



Single Quadrupole LC/MS Acquisition
Release Notes

Notices

Document Identification

D0133350 Rev. A.00

July 2025

Copyright

© Agilent Technologies, Inc. 2025

No part of this manual may be reproduced in any form or by any means (including electronic storage and retrieval or translation into a foreign language) without prior agreement and written consent from Agilent Technologies, Inc. as governed by United States and international copyright laws.

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051

Software Revision

This guide is valid up to the 3.2 revision of Agilent Single Quadrupole LC/MS Acquisition.

Warranty

The material contained in this document is provided "as is," and is subject to being changed, without notice, in future editions. Further, to the maximum extent permitted by applicable law, Agilent disclaims all warranties, either express or implied, with regard to this manual and any information contained herein, including but not limited to the implied warranties of merchantability and fitness for a particular purpose. Agilent shall not be liable for errors or for incidental or consequential damages in connection with the furnishing, use, or performance of this document or of any information contained herein. Should Agilent and the user have a separate written agreement with warranty terms covering the material in this document that conflict with these terms, the warranty terms in the separate agreement shall control.

Technology Licenses

The hardware and/or software described in this document are furnished under a license and may be used or copied only in accordance with the terms of such license.

Restricted Rights Legend

U.S. Government Restricted Rights. Software and technical data rights granted to the federal government include only those rights customarily provided to end user customers. Agilent provides this customary commercial license in Software and technical data pursuant to FAR 12.211 (Technical Data) and 12.212 (Computer Software) and, for the Department of Defense, DFARS 252.227-7015 (Technical Data - Commercial Items) and DFARS 227.7202-3 (Rights in Commercial Computer Software or Computer Software Documentation).

Safety Notices

CAUTION

A **CAUTION** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in damage to the product or loss of important data. Do not proceed beyond a **CAUTION** notice until the indicated conditions are fully understood and met.

WARNING

A **WARNING** notice denotes a hazard. It calls attention to an operating procedure, practice, or the like that, if not correctly performed or adhered to, could result in personal injury or death. Do not proceed beyond a **WARNING** notice until the indicated conditions are fully understood and met.

Table of Contents

1	Introduction	4
	For our Regulated Customers	4
2	Version History	5
	Version 3.2	6
	Features	6
	Version 3.0	7
	Features	7
3	Important Support Information	8

1

Introduction

This document provides a listing of the major feature modifications made in each release of actively supported Single Quadrupole LC/MS Acquisition.



Documentation is found by scanning the code or navigating to <https://aglt.co/LCMSUserDocs>.

For our Regulated Customers

When any change is made to Agilent software, the validation status of the software needs to be re-established by the user/customer.

Whenever software is changed, a validation analysis should be conducted not just for the validation of an individual change, but also to determine the extent and impact of that change on the entire software system.



2

Version History

Version 3.2 6

Features 6

Version 3.0 7

Features 7

Version 3.2

After upgrading LC/MSD iQ systems with LC/MS SQ driver v3.2, the instrument must be reconfigured, and methods must be manually resolved to set the correct model name and instrument mode.

Features

- Support for new instruments: InfinityLab Pro iQ and InfinityLab Pro iQ Plus
- Mass-based fraction collection (LC/MSD iQ, Pro iQ, and Pro iQ Plus): Mass-based fraction collection support allows users to set up their methods to rapidly purify and characterize targets from complex mixtures, even at the smallest quantities, with high purity.
- Download MS instrument information from the instrument to your local computer (Pro iQ and Pro iQ Plus): Information includes factory autotune files, instrument performance reports (IDL Report), and regulatory conformity documentation.
- Calibrant valve switching during acquisition (CERNO MassWorks compatibility)
- Tune report enhancements
 - "Last tuned by" added to tune report
 - Tune ion abundances and ion profiles added to checktune report
 - Autotune and checktune reports saved in the data file

Version 3.0

Features

- Ramp the fragmentor voltage for Scan segments in the Acquisition Parameters table
- Calculation of Actual targeted points per second values. Migration of methods from previous versions is updated to change the actual points per second value.
- Lower ion source N2 consumption mode during standby
- Automatically launch an Autotune upon Checktune Out of Tolerance (LC/MSD iQ Only)
- Single polarity tuning (LC/MSD iQ Only)
- Extended detailed tune reports (LC/MSD iQ Only)

3

Important Support Information

For known issues and workarounds in the OpenLab software at the time of release see the Software Status Bulletins (SSB), which is found online at:

<https://www.agilent.com/cs/library/support/Patches/SSBs/M84xx.html>

Agilent Community

To get answers to your questions, join over 10,000 users in the Agilent Community. Review curated support materials organized by platform technology. Ask questions to industry colleagues and collaborators. Get notifications on new videos, documents, tools, and webinars relevant to your work.

<https://community.agilent.com/>

Agilent on the web

Visit Agilent's website [Lab Informatics Software - OpenLab Software Suite | Agilent](#) for up-to-date information.

www.agilent.com

© Agilent Technologies, Inc. 2025

D0133350 Rev. A.00

July 2025

