

Flexible Automation Enablement

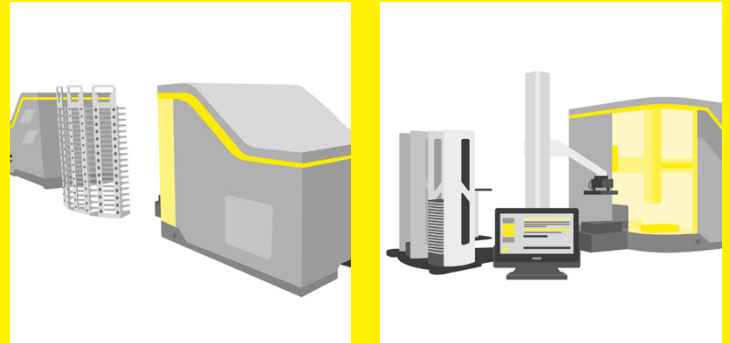
On the iQue® High-Throughput Screening (HTS) Cytometry Platform

Maximize Plate-Processing Power

Utilize microplate loading robotic arms to allow unattended analysis of samples (96-well or 384-well) on the iQue® Platform.

- Reduce manual intervention
- Automated cleaning and shut-down
- On average, an iQue® user runs 45 plates | experiment with a variety of robotic integrations

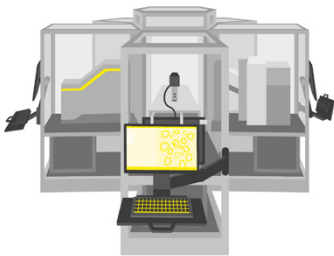
The iQue® Platform is an automation agnostic system by design and pairs well with a variety of established liquid-handling robotics like Hudson Robotics PlateCrate EX™ Microplate Handler, HighRes® Biosolutions Acell, Thermo Scientific™ Spinnaker™ Microplate robot.



Integrate an iQue® Into an Automated Workflow

Create multifaceted workflows using the iQue® Platform alongside multiple systems and detectors to amplify data analysis capabilities.

- Reduce assay variability
- Improve unattended run-times
- Remove manual intervention

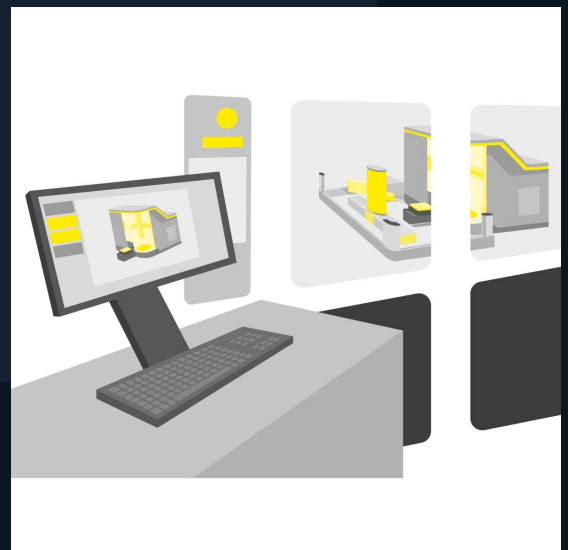


For example, use the iQue® Platform with an autosampler to amplify throughput, automated incubator to control experimental conditions, plate readers for powerful secondary analysis, a dispenser system to automate sample administration, a transfer system to remove manual transfers between systems, a hotel rack microplate system for additional throughput, and many more.

Implement Environmental Control

iQue® systems were built from the beginning to support automation compatibility and integrate the iQue® Platform into a workstation to enable tighter control of broader experiments.

- Control temperature, humidity and more
- Centralize or remote monitoring
- Safety for samples and personnel



Speed.

Unparalleled speed allows library screening in days, not months

Scalability.

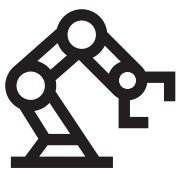
Automation enablement to greater throughput and less manual intervention

Simplicity.

One centralized control of a multi-user environment



Connecting iQue® to Automation



3rd Party Automation Controller



iQue Forecyt® Controller

1. Connect iQue® using the **Robotic Integration Automation Programming Interface (API) License** via a Serial | TCP connection

- License includes all the tools necessary including Screen Development Kit (SDK)
- Compatible with 21 CFR Part 11 compliance

2. Remotely control the iQue® functionality

- Continue to utilize iQue® stand-alone if needed



For further information, visit

www.sartorius.com/en/products/flow-cytometry/flow-cytometry-software

Specifications subject to change without notice.

©2024. All rights reserved. All names of Sartorius products are registered trademarks and the property of Sartorius AG and | or one of its affiliated companies.

Status: 04 | 2024



Supporting Additional Run-Time on the iQue®

Extend uninterrupted instrument operation with the Qmax™ Refill Module.

- 5x the run time (extend sheath and waste levels from 7 hours to 48 hours)
- Optimize iQue® location with remote placement of carboys up to 6 meters away
- Never stop an experiment with hot-swapping capabilities

Catalog number 90208 Robotic Integration Automation Programming Interface (API) License

Catalog number 91309 QMax Refill Module (all configurations)

Catalog number 91310 Bacteriostatic Concentrate (100 mL)

Catalog number 91311 QMax Refill Module Sheath Carboy Filter