



# GenePix Pro

Microarray Acquisition and Analysis Software

Software Release Notes

## GenePix Pro Microarray Acquisition and Analysis Software Release Notes

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

### **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. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/patents>

Product manufactured by Molecular Devices, LLC.  
3860 N. First Street, San Jose, California, 95134, United States of America.  
Molecular Devices, LLC is ISO 9001 registered.  
©2023 Molecular Devices, LLC.  
All rights reserved.



# Contents

- Chapter 1: GenePix Pro Software General Information ..... 5**
  - Introduction ..... 5
  - Computer System Requirements ..... 5
  - Software Installation ..... 5
  - License Key Installation ..... 6
  - Troubleshooting the Installation ..... 7
  - Obtaining Support ..... 7
- Chapter 2: Version 7.4 Release Notes ..... 9**
  - New in Version 7.4 ..... 9
- Chapter 3: Version 7.3.1 Release Notes ..... 11**
  - Issues Addressed in Version 7.3.1 ..... 11
- Chapter 4: Version 7.3 Release Notes ..... 13**
  - New in Version 7.3 ..... 13



# Chapter 1: GenePix Pro Software General Information

# 1

## Introduction

The GenePix® Pro Microarray Acquisition and Analysis Software version 7.4 is a fully featured array acquisition and first-pass analysis solution, providing advanced imaging and data manipulation capabilities.

## Computer System Requirements

GenePix Pro Microarray Acquisition and Analysis Software can be installed on a computer with the following minimum system specifications:

### Supported Operating Systems

- Windows 7, 32-bit and 64-bit
- Windows 10, 32-bit and 64-bit
- Windows 11, 64-bit

### Hardware

- 1.2 GHz or faster processor
- 4 GB RAM or more
- 250 GB hard drive (for image storage)
- USB 2.0 ports for:
  - USB software protection key
  - Scanner
  - SL50 slide loader (Discontinued)

## Software Installation

Use the following procedure to install the GenePix Pro Microarray Acquisition and Analysis Software.

1. Double-click the **GenePixPro\_7\_4.exe** file to start the installation. Install Wizard dialog is displayed.
2. Click **Next**.
3. In the **License Agreement** dialog, select **I accept the terms of the license agreement**, and click **Next**.
4. In the **Destination Folder** dialog, the **Install GenePix Pro 7 to** field displays the default installation directory. To change the installation directory, click **Change**, navigate to the desired directory, and then click **OK**.
5. Click **Next**.
6. In the **Select Program Folder** dialog, leave the displayed default **Program Folder** settings. Select **Anyone who uses this computer**, then click **Next**.
7. In the **Simulation Images Setup** dialog, click **OK**.
8. In the **Software License Setup** dialog, click **OK**.
9. In the **Sentinel Runtime Setup** dialog, click **Next**.
10. When the installation is complete, click **Finish** to exit the wizard.

## License Key Installation

GenePix Pro Software requires a license key (sometimes called a dongle) to operate fully. Without the key, you can open images, but you cannot save settings files or analysis results, and you can analyze only up to 1000 features.

The license key is a USB device that provides security authorization for a single-user or for multiple users across a network, which avoids the inconvenience of installing individual devices for each computer on your network. The license keys can also be used across a network for remote access. Whether you have a single-user or multi-user license, your key is programmed with the number of licenses purchased.

### Single-User License Keys

For single-user licenses, a single physical license key device is provided. To install the single-user license key, insert it into a USB port on your computer.

### Multi-User Network License Keys

If you have a multi-user license, a network USB license key and a HASP Admin Control Center software package is provided. To take advantage of multiple-site licenses, use the USB license key on a local area network. You must choose a single computer on your network to act as your GenePix Pro system server. On this computer, install the physical network key, the HASP Admin Control Center software, and the related device drivers.

### Installing the License Key on a Network

If you have purchased a multi-site license for your software, you will be provided with a USB license key that you need to install to access the full functionality of the software. The USB license key can be attached to any computer on the network that has an available USB port, with the following condition:

- The chosen computer should be one that is always switched on and rarely rebooted, such as a mail, print, or other software server.

The GenePix Pro Software license key can be used in conjunction with other HASP USB license keys; in particular, if you use two or more Molecular Devices software products on your network, each with a multi-site license and USB license key, you can install all the network keys on the same computer, as long as they are all HASP Version 3.0 or greater. To install the network key, attach it to an available USB port on your chosen computer.

To see your HASP version, connect the key to your computer and open the HASP Admin Control Center. In the HASP Keys Options section, the connected keys are listed together with their properties and version number.



**Note:** If you have a previous version of the GenePix Pro Software, you might be using HASP 2.x version keys. These require the HASP License Manager for network configuration. You can install and run this utility on the same computer as your HASP Admin Control Center. See your original GenePix Pro user guide for installation instructions.

---

## HASP Admin Control Center

To provide centralized control and reporting of license and user information within your network, the GenePix Pro software package automatically installs the HASP Admin Control Center and HASP protection key drivers. The HASP Admin Control Center runs on your Internet browser. It serves two purposes:

- For license servers, to manage licenses of any attached USB License keys.
- For client computers, to view available USB License keys and licenses on the network.

After you install the GenePix Pro Software and attach the license key to your computer, you can view key properties and manage licenses in the HASP Admin Control Center.

You can access the HASP Admin Control Center from any computer on your network where the GenePix Pro Software has been installed. Type **HTTP://LOCALHOST:1947/** in the address bar of your Internet browser. You can see all the GenePix Pro Software license keys installed on your local network.

If your multi-site license software came with a bundled computer package, that computer already has the GenePix Pro Software and the HASP Admin Control Center installed. To use a different computer as the license server, install the GenePix Pro Software on that computer to get the HASP Admin Control Center. To install the HASP Admin Control Center alone, download the HASP SRM Runtime Setup files from <http://www.Aladdin.com/support/hasp-srm/enduser.aspx>.

## Troubleshooting the Installation

If you experience difficulty running the GenePix Pro Software, verify the following:

- Your computer has the software required minimum 4 GB of RAM to acquire images.
- The license key is installed. The software requires a license key to save results and settings files, and to analyze images. Without the key, the software runs with limited analysis and file-saving functionality.

## Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest level of technical service.

Our Support website—[support.moleculardevices.com/](http://support.moleculardevices.com/)—describes the support options offered by Molecular Devices, including service plans and professional services. It also has a link to the Molecular Devices Knowledge Base, which contains documentation, technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance, you can submit a request to Molecular Devices Technical Support.

Please have your instrument serial number or Work Order number and your software version number available when you call.





## Chapter 2: Version 7.4 Release Notes



# 2

The GenePix Pro Microarray Acquisition and Analysis Software version 7.4 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 7.3.1, the last general release of the GenePix Pro Microarray Acquisition and Analysis Software.

### New in Version 7.4

The following new features are included in GenePix Pro Microarray Acquisition and Analysis Software version 7.4.

#### **Windows 11 Operating System Support**

Adds software support for the Windows 11 operating system.



## Chapter 3: Version 7.3.1 Release Notes



# 3

The GenePix Pro Microarray Acquisition and Analysis Software version 7.3.1 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 7.3, the last general release of the GenePix Pro Microarray Acquisition and Analysis Software.

### Issues Addressed in Version 7.3.1

The following issue was addressed in GenePix Pro Microarray Acquisition and Analysis Software version 7.3.1.

#### **Need Support for the GenePix 4000B Instrument on Windows 10**

Tracking ID: 5268

GenePix Pro software is missing the Windows signed driver for the GenePix 4000B instrument.

#### **Resolution:**

This issue has been fixed in the software.

#### **Impact of fix:**

This fix has no impact on current workflow or data.



## Chapter 4: Version 7.3 Release Notes



# 4

The GenePix Pro Microarray Acquisition and Analysis Software version 7.3 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 7.2, the last general release of the GenePix Pro Microarray Acquisition and Analysis Software.

### New in Version 7.3

The following new features are included in GenePix Pro Microarray Acquisition and Analysis Software version 7.3.

#### Windows 10 Operating System Support

Adds software support for the Windows 10 (32-bit and 64-bit) operating system.

## Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)  
Web: [moleculardevices.com](http://moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)

Visit our website for a current listing of worldwide distributors.