



MassHunter Acquisition for LC/TQ
Release Notes

Notices

Document Identification

D0131337 Revision A.00

May 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 for the 12.3 revision or higher of Agilent MassHunter Acquisition, for LC/TQ

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 12.3 6

Version 12.2 7

Restored Worklist Editing Capabilities 7

Real-Time Worklist Management 7

Enhanced Excel Integration 7

Performance Improvements 7

New Hardware Support 8

Version 12.1 9

Manual Tune File Support 9

Compound-by-Compound Optimization 9

Skyline Integration Enhancements 9

Real-Time Worklist Editing 9

Version 12.0 10

Updated Worklist Interface 10

Dual Configuration Support for Data Acquisition 10

Intelligent Reinjection Logic (iReflex) 11

Enforced Licensing Model 11

Role-Based Access Control 11

Comprehensive Audit Trails 11

3 Important Support Information 12

1

Introduction



This document provides a listing of the major feature modifications made in each release of actively supported MassHunter Acquisition Software including selected add-ons.

Instrument documentation, installation instructions, step by step videos, and more can be 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 12.3

Supported Software

- MassHunter Qualitative Analysis 13.0
- MassHunter Quantitative Analysis 12.1 Update 3
- Embedded Software 11.2.30; SC4 Firmware 9.12

Features

OpenLab ECM Support

- Compatible with OpenLab ECM 3.6 (with OpenLab Shared Services Server 2.7 Update 11 and later)
- Compatible with OpenLab Server / ECM XT 2.7 Update 11 and later

Version 12.2

Restored Worklist Editing Capabilities

- Reintroduces the comprehensive editing functionality from version 10.1, allowing users to modify a running worklist without interruption.

Real-Time Worklist Management

- Users can now add, delete, reorder entries, and import data from CSV files or existing worklists—all while the worklist is actively running.

Enhanced Excel Integration

- Improved copy/paste functionality enables seamless data exchange between the application and Microsoft Excel.

Performance Improvements

- Worklist editing operations are now significantly faster, providing a more responsive and efficient user experience.

New Hardware Support

- Adds compatibility with the InfinityLab Sample ID Reader, expanding the system's integration with laboratory hardware.

Version 12.1

Manual Tune File Support

- Users can now create and utilize manual tune files. This functionality is currently available for the 6495D instrument model and will be extended to additional models in the next release.

Compound-by-Compound Optimization

- In addition to existing compound and source optimization features, compound-by-compound optimization is now supported across all instrument models, enabling more granular control and improved performance.

Skyline Integration Enhancements

- Expanded support for Skyline automation workflows is now available for all instrument models, streamlining data analysis and reporting processes.

Real-Time Worklist Editing

- Users can now edit the worklist while it is running, including adding, deleting, and reordering entries, as well as importing from CSV or other worklists—enhancing flexibility and operational efficiency.

Version 12.0

Updated Worklist Interface

- The worklist user interface has been modernized with a ribbon-based layout, replacing the traditional menu structure for improved usability and navigation.

Dual Configuration Support for Data Acquisition

- Data Acquisition now supports two distinct configurations:

MassHunter Data Acquisition Workstation

- Data is stored locally with no content management capabilities.
- Authentication Mode can be toggled on or off via the Agilent Control Panel, with options for Internal or Domain authentication.

MassHunter Data Acquisition Networked Workstation

- Data is stored and managed within the Content Management component of OpenLab Server or ECM XT.
- ECM 3.x is also supported when used with OpenLab Shared Services Server.
- Authentication Mode is always enabled.
- Users can import existing Methods, Worklists, Worklist Templates, and Studies into projects managed by Content Management.

Intelligent Reinjection Logic (iReflex)

- Introduces iReflex, a data-dependent reinjection feature that uses real-time processing to automatically reinject samples based on predefined scenarios:
 - Carryover
 - Above Calibration Range
 - Fast Screening

Enforced Licensing Model

- Data Acquisition now includes enforced licensing. A 60-day startup license is provided for full-featured use. Continued operation beyond this period requires license registration.

Role-Based Access Control

- This release introduces user roles and privileges. Access to certain features in Data Acquisition is restricted based on assigned roles.

Comprehensive Audit Trails

- Audit trails are now maintained for Methods, Worklists, Data Files, and Studies in both Workstation and Networked Workstation configurations.
- A new Audit Trail Viewer provides a streamlined interface for reviewing entries and managing audit review workflows.

3

Important Support Information

For a list of defects fixed with this release, please see the Software Release Bulletin (SRB) included in the \Support\SSB and SRB\ folder.

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

<https://www.agilent.com/cs/library/support/Patches/SSBs/MHAcqLCTQ.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

<https://www.agilent.com/en/product/software-informatics/mass-spectrometry-software/data-acquisition/acquisition-for-lc-ms> for up-to-date information on Data Acquisition for LC/MS for up-to-date issue information on Agilent's Analytical Software suite.

www.agilent.com

© Agilent Technologies, Inc. 2025

D0131337

May 2025

