



MassHunter BioConfirm Software

Release Notes

Notices

Document Identification

D0145963 A.00

May 2026

Copyright

© Agilent Technologies, Inc. 2026

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
www.agilent.com

Software Revision

This guide is valid for the 14.0 revision or higher of Agilent MassHunter BioConfirm until superseded.

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.

Contents

1	Introduction	4
	For our Regulated Customers	6
2	Version History	7
	Version 14.0	8
	Supported Instruments	8
	Supported Software	8
	Supported Operating System	8
	Features	9
	Version 12.0	11
	Features	11
3	Important Support Information	12
	Overview	13
	Agilent Community	13
	Agilent on the web	13



1

Introduction

For our Regulated Customers

6

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



Instrument documentation, installation instructions, step-by-step videos, and more can be found by scanning the QR 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 14.0	8
Supported Instruments	8
Supported Software	8
Supported Operating System	8
Features	9
Version 12.0	11
Features	11

Version 14.0

MassHunter BioConfirm 14.0 introduces compatibility with OpenLab CDS 3.0 to read and process (Q-)TOF data and supports e-signature functionality. Users can open and process data files acquired on a (Q-)TOF using OpenLab CDS 3.0, use, modify, and generate methods that are compatible with OpenLab CDS 3.0, and operate BioConfirm in a Client/Server configuration with OpenLab CDS 3.0.

Supported Instruments

- Agilent LC/TOF
- Agilent LC/Q-TOF

Supported Software

- OpenLab CDS 3.0
- MassHunter Qualitative Analysis 14.0
- ChemVista

Supported Operating System

Support for Windows Server operating systems is limited to environments in which OpenLab CDS is installed first and provides the acquisition framework. In this configuration, the software operates as data analysis components within an OpenLab CDS Client/Server deployment.

MassHunter BioConfirm 14.0 is supported on Windows Server 2022 and Windows Server 2025 only when deployed as part of an OpenLab CDS Client/Server configuration.

Windows Server 2022 and Windows Server 2025 are not supported for MassHunter BioConfirm 14.0 in a MassHunter Acquisition-only (non-CDS)

environment. For standalone or non-CDS MassHunter Acquisition workflows, supported client operating systems must be used.

Features

- **Revident Data Processing Support:**
 - Enables processing of data collected using the Agilent Revident platform.
- **Distributed Database Architecture:**
 - A distributed database for Sequence Manager and the Chemical Data Dictionary allows all users in a Client/Server configuration to access the same sequence and chemical content (CDS Client/Server topology only).
- **OpenLab Reprocessor Tool:**
 - A reprocessing tool that supports batch processing of data collected using OpenLab CDS 3.0.
- **Regulatory Compliance Controls:**
 - Provides technical controls to securely acquire, process, report, and store data in compliance with FDA 21 CFR Part 11, EU Annex 11, GAMP 5, ISO/IEC 17025, and EPA 40 CFR Part 160.
- **PCDL (.cdbx) Format:**
 - A redesigned PCDL format that reduces database and library size and addresses software security concerns, with a standalone PCDL Converter tool supporting batch conversion of .cdb-format PCDL files.
- **Algorithm Changes:**
 - The integration algorithm used in BioConfirm has been updated in version 14.0. As a result, customers may observe differences in reported results for the following workflows:
 - Intact Protein
 - Oligos – Target Plus Impurities (Deconvolution)

Due to the updated integration behavior, the number or properties of detected biomolecules may differ from previous releases.

- **Ribbon UI Updates:**

- Home > Print includes Tune Report and Instrument Configuration Report.
- View tab includes a Sample Audit Trail window (visible with .sirslit data) and Bar Code Reader fields in the Sample Table.
- Processing > Actions: Find by tools (except Molecular Feature) include a drop-down for Entire Chromatogram or Over Selected Range.
- Tools > Export includes Acq (acquisition) Compound Info CSV and .d Conversion (conversion shown only with .sirslit data).

- **Method Editor Updates:**

- Method Automation > Report > Templates includes a Refresh button.

- **Data and Method Format Compatibility:**

- Supports both MassHunter (.d) and OpenLab CDS (.sirslit) data files. While these formats differ in structure and metadata, the underlying chromatographic and spectral data are equivalent.
- The application continues to support the traditional .m method format and introduces a new CDS native format— .bcpmx —which is fully interchangeable within the application.

Version 12.0

Features

- **Oligonucleotide Analysis Support:** First-time support for full-length product (FLP) confirmation and impurity profiling. MS/MS-based sequence confirmation dramatically reduces analysis time.
- **Workflow Automation:** Integration with MassHunter Acquisition and WalkUp for seamless data processing.
- **Unified Interface:** Consistent UI across workflows for proteins, peptides, glycans, and oligonucleotides.
- **Regulatory Compliance:** GxP-ready with audit trails, access control, and user roles.
- **Productivity Gains:** Reduced manual steps and training time; faster turnaround for results.



3 **Important Support Information**

Overview	13
Agilent Community	13
Agilent on the web	13

Overview

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 MassHunter BioConfirm 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/BioConfirm.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 web <https://www.agilent.com/en/product/software-informatics/mass-spectrometry-software/data-acquisition> for up-to-date information on Data Acquisition for LC/MS and GC/MS and for up-to-date issue information.

www.agilent.com

© Agilent Technologies, Inc. 2026

D0145963 A.00

May 2026

