

April, 2026

Keywords or phrases:

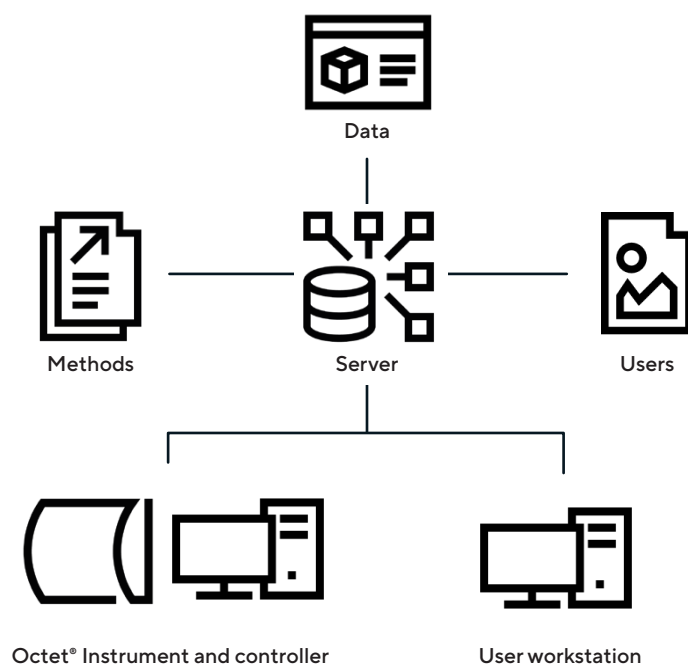
Octet® Hub; Octet® Software v14; Octet® Research Suite; Octet® Compliance Suite; Research Hub; Compliance Hub; Octet® BLI software; Bio Layer Interferometry; BLI data management; project based data management

Octet® Software Version 14 with Octet® Hub

Correspondence
E-Mail: octet@sartorius.com

Introduction

Octet® Software v14 introduces Octet® Hub, a project-based workspace backed by a central database. It brings methods, experiments, analysis settings, and results together in one place. Instead of managing work across local folders and duplicate file copies, teams organize their work as Hub projects and access shared assets from authorized workstations.



Octet® Hub connects methods, data, systems, and users in one shared workspace. Octet® BLI systems and user workstations connect to a central Hub, providing teams a single source of truth to store, find, and reuse methods and results.

Two Suites. Two Outcomes.

- **Research Suite + Research Hub:** Project-based collaboration, fast search, version history, and organized data for research teams.
- **Compliance Suite + Compliance Hub:** Includes all Research Suite capabilities and adds governance such as custom roles, audit-ready reporting, and workflow traceability) to support regulated workflows when configured and validated by the customer.

Capability	Octet® Research Hub	Octet® Compliance Hub
Centralized storage (methods, results, analyses)	Yes	Yes
Search across metadata and projects	Yes	Yes
Project-based organization (hierarchy)	Yes	Yes
Version history	Yes	Yes
Backup / restore	Yes	Yes
User groups and role-based access	Yes	Yes
Configurable security policies	Yes	Yes
Identity integration (Active Directory / SSO, where enabled)	Yes	Yes
Custom roles	No	Yes
Workflow traceability (method → experiment → analysis → report)	No	Yes
Audit-ready reporting (user activity, logs)	No	Yes
Access and role reports	No	Yes

Frequently Asked Questions

1. What problem does Octet® Software Version 14 with Octet® Hub solve?

Octet® Hub eliminates file and version confusion by replacing scattered local folders and duplicate copies with project-based, database-backed storage. Octet® Hub makes methods and results easy to find and reuse through project organization and search, and it scales from a single team to multi-user and multi-instrument environments.

2. What Changes with the Octet® Hub architecture?

Earlier Octet® workflows managed methods, experiments, and analysis data were managed as files, stored locally, and shared by copying. With the introduction of the Octet® Hub, work is organized as projects. Each project contains a managed collection of methods, runs, analyses, and results, so teams work from the same project context.

3. Can previous Octet® software versions support connect to the Octet® Hub database?

No. Octet® Hub works exclusively with Octet® Discovery v14 and Octet® Analysis Studio v14.

4. Is Octet® Hub cloud-based?

No. Octet® Hub is deployed on premises within the customer's environment (on-premise), either on a dedicated server or a designated workstation.

5. What does “project-based” mean in Octet® Software v14 and with the Octet® Hub?

In Octet® Hub projects are the top-level structure. Each project contains its own collection of records such as methods (and versions), experiments/runs, analysis settings, and results.

6. Can I migrate methods/experiments from earlier Octet® software versions?

Yes. Users can import existing methods and experiments into Hub projects using the Octet® Hub migration tool that were generated by Octet® Discovery Software and analyzed with Octet® Analysis Studio software.

Octet® Research Suite with the Octet® Research Hub

7. Who should use Octet® Research Suite?

Octet® Research Suite is ideal for discovery and assay development teams, startups, academic labs, and core facilities that want faster reuse and organization without the overhead of regulated governance.

8. Does Octet® Research Suite support identity integration and session timeout?

Yes. Octet® Research Suite includes identity integration options (e.g., Active Directory/Single Sign On (SSO), where enabled) and configurable session timeout controls.

9. Can multiple users share methods and results without making copies?

Yes. Octet® Research Suite users can access shared Octet® Hub projects and standardized assets based on permissions, reducing uncontrolled file copies.

Octet® Compliance Suite with the Octet® Compliance Hub

10. What does Octet® Compliance Suite with Octet® Compliance Hub add beyond Octet® Research Suite?

Octet® Compliance Suite includes all Research capabilities and adds governance features for regulated environments: custom roles in addition to predefined roles; audit logs and audit ready reports; end to end workflow traceability from method to experiment to analysis to reporting; and controls that support regulated workflows (for example, 21 CFR Part 11) when configured and validated by the customer.

11. Does using Octet® Compliance Suite automatically make my lab compliant?

No. Octet® Compliance Suite provides technical controls, including role-based access, audit-ready activity logging, workflow traceability, and controlled record handling, to support compliance with regulations such as FDA 21 CFR Part 11. Achieving compliance depends on how the system is configured and used within your organization's quality processes, including SOPs, training, change control, and validation.

12. What does Sartorius provide to help customers labs validate Octet® BLI System performance?

Sartorius provides documentation, qualification materials, and optional services that customers can use as part of their validation approach. Depending on the package and system, this may include:

- Instrument documentation (e.g., user guides)
- IQ/OQ kits and manuals (Installation/Operational Qualification)
- Software validation package (validation documentation to support software qualification activities)
- IQ/OQ service (optional service support for executing qualification)
- PQ kits and manuals (Performance Qualification), including PQ-K and PQ-Q kits where offered
- PQ services (e.g., PQ-Kinetics and PQ-Quantitation service) where offered
- PMOQ service (where offered)

Deployment and IT

13. What are the common deployment options?

- Option 1 (recommended): Shared Hub on an IT-managed server. Ideal for multi-user, multi-instrument, and regulated environments.
- Option 2: Local Hub on a workstation. Best for small teams or pilots with minimal infrastructure.

14. Do we need a dedicated server to run Octet® Hub?

A dedicated server (physical or virtual) is recommended for multi-user environments. Smaller groups may host on a designated workstation, but IT should align uptime expectations and backups.

15. What are the basic system requirements to host Octet® Hub?

Octet® Hub server requirements: Windows Server 2019, 2022, or 2025; 4 dedicated CPU cores; 8 GB RAM minimum/16 GB recommended; 250 GB storage minimum/512 GB recommended; customer network connection; customer-managed backup. Compliance Suite deployments require time synchronization, preferably using an NTP or domain time source.

16. Does Octet® Hub support HTTPS (TLS encryption)?

Yes. Octet® Hub supports HTTPS. Sartorius recommends configuring an SSL certificate, so communication uses HTTPS as a best practice.

17. Can we use a self-signed certificate for HTTPS?

Yes. The self-signed certificate must be trusted on the server and on each client computer that will connect to Octet® Hub.

18. What database does Octet® Hub use?

The Octet® Hub installer installs and configures a dedicated Microsoft SQL Express instance.

19. Can Octet® Hub integrate with Active Directory?

Yes. Octet® Hub supports Active Directory integration for authentication and user addition. If you plan to use AD integration, the Octet® Hub server must be joined to your domain.

What Else has Changed Since Version 13 of the Software?

20. What OS and upgrade prerequisites were introduced in v13.1?

Octet® Software v13.1 added support for Windows 10 (64-bit) and Windows 11; older Windows versions (e.g., Windows 7) are not supported. .NET Framework 4.8 is required for v13.1; offline PCs may need IT/support-assisted installation.

21. What were the key Octet® BLI Discovery improvements since v13?

Octet® Discovery added Octet® R8e instrument support, including expanded sampling rate options, increased the step recording time limit to 24 hours in 5 Hz mode for supported instruments, captured the time zone offset of the computer that created the file, and delivered additional templates (e.g., epitope binning templates and 384-well workflows) along with stability fixes.

22. What were the key Octet® Analysis Studio enhancements since v13?

Octet® Studio expanded sample and sensor information editing and CSV import capabilities, redesigned tables so exports match on-screen column order and sorting, introduced new kinetics capabilities (including the Two-State model), and delivered broad usability improvements.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
3874 Research Park Drive
Ann Arbor, MI 48108
Phone +1 734 769 1600

 For further contacts, visit
www.sartorius.com