



LC/(Q-)TOF Driver **Release Notes**

For OpenLab CDS Acquisition Software

Notices

Manual Part Number

D0147608

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This guide is valid for the 1.0 release or higher of the OpenLab CDS LC/ (Q-)TOF driver until superseded.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	Features	7
	Version 1.0	8
3	Important Support Information	10



1

Introduction

For our Regulated Customers

6

This document provides a listing of the major features included in release 1.0 of the OpenLab CDS LC/(Q-)TOF driver.



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 **Features**

Version 1.0

8

Version 1.0

Integration with the OpenLab CDS 3.0 Acquisition platform, including support for the following time-of-flight and quadrupole time-of-flight instruments, ion sources, and features:

- Supported instruments:
 - 6230C LC/MS (TOF)
 - G6545XT LC/MS (Q-TOF)
 - 6575 Revident LC/MS (Q-TOF)
- Supported ion sources:
 - ESI
 - Dual ESI
 - Dual AJS
 - Nanospray (not supported on Revident)
- Supported acquisition modes:
 - 6230C: MS, FPS
 - G6545XT: MS, FPS, All Ions, Targeted MS/MS, Directed MS/MS, Auto MS/MS
 - Revident: MS, FPS, All Ions, Q-RAI, Targeted MS/MS, Directed MS/MS, Auto MS/MS
- Application-based tune architecture
 - Simplified UI for routine users
 - Select and run multiple tunes at a time
 - Estimation of time to run multiple tunes
 - Scheduled tune (Revident only)
 - Editable description for application tunes
 - Diagnostic tune
- Early Maintenance Feedback (EMF)
 - Provides an easy-to-read dashboard to help visualize LC/MS instrument health and anticipate upcoming maintenance events.

- Scheduled Tune (Revident only)
 - Scheduled tuning alleviates the burden of routine maintenance and ensures that the mass spectrometer is always at peak performance.
- Iterative MS/MS support (Q-TOF only)
- Method Resolution support
- Single Sample Results data file (.sirsIt) format
- Distributed (Client/Server) environments
- Cloud and virtualized environments



3 Important Support Information

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 at <https://www.agilent.com/en/product/software-informatics/analytical-software-suite/chromatography-data-systems/openlab-cds> for up-to-date information on OpenLab CDS and the rest of Agilent's Analytical Software suite.

www.agilent.com

© Agilent Technologies, Inc. 2026

D0147608

May 2026

