



Agilent RapidFire Acquisition and Integrator Software

Version 6.3

Release Notes



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2010-2024

No part of this document 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.

Software Revision

Agilent RapidFire Acquisition and Integrator Software version 6.3

March, 2024

Acknowledgements

Microsoft ® and Windows are registered trademarks of Microsoft Corporation in the United States and other countries.

This product is intended for research use only. Not for use in diagnostics procedures.

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.

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.

Introduction

This document describes the functions, features and changes introduced with this release of the Agilent RapidFire Acquisition and Integrator Software. References to product documentation regarding known issues and workarounds are also provided. Software compatibility with Agilent MassHunter Acquisition Software versions are listed. Licensing information for programs used by RapidFire is included at the end of this document.

RapidFire Acquisition 6.3

Updates

- Added support for MassHunter Acquisition Software 12.0 for Revident LC/QTOF
- Added support for MassHunter Acquisition Software 12.1 for LC/TQ
- Added support for Windows 11 for RF 400 System

Supported RapidFire Models

| Product Number | Product Description |
|----------------|---|
| G9532AA | RapidFire 400 LC/MS System |
| G9531AA | RapidFire 365 LC/MS System (RoHS compliant) |
| G9530AA | RapidFire 365 LC/MS System (not RoHS compliant) |

MassHunter and Mass Spec Compatibility

| MassHunter Acquisition Software | LC/MS Instrument |
|---------------------------------|---|
| 12.1 for LC/TQ | See MassHunter Acquisition release for supported mass specs |
| 11.0 for LC/TOF and LC/QTOF | See MassHunter Acquisition release for supported mass specs |
| 12.0 for Revident LC/QTOF | See MassHunter Acquisition release for supported mass specs |

Operating System Compatibility

| Instrument | OS |
|---------------|--------------------------|
| RapidFire 365 | Windows 10 |
| RapidFire 400 | Windows 10 Windows 11 |

Software Dependency

The RapidFire Acquisition 6.3 Software requires Microsoft Excel (version >2013) to be fully functional. It is the customer's responsibility to procure licenses for Microsoft Excel prior to system installation.

Known Issues

RapidFire Acquisition 6.3

| Issue Number | Summary |
|--------------|--|
| 227069 | <p>Description: Software did not prompt user to save changes to the current method file when loading a different method file.</p> <p>Workaround: Save the changes to the current method file before opening another method file.</p> |
| 227641 | <p>Description: There was no “communication error” message when attempting to connect to a pump that was powered off.</p> <p>Workaround: Check to ensure the pump is turned on before establishing communication.</p> |
| 227650 | <p>Description: When auto scrolling down the Injection List Window, the highlight and cell content in the window were scrambled.</p> <p>Workaround: Minimize the window, and open up the window again. This will refresh the list.</p> |
| 243621 | <p>Description: Minimum gripper offset in the teach dialog did not take effect.</p> <p>Workaround: This parameter is not used to teach the Benchbot.</p> |
| 243633 | <p>Description: RapidFire remote connect was unable to see error messages (log entries) from the system.</p> <p>Workaround: Remotely access the desktop of the RapidFire computer to see the error message or map the directory of where the log entries are located.</p> |
| 471943 | <p>Description: When a run has been initiated and the system is waiting for MassSpec to warmup, user won't be able to abort the run for 10 minutes.</p> <p>Workaround: Shutdown RapidFireUI, click on Stop button on RapidFire Control Panel and start again.</p> |
| 539646 | <p>Description: RapidFire UI displays the currently loaded batch file from the <FileName> xml tag and not the actual file name</p> <p>Workaround: If a .bat file name is changed outside of RapidFire software, then perform a save As on the batch file using batch editor.</p> |
| 738112 | <p>Description: Default data directory in RapidFire.cfg file is incorrect.</p> <p>Workaround: Edit directory in RapidFire.cfg to preference.</p> |

| | |
|---------|---|
| 731472 | <p>Description: User Account Control window shows Unknown Publisher when uninstalling RapidFire Acquisition, Integrator, and Communicator.</p> <p>Workaround: None, continue uninstall.</p> |
| 1050523 | <p>Description If the Mass Spec is not the first module in the OpenLab Instrument (for example, following an LC), the first batch run of a TQ or QTOF will fail and display the message “Invalid MassSpec Detected”.</p> <p>Workaround: Delete the LC from the Instrument, or create a new Instrument containing only the MS.</p> |
| 1057056 | <p>Description Misleading error message when an unrelated axis has an error during a move.</p> <p>Workaround: If you see the message “Error : GeminiPci7344::check_move_complete_status on axis” followed by an axis number, ignore the axis number given in the error message, and inspect the Gemini logs to determine which axis actually had the error.</p> |

RapidFire Integrator 6.3

| Issue Number | Summary |
|--------------|--|
| 261366 | <p>Description: Software did not validate user input in the field “Metabolite Analysis”.</p> <p>Workaround: Enter a valid mass or formula.</p> |

Computer Requirements

RapidFire Acquisition software has specific computer requirements. Please contact Agilent Technologies Inc. for any questions regarding a computer for your RapidFire. RapidFire acquisition software requires Microsoft Excel 2013 or greater for some of its functionality to work.

Installing the RapidFire Software

Refer to the Installation Guide (part number D0027707) for detailed instructions on how to install and configure RapidFire software for RapidFire 365.

Refer to the Installation Guide (part number D0027707) for detailed instructions on how to install and configure RapidFire software for RapidFire 400.

Where To Find User Guides

The user guides for Agilent RapidFire products can be found on the RapidFire Resource App. This is a separate installer that comes with the instrument and software.

Once installed, the Resource App can be opened from the desktop. The Resource App gives you access to these guides:

- Quick Start Guide
- User Guide
- Data Analysis Guide
- Safety Guide

These guides are provided in PDF format and requires a PDF viewer to open. These guides are also available on www.agilent.com.

AGILENT TECHNOLOGIES, INC. LICENSE TERMS.

LEGAL NOTICE: PLEASE READ THESE TERMS BEFORE INSTALLING OR OTHERWISE USING THE SOFTWARE. ALL USE OF THE SOFTWARE IS SUBJECT TO THE LICENSE TERMS SET FORTH BELOW. "SOFTWARE" MEANS ONE OR MORE COMPUTER PROGRAMS IN EXECUTABLE FORMAT, WHETHER STAND-ALONE OR BUNDLED WITH OTHER PRODUCTS, AND RELATED DOCUMENTATION.

NOTWITHSTANDING ANYTHING TO THE CONTRARY IN THIS NOTICE, INSTALLING OR OTHERWISE USING THE SOFTWARE INDICATES YOUR ACCEPTANCE OF THESE TERMS. IF YOU DO NOT AGREE TO THESE LICENSE TERMS, DO NOT INSTALL THE SOFTWARE, OR DE-INSTALL THE SOFTWARE, AND YOU MAY RETURN THE SOFTWARE FOR A FULL REFUND.

AGILENT SOFTWARE LICENSE TERMS

The following License Terms govern your use of the Software unless you have a separate written agreement with Agilent, in which case, that written agreement will control and take precedence.

Software. "Software" means one or more computer programs in executable format, whether stand-alone or bundled with other products, and related documentation.

License Grant. Agilent grants you a non-exclusive license to use one copy of the Software for internal purposes in accordance with these license terms and the documentation provided with the Software. Such documentation may include license terms provided by Agilent's third party suppliers, which will apply to the use of the Software and take precedence over these license terms. In the absence of documentation specifying the applicable license, you may use one copy of the Software on one machine or instrument. If the software is licensed for concurrent or network use, you may not allow more than the maximum number of authorized users to access and use the software concurrently.

License Restrictions. You may make copies or adaptations of the Software only for archival purposes or only when copying or adaptation is an essential step in the authorized use of the Software. You must reproduce all copyright notices in the original Software on all permitted copies or adaptations. You may not copy the Software onto any public or distributed network.

Upgrades. This license does not entitle you to receive upgrades, updates or technical support. Such services may be purchased separately.

Ownership. The Software is owned and copyrighted by Agilent or its third party suppliers. Agilent and its third party suppliers retain all right, title and interest in the Software. Agilent's third party suppliers may protect their rights in the Software in the event of any violation of these License Terms.

No Disassembly. You may not disassemble or otherwise modify the Software without written authorization from Agilent, except as permitted by law. Upon request, you will provide Agilent with reasonably detailed information regarding any permitted disassembly or modification.

High Risk Activities. The Software is not specifically designed, manufactured or intended for use in the planning, construction, maintenance or direct operation of a nuclear facility, nor for use in on line control or fail safe operation of aircraft navigation, control or communication systems, weapon systems or direct life support systems.

Transfer. You may transfer the license granted to you here provided that you deliver all copies of the Software to the transferee along with these License Terms. The transferee must accept these License Terms as a condition to any transfer. Your license to use the Software will terminate upon transfer.

Termination. Agilent may terminate your license upon notice for breach of these License Terms. Upon termination, you must immediately destroy all copies of the Software.

Export Requirements. If you export, re-export or import Software, technology or technical data licensed hereunder, you assume responsibility for complying with applicable laws and regulations and for obtaining required export and import authorizations. Agilent may terminate this license immediately if you are in violation of any applicable laws or regulations.

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).

Additional program licensing information:

RapidFire uses ZedGraph Class Library and Visual Leak Detector (Vld) Library that are covered by the GNU General Public License v2 found below.

RapidFire also used Antlr software for development.

GNU GENERAL PUBLIC LICENSE

Version 2, June 1991

Copyright (C) 1989, 1991 Free Software Foundation, Inc.
51 Franklin Street, Fifth Floor, Boston, MA 02110-1301, USA

Everyone is permitted to copy and distribute verbatim copies of this license document, but changing it is not allowed.

Preamble

The licenses for most software are designed to take away your freedom to share and change it. By contrast, the GNU General Public License is intended to guarantee your freedom to share and change free software--to make sure the software is free for all its users. This General Public License applies to most of the Free Software Foundation's software and to any other program whose authors commit to using it. (Some other Free Software Foundation software is covered by the GNU Lesser General Public License instead.) You can apply it to your programs, too.

When we speak of free software, we are referring to freedom, not price. Our General Public Licenses are designed to make sure that you have the freedom to distribute copies of free software (and charge for this service if you wish), that you receive source code or can get it if you want it, that you can change the software or use pieces of it in new free programs; and that you know you can do these things.

To protect your rights, we need to make restrictions that forbid anyone to deny you these rights or to ask you to surrender the rights. These restrictions translate to certain responsibilities for you if you distribute copies of the software, or if you modify it.

For example, if you distribute copies of such a program, whether gratis or for a fee, you must give the recipients all the rights that you have. You must make sure that they, too, receive or can get the source code. And you must show them these terms so they know their rights.

We protect your rights with two steps: (1) copyright the software, and (2) offer you this license which gives you legal permission to copy, distribute and/or modify the software.

Also, for each author's protection and ours, we want to make certain that everyone understands that there is no warranty for this free software. If the software is modified by someone else and passed on, we want its recipients to know that what they have is not the original, so that any problems introduced by others will not reflect on the original authors' reputations.

Finally, any free program is threatened constantly by software patents. We wish to avoid the danger that redistributors of a free program will individually obtain patent licenses, in effect making the program proprietary. To prevent this, we have made it clear that any patent must be licensed for everyone's free use or not licensed at all.

The precise terms and conditions for copying, distribution and modification follow.

TERMS AND CONDITIONS FOR COPYING, DISTRIBUTION AND MODIFICATION

0. This License applies to any program or other work which contains a notice placed by the copyright holder saying it may be distributed under the terms of this General Public License. The "Program", below, refers to any such program or work, and a "work based on the Program" means either the Program or any derivative work under copyright law: that is to say, a work containing the Program or a portion of it, either verbatim or with modifications

and/or translated into another language. (Hereinafter, translation is included without limitation in the term "modification".) Each licensee is addressed as "you".

Activities other than copying, distribution and modification are not covered by this License; they are outside its scope. The act of running the Program is not restricted, and the output from the Program is covered only if its contents constitute a work based on the Program (independent of having been made by running the Program). Whether that is true depends on what the Program does.

1. You may copy and distribute verbatim copies of the Program's source code as you receive it, in any medium, provided that you conspicuously and appropriately publish on each copy an appropriate copyright notice and disclaimer of warranty; keep intact all the notices that refer to this License and to the absence of any warranty; and give any other recipients of the Program a copy of this License along with the Program.

You may charge a fee for the physical act of transferring a copy, and you may at your option offer warranty protection in exchange for a fee.

2. You may modify your copy or copies of the Program or any portion of it, thus forming a work based on the Program, and copy and distribute such modifications or work under the terms of Section 1 above, provided that you also meet all of these conditions:

a) You must cause the modified files to carry prominent notices stating that you changed the files and the date of any change.

b) You must cause any work that you distribute or publish, that in whole or in part contains or is derived from the Program or any part thereof, to be licensed as a whole at no charge to all third parties under the terms of this License.

c) If the modified program normally reads commands interactively when run, you must cause it, when started running for such interactive use in the most ordinary way, to print or display an announcement including an appropriate copyright notice and a notice that there is no warranty (or else, saying that you provide a warranty) and that users may redistribute the program under these conditions, and telling the user how to view a copy of this License. (Exception: if the Program itself is interactive but does not normally print such an announcement, your work based on the Program is not required to print an announcement.)

These requirements apply to the modified work as a whole. If identifiable sections of that work are not derived from the Program, and can be reasonably considered independent and separate works in themselves, then this License, and its terms, do not apply to those sections when you distribute them as separate works. But when you distribute the same sections as part of a whole which is a work based on the Program, the distribution of the whole must be on the terms of this License, whose permissions for other licensees extend to the entire whole, and thus to each and every part regardless of who wrote it. Thus, it is not the intent of this section to claim rights or contest your rights to work written entirely by you; rather, the intent is to exercise the right to control the distribution of derivative or collective works based on the Program. In addition, mere aggregation of another work not based on the Program with the Program (or with a work based on the Program) on a volume of a storage or distribution medium does not bring the other work under the scope of this License.

3. You may copy and distribute the Program (or a work based on it, under Section 2) in object code or executable form under the terms of Sections 1 and 2 above provided that you also do one of the following:

a) Accompany it with the complete corresponding machine-readable source code, which must be distributed under the terms of Sections 1 and 2 above on a medium customarily used for software interchange; or,

b) Accompany it with a written offer, valid for at least three years, to give any third party, for a charge no more than your cost of physically performing source distribution, a complete machine-readable copy of the corresponding source code, to be distributed under the terms of Sections 1 and 2 above on a medium customarily used for software interchange; or,

c) Accompany it with the information you received as to the offer to distribute corresponding source code. (This alternative is allowed only for noncommercial distribution and only if you received the program in object code or executable form with such an offer, in accord with Subsection b above.)

The source code for a work means the preferred form of the work for making modifications to it. For an executable work, complete source code means all the source code for all modules it contains, plus any associated interface definition files, plus the scripts used to control compilation and installation of the executable. However, as a

special exception, the source code distributed need not include anything that is normally distributed (in either source or binary form) with the major components (compiler, kernel, and so on) of the operating system on which the executable runs, unless that component itself accompanies the executable.

If distribution of executable or object code is made by offering access to copy from a designated place, then offering equivalent access to copy the source code from the same place counts as distribution of the source code, even though third parties are not compelled to copy the source along with the object code.

4. You may not copy, modify, sublicense, or distribute the Program except as expressly provided under this License. Any attempt otherwise to copy, modify, sublicense or distribute the Program is void, and will automatically terminate your rights under this License. However, parties who have received copies, or rights, from you under this License will not have their licenses terminated so long as such parties remain in full compliance.

5. You are not required to accept this License, since you have not signed it. However, nothing else grants you permission to modify or distribute the Program or its derivative works. These actions are prohibited by law if you do not accept this License. Therefore, by modifying or distributing the Program (or any work based on the Program), you indicate your acceptance of this License to do so, and all its terms and conditions for copying, distributing or modifying the Program or works based on it.

6. Each time you redistribute the Program (or any work based on the Program), the recipient automatically receives a license from the original licensor to copy, distribute or modify the Program subject to these terms and conditions. You may not impose any further restrictions on the recipients' exercise of the rights granted herein. You are not responsible for enforcing compliance by third parties to this License.

7. If, as a consequence of a court judgment or allegation of patent infringement or for any other reason (not limited to patent issues), conditions are imposed on you (whether by court order, agreement or otherwise) that contradict the conditions of this License, they do not excuse you from the conditions of this License. If you cannot distribute so as to satisfy simultaneously your obligations under this License and any other pertinent obligations, then as a consequence you may not distribute the Program at all. For example, if a patent license would not permit royalty-free redistribution of the Program by all those who receive copies directly or indirectly through you, then the only way you could satisfy both it and this License would be to refrain entirely from distribution of the Program.

If any portion of this section is held invalid or unenforceable under any particular circumstance, the balance of the section is intended to apply and the section as a whole is intended to apply in other circumstances.

It is not the purpose of this section to induce you to infringe any patents or other property right claims or to contest validity of any such claims; this section has the sole purpose of protecting the integrity of the free software distribution system, which is implemented by public license practices. Many people have made generous contributions to the wide range of software distributed through that system in reliance on consistent application of that system; it is up to the author/donor to decide if he or she is willing to distribute software through any other system and a licensee cannot impose that choice.

This section is intended to make thoroughly clear what is believed to be a consequence of the rest of this License.

8. If the distribution and/or use of the Program is restricted in certain countries either by patents or by copyrighted interfaces, the original copyright holder who places the Program under this License may add an explicit geographical distribution limitation excluding those countries, so that distribution is permitted only in or among countries not thus excluded. In such case, this License incorporates the limitation as if written in the body of this License.

9. The Free Software Foundation may publish revised and/or new versions of the General Public License from time to time. Such new versions will be similar in spirit to the present version, but may differ in detail to address new problems or concerns.

Each version is given a distinguishing version number. If the Program specifies a version number of this License which applies to it and "any later version", you have the option of following the terms and conditions either of that version or of any later version published by the Free Software Foundation. If the Program does not specify a version number of this License, you may choose any version ever published by the Free Software Foundation.

10. If you wish to incorporate parts of the Program into other free programs whose distribution conditions are different, write to the author to ask for permission. For software which is copyrighted by the Free Software Foundation, write to the Free Software Foundation; we sometimes make exceptions for this. Our decision will be guided by the two goals of preserving the free status of all derivatives of our free software and of promoting the sharing and reuse of software generally.

NO WARRANTY

11. BECAUSE THE PROGRAM IS LICENSED FREE OF CHARGE, THERE IS NO WARRANTY FOR THE PROGRAM, TO THE EXTENT PERMITTED BY APPLICABLE LAW. EXCEPT WHEN OTHERWISE STATED IN WRITING THE COPYRIGHT HOLDERS AND/OR OTHER PARTIES PROVIDE THE PROGRAM "AS IS" WITHOUT WARRANTY OF ANY KIND, EITHER EXPRESSED OR IMPLIED, INCLUDING, BUT NOT LIMITED TO, THE IMPLIED WARRANTIES OF MERCHANTABILITY AND FITNESS FOR A PARTICULAR PURPOSE. THE ENTIRE RISK AS TO THE QUALITY AND PERFORMANCE OF THE PROGRAM IS WITH YOU. SHOULD THE PROGRAM PROVE DEFECTIVE, YOU ASSUME THE COST OF ALL NECESSARY SERVICING, REPAIR OR CORRECTION.

12. IN NO EVENT UNLESS REQUIRED BY APPLICABLE LAW OR AGREED TO IN WRITING WILL ANY COPYRIGHT HOLDER, OR ANY OTHER PARTY WHO MAY MODIFY AND/OR REDISTRIBUTE THE PROGRAM AS PERMITTED ABOVE, BE LIABLE TO YOU FOR DAMAGES, INCLUDING ANY GENERAL, SPECIAL, INCIDENTAL OR CONSEQUENTIAL DAMAGES ARISING OUT OF THE USE OR INABILITY TO USE THE PROGRAM (INCLUDING BUT NOT LIMITED TO LOSS OF DATA OR DATA BEING RENDERED INACCURATE OR LOSSES SUSTAINED BY YOU OR THIRD PARTIES OR A FAILURE OF THE PROGRAM TO OPERATE WITH ANY OTHER PROGRAMS), EVEN IF SUCH HOLDER OR OTHER PARTY HAS BEEN ADVISED OF THE POSSIBILITY OF SUCH DAMAGES.

END OF TERMS AND CONDITIONS