

Octet® Hub

Centralized Data Management for Octet® Workflows



Octet® Software Version 14 | Research and Compliance Suites

Octet® Software Version 14 introduces Octet® Hub, a project-based workspace backed by a central database. It brings methods, projects, analysis settings, and results in a single, governed workspace, so teams can find what they need faster, collaborate without duplicating work, and reduce the risk of using the wrong version of a file.

Centralization of projects in the Octet® Hub removes the potential for work to be spread across several folders, shared drives, and email keeps, which helps teams align on a single project.

Octet® Software Version 14 is available in two suites:

- Octet® Research Suite with Research Hub
- Octet® Compliance Suite with Compliance Hub

Together, they provide a modern platform for both discovery research and regulated laboratory environments.

Key Features

- Hub-based data management: Centralize methods, experiments, analysis settings, and results in one source of truth.
- Project-based organization: Structure work by lab, team, or program for clearer ownership and collaboration.
- Version history: Track changes to methods and analysis settings and revert when needed to support reproducibility.
- Collaboration with control: Share within projects without creating uncontrolled copies; prevent simultaneous editing conflicts.
- Backup and restore: Admin-friendly workflows to protect data and simplify recovery.
- Two suites for two workflows:
 - Research Suite + Research Hub for productivity and collaboration.
 - Compliance Suite + Compliance Hub for governed audit-ready workflows in regulated environments.

What Teams Will Gain

- Spend less time searching for and recreating methods, runs, and results for each project.
- Reduce duplicate files and eliminate version confusion by reusing validated methods and data files, allowing for more confident work.
- Enhance consistency across teams by standardizing how work is organized and shared.
- Benefit from improved inspection readiness through streamlined traceability and audit reporting provided by the Octet® Compliance Suite in regulated environments.

Advantages of the Octet® Hub

Centralize & Find

With Octet® Hub, your methods, projects, and results live in one trusted place, so everyone works from the same, up-to-date information. Instead of navigating folders and shared drives, teams can quickly search across projects to which they have access finding exactly what they need. Validated methods are easy to reuse, results are simple to locate, and workflows stay consistent, helping scientists spend less time managing data and more time advancing their work.

Collaborate & Control

Collaborate efficiently while preserving data integrity. Octet® Hub can be used to organize data by project, giving teams a shared workspace where methods and analyses can be accessed securely, without uncontrolled copies. User groups and role-based access ensure the right people can contribute, while permission-based archiving and restoring support the full lifecycle of your projects. Built-in protections prevent accidental overwrites, so teams collaborate with confidence, clarity, and accountability.

Govern & Prove

Turn compliance into a natural part of your workflow. The Octet® Compliance Suite adds governance and traceability to the Octet® Compliance Hub, connecting every step from method to experiment to analysis and instrument. Automated activity tracking, role-based access, and session controls help protect sensitive data while supporting audit-ready reporting. These capabilities simplify inspection preparation and help teams confidently demonstrate data integrity.



Two Suites. Two Outcomes: Octet® Research Suite with the Octet® Research Hub

Move faster

Stay organized. Collaborate with confidence. Designed for research teams who want quicker access to data, clearer organization, and smoother collaboration, without the need for regulated controls. With Research Hub, methods and results live in one shared workspace, making it easy to find past work, reuse validated procedures, and keeping teams aligned.

Ideal for discovery labs, startups, academia, and core facilities.

What You Gain

- A centralized home for methods and results; no more scattered folders.
- Searchable projects and assets for faster reuse.
- Version history to support reproducibility.
- Simple sharing across teams without duplicate copies
- Identity integration options (e.g., Active Directory / SSO, where enabled).
- Backup and restore tools to help protect valuable work.

Octet® Compliance Suite with Octet® Compliance Hub

Bring governance into everyday workflows—and be ready when it matters.

Designed for laboratories operating in regulated environments, the Octet® Compliance Suite builds on the same Octet® Hub foundation by adding governance and traceability, supporting compliance with regulations such as FDA 21 CFR Part 11 when configured and validated by the customer.

Ideal for QC, QA, and regulated development environments.

Additional Capabilities Include

- Role-based access and project-level permissions.
- End-to-end workflow traceability and audit reporting.
- Access and role reports for accountability.
- Identity integration options (e.g., Active Directory / SSO, where enabled).
- Session management and configurable security policies to help protect sensitive work.
- Ability to create custom roles (Compliance Suite) for tailored governance needs.

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

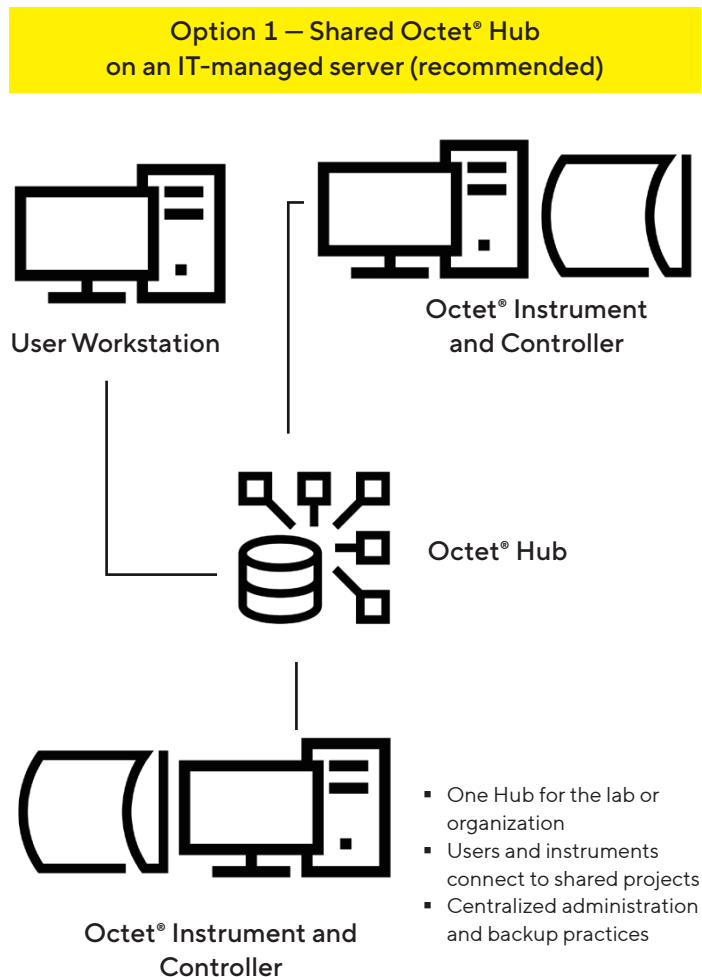
Deployment Options

Octet® Hub is deployed on customer-managed, on premise infrastructure and can support multiple deployment models, from a single user desktop workstation to a centralized server handling multiple work groups. The best option depends on how users collaborate, how IT manages infrastructure and security, and user needs. Two deployment options are discussed below.

Recommended guidance: Option 1 is best for multi-user teams and regulated environments; Option 2 is best for small teams and pilots with minimal infrastructure.

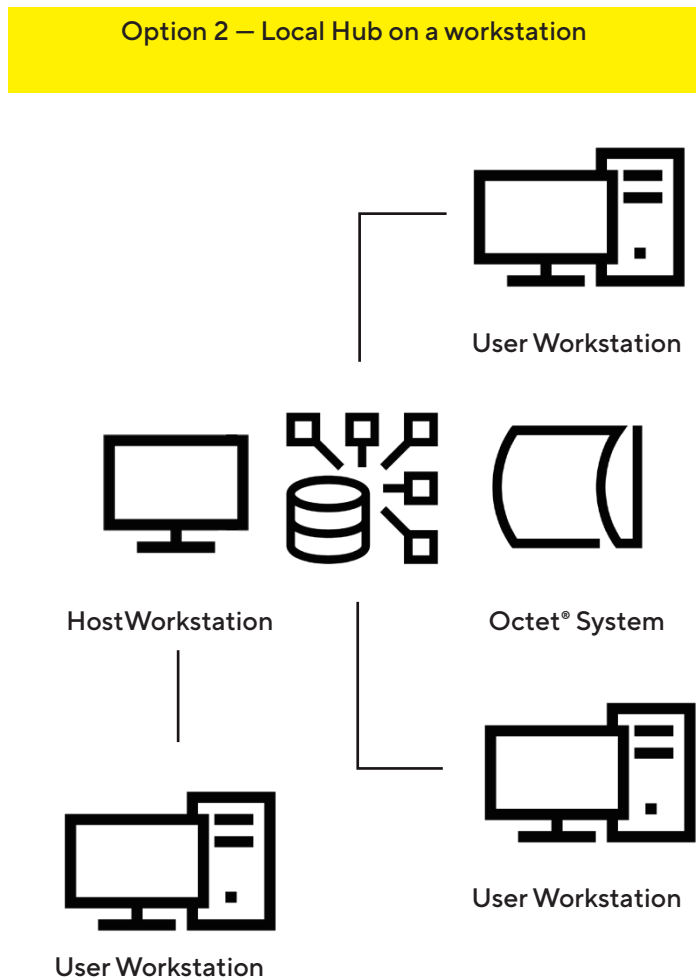
Option 1 – Shared Octet® Hub on an IT-managed server (recommended)

In this model, Octet® Hub is hosted on a dedicated server environment managed by the customer's IT team. Scientists and teams connect to the same Hub to access shared projects, standardized methods, and centralized results from any authorized workstation. For IT, this approach supports central administration, user access management, and the establishment of backup and security practices.



Option 2 – Local Hub on a workstation

For small teams or pilot deployments, Octet® Hub can run on a designated “host” workstation without standing up a separate server. Users in the same workgroup can still benefit from project-based organization and shared access with minimal infrastructure. Because the Hub resides on a workstation, IT and the lab should align on host PC resources, uptime expectations, and backup practices.



Technical Specifications

Server Computer Requirements – Octet® Hub		
Item	Minimum	Recommended
Operating system	Windows Server 2019, 2022 or 2025	Windows Server 2025
Processor	Dedicated 4 cores	Dedicated 4 cores
Memory	8 GB	16 GB
Storage	250 GB	512 GB
Network	Customer network with 100 Mbps required dependent on the deployment scenario	
Backup	Customer-managed	Standard IT backup policy
Time synchronization	Required for Compliance	NTP/domain time source recommended
Identity integration		Domain joined server for Active Directory/SSO login
Client Workstation Requirements		
Item	Minimum	Recommended
Operating system	Windows 10 64-bit Professional/ Enterprise or Windows 11	Windows 11
Processor	Dedicated 2 cores	Dedicated 4 cores
Memory	4 GB	16 GB
Display	1920 X 1080	1920 x 1080
Network connection	Required for Octet® Hub access	Stable LAN/VPN connection with 100 Mbps or higher
Additional Software		Microsoft® Excel
Octet® Instrument Controller Requirements		
Item	Requirement	
Operating system	Windows 10 Professional/ Enterprise or Windows 11	
Processor	Dedicated 4 cores	
Memory	16 GB	
Storage	512 GB	
Instrument connection	USB 3.0/RS-232	
Display	1920 X 1080 or better	
Network	Required for Octet® Hub access	
Laptop use	Not recommended	

Ordering Information

Octet® Research Suite

Order No.	Description
50-5030	Octet® Research Suite v14 Base – 5 Users. Octet® Research Suite v14 provides the complete Octet® software environment for data acquisition, analysis, and centralized data management, including Octet® BLI Discovery, Octet® Analysis Studio, and the Octet® Research Hub. - This package includes five user licenses ideal for collaborative research teams needing shared access to methods, experiments, and analysis results within a centralized Hub workspace. This product is also shipped with any new Octet® automated instrument. - Additional users can be added later using license expansion packages.
50-5200	Octet® Research Suite v14 Base – 1 User Octet® Research Suite v14 provides the complete Octet® software environment for data acquisition, analysis, and centralized data management, including Octet® BLI Discovery, Octet® Analysis Studio, and the Octet® Research Hub. - This package includes one user licenses designed for individual researchers starting on Octet® workflows. It establishes a Hub-based workspace for storing, searching, and organizing methods, experiments, and results. - Additional users can be added later using license expansion packages.
50-0400	Octet® Research Suite v14 – Add 5 Users Provides five additional user licenses for an existing Octet® Research Suite v14 installation. - This expansion package increases the number of authorized users who can access the Research Hub and associated Octet® software. This package can be purchased in multiples to support growing teams. - Requires an existing Octet® V14 Research Suite base license.

Octet® Compliance Suite

Order No.	Description
50-5035	Octet® Compliance Suite v14 Base – 5 Users Octet® Compliance Suite v14 includes Octet® BLI Discovery, Octet® Analysis Studio, and the Octet® Compliance Hub with governance, traceability, and access-control capabilities designed for regulated laboratory workflows. - This package includes five user licenses and provides centralized controlled data storage, role-based access, workflow traceability, and audit-ready reporting to support compliance with regulations such as 21 CFR Part 11 when configured and validated by the customer. - Additional users can be added later using license expansion packages.
50-5300	Octet® Compliance Suite v14 Base – 10 Users Octet® Compliance Suite v14 includes Octet® BLI Discovery, Octet® Analysis Studio, and the Octet® Compliance Hub with governance and traceability features for regulated environments. - This package includes ten user licenses, providing a larger initial deployment option for regulated laboratories or multi-user teams requiring centralized controlled workflows from the outset. - Supports compliance with regulations such as 21 CFR Part 11 when configured and validated by the customer. - Additional users can be added later using license expansion packages.
50-0401	Octet® Compliance Suite v14 – Add 5 Users Provides five additional user licenses for an existing Octet® Compliance Suite v14 installation. - This expansion package increases the number of authorized users while maintaining governance controls, traceability, and auditability required in regulated environments. - Requires an existing Octet® v14 Compliance Suite base license.

Germany

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

USA

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

 **For further information, visit**
[sartorius.com](https://www.sartorius.com)