

MOLECULAR DEVICES, LLC
END USER SOFTWARE LICENSE AGREEMENT
and
LIMITED PRODUCT WARRANTY
GenePix Pro Software v7.3

NOTICE TO USER: PLEASE READ THIS DOCUMENT CAREFULLY. THIS IS THE CONTRACT BETWEEN YOU AND MOLECULAR DEVICES REGARDING THIS SOFTWARE PRODUCT. THIS AGREEMENT INCLUDES LICENSE TERMS AND RESTRICTIONS, WARRANTY AND LIABILITY DISCLAIMERS, AND LIMITATIONS.

YOUR INSTALLATION AND USE OF THIS MOLECULAR DEVICES SOFTWARE IS SUBJECT TO THE TERMS AND CONDITIONS CONTAINED IN THIS END USER SOFTWARE LICENSE AGREEMENT, AND YOU WILL BE BOUND BY THESE TERMS AND CONDITIONS IF YOU INSTALL OR USE THE SOFTWARE.

IF YOU AGREE TO THESE TERMS AND CONDITIONS, SELECT THE "I AGREE" OR "ACCEPT" BUTTON BELOW. IF YOU DO NOT AGREE TO THESE TERMS AND CONDITIONS, YOU SHOULD SELECT THE "I DO NOT AGREE" OR "DECLINE BUTTON" BELOW, WHEREUPON INSTALLATION WILL STOP, AND THEN PROMPTLY RETURN THIS SOFTWARE, TOGETHER WITH ALL PACKAGING, TO MOLECULAR DEVICES AND YOUR LICENSE PRICE FOR THIS SOFTWARE WILL BE REFUNDED.

IF THIS SOFTWARE PRODUCT IS INSTRUMENT OPERATING SOFTWARE OR SOFTWARE BUNDLED AND LOADED WITH INSTRUMENT OPERATING SOFTWARE, YOUR QUOTATION FROM MOLECULAR DEVICES WOULD HAVE REFERENCED THIS END USER SOFTWARE LICENSE AGREEMENT OR IT WOULD HAVE BEEN REFERENCED ON MOLECULAR DEVICES' WEBSITE (ERROR! HYPERLINK REFERENCE NOT VALID., "SALE AND LICENSE TERMS"), FOR YOUR INFORMATION AND AGREEMENT TO ITS TERMS AS A PART OF YOUR PURCHASE OF THE INSTRUMENT. IF YOU HAVE ANY QUESTIONS ABOUT THIS AGREEMENT, PLEASE CONTACT YOUR MOLECULAR DEVICES SALES OR SERVICE REPRESENTATIVE PROMPTLY.

This Molecular Devices End User Software License Agreement (EULA) accompanies a Molecular Devices software product ("Software") and related explanatory materials ("Documentation"). The term "Software" also includes any upgrades, modified versions, updates, additions and copies of the Software licensed to you by Molecular Devices. The term "Molecular Devices," as used in this License, means Molecular Devices, LLC. and its subsidiaries and affiliates. The term "License" or "Agreement" means this EULA including any and all Appendixes. The term "you" or "Licensee" means the purchaser of this license to the Software.

A copy of this End User Software License Agreement is available as either (i) an .rtf file on the installation media or (ii) a printable text file. To print the text file, select the "Print" button at the end of this Agreement.

THIRD PARTY PRODUCTS

This Software may include software products licensed by third party provider(s). In addition, the Software may enable you to run multiple instances of third-party software and application programs. You are responsible for reviewing and complying with any licenses necessary to use any such third-party software. Such licenses may be included in your Molecular Devices installation package.

TITLE

Title, ownership rights and intellectual property rights in and to the Software and Documentation shall at all times remain with Molecular Devices and its subsidiaries, and their suppliers. All rights not specifically granted by this License, including Federal and international copyrights, are reserved by Molecular Devices or their respective owners.

COPYRIGHT

The Software, including its structure, organization, code, user interface and associated Documentation, is a proprietary product of Molecular Devices or its suppliers, and is protected by international laws of copyright. The law provides for civil and criminal penalties for anyone in violation of the laws of copyright.

LICENSE

Use of the Software

1. If the Software is not instrument operating software, subject to the terms and conditions of this Agreement, Molecular Devices grants the purchaser of this product a non-exclusive license only to install and use the Software on a single computer (or in a single network, if your Software is designated as a network version). You may transfer the Software to another single computer (or network, if a network version), but the Software may never be installed on more than one computer (or more than one network, if a network version) at any one time.
2. If the Software is instrument operating software, you may install the Software on a single computer (or in a single network, if your Software is designated as a network version) and use the Software only for the operation of the instrument on which the Software is installed.
3. If the Software uses registration codes or product keys, access to the number of licensed copies of Software is controlled by a registration code or product key. For example, if you have a registration code or product key that enables you to use three copies of the Software simultaneously, you cannot install the Software on more than three separate computers.
4. You may make one copy of the Software in machine-readable form solely for backup or archival purposes. You must reproduce on any such copy all copyright notices and any other proprietary legends found on the original. You may not make any other copies of the Software.
5. From time to time, Molecular Devices may change the terms of this Agreement. Molecular Devices will notify you of such change. Your continued use of the Software will indicate your agreement to the change.

Restrictions

1. You may not use the Software, Documentation, or any associated documentation, in whole or in part, except as expressly permitted in this Agreement.
2. You may not sublicense, copy, merge, modify, sell, resell, rent, lease, assign, transfer (except temporarily in the event of a computer malfunction or as expressly permitted under Section 1 of "Use of the Software"), or distribute this license or the Software, or create derivative works based on the Software, or any part thereof or any interest therein.
3. You may not reverse assemble, decompile, modify, create any derivative works, or otherwise attempt, cause, or permit any reverse engineering of the Software or any part thereof.
4. You acknowledge that the Software in human-readable (source code) form, including its structure, sequence, and organization, is and remains the confidential trade secret of Molecular Devices. You agree that you shall not attempt, cause or permit the reverse engineering, modification, decompilation, disassembly of the Software to any human readable or perceivable form, or creation of any derivative works based on the Software or any part thereof.
5. You may not remove any proprietary, copyright, trade secret or warning legend from the Software or any Documentation and you acknowledge that the Software in human-readable (source code) form, including its structure, sequence, and organization, is and remains the confidential trade secret of Molecular Devices.
6. You agree to comply fully with all export laws and restrictions and regulations of the United States or applicable foreign agencies or authorities. You agree that you will not export or reexport, directly or indirectly, the Software into any country prohibited by the United States Export Administration Act and the regulations thereunder or other applicable United States law.

7. Oracle Corporation prohibits any transfer of Oracle Corporation products embedded in the Software. If this Software includes Oracle Corporation products, please contact Molecular Devices to obtain replacement operating software in connection with any sale of the instrument. A re-licensing fee may be charged for any such replacement operating software. (See "Third Party Products" above to determine if the Software includes Oracle Corporation products. If Oracle Corporation is not listed, the Software does not include Oracle Corporation products.)

Trial

If this license is granted on a trial basis, you are hereby notified that license management software may be included to automatically cause the Software to cease functioning at the end of the trial period.

Termination

You may terminate this Agreement by discontinuing use of the Software, removing all copies from your computers and storage media, and returning the Software and Documentation, and all copies thereof, to Molecular Devices. Molecular Devices may terminate this Agreement if you fail to comply with all of its terms, in which case you agree to discontinue using the Software, remove all copies from your computers and storage media, and return the Software and Documentation, and all copies thereof, to Molecular Devices.

U.S. Government End Users

The Software is a "commercial item," as that term is defined in 48 C.F.R. 2.101 (Oct. 1995), consisting of "commercial computer software" and "commercial computer software documentation," as such terms are used in 48 C.F.R. 12.212 (Sept. 1995). Consistent with 48 C.F.R. 12.212 and 48 C.F.R. 227.7202-1 through 227.7202-4 (June 1995), all U.S. Government End Users acquire the Software with only those rights set forth herein.

European Community End Users

If this Software is used within a country of the European Community, nothing in this Agreement shall be construed as restricting any rights available under the European Community Software Directive, O.J. Eur. Comm. (No. L. 122) 42 (1991).

Regulated Uses

You acknowledge that the Software has not been cleared, approved, registered or otherwise qualified (collectively, "Approval") by Molecular Devices with any regulatory agency, including without limitation for use in diagnostic or therapeutic procedures or any other use requiring compliance with any federal or state law regulating diagnostic or therapeutic products, blood products, medical devices or any similar product (hereafter collectively referred to as "federal or state drug laws"). The Software may not be used for any purpose that would require any such Approval unless proper Approval is obtained. You agree that if you elect to use the Software for a purpose that would subject you or the Software to the jurisdiction of any federal or state drug laws, you will be solely responsible for obtaining any required Approvals and otherwise ensuring that your use of the Software complies with such laws.

LIMITED WARRANTY and LIMITATION OF REMEDIES

Limited Warranty. Molecular Devices warrants that for a period of ninety (90) days from the beginning of the applicable warranty period (as described below), or for the designated warranty period if a different warranty period is designated as the warranty period for the Software in the current version of an instrument operating manual or catalog or in a specific written warranty including with and covering the Software, the Software will function substantially in accordance with the functions and features described in the Documentation delivered with the Software when properly installed, and that for a period of ninety days from the beginning of the applicable warranty period (as described below) the tapes, CDs, diskettes or other media bearing the Software will be free of defects in materials and workmanship under normal use.

The above warranties do not apply to defects resulting from misuse, neglect, or accident, including without limitation: operation outside of the environmental or use specifications, or not in conformance with the instructions for any instrument system, software, or accessories; improper or inadequate maintenance

by the user; installation of software or interfacing, or use in combination with software or products not supplied or authorized by Molecular Devices; and modification or repair of the products not authorized by Molecular Devices.

Warranty Period Commencement Date. The applicable warranty period for software begins on the earlier of the date of installation or three (3) months from the date of shipment for software installed by Molecular Devices personnel. For software installed by the purchaser or anyone other than Molecular Devices, the warranty period begins on the date the software is delivered to you. The applicable warranty period for media begins on the date the media is delivered to the purchaser.

MOLECULAR DEVICES MAKES NO OTHER WARRANTIES OF ANY KIND WHATSOEVER, EXPRESS OR IMPLIED, WITH RESPECT TO THE SOFTWARE OR DOCUMENTATION, INCLUDING BUT NOT LIMITED TO WARRANTIES OF FITNESS FOR A PARTICULAR PURPOSE OR MERCHANTABILITY OR THAT THE SOFTWARE OR DOCUMENTATION IS NON-INFRINGEMENT. ALL OTHER WARRANTIES ARE EXPRESSLY DISCLAIMED. WITHOUT LIMITING THE GENERALITY OF THE FOREGOING, MOLECULAR DEVICES MAKES NO WARRANTIES THAT THE SOFTWARE WILL MEET YOUR REQUIREMENTS, THAT OPERATION OF THE LICENSED SOFTWARE WILL BE UNINTERRUPTED OR ERROR FREE OR WILL CONFORM EXACTLY TO THE DOCUMENTATION, OR THAT MOLECULAR DEVICES WILL CORRECT ALL PROGRAM ERRORS. MOLECULAR DEVICES' SOLE LIABILITY AND RESPONSIBILITY FOR BREACH OF WARRANTY RELATING TO THE SOFTWARE OR DOCUMENTATION SHALL BE LIMITED, AT MOLECULAR DEVICES' SOLE OPTION, TO (1) CORRECTION OF ANY ERROR IDENTIFIED TO MOLECULAR DEVICES IN A WRITING FROM YOU IN A SUBSEQUENT RELEASE OF THE SOFTWARE, WHICH SHALL BE SUPPLIED TO YOU FREE OF CHARGE, (2) ACCEPTING A RETURN OF THE PRODUCT AND REFUNDING THE PURCHASE PRICE UPON RETURN OF THE PRODUCT AND REMOVAL OF ALL COPIES OF THE SOFTWARE FROM YOUR COMPUTERS AND STORAGE DEVICES, (3) REPLACEMENT OF THE DEFECTIVE SOFTWARE WITH A FUNCTIONALLY EQUIVALENT PROGRAM AT NO CHARGE TO YOU, OR (4) PROVIDING A REASONABLE WORK AROUND WITHIN A REASONABLE TIME. MOLECULAR DEVICES SOLE LIABILITY AND RESPONSIBILITY UNDER THIS AGREEMENT FOR BREACH OF WARRANTY RELATING TO MEDIA IS THE REPLACEMENT OF DEFECTIVE MEDIA RETURNED WITHIN 90 DAYS OF THE DELIVERY DATE. THESE ARE YOUR SOLE AND EXCLUSIVE REMEDIES FOR ANY BREACH OF WARRANTY. WARRANTY CLAIMS MUST BE MADE WITHIN THE APPLICABLE WARRANTY PERIOD.

LIMITATION OF LIABILITY

IN NO EVENT SHALL MOLECULAR DEVICES OR ITS AGENTS, CONSULTANTS, CONTRACTORS OR SUPPLIERS BE RESPONSIBLE OR LIABLE, WHETHER IN CONTRACT, TORT, WARRANTY OR UNDER ANY STATUTE (INCLUDING WITHOUT LIMITATION ANY TRADE PRACTICE, UNFAIR COMPETITION OR OTHER STATUTE OF SIMILAR IMPORT) OR ON ANY OTHER BASIS FOR SPECIAL, INDIRECT, INCIDENTAL, MULTIPLE, PUNITIVE, OR CONSEQUENTIAL DAMAGES ARISING OUT OF THE POSSESSION OR USE OF, OR THE INABILITY TO USE, THE SOFTWARE OR DOCUMENTATION, EVEN IF MOLECULAR DEVICES IS ADVISED IN ADVANCE OF THE POSSIBILITY OF SUCH DAMAGES, INCLUDING WITHOUT LIMITATION DAMAGES ARISING FROM OR RELATED TO LOSS OF USE, LOSS OF DATA, DOWNTIME, OR FOR LOSS OF REVENUE, PROFITS, GOODWILL OR BUSINESS OR OTHER FINANCIAL LOSS. IN ANY CASE, THE ENTIRE LIABILITY OF MOLECULAR DEVICES AND ITS AGENTS, CONSULTANTS, CONTRACTORS AND SUPPLIERS UNDER THIS LICENSE, OR ARISING OUT OF THE USE OF THE SOFTWARE, SHALL NOT EXCEED IN THE AGGREGATE THE PURCHASE PRICE OF THE LICENSE OR LICENSES TO THE SOFTWARE.

SOME STATES, COUNTRIES OR JURISDICTIONS LIMIT THE SCOPE OF OR PRECLUDE LIMITATIONS OR EXCLUSION OF REMEDIES OR DAMAGES, OR OF LIABILITY, SUCH AS LIABILITY FOR GROSS NEGLIGENCE OR WILLFUL MISCONDUCT, AS OR TO THE EXTENT SET FORTH ABOVE, OR DO NOT ALLOW IMPLIED WARRANTIES TO BE EXCLUDED. IN SUCH STATES, COUNTRIES OR JURISDICTIONS, THE LIMITATION OR EXCLUSION OF WARRANTIES, REMEDIES, DAMAGES OR LIABILITY SET FORTH ABOVE MAY NOT APPLY TO YOU. HOWEVER, ALTHOUGH

THEY SHALL NOT APPLY TO THE EXTENT PROHIBITED BY LAW, THEY SHALL APPLY TO THE FULLEST EXTENT PERMITTED BY LAW. YOU MAY ALSO HAVE OTHER RIGHTS THAT VARY BY STATE, COUNTRY OR OTHER JURISDICTION.

GENERAL

This Agreement shall be governed by laws of the Commonwealth of Massachusetts, U.S.A., exclusive of its conflict of laws provisions. This Agreement shall not be governed by the United Nations Convention on Contracts for the International Sale of Goods. This Agreement contains the complete agreement between the parties with respect to the subject matter hereof, and supersedes all prior or contemporaneous agreements or understandings, whether oral or written. If any provision of this Agreement is held by a court of competent jurisdiction to be contrary to law, that provision will be enforced to the maximum extent permissible, and the remaining provisions of this Agreement will remain in full force and effect. The controlling language of this Agreement, and any proceedings relating to this Agreement, shall be English. You agree to bear any and all costs of translation, if necessary. The headings to the sections of this Agreement are used for convenience only and shall have no substantive meaning. All questions concerning this Agreement shall be directed to: Molecular Devices, LLC c/o AB Sciex, 500 Old Connecticut Path, Framingham MA U.S.A., Attention: Legal Department.

Unpublished rights reserved under the copyright laws of the United States.
Molecular Devices, LLC, 1311 Orleans Drive, Sunnyvale CA, 94089 U.S.A.

For Research Use Only. Not for use in diagnostic procedures. The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. © 2016 Molecular Devices, LLC
V9/2016