



Agilent Technologies, Inc. Software License Agreement

ATTENTION: DOWNLOADING, COPYING, PUBLICLY DISTRIBUTING, OR USING THIS SOFTWARE IS SUBJECT TO THE AGREEMENT SET FORTH BELOW.

Software: "Software" means one or more computer programs in object code format, whether stand-alone or bundled with other products, and related documentation. It does NOT include programs in source code format.

License grant: Agilent grants you a non-exclusive license to download one copy of the Software, and to store or run that copy of the Software for internal use and purposes in accordance with this Agreement and the documentation provided with the Software. Such documentation may include license terms provided by Agilent 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 store or run one copy of the Software on one machine or instrument. If the Software is otherwise 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 for archival purposes or when copying or adaptation is an essential step in the authorized use of the Software, but for no other purpose. 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 and all copies thereof are 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 and its third party suppliers may protect their respective rights in the Software in the event of any violation of this Agreement.

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 written, 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 this Agreement, and pay any applicable fees to the extent permissible under local laws. The transferee must accept this Agreement as a condition to any transfer. Your license to use the Software will terminate upon transfer.

Termination: Agilent may terminate this license upon notice for breach of this Agreement. 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).

NO WARRANTY: TO THE EXTENT ALLOWED BY LOCAL LAW, AND EXCEPT TO THE EXTENT AGILENT HAS PROVIDED A SPECIFIC WRITTEN WARRANTY APPLICABLE TO THIS PRODUCT, THIS SOFTWARE IS PROVIDED TO YOU "AS IS" WITHOUT WARRANTIES OR CONDITIONS OF ANY KIND, WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED. AGILENT SPECIFICALLY DISCLAIMS ANY IMPLIED WARRANTIES OR CONDITIONS OF MERCHANTABILITY, SATISFACTORY QUALITY, NON-INFRINGEMENT AND FITNESS FOR A PARTICULAR PURPOSE. SHOULD THE SOFTWARE PROVE DEFECTIVE, YOU ASSUME THE ENTIRE RISK AND COST RESULTING FROM OR RELATING TO THE DEFECT. SOME JURISDICTIONS DO NOT ALLOW EXCLUSIONS OF IMPLIED WARRANTIES OR CONDITIONS, SO THE ABOVE EXCLUSION MAY NOT APPLY TO YOU. YOU MAY HAVE OTHER RIGHTS THAT VARY ACCORDING TO LOCAL LAW.

LIMITATION OF LIABILITY: TO THE EXTENT ALLOWED BY LOCAL LAW, IN NO EVENT WILL AGILENT OR ITS SUBSIDIARIES, AFFILIATES OR SUPPLIERS BE LIABLE FOR DIRECT, SPECIAL, INCIDENTAL, CONSEQUENTIAL, OR OTHER DAMAGES (INCLUDING LOST PROFIT, LOST DATA, OR DOWNTIME COSTS), ARISING OUT OF THE USE, INABILITY TO USE, OR THE RESULTS OF USE OF THE SOFTWARE, WHETHER BASED IN WARRANTY, CONTRACT, TORT OR OTHER LEGAL THEORY, AND WHETHER OR NOT ADVISED OF THE POSSIBILITY OF SUCH DAMAGES. YOUR USE OF THE SOFTWARE IS ENTIRELY AT YOUR OWN RISK. SOME JURISDICTIONS DO NOT ALLOW THE EXCLUSION OR LIMITATION OF LIABILITY FOR DAMAGES, SO THE ABOVE LIMITATION MAY NOT APPLY TO YOU.

Applicable law: Disputes arising in connection with this Agreement will be governed by the laws of the United States and of the State of New York, without regard to choice of law provisions. The United Nations Convention for Contracts for the International Sale of Goods will not apply to this Agreement.

Unenforceability: To the extent that any provision of this Agreement is determined to be illegal or unenforceable, the remainder of this Agreement will remain in full force and effect.

Entire agreement: This Agreement constitutes the entire agreement between you and Agilent, and supersedes any previous communications, representations, or agreements between the parties, whether oral or written, regarding transactions hereunder except for a specific warranty issued by Agilent with regard to this product. Your additional or different terms and conditions will not apply. This Agreement may not be changed except by an amendment signed by an authorized representative of each party.



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.

© Agilent Technologies, Inc. 2017
Printed in USA
Agilent Technologies, Inc.
2850 Centerville Road
Wilmington, DE 19808-1610 USA
For Research Use Only
Not for use in diagnostic procedures.