQue Forecyt® Enterprise Edition – Multi-User Access and Data Sharing



Sync iQue® data safely and seamlessly across a network

A scalable solution for multi-user or multi-instrument environments. Multiple users connect directly to an iQue Forecyt® database over the network using floating licenses, eliminating the need to import | export data and allowing multiple users to set up, run and analyze experiments simultaneously from any PC connected to the network. Additional floating licenses can be added as your demand increases.

Optional Licensable Features

iQue® Automation API License – Robotic Integration



Automation

Enable continuous plate loading with the Robotic Integration API to relieve capacity bottlenecks throughout your entire experiment with the fastest plate sampling, integrated analysis, and novel data reduction tools.

iQue® 21 CFR Part 11 Software Module – Traceability and Accountability



Documentation

Audit trail, electronic records and electronic signature features to support 21 CFR Part 11 workflows. Every step from acquisition through to exported results will be tracked within a single Audit Log. Custom standardized acceptance criteria and workflows to create the best fit for your organizational day to day needs.

Acquire

Intuitive and user-friendly acquisition panels facilitate quick-start experiment set-up. Define wells to be sampled and additional annotation layers such as dilution series, cell types and other parameters that can be used in downstream analysis. View data as it is being acquired.

Analyze

Use iQue Forecyt®’s gating tools to identify cell populations, including “smart gates” which automatically adjust to different data sets. Changing gate position results in real-time updates, allowing gating strategy effects to be visualized immediately across the entire plate.

Quickly generate Dose Response graphs to fit concentration curves to well metrics, across an entire plate.

Visualize

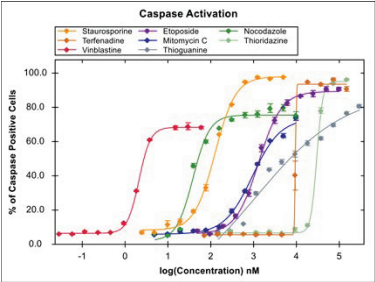
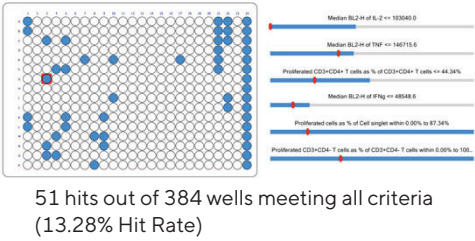
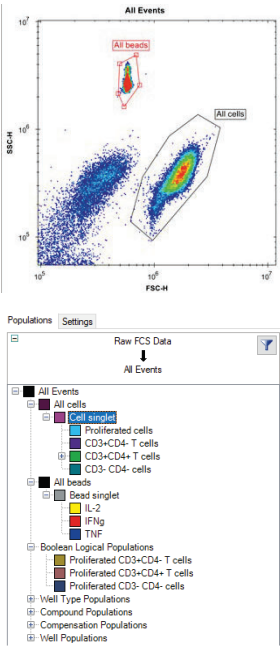
Profile maps provide dynamic adjustment of activity ranges within the selected metrics, enabling real-time hypothesis testing or optimization of activity ranges to align unknown samples with positive and negative controls.

With Panorama, view multiple plates simultaneously to compare heat maps, identify hits across profile maps, and rank wells by any metric with user definable line graphs.

Interpret

Generate the data you need with custom metrics derived from basic metric building blocks using a simple wizard. All metrics become available for additional analysis, visualization, and/or exporting. Go from sample acquisition to data interpretation with publication ready images.

Ordering Information



Description	Order No.
Forecyt® Enterprise Edition	90212
Forecyt® Enterprise Lite	60036
Forecyt® Enterprise Edition Floating Client License (1)	60028
Forecyt® Enterprise Edition Floating Client License (5)	60029
Forecyt® Offline Edition for single PC	90272
Robotic Integration API License (1)	90208
iQue® 21 CFR Part 11 Software Module	60035
iQue® Signal Reduction Software Module	BA-60037



To Place an Order or Request Additional Information Info:
info.intellicyt@sartorius.com