



Agilent Genomic Workbench Feature Extraction 10.10

Quick Start Guide

Starting Feature Extraction on PCs with Agilent Scanners 2

Getting Help 6

What's new in version 10.10? 9

Starting the program and setting up eArray 13

Setting up a Standard Project to extract Agilent and non-Agilent image files 21

Setting up an On-Time project to extract Agilent image files in real time 47

This Quick Start Guide provides an overview of how to use the Feature Extraction 10.10 software, with step-by-step instructions on how to start the program and run an extraction. Exercises are also provided to help you, through the use of example images provided with the software.

This guide assumes you have already installed Feature Extraction 10.10 and its license on your computer. Feature Extraction 10.10 is installed automatically as part of the Agilent Genomic Workbench Standard Edition. If you need to install the software, see the installation guide for the Agilent Genomic Workbench Standard Edition.

For a complete description of the Feature Extraction main window and its components, see the *Feature Extraction 10.10 Users Guide*.

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



Agilent Technologies

Starting Feature Extraction on PCs with Agilent Scanners

Agilent Genomic Workbench software (including Feature Extraction) is pre-installed on PCs that are shipped with Agilent scanners. When you start the Feature Extraction or Agilent Genomic Workbench software for the first time on the PC that was shipped with your Agilent scanner, you will receive an error indicating that the software cannot find the database. Before you can use the software, you must set the database host name to the name of the PC.

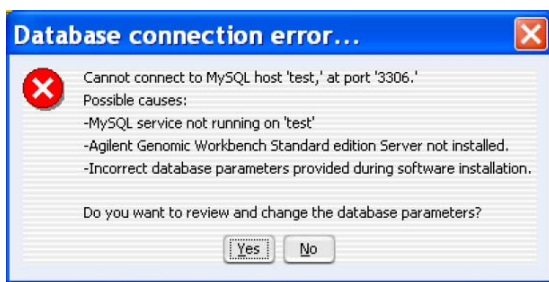


Figure 1 Database connection error for Agilent Genomic Workbench

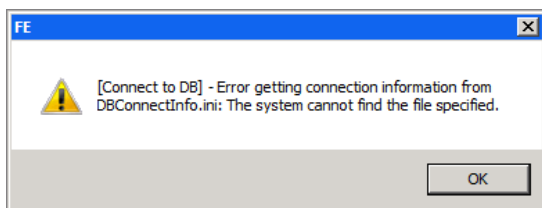


Figure 2 Database connection error for Feature Extraction

Use the following procedure to set the database host computer name.

To set the database host name for Agilent Genomic Workbench and Feature Extraction

- 1 Start the Agilent Genomic Workbench software.
- 2 When the Database connection error message appears, click **Yes**.

To set the database host name for Agilent Genomic Workbench and Feature Extraction

The Enter the database parameters dialog box appears.

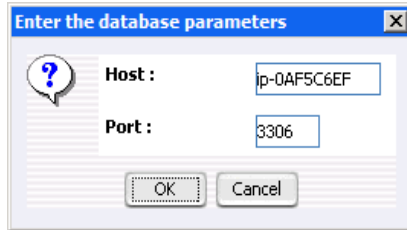


Figure 3 Enter the database parameters dialog box

- 3** In the box next to Host, type the name of your PC. You can find the name of the computer by opening the Windows Control Panel, and selecting System. Leave the Port set to 3306.
- 4** Click **OK** to close the dialog box.
- 5** You can now log into Agilent Genomic Workbench, or close it and start the Feature Extraction software.

What is Agilent Feature Extraction software?

Agilent Feature Extraction 10.10 software automatically reads and processes up to 100 raw microarray image files in the batch mode, or real-time extraction as soon as the .tif image is produced by the Agilent scanner. It finds and adjusts microarray grids, rejects outlier pixels, accurately calculates feature intensities and ratios, flags outlier features, and calculates statistical confidences. Application-specific QC reports summarize multiple analyses.

Feature Extraction is a key component of Agilent Genomics Workbench, a comprehensive informatics platform that integrates complementary technologies to provide a complete genomic solution, from design of microarray and target enrichment libraries through analysis and reporting.

Agilent Feature Extraction software extracts data from microarray images produced in two different situations:

- Agilent microarrays scanned on an Agilent scanner
- Non-Agilent microarrays scanned on an Agilent scanner

Feature Extraction handles each of these images differently. These differences are described throughout the documentation.

NOTE

Feature Extraction versions 10.7 and higher do not support microarray images generated by the GenePix/Axon scanner.

If you are a new user

Read [“Getting Help”](#) on page 6.

[“Starting the program and setting up eArray”](#) on page 13.

If you intend to analyze...

- Saved Agilent images, read the [“Instructional overview”](#) on page 21, and go to [“Example Exercise 1: Set up and run batch extractions on Agilent images”](#) on page 31.
- Saved non-Agilent images, read the [“Instructional overview”](#) on page 21, and go to [“Example Exercise 2: Set up and run batch extractions on non-Agilent images”](#) on page 43.

Agilent images in real time, read the “[Instructional overview](#)” on page 47, and go to “[General Instructions](#)” on page 48.

If you are an experienced user

Read “[What’s new in version 10.10?](#)” on page 9.

Do the exercises again if you want to sharpen your skills.

To manually download updates

When Feature Extraction is installed, Feature Extraction protocols and QC metric sets are also installed. Grid templates are not included with the installation. But if you set up an eArray login in Feature Extraction, the software will automatically download and update the necessary grid templates and Feature Extraction protocols from the eArray server. You can also go to the following web sites to download updates manually.

Listed below are some useful Agilent Web sites:

www.agilent.com/chem/earray - To register for eArray account

www.agilent.com/chem/feprotocols - To download Feature Extraction protocols

www.agilent.com/chem/designfile - To download grid templates

www.agilent.com/chem/feqcmetrics - To download QC metrics

www.agilent.com/chem/registration - To subscribe to E-Notes for monthly updates on Agilent products

Getting Help

To get help for Feature Extraction

The following tables list the resources available for getting help. Guides and manuals are opened using Adobe® Reader® software.

Help Resource	Description/Instructions
Feature Extraction Quick Start Guide	This guide, which you are reading, gives you an overview of the Feature Extraction software, including tutorial exercises. To open this guide from within Feature Extraction, click Help > Quick Start Guide .
Feature Extraction User Guide	A comprehensive guide for using the Feature Extraction software, that includes the following chapters: <ul style="list-style-type: none"> • Chapter 1– Working with Feature Extraction • Chapter 2– Extracting Microarrays Automatically • Chapter 3– Creating Grid Files and Templates • Chapter 4– Changing Protocol Settings • Chapter 5– Changing Image Displays To open this guide from within Feature Extraction, click Help > User Guide .
Feature Extraction Reference Guide	A reference guide that contains information about protocols, QC reports, results, and algorithms, that includes the following chapters: <ul style="list-style-type: none"> • Chapter 1 – Default Protocol Settings • Chapter 2 – QC Report Results • Chapter 3 – Text File Parameters and Results • Chapter 4 – MAGE-ML Results • Chapter 5 – How Algorithms Calculate Results • Chapter 6 – Command Line Feature Extraction To open this guide from within Feature Extraction, click Help > Reference Guide .
Technical Support	Opens a link to the Agilent Technical Support Web site where you can find literature and manuals, view Frequently Asked Questions, and download updates and revision information. There is also a Contact Us link where you can locate the technical support phone number or send an e-mail to technical support for your country.

Help Resource	Description/Instructions
Feature Extraction Training	<p>Opens a link to the Feature Extraction Tutorials Web Site, where you can view a number of self-paced instructional tutorials.</p> <p>To open this Web site from within Feature Extraction, click Help > Feature Extraction Training.</p>
Feature Extraction on the Web	<p>Opens a menu that provides links to the following:</p> <ul style="list-style-type: none"> • Array annotation – Find available design files by barcode • Feature Extraction Software – Feature Extraction software downloads and information • Feature Extraction Protocols – Download current and previous versions of Feature Extraction protocols
Agilent License	<p>Opens the Agilent software license redemption Web site, where you use your entitlement certificate to redeem your software license.</p>
Memory Status	<p>Opens an information window that shows how much computer memory is currently available and how much memory is in use.</p>
About	<p>Opens an information window that shows your Feature Extraction version, license, and database information. This window also displays your computer Host ID, which you need in order to redeem your license.</p>

To get help within Agilent Genomic Workbench

Help Resource	Description/Instructions
Feature Extraction Guides	<p>From the Help tab in Agilent Genomic Workbench, a Feature Extraction button appears if the selected application can use Feature Extraction. Click Feature Extraction to open a menu that lets you open the following guides:</p> <ul style="list-style-type: none"> • Feature Extraction Quick Start Guide • Feature Extraction User Guide • Feature Extraction Reference Guide

To learn about Agilent products and services

Help Resource	Description/Instructions
Other User Guides	<p>The Help tab in Agilent Genomic Workbench lets you view any of the available user guides that apply to the currently selected application type.</p> <ol style="list-style-type: none">1 Set the desired application type from the Switch Application menu.2 In the Agilent Genomic Workbench tab bar, click Help. The names of the available user guides appear in the command ribbon.3 Click the desired user guide. The selected user guide opens in Adobe Reader.

To learn about Agilent products and services

To view information about the Life Sciences and Chemical Analysis products and services that are available from Agilent, go to www.chem.agilent.com.

What's new in version 10.10?

Feature Extraction 10.10 includes the following new and changed functions:

- Perform Feature Extraction on CGH+SNP (Comparative Genomic Hybridization with Single Nucleotide Polymorphism) microarrays containing new SNP control probes.
- Import Dye Normalization lists from eArray and store them in the database.
- For manual gridding, associate Dye Normalization list (if present) automatically to grid file from the corresponding grid template.
- Use a Dye Normalization list associated with a grid template (if present) instead of the method specified in the protocol.
- Details of the Dye Normalization list used during extraction are included in the XML, QC Report, and Batch Summary output, and in the Run Monitor Window.
- External Dye Normalization lists are only used for non-Agilent arrays.
- Automatically create zip file of text output package.
- Changes to the Array Format in the Place Grid protocol step.
- Addition of two parameters to all array formats for the Place Grid protocol step.
- Ability to report population outliers as Failed in MAGEML file.
- The CGH report has been updated and renamed CGH_ChIP. It is now the default report for the ChIP protocol.
- New QC Report for Agilent test slides.
- Parameters (params) are now included in the output XML.
- New Corner Pattern Detection (CPD) protocol (available by request) for fully automatic grid placement of dim arrays.

See the Release Notes installed with the Feature Extraction software for more information about Feature Extraction 10.10 (**Start > Programs > Agilent > Agilent Genomic Workbench <version>\FeatureExtraction > Release Notes**).

Perform Feature Extraction on CGH+SNP (Comparative Genomic Hybridization with Single Nucleotide Polymorphism) microarrays containing new SNP control probes

The SurePrint G3 Human Comparative Genomic Hybridization (CGH) platform has been extended to include a set of Single Nucleotide Polymorphism (SNP) probes on the same microarray. This allows for the simultaneous detection of high resolution copy number and copy neutral variations and eliminates the need to run two separate microarrays. Agilent Feature Extraction software now supports feature extraction of new CGH+SNP microarrays that contain SNP control probes.

Import Dye Normalization lists from eArray and store them in the database

When you download a grid template from eArray.com Feature Extraction also downloads the corresponding DyeNormList file, if present. Feature Extraction automatically imports and stores the DyeNormList in the database and associates it with the grid template.

Associate Dye Normalization list automatically to grid file from the corresponding grid template

Grid files, used to help locate the spot centroids (geometric centers) on the microarrays, now are automatically associated with the dye normalization list (if present) from the grid template on which the grid file is based.

Use a Dye Normalization list associated with a grid template (if present) instead of the method specified in the protocol

You now have the option to use a dye normalization list associated with the grid template instead of the using the Dye Norm List setting from the protocol

Details of the Dye Normalization list used during extraction are included in the XML, QC Report, Batch Summary output, and in the Run Monitor Window

The name of the dye normalization list and the number of probes it contains are now displayed in the QC report header, Batch Summary Report, the XML output file, and in the Run Monitor Window.

External Dye Normalization lists are only used for non-Agilent arrays

External dye normalization lists are no longer used for Agilent arrays, which use the dye normalization lists associated with the array designs. You can specify an external dye normalization list for non-Agilent arrays, however.

Automatically create .zip file of text output package

You can now select to automatically create a .zip file containing the text output from an extraction set. This option is turned on or off in the Project Properties.

Changes to the Place Grid>Array Format choices

The 30 micron feature size choice was removed. Two new choices, 30 micron single pack, and 30 micron multi-pack are now available.

New options in the Place Grid step for all array formats

“Use central part of pack for peak and skew calculation” and “Use the correlation method to obtain origin x of subgrids” options are now available in the Place Grid step for all array formats. These options are useful particularly in cases where pack edges have dim spots and are failing to grid.

Report Population outliers as Failed in MAGEML file

With this new option turned on, any feature that is flagged as a population outlier is reported as Failed in the MAGE-ML output file.

The CGH report has been updated and renamed CGH_ChIP

The CGH report now includes a plot of the log ratio distributions, and has been renamed CGH_ChIP. This report is now the default report for the ChIP protocol.

New QC Report for Agilent test slides

A new QC Report, “Test Chip” is now available that is designed for use with Agilent test slides. This report is used by Agilent service personnel for scanner performance verification.

Parameters (params) are included in the output XML

The XML output file now contains Feature Extraction parameters.

New Corner Pattern Detection (CPD) protocol (available by request) for fully automatic grid placement of dim arrays

The Corner Pattern Detection (CPD) is a technique for fully automatic grid placement using a special pattern placed at the four corners of each array on a slide.

A template is formed by a unique arrangement of control probes. The template is printed and hybridized at four corners of each array along with other experimental probes. The grid placement step searches for these templates using image correlation technique. The locations obtained are used to calculate the origin, slope and skew. The grid geometry information is further used to place the grid on the scan.

Since the grid estimation is based on the locations obtained for the templates and the templates are placed at the corners of arrays, the irregularity in the spot grid due to missing rows/columns and the dust particles spread over image do not affect the grid placement. Further, the template size (10X10 spots) and the particular arrangement of spots within the template help to obtain correct locations even in the presence of noise and presence/absence of few probes and dust particles in the template area. Thus the technique is very robust to all common types of artifacts present in the slide image. In addition, the templates aid in accurate calculation of rotation and skew even if they are large in values. Above all, the technique has delivered very good grid irrespective of number of bright spots present on the array.

The prerequisite of the technique is presence of the corner patterns in the assays. For further information on this please contact Agilent Technical Support.

Starting the program and setting up eArray


The Feature Extraction program can be started in one of two ways:

- From Microsoft® Windows®
- From the Agilent Genomic Workbench Open Application tab

To start the program from Windows

- Double-click the **Feature Extraction** shortcut on your desktop.
This shortcut was created when the Feature Extraction program was installed.
OR
- From Windows, click **Start > All Programs > Agilent Genomic Workbench Standard Edition<revision> AGW Feature Extraction <revision>**.

To start the program from Agilent Genomic Workbench

- 1 Open Agilent Genomic Workbench, if you have not already done so.
- 2 In the Agilent Genomic Workbench main window, click the **Open Application** tab.
- 3 In the Open Application tab, click the Feature Extraction icon .

NOTE

You must have a valid license installed in order to use Feature Extraction. If you do not have a license installed, you will see an error message.

To set up a Feature Extraction license

Redeem your software licenses

To obtain your Feature Extraction license:

- 1 Start the Feature Extraction program and select **Help > Agilent License**.

Have the following items available:

- **Software Entitlement Certificate (SEC)** - This PDF contains the Order Number and Certificate Number that you need to provide in order to redeem your purchased license.
- **Host ID** - This can be found under the menu Help > About. You need to provide the host ID to redeem your purchased license.

- 2 Follow the instructions on the Agilent Software License Redemption Web site to redeem your purchased license.

After redeeming the license online, you will receive an e-mail with the license attached.


- 3 Save the unzipped license file to the folder on the drive where you installed the software. By default, Feature Extraction is installed in the Agilent Genomic Workbench installation folder:

C:\Program Files\Agilent\Agilent Genomic Workbench Standard Edition <version>\FeatureExtraction

Set up the license from Agilent Genomic Workbench

- 1 Locate your license file and double-click it to open the file in Notepad.
- 2 Highlight the license text, and copy it to the Clipboard.
- 3 Start the Agilent Genomic Workbench software. See [“To start the program from Agilent Genomic Workbench”](#) on page 13.
- 4 In the Open Application tab in Agilent Genomic Workbench, next to Feature Extraction, click **License**.
- 5 Click **License** and then paste the license in the blank license field. (To paste the text, press **Ctrl V** on your keyboard.)

Set up the license from Feature Extraction

- 1 Start the Feature Extraction software, if it is not already running.
If the Feature Extraction program is already running (unlicensed), click the **New Project** icon .
- 2 A message appears that informs you that you do not have a valid license installed. Click **Retry**.
The Select Agilent Feature Extraction License File dialog box opens.
- 3 Find the folder that contains the license file.
- 4 Select the license file (.lic), and click **Open**.

To set up eArray login

When you first open Feature Extraction, the **eArray Login Setting** window appears on top of the Feature Extraction main window. If you do not see the eArray Login Setting window, you can access it using Tools > eArray settings.

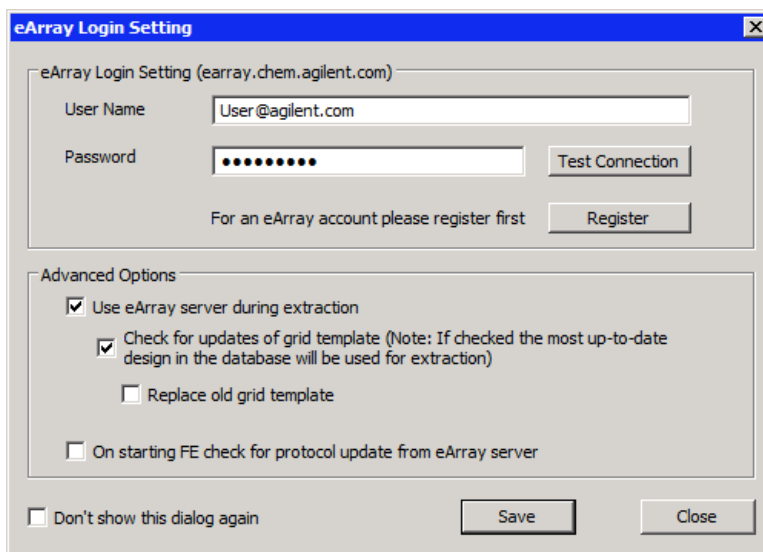


Figure 4 eArray Login Setting dialog box

To set up eArray login

Feature Extraction connects to the Agilent eArray server to automatically download, install, and use the required grid templates, Feature Extraction protocols, and QC metric sets during extraction.

For information on the eArray server and how to use the eArray_{XD} module in Agilent Genomic Workbench to create and manage custom microarray content for CGH, ChIP, methylation, gene expression, and microRNA applications, see the *eArray_{XD} User Guide*.

It is highly recommended that you set up an eArray login to take advantage of this feature *before* you run the extraction project.

NOTE

An internet connection is required in order to log into eArray.

- 1 Type your eArray **User Name** and **Password** in the eArray Login Setting window, if you already have an existing eArray account. Hint: The user name is the e-mail address you provided when you created an eArray account.

If you do not have an eArray account, you must first register for one. To start the registration process, click **Register**. This will take you to the Agilent eArray registration Web site where you can request a free eArray account. Follow the instructions provided on the eArray Web site to complete the registration.

- 2 Click **Test Connection**. The message “Login Successful!!!” is displayed if the connection to the eArray server is successful. If you get the message “Login denied”, check that the user name and password you typed are correct.
- 3 Under Advanced Options, select the appropriate boxes:
 - **Use eArray server during extraction**

Select this check box to log into eArray server to automatically download, install, and use grid templates (design files), Feature Extraction protocols, and QC metric sets during extraction.

- **Check for updates of grid template**

Select this check box to download updates to existing grid templates already in the Grid Template Browser. Updated grid templates are added to the database, leaving existing grid templates, unless you mark **Replace old grid template**. When this is marked, the newest grid template in the database will be used for feature extraction.

If several versions of a design are in the database and you want to do an extraction using an older design, clear this box. Otherwise, the extraction will automatically use the most recent version of the design (whether or not you are using the eArray server during extraction.)

- **Replace old grid template**

When you select this check box, Feature Extraction will replace old grid templates in the Grid Template Browser with the latest ones from the eArray server.

- **On starting Feature Extraction check for protocol update from eArray server**

Select this box to download updates from the eArray server for existing default protocols and QC metric sets in the FE Protocol Browser.

- 4 To save the eArray login settings, click **Save**. You will get a message "Setting updated successfully". Click **OK** to close the message, and click **Close** to exit the eArray Login Setting window.

Now you see the Feature Extraction main window.

To set up eArray login

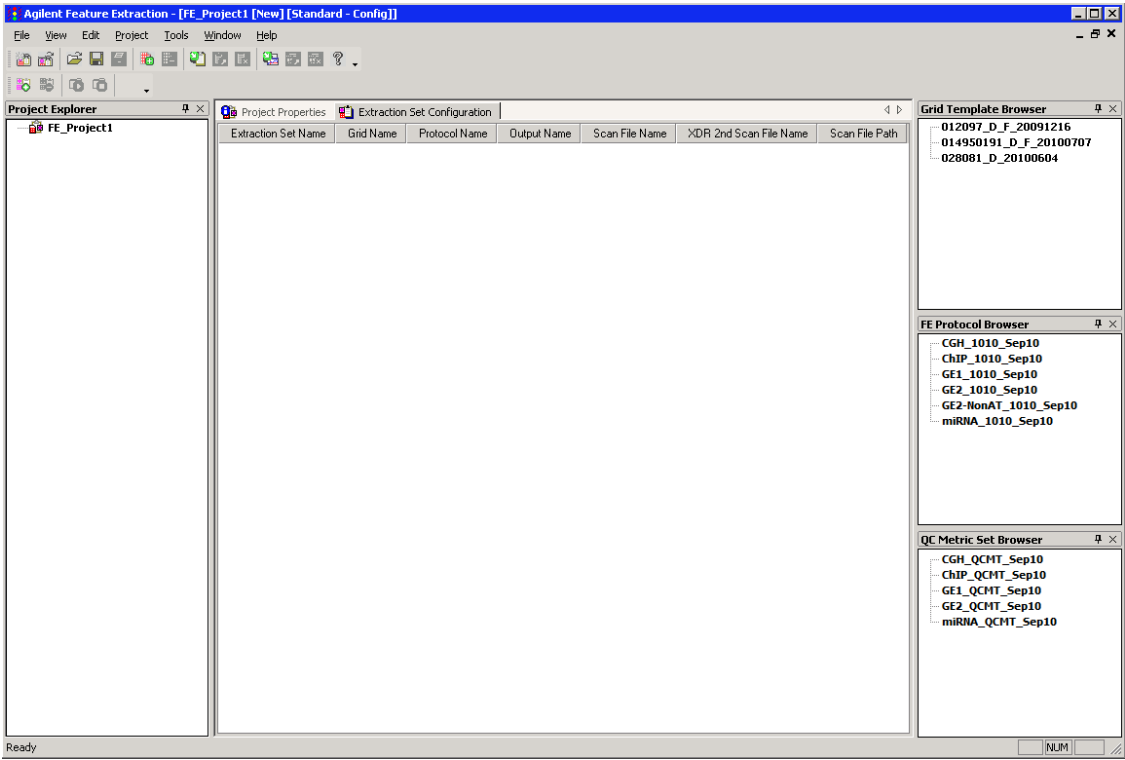


Figure 5 Feature Extraction Main Window

To open an image file

- 1 With the Feature Extraction main window open, click **File > Open > Image**, and browse to the folder that contains the image file. Double-click the image file (.tif) that you want to open.

If the Feature Extraction program is running, you can also “drag and drop” the image from your folder to the open Feature Application window.

If you do not have the Feature Extraction program running, you can also open an image file by dragging and dropping an image file to the Feature Extraction shortcut on your desktop.

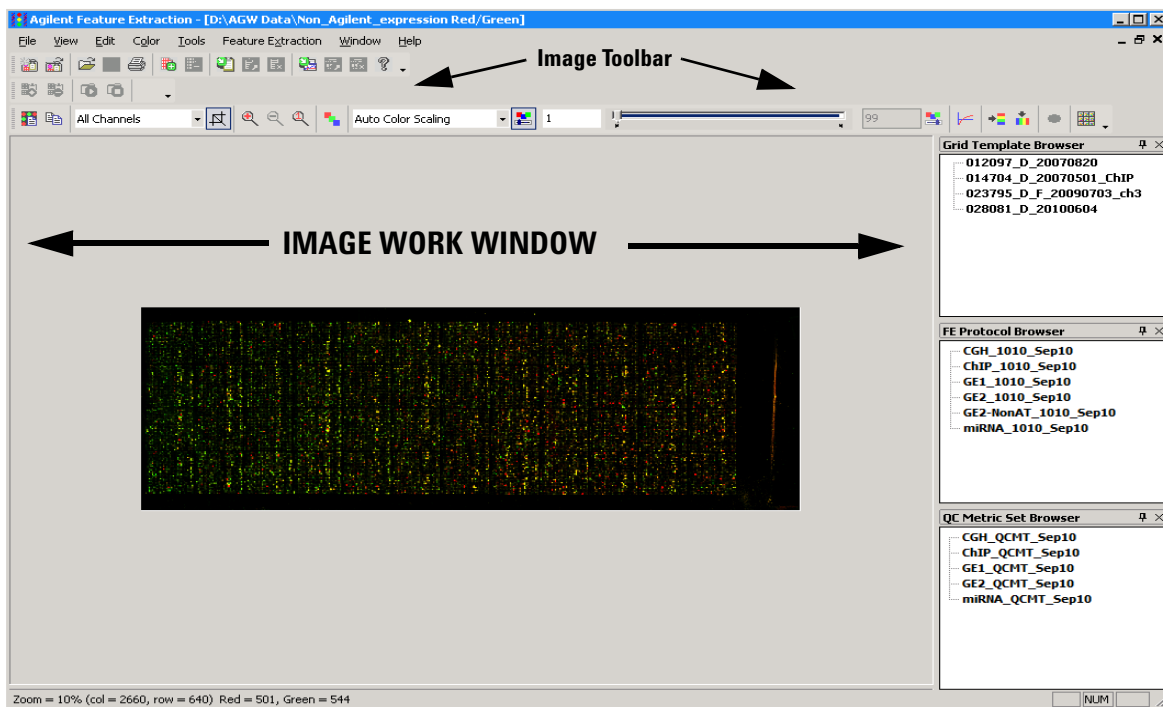


Figure 6 Feature Extraction main window with open image

NOTE

For high resolution images (2 or 3 micron or 20-bit), you see a low-resolution preview image until you crop to roughly 1/10 of the preview image size. Then you see the original high-resolution image. [The Crop Cursor turns red when you are outside the high-resolution (High Fidelity) region.]

To learn more about interactively finding spots and creating grid files for non-Agilent images, see Chapter 3, “Creating Grid Files”, in the *Feature Extraction 10.10 User Guide*.

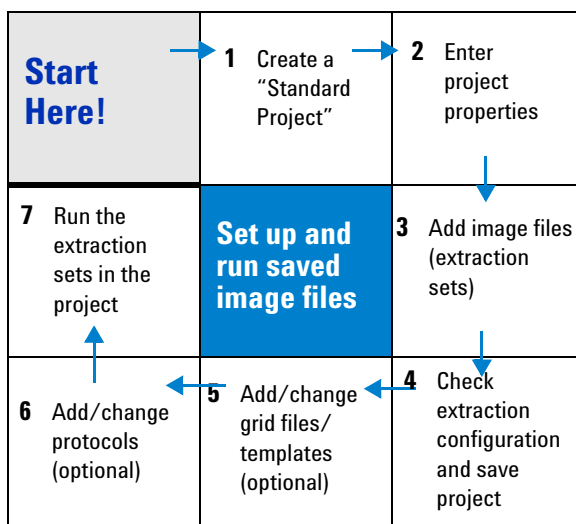
To learn more about changing the image display, see Chapter 5, “Changing Image Displays”, in the *Feature Extraction 10.10 User Guide*.

Setting up a Standard Project to extract Agilent and non-Agilent image files

Instructional overview

Follow the road map below and the instructions on the next pages to get started.

See Chapter 2 in the *Feature Extraction 10.10 User Guide* for details on this task.



The Agilent Feature Extraction program works with Agilent or non-Agilent microarray images produced on Agilent scanners using projects, extraction sets, grid templates or files, and protocols. The images are organized into extraction sets that contain the grid and protocol information to run the extraction. The extraction sets, in turn, are organized into projects.

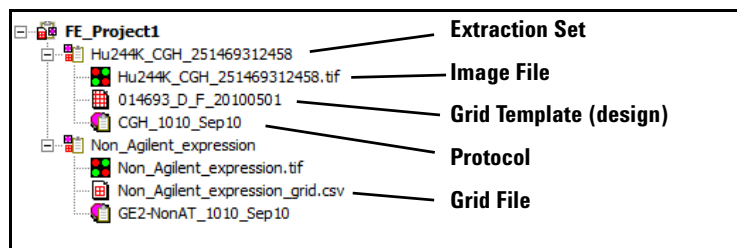


Figure 7 Project Explorer with expanded extraction sets

- project** A grouping of information that contains one or more *extraction sets* that define the files to use in the extraction.
- extraction set** Each group that contains a microarray image file, grid template (or grid file) and protocol within a project. This term is used both before and after extraction.
- grid template** Grid information from Agilent design files or grid files stored in the database. A grid template includes feature annotation and general geometry about the microarray (number of rows, columns, subgrids, feature spacings), which is used to find spots before data analysis takes place. Although not specific to any microarray image, grid template information is usually applied to the image for which the template was designed.
- grid file** Grid files are required in order to do feature extraction for non-Agilent images. Rarely, the automatic grid placement of Agilent images using a grid template may produce an incorrect grid, which will also require the use of a user-generated grid file. When using a grid file, extraction uses the nominal and “found” spot locations specific to the image from which the grid file was created. Grid files are created in the Feature Extraction program using a microarray image file and a microarray layout file (such as an Agilent Design file or a .GAL [GenePix Array List] file). The grid file contains the locations of the “found” spots if the “Calculate spot size and centroids” option is run before the file is saved. You can create grid templates from grid files, although the templates have only geometry information, no found spots.
- protocol** A list of steps and parameter values that define the data analysis algorithms and calculations used for feature extraction of a microarray image or images.

General Instructions

Step 1. Create a Standard Project

- If the Feature Extraction program is not already running, double-click the **Feature Extraction** shortcut on your desktop to start, or click the **Feature Extraction** icon in the Open Application tab of Agilent Genomic Workbench.

Feature Extraction automatically creates a standard project (FE_Project1) every time you start the program. The Feature Extraction main window appears with the Project1 work window and an empty Extraction Set Configuration tab window.

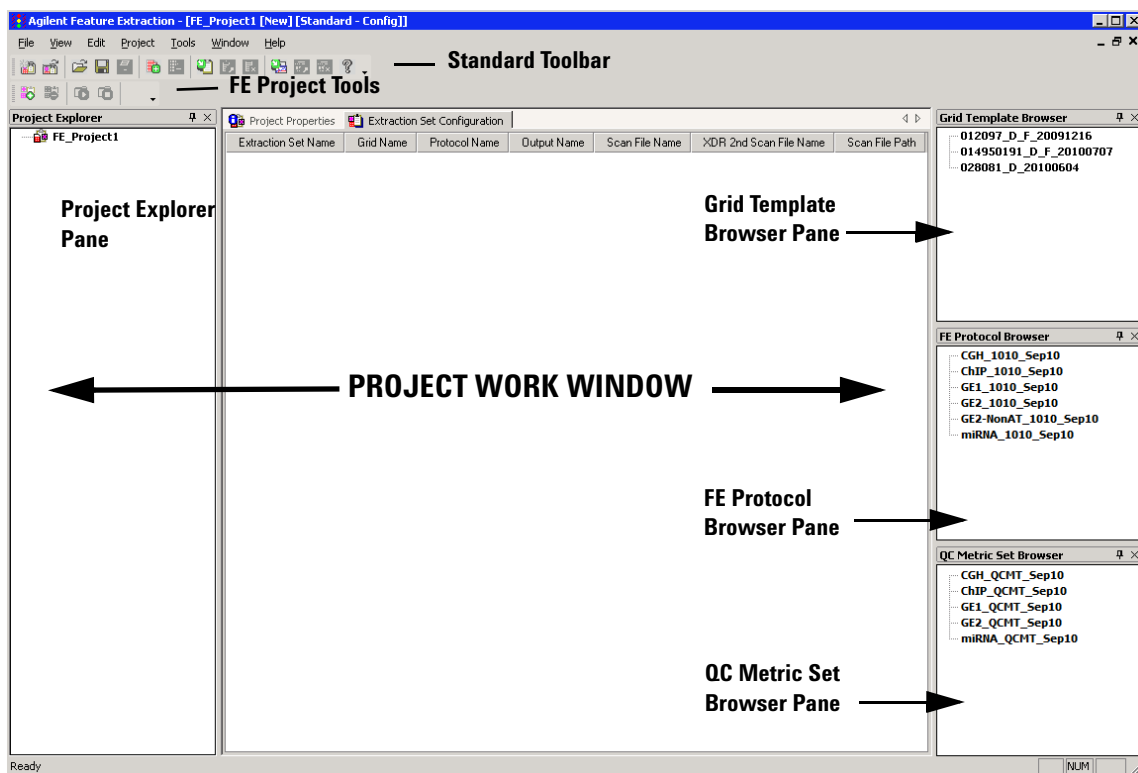


Figure 8 Feature Extraction main window

You can change where and how the project work window and the browser panes dock when you start the program. To learn more about working with this interface, see the *Feature Extraction 10.10 User Guide*.

Step 2. Enter the project properties

- 1 Click the **Project Properties** tab to view the default project properties.
- 2 To change a default setting, select a property and click the down arrow that appears on the right column. Choose from the available options.

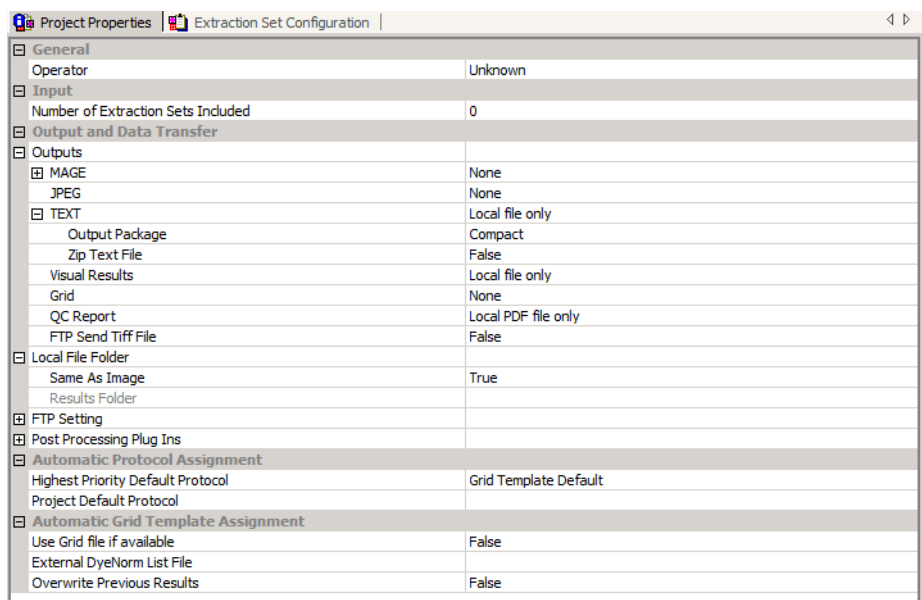


Figure 9 Default project properties for a standard project

You can run Feature Extraction for one or more Agilent images generated by the Agilent scanner with these default project properties. You can also set up and run multiple extractions with a combination of Agilent and non-Agilent images.

To learn more about setting up and running batch extractions with the Project Work Window, see the exercises in this guide:

- “[Example Exercise 1: Set up and run batch extractions on Agilent images](#)” on page 31.
- “[Example Exercise 2: Set up and run batch extractions on non-Agilent images](#)” on page 43.

Step 3. Add image files (extraction sets) to the project

- 1 In the Project Explorer, right-click **FE_Project1**, and click **Add Extraction**.
- 2 Select one or more images, and click **Open**.
 - To select multiple images at a time, hold down the **Ctrl** key and select the images.
 - You can also add one or more image files by dragging and dropping them from the folder to the open project in the Feature Extraction window.

Any extraction set that is missing the required components, such as grid template or protocol, shows a lightning bolt in Project Explorer and a “?” in the icon of the missing component. (Provided eArray is enabled and the image is an Agilent image.) In versions prior to 10.5, if any of the required components were not attached to the tiff image, then the extraction set would not be extracted. This changed beginning with version 10.5.

You do not need to assign a grid template/grid file and a protocol to an image file in order to run the extraction. Extraction sets without a grid template and/or a protocol attached to the project can be extracted, by setting the program to automatically determine the correct template and protocol to use. For this functionality to work, **one** of the following must be true:

- Prior to running the extraction, you set up to use the eArray server during extraction, as discussed in the section “[To set up eArray login](#)” on page 15.
- Prior to running the extraction, the required grid templates and protocols already exists in the database and are shown in the Grid Template Browser and FE Protocol Browser.

For custom grid templates and protocols that cannot be found on the eArray server, you must add them manually to the database before you run the extraction. See the *Feature Extraction 10.10 User Guide*.

If grid templates are not loaded automatically, or you want to use a different grid template or protocol, you can assign them manually. See Step 4 “[Assign grid template/grid file to the extraction set \(depending on type of images\)](#)” on page 26 and Step 5 “[Assign a protocol to the extraction set \(optional\)](#)” on page 27.

Step 3. Assign grid template/grid file to the extraction set (depending on type of images)

Agilent images By default, if the required grid templates are in the database, when you add Agilent tiff images to the project, Feature Extraction will automatically associate the appropriate grid templates to the extraction sets. You can also assign a grid template or grid file to the Agilent image.

To assign a grid template for Agilent images:

- 1 Expand the extraction set.
- 2 Select a grid template from the Grid Template Browser, hold down the mouse button and drag it to the second position of the extraction set in Project Explorer.
- 3 Let go of the mouse button to drop it at this position.

To assign a grid file for Agilent images:

- 1 Right-click the second position of the expanded extraction set and click **Select Grid File**.
- 2 Select the grid file, and click **Open**.

The grid file name will appear in the extraction set, along with the grid template on which it is based.

Non-Agilent images You must assign a grid file to non-Agilent images in the project. The extraction automation feature is not supported for non-Agilent images. Feature Extraction will not extract non-Agilent images that are missing grid files or protocols.

To assign a grid file for non-Agilent images:

- 1 Right-click the second position of the expanded extraction set and click **Select Grid File**.
- 2 Select the grid file, and click **Open**.

AMADID is the
Agilent MicroArray
Design Identifier.

Agilent microarray images scanned with an Agilent scanner For Agilent microarray images with barcode information included in the image, the program selects a grid template from the database using the AMADID portion of the barcode associated with the image file. If there are grid templates for that AMADID that come from multiple design files, the program selects the grid template from the most recently created Agilent design file. To add a barcode if the image has none, see the *Feature Extraction 10.10 User Guide*.

For this automated assignment to work, the grid template must exist in the database. If it does not, you can let the Feature Extraction try to automatically download the required grid template from the eArray server or you can interactively associate a grid template or grid file with the extraction set. See [page 26](#) for details on how to assign a grid template or grid file.

Agilent uses barcode information for many workflow automation and data tracking steps. Inclusion of the Agilent barcode information in the TIFF file helps ensure that the data is quickly, easily and accurately processed and that informatics programs know how to analyze the results.

Non-Agilent images scanned with an Agilent scanner Non-Agilent image files from an Agilent scanner do not point to grid information that the Agilent Feature Extraction program can read. Therefore, when you load the non-Agilent image file, you must add a grid file that you have already created.

See Chapter 3 in the *Feature Extraction 10.10 User Guide* to learn how to create grid files.

Step 3. Assign a protocol to the extraction set (optional)

Automatically assigned protocols come from two sources: a default protocol associated with the grid template, or a project level default protocol. You select which of these options has the higher priority. (See Chapter 2 of the *Feature Extraction 10.10 User Guide* to learn how to associate a protocol with a grid template and where to enter the project level default protocol.) You can also add a protocol to the extraction set manually.

To add a protocol manually:

- 1 From the Protocol Browser select a protocol, hold down the mouse button and drag it to the third position of the extraction set.
- 2 Release the mouse button to drop the protocol onto the third position of the set.

Feature Extraction can run analyses on Agilent 1-color gene expression (GE1), 2-color gene expression (GE2), CGH, CGH+SNP, ChIP, miRNA, and non-Agilent 2-color GE (GE2-nonAT) microarray images scanned on the Agilent scanner. Agilent includes Feature Extraction protocols for extracting each of these image types.

CAUTION

These Feature Extraction Protocols were optimized using data from Agilent catalog arrays, which have many replicated probes and validated Negative Control probes. If custom arrays without enough replicated probes are used, or arrays with custom probes designated as Negative Control probes, the default Feature Extraction Protocols may not be optimal. The Feature Extraction Protocol optimization also uses Agilent's default lab protocols; that is, label, hybridization, wash, and scanner procedures.

Step 3. Add or change grid templates from design files (optional)

For the Agilent images you intend to extract, you add grid templates based on the images' design files to the Grid Template Browser.

- 1 Right-click the Grid Template Browser pane, and click **Add**.
- 2 Select a design file or files, and click **Open**.

Every extraction set must have an image file, assigned grid template/grid file and protocol. Grid templates and protocols for Agilent images are loaded to the extraction set automatically by the program, if the grid template associated with the loaded image's design file is available in the database. If a grid template is in the database, it is displayed automatically in the Grid Template Browser. If a grid is not in the database, the program will automatically get the grid template from eArray when the extraction starts. For details on how to update or add grid templates using Agilent eArray, see the *Feature Extraction 10.10 User Guide*.

Step 3. Change QC metric set associated with the protocol to monitor QC metrics

At times the Feature Extraction program is used in a production environment, where the microarray processing protocols are similar and monitoring run-to-run consistency is an important goal. Feature Extraction can help monitor this consistency with the QC metric sets.

In earlier versions of Feature Extraction, you choose one QC metric set in the Project Properties tab that was used for all extractions in a project.

In version 10.5 and higher, the QC metric sets and protocols are linked. The Feature Extraction protocol defines which metric set is used with it.

To change the QC metric set associated to the Feature Extraction protocol:

- 1 In the FE Protocol Browser, double-click a protocol to open the FE Protocol Editor.
- 2 Select **Protocol Properties** to see the protocol properties.
- 3 Under General, click the cell next to **Metricset**.
- 4 Click the down arrow, and select the QC metric set appropriate for your microarray application.

If a metric set you want to use is not available in the metric set list, you must first import it to the database.

To import a QC metric set into the database:

- 1 Select **View > QC Metric Set Browser** from Feature Extraction main window, if it is not already visible.
- 2 The QC Metric Set Browser displays all of the default metric sets and custom metric sets in the database. Custom metric sets are created using the Quality module in Agilent Genomic Workbench. Right-click in the QC Metric Set Browser and select **Import** to import metric sets that are not shown in the QC Metric Set Browser.

The default QC metric sets installed with the Feature Extraction software are read-only and cannot be removed from the database. However, you can remove or change a QC metric set associated with the protocol and then save it as a new protocol with a different name.

To remove a QC metric set from the protocol:

- Select **<Set empty>** from the metric set list in the Protocol Properties.
- 5 Click **Save As** to save the new protocol.

QC metric sets Metric sets are loaded into the database when you install the Feature Extraction software. If you set up the eArray login as discussed in the section “[To set up eArray login](#)” on page 15, then Feature Extraction will automatically download and install the latest QC metric sets from the eArray server.

If you remove a QC metric set, updates of QC metric sets are also available at www.agilent.com/chem/feqcmetrics, if you need to manually download and install them in Feature Extraction.

CAUTION

Agilent-supplied QC metric sets are intended to assist users in monitoring microarray processing issues. They were not developed to detect microarray manufacturing issues.

When you create QC metric sets using the Agilent Genomic Workbench Quality tools, you can associate them to protocols in Feature Extraction. When that project is run, the Feature Extraction program summarizes the metric statistics on each microarray's QC report and shows how the results compare to the thresholds (if any) defined in the metric set. It also reports this information in the output files in the statistics table. For more information on using the Quality module in Agilent Genomic Workbench, see the Agilent Genomic Workbench *Quality Tools User Guide*.

Step 6. Save the project

- 1 Select **File > Save As**.

By default, your projects are saved in **C:\Program Files\Agilent\Agilent Genomic Workbench Standard Edition <version>\FeatureExtraction\FEProjects**.

- 2 To save in a different location, go to or create the folder for your projects.
- 3 Type a project name, and click **Save**.

Step 4. Run Feature Extraction

- 1 Click the **Project Run mode On/Off** button on the Feature Extraction Project toolbar , or select **Project > Start Extracting**.

Example Exercise 1: Set up and run batch extractions on Agilent images

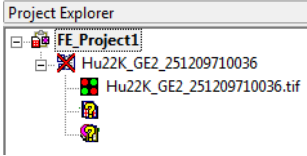
The following exercise takes you through the setup, analysis, and reporting steps of Feature Extraction with example images provided by Agilent. These images are provided on the Agilent Genomic Workbench installation disk.

With this exercise you learn how to:

- Set up and extract a 2-color GE microarray image and produce a 2-color GE QC report with spike-ins.
- Set up and extract four 1-color GE images on a 4-multiplex (multi pack) microarray and produce a 1-color GE QC report.

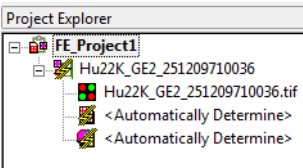
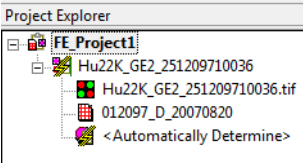
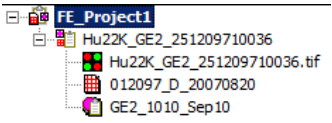
Try the Steps on the left before reading the Detailed Instructions.

Task 1: Add an extraction set for a 2-color Gene Expression microarray image file

Steps	Detailed Instructions	Comments
1 Add the 2-color gene expression image file that was installed with the Feature Extraction software to Project 1.	<p>a Start the Feature Extraction program.</p> <p>b Click the Add Extraction Set(s) button on the toolbar, or right-click Project Explorer, and select Add Extraction.</p> <p>c The example images are located in the installation folder for the Agilent Genomic Workbench client. By default, this is C:\Program Files\Agilent\Agilent Genomic Workbench Standard Edition <version>\FeatureExtraction\ExampleImages. Select the Hu22K_GE2_251209710036.tif file, and click Open.</p> <p>d Expand the extraction set that appears in Project Explorer.</p>	<ul style="list-style-type: none">• There may be times when you need to manually assign a grid template and/or a protocol to the extraction set. Task 1 is intended to show you how to do this.• The grid template associated to this image is automatically assigned to the extraction set when the grid template is available in the Grid Template Browser.
		

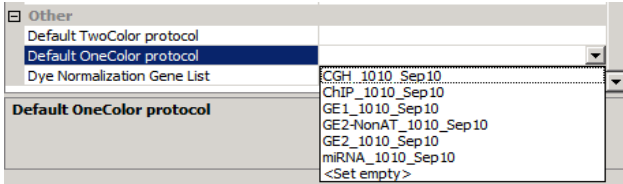
Example Exercise 1: Set up and run batch extractions on Agilent images

Task 1: Add an extraction set for a 2-color Gene Expression microarray image file (continued)

Steps	Detailed Instructions	Comments
<p>2 Add a grid template to the extraction set, if one has not been automatically assigned.</p> <p>Note: The grid template for this image is a design file that begins with 012097_D...</p>	<p>a Right-click the Grid Template Browser, and click Add.</p> <p>b Go to the ExampleImages folder, or the folder where the image file is stored.</p> <p>c Select the design file (012097_D_20070820.xml), and click Open.</p> <p>d When the message appears that tells you the grid template was successfully added, click OK.</p>  <p>e Drag and drop the new grid template to the second row, after the image name.</p> 	<ul style="list-style-type: none">• The program will automatically assign the appropriate grid template and protocol for the image if <Automatically Determine> is displayed in the extraction set, and you have selected in the eArray settings to use the eArray server during extraction.
<p>3 Assign the gene expression 2-color protocol to the extraction set.</p> <p>Note: The 2-color GE protocol for this extraction set has spike-ins turned on as the default setting. The QC report will show all the spike-in results.</p>	<ul style="list-style-type: none">• From the FE Protocol Browser, drag and drop the GE2_1010_Sep10 Feature Extraction protocol to the third row of the gene expression extraction set. 	<ul style="list-style-type: none">• You can add a GE2_105 (version 10.5) Feature Extraction protocol to the Grid Template Browser if you want to compare data run on v10.5 vs. v10.10. The extraction set using v10.5 Feature Extraction protocol is analyzed using v10.5 data processing, but with v10.10 for image processing.• The latest protocols can be automatically downloaded from eArray server and installed on Feature Extraction, if you select this in the eArray settings.

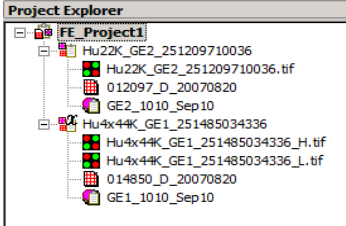
Task 2: Add an extraction set for a 1-color Gene Expression image file of 4-multiplex microarray

Steps	Detailed Instructions	Comments
1 Add a grid template from the design file. Note: The design file for this image is 014850_D_20070207.xml.	a Right-click the Grid Template Browser, and click Add . b Go to the directory where the image file is stored. c Select the design file (014850_D_20070820.xml), and click Open . d When the message appears that tells you the design was added successfully, click OK .	<ul style="list-style-type: none">Task 2 shows you how to associate a default Feature Extraction protocol with a grid template. When the image and grid template are loaded in the project, Feature Extraction automatically attaches the default protocol to the extraction set.
2 Assign a default 1-color GE protocol to the grid template to use for the microarray image with barcode 25148503446.	a Right-click grid template 14850_D_20070820 in the Grid Template Browser and select Properties . b In the Grid Template Properties page, click the Default OneColor protocol row. c From the Default OneColor protocol list select GE1_11010_Jun10 , and click OK .	<ul style="list-style-type: none">The software recognizes to use the default 1-color protocol when the 1-color scan file is loaded.



Example Exercise 1: Set up and run batch extractions on Agilent images

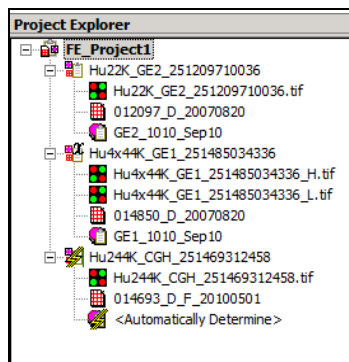
Task 2: Add an extraction set for a 1-color Gene Expression image file of 4-multiplex microarray (continued)

Steps	Detailed Instructions	Comments
<p>3 Add the 1-color GE example image file installed with the Feature Extraction software to Project 1.</p> <p>Note: This image has been renamed from the original scan name. The “14850” in the name corresponds to the AMADID.</p>	<p>a Right-click Project Explorer, and select Add Extraction.</p> <p>b Select the Hu4x44K_GE1_251485034336_H.tif file, and click Open.</p> <p>c Expand the extraction set that appears in Project Explorer.</p> 	<ul style="list-style-type: none">• Feature Extraction lets you extract the multiplex images as one image file and produces one .shp shape file for the four microarrays in the multiplex.• This is an XDR image scanned at high and low PMTs. Adding a high PMT scan (_H) automatically loads the corresponding low PMT scan (_L) that is in the same directory.• Feature Extraction automatically loads the grid template which has the default 1-color GE protocol assigned to it. As a result, the protocol is also loaded to the extraction set.

Example Exercise 1: Set up and run batch extractions on Agilent images

Task 3: Add an extraction set for a CGH microarray image file

Steps	Detailed Instructions	Comments
1 Add the 20-bit CGH image file installed with the Feature Extraction software to Project 1. Note: The grid template for this image is not provided as part of the Example Images installation.	a Click the Add Extraction Set(s) button on the toolbar, or right-click Project Explorer, and select Add Extraction . b Select the file Hu244K_CGH_251469312458.tif , and click Open . c Expand the extraction set that appears in Project Explorer.	<ul style="list-style-type: none"> Task 3 shows you how to extract a tiff image without a grid template and protocol attached in the project when you set up the eArray login.



2 Check the Extraction Set Configuration tab to make sure that the assignments are OK for the GE 1-color and 2-color extraction sets.	a Expand the columns in the Extraction Set Configuration tab to see all the entries for the three extraction sets.	<ul style="list-style-type: none"> Note: The output file name is the name of the extraction set plus the protocol name. Note: The CGH extraction set does not have a protocol assigned to it.
--	---	---

Extraction Set Name	Grid Name	Protocol Name	Output Name	Scan File Name	XDR 2nd Scan File Name	Scan File Path
Hu22K_GE2_25120971_D_20070820		GE2_1010_Sep10	Hu22K_GE2_25120971	Hu22K_GE2_25120971	<None>	D:\AGW Data\
Hu244K_CGH_25146931_F_20100501			Hu244K_CGH_25146931	Hu244K_CGH_25146931	<None>	D:\AGW Data\
Hu4x44K_GE1_25148503_D_20070820		GE1_1010_Sep10	Hu4x44K_GE1_25148503	Hu4x44K_GE1_25148503	Hu4x44K_GE1_25148503	D:\AGW Data\

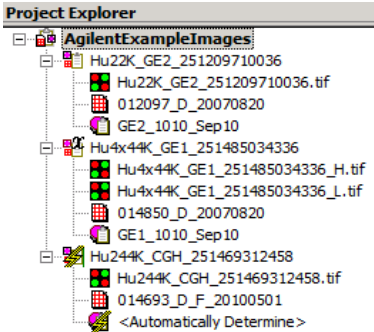
Example Exercise 1: Set up and run batch extractions on Agilent images

Task 3: Add an extraction set for a CGH microarray image file (continued)

Steps	Detailed Instructions	Comments
3 Enter eArray login settings, if you have not yet done so. You must select the Use eArray server during extraction in order for this exercise to work. Note: You must have an eArray account before you can set up the login settings in Feature Extraction.	a Select Tools > eArray Settings . b In the eArray Login Setting window, if you do not have an eArray account, click the Register button to sign up for one. c Enter your eArray login information, and click the Test Connection button. You should get a message indicating successful login. d In Advanced Options , select Use eArray server during extraction . Click Save and then close the eArray Settings dialog box.	<ul style="list-style-type: none">The settings let Feature Extraction automatically log into the eArray server to check for updates of grid templates, protocols, and QC metric sets.

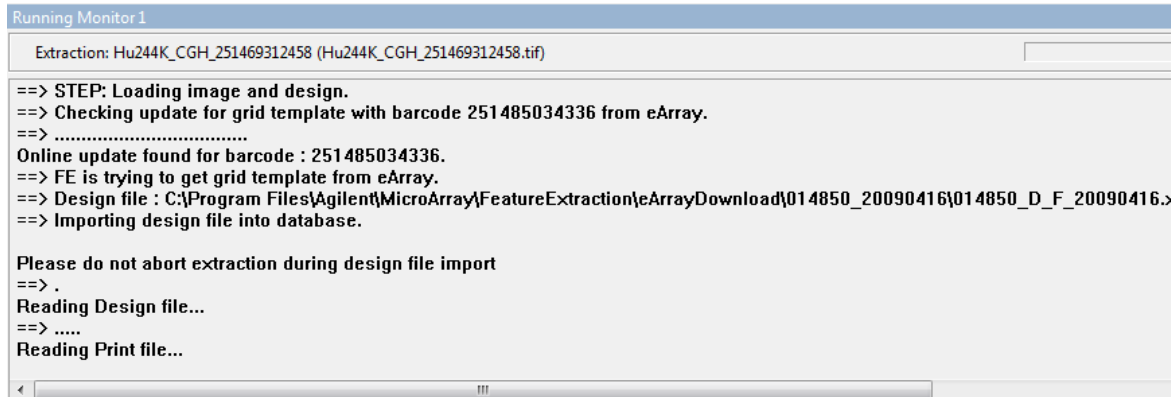
Task 4: Save the project and start the Feature Extraction batch run

Steps	Detailed Instructions	Comments
1 Save the project with a new name.	a Select File > Save As . b Go to or create the folder for your projects, if you want to save in a location different from the default FEProjects directory. c Enter AgilentExampleImages (or anything you want) in the File Name box, and click Save .	<ul style="list-style-type: none">You must save the project before the program will start extracting.After you save the project, the project name changes in both the title bar and Project Explorer.



Task 4: Save the project and start the Feature Extraction batch run (continued)

Steps	Detailed Instructions	Comments
2 Run Feature Extraction on AgilentExampleImages project. Note: The Running Monitor window shows the progress of each extraction set.	a Click the Project Run mode On/Off button, or select Project > Start Extracting . The project extraction starts to run.	<ul style="list-style-type: none">• In Feature Extraction 10.10 only one extraction set is extracted at a time.



Task 5: View project run summary and QC report

Steps	Detailed Instructions	Comments
1 View the Summary Report tab for the project run summary.	<ul style="list-style-type: none">• The Summary Report tab appears next to the Extraction Set Configuration tab when the project run starts.• See Figure 10 on page 39 for details.	<ul style="list-style-type: none">• Once extraction is done, the Project Run Summary shows information, warnings, errors, and QC reports associated to each extraction set.• For multiplex microarray images, a QC report is generated for each microarray.• Note: The 4-multiplex 1-color GE microarray image will have four QC reports, one for each microarray.

Example Exercise 1: Set up and run batch extractions on Agilent images

Task 5: View project run summary and QC report (continued)

Steps	Detailed Instructions	Comments
<p>2 Open the QC report for the 2-color GE extraction set, and see whether the results seem reasonable or not.</p> <p>See Chapter 2 in the <i>Feature Extraction 10.10 Reference Guide</i> for information on how to interpret the QC Report.</p>	<p>a In the Project Run Summary on-screen report, double-click the link to the QC report for Hu22K_GE2.</p> <p>Or do the following:</p> <ul style="list-style-type: none">Go to the folder where the .tif image file is saved.Double-click the .pdf file for the QC report. <p>b Display the grid, spatial distribution plot and plot of median % CV to see any anomalies.</p>	<ul style="list-style-type: none">The QC report (.pdf file) by default is put in the same folder as the image file. You can change the default folder setting by selecting Tools>Preferences from the toolbar, and then opening the Output & Data Transfer folder. The default results folder is defined in the Local File Folder setting.If you are running Feature Extraction with a 30-day demo license, the QC report is an .html file. See Chapter 2 of the <i>Feature Extraction User Guide</i> to learn to how print this .html file.

Task 5: View project run summary and QC report (continued)

Steps	Detailed Instructions	Comments
<div> Project Properties Extraction Set Configuration Summary Report </div> <h2>Project Run Summary: <i>AgilentExampleImages</i></h2> <p>Project started on Fri, Sep 10, 2010 at 16:32:35.</p> <p>D:\AGW Data\Hu22K_GE2_251209710036.tif Fri Sep 10 16:32:35 2010</p> <p>INFO: Grid Template in use : 012097_D_20070820</p> <p>INFO: Protocol in use: GE2_1010_Sep10</p> <p>INFO: MetricSet in use: GE2_QCMT_Sep10.</p> <p>INFO: Agilent supports but has not tested this configuration, please consult the support matrix in the Feature Extrac</p> <p>INFO: FE automatically determined that ArrayFormat is Single Density</p> <p>INFO: FE automatically determined that GridFormat is 65 micron feature size</p> <p>INFO: FE automatically determined that SpotFormat is Single Density</p> <p>INFO: FE automatically determined that Scanner is Agilent Scanner</p> <p>INFO: FE automatically determined that GridTestFormat is 65 micron feature size</p> <p>INFO: There are 1829 (Red) negative features (Non-Controls).</p> <p>INFO: There are 5445 (Green) negative features (Non-Controls).</p> <p>INFO: 16 (Red) and 1 (Green) saturated features</p> <p>INFO: 4 (Red) and 8 (Green) feature non-uniformity outliers</p> <p>INFO: 98 (Red) and 73 (Green) feature population outliers</p> <p>INFO: 9 (Red) and 0 (Green) background non-uniformity outliers</p> <p>INFO: 48 (Red) and 0 (Green) background population outliers</p> <p>INFO: QCMetrics Totals: Found 12 Metrics In Good Range</p> <p>INFO: Extraction completed successfully.</p> <p>QCREPORT: <D:\AGW Data\Hu22K_GE2_251209710036_GE2_1010_Sep10.pdf></p> <p>D:\AGW Data\Hu4x44K_GE1_251485034336_H.tif Fri Sep 10 16:33:31 2010</p> <p>Processing 4 arrays in single scan.</p> <p>INFO: Grid Template in use : 014850_D_20070820</p> <p>INFO: Protocol in use: GE1_1010_Sep10</p> <p>INFO: multiple extractions</p> <p>INFO: MetricSet in use: GE1_QCMT_Sep10.</p> <p>INFO: FE automatically determined that ArrayFormat is 65 micron feature size</p> <p>INFO: FE automatically determined that GridFormat is 65 micron feature size</p>		

Figure 10 Summary report for AgilentExampleImages project

Example Exercise 1: Set up and run batch extractions on Agilent images

Task 5: View project run summary and QC report (continued)

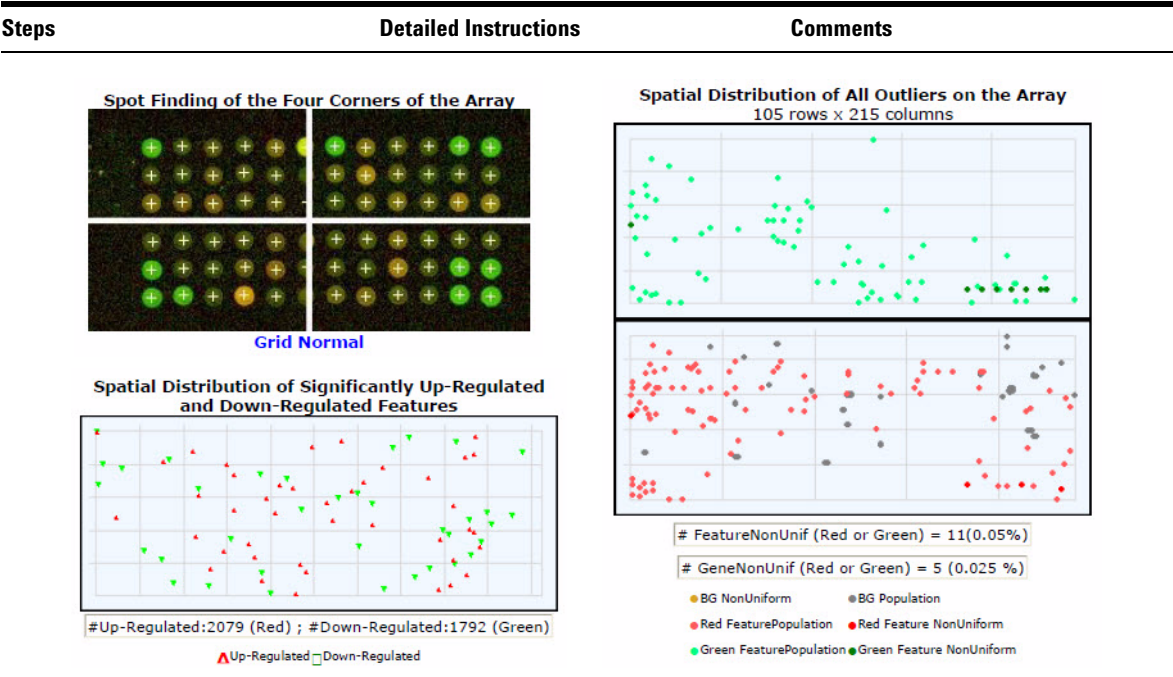


Figure 11 Grid placement at four corners and spatial distribution plots for Hu22K_GE2 QC report

Task 5: View project run summary and QC report (continued)

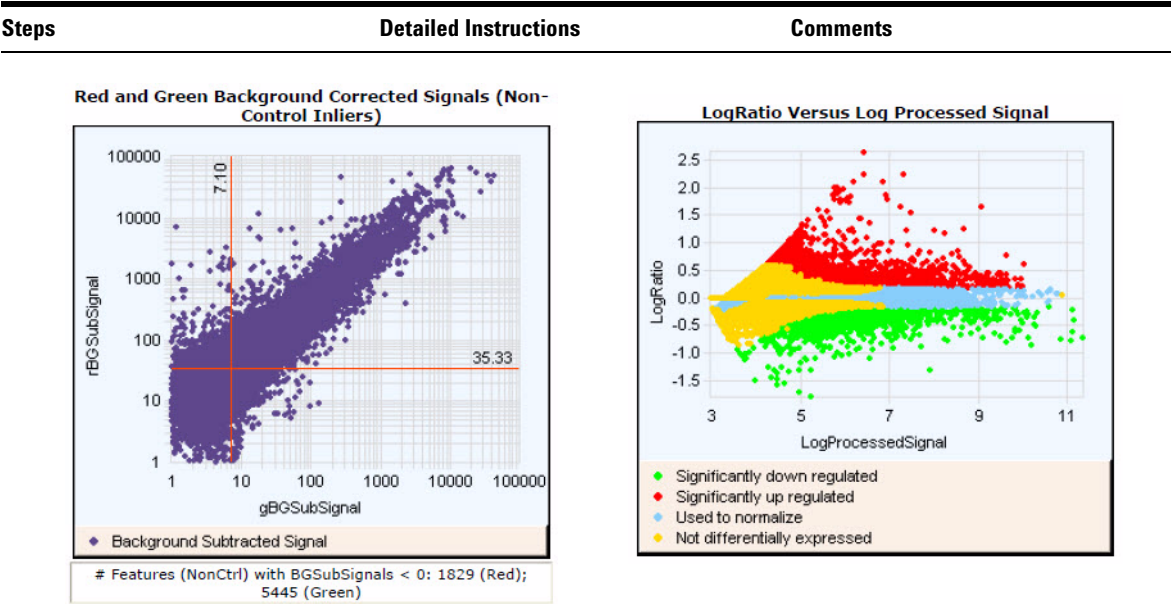


Figure 12 Background corrected signals and log ratio vs. processed signals plots for Hu22K_GE2 QC report

3 Print the latest Project Run Summary report for the project.

- Go to the results folder where the **AgilentExampleImages** project is saved.
- Open and print the **_LastBatchReport.rtf** with Microsoft Word.

- All Summary Reports are saved to the folder containing the project associated with the report.
- Note that there are two reports with the .rtf extension. One is always a specifically-named rtf (ProjectName_Date). The Computer Name_LastBatch file is used for the last extraction run and is overwritten by the next project, if same folder is used.
- If an image was extracted previously, the previous .rtf files are named as Project Name_Date and Time.rtf.

Name	Size	Type	Date Modified
AgilentExampleImages.fep	3 KB	FEP File	12/13/2008 6:15 PM
AgilentExampleImages_200812131841.rtf	12 KB	Rich Text Format	12/13/2008 6:41 PM
T43-2687D3U_LastBatchReport.rtf	12 KB	Rich Text Format	12/13/2008 6:41 PM

Example Exercise 1: Set up and run batch extractions on Agilent images

CAUTION

Agilent-supplied QC metric sets are intended to assist users in monitoring microarray processing issues. They were not developed to detect microarray manufacturing issues.



Example Exercise 2: Set up and run batch extractions on non-Agilent images

This exercise shows you how to create a new project, change project properties and run Feature Extraction for the non-Agilent image shipped with the software.

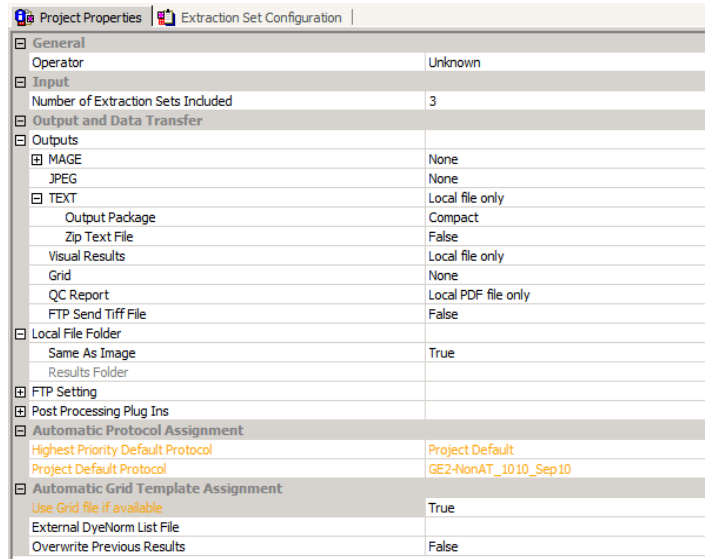
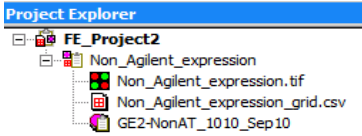
Repeat these instructions to add non-Agilent images of your own to create a batch. You can also set up and run a combination of Agilent and non-Agilent images in one project.

Scans must be from Agilent scanner. Other scanners are not supported.

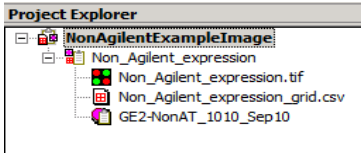
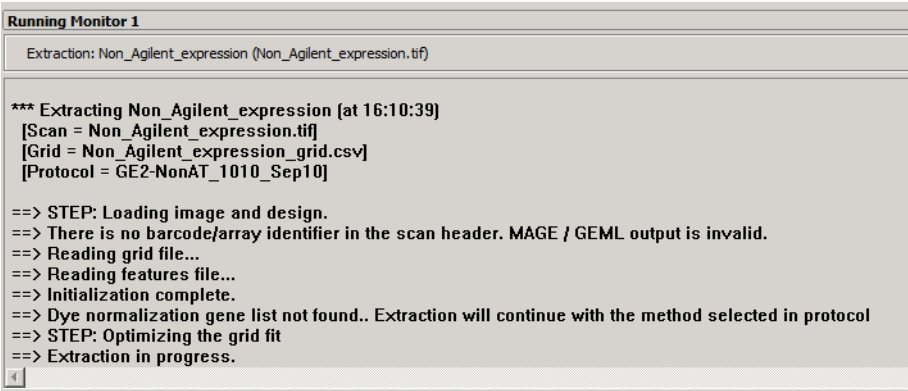
Try the Steps on the left before reading the Detailed Instructions.

Steps	Detailed Instructions	Comments
1 Create a new standard project. Note: After FE_Project1, the next project you open is FE_Project2. The name of every project that you open after that follows in numerical order, whether you save FE Project N or not or close FE Project N or not.	<ul style="list-style-type: none"> Select File > New > Standard Project, OR Click the New Project icon,  	<ul style="list-style-type: none"> A standard project is one where extraction sets are set up from existing scan files to run automatically as a batch. Each time you restart the software, the first new project is FE_Project1.
2 Change the following Project Properties: <ul style="list-style-type: none"> Highest Priority Default Protocol: Project Default Project Default Protocol: GE2-nonAT_105_Sep09 Use Grid file if available: True 	<ol style="list-style-type: none"> Click the Project Properties tab. Click the cell next to Highest Priority Default Protocol, and select Project Default. Click the cell next to Project Default Protocol, and select GE2-nonAT_1010_Sep10. Click the cell next to Use Grid file if available, and select True. 	<ul style="list-style-type: none"> In the Project Properties tab screen, you can select a number of variables that pertain to the project as a whole, rather than to individual extraction sets. Because some of these properties, such as Project Default Protocol, affect how the extraction set is constructed, you may want to set these before you add the image files (extraction sets).

Example Exercise 2: Set up and run batch extractions on non-Agilent images

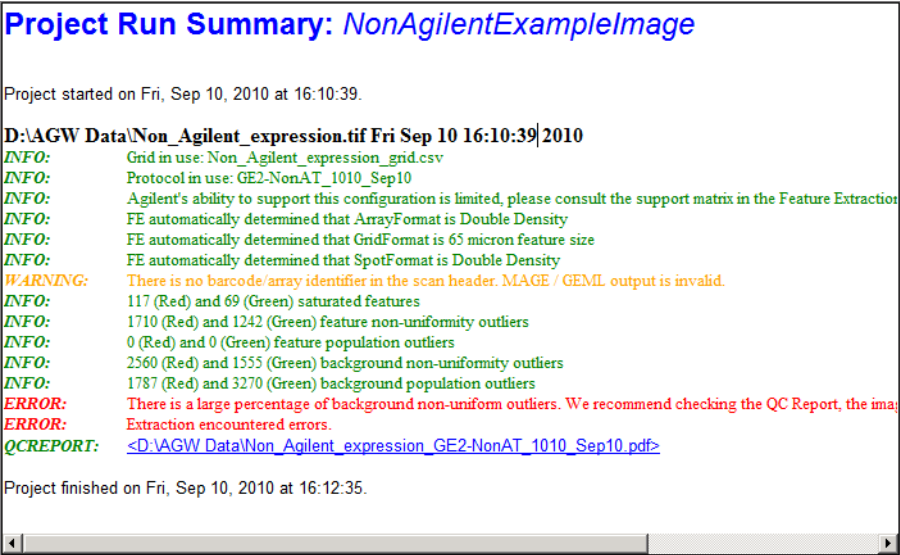
Steps	Detailed Instructions	Comments
		<ul style="list-style-type: none">• See Chapter 2 in the <i>Feature Extraction 10.10 User Guide</i> for instructions on how to change other project properties• Note that any changes from the default settings are shown in orange.
3	<p>Add the non-Agilent image shipped with the software to Project 2.</p> <ul style="list-style-type: none">• Add any other of your non-Agilent images that you want in the batch. <p>a Click the Add Extraction Set(s) button, or right-click anywhere in Project Explorer, and select Add Extraction.</p> <p>b Go to the directory containing the non-Agilent image, and click Open.</p> <p>c Expand this extraction set.</p> <p>d Right-click grid icon (below the tif file), and click Select Grid file.</p> <p>e Browse to the folder containing the grid file. For this example, it is located in the same folder as the image.</p> <p>f Select the grid file (_grid.csv) and click Open.</p>	<ul style="list-style-type: none">• The grid is automatically assigned, if its name is the same as the image name and you set the Project Properties parameter "Use Grid File if available" to "True".• Note: In this example, because the image file name "Non_Agilent_expression" is NOT the same as the grid file name "Non-Agilent", the grid will not be automatically assigned to this image. You must manually select a grid file to use.• The protocol was automatically assigned to the extraction set because you set the highest priority default protocol to be the one for the project and then selected the non-Agilent protocol as the default project protocol.
		

Example Exercise 2: Set up and run batch extractions on non-Agilent images

Steps	Detailed Instructions	Comments
4 Save the project with a new name.	<p>a Select File > Save As.</p> <p>b Go to or create the directory for your projects.</p> <p>a Enter nonAgilentExampleImage, and click Save.</p>	<ul style="list-style-type: none"> For non-Agilent images you must create a grid file (grid.csv) for use in the extraction set. For more information see Chapter 3 in the <i>Feature Extraction User Guide</i>.
		
5 Run Feature Extraction on nonAgilentExampleImage project. Note: The Running Monitor window shows the progress of each extraction set.	<p>a Click the Project Run mode On/Off button, or select Project > Start Extracting.</p> <p>The project extraction starts to run.</p>	<ul style="list-style-type: none"> In Feature Extraction 10.10 only one extraction set is extracted at a time.
		

Example Exercise 2: Set up and run batch extractions on non-Agilent images

Steps	Detailed Instructions	Comments
6 View the Summary Report tab for the project run summary.	<ul style="list-style-type: none">Summary Report tab appears next to the Extraction Set Configuration tab when the project run starts.	The Project Run Summary report shows a red Error message that there were a large percentage of background non-uniform outliers. User should check the QC Report and may need to make a grid file to re-extract this image. See the <i>Feature Extraction User Guide</i> for information on how to create a grid file.



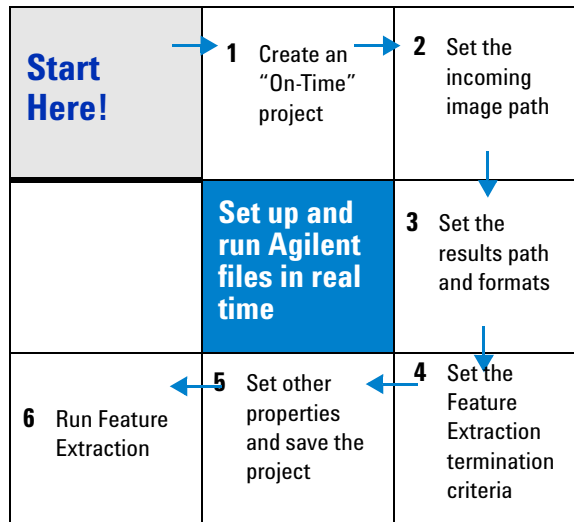
7 View the QC Report	b Double-click the blue line containing the link to the QC report.	On the QC report, note the red Evaluate Grid message under Spot Finding of the Four Corners of the Array.
----------------------	---	--

Setting up an On-Time project to extract Agilent image files in real time

Instructional overview

An On-Time project lets you set up the Feature Extraction software to automatically extract Agilent microarray image files as they are transferred from the Agilent scanner to a folder of your choosing.

Follow the workflow below and the instructions on the next pages to get started.



NOTE

On-Time projects work only with Agilent microarrays scanned on the Agilent scanner.

General Instructions

These instructions help you set up the program to run a batch of Agilent microarray images in real time as they arrive from the Agilent scanner to a folder of your choosing. These instructions do not include running Feature Extraction.

Step 1. Create an on-time project

- Select **File > New > On-Time Project**.
Note that you do not add extraction sets to on-time projects.

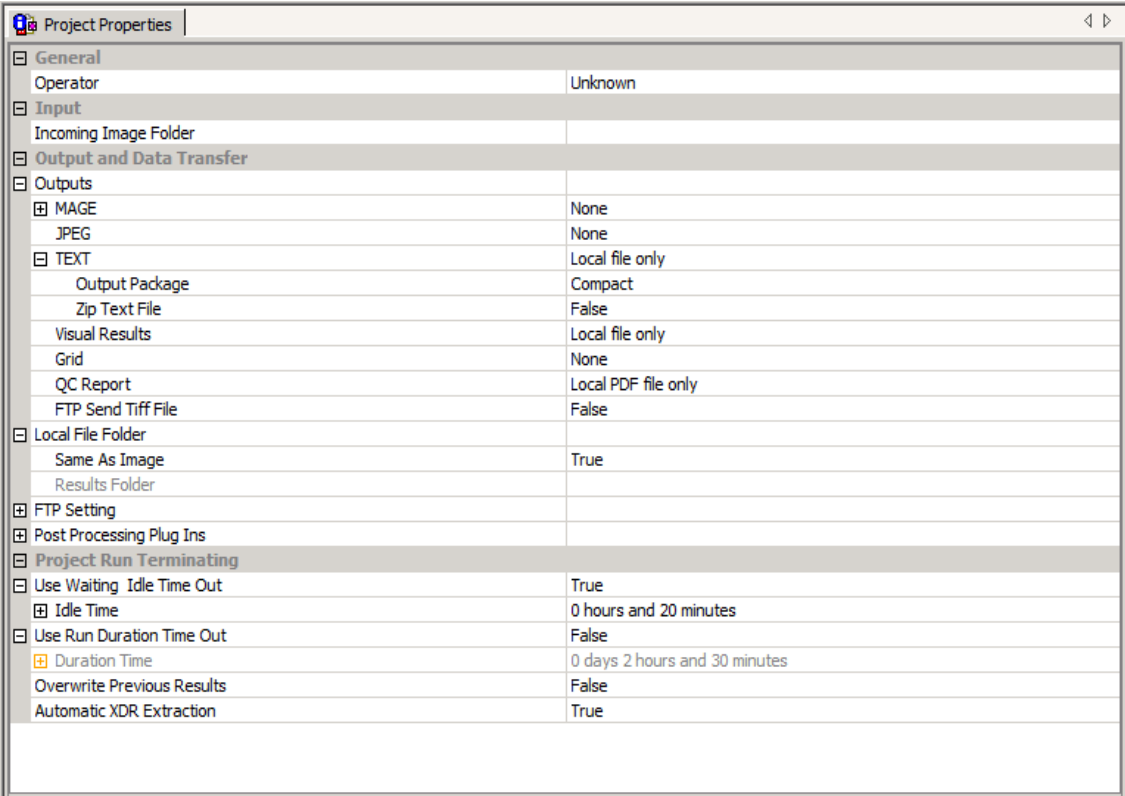


Figure 13 On-Time Project Properties window

Step 2. Enter the incoming image path

- 1** Click the cell next to Incoming Image Folder, and select the path for your scanned images.
- 2** Click **OK**.

The software assigns the appropriate grid template to each of the Agilent microarrays and the appropriate protocol included in the Grid Template Properties sheet.

If you want the On-Time project to process microarrays whose grid templates are missing from the database, you must set up the software to use the eArray server during extraction. The required grid templates, protocols, and QC metric sets are automatically downloaded and installed in the database. See [“To set up eArray login”](#) on page 15 for more details.

Step 3. Enter termination criteria

- 1** Click the cell next to Use Run Duration Time Out, and select **True**.
- 2** If not expanded, expand the Duration Time folder.
- 3** Use the sliders to enter 0 for **Days**, 2 for **Hours** and 30 for **Minutes**.

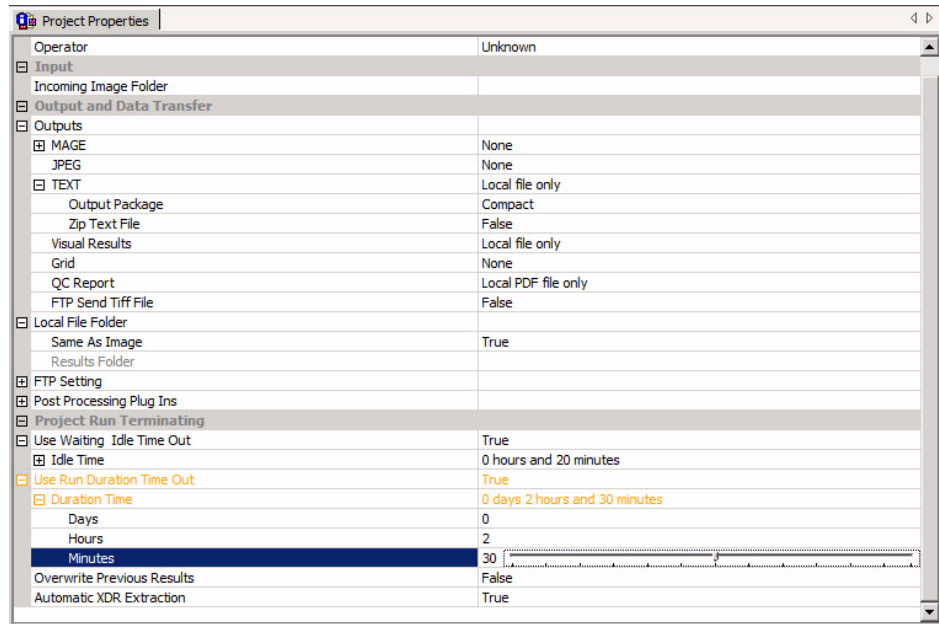


Figure 14 On-Time Project run terminating

Step 4. Change QC metric set associated with a protocol to monitor QC metrics

To change the QC metric set associated to the Feature Extraction protocol:

- 1 In the FE Protocol Browser, double-click a protocol to open the FE Protocol Editor.
- 2 Select the **Protocol Properties** tab to see the protocol properties.
- 3 Under General, click the cell next to **Metricset**.
- 4 Click the down arrow, and select the QC metric set appropriate for your microarray application.

If a metric set you want to use is not available in the metric set list, you must first import it to the database.

To import a QC metric set into the database:

- 1 Select **View > QC Metric Set Browser** from Feature Extraction main window, if it is not already visible.

- 2 Right-click in the QC Metric Set Browser and select **Import** to import either Agilent metric sets or metric sets created with the QC Chart Tool.

The default QC metric sets installed with the Feature Extraction software are read-only and can not be removed from the database. But you can remove a QC metric set from the protocol.

To remove a QC metric set associated with the protocol:

- 1 Select **<Set empty>** from the metric set list in the Protocol Properties.
- 5 Click **Save As** to save the new protocol.

Step 6. Save the project

- 1 Select **File > Save As**.
By default, your projects are saved in **C:\Program Files\Agilent\Agilent Genomic Workbench Standard Edition <version>\FeatureExtraction\FEProjects**.
- 2 To save in a different location, go to or create the folder for your projects.
- 3 Type a project name, and click **Save**.

Step 4. Run Feature Extraction

- 1 Make sure that the scanner is ready.
- 2 Click the **Project Run mode On/Off** icon in the Feature Extraction Project Tools, or select **Project > Start Extracting**.

The image files are saved to the Incoming Image Folder as they complete scanning and are then extracted automatically.

You can display the progress of the extractions, the Summary Report, and the QC Report as you did for the Standard Project extractions.

The Summary Report notes how many metrics are within range for each extraction. The QC report for each image contains the values for the metrics and an evaluation of whether they were within or exceeded the threshold limits.

In this book

This book contains brief instructions to help you get started with your Agilent Feature Extraction 10.10 software. This book shows you:

- What is new for version 10.10
- How to set up and run batch extractions of existing Agilent and non-Agilent image files
- How to set up and run batch extractions of Agilent image files in real time

© Agilent Technologies, Inc. 2015

Revision A1, October 2015



G4460-90035



Agilent Technologies