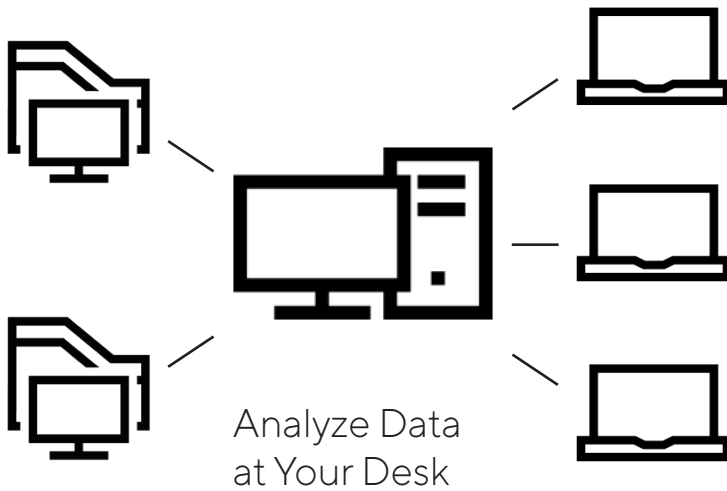


# SARTORIUS

## iQue Forecyt<sup>®</sup> Enterprise Unlimited



The Forecyt<sup>®</sup> Enterprise Unlimited solution syncs iQue<sup>®</sup> data from multiple instruments safely and seamlessly across a network. Secure remote access allows users and administrators to access their research without having to leave their desks. This access includes setting up experiment conditions before heading to the lab, performing analysis, and exporting results for presentations and papers. No longer will you have to wait for someone else to be done with the machine to review your data. Remove the need to be at the instrument itself when setting up experiments or to complete analysis.

## Enterprise Scales with Your Research

- Unlimited User Licences for everyone in your laboratory
- Collaboration via data sharing among multiple users
- Throughput enhancement by allowing experiments to be run on multiple instruments
- Scalability and accessibility via an enterprise-caliber software architecture
- Ease of deployment via simple installers and flexible physical IT layout
- Security and reliability via centralized data storage and encrypted network communication
- User Security through controlled project access and user level permissions



Forecyt<sup>®</sup> Enterprise supports iQue<sup>®</sup> usage across your organization

Researchers — everyday ease of use and collaboration

IT — provides data security and deployment flexibility

Cores — whether you are running a core or using a core, access can be granted between researcher and data directly

Screeners — scalability when using multiple instruments generating high volumes of data

## Architecture

Forecyt® Enterprise Unlimited uses a three-tier architecture, providing storage, access, and ensuring data integrity, while connecting directly with Clients for a seamless experience using Forecyt®. Users won't even know they aren't in front of the equipment.

Tier	Name	Purpose
Server	Data Storage	Physical storage of experimental data
	Data Access	Provides validation, security, and ensures data integrity
Client	Application	The user interface that controls data acquisition and analysis

## Deployment Options

Enterprise offers flexible options to support your organizational needs as the number of users and instruments grows. Enterprise removes the stress of deployment by providing multiple options accounting for cost, complexity, timing, and infrastructure.

- Dedicated Server: real or virtual server hosts the Forecyt® Enterprise Server components
- Dedicated Server + Network File Storage: FCS files are moved to a shared server or "drive"

## Windows Authentication Mode

Users are automatically logged onto the system with their companywide login – no additional passwords.

## Unlimited Access

Provide access to Forecyt® for everyone in your laboratory.

## Nightly Database Backup

By default, the Forecyt® Enterprise server performs a nightly database backup complimenting existing redundancy solutions - no more lost files.

## Distributed Licensing

Additional instrument modules for automation, signal reduction, and 21 CFR Part 11 compliance make Enterprise even more powerful.

## Forecyt® Tools and Software

Description	Order No.
iQueForecyt® Enterprise Unlimited	90212
Forecyt® Enterprise Lite	60036
iQue Forecyt® Enterprise Upgrade - Lite to Unlimited	BA-97136
Robotic Integration API License	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](mailto:info.intellicyt@sartorius.com)