



Declaration of Software Quality

We herewith inform you that the software product/systems

Product Name	Revision Number
Sample Scheduler for OpenLab	2.8

were developed according to the quality process and product development lifecycle established by the Life Sciences and Applied Markets Group (LSAG) of Agilent Technologies. Life cycle check-point details were reviewed and approved by the responsible management. The products were verified and validated to meet their functional and performance specifications and release criteria prior to ship release.

Users of this product may need to fulfill additional regulatory requirements according to applicable regulations and quality standards as defined by the User organizations. In support of our customers, Agilent Technologies will make design and supporting documentation of quality management system available to customer audit upon request at its facilities in Santa Clara, California (terms and conditions to be negotiated).

Agilent Technologies will maintain possession of all documents and their reproductions and may require a confidential disclosure agreement to be provided by those requiring access to these documents.

Date: February 2024

Quality Manager *Iris Feng*
Iris Feng (Feb 14, 2024 12:38 PST)
Iris Feng



软件质量声明

我们在此通知您软件产品/系统

产品名称	版本号
Sample Scheduler for OpenLab	2.8

是根据 Agilent Technologies 的 Life Science and Applied Markets Group (LSAG) 部门制定的软件质量管理流程和软件产品开发生命周期管理规范开发的，并且生命周期检查点的详细信息已由管理部门进行审核和批准。产品在发货前已经通过验证和确认，符合相关的功能和性能规格标准。

根据用户组织定义的适用法规和质量标准，本产品的用户可能需要满足其他合规要求。为支持客户，Agilent Technologies 可根据客户要求，在位于加利福尼亚州圣克拉拉的工厂提供相关质量管理体系的设计和支持文档供客户审核（条款和条件待商定）。

Agilent Technologies 将保留所有文档及其复制品的所有权，并可能需要提供保密协议才能访问这些文档。最终解释权以英文为准。

日期： 2024 年 2 月

质量经理：

[Iris Feng \(Feb 14, 2024 12:38 PST\)](#)

Iris Feng



Sample Scheduler for OpenLab_v2.8_DeclarationSoftwareQuality

Final Audit Report

2024-02-14

Created:	2024-02-14
By:	Katja Bertz (katja_bertz@agilent.com)
Status:	Signed
Transaction ID:	CBJCHBCAABAAyTg-9q9NG4lw79A7Cqja0lmlApjSQrj

"Sample Scheduler for OpenLab_v2.8_DeclarationSoftwareQuality" History

-  Document created by Katja Bertz (katja_bertz@agilent.com)
2024-02-14 - 1:40:18 PM GMT- IP address: 18.185.66.42
-  Document emailed to iris-nk_feng@agilent.com for signature
2024-02-14 - 1:41:14 PM GMT
-  Email viewed by iris-nk_feng@agilent.com
2024-02-14 - 8:36:28 PM GMT- IP address: 104.47.55.254
-  Signer iris-nk_feng@agilent.com entered name at signing as Iris Feng
2024-02-14 - 8:38:10 PM GMT- IP address: 4.53.28.20
-  Document e-signed by Iris Feng (iris-nk_feng@agilent.com)
Signature Date: 2024-02-14 - 8:38:12 PM GMT - Time Source: server- IP address: 4.53.28.20
-  Agreement completed.
2024-02-14 - 8:38:12 PM GMT