



Agilent CytoGenomics 5.0

Product Overview and Installation Guide

**For Research Use Only. Not for use in diagnostic
procedures.**



Agilent Technologies

Notices

© Agilent Technologies, Inc. 2018

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.

Manual Part Number

G1662-90057

Edition

Revision C0, December 2018

Agilent Technologies, Inc.
5301 Stevens Creek Blvd.
Santa Clara, CA 95051

Technical Support

For US and Canada

Call 800-227-9770 (option 3, 4, 4)
Or send an email to
informatics_support@agilent.com

For Europe, Middle East, Africa, and India

Call 00800 345 600 (toll free) or
+49 69 8679 7730
Or send an email to
genomics_tech_europe@agilent.com

For all other regions

Agilent's world-wide sales and support center telephone numbers can be obtained at www.agilent.com/genomics under Contact Us.
Or send an email to genomics@agilent.com.

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.

In This Guide...

This *Product Overview and Installation Guide* provides an overview of Agilent CytoGenomics 5.0 and instructions for installing and administering the software.

1 Product Overview

This chapter provides an overview of Agilent CytoGenomics. It explains the overall design of the program and how different users can work with it to fulfill their roles in the laboratory.

2 Installation Instructions for Windows

This chapter shows you how to install Agilent CytoGenomics 5.0 and migrate data from a previous version of CytoGenomics on computers running a compatible Microsoft Windows operating system.

3 Installation Instructions for Macintosh

This chapter shows you how to install Agilent CytoGenomics 5.0 and migrate data from a previous version of CytoGenomics on systems running a Macintosh OS.

4 Administering the System

This chapter shows you how to add users, assign roles and manage the database.

Contents

1 Product Overview 9

What is Agilent CytoGenomics? 10

Getting Help 12

To get help within Agilent CytoGenomics 5.0 12

To contact Agilent Technical Support 12

To learn about Agilent Genomics products and services 12

Key Features 13

The Home tab - getting started with CytoGenomics 13

The Analysis Workflow tab – selecting samples and running analyses 13

The Sample Review tab – reviewing analyses 14

The Configure Settings tab – setting up system defaults and analysis methods 14

The Auto Processing tab – setting up Auto-Processing workflows 15

The Supporting Files tab – managing samples, designs, gene lists, tracks, and other supporting files 15

The Admin tab – managing users and database 15

User Roles and Capabilities 16

2 Installation Instructions for Windows 17

Installing Agilent CytoGenomics 5.0 – Windows 18

Step 1. Prepare for installation 18

Step 2. Install the database server and client on a single Windows computer 20

Step 3. Install the client software on additional Windows computers 36

Step 4. Start the Agilent CytoGenomics 5.0 client software 39

Step 5. Log in to Agilent CytoGenomics 39

Step 6. Add your software license 41

Step 7. Add users and assign roles 43

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows 44

Step 1. Install CytoGenomics 5.0 on the server computer	44
Step 2. Migrate data to CytoGenomics 5.0 server	47
Step 3. Upgrade to CytoGenomics 5.0 on the client computers	52
Step 4. Start the Agilent CytoGenomics 5.0 client software, log in, and add new software license	54
Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2	55
Step 1. Install CytoGenomics 5.0.2 client and uninstall CytoGenomics 5.0.0 or 5.0.1 client	55
Step 2. Start the Agilent CytoGenomics 5.0.2 client software, log in, and add new software license	60
Uninstalling Agilent CytoGenomics on Windows computers	61
Troubleshooting	63
If the Agilent CytoGenomics program does not open	63
3 Installation Instructions for Macintosh	65
Installing Agilent CytoGenomics 5.0 – Macintosh	66
Step 1. Prepare for installation	66
Step 2. Install the database server and client on a single Macintosh computer	68
Step 3. Install the client software on additional Macintosh computers	84
Step 4. Start the Agilent CytoGenomics 5.0 client software	87
Step 5. Log in to Agilent CytoGenomics	87
Step 6. Add your software license	89
Step 7. Add users and assign roles	90
Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh	91
Step 1. Install CytoGenomics 5.0 on the server computer	91
Step 2. Migrate data to CytoGenomics 5.0 server	94
Step 3. Upgrade to CytoGenomics 5.0 on the client computers	98
Step 4. Start the Agilent CytoGenomics 5.0 client software, log in, and add new software license	98
Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2	99

Step 1. Uninstall CytoGenomics 5.0.0 or 5.0.1 client	99
Step 2. Install CytoGenomics 5.0.2 client	102
Step 3. Start the Agilent CytoGenomics 5.0.2 client software, log in, and add new software license	104
Uninstalling Agilent CytoGenomics on Macintosh computers	105
4 Administering the System	109
Administering the System	110
To manage users and roles	110
To manage the database	111



1

Product Overview

What is Agilent CytoGenomics? [10](#)

Getting Help [12](#)

To get help within Agilent CytoGenomics 5.0 [12](#)

To contact Agilent Technical Support [12](#)

To learn about Agilent Genomics products and services [12](#)

Key Features [13](#)

The Analysis Workflow tab – selecting samples and running analyses [13](#)

The Sample Review tab – reviewing analyses [14](#)

The Configure Settings tab – setting up system defaults and analysis methods [14](#)

The Supporting Files tab – managing samples, designs, gene lists, tracks, and other supporting files [15](#)

The Admin tab – managing users and database [15](#)

User Roles and Capabilities [16](#)

This chapter provides an overview of Agilent CytoGenomics. It explains the overall design of the program and how different users can work with it to fulfill their roles in the laboratory.



Agilent Technologies

What is Agilent CytoGenomics?

Cytogenetic research labs are increasingly turning to array CGH to improve detection of copy number changes and increase sample throughput. Alongside Agilent's gold standard CGH+SNP and CGH microarrays, Agilent CytoGenomics software completes Agilent's array-based cytogenetic research solution. Agilent CytoGenomics provides a statistically and visually powerful tool to streamline the day-to-day cytogenetic sample analysis research workflow. It offers full support for efficient analysis of cytogenetic samples and easy report generation, based on information from both external databases and an internal database. Key automation features enable a seamless flow through the experimental process, starting with full automation of data upload and analysis, and ending with customized report generation.

With Agilent CytoGenomics, cytogenetic research labs can analyze data from a microarray image in 5–30 minutes. This increases efficiency and throughput, while enabling cytogenetic researchers to better detect copy number and copy-neutral aberrations across the genome.

Agilent CytoGenomics supports CGH analysis for both constitutional and tumor samples from either fresh or FFPE sample sources. The SNP-based LOH (Loss or Lack of Heterozygosity) and UPD (Uniparental Disomy) analysis is supported for both constitutional and liquid tumor samples from fresh sample sources only.

Agilent CytoGenomics provides cytogenetic laboratories security data access through three types of user roles: Technician, Scientist, and Administrator. Each of the three user roles has unique permissions to access a selected set of features and data in the software. (See [Table 1.](#)) In addition to these three default user roles, Administrators can create custom user roles with a defined set of user privileges.

An Agilent CytoGenomics workflow ([Figure 1](#)) automatically feature extracts images generated by a microarray scanner. The extracted data is then filtered and analyzed, and aberrant intervals are identified. You can edit, suppress, add, or classify identified aberration calls during sample triage. A set of reports is generated at the end of the workflow, and a Cyto report is generated upon sign-off of the record. Audit trails of changes are available for display or inclusion in the Cyto report.

* The exact time required to process a workflow job is dependent on the format of the microarray and the specifications of the computer running the software.

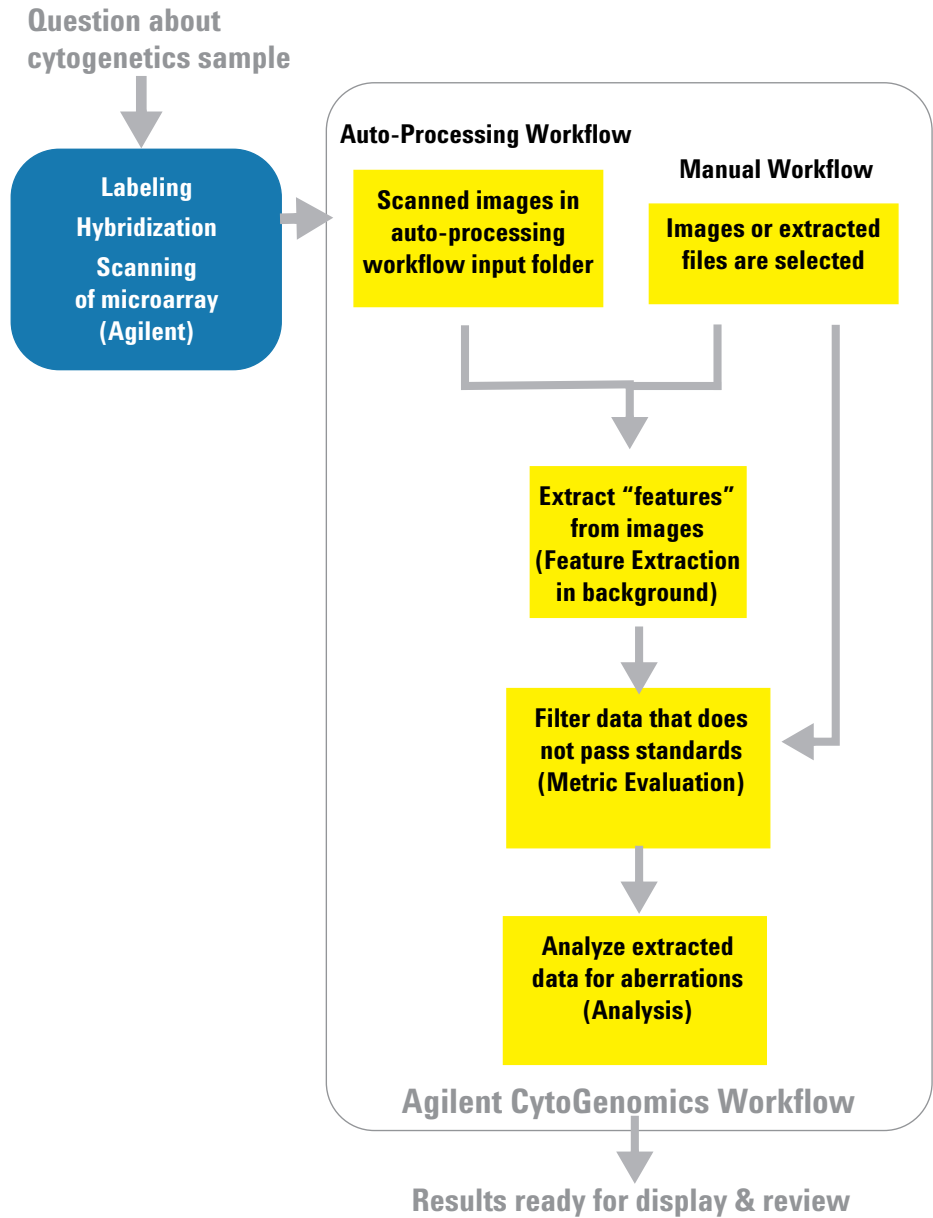


Figure 1 Agilent CytoGenomics Workflow — from image to result

Getting Help

To get help within Agilent CytoGenomics 5.0

Agilent CytoGenomics 5.0 has an online help system that provides detailed instructions on using the software and descriptions of the user interface. To open the help, click **Help** near the top right corner of the program window.

Help videos are also available from within the Agilent CytoGenomics program. These short videos give you instructions for doing basic tasks within the program. You can access the videos from the CytoGenomics Home screen.

To contact Agilent Technical Support

Technical support is available by phone and/or e-mail.

Resource	To find technical support contact information
Contact Agilent Technical Support by telephone or e-mail message (United States and Canada)	Telephone: 800-227-9770 (option 3, 4, 4) E-mail message: informatics_support@agilent.com
Contact Agilent Technical Support by telephone or e-mail message (for your country)	1 Go to www.agilent.com/en/contact-us/page. 2 Under Worldwide click to select a country or region. Complete e-mail message and telephone contact information for your country is displayed.

To learn about Agilent Genomics products and services

To view information about the Genomics products and services that are available from Agilent, go to www.agilent.com.

Key Features

This section provides an overview of the features and capabilities available in Agilent CytoGenomics 5.0. The “tabs” at the top of the program window provide access to related tasks and commands. Users see only the tabs for which they have access, according to their assigned user role. (See “[User Roles and Capabilities](#)” on page 16.)

The Home tab - getting started with CytoGenomics

Capabilities within the Home tab include:

- Entering your user license for the software
- Watching help videos for CytoGenomics
- Viewing news and updates from Agilent
- Opening links to other Agilent Genomics software programs
- Opening links to Agilent social media sites

The Analysis Workflow tab – selecting samples and running analyses

Capabilities within the Analysis Workflow tab include:

- Running manual workflows
- Adding sample information
- Importing design files
- Viewing array images by launching the Feature Extraction program

The Sample Review tab – reviewing analyses

Capabilities within the Sample Review tab include:

- Monitoring and searching for workflow jobs
- Displaying Cyto reports and aberration results
- Selecting sample results for triaging tasks:
 - Check in/out samples
 - Compare results from multiple samples
 - Add sample and interval notes
 - Modify, add, or suppress calls
 - Add/change interval classifications
 - Search database for similar classified intervals
 - Sign off samples and generate sign off reports
- Reanalyzing samples
- Reviewing QC metrics and SNP fit, and changing copy number peak assignments
- Transferring results to Cartagenia BENCH

The Configure Settings tab – setting up system defaults and analysis methods

Capabilities within the Configure Settings tab include:

- Setting up analysis method parameters
- Creating customized QC metric sets
- Creating customized Cyto report templates
- Viewing, creating, and updating dynamic and static tracks
- Configuring system preferences, including:
 - Default view preferences
 - Default input/output folders
 - Default genome build
 - Standard notes and classifications
 - Cartagenia BENCH data transfer settings
- Entering Cartagenia BENCH and Agilent SureDesign accounts

The Auto Processing tab – setting up Auto-Processing workflows

Capabilities within the Auto Processing tab include:

- Designating input folders and archive folders for tiff images and sample attribute files
- Setting up criteria for stopping or failing Auto-Processing
- Viewing Auto-Processing workflow logs

Note that Auto-Processing is not available in the Macintosh version of CytoGenomics.

The Supporting Files tab – managing samples, designs, gene lists, tracks, and other supporting files

Capabilities within the Supporting Files tab include:

- Viewing sample attributes and data
- Downloading new builds of a design from eArray
- Exporting sample attribute files
- Configuring spike-in identifiers
- Importing design files
- Importing genotype reference files
- Importing and exporting probe ID lists
- Importing and exporting tracks
- Opening the workflow log
- Displaying list of samples in the database, organized by design/build

The Admin tab – managing users and database

Capabilities within the Admin tab include:

- Adding new users and creating user roles
- Changing roles for existing users
- Enabling/disabling user access to the program
- Displaying or changing database connection parameters for the client computer
- Backing up the CytoGenomics database
- Clearing cached data

User Roles and Capabilities

The following table lists the capabilities allowed for the three different default user roles within Agilent CytoGenomics 5.0.

Table 1 User roles and capabilities

Role	Tasks and capabilities
Technician	• Run manual or auto-processing workflows • Add sample information • Monitor workflow jobs • Display reports and aberration results • View sample results in Triage View • Display auto-processing logs (Technicians cannot sign off samples.)
Scientist	Technician tasks, plus: • Configure system preferences, including • Default input/output folders • Auto-processing workflow settings • Default view preferences • Default genome build • Standard notes, classifications, and queries • Dynamic and static tracks • Cartagenia BENCH data transfer settings • Set analysis and workflow settings • Create standard notes, classifications, and queries • Create customized Cyto report templates • Manage samples • Evaluate and chart QC trends • Sign off and unlock results
Administrator	Complete system access, including all Technician and Scientist tasks, plus • Add users and create user roles • Change database connection settings for client systems • Backup and restore database



2

Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows 18

Step 1. Prepare for installation 18

Step 2. Install the database server and client on a single Windows computer 20

Step 3. Install the client software on additional Windows computers 36

Step 4. Start the Agilent CytoGenomics 5.0 client software 39

Step 5. Log in to Agilent CytoGenomics 39

Step 6. Add your software license 41

Step 7. Add users and assign roles 43

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows 44

Step 1. Install CytoGenomics 5.0 on the server computer 44

Step 2. Migrate data to CytoGenomics 5.0 server 47

Step 3. Upgrade to CytoGenomics 5.0 on the client computers 52

Step 4. Start the Agilent CytoGenomics 5.0 client software, log in, and add new software license 54

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2 55

Uninstalling Agilent CytoGenomics on Windows computers 61

Troubleshooting 63

This chapter shows you how to install Agilent CytoGenomics 5.0 and migrate data from a previous version of CytoGenomics on computers running a compatible Microsoft Windows operating system.



Agilent Technologies

Installing Agilent CytoGenomics 5.0 – Windows

This section includes detailed instructions and important notes for installing Agilent CytoGenomics 5.0 on a Windows system that does not have a previous version of CytoGenomics installed.

Agilent CytoGenomics consists of two parts: a PostgreSQL database server and the client software. Once you have installed the database server and client software on one computer, you can install additional copies of the client software on client computers that have network access to the database computer.

For instructions on how to upgrade an existing Agilent CytoGenomics installation, see “[Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows](#)” on page 44.

Step 1. Prepare for installation

a. Download the Windows version of the software and release notes

- 1 Go to the CytoGenomics download page on the Agilent website at: www.agilent.com/en/download-agilent-cyto-genomics-software.
- 2 Click the **Download** link for the Windows version of the software.
- 3 Save the Installer on your computer. (Do not start the installation at this time.)
- 4 Click the **Release Notes** link to download the software release notes.
- 5 Save the release notes to your computer.

b. Check system requirements before you start

Check to make sure that your computer systems meet the following requirements. All client installations point to the computer on which the server software is installed. Requirements are listed in [Table 2](#).

NOTE

If an operating system is not listed in [Table 2](#), it is not supported.

Table 2 Requirements for Agilent CytoGenomics computers – Windows

Operating system	64-bit Windows 7 Enterprise, or 64-bit Windows 10 Enterprise or Professional, or Windows Server 2012
	Note: For all of these operating systems, the regional setting must be set to English
Programs	Any program that enables you to open PDF files (for example, Adobe Reader)
Processor	Minimum: > 2 GHz Recommended: > 3 GHz
Working memory (RAM)	Minimum: 4 GB Recommended: 8 GB
Hard disk space	500 GB
Display Resolution	1280 x 768 or higher

c. Set up permissions for Installer and users

- Make sure that the user account used to install the Agilent CytoGenomics 5.0 software has administrator permissions on the installation computer.
- Make sure that every user account running the Agilent CytoGenomics 5.0 software has read/write permissions in the computer where the client software is installed. In addition, users must have read/write permissions in the installation folder of the computer where the database server is installed.

The way you set permissions varies between operating systems. Contact your network administrator for help in checking or changing user account permissions. If the user account permissions are not set correctly, the software will fail to install or run properly.

d. Adjust User Account Control settings for Windows 7 and Windows 10

If you are using the Windows 7 or Windows 10 operating system, change the User Account Control settings from the default setting to “Never Notify.”

- 1** Go to Control Panel.
- 2** Click **System and Security**.

- 3 Click **Change User Account Control setting** (found under **Action Center** for Windows 7 and under **Security and Maintenance** for Windows 10).
- 4 Change the setting to “Never Notify” and click **OK**.
- 5 Restart the machine.

Step 2. Install the database server and client on a single Windows computer

NOTE

Whether you plan on having a centralized, shared database, or plan on having one standalone installation, you must first install both the CytoGenomics client and server software on a machine with adequate disk space. (Available and required disk space are provided during installation at the Pre-Installation Summary screen. See [Figure 14](#) on page 32.)

The server is the program that hosts the CytoGenomics database. The client is the program you run when you want to use CytoGenomics. You can run the client program on the same computer as the server, or from any other computer with the client program installed. After installation, the database server software runs silently in the background and manages the shared database for all Agilent CytoGenomics 5.0 users. All client computers must have network access to the server computer and users must have read/write permissions.

- 1 Log on to the Windows system where you want to install Agilent CytoGenomics 5.0 server. Use an account that has administrator permissions.
- 2 Locate the installation program that you downloaded in Step 1, “a. [Download the Windows version of the software and release notes](#)” on page 18. Double-click the Agilent CytoGenomics 5.0.X.X application file to start the installation wizard.

The following message is displayed.

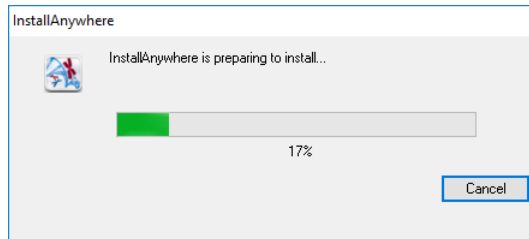


Figure 2 InstallAnywhere dialog box - preparing to install

The installation wizard opens to the Introduction screen.

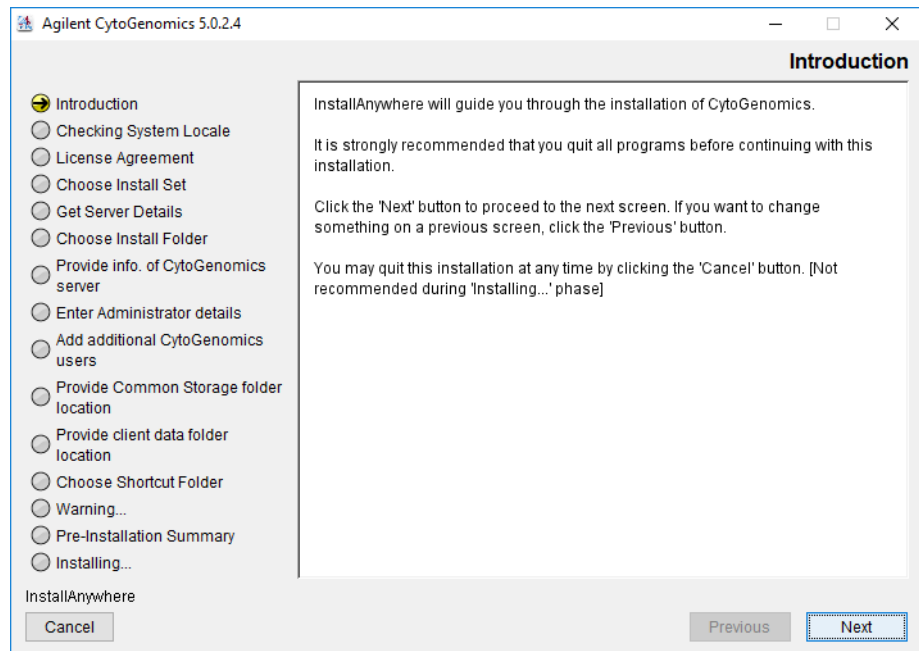


Figure 3 Introduction screen for the installation wizard – Windows

3 Read the introductory information.

4 Click **Next**.

The License Agreement screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

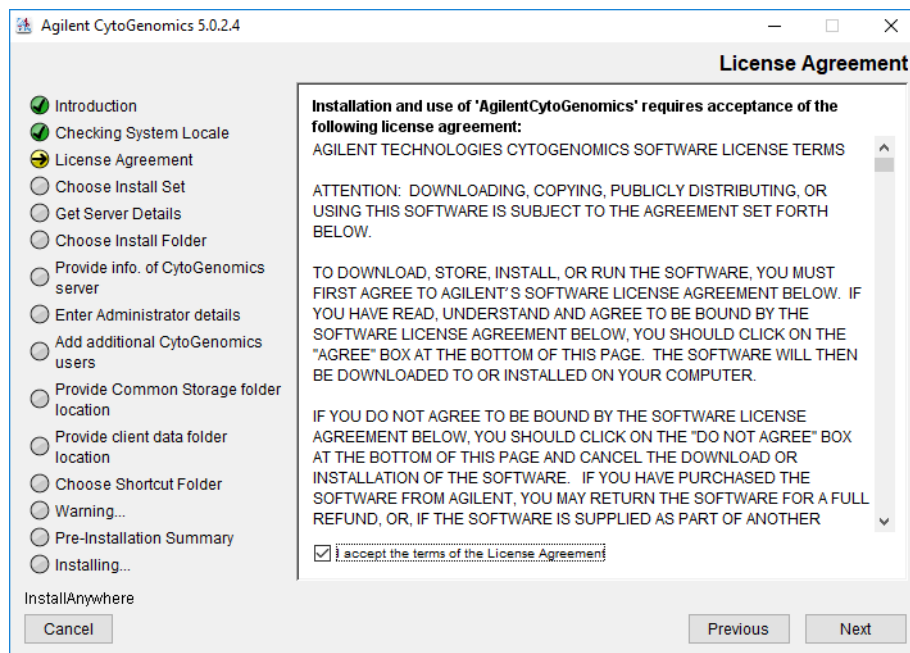


Figure 4 License Agreement screen – Windows

- 5 Read the license agreement. Click **I accept the terms of the License Agreement** and then click **Next**. (Click **Previous** to review information in a previous section.)

The Choose Install Set screen opens.

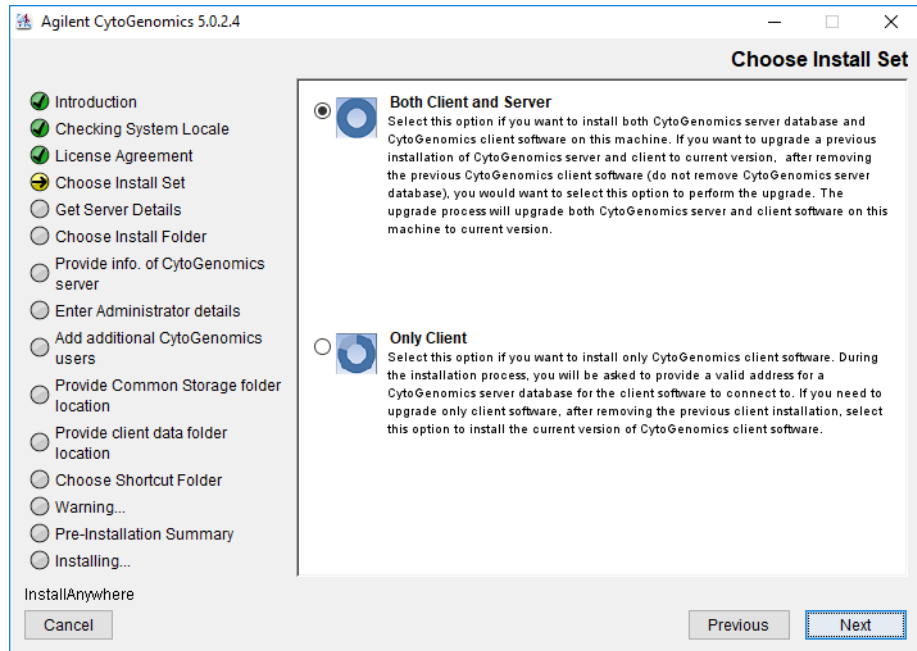


Figure 5 Choose Install Set screen – Windows

- 6** Click the icon next to **Both Client and Server** to install both the database server and client software on this computer.
- 7** Click **Next**.

An Information message opens. Click **OK** to continue. The Choose Install Folder screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

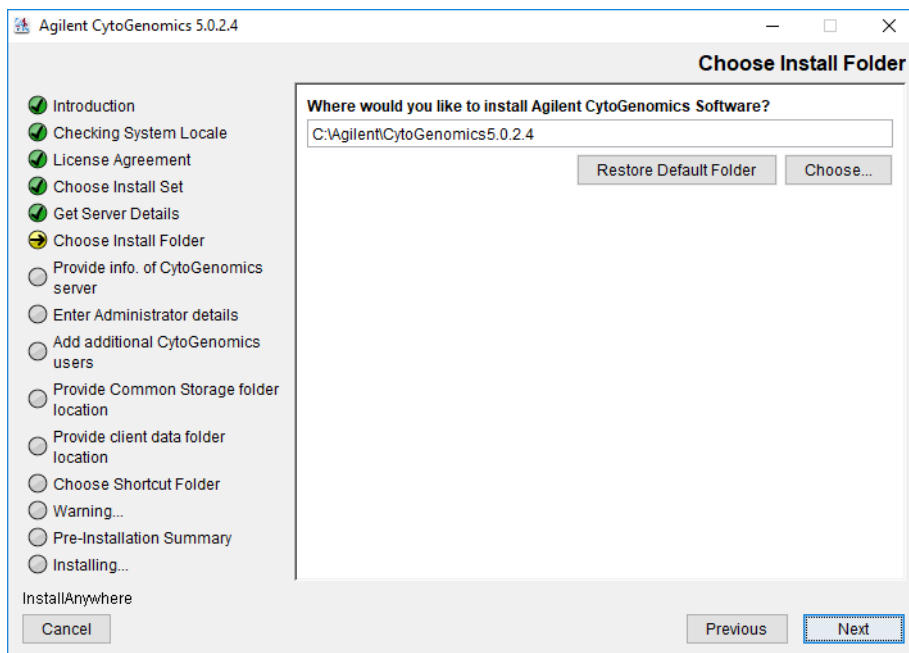


Figure 6 Choose Install Folder screen – Windows

- 8 By default, the program is installed in:
C:\Agilent\AgilentCytoGenomics5.0.X.X (where X.X is the revision number). Leave this folder as-is, or type the location and name of the folder where you want to install Agilent CytoGenomics 5.0. Within this folder, one folder is created for the server software and one folder is created for the client software.

OR

Click **Choose**, browse to a location where you want to install the program, and then click **Save**.

NOTE

For installations on Windows 7, the installation program folder must not be in the C:\Program Files folder.

- 9 Click **Next**.

The Enter Administrator details screen opens.

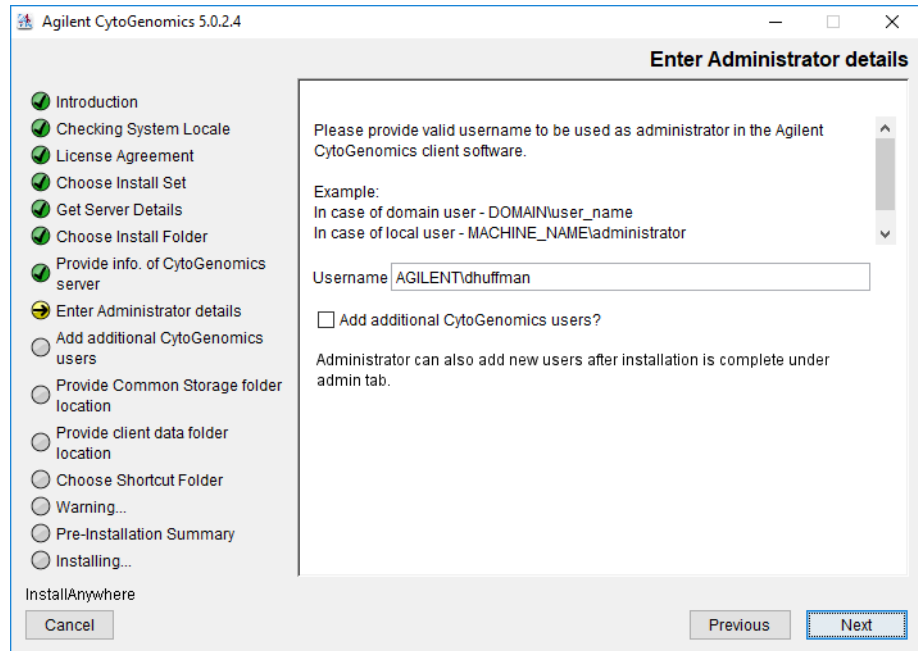


Figure 7 Enter Administrator details screen – Windows

- 10** The domain and user name of the person currently logged on to the computer appear in Username. If this Username is not the administrator for the Agilent CytoGenomics 5.0 software, type the correct user, using the format example shown.

For domain user: DOMAIN\username

For local machine: MACHINENAME\username

- 11** To set up additional CytoGenomics users, mark the check box labeled **Add additional CytoGenomics users?**. In the Password field that appears, type the password for the administrator account.

If you do not want to set up new users at this time, proceed directly to [step 16 on page 27](#). Note that the administrator can set up new users from within CytoGenomics once installation is complete (see “[To manage users and roles](#)” on page 110).

- 12** Click **Next**.

The Add additional CytoGenomics users screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

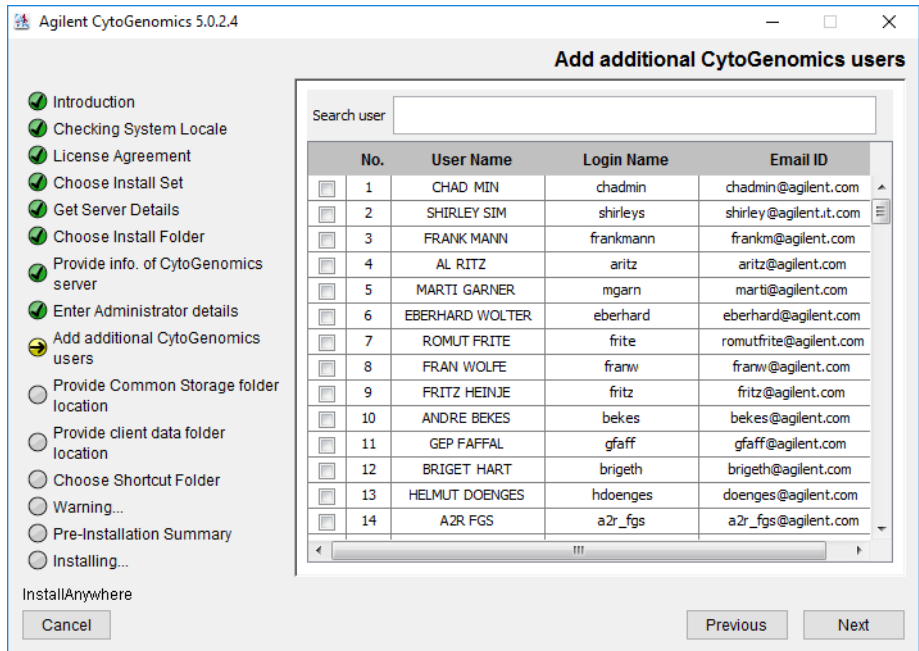


Figure 8 Add additional CytoGenomics users screen – Windows

13 The Add additional CytoGenomics users screen lists all of the users in the domain or on the local machine. Mark the check box next to the users that you want to add to the CytoGenomics database.

To search for a user, type the user name into the search box at the top of the screen.

14 Click **Next**.

The Add additional CytoGenomics users screen advances to the step for configuring user roles.

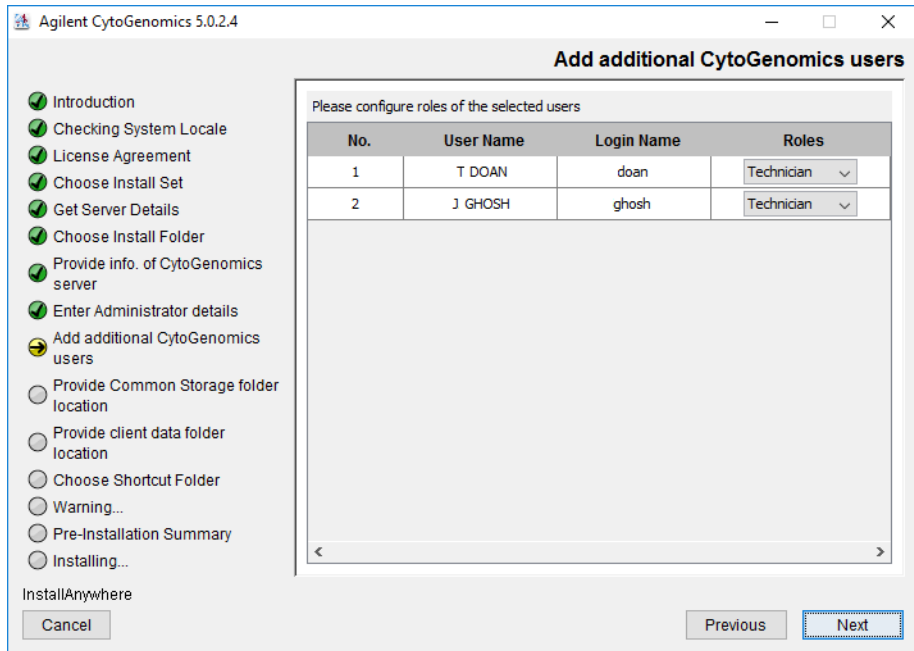


Figure 9 Add additional CytoGenomics users screen – Configure roles step – Windows

15 For each user listed on the screen, select a role in the Roles drop-down list. The default selection is *Technician*.

See “[User Roles and Capabilities](#)” on page 16 for a description of each role.

16 Click **Next**.

The Choose common storage folder screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

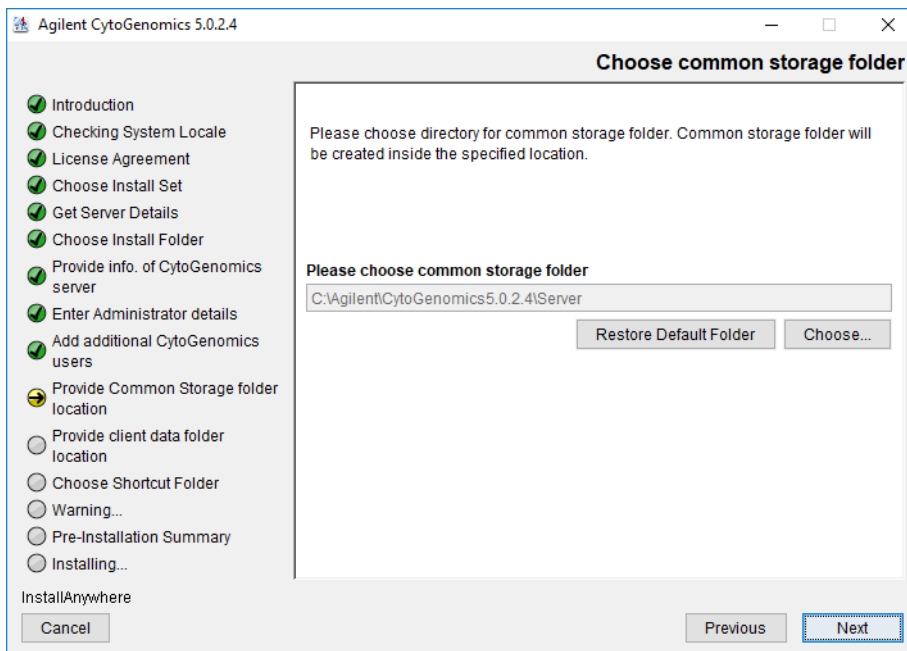


Figure 10 Choose common storage folder screen – Windows

17 By default, the common storage folder is created inside the folder C:\Agilent\AgilentCytoGenomics5.0.X.X\Server (where X.X is the revision number). Leave this folder as-is, or Click **Choose**, browse to the desired folder, and then click **Save**.

The common storage folder is where the program saves files that are used at the server level.

18 Click **Next**.

The Choose client data folder screen opens.

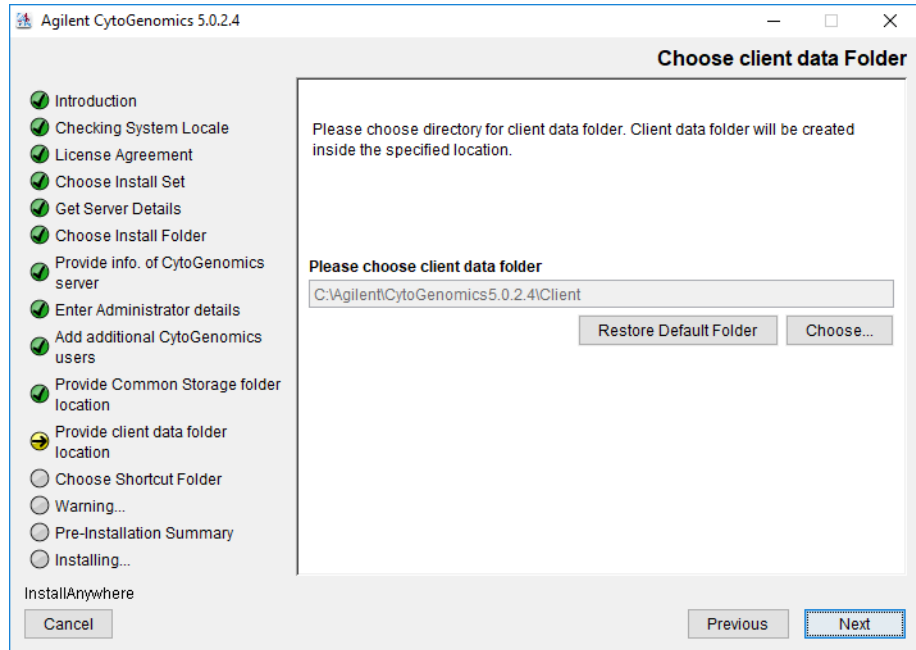


Figure 11 Choose client data folder screen – Windows

19 By default, the client data folder is created inside the folder C:\Agilent\AgilentCytoGenomics5.0.X.X\Client (where X.X is the revision number). Leave this folder as-is, or Click **Choose**, browse to the desired folder, and then click **Save**.

The client data folder is where the program installs the files for the genome build and reference genomes.

20 Click **Next**.

The Choose Shortcut Folder screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

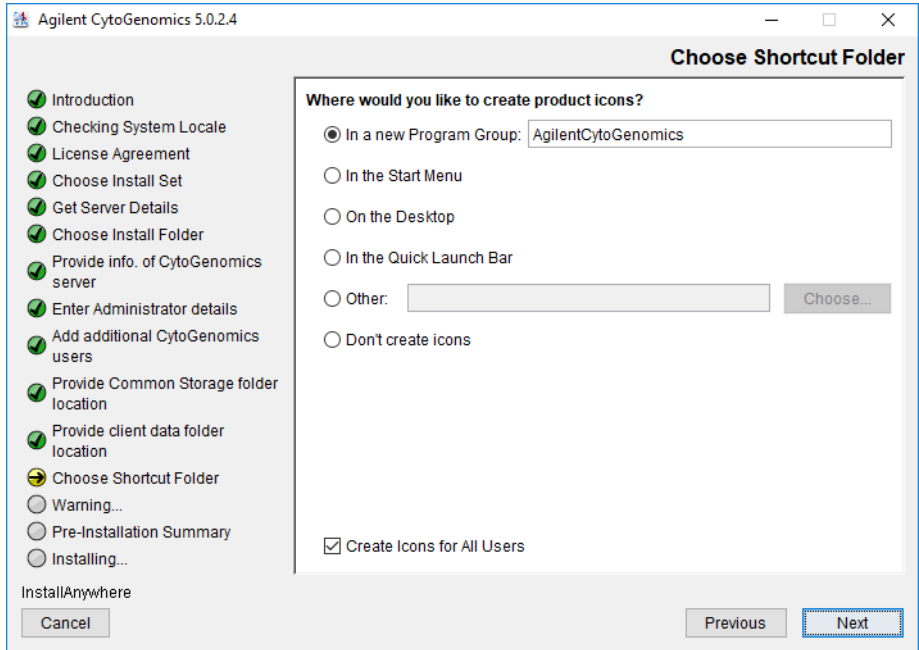


Figure 12 Choose Shortcut Folder screen – Windows

21 By default, program icons are installed in the Agilent CytoGenomics 5.0.X.X program group on your computer (where X.X is the revision number). To select another location for the program icons, select the desired location (on the desktop, for example) and then click **Next**.

NOTE

If more than one user plans to use Agilent CytoGenomics 5.0 on this computer, select **Create Icons for All Users**. Otherwise, other users do not see the program icons.

The Warning screen opens.

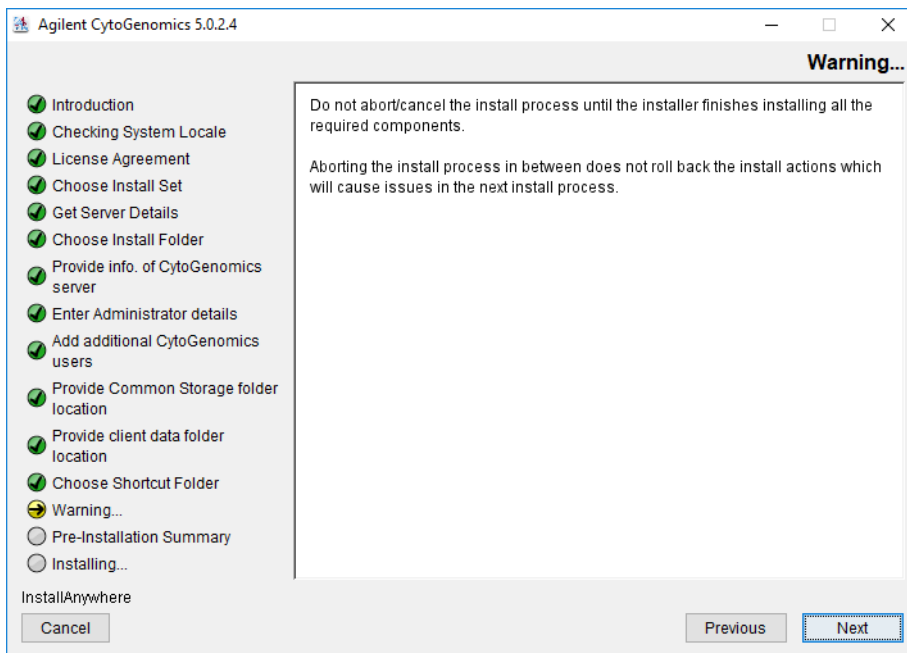


Figure 13 Warning screen – Windows

22 Read the message that cautions you not to stop the installation once it has started, and then click **Next**.

The Pre-Installation Summary screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

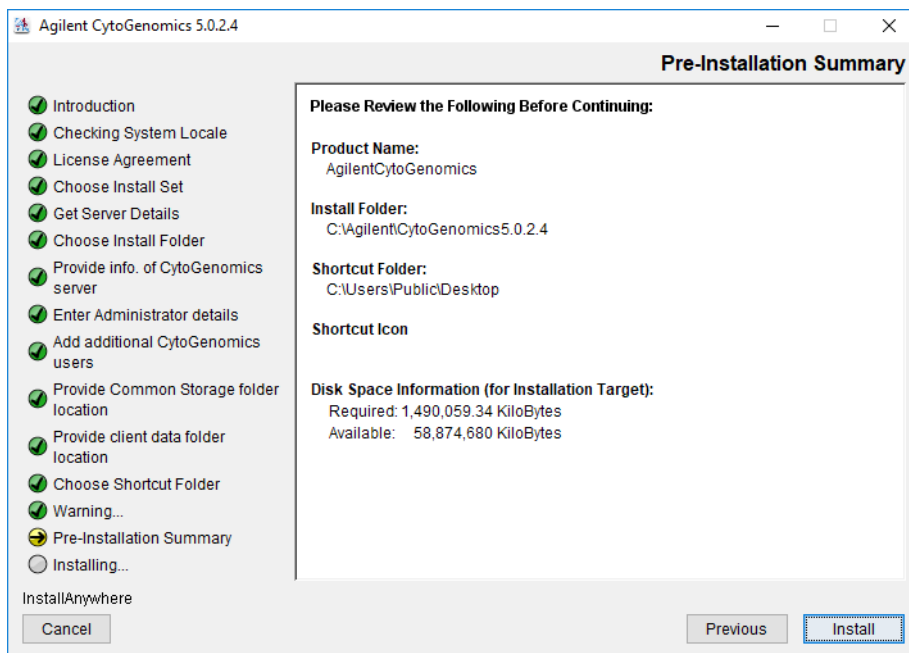


Figure 14 Pre-Installation Summary screen – Windows

23 Review the installation setup information. If you want to change any of the settings, click **Previous** and go back and change the settings. To start the installation, click **Install**.

The Installing Agilent CytoGenomics screen opens and is displayed until the installation is complete.

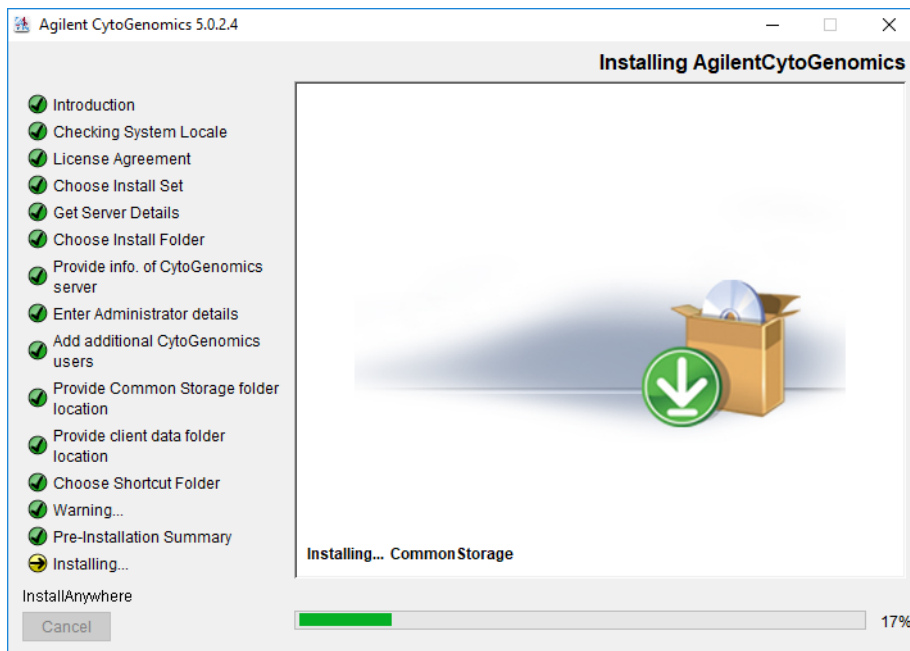


Figure 15 Installing Agilent CytoGenomics screen – Windows

When the installation is complete, the Installation Complete screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

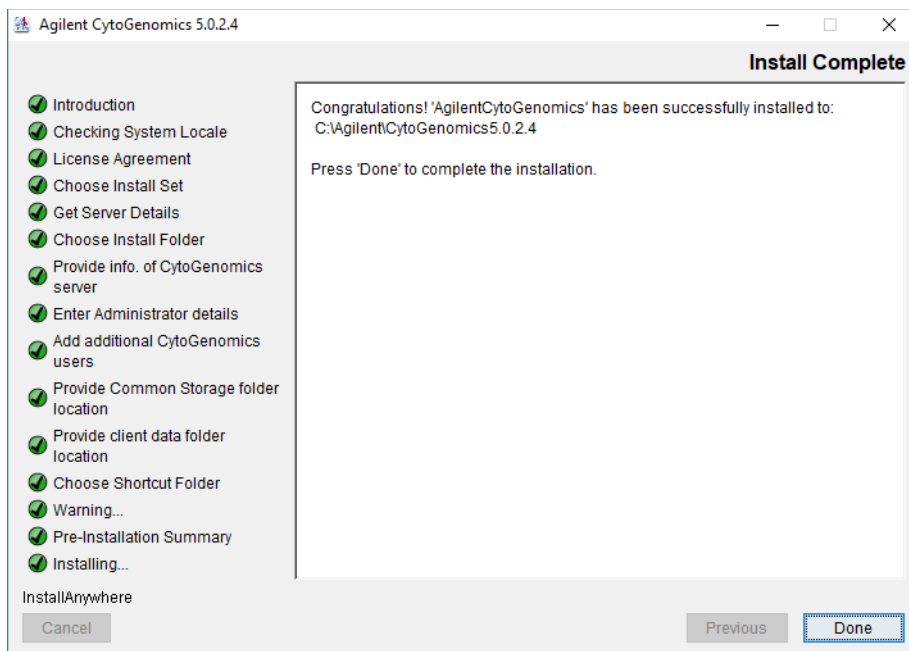


Figure 16 Installation Complete screen – Windows

24 Click **Done**.

The Launch Migration Utility dialog box opens asking if you want to launch Agilent’s migration utility tool designed to transfer data from a previous version of CytoGenomics server to the CytoGenomics 5.0 server.

NOTE

If the installer was unable to complete installation of CytoGenomics, then the Install Complete screen will note the installation errors and provide instructions on how to proceed.

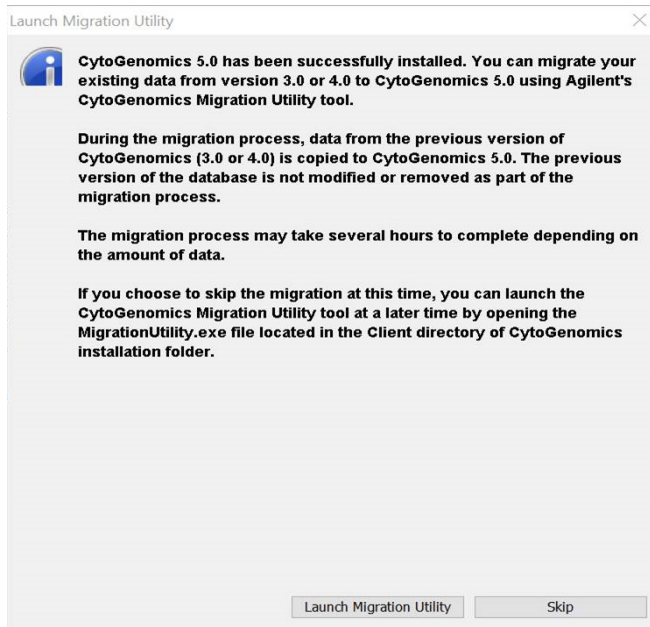


Figure 17 Launch Migration Utility dialog box – Windows

- 25** Click **Skip** to close the dialog box without launching the migration utility tool.

Because these instructions are for installing Agilent CytoGenomics 5.0 on a Windows system that does not have a previous version of CytoGenomics installed, there is no need to use the migration utility tool to transfer data.

- 26** Make sure the folder where you installed the software is shared with read/write permission for all client computers and for all users who run Agilent CytoGenomics.

Step 3. Install the client software on additional Windows computers

To use other computers to run Agilent CytoGenomics 5.0 client software (with a shared database), install the client software on each of the computers.

- 1 On each computer you want to run the Agilent CytoGenomics 5.0 program, run the installation wizard, as described in “[Step 2. Install the database server and client on a single Windows computer](#)” on page 20.
- 2 When the Choose Install Set screen opens, select **Only Client**.

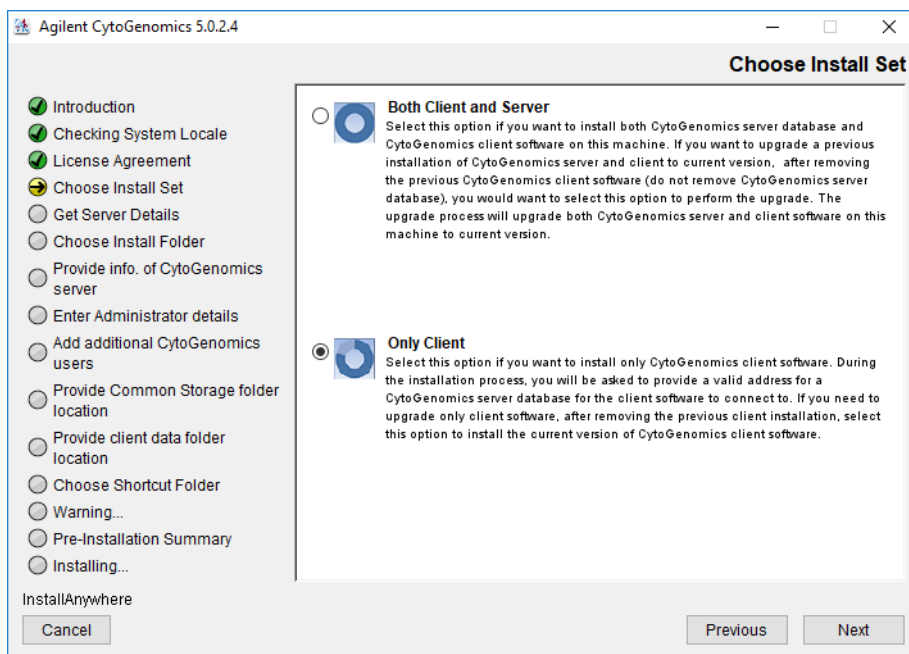


Figure 18 Choose Install Set screen – Windows – select **Only Client**

- 3 Click **Next**.

The Choose Install Folder screen opens.

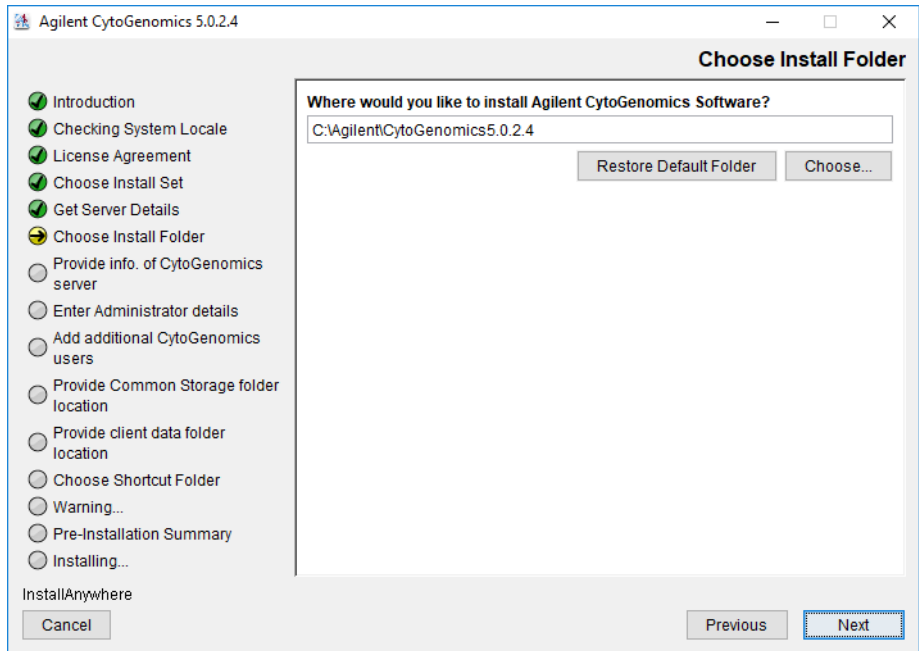


Figure 19 Choose Install Folder screen – Windows

- 4 By default, the program is installed in C:\Agilent\AgilentCytoGenomics5.0.X.X (where X.X is the revision number). Leave this folder as-is, or type the location and name of the folder where you want to install Agilent CytoGenomics 5.0.

OR

Click **Choose** and browse to a location where you want to install the program, and then click **Save**.

- 5 Click **Next**.

The Provide Info. of CytoGenomics server screen opens.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

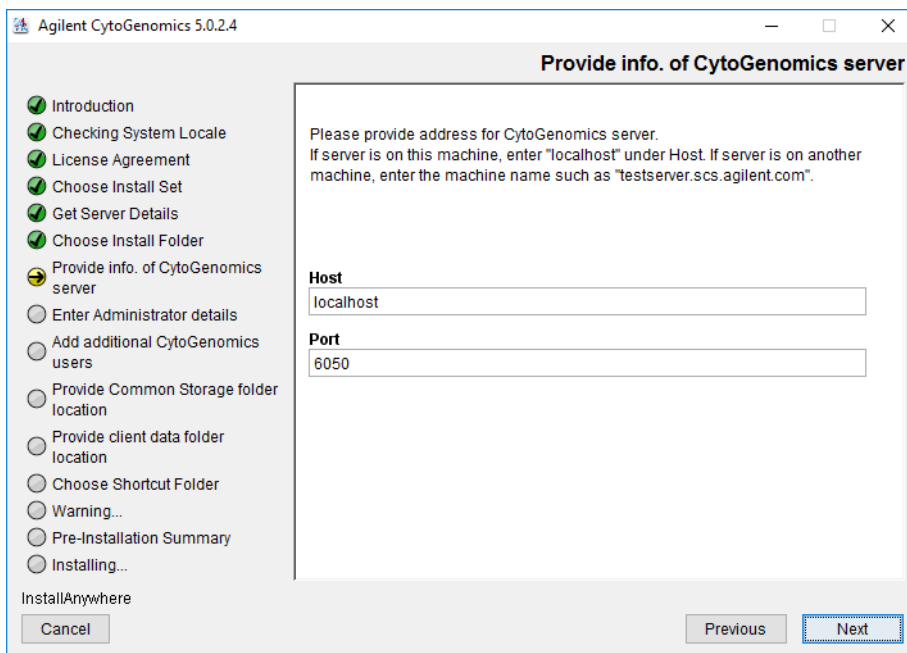


Figure 20 Provide Info. of CytoGenomics server screen – Windows

- 6 Under **Host**, type the assigned computer name or IP address of the computer (from [step 2](#)) where you installed the Agilent CytoGenomics 5.0 database server. (To find the computer name of your current computer, click **Start > Control Panel > System**.)
- 7 Accept the default **Port**, unless your system administrator tells you otherwise.
- 8 Click **Next**.

Installer immediately tests the connection to the specified database. If incorrect database parameters are provided, or the PostgreSQL database service is not running, a message is displayed.

You can choose to complete the installation and change database connection parameters when launching the client software. However, Agilent recommends that you provide the correct database information during installation.

Continue the installation wizard. The next screen is the Choose client data folder screen. Follow the instructions for the installation wizard to enter the client data folder location, choose a shortcut location, and start the installation. For information, see [step 19](#) through [step 26](#) in the section “[Step 2. Install the database server and client on a single Windows computer](#)” on page 20.

Step 4. Start the Agilent CytoGenomics 5.0 client software

After you install the Agilent CytoGenomics database server and client programs as described in this manual, start the program.

- Double-click the Agilent CytoGenomics 5.0 icon, or select Agilent CytoGenomics 5.0 from the Windows Start menu.

NOTE

How you start the program depends on the choice you made for Shortcut Folder during installation of the Agilent CytoGenomics client.

Step 5. Log in to Agilent CytoGenomics

After you start the program, the Login screen opens. The username and domain for the administrator set during installation appears in the Sign In area of the screen.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

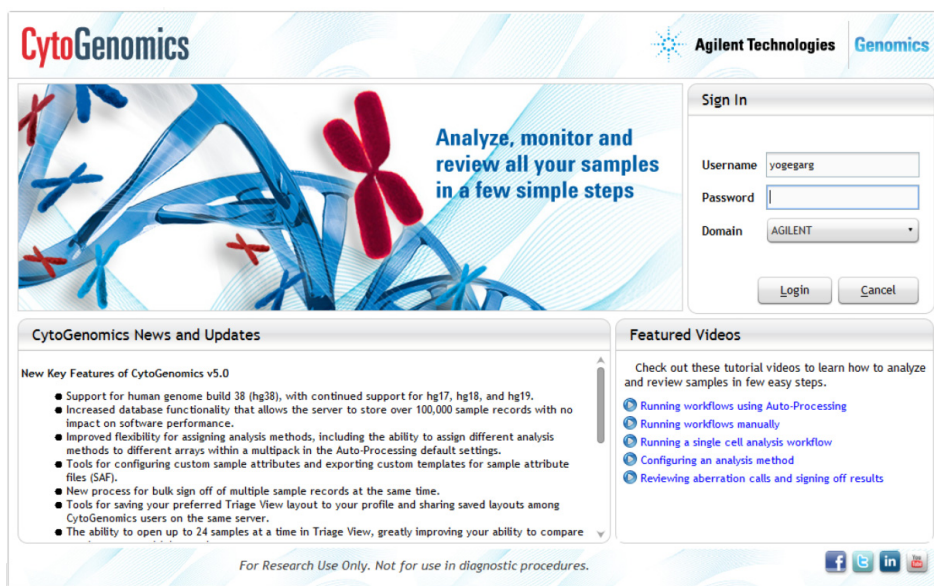


Figure 21 Login screen

1 Enter your username and password.

To change the domain, click the arrow next to **Domain** and select the correct domain. Otherwise, leave it as-is.

2 Click **Login**.

After the software is started, users with administrator privileges can add additional users and assign their roles.

NOTE

If you are using a local machine that is not part of a domain, you will see a message that the software cannot find the domain. Click **OK** to ignore this message.

If the installer did not properly share the common storage folder during installation, then an Application Initialization Error message will open notifying you that you do not have access to the common storage folder. This error message provides instructions on how to share the folder with all users. Follow the steps in the error message, then click **OK** to close the message and restart the application.

Step 6. Add your software license

When you log in for the first time, the Enter License dialog box opens (Figure 22).

- 1 If you do not already have the license file saved to your computer, click **Request for license** to open the Agilent website to the CytoGenomics license request page. Complete the fields in the request form to obtain a CytoGenomics software license. The license is contained in a text file and sent by email.
- 2 At the bottom of the Enter License dialog box, click **Browse**.
The Open dialog box opens.

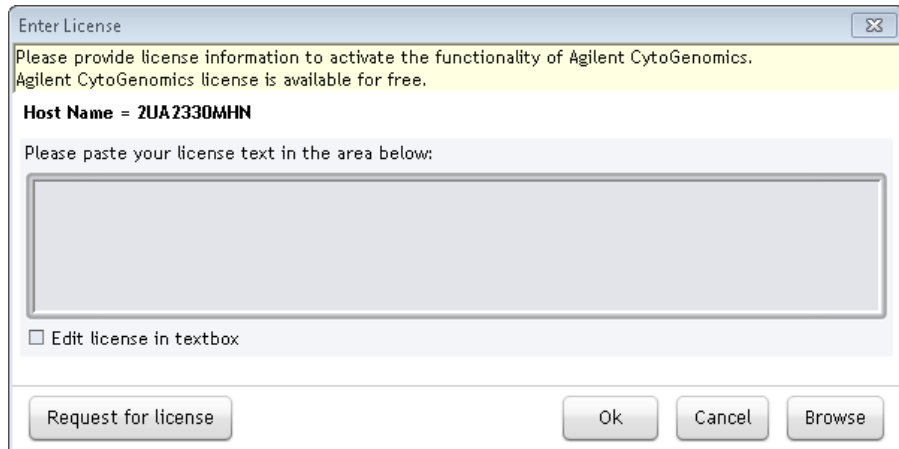


Figure 22 Enter License dialog box

- 3 Browse to the saved license text file. Select the file and click **Open**.
The program loads the contents of the file into the text box on the Enter License dialog box.
- 4 Click **OK**.

The Agilent CytoGenomics program opens to the Home screen (Figure 23).

Upon opening the program, you may see a message notifying you that updates are available for the software.

2 Installation Instructions for Windows

Installing Agilent CytoGenomics 5.0 – Windows

- If you want to download the available updates at this time, mark the check box next to **Software update** then click **Download**. During the download process, you are blocked from using CytoGenomics. Download time is dependent on internet speed.
- If you do not want to download the available update at this time, click **Cancel**.
- To stop being automatically notified of available updates, clear the check box labeled **Check for updates at application start up**. This check box is also accessible from Local Settings tab of the Configure Settings > Settings screen of CytoGenomics.

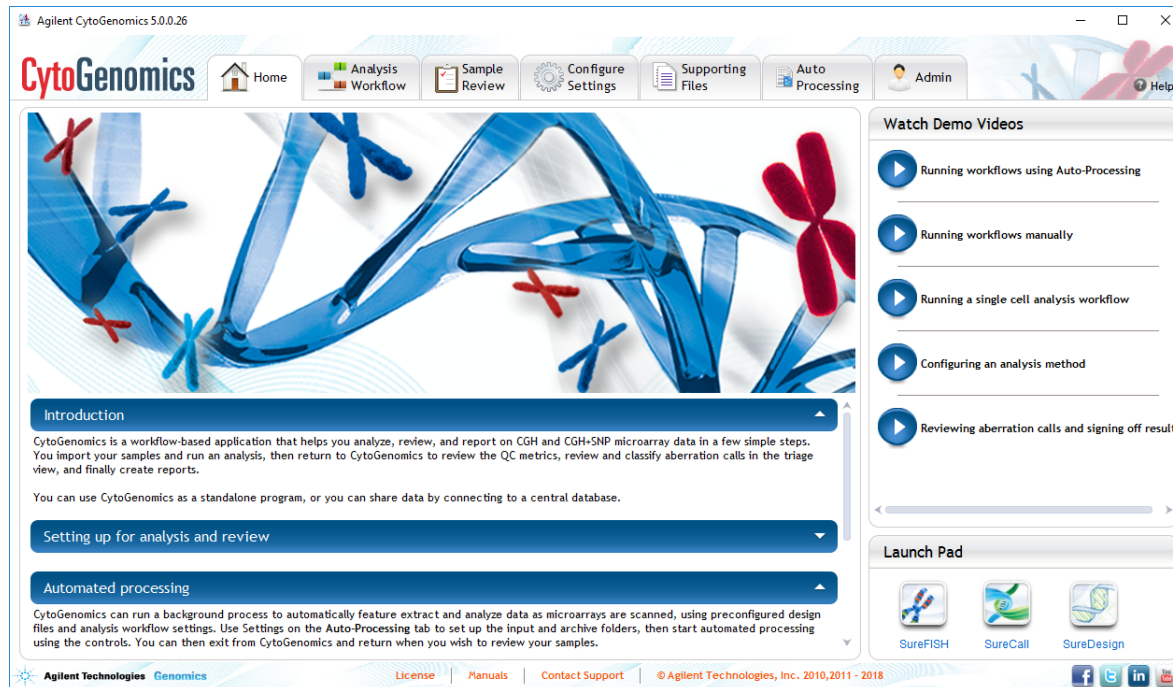


Figure 23 Agilent CytoGenomics program Home screen

Step 7. Add users and assign roles

Before other users can log in to Agilent CytoGenomics, a user with administrator privileges must add the users and assign user roles (if the users were not added during installation of the server). The user role determines what capabilities the user has within Agilent CytoGenomics. For information on what capabilities each user role has, see *“User Roles and Capabilities”* on page 16. For information on how to add users and assign roles, see *“To manage users and roles”* on page 110.

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows

These instructions are for users of CytoGenomics 3.0 or 4.0 who want to upgrade to CytoGenomics 5.0 and migrate all sample records from the previous database to the new CytoGenomics 5.0 database.

NOTE

CytoGenomics 5.0 is compatible with and can co-exist on the same machine as CytoGenomics 3.0 or 4.0. It is not necessary to uninstall the previous version of CytoGenomics (i.e., v3.0 or 4.0) in order to install v5.0.

Step 1. Install CytoGenomics 5.0 on the server computer

Follow these steps to install Agilent CytoGenomics 5.0 on the computer where the CytoGenomics server software is installed.

- 1 Download Agilent CytoGenomics 5.0.X.X from Agilent Technologies and check the system requirements.

See “a. [Download the Windows version of the software and release notes](#)” on page 18 for detailed instructions on this step.

- 2 Double-click the Agilent CytoGenomics 5.0 application file that you downloaded from Agilent.
- 3 On the Introduction screen, read the introductory information and then click **Next**.

The License Agreement screen opens.

- 4 Read the license agreement. Click **I accept the terms of the License Agreement** and then click **Next**.

The Choose Install Set screen opens.

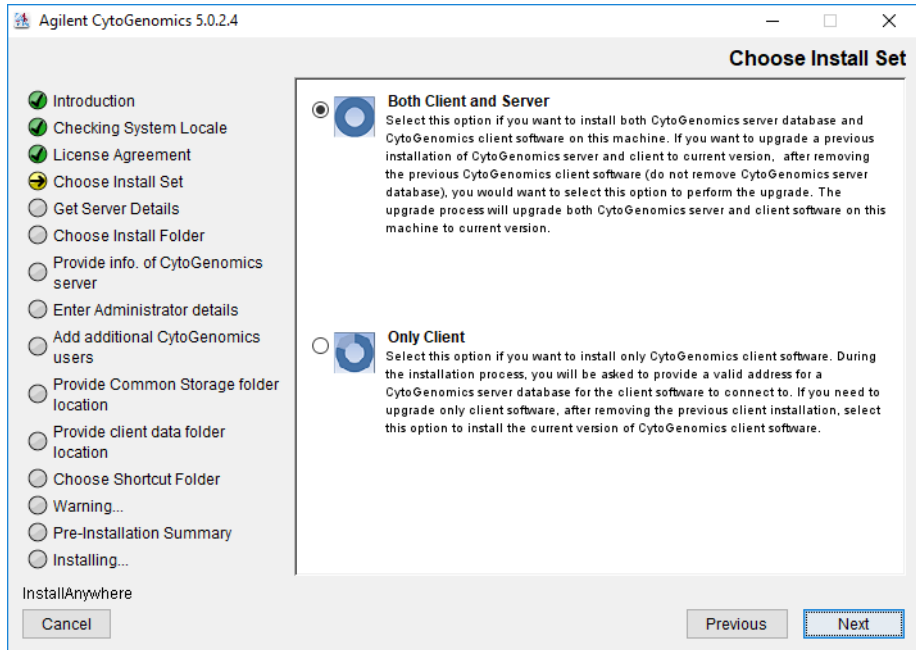


Figure 24 Choose Install Set screen – Windows – select **Both Client and Server**

- 5** On the Choose Install Set screen, select **Both Client and Server**. Click **Next**.

The following dialog box opens to inform you that you have selected to install the CytoGenomics server and client.

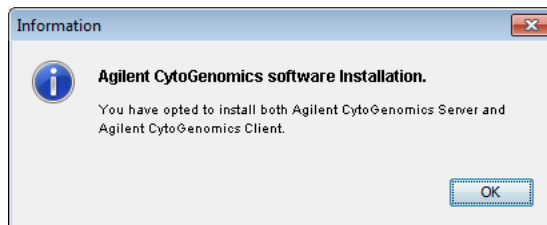


Figure 25 Dialog box – Agilent CytoGenomics software installation

- 6** Click **OK** in the dialog box to continue.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows

- 7 Continue the installation wizard. The next screen is the Choose Install Folder screen. Follow the instructions for the installation wizard to enter the client data folder location, choose a shortcut location, and start the installation. For information, see [step 8](#) through [step 24](#) in the section “[Step 2. Install the database server and client on a single Windows computer](#)” on page 20.

At the completion of installation, the Launch Migration Utility dialog box opens asking if you want to launch Agilent’s migration utility tool designed to transfer data from a previous version of CytoGenomics server to the CytoGenomics 5.0. server.

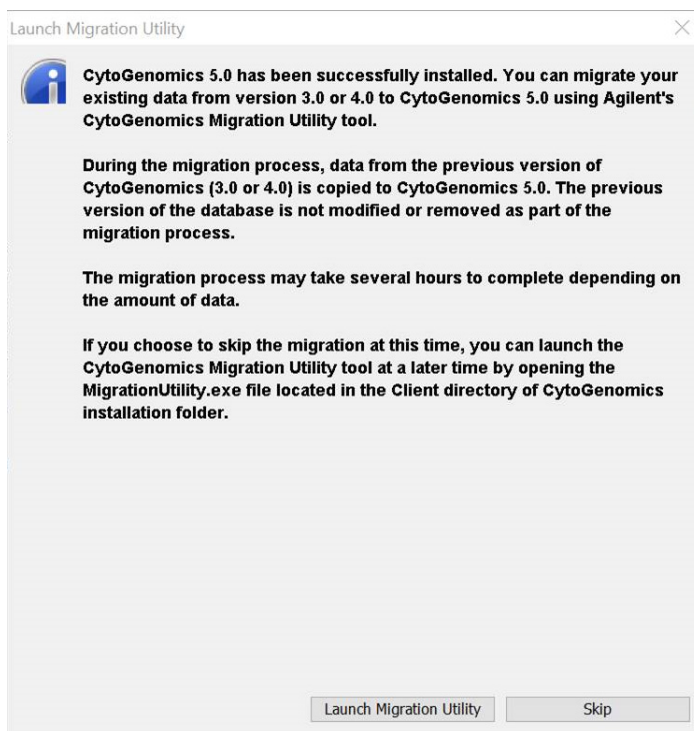


Figure 26 Launch Migration Utility dialog box – Windows

- 8 Depending on your needs, click either **Launch Migration Utility** or **Skip**.

- Click **Launch Migration Utility** if you have data saved to the previous version of CytoGenomics server (e.g., sample results, track files, and other supporting files) and you want to transfer that data to the CytoGenomics 5.0 server. The migration utility tool launches on your PC. Proceed to [“Step 2. Migrate data to CytoGenomics 5.0 server”](#) on page 47.
- Click **Skip** if you do not want to transfer any data (e.g., sample results, track files, and other supporting files) from the previous version of CytoGenomics server to the CytoGenomics 5.0 server at this time. The dialog box will close and that data will remain on the previous version of the server. Proceed to [“Step 3. Upgrade to CytoGenomics 5.0 on the client computers”](#) on page 52.

NOTE

To protect against loss of data and ensure migration of all analyzed samples, Agilent strongly recommends migrating data immediately upon completion of the CytoGenomics 5.0 installation.

If you click **Skip** with the intention of migrating your data at a later time, be aware that if CytoGenomics 5.0 already contains data at the time of migration, the Migration Utility application will delete that data and overwrite it with data from the previous version of CytoGenomics.

Step 2. Migrate data to CytoGenomics 5.0 server

If you selected to launch the migration utility tool, follow these step to use the tool to transfer data from the previous version of CytoGenomics server to CytoGenomics server 5.0. If you do not want to transfer data, bypass

2 Installation Instructions for Windows

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows

this step and proceed to “Step 3. Upgrade to CytoGenomics 5.0 on the client computers” on page 52.

- 1 On the Sign In screen of the migration utility application (Figure 27), review the list of migration prerequisites and make sure they are all met before you continue.

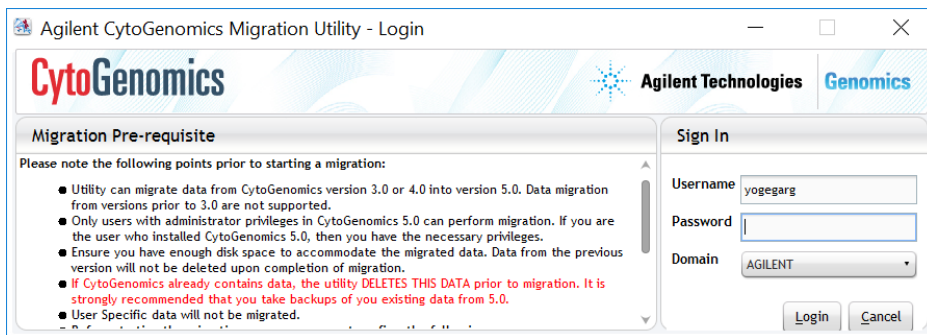


Figure 27 CytoGenomics Migration Utility application login screen

- 2 Log in with your username and password. The username and domain for the administrator set during installation appears in the Sign In area of the screen. (Only accounts with administrator privileges can migrate data.)

The CytoGenomics Migration Utility application opens.

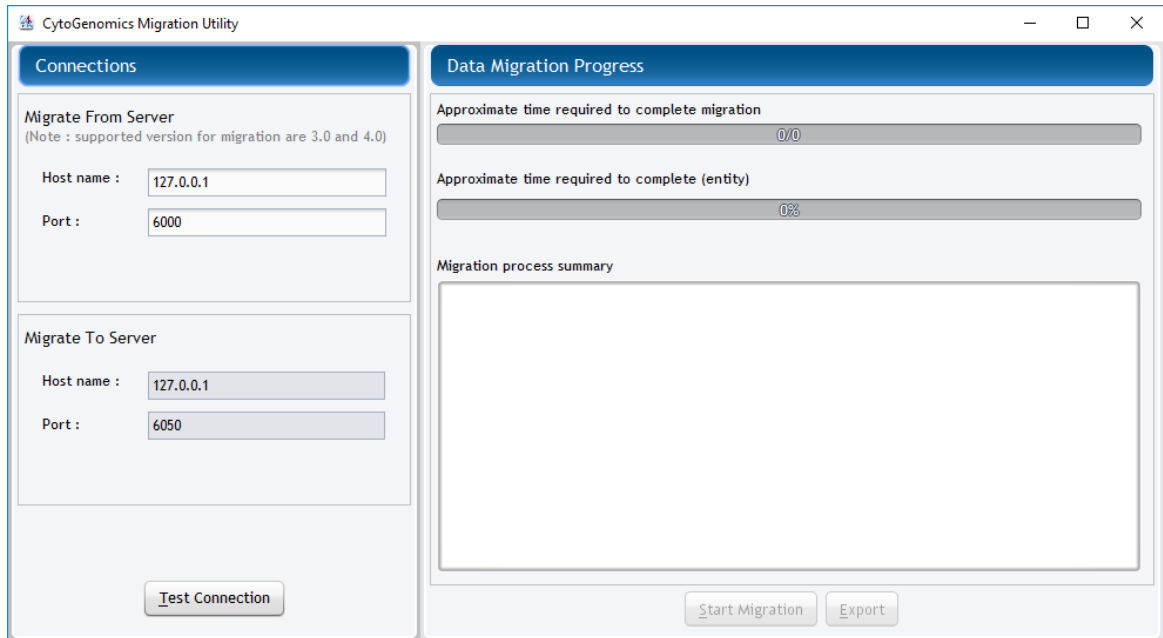


Figure 28 CytoGenomics Migration Utility application

- 3 Under **Migrate From Server**, enter the host name (i.e., IP address) and port number where the previous version of CytoGenomics server is installed.
- 4 Under **Migrate To Server**, enter the host name (i.e., IP address) and port number where the CytoGenomics 5.0 server is installed. If CytoGenomics 5.0 server is installed on the same PC as the previous version of CytoGenomics server, then the host names are the same.

NOTE

If CytoGenomics 5.0 server is installed on the same PC as the previous version of CytoGenomics server, then the host names of the two servers are the same.

- 5 Click **Test Connection**.

The screen displays the results of connection test with each server.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows

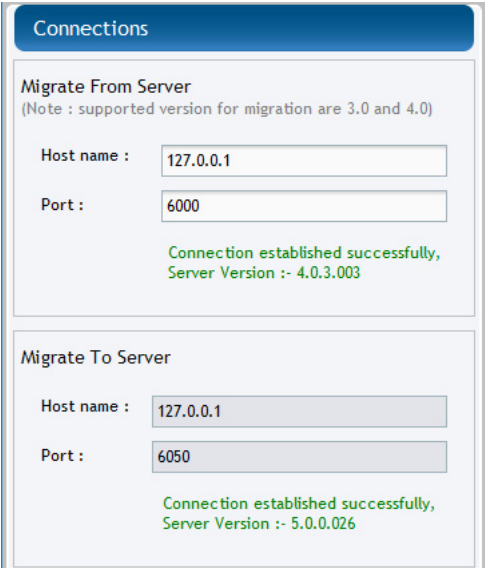


Figure 29 CytoGenomics Migration Utility application – Connections established

NOTE

If the application was unable to successfully establish a connection to one or both servers, verify that the correct host name and port number are entered and test the connection again. You cannot proceed to data migration until the application connects to the servers.

6 Once the application has successfully connected to the servers, click **Start Migration**.

The Warning message box opens.

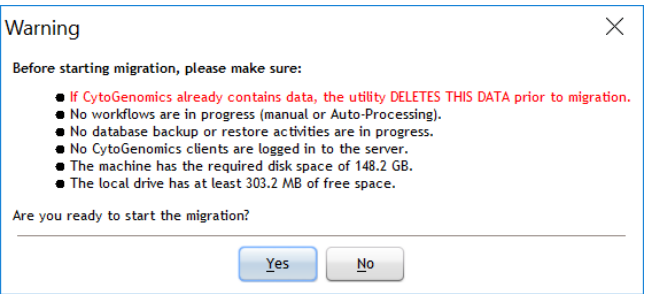


Figure 30 Warning message box

- 7 As instructed in the Warning message box, make sure that CytoGenomics is not actively running any workflows or performing any database backup or restore activities and that all clients are logged out. If your CytoGenomics 5.0 database already contains data, make sure that data has been backed up. Then, click **Yes** to continue with the migration.

The right side of the screen (under **Data Migration Process**) displays the progress of the migration and the migration process summary.

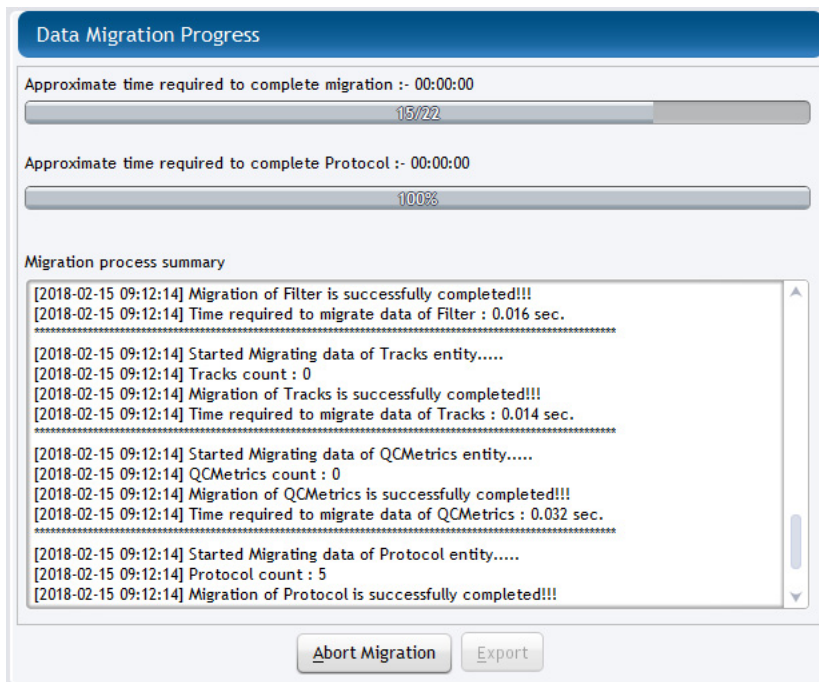


Figure 31 CytoGenomics Migration Utility application – Migration in progress

Upon completion of the migration process, the following message box opens.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows

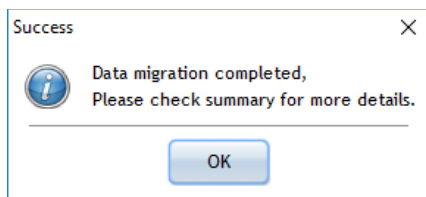


Figure 32 CytoGenomics Migration Utility application – Data migration complete

- 8 Click **OK** to close the message box.
- 9 If desired, click **Export** to export a summary of the migration job.
- 10 Click the X in the top right corner of the Migration Utility application window to close the application.

Step 3. Upgrade to CytoGenomics 5.0 on the client computers

Follow these steps to install Agilent CytoGenomics 5.0 on each client computer.

- 1 Double-click the Agilent CytoGenomics 5.0.X.X application file that you downloaded from Agilent.
- 2 On the Introduction screen, read the introductory information then click **Next**.
The License Agreement screen opens.
- 3 Read the license agreement. Click **I accept the terms of the License Agreement** and then click **Next**.
The Choose Install Set screen opens.

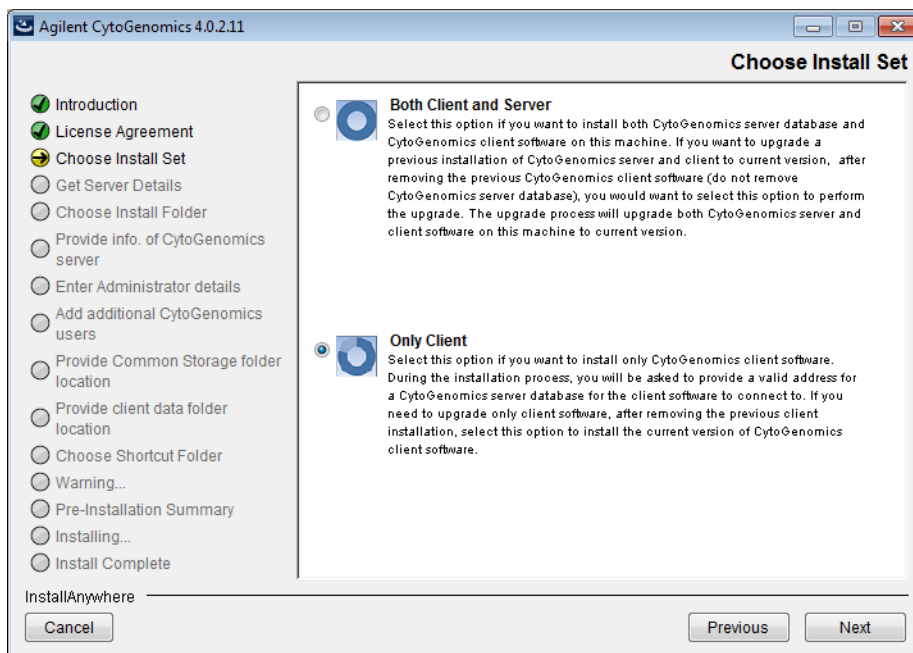


Figure 33 Choose Install Set screen – Windows – select **Only Client**

4 On the Choose Install Set screen, select **Only Client**. Click **Next**.

The following dialog box opens to inform you that you have selected to install only the Agilent CytoGenomics 5.0 client.

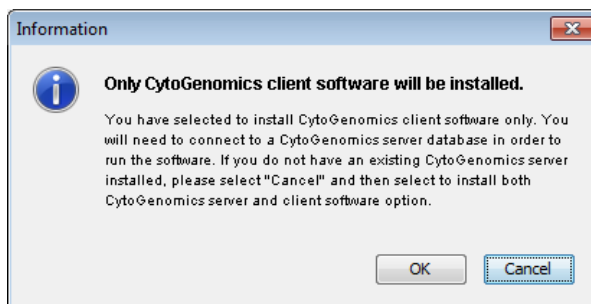


Figure 34 Dialog box – Only CytoGenomics client software will be installed

5 Click **OK** in the dialog box to continue.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Windows

The Choose Install Folder screen opens.

- 6 At the Choose Install Folder screen, the default folder for installation of the program is C:\Agilent\AgilentCytoGenomics5.0.X.X (where X.X is the revision number). Leave this folder as-is, or click **Choose** and browse to a location where you want to install the program, and then click **Save**.
- 7 Click **Next**.
- 8 Continue the installation wizard. Follow the instructions to provide server information, choose a shortcut location, and start the installation. See [step 6](#) through [step 8](#) in the section “[Step 3. Install the client software on additional Windows computers](#)” on page 36.

Step 4. Start the Agilent CytoGenomics 5.0 client software, log in, and add new software license

- 1 On the server and client computers, start the Agilent CytoGenomics program.
See “[Step 4. Start the Agilent CytoGenomics 5.0 client software](#)” on page 39.
- 2 Log in.
See “[Step 5. Log in to Agilent CytoGenomics](#)” on page 39.
- 3 Enter your updated license information.
See “[Step 6. Add your software license](#)” on page 41.

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

If you are using CytoGenomics 5.0.0 or 5.0.1 and want to upgrade to CytoGenomics 5.0.2, you only need to upgrade the CytoGenomics client application. You do not need to upgrade to the CytoGenomics server application or migrate data from the previous version to the new version.

Step 1. Install CytoGenomics 5.0.2 client and uninstall CytoGenomics 5.0.0 or 5.0.1 client

- 1 Download Agilent CytoGenomics 5.0.2 from Agilent Technologies and check the system requirements.

See “a. [Download the Windows version of the software and release notes](#)” on page 18 for detailed instructions on this step.

- 2 Double-click the Agilent CytoGenomics 5.0.2 application file that you downloaded from Agilent.

The InstallAnywhere dialog box is displayed. Then, the installation wizard opens to the Introduction screen.

- 3 Read the introductory information, then click **Next**.

The following message box opens notifying you that the existing version of CytoGenomics client must be uninstalled before installing CytoGenomics 5.0.2 client.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

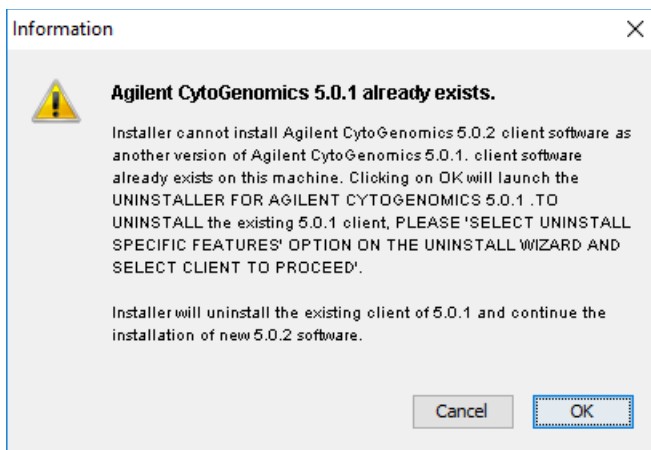


Figure 35 “Agilent CytoGenomics 5.0.1 already exists” message box

4 Click **OK** in the message box.

The CytoGenomics uninstaller opens.

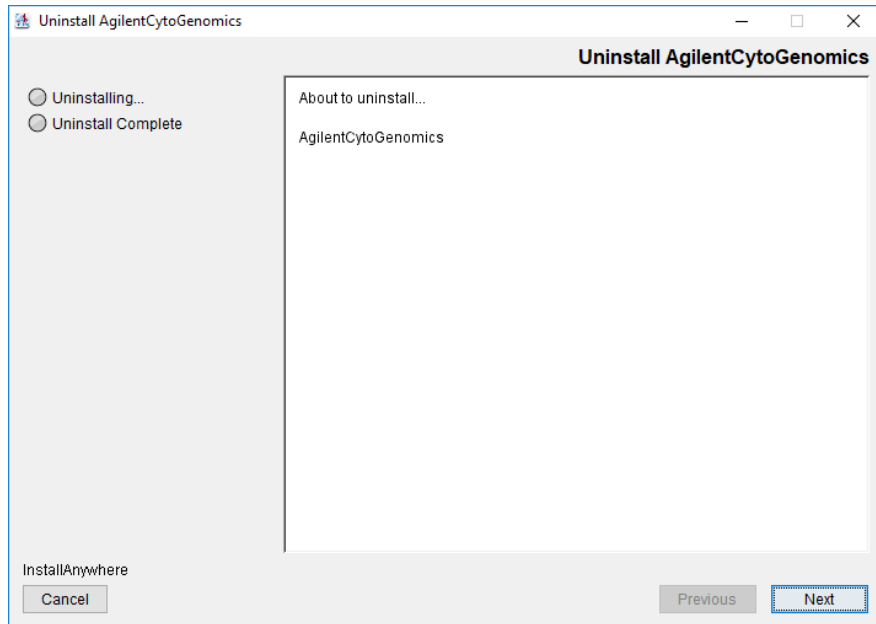


Figure 36 CytoGenomics Uninstaller – Windows

- 5 Click **Next**.
- 6 Read the warning information. You are cautioned not to stop the Uninstaller once it has started to remove the programs. Click **Next**. The Uninstall Options screen opens.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

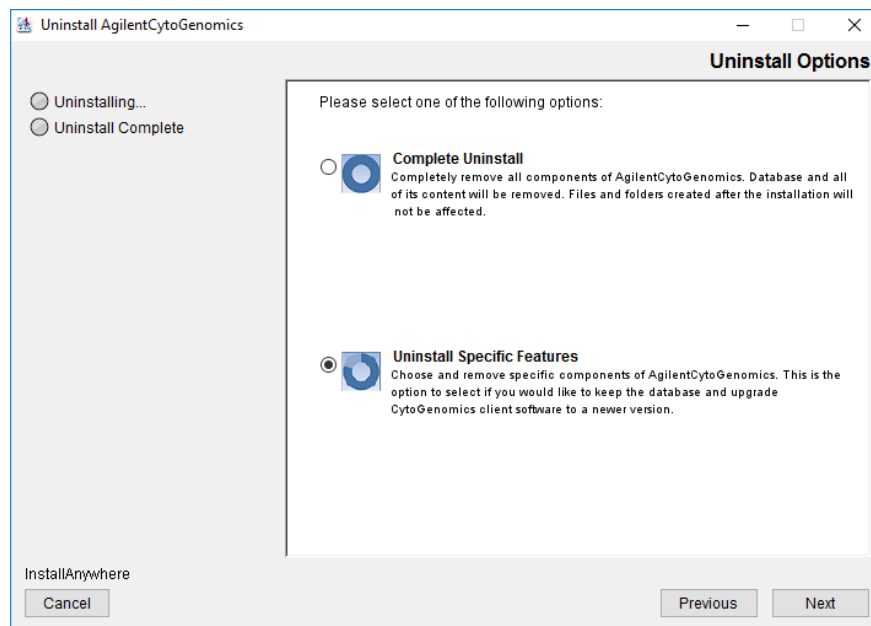


Figure 37 Uninstall Options screen – Windows – select **Uninstall Specific Features**

7 Select **Uninstall Specific Features**.

The Choose Product Features screen opens.

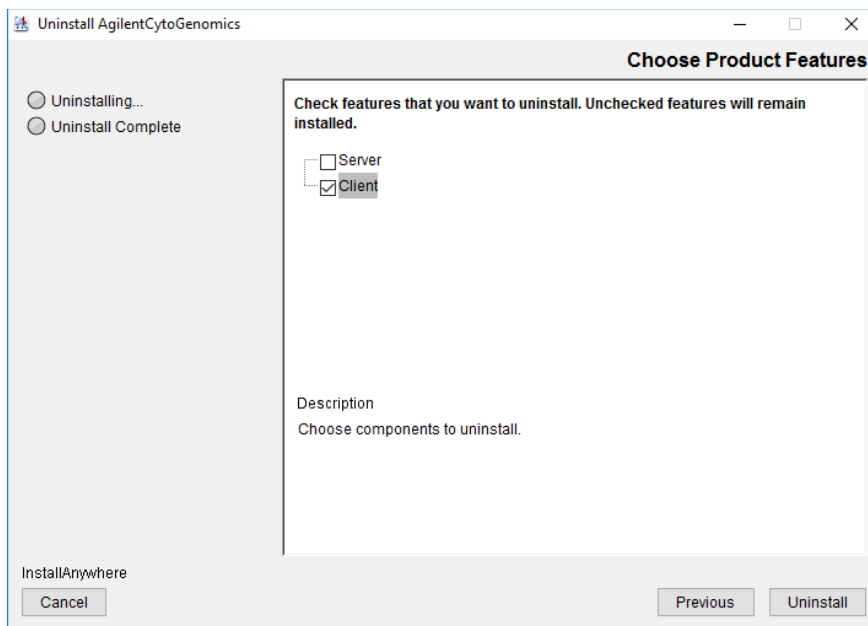


Figure 38 Choose Product Features screen – Windows – select **Client**

8 Mark **Client**, then click **Uninstall**.

NOTE

Do not mark **Server** on the Choose Product Features screen. Doing so will result in removal of all previously analyzed data from your existing CytoGenomics server.

- 9** In the confirmation message box that opens, click **Yes** to continue uninstalling CytoGenomics client for version 5.0.0 or 5.0.1.
- 10** When the uninstallation is complete, a message opens. Click **Done**.
The uninstaller closes and you are directed to the License Agreement screen of the CytoGenomics 5.0.2 installer.
- 11** Continue the installation wizard, as described in “[Step 2. Install the database server and client on a single Windows computer](#)” on page 20.
- 12** When the Choose Install Set screen opens, select **Only Client**.

2 Installation Instructions for Windows

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

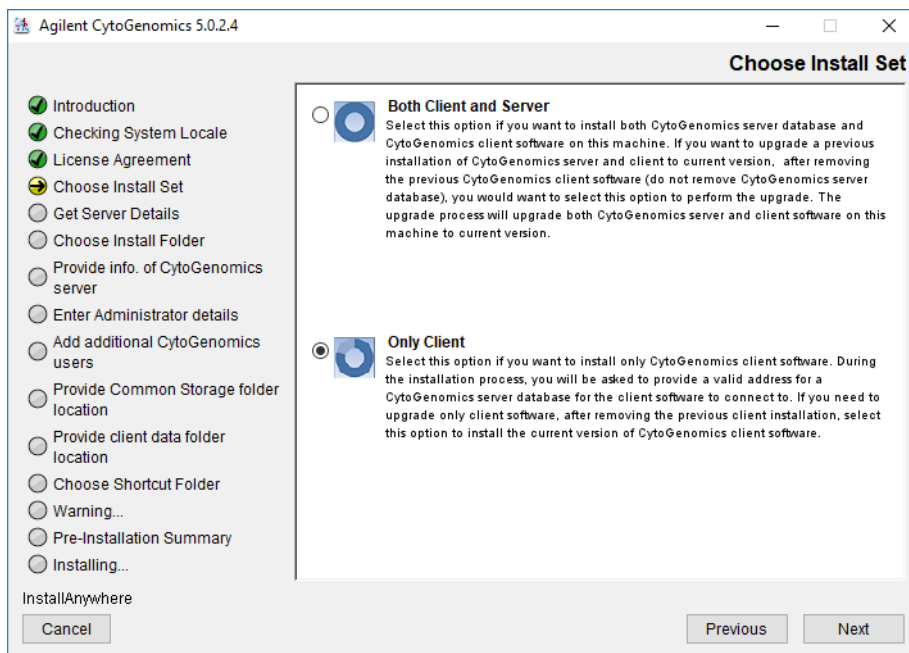


Figure 39 Choose Install Set screen – Windows – select **Only Client**

13 Click **Next**.

14 Continue the installation wizard. For information, refer to “[Step 3. Install the client software on additional Windows computers](#)” on page 36.

Step 2. Start the Agilent CytoGenomics 5.0.2 client software, log in, and add new software license

- 1** Start the Agilent CytoGenomics program.
See “[Step 4. Start the Agilent CytoGenomics 5.0 client software](#)” on page 39.
- 2** Log in.
See “[Step 5. Log in to Agilent CytoGenomics](#)” on page 39.
- 3** Enter your updated license information.
See “[Step 6. Add your software license](#)” on page 41.

Uninstalling Agilent CytoGenomics on Windows computers

The Uninstall program for Agilent CytoGenomics provides an easy way to remove the Agilent CytoGenomics client, server, and PostgreSQL programs from your computer.

NOTE

CytoGenomics 5.0 is compatible with and can co-exist on the same machine as CytoGenomics 3.0 or 4.0. It is not necessary to uninstall the previous version of CytoGenomics (i.e., v3.0 or 4.0) in order to install v5.0.

When you completely uninstall Agilent CytoGenomics, the database and the data contained in it are completely removed. Do not uninstall the entire program unless you have backed up your database or no longer wish to use it.

- 1 Start the **Uninstall AgilentCytoGenomics.exe** file, located in the \Uninstall_AgilentCytoGenomics folder of your program folder.
The Uninstaller program starts.
- 2 Read the introductory information, and then click **Next**.
- 3 Read the warning information. You are cautioned not to stop the Uninstaller once it has started to remove the programs. Click **Next**.
The Uninstall Options screen opens.

2 Installation Instructions for Windows

Uninstalling Agilent CytoGenomics on Windows computers

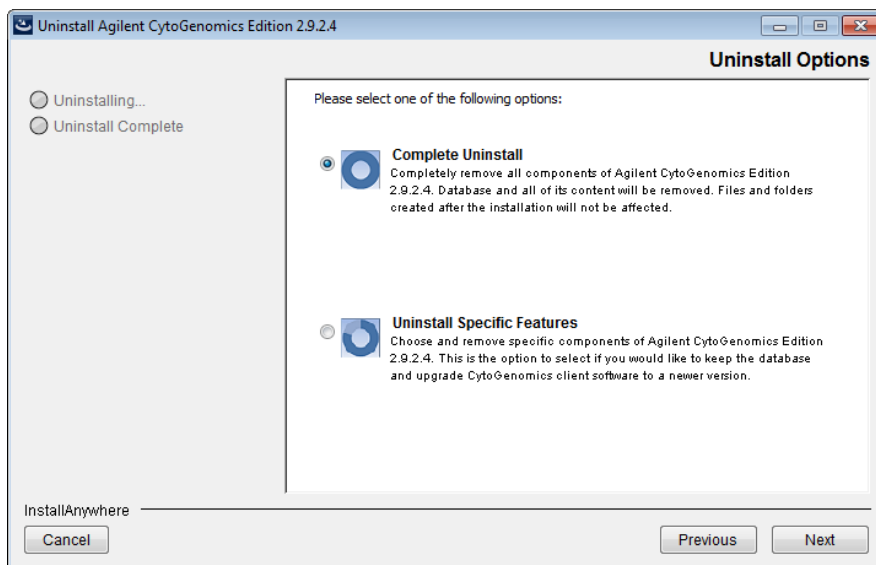


Figure 40 Uninstall Options screen – Windows – select **Complete Uninstall**

4 Select **Complete Uninstall**.

Depending on which version of CytoGenomics you are uninstalling, you may see a warning message advising you to backup the database before proceeding with the uninstall.

5 When the uninstallation is complete, a message opens. Click **Done**.

The Agilent CytoGenomics client program, server program and database, and PostgreSQL program were removed from your computer. Program folders and data folders (such as data output folders) created after installation were not removed.

Troubleshooting

If the Agilent CytoGenomics program does not open

- ✓ Try opening the program by running the *run_CytoGenomics.bat* file.

- a Open the installation folder AgilentCytoGenomics5.0.X.X.

- b Double-click *run_CytoGenomics.bat*.

Agilent CytoGenomics opens and you can operate the program normally.

If the program still does not open, try adjusting the amount of RAM allocated to the program using the approach provided below.

- ✓ Adjust the amount of RAM that is allocated to the program.

- a Open the Client folder inside the AgilentCytoGenomics5.0.X.X installation folder.

- b Open the file *Agilent CytoGenomics.lax* in Notepad.

- c In this file, locate the following text.

- lax.nl.java.option.additional=Xms1000m

- d Change Xms1000m to Xms900m.

- e Save and close the file.

- f Open the Agilent CytoGenomics client program.

- Double-click the Agilent CytoGenomics 5.0 icon, or select Agilent CytoGenomics 5.0 from the Windows Start menu.

If Agilent CytoGenomics opens, you can operate the program normally.

If Agilent CytoGenomics still does not open, repeat [step a](#) through [step f](#), and reduce the Xms value by 100 again. Continue to reduce this value in increments of 100 until the program opens successfully.

2 Installation Instructions for Windows

Troubleshooting



3

Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh 66

Step 1. Prepare for installation 66

Step 2. Install the database server and client on a single Macintosh computer 68

Step 3. Install the client software on additional Macintosh computers 84

Step 4. Start the Agilent CytoGenomics 5.0 client software 87

Step 5. Log in to Agilent CytoGenomics 87

Step 6. Add your software license 89

Step 7. Add users and assign roles 90

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh 91

Step 1. Install CytoGenomics 5.0 on the server computer 91

Step 2. Migrate data to CytoGenomics 5.0 server 94

Step 3. Upgrade to CytoGenomics 5.0 on the client computers 98

Step 4. Start the Agilent CytoGenomics 5.0 client software, log in, and add new software license 98

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2 99

Uninstalling Agilent CytoGenomics on Macintosh computers 105

This chapter shows you how to install Agilent CytoGenomics 5.0 and migrate data from a previous version of CytoGenomics on systems running a Macintosh OS.

NOTE

The Macintosh version of CytoGenomics does not permit analysis of scanner image files. To run an analysis workflow, you must first extract the image file using the Agilent Feature Extraction program, then use the extracted FE file in the workflow.



Agilent Technologies

Installing Agilent CytoGenomics 5.0 – Macintosh

This section includes detailed instructions and important notes for installing Agilent CytoGenomics 5.0 on a Macintosh system that does not have a previous version of CytoGenomics installed.

Agilent CytoGenomics consists of two parts: a PostgreSQL database server and the client software. Once you have installed the database server and client software on one computer, you can install additional copies of the client software on client computers that have network access to the database computer.

For instructions on how to upgrade an existing Agilent CytoGenomics installation, see “[Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh](#)” on page 91.

Step 1. Prepare for installation

a. Download the Macintosh version of the software and release notes

- 1 Go to the CytoGenomics download page on the Agilent website at: www.agilent.com/en/download-agilent-cyto-genomics-software.
- 2 Click the **Download** link for the Macintosh version of the software.
- 3 Save the Installer on your computer. (Do not start the installation at this time.)

The Mac installer is downloaded as a compressed folder (Agilent_CytoGenomics_5_0_X_X_MAC.zip). Do not start the installation at this time.
- 4 Click the **Release Notes** link to download the software release notes.
- 5 Save the release notes to your computer.

b. Check Macintosh system requirements before you start

Check to make sure that your computer systems meet the following requirements. All client installations point to the computer on which the server software is installed. Requirements are listed in [Table 3](#).

NOTE

If an operating system is not listed in [Table 3](#), it is not supported.

Table 3 Requirements for Agilent CytoGenomics on Macintosh computers

Operating System	Macintosh OS Sierra, or Macintosh OS X El Capitan
Programs	Java 1.8 Any program that enables you to open PDF files (for example, Adobe Reader)
Processor	3 GHz Intel Core 2 Duo CPU or better
Working memory (RAM)	Minimum: 4 GB Recommended: 8 GB
Hard disk space	40 GB (For analysis of large datasets, more space is required)
Display resolution	1280 x 768 or higher

c. Set up permissions for installer and users

- Make sure that the user account used to install the Agilent CytoGenomics 5.0 software has administrator permissions on the installation computer.
- Make sure that every user account running the Agilent CytoGenomics 5.0 software has read/write permissions in the computer where the client software is installed. In addition, users must have read/write permissions in the installation folder of the computer where the database server is installed.

The way you set permissions varies between operating systems. Contact your network administrator for help in checking or changing user account permissions. If the user account permissions are not set correctly, the software will fail to install or run properly.

Step 2. Install the database server and client on a single Macintosh computer

NOTE

Only one copy of the server software is installed. To run the client software, the database server must be installed. The client is the program you run when you want to use Agilent CytoGenomics. You can run the client program on the same computer as the server, or from any other computer with the client program installed. After installation, the database server software runs silently in the background and manages the shared database for all Agilent CytoGenomics 5.0 users. All client computers must have network access to the server computer and users must have read/write permissions.

- 1** Log on to the Macintosh system where you want to install Agilent CytoGenomics 5.0 server. Use an account that has administrator permissions.
- 2** Locate the compressed folder (Agilent_CytoGenomics_5_0_X_X_MAC.zip) that you downloaded in Step 1, “[a. Download the Macintosh version of the software and release notes](#)” on page 66. Double-click the folder to expand the contents of the folder and save the installer on your computer.
- 3** Double-click the Mac installer.
The following window opens.

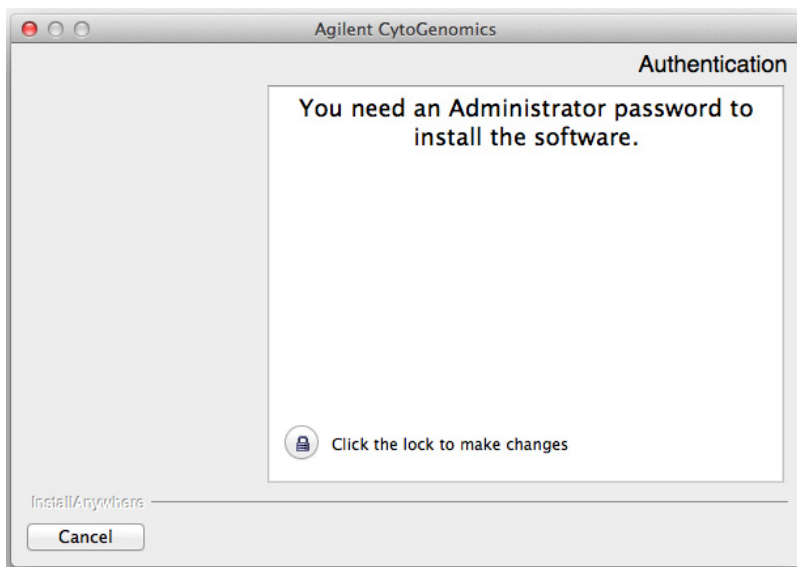


Figure 41 Authentication dialog box - preparing to install

- 4 Click directly on the lock icon near bottom of the dialog box.
The following dialog box opens.

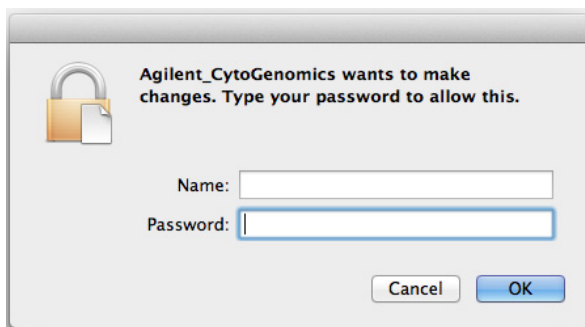


Figure 42 Administrator credentials dialog box - preparing to install

- 5 In this dialog box, enter the Name and Password for a user with administrator rights to the computer, then click **OK**.
The installation wizard starts.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

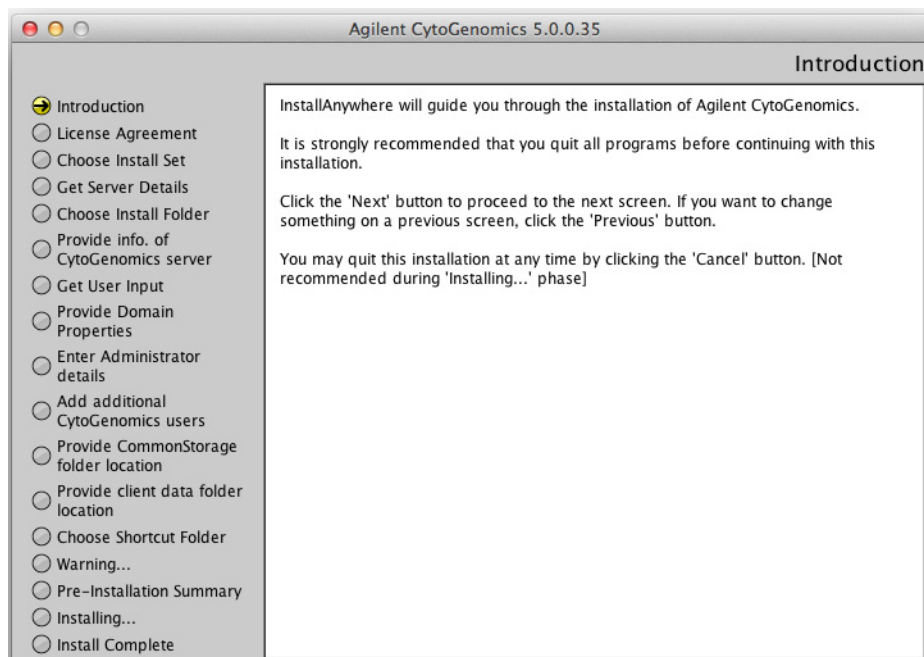


Figure 43 Introduction screen for the installation wizard – Macintosh

- 6 Read the introductory information.
- 7 Click **Next** to display the next section.
The License Agreement screen opens.

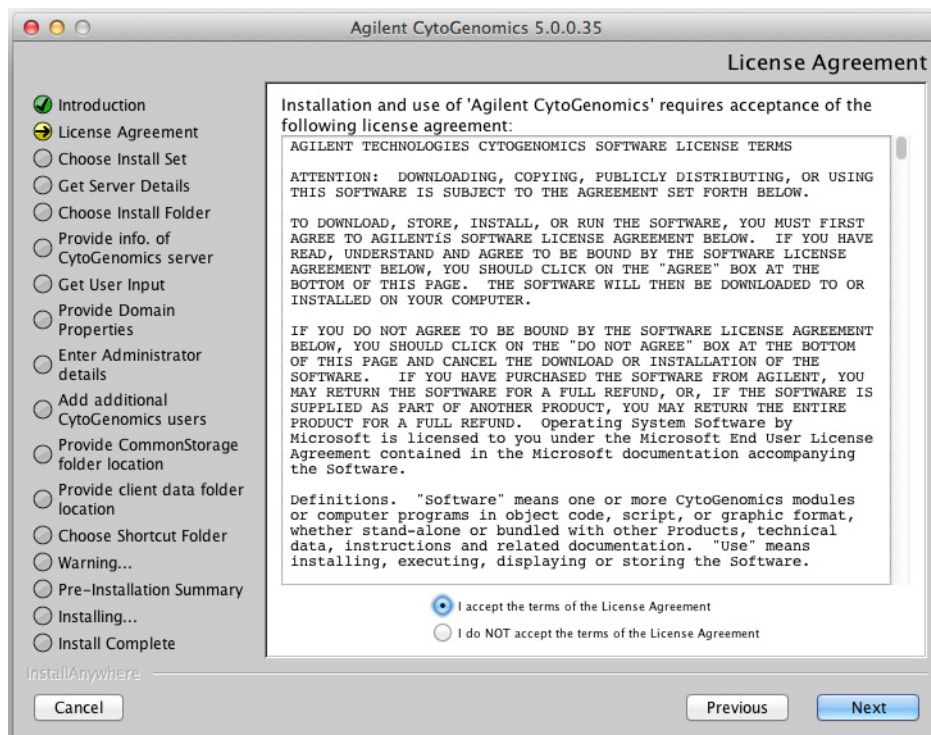


Figure 44 License Agreement screen – Macintosh

- 8 Read the license agreement. Click **I accept the terms of the License Agreement** and then click **Next**. (Click **Previous** to review information in a previous section.)

The Choose Install Set screen opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

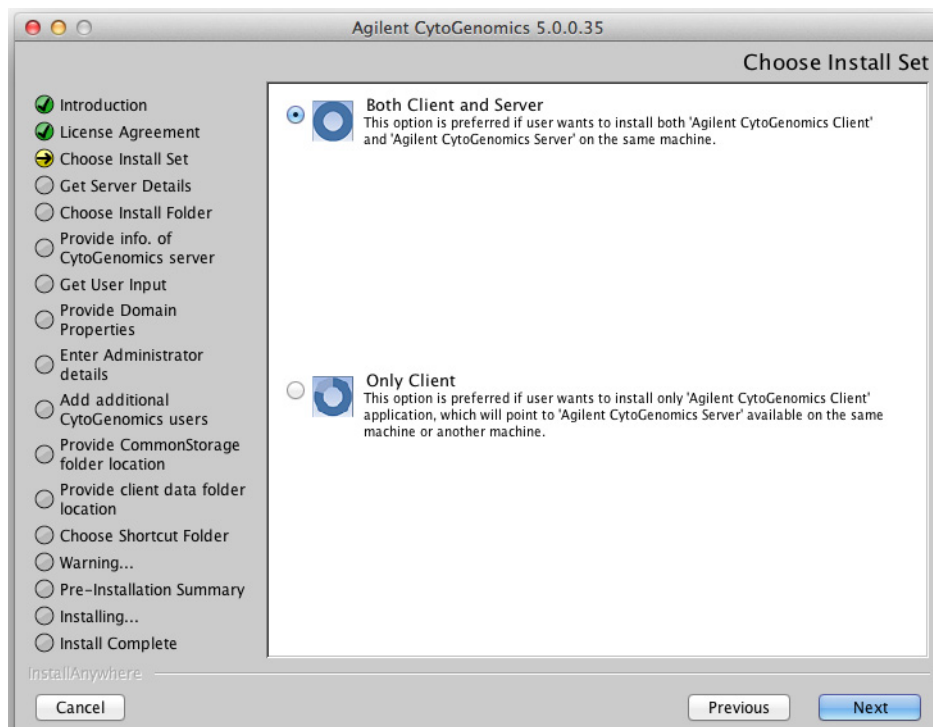


Figure 45 Choose Install Set screen – Macintosh

9 Click the icon next to **Both Client and Server** to install both the database server and client software on this computer.

10 Click **Next**.

11 An Information message opens.

Click **OK**. The Choose Install Folder screen opens.

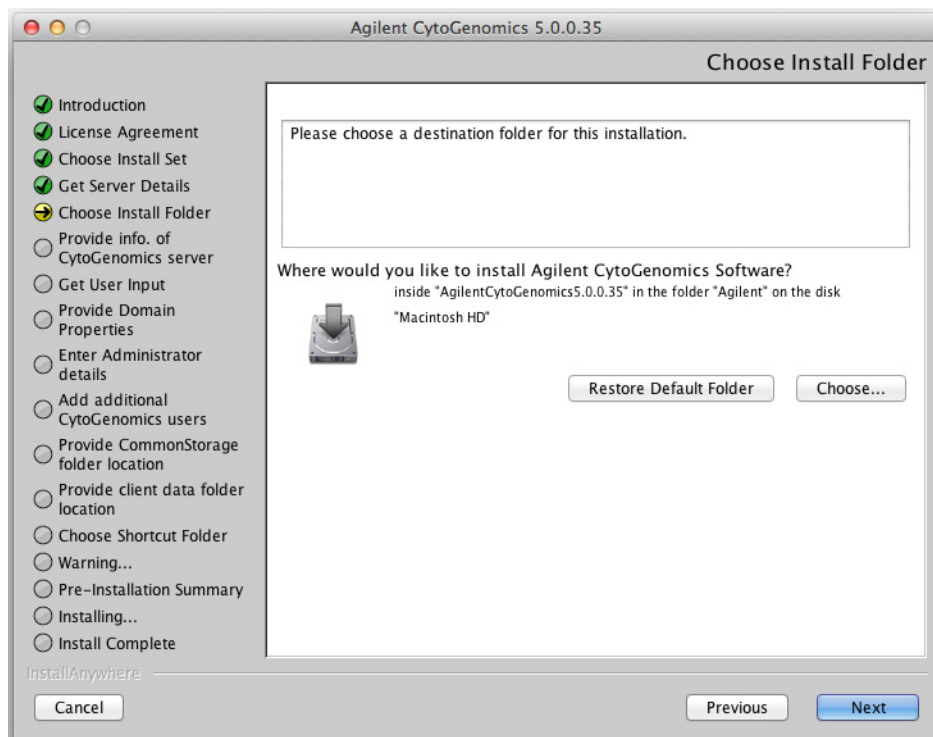


Figure 46 Choose Install Folder screen – Macintosh

12 By default, the program is installed in the AgilentCytoGenomics5.0.X.X folder (where X.X is the revision number) in the Agilent folder on the Macintosh HD disk. Leave this folder as-is, or click **Choose** to select a new location where you want to install Agilent CytoGenomics 5.0 folder. Within this folder, one folder is created for the server software and one folder is created for the client software.

13 Click **Next**.

The Provide Domain Properties screen opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

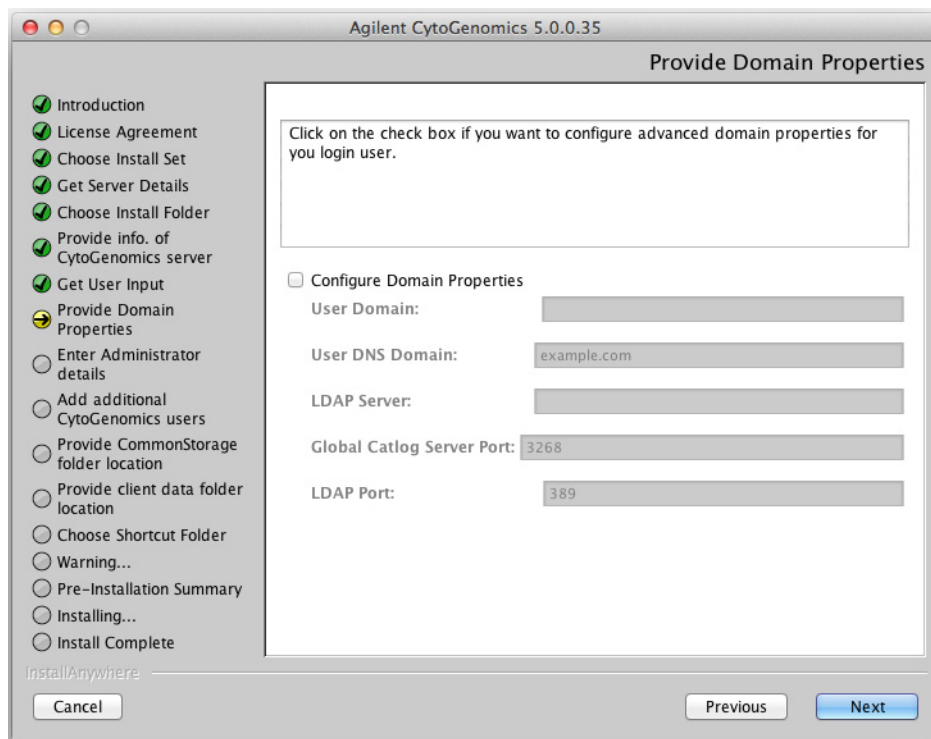


Figure 47 Provide Domain Properties screen – Macintosh

14 If only the administrator will be accessing Agilent CytoGenomics 5.0 from this computer, click **Next**. The Enter Administrator details screen opens, and you can proceed to [step 26](#).

If more than one user plans to use Agilent CytoGenomics 5.0 on this computer, configure the domain properties using the instructions in the following substeps.

- a** Mark **Configure Domain Properties**.
- b** Using the domain properties for your system, complete the fields for User Domain, User DNS Domain, LDAP Server, Global Catalog Server Port, and LDAP Port.
- c** Click **Next**.
The Enter Administrator details screen opens.

NOTE

If you see a warning message titled **Domain validation failed**, click **OK** in the message box, and then click **Previous** in the install window to return to the Provide Domain Properties screen. Repeat [step b](#) above with the correct domain properties. Do not proceed with the installation until you enter the domain properties and do not receive the **Domain validation failed** warning message.

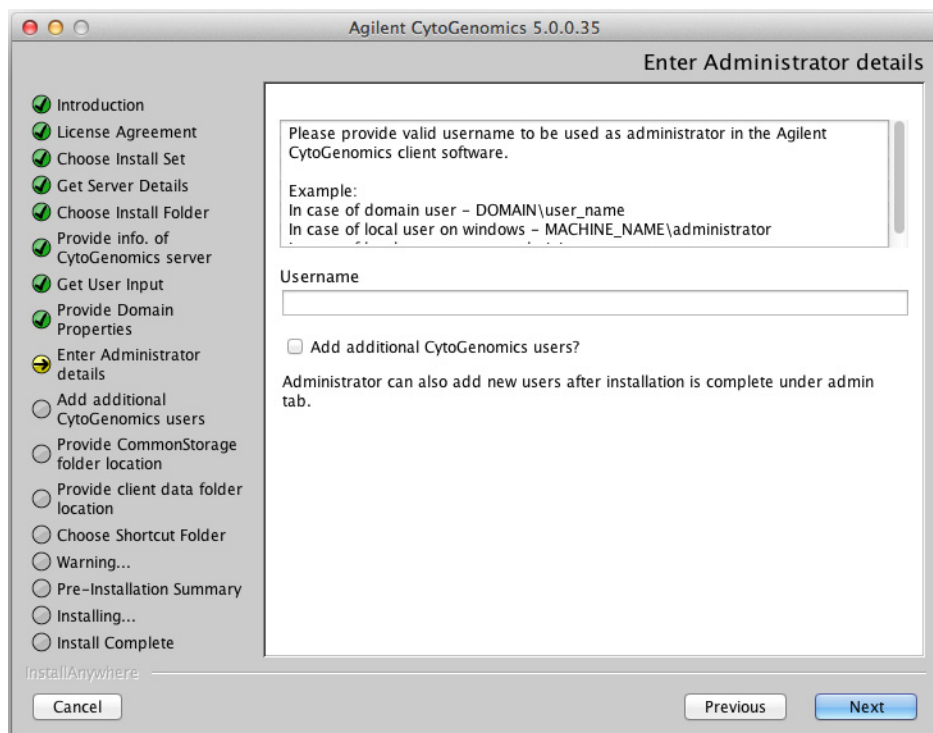


Figure 48 Enter Administrator details screen – Macintosh

15 In the Username field, type the user name of the person who will be the administrator for the Agilent CytoGenomics 5.0 software, using the format example shown.

For domain user: DOMAIN\username

For local machine: \username

16 To set up additional CytoGenomics users, mark the check box labeled **Add CytoGenomics users?**. In the Password field that appears, type the password for the administrator account.

If you do not want to set up new users at this time, proceed directly to [step 21](#) on [page 77](#). Note that the administrator can set up new users from within CytoGenomics once installation is complete (see “[To manage users and roles](#)” on [page 110](#)).

17 Click **Next**.

The Add CytoGenomics Users screen opens.

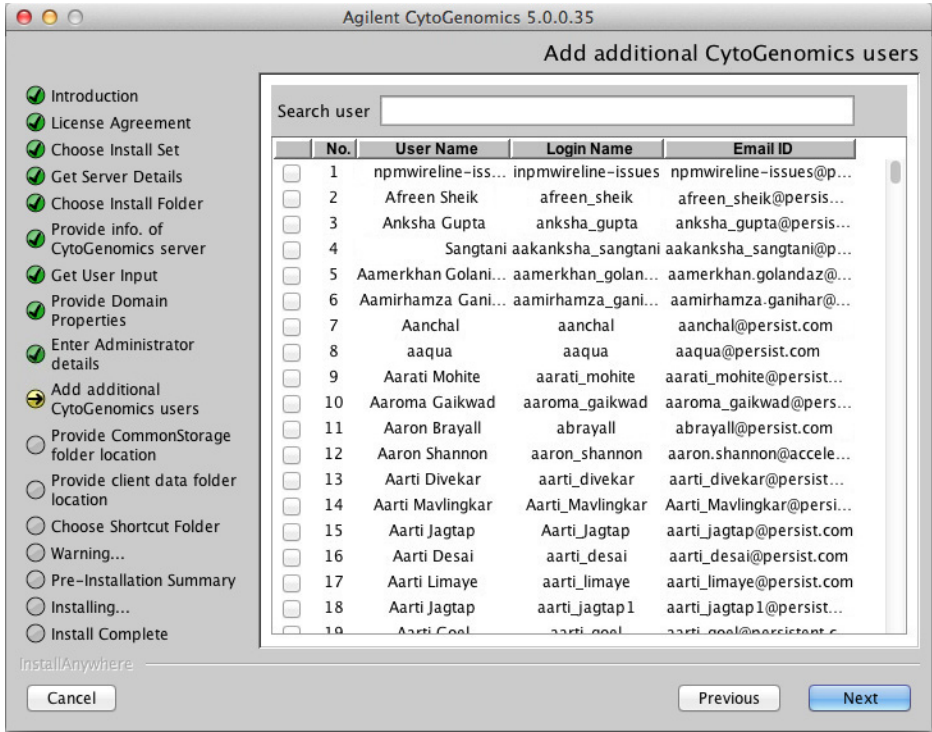


Figure 49 Add CytoGenomics Users screen – Macintosh

18 The Add CytoGenomics Users screen lists all of the users in the domain or on the local machine. Mark the check box next to the users that you want to add to the CytoGenomics database.

To search for a user, type the user name into the search box at the top of the screen.

19 Click Next.

The Add CytoGenomics Users screen advances to the step for configuring user roles.

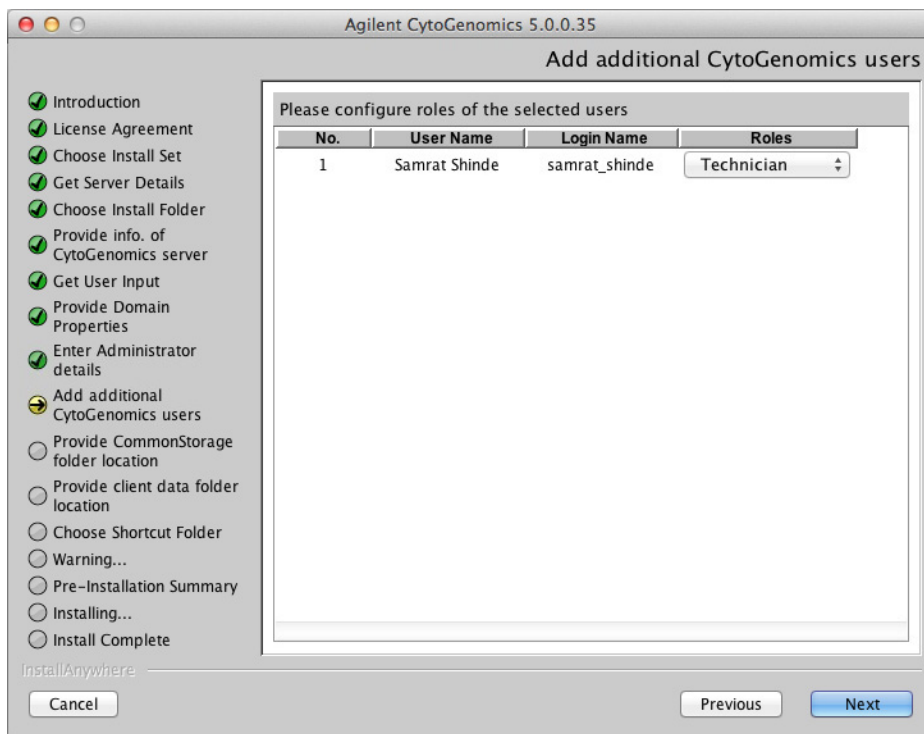


Figure 50 Add CytoGenomics Users screen – Configure roles step – Macintosh

20 For each user listed on the screen, select a role in the Roles drop-down list. The default selection is *Technician*.

See “[User Roles and Capabilities](#)” on page 16 for a description of each role.

21 Click Next.

The Choose common storage folder screen opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

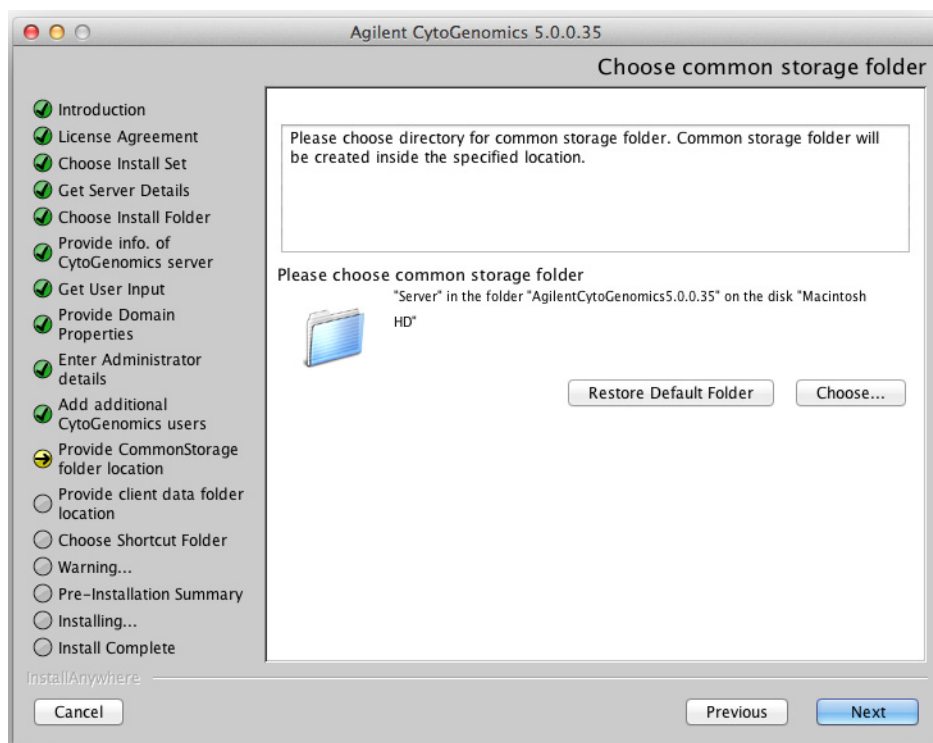


Figure 51 Choose common storage folder screen – Macintosh

22 By default, the common storage folder is created inside the Server folder in the AgilentCytoGenomics5.0.X.X folder (where X.X is the revision number) on the Macintosh HD disk. Leave this folder as-is, or Click **Choose**, browse to the desired folder, and then click **Save**.

The common storage folder is where the program saves files that are used at the server level.

23 Click **Next**.

The Choose client data folder screen opens.

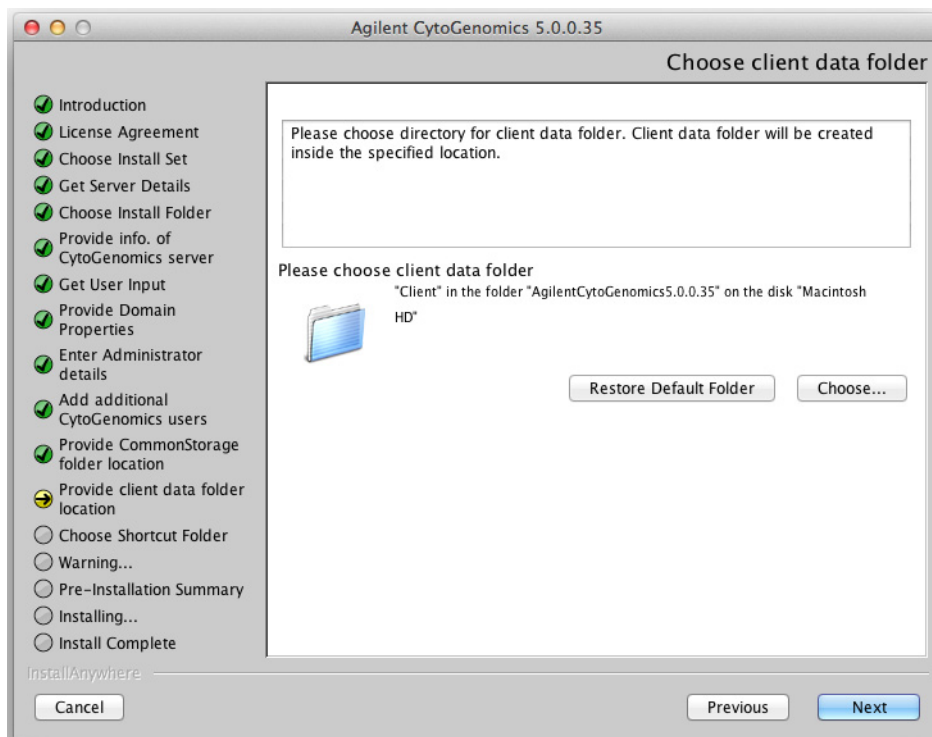


Figure 52 Choose client data folder screen – Macintosh

24 By default, the client data folder is created inside the Client folder in the AgilentCytoGenomics5.0.X.X folder (where X.X is the revision number) on the Macintosh HD disk. Leave this folder as-is, or Click **Choose**, browse to the desired folder, and then click **Save**.

The client data folder is where the program installs the files for the genome build and reference genomes.

25 Click **Next**.

The Choose Alias Folder screen opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

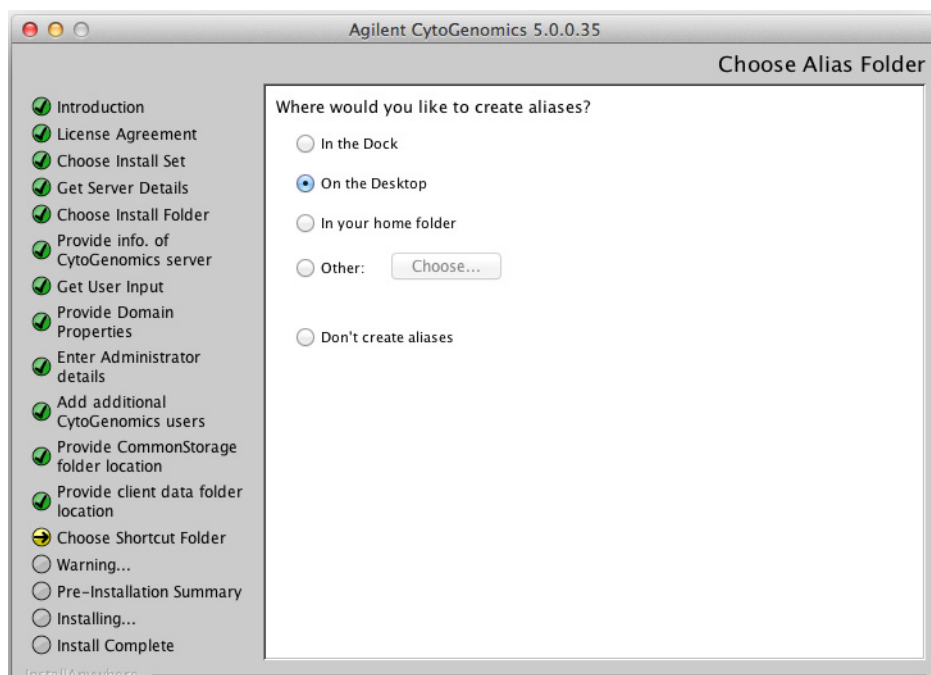


Figure 53 Choose Alias Folder screen – Macintosh

26 By default, program icons are installed on the Desktop of your computer. To select another location for the program icons, select the desired location (In the Dock, for example) and then click **Next**.

The Warning screen opens.

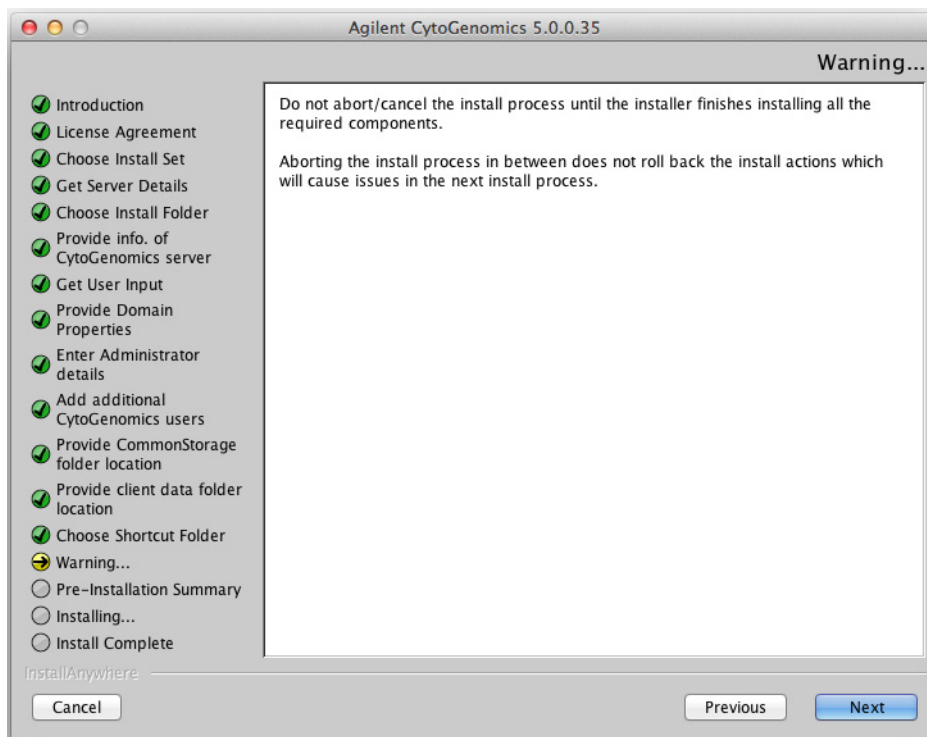


Figure 54 Warning screen – Macintosh

27 Read the message that cautions you not to stop the installation once it has started, and then click **Next**.

The Pre-Installation Summary screen opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

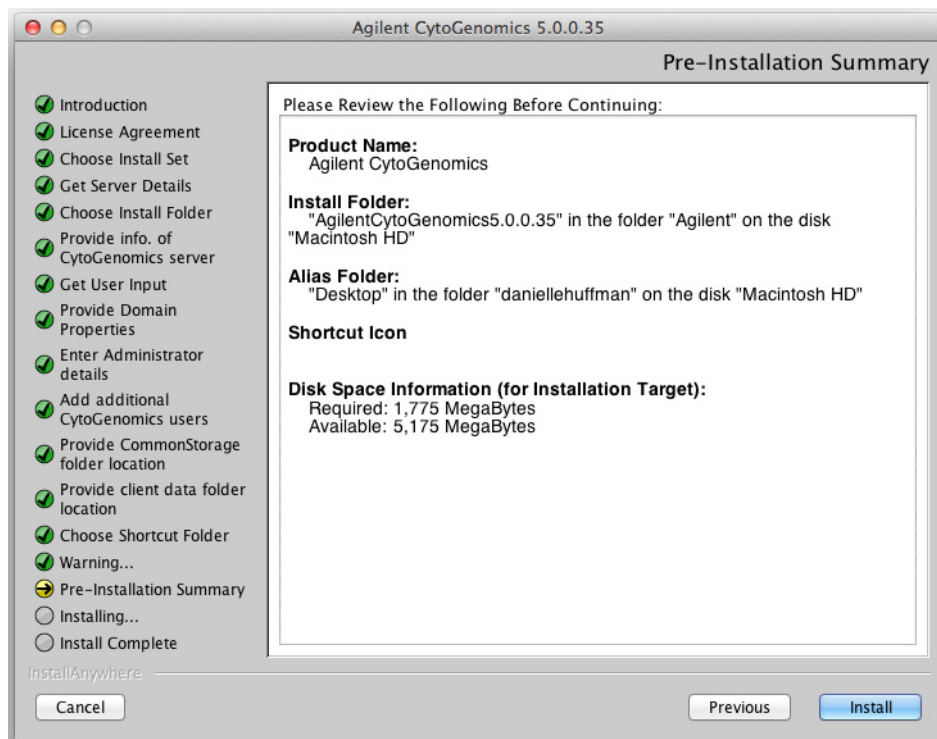


Figure 55 Pre-Installation Summary screen – Macintosh

28 Review the installation setup information. If you want to change any of the settings, click **Previous** and go back and change the settings. To start the installation, click **Install**.

The Installing Agilent CytoGenomics 5.0.X.X screen opens and is displayed until the installation is complete.

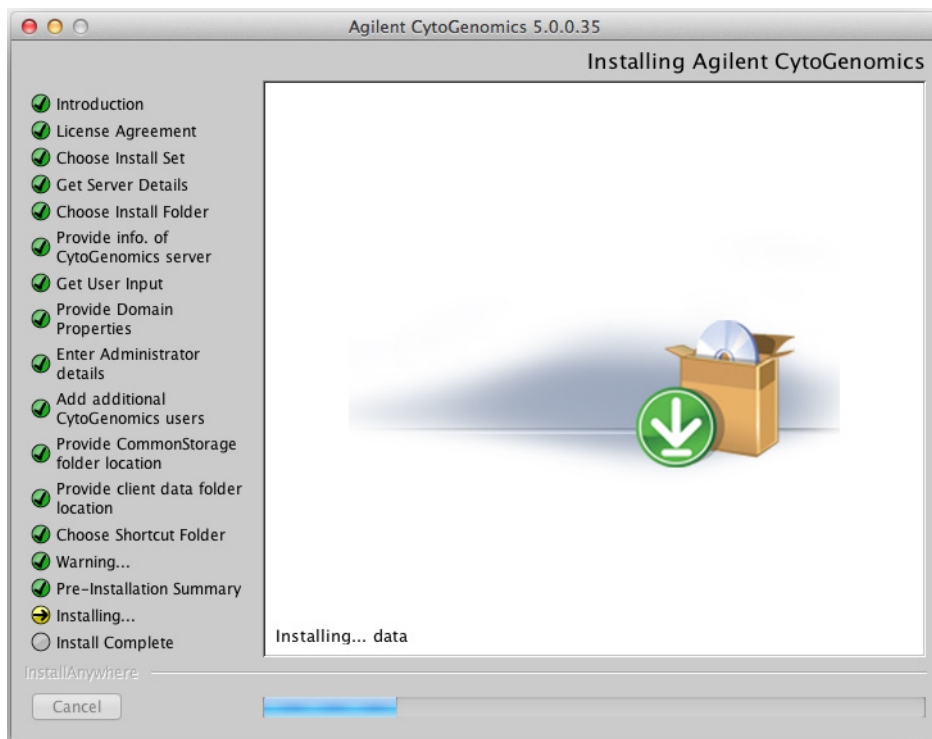


Figure 56 Installing Agilent CytoGenomics 5.0.X.X screen – Macintosh

When the installation is complete, the Installation Complete screen is displayed.

29 Click **Done**.

30 Make sure the folder where you installed the software is shared with read/write permission for all client computers and for all users who run Agilent CytoGenomics.

Step 3. Install the client software on additional Macintosh computers

To use other computers to run Agilent CytoGenomics 5.0 client software (with a shared database), install the client software on each of the computers.

- 1 On each computer you want to run the Agilent CytoGenomics 5.0 program, run the installation wizard, as described in “[Step 2. Install the database server and client on a single Macintosh computer](#)” on page 68.
- 2 When the Choose Install Set screen appears, select **Only Client**.

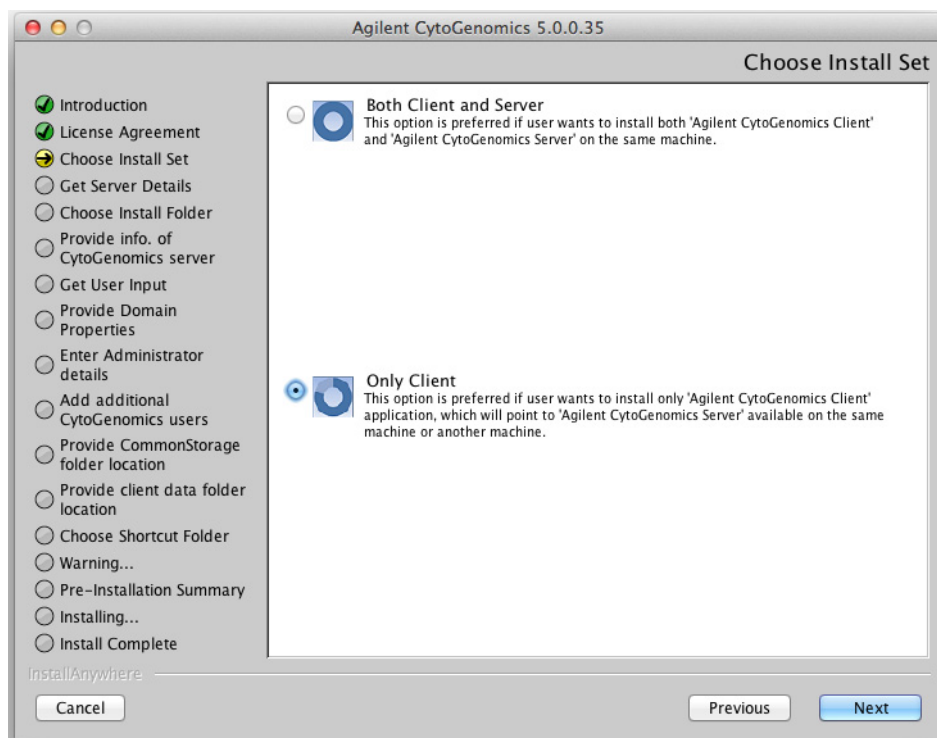


Figure 57 Choose Install Set screen – Macintosh

- 3 Click **Next**.

The Choose Installation Folder opens.

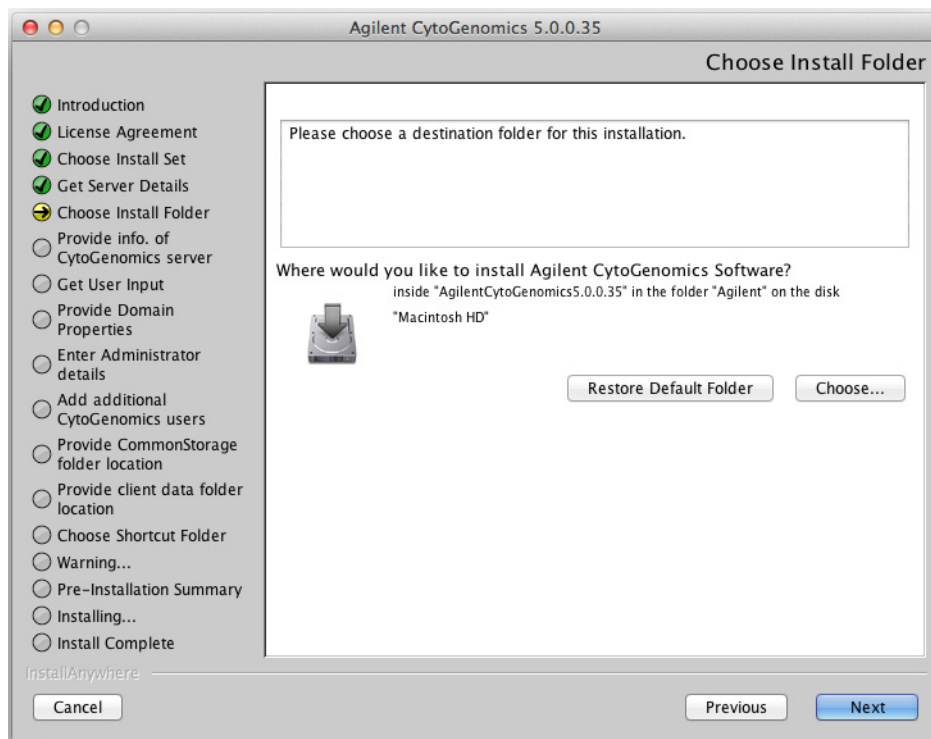


Figure 58 Choose Install Folder screen – Macintosh

- 4 By default, the program is installed in the AgilentCytoGenomics5.0.X.X folder (where X.X is the revision number) in the Agilent folder on the Macintosh HD disk. Leave this folder as-is, or type the location and name of the folder where you want to install Agilent CytoGenomics 5.0.

OR

Click **Choose** and browse to a location where you want to install the program, and then click **Save**.

- 5 Click **Next**.

The Provide Information of CytoGenomics server screen opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

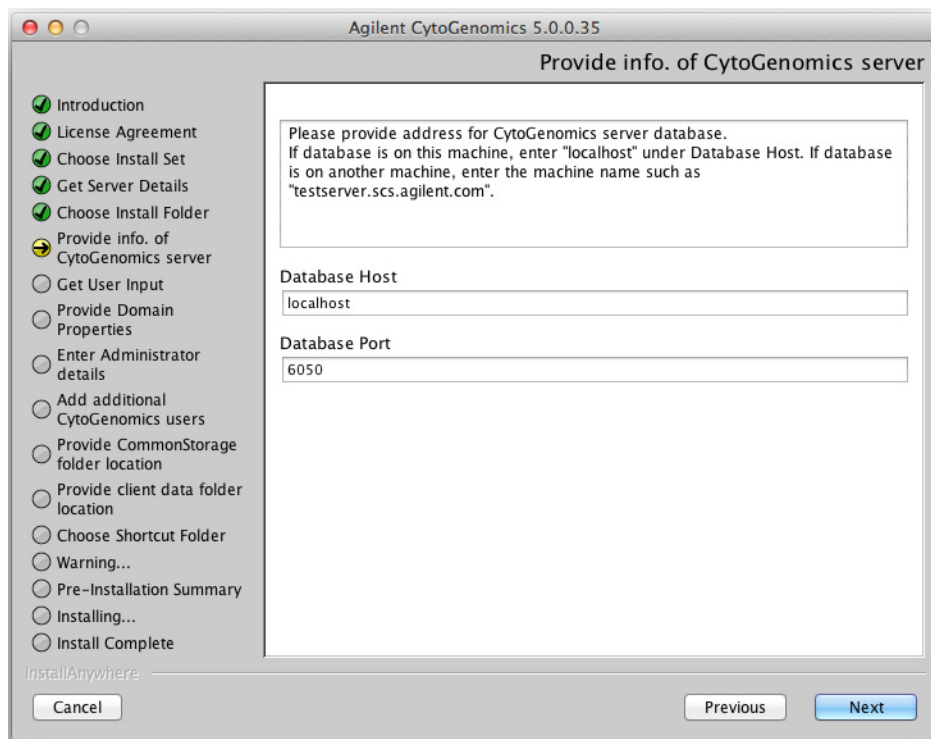


Figure 59 Provide information of CytoGenomics server screen – Macintosh

- 6 Under Database Host, type the assigned computer name or IP address of the computer (from [step 2](#)) where you installed the Agilent CytoGenomics 5.0 database server.
- 7 Accept the default **Database Port**, unless your system administrator tells you otherwise.
- 8 Click **Next**.

Installer immediately tests the connection to the specified database. If incorrect database parameters are provided, or the PostgreSQL database service is not running, a message is displayed.

You can choose to complete the installation and change database connection parameters when launching the client software. However, Agilent recommends that you provide the correct database information during installation.

Continue the installation wizard. Follow the instructions for the installation wizard to provide domain properties, enter administrator information, choose an alias location, and start the installation. For information, see [step 15](#) through [step 30](#) in the section “[Step 2. Install the database server and client on a single Macintosh computer](#)” on page 68.

Step 4. Start the Agilent CytoGenomics 5.0 client software

After you install the Agilent CytoGenomics database server and client programs as described in this manual, start the program.

- Double-click the Agilent CytoGenomics 5.0 icon to open the program.

NOTE

How you start the program depends on the choice you made for the Alias or Install Folder during installation of the Agilent CytoGenomics client.

Step 5. Log in to Agilent CytoGenomics

After you start the program, the Login screen appears. The name and domain for the administrator set during installation appears in the dialog box.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

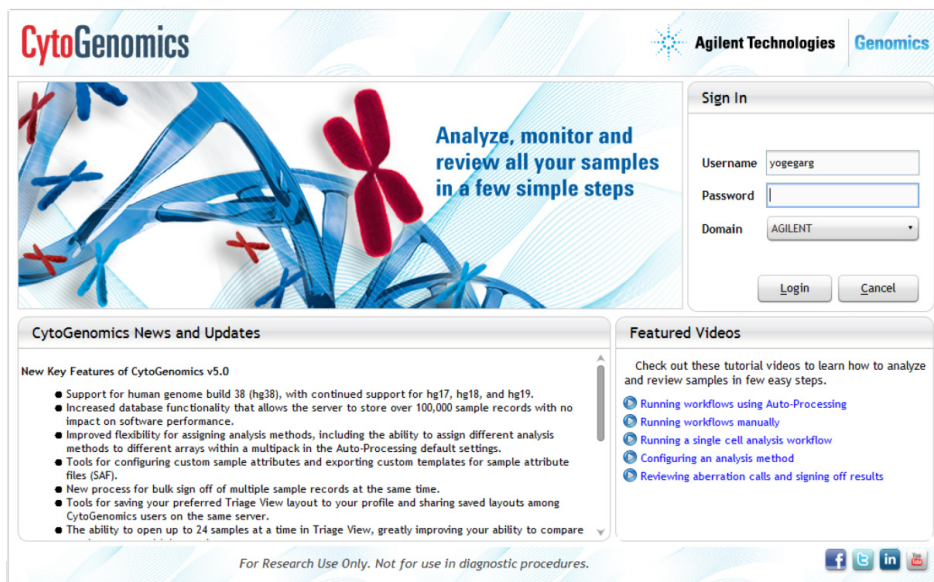


Figure 60 Login screen

- 1 Log in with your username and password. After the software is started, administrators can add additional users and assign their roles.
- 2 Next to **Password** type your password.
- 3 To change the domain, click the arrow next to **Domain** and select the correct domain. Otherwise, leave it as-is.
- 4 Click **Ok**.

NOTE

If you are using a local machine that is not part of a domain, you see a message that the software cannot find the domain. Click **OK** to ignore this message.

If the installer did not properly share the common storage folder during installation, then an Application Initialization Error message will open notifying you that you do not have access to the common storage folder. This error message provides instructions on how to share the folder with all users. Follow the steps in the error message, then click **OK** to close the message and restart the application.

Step 6. Add your software license

After you log in for the first time, the following dialog box opens.

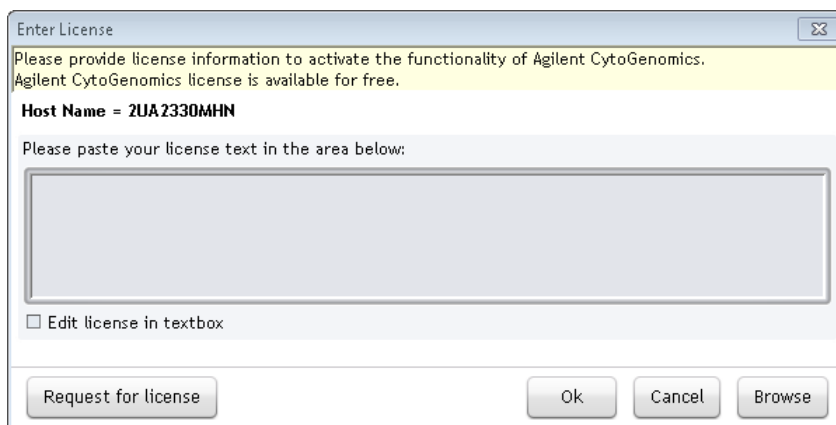


Figure 61 Enter License dialog box

- 1 If you do not already have the license file saved to your computer, click **Request for license** to open the Agilent website to the CytoGenomics license request page. Complete the fields in the request form to obtain a CytoGenomics software license. The license is contained in a text file and sent by email.
- 2 At the bottom of the Enter License dialog box, click **Browse**.
The Open dialog box opens.
- 3 Browse to the saved license text file. Select the file and click **Open**.
The program loads the contents of the file into the text box on the Enter License dialog box.
- 4 Click **OK**.
The Agilent CytoGenomics program opens.

3 Installation Instructions for Macintosh

Installing Agilent CytoGenomics 5.0 – Macintosh

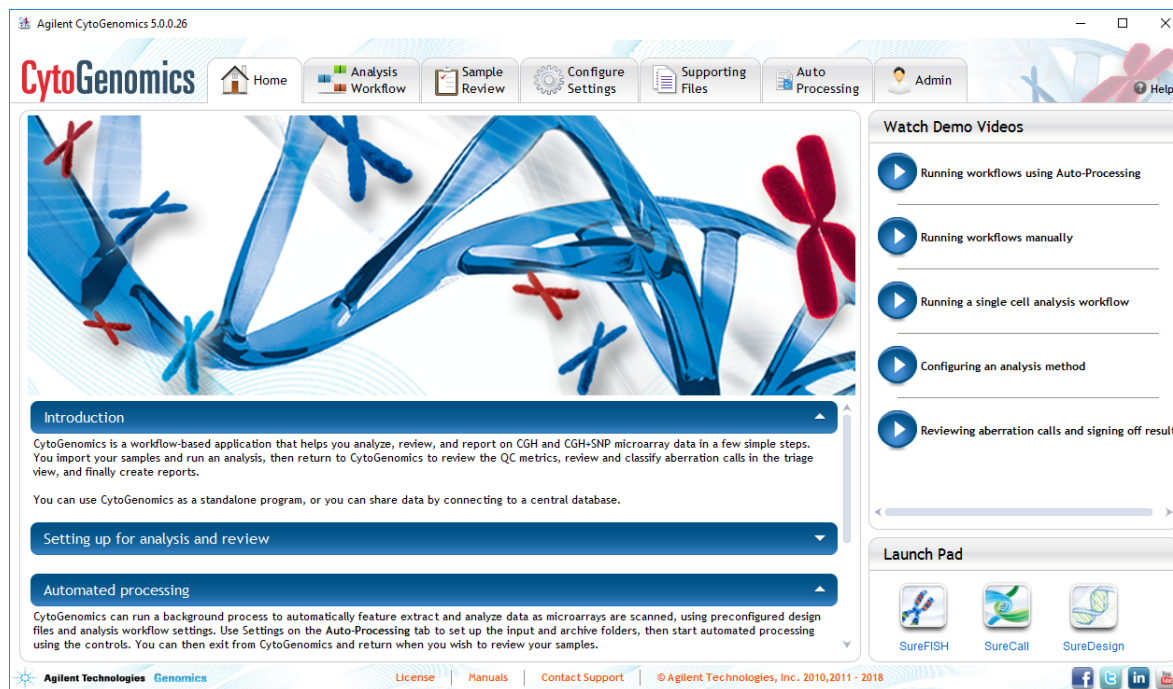


Figure 62 Agilent CytoGenomics program Home screen

Step 7. Add users and assign roles

Before other users can log in to Agilent CytoGenomics, the administrator must add the users and assign user roles. The user role determines what capabilities the user has within Agilent CytoGenomics. For information on what capabilities each user role has, see [“User Roles and Capabilities”](#) on page 16. For information on how to add users and assign roles, see [“To manage users and roles”](#) on page 110.

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh

These instructions are for users of a previous version of CytoGenomics (specifically, version 3.0 or 4.0) who want to upgrade to CytoGenomics 5.0 and migrate all sample records from the previous database to the new CytoGenomics 5.0 database.

NOTE

CytoGenomics 5.0 is compatible with and can co-exist on the same machine as CytoGenomics 3.0 or 4.0. It is not necessary to uninstall the previous version of CytoGenomics (i.e., v3.0 or 4.0) in order to install v5.0.

Step 1. Install CytoGenomics 5.0 on the server computer

Follow these steps to install the CytoGenomics 5.0 server and client software.

- 1 Download the new software revision from Agilent Technologies and check the system requirements.
See “a. [Download the Macintosh version of the software and release notes](#)” on page 66 for detailed instructions on this step.
- 2 Start the installation program you downloaded from Agilent.
- 3 Read the Introduction and then click **Next**.
- 4 Click **Accept** and then click **Next** to accept the license agreement.
- 5 On the Choose Install Set screen, select **Both Client and Server**.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh

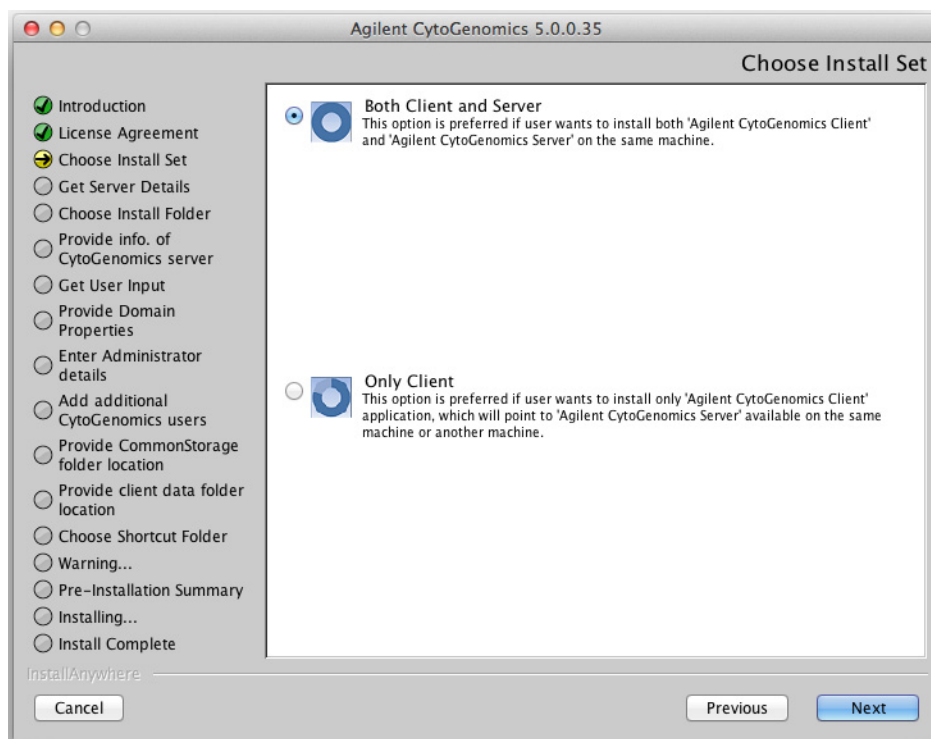


Figure 63 Choose Install Set screen – Macintosh - select **Both Client and Server**

6 Click **Next**.

7 Continue the installation wizard. Follow the instructions to enter administrator information, choose a shortcut location, and start the installation. For more information, see [step 12](#) through [step 29](#) in the section “[Step 2. Install the database server and client on a single Macintosh computer](#)” on page 68.

At the completion of installation, the Launch Migration Utility dialog box opens asking if you want to launch Agilent’s migration utility tool designed to transfer data from a previous version of CytoGenomics server to the CytoGenomics 5.0 server.

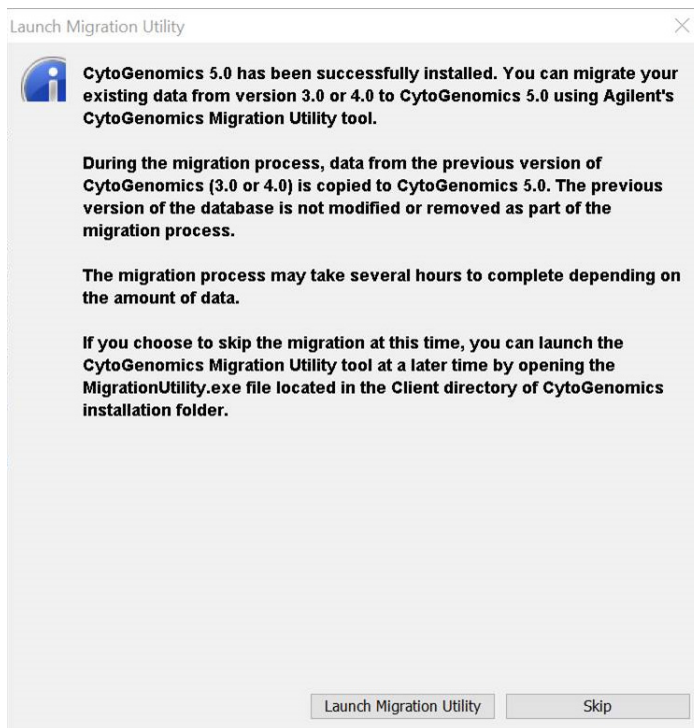


Figure 64 Launch Migration Utility dialog box – Macintosh

8 Depending on your needs, click either **Launch Migration Utility** or **Skip**.

- Click **Launch Migration Utility** if you have data saved to the previous version of CytoGenomics server (e.g., sample results, track files, and other supporting files) and you want to transfer that data to the CytoGenomics 5.0 server. The migration utility tool launches on your PC. Proceed to [“Step 2. Migrate data to CytoGenomics 5.0 server”](#) on page 94.
- Click **Skip** if you do not want to transfer any data (e.g., sample results, track files, and other supporting files) from the previous version of CytoGenomics server to the CytoGenomics 5.0 server at this time. The dialog box will close and that data will remain on the previous version of the server. Proceed to [“Step 3. Upgrade to CytoGenomics 5.0 on the client computers”](#) on page 98.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh

NOTE

To protect against loss of data and ensure migration of all analyzed samples, Agilent strongly recommends migrating data immediately upon completion of the CytoGenomics 5.0 installation.

If you click **Skip** with the intention of migrating your data at a later time, be aware that if CytoGenomics 5.0 already contains data at the time of migration, the Migration Utility application will delete that data and overwrite it with data from the previous version of CytoGenomics.

Step 2. Migrate data to CytoGenomics 5.0 server

If you selected to launch the migration utility tool, follow these step to use the tool to transfer data from the previous version of CytoGenomics server to CytoGenomics server 5.0. If you do not want to transfer data, bypass this step and proceed to “[Step 3. Upgrade to CytoGenomics 5.0 on the client computers](#)” on page 98.

- 1 On the Sign In screen of the migration utility application ([Figure 65](#)), review the list of migration prerequisites and make sure they are all met before you continue.

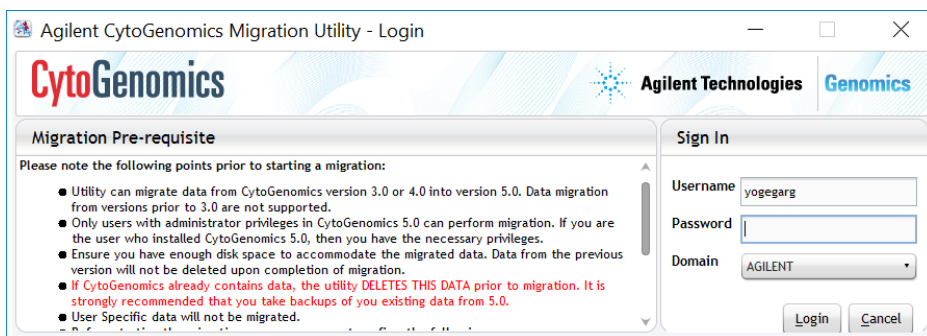


Figure 65 CytoGenomics Migration Utility application login screen

- 2 Log in with your username and password. The username and domain for the administrator set during installation appears in the Sign In area of the screen. (Only accounts with administrator privileges can migrate data.)

The CytoGenomics Migration Utility application opens.

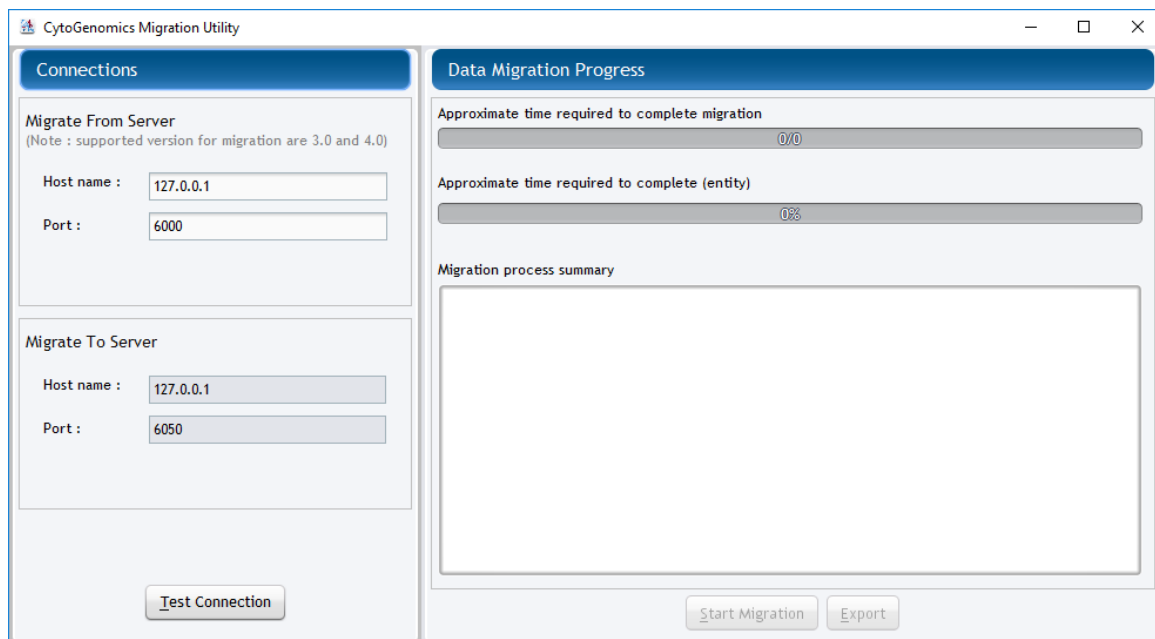


Figure 66 CytoGenomics Migration Utility application

- 3 Under **Migrate From Server**, enter the host name (i.e., IP address) and port number where the previous version of CytoGenomics server is installed.
- 4 Under **Migrate To Server**, enter the host name (i.e., IP address) and port number where the CytoGenomics 5.0 server is installed. If CytoGenomics 5.0 server is installed on the same PC as the previous version of CytoGenomics server, then the host names are the same.

NOTE

If CytoGenomics 5.0 server is installed on the same PC as the previous version of CytoGenomics server, then the host names of the two servers are the same.

- 5 Click **Test Connection**.

The screen displays the results of connection test with each server.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh

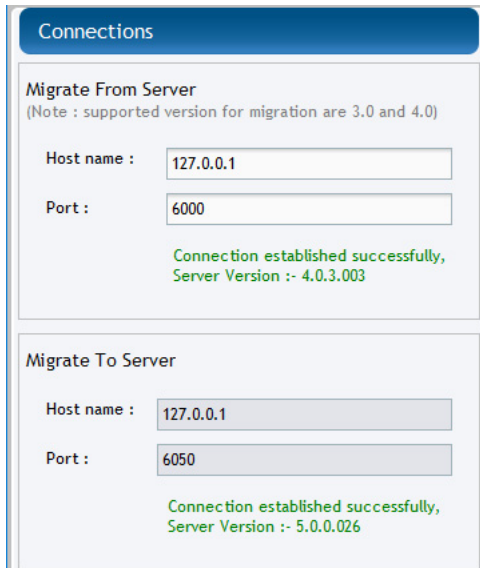


Figure 67 CytoGenomics Migration Utility application – Connections established

NOTE

If the application was unable to successfully establish a connection to one or both servers, verify that the correct host name and port number are entered and test the connection again. You cannot proceed to data migration until the application connects to the servers.

- 6 Once the application has successfully connected to the servers, click **Start Migration**.

The Warning message box opens.

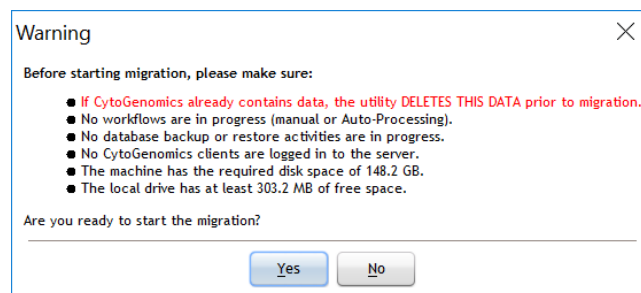


Figure 68 Warning message box

- 7 As instructed in the Warning message box, make sure that CytoGenomics is not actively running any workflows or performing any database backup or restore activities and that all clients are logged out. If your CytoGenomics 5.0 database already contains data, make sure that data has been backed up. Then, click **Yes** to continue with the migration.

The right side of the screen (under **Data Migration Process**) displays the progress of the migration and the migration process summary.

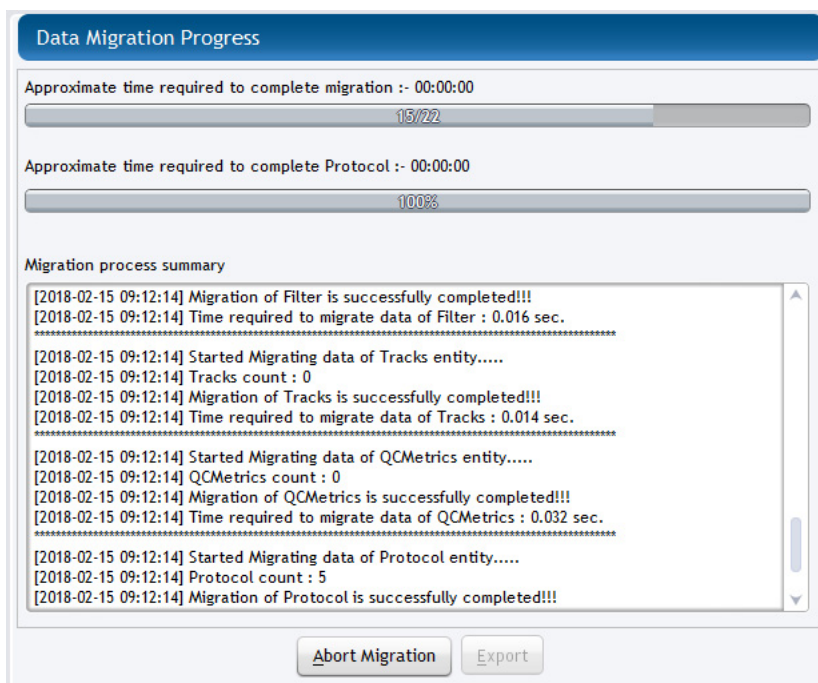


Figure 69 CytoGenomics Migration Utility application – Migration in progress

Upon completion of the migration process, the following message box opens.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 3.0 or 4.0 to CytoGenomics 5.0 and Migrating Data – Macintosh

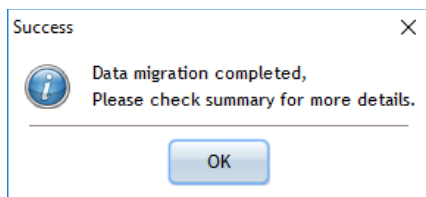


Figure 70 CytoGenomics Migration Utility application – Data migration complete

- 8 Click **OK** to close the message box.
- 9 If desired, click **Export** to export a summary of the migration job.
- 10 Click the X in the top right corner of the Migration Utility application window to close the application.

Step 3. Upgrade to CytoGenomics 5.0 on the client computers

Follow these steps to install the CytoGenomics 5.0 client software.

- This procedure is the same as installing on client computers for a new installation.

For more information, see [“Step 3. Install the client software on additional Macintosh computers”](#) on page 84.

Step 4. Start the Agilent CytoGenomics 5.0 client software, log in, and add new software license

- 1 Start the Agilent CytoGenomics program.
See [“Step 4. Start the Agilent CytoGenomics 5.0 client software”](#) on page 87.
- 2 Log in.
See [“Step 5. Log in to Agilent CytoGenomics”](#) on page 87.
- 3 Enter your updated license information.
See [“Step 6. Add your software license”](#) on page 89.

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

If you are using CytoGenomics 5.0.0 or 5.0.1 and want to upgrade to CytoGenomics 5.0.2, you only need to upgrade the CytoGenomics client application. You do not need to upgrade to the CytoGenomics server application or migrate data from the previous version to the new version.

Step 1. Uninstall CytoGenomics 5.0.0 or 5.0.1 client

- 1 Double-click the Uninstaller file, located in the \Uninstall_Agilent CytoGenomics folder of your program folder.

The following window opens.

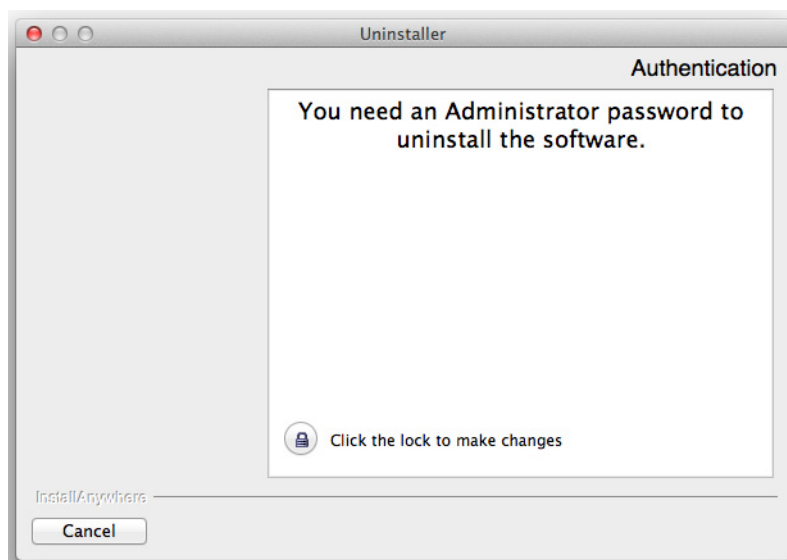


Figure 71 Authentication dialog box - preparing to uninstall

- 2 Click directly on the lock icon near bottom of the dialog box. The following dialog box opens.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

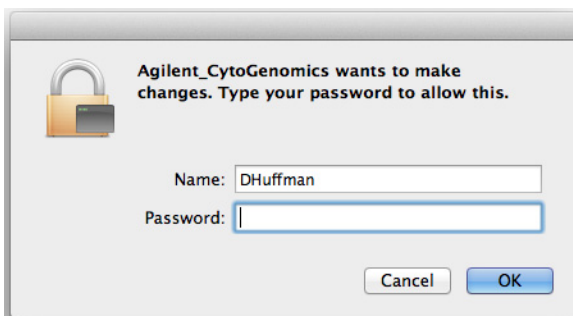


Figure 72 Administrator credentials dialog box - preparing to uninstall

- 3 In this dialog box, enter the Name and Password for a user with administrator rights to the computer, then click **OK**.
The uninstaller starts.
- 4 Read the introductory information, and then click **Next**.
- 5 Read the warning information. You are cautioned not to stop the uninstaller once it has started to remove the programs. Click **Next**.
The Uninstall Options screen opens.

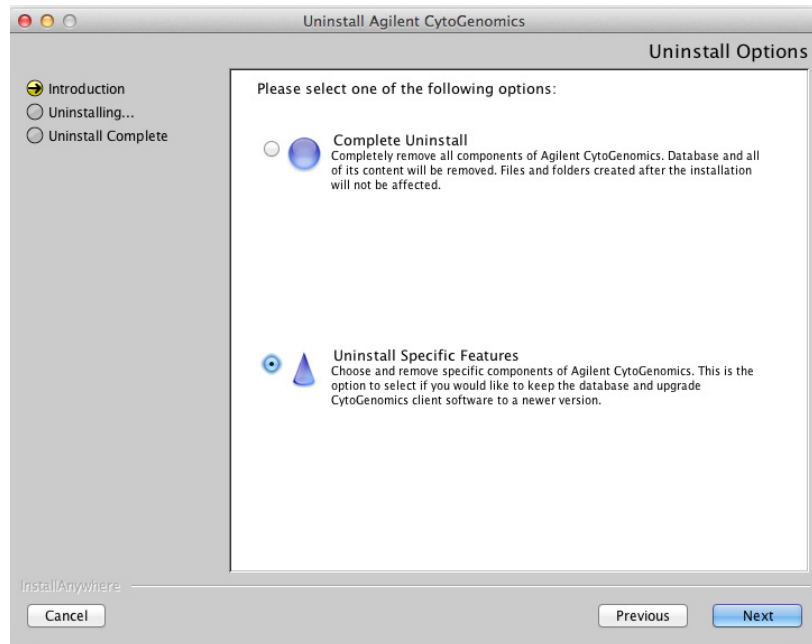


Figure 73 Uninstall Options screen – Macintosh – select **Uninstall Specific Features**

6 Select **Uninstall Specific Features** and click **Next**.

The Choose Product Features screen opens.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

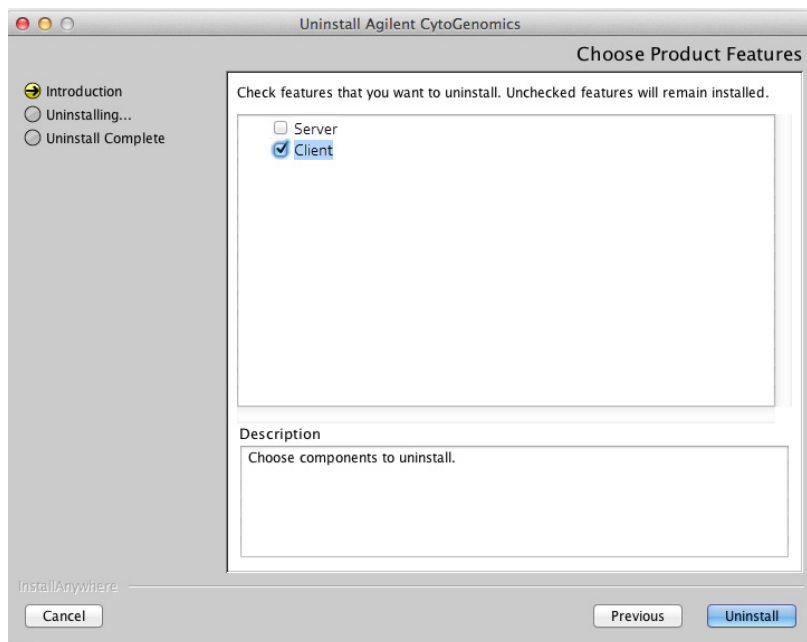


Figure 74 Choose Product Features screen – Windows – mark **Client**

7 Mark **Client** and click **Uninstall**.

NOTE

Do not mark **Server** on the Choose Product Features screen. Doing so will result in removal of all previously analyzed data from your existing CytoGenomics server.

- 8** In the message box that opens click **Yes** to confirm that you want to uninstall the client software.
- 9** When the uninstallation is complete, click **Done** to close the uninstaller.

Step 2. Install CytoGenomics 5.0.2 client

- 1** Download Agilent CytoGenomics 5.0.2 from Agilent Technologies and check the system requirements.
See [“a. Download the Macintosh version of the software and release notes”](#) on page 66 for detailed instructions on this step.

- 2 Double-click the compressed folder (Agilent_CytoGenomics_5_0_2_X_MAC.zip) that you downloaded and save the installer on your computer.
- 3 Double-click the Mac installer.
- 4 Run the installation wizard, as described in “Step 2. Install the database server and client on a single Macintosh computer” on page 68.
- 5 When the Choose Install Set screen opens, select **Only Client**.

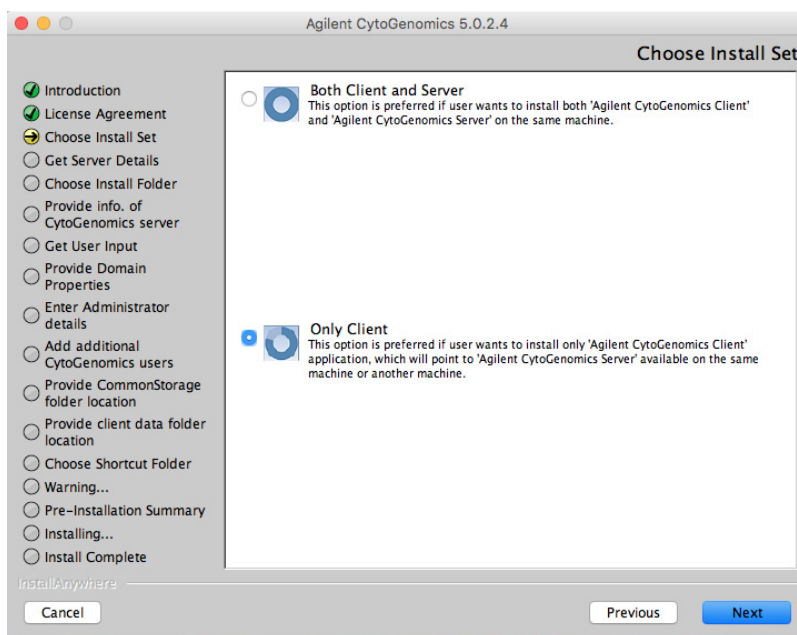


Figure 75 Choose Install Set screen – Windows – select **Only Client**

- 6 Click **Next**.
- 7 Continue the installation wizard. For information, refer to “Step 3. Install the client software on additional Macintosh computers” on page 84.

3 Installation Instructions for Macintosh

Upgrading from CytoGenomics 5.0.0 or 5.0.1 to CytoGenomics 5.0.2

Step 3. Start the Agilent CytoGenomics 5.0.2 client software, log in, and add new software license

- 1** Start the Agilent CytoGenomics program.

See “[Step 4. Start the Agilent CytoGenomics 5.0 client software](#)” on page 87.

- 2** Log in.

See “[Step 5. Log in to Agilent CytoGenomics](#)” on page 87.

- 3** Enter your updated license information.

See “[Step 6. Add your software license](#)” on page 89.

Uninstalling Agilent CytoGenomics on Macintosh computers

The Uninstall program for Agilent CytoGenomics provides an easy way to remove the CytoGenomics client, server, and PostgreSQL programs from your computer.

NOTE

CytoGenomics 5.0 is compatible with and can co-exist on the same machine as CytoGenomics 3.0 or 4.0. It is not necessary to uninstall the previous version of CytoGenomics (i.e., v3.0 or 4.0) in order to install v5.0.

When you completely uninstall Agilent CytoGenomics, the database and the data contained in it are completely removed. Do not uninstall the entire program unless you have backed up your database or no longer wish to use it.

- 1 Double-click the Uninstaller file, located in the **\Uninstall_Agilent CytoGenomics 3.0.X.X** or **\Uninstall_Agilent CytoGenomics** folder of your program folder.

The following window opens.

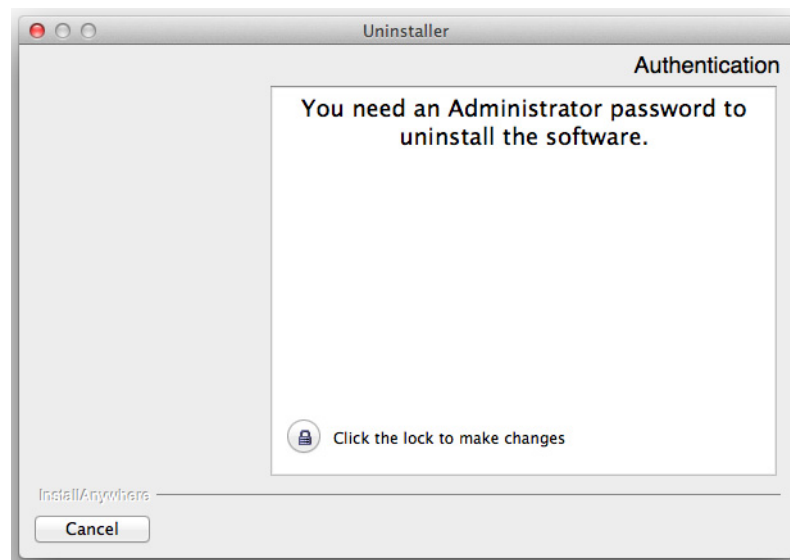


Figure 76 Authentication dialog box - preparing to uninstall

3 Installation Instructions for Macintosh

Uninstalling Agilent CytoGenomics on Macintosh computers

- 2 Click directly on the lock icon near bottom of the dialog box.
The following dialog box opens.

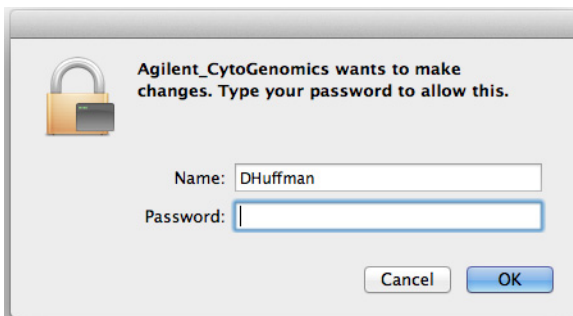


Figure 77 Administrator credentials dialog box - preparing to uninstall

- 3 In this dialog box, enter the Name and Password for a user with administrator rights to the computer, then click **OK**.
The uninstaller starts.
- 4 Read the introductory information, and then click **Next**.
- 5 Read the warning information. You are cautioned not to stop the uninstaller once it has started to remove the programs. Click **Next**.
The Uninstall Options screen opens.

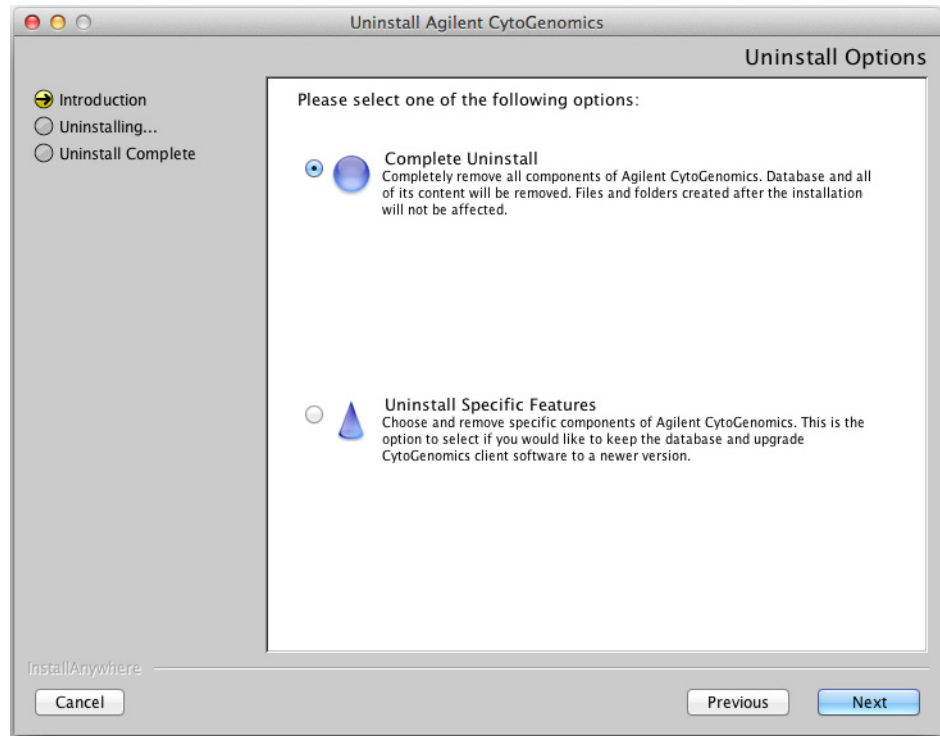


Figure 78 Uninstall Options screen – Macintosh – select **Complete Uninstall**

6 Select **Complete Uninstall**.

7 Click **Next**.

8 When the uninstallation is complete, a message appears. Click **Done** and then restart the system.

The Agilent CytoGenomics client program, server program and database, and PostgreSQL program were removed from your computer. Program folders and data folders (such as data output folders) created after installation were not removed.

3 Installation Instructions for Macintosh

Uninstalling Agilent CytoGenomics on Macintosh computers



4

Administering the System

Administering the System 110

To manage users and roles 110

To manage the database 111

This chapter shows you how to add users, assign roles and manage the database.

See the *Product Overview Guide* for an overview of the program and organization of the tasks within it.



Agilent Technologies

Administering the System

As an Administrator for Agilent CytoGenomics, you have access to everything. That is, you can perform any task available in the program. Your primary responsibility is to add users and their roles to the system, and you also manage the database.

See “User Roles and Capabilities” on page 16 for an explanation of the tasks that each role can perform in the program. Users who are assigned the role of Technician can only run workflows and view the results. They cannot set up workflows, nor can they sign off on results. Those are tasks for users assigned the role of Scientist.

To manage users and roles

Only the Administrator role can perform these tasks.

Table 4 Tasks for managing users and roles

To do this task	Follow these instructions	Comments
Add new users	1 Click Admin. 2 Click Manage Users & Roles. 3 Click Add New User. A list of all the users in the domain or on the local machine appears. 4 Next to Add New User, type the user name, and then click Find. (See comments for information about adding users for local machines.) 5 At the bottom of the Add User Dialog dialog box, click Add User. 6 When the Add User dialog box opens asking you to confirm, click Yes. 7 When the notice of success opens, click OK. 8 When you are finished adding users, click Close. The user now appears in the list of users with the role of Technician, with a status of <i>Enabled</i>.	• A user cannot be deleted but can be disabled. a Click Edit for the user you intend to disable. b Clear the Enabled check box, and click Save. • For local machines with no Domain, a Next to User, type the computer name and user in the format <code>machinename\username</code>, or <code>\username</code> for Macintosh systems. (Do NOT click Find.) b Click Add User. A warning dialog box opens stating that there was a problem fetching user details from the LDAP server. You can ignore this warning. Click OK to close the dialog box.

Table 4 Tasks for managing users and roles (continued)

To do this task	Follow these instructions	Comments
Change user roles	1 Click Admin. 2 Click Manage Users & Roles. 3 On the line for the user whose role you intend to change, click View/Edit. 4 Highlight the role on the right side of the panel, and then click < to remove the role on the right. 5 Highlight a role on the left side of the panel, and then click > to add the role to the right. 6 Click Save.	• See “User Roles and Capabilities” on page 16 for descriptions of each user role. • You can create custom roles that allow for a desired set of privileges. See the CytoGenomics help system for instructions.

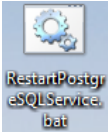
To manage the database

Although only the Administrator role can change the location of the database or the common storage folder, any user can restart the PostgreSQL service.

Table 5 Tasks for managing the database

To do this task	Follow these instructions	Comments
Change database locations and information	1 Click Admin. 2 Click Database Settings. 3 Under Database Settings, click Change. 4 Type: • Common Storage Location • Database Host • Database Port Click Apply to save your changes.	• If you change the location of your common storage, copy the contents of the previous location to the new location manually. Failure to do so can cause unexpected behavior in the program. • Enter the location of the common storage using UNC standard notation. (\\machine2\CommonStorage)

Table 5 Tasks for managing the database

To do this task	Follow these instructions	Comments
	If the PostgreSQL service stops, to restart it, click Restart PostgreSQL Service on the desktop. 	

For instructions on backing up and restoring the database, see the CytoGenomics help system.

www.agilent.com

In this book

This book provides an overview of Agilent CytoGenomics 5.0 and gives you instructions for installing and administering the software.

© Agilent Technologies, Inc. 2018

Revision C0, December 2018



G1662-90057



Agilent Technologies