

Security and Architecture Policy for Agilent SLIMS 6.7

Overview

Data security, privacy, and their various aspects, such as availability, integrity, and confidentiality, are of critical importance to our customers. Agilent commits to delivering the highest quality services to our customers, addressing security at all stages of the product development life cycle.

This document explains how Agilent SLIMS implements security measures that comply with assorted U.S., European and other regulations and standards. We have defined a set of policies, processes, and controls for security and privacy in accordance with the internationally accepted ISO 27001 and 27002 security standards.

The document is divided into three sections:

1. The platform architecture and implementation that enable us to deliver the highest level of security and to comply with the regulations
2. The security measures we have deployed to ensure data availability, data integrity, and data confidentiality
3. The applicable data privacy regulations and how we comply

Platform Architecture

This section describes the hardware, software, and technologies used to support the SLIMS platform and our customers.

Agilent SLIMS offers flexible installation options:

- SLIMS Customer Hosted: hosted by customer, either on-premises or on their own cloud solution
- SLIMS Agilent Hosted: hosted by Agilent as Software-As-A-Service on a scalable, powerful, secure server.

Technology Architecture

SLIMS relies on tools that provide a scalable, web-oriented architecture. It is highly configurable and is extensible through various APIs. Java-based plugins can be created and activated from several points within the system. Other interfaces such as a REST API allow for other applications to integrate with SLIMS. The system is suitable for demanding and high-profile customers that appreciate the system's stability and flexibility.

SLIMS team operations are furthermore based on best practices and a commitment to excellence. As documented in its corporate Quality Management Policy, the company promotes and supports a methodology that has achieved ISO 9001:2015 certification. For more details about Quality at Agilent, please visit www.agilent.com/quality.

SLIMS Platform

The SLIMS architecture is layered in the following way:

The **data** is stored in a relational database and files are stored in a dedicated file store. The data is exposed to the rest of the framework by the Platform API. This layer provides low-level data access while guaranteeing business logic such as traceability and authorization.

SLIMS Server is the server side of the main UI (User Interface). It is hosted on an application server (Apache Tomcat) and presents the Web UI to the user.

The SLIMS UI can be extended by a plugin framework based on **Vaadin**.

Plugins can be useful for creating dedicated data entry forms or portals or to visualize sample-specific data using pluggable external web tools.

SLIMS GATE is an integration layer based on Apache Camel, an Enterprise Service Bus. This technology is built for integration with a wide variety of systems using a variety of protocols (e.g., direct database access, file exchange, HL7, ftp) and can be easily extended.

SLIMS GATE furthermore includes XML configuration for defining small UI components that interact directly with the user from the SLIMS interface. Plugins can be installed and deployed from the user interface and there is a mechanism to pass parameters to the plugins. Another functionality plugins provide are file

upload integration, enabling instrument uploads to file sharing. In addition to user-triggered actions, programmed actions such as nightly dumps or imports are also achieved.

The **REST API** is a language-agnostic API that provides read/write access to all system tables (location, content, result, etc.) and the file repository. This can be used to perform tasks such as submitting orders from a customer portal and requesting sample information or order statuses.

A **Python API** is also available. This API makes it possible to achieve complex tasks using concise scripts. An open cookbook with examples is available at python cookbook on [GitHub](https://github.com) and provides examples on data manipulation, defining a multi-step flow interacting with the user, a live report, an example portal for order submission, and an example of data plotting.

Connections

SLIMS connects to instruments and equipment in the laboratory through industry standard communications protocols and interfaces. The SLIMS connection library can be used with a valid SLIMS subscription license.

Integrations

Integration with external systems such as EMR/EHR, environmental monitoring, billing, and CRM systems can be implemented via either file transfers (csv, xml, HL7) or APIs. Controlled vocabularies and industry standard ontologies are supported. The SLIMS team have encountered and integrated many instruments and a plugin library is available for easier implementations.

Barcode use

SLIMS use is optimized with the use of barcode scanners and printers. Scanners can facilitate data entry and automate repetitive tasks when combined with Macros. Label Templates can be edited in SLIMS so print operations can be automatically run at specific steps of the workflow.

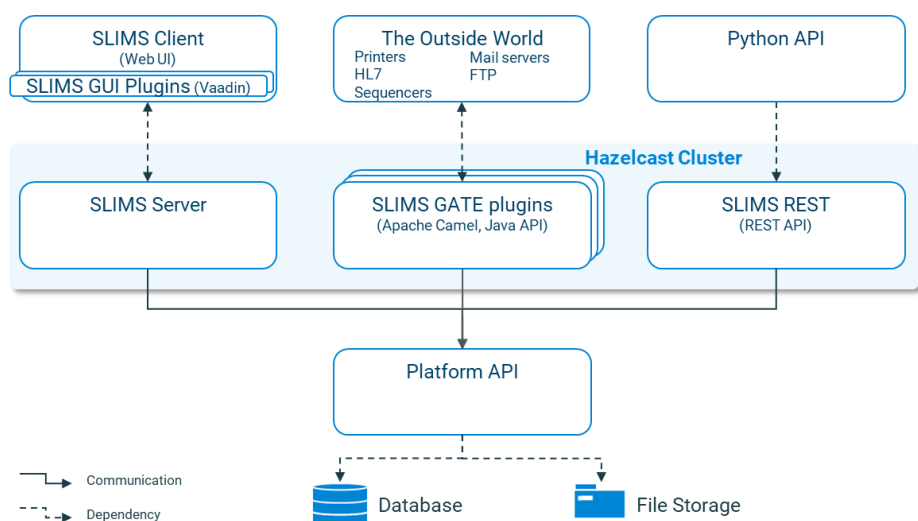


Figure 1. SLIMS architecture

Manual files import

External documents can be added to SLIMS in a variety of ways. Digital media in common formats can be associated in the various phases of the process via drag and drop. Scanned laboratory information such as equipment certification or results can be uploaded independently of a sample or order.

Generic Excel import

SLIMS supports importing from and exporting to XLS(X) files out of the box. This approach is used when the integration is not frequent, or the third-party system is flexible in accepting or generating flat files.

Configuration of flows for CSV or XML files

Common formats are supported with predefined scripts available for importing from or exporting to files that can be configured by simple field mapping. Fields from the CSV/XML are mapped to fields in SLIMS and vice versa. This approach is used when the third-party software is less flexible or frequent interaction is needed and the customer wants to avoid manual manipulation of files.

Dedicated import/export flows

Dedicated scripts can be created in the case of more complex operations. An example is importing a sample taken from a subject; checking if the subject exists, and either adding the sample to the existing one or creating a new one.

Create dedicated API integrations

This type of interaction is used when deeper integration is needed (e.g., to query the live status of an instrument).

Platform Implementations

Customer Hosted

This section applies to customers using the Customer hosted solution. The minimum server requirements for running SLIMS are listed in the *Agilent SLIMS 6.7 - Technical Requirements* document though the hardware specifications should be adjusted based on the actual usage patterns in the lab.

SLIMS runs on a web server that uses the HTTP(S) protocol and TCP/IP as its network transport. SLIMS needs to be installed within the same network range as the connected client computers, lab printers, instruments, and devices to connect with (either through physical network connections or via VPN bridges).

Agilent Hosted

This section applies to customers using SLIMS Agilent Hosted. Also, for customers that decide to use SLIMS Customer Hosted with SLIMS Remote, some aspects might be relevant.

The cloud infrastructure is a hybrid implementation between Single-Tenant and Multi-Tenant Cloud Hosting supported by Amazon Web Services (AWS). Each customer is assigned a dedicated database/file storage with platform HA (high availability) features. The SLIMS application layer, while dedicated to a single customer, is hosted as a containerized application to provide resource optimization and high availability.

The use of containers provides a

generic build for all customers which ensures they are replaceable and mobile across hardware with the previously stated advantages. Deployments are faster, environments are standardized, and are easily accessed by SLIMS personnel for maintenance. Customers can rely on our DevOps experts for hosting and maintaining the system.

For connectivity to local instruments, the cloud deployment requires the installation of a small service called SLIMS Remote that is downloaded from SLIMS. This service can be installed on Windows or Linux. It runs as a Java service and the workstation or server it is installed on must remain connected constantly. The SLIMS Remote service acts as a proxy for SLIMS GATE, ensuring connectivity between SLIMS and the local equipment by initiating REST calls to the server. This eliminates firewall configuration issues.

SLIMS Remote uses a plugin system to download functionality from its server. This way, downloading new functionality or updating existing functionality does not require reinstalling the service. This limits maintenance of the component.

Some limitations to the possible plugin integrations may apply for Agilent hosted versions of SLIMS. Customer hosted instances can also use the SLIMS Remote plugin system.

See also the full document describing the [SLIMS Cloud Technical Security Policy](#) with further specifications about our security measures and the [SLIMS Cloud Acceptable Use Policy](#).

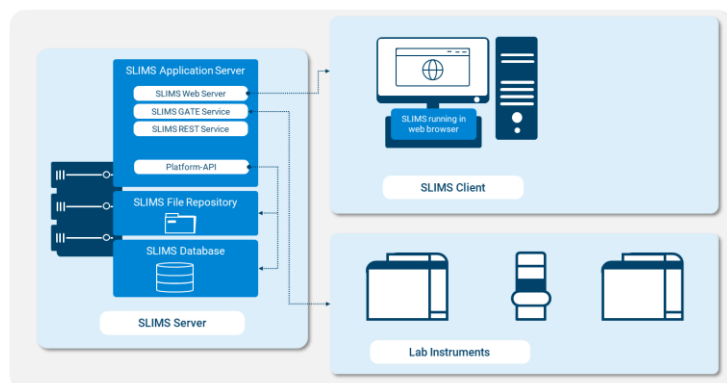


Figure 2. Customer Hosted SLIMS (One-Server Configuration)

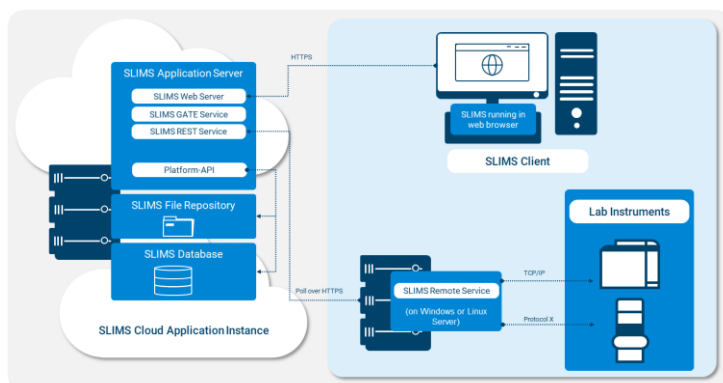


Figure 3. Agilent Hosted SLIMS (with Remote Service)

Security Measures

SLIMS Deployment Types and Responsibilities

SLIMS Agilent Hosted and SLIMS Customer Hosted each have different security responsibilities and implications.

| SLIMS Deployment Type | | |
|---|--|-----------------------|
| Responsibility | SLIMS Agilent Hosted | SLIMS Customer Hosted |
| Infrastructure provider | Agilent managed AWS infrastructure | Customer |
| Infrastructure managed by | Shared responsibility model between Agilent & AWS | Customer |
| Infrastructure security measures managed by | Shared responsibility model between Agilent & AWS | Customer |
| Secure SLIMS configuration & customization | Shared responsibility model between Agilent & Customer | |
| SLIMS customizations | Shared responsibility model between Agilent & Customer | |

Security Measures Overview

Availability

| Feature | SLIMS Agilent Hosted | SLIMS Customer Hosted |
|---------------------------------------|---|-------------------------|
| Facilities | AWS data centers demonstrate a strong physical security process, as acknowledged by their ISO 27001, ISO 27018, SOC1, SOC2, and SOC3 certifications. | Customer responsibility |
| Backup / restore | All AWS features used to store data (S3, EC2, EBS, RDS) are backed up and replicated in different physical availability zones; restore tests are performed at least once a year. | Customer responsibility |
| High Availability & Disaster Recovery | In case of a major incident in one of the physical availability zones (AZ) your SLIMS cloud instance will start in one of the other 2 physical AZs in the region. Application instance data is available in all physical AZs. | Customer responsibility |

Integrity

| Feature | SLIMS Agilent Hosted | SLIMS Customer Hosted |
|---------------------------|---|-------------------------|
| Public Key Infrastructure | All communications between client and application are protected by encryption provided by certificates issued by AWS CA. AWS PKI is used for encryption at rest of the customer data. | Customer responsibility |
| Data durability | Replicated and fault tolerant storage of database volume and backups in 3 separate physical availability zones (within one geographical region). | Customer responsibility |
| Hardware checks | All hardware underlying AWS services are permanently checked for failures and proactive migrations are performed. | Customer responsibility |
| Intrusion detection | Continuous scanning of incoming and outgoing traffic is performed with AWS Guard Duty, which triggers alarms in case of suspicious events. | Customer responsibility |

Confidentiality

| Feature | SLIMS Agilent Hosted | SLIMS Customer Hosted |
|---------------------|---|-------------------------|
| Authentication | SLIMS authentication options <ul style="list-style-type: none"> Local authentication Single Sign-On (SSO) with OpenID Connect (Okta, Google, Identity providers supporting OpenID Connect Discovery) OAuth 2.0 for API Clients SLIMS security options <ul style="list-style-type: none"> Configurable password complexity and lifetime Configurable failed login attempts before user lockout and session time-out | |
| | | LDAP integration |
| Authorization | SLIMS Role Based Access Control (RBAC) | |
| Isolation | Dedicated SLIMS database instance & dedicated SLIMS File Storage. Containerized SLIMS application on shared compute cluster with network policies in place to isolate traffic. | Customer responsibility |
| Encryption | <ul style="list-style-type: none"> Data at rest encrypted with AES-256 (database, file storage, backups) Encrypted database connection (TLS) Encrypted application endpoint (TLS), terminated at Application Load Balancer | Customer responsibility |
| Firewall | Packet filtering firewall for ingress HTTP(S) traffic. Network best practices (data storage on private network segments). | Customer responsibility |
| Intrusion detection | Automated anomaly detection is performed on infrastructure logs for suspicious events. | Customer responsibility |

Auditability

| Feature | SLIMS Agilent Hosted | SLIMS Customer Hosted |
|-------------------------------|---|--|
| Application data traceability | SLIMS keeps a detailed history of records on both data and configuration tables. The behavior of electronic signatures can be configured to regulate record changes. | |
| Logging | <ul style="list-style-type: none"> Application logging Access logging Infrastructure and security logging | <ul style="list-style-type: none"> Application logging Access logging Infrastructure and security logging are customer's responsibility |
| Change management | SLIMS development uses a Secure SDLC (Software Development Life Cycle). | |
| | Infrastructure changes are thoroughly tested, reviewed and version controlled. | Infrastructure managed by customer. |
| Incident management | SLIMS customer support and incidents managed in online Agilent Service Desk portal. | |
| | SLIMS infrastructure incidents are managed by Agilent and logged and managed in an internal and central system. An RCA (root cause analysis) is performed when the cause of the incident is unknown, to guarantee continuous improvement. | SLIMS infrastructure incidents managed by customer. |

Applicable security & data

privacy regulations

Agilent security team works with security experts to ensure that our platform is compliant with international regulations, and that it helps our customers comply with international standards.

GDPR

In May 2018, the European Union (EU) General Data Protection Regulation (GDPR) replaced the 1995 EU Data Protection Directive (European Directive 95/46/EC).

GDPR at Agilent is managed at the corporate level. Agilent has a comprehensive GDPR compliance program and provides a processing solution that incorporates the relevant GDPR requirements and that allows a customer to be assured that in choosing Agilent they are making a GDPR compliant choice. To facilitate our global business, we adhere to the requirements for safeguarding transfers of personal data internationally, including using Standard Contractual Clauses. Agilent personnel receive training on the GDPR and Agilent's obligations as both a data controller and a data processor to our customers.

HIPAA

The Health Insurance Portability and Accountability Act (HIPAA), enacted in 1996, established requirements for the protection and security of patient health information held by Covered Entities and Business Associates in the United States. HIPAA was expanded by the Health Information Technology for Economic and Clinical Health (HITECH) Act, as incorporated in the American Recovery and Reinvestment Act of 2009, to address increasing reliance on electronic maintenance and storage of patient health information.

The requirements of HIPAA/HITECH are contained in rules that include the:

- Privacy Rule: Protects the privacy of Protected Health Information (PHI) in any form (that is, written, recorded,

spoken orally, or electronic).

- Security Rule: Sets forth standards for the security—that is, the confidentiality, integrity, and availability—of PHI maintained in electronic form (known as ePHI) only.
- Breach Notification Rule: Requires Covered Entities and Business Associates to provide certain notifications following breaches of unsecured PHI.

Where Agilent is a Business Associate, we partner with our Covered Entity customers to ensure that appropriate HIPAA-compliant agreements and controls are in place.

ISO 27001

ISO/IEC 27001:2013 is an information-security standard that controls the following aspects of the security-management system of a company:

- Information-security policies
- Organization of information security
- Human resource security
- Asset management
- Access control
- Cryptography
- Physical and environmental security
- Operations security
- Communications security
- System acquisition, development, and maintenance
- Supplier relationships
- Information-security incident management
- Information-security aspects of business continuity management
- Compliance with internal requirements such as policies, and with external requirements such as laws

The SLIMS platform has been ISO 27001 certified by an independent auditor for its development, management, and support for its cloud deployment.

21 CFR Part 11

US FDA Part 11 in Title 21 of the Code of Federal Regulations (CFR), and its EU analog, Eudralex Chapter 4, Annex 11, describe the requirements for electronic records and electronic signatures for regulated pharmaceutical organizations. Released in 1997, 21 CFR Part 11 has been enforced since 1999. The intent of these guidelines is to ensure that all appropriate electronic records are attributable, legible, contemporaneous, original, accurate, and maintained with integrity.

A White Paper issued for each version of SLIMS provides a detailed description of how Agilent supports users and their organizations in achieving the requirements of each section of those regulations. The descriptions assume that system access, including instrument hardware and software, is controlled by the staff responsible for the electronic records contained on the system. Thus, the system is designed as a "closed system" as defined in 21 CFR Part 11.3(b)(4).

For more details, please refer to the resource *Support for Title 21 CFR Part 11 and Annex 11 compliance: Agilent SLIMS 6.7*.

Conclusion

Digital solutions such as SLIMS are an integral part of the laboratory. As described in this document, an extensive set of security considerations are put in place by Agilent during development, implementation, deployment, and support of SLIMS.

www.agilent.com/chem/agilentslims

DE44431.3478125

This information is subject to change without notice.

© Agilent Technologies, Inc. 2021
Printed in the USA, Aug 30, 2021
5994-4844EN

