

GeneChip™ miRNA 2.0 Assay Software Module Version 1.0

User Guide

Information in this document is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information

These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

Corporate entity

Life Technologies | Carlsbad, CA 92008 USA | Toll free in USA 1.800.955.6288

Trademarks

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified. All other trademarks are the property of their respective owners.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.

P/N 703073

Contents

Chapter 1 GeneChip™ miRNA 2.0 Assay Software Module, Version 1.0	3
Introduction	3
Nomenclature	3
Major Features of miRNA 2.0 ASM v1.0	3
Reagents	3
Installing the Assay Software Module	4
Installing a Certificate	4
Installing the miRNA 2.0 ASM v1.0	5
Creating the Test Request	12
Adding Additional Information to the Specimen	13
The Assay Information Screen—Adding Reagent Information	15
Batching Additional Test Request Information	20
Gridding Manually	22
Transferring Data	23
Reviewing the Test Report	24
Ordering Information	27

Chapter 1 GeneChip™ miRNA 2.0 Assay Software Module, Version 1.0

Introduction

This user's guide shows you how to create the GeneChip™ miRNA 2.0 Assay Software Module v1.0 RUO assay test request. This module is for the GeneChip™ miRNA 2.0 Assay that is run on the GCS3000Dx v.2 system using the Molecular Diagnostic Software (AMDS).

This assay software is compatible with the GCS3000Dx v.2 system and the GeneChip™ miRNA 2.0 Array Reagent Kits.

Nomenclature

The assay name is "GeneChip™ miRNA 2.0 Assay" The assay display name is "miRNA 2.0 ASM v1.0" The term ASM refers to Assay Software Module.

Major Features of miRNA 2.0 ASM v1.0

The major features of the miRNA 2.0 ASM v1.0 are four screens that augment assay record keeping. This user guide will also discuss the use of these screens in setting up the assay. These screens include:

1. Additional Information screen
2. Assay Home screen (or Assay Landing screen)
3. Batch Edit screen
4. Report screen

In addition, the miRNA 2.0 ASM v1.0, under the control of the AMDS application, can transfer all specimen information to the server.

It is not the purpose of this short user's guide to instruct you on how to run the assay. This user's guide will show you how to create the RUO test request. To complete the assay run, you will still need to follow the standard AMDS assay protocols and process it through the workflow to register, hybridize, wash/ stain, and scan the assay (as part of the array cartridge). The details of which are discussed in the *Molecular Diagnostic Software User Guide* (P/N 08-0261).

You must be thoroughly familiar with the information contained in the *Molecular Diagnostic Software User Guide*, the *Molecular Diagnostic Software Quick Reference Card* (P/N 08-0262) and the *FlashTag™ Biotin HSR Labeling Protocol* before you can use the information contained in this supplement to run the assay.

Reagents

Reagents for the GeneChip™ miRNA 2.0 Array (single array P/N 901752, 2 arrays P/N 901753, Box of 6 arrays P/N 701754, Box of 30 arrays P/N 901755) comprise the following reagent sub-kits and associated part numbers:

- FlashTag™ Biotin HSR RNA Labeling Kit (10 rxn) - P/N: HSR10FTA
- FlashTag™ Biotin HSR RNA Labeling Kit (30 rxn) - P/N: HSR30FTA
- GeneChip™ Hybridization, Wash, and Stain Kit - P/N: 900720
- GeneChip™ Eukaryotic Hybridization Control Kit - P/N: 900454

Installing the Assay Software Module

To create a miRNA 2.0 ASM v1.0 test request, you must first install a certificate, second, install the miRNA 2.0 ASM v1.0, and finally enter the test request.



NOTE: In most cases, a field service technician will install the certificate and assay software module.

Installing a Certificate

As a security measure, AMDS requires that you or the field service technician install a SSL server certificate on your local workstation in order for your workstation to communicate with the server. This is required for the transfer of the miRNA 2.0 ASM v1.0 test request data to the server.

You should have the certificate installed before installing the assay.

The service technician should have installed a certificate at the time of the system's installation. If for some reason the service technician did not install a certificate or if the server has changed after the system's initial installation, you must install or reinstall a certificate.

If a proper certificate has not been installed, you can still install the assay and process test requests; however, you will have no permission to access the server, and you cannot transfer test request data to the server.

If you cannot access the server and have already installed a certificate, contact Thermo Fisher technical support.

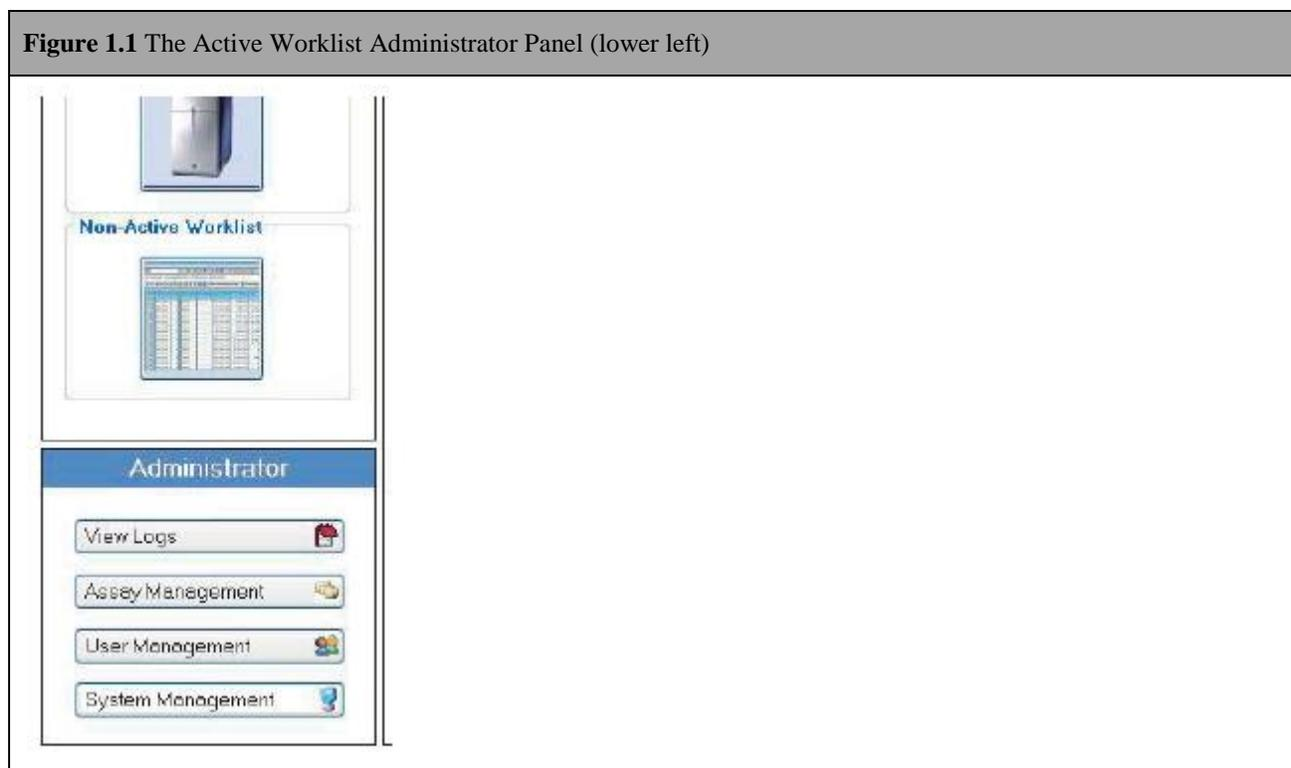


NOTE: You must be logged in with either Laboratory Supervisor or System Maintainer privileges to install a certificate.

Procedure for Installing a Certificate

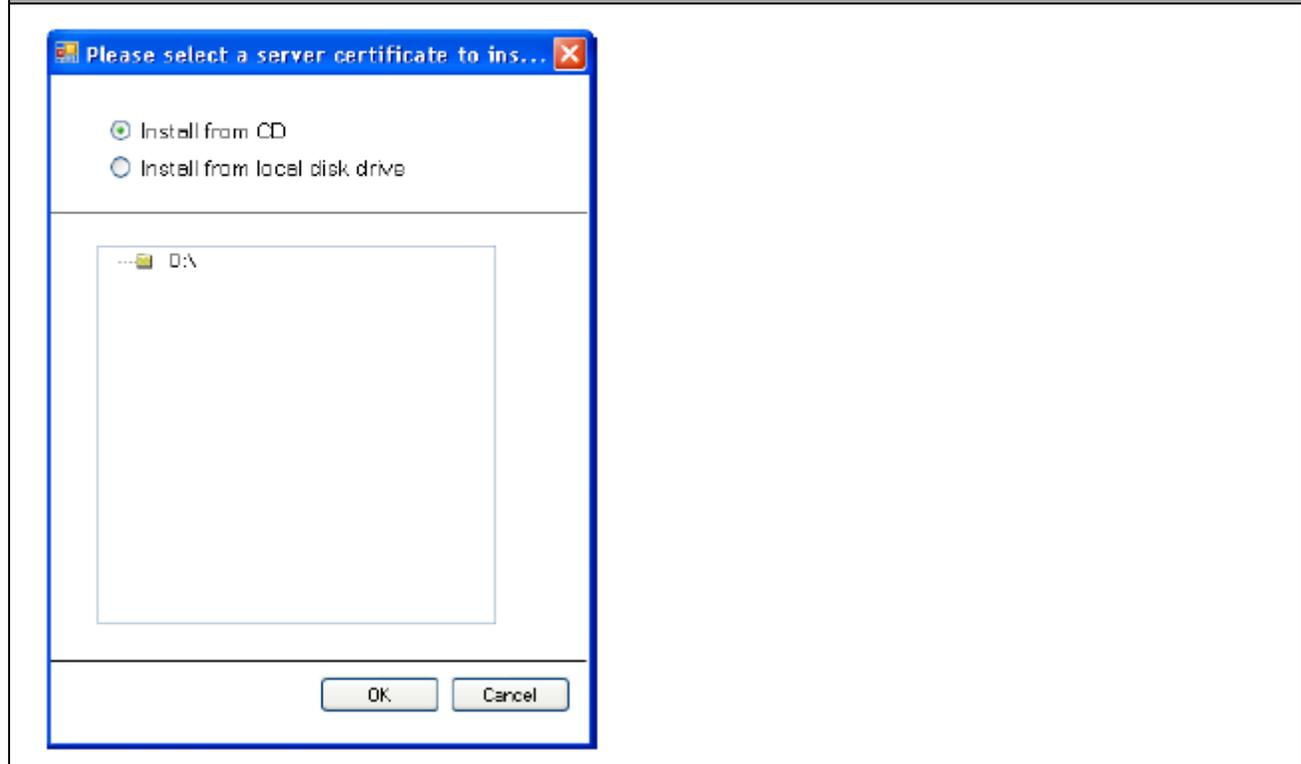
The Active Worklist Administrator panel contains the **Assay Management** button and provides the starting point for installing a certificate (Figure 1.1).

Figure 1.1 The Active Worklist Administrator Panel (lower left)



1. Click on the **Assay Management** button. The Assay Management window appears.
2. Click the **Install Certificate** button. The server certificate browse dialog box appears (Figure 1.2).

Figure 1.2 The Server Certificate browse dialog box: note that the CD/DVD drive is the default drive.



3. Select either the default drive, **Install from CD** (i.e., the internal drive D), or **Install from local disk drive**.
4. Browse for your desired xxxxx.cer file; where xxxxx is the name of the certificate that the installer of the RUO server provided to you.
5. Click the **OK** button or the **Cancel** button.
6. If successful, you are notified by a message “You have successfully added the server certificate “xxxxx.cer” to the AMDS trusted certificate store.”

If the certificate installation fails, AMDS notifies you with the particular algorithm error message. The software prevents you from accessing the server and from transferring data without the proper certificate installed.

7. Click the **OK** button on the message to return to the Assay Management screen.

Installing the miRNA 2.0 ASM v1.0

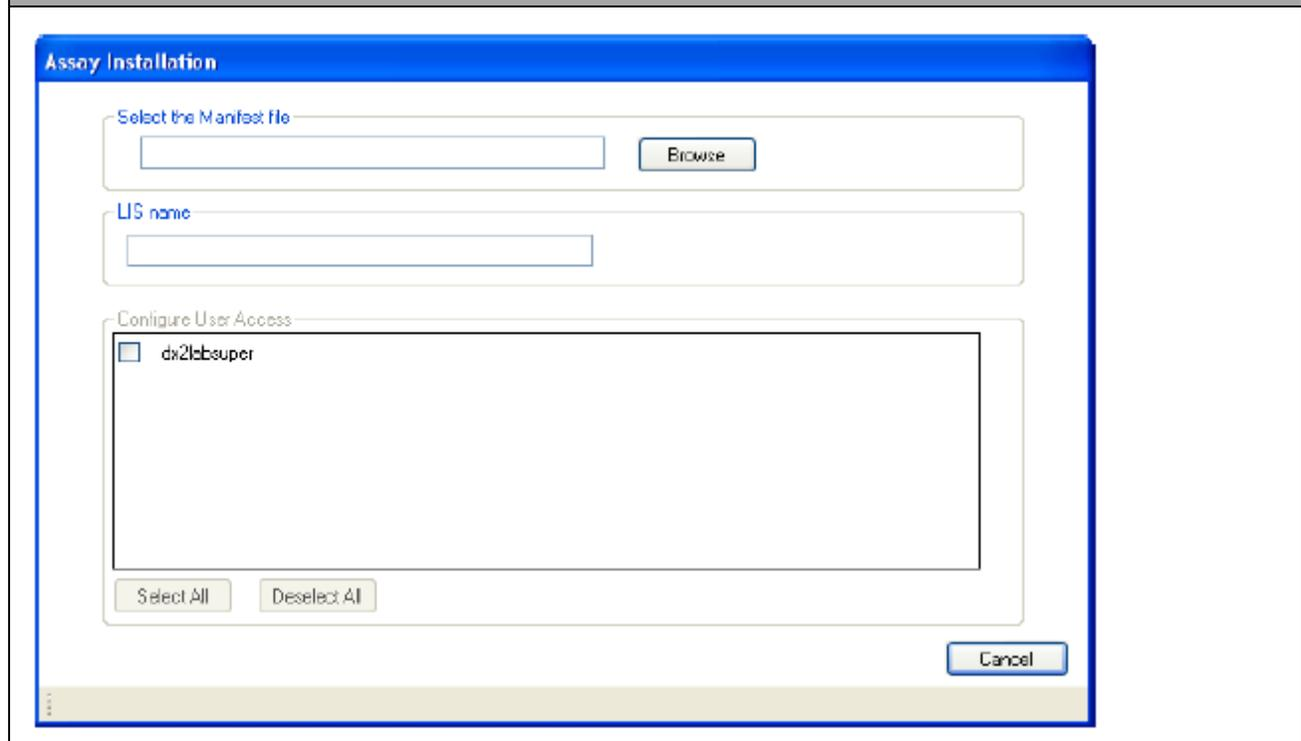
The assay installation process is relatively simple. It requires the selection of an assay manifest file and the selection of approved user access.



NOTE: You must be logged in with either **Laboratory Supervisor** or with **System Maintainer** privileges to install an Assay Software Module.

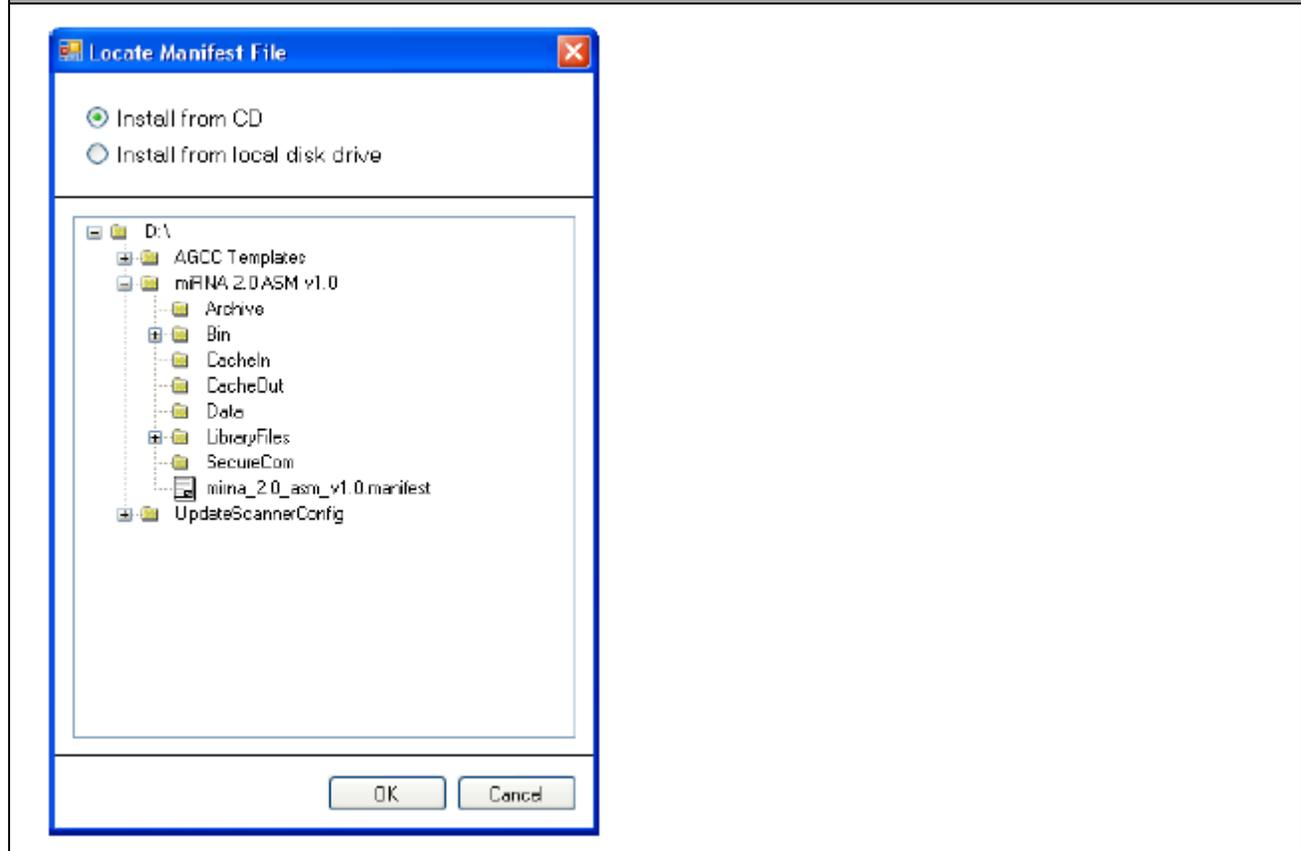
1. Click on the **Assay Management** button. The Assay Management window appears.
2. Click the **Install Assay** button. The Assay Installation browse dialog box appears (Figure 1.3).

Figure 1.3 The Assay Installation browse dialog box



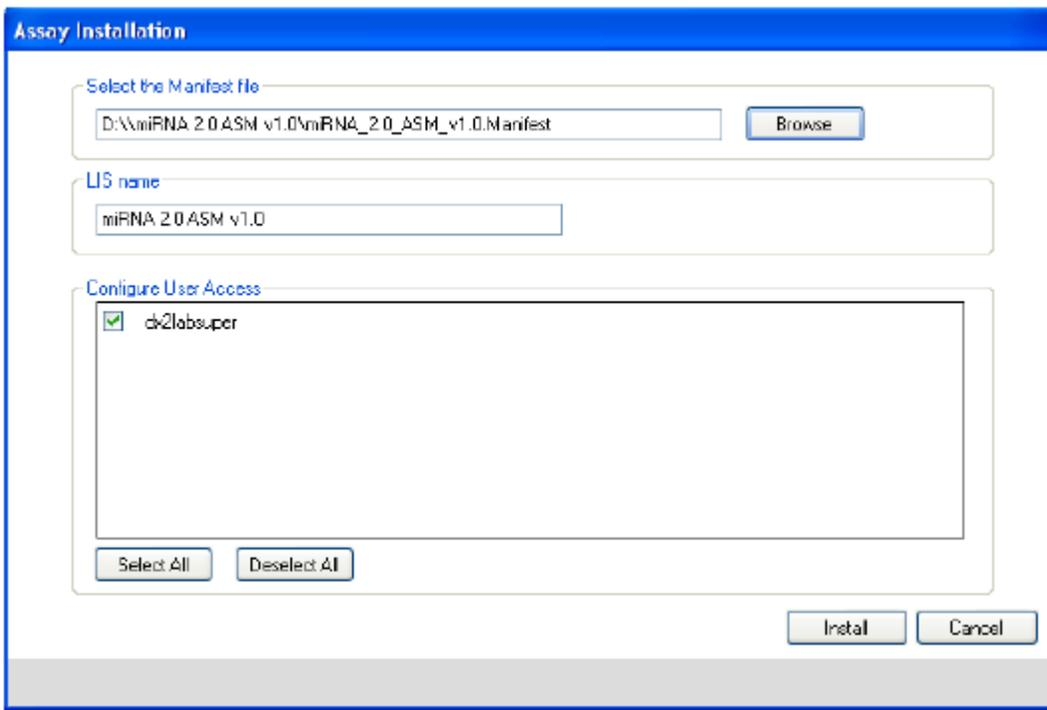
3. Click the **Browse** button. The Assay Installation browse dialog box appears (Figure 1.4).

Figure 1.4 The Assay Installation browse dialog box: note this figure shows internal CD drive D:



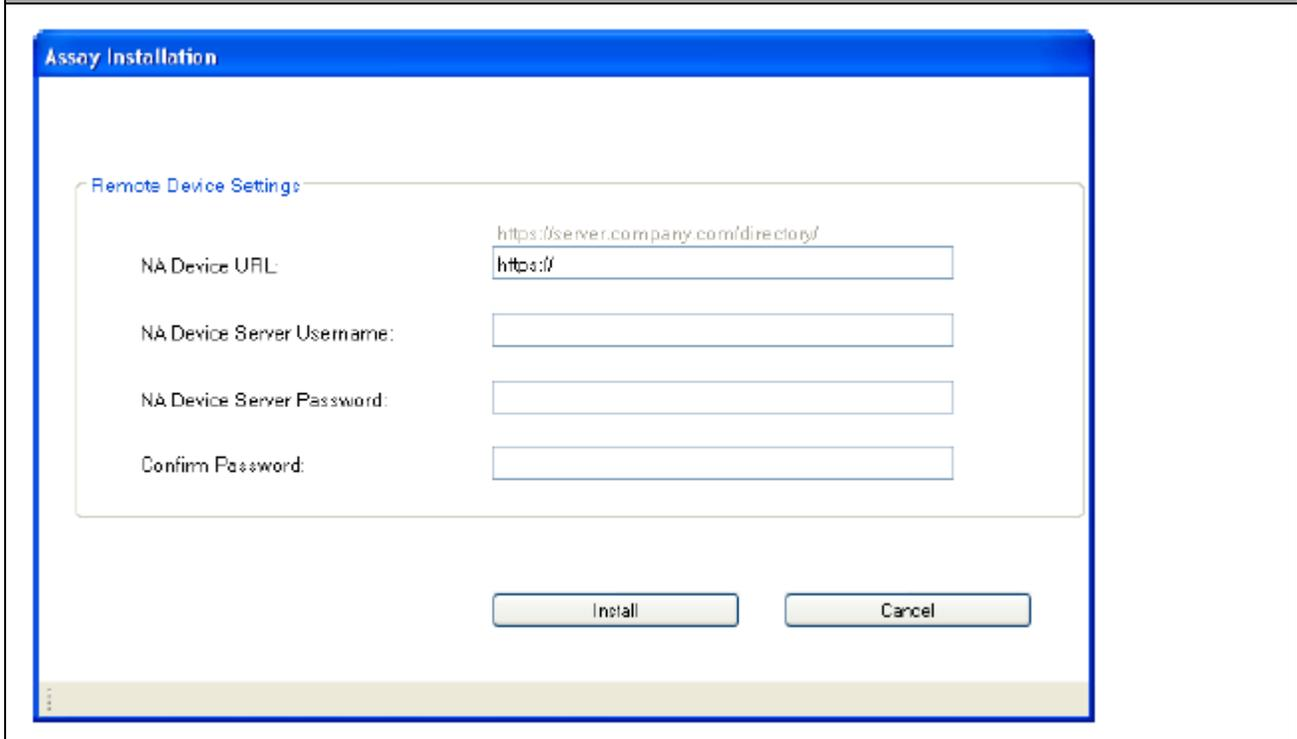
4. Select either the default drive, **Install from CD** (i.e., an internal CD drive D), or **Install from local disk drive** (i.e., the C: drive).
5. Browse for your desired xxxxx.Manifest file. In the present case, it is the miRNA_2.0_ASM_v1.0.Manifest file. Typically you will access the manifest file from the D: drive. It should be in the folder similar to D:\miRNA 2.0 ASM v1.0. Browse for it and select it.
6. Click the **OK** button. The Assay Installation browse screen appears (Figure 1.5).

Figure 1.5 The Assay Installation browse screen for the miRNA 2.0 ASM v1.0



7. Choose those users who will have access to this assay either by selecting individual users, or by clicking the **Select All** button, or if you desire to start the selection over, click the **Deselect All** button.
8. Click the **Install** button or the **Cancel** button.
9. If you clicked the Install button, the Data Transfer Server credentials dialog box window appears (Figure 1.6).

Figure 1.6 The Data Transfer Server Credentials dialog box



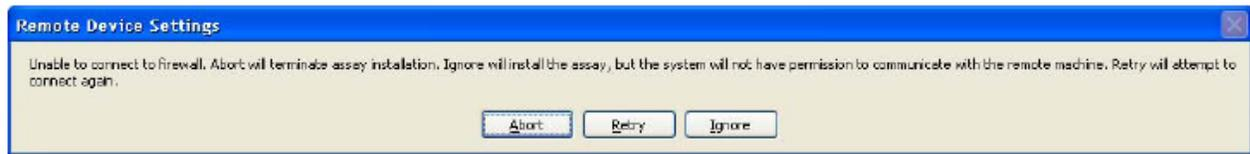
10. Locate and enter the **NA Device URL**, i.e., the secure URL server where AMDS will send the completed assay data (as an example: <https://dx2webdavserver/RUO>).
11. Enter the **NA Device Server Username**. This is your user name for the server, usually your workstation user name or network user credentials.
12. Enter the **NA Device Server Password**. This is your password for the server, usually your workstation user password or network user credentials.
13. Confirm your password.



NOTE: Your NA Device Server Username and Password are not the user name and password that you used to log into AMDS. You must enter your user name and password that you use to access the server.

14. Confirm that the firewall is on and that your workstation is connected to the network.
15. Click the **Install** button or the **Cancel** button.
 - If you clicked Install and if AMDS accepts your user ID and password, the system checks for the appropriate firewall parameters. The system also checks the proper server credentials, and displays an error message if it detects the wrong credentials. If the system detects the proper firewall setup, it accepts the assay installation and returns to the Assay Management screen. Setup is done. You may now begin to use the assay software module.
16. If the system does not pass your system's firewall settings, the following error message is displayed (Figure 1.7).

Figure 1.7 Error Message



17. Click the **Abort** button to exit the assay installation procedure.

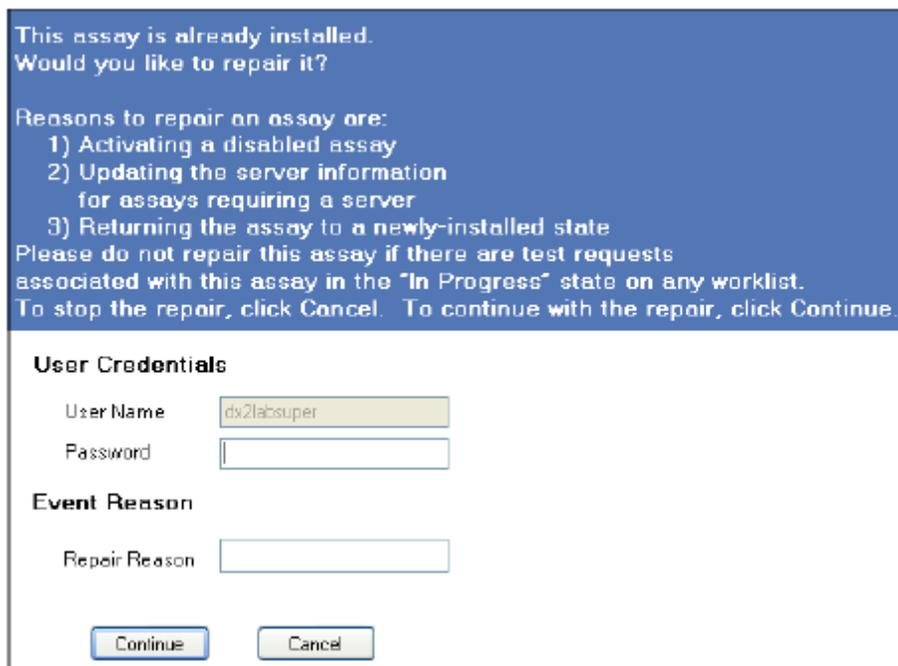
- or click the **Retry** button to attempt another try.
- or click the **Ignore** button to continue.



IMPORTANT: If you select **Ignore**, the AMDS will still install the assay, but you will not have permission to communicate with any remote server.

18. If the same assay already exists in the system, a message appears asking if you would like to repair the assay (Figure 1.8).

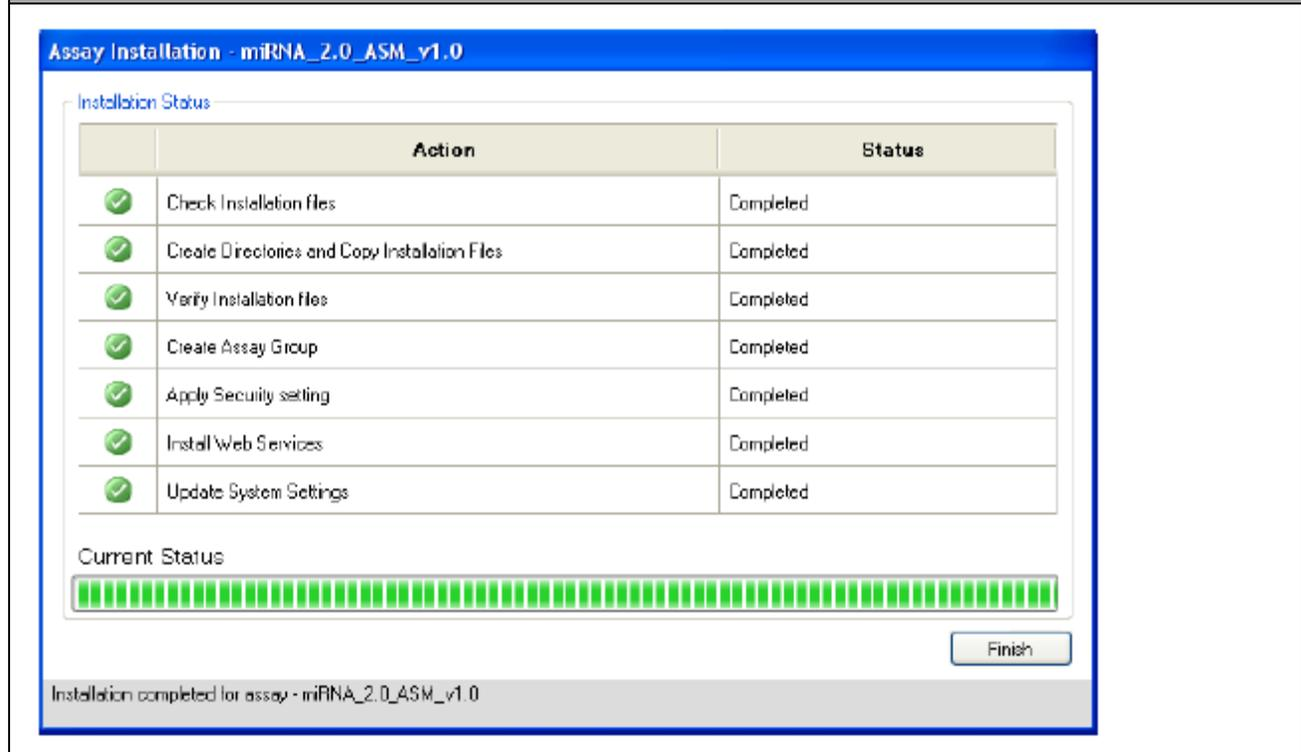
Figure 1.8 The Repair Assay dialog box



- Any AMDS user with valid AMDS credentials and with the right permissions can repair, or reinstall, the assay. You can include a reason, but this is optional.

19. Click the **Continue** button or the **Cancel** button. If you clicked **Continue**, an Assay Installation status window appears and displays the progress of the assay installation (Figure 1.9).

Figure 1.9 The Assay Installation status window



- If AMDS detects no errors, it enables the **Finish** button.

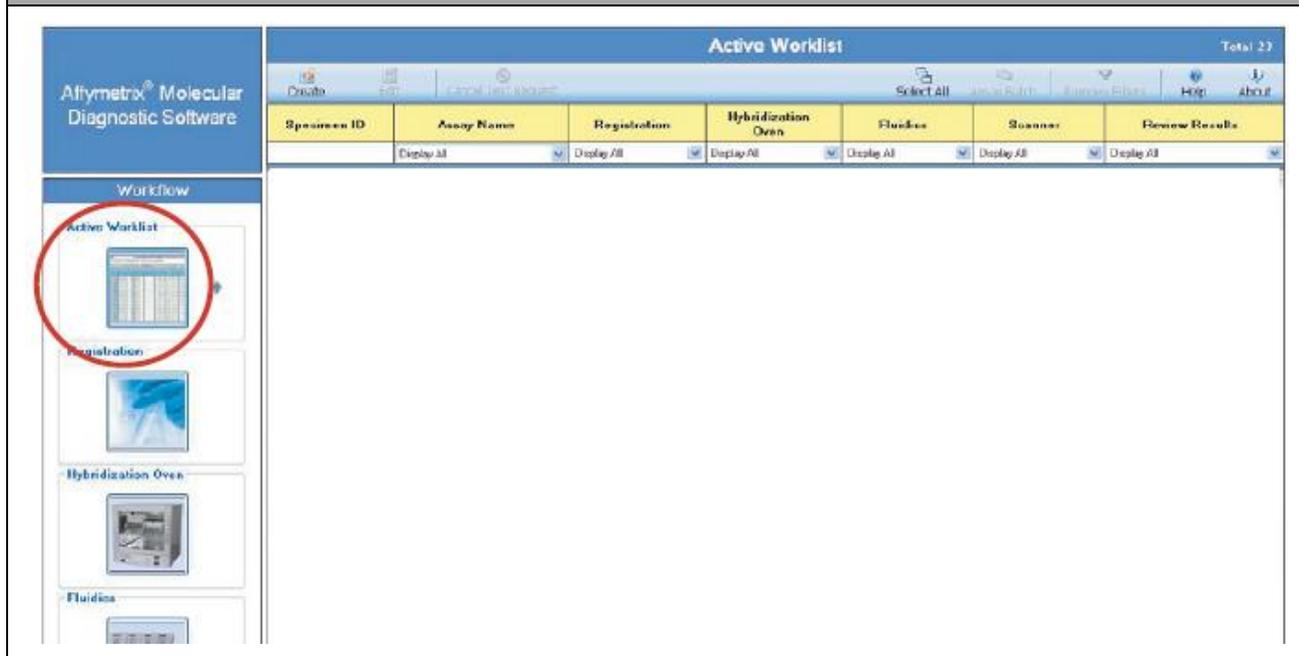
20. Click the **Finish** button to conclude the assay installation procedure and return to the Assay Management screen.

Creating the Test Request

You must be a Laboratory Supervisor, Technician or a Technologist in order to create a test request.

1. If you are not already in the Active Worklist window, select **Active Worklist** from the Workflow panel on the left (Figure 1.10). The Active Worklist panel opens.

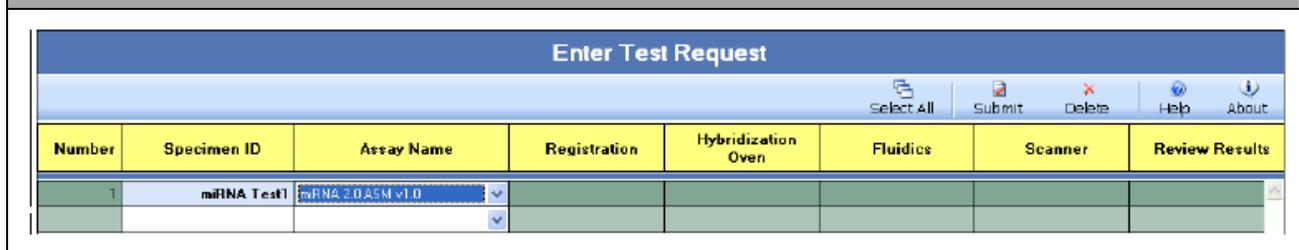
Figure 1.10 The Active Worklist window



2. Click the **Create** button on the toolbar of the Active Worklist.

The Enter Test Request Screen appears (Figure 1.11).

Figure 1.11 The Enter Test Request window



3. Enter one or more **Specimen ID** and **Assay Name** combinations. The current Assay Name for the GeneChip miRNA 2.0 ASM v1.0 is miRNA 2.0 ASM v1.0.
4. Once you have made the entries, click the **Submit** button and the Enter Test Request screen closes. The software displays the Active Worklist window with the newly created test request(s) on the screen.



NOTE: After you have submitted the test request, the procedures for Registration, Hybridization, Fluidics Station processing and Scanning are the same as those steps outlined in the *Molecular Diagnostic Software User's Guide* (P/N 08-0261). Refer to that document for further instructions.

Adding Additional Information to the Specimen

After you add a specimen, and return to the Active Worklist, you can add certain types of pertinent information about the specimen. You can also add this information later here, or in other worklists. This information is not necessary in order to run the assay.



IMPORTANT: You cannot associate any additional information with test requests after scanning has started. You will not be able to save your edits. If you attempt to save, the software displays an error message.

Saving and editing specimen information requires an e-signature upon saving.

1. Click on the desired **Specimen ID** field. The Additional Information dialog box opens (Figure 1.12).

Figure 1.12 The Specimen ID Additional Information dialog box

Field Name	Field Values
Species	Human
Patient ID	G55507190
Patient Last Name	Lynn
Patient First Name	Wilson
Sex	Male
Date of Birth	2011-04-07
Specimen Type	Blood
Specimen Size Unit	mL
Specimen Size Value	3
Method of Collection	venipuncture
Collection Date	2011-05-11
Collection Time	4:33:00 PM
Requestor Last Name	Clark
Requestor First Name	Sydney
Requesting Institution	Stanford Hospital
Request Date	2011-05-11
Additional info 1	
Additional info 2	
Additional info 3	
Additional info 4	
Additional info 5	

2. Add the following information. This information includes:
 - Species - manually entered as freeform text.
 - Patient ID - manually entered as freeform text or read from a barcode.
 - Patient Last Name - manually entered as freeform text.

- Patient First Name - manually entered as freeform text.
- Sex-chosen from a drop-down menu. The menu includes the following choices:
 - Male
 - Female
 - Unknown
 - blank field
- Date of Birth - manually entered with defined format. The format must be entered in an ISO 8601 format: that is with 4-digit year first, then month, then day. (e.g., 2001-11-09 for November 9, 2001) No other format is allowed.
- Specimen Type - chosen from a drop-down menu
 - Blood
 - Solid Tissue
 - Soft Tissue
 - Saliva
 - Buccal Swab
 - Plasma
 - blank field
- Specimen Size Unit - manually entered or from a drop-down menu
 - mL
 - μ L
 - oz
 - mg
 - g
 - inch
 - mm
 - cm
 - blank field
- Specimen Size Value - manually entered floating-point numeric value.
- Method of Collection - manually entered as freeform text
- Collection Date - manually entered as an ISO 8601 format date
- Collection Time - manually entered using either a 12- or 24-hour time scale. (e.g., 1:15 pm or 13:15). The software will display the time using 12-hour time scale (e.g., 1:15 pm).
- Requestor Last Name - manually entered as freeform text
- Requestor First Name - manually entered as freeform text
- Requesting Institution - manually entered as freeform text
- Request Date - manually entered with defined format manually entered as an ISO 8601 format date
- Additional Info 1-5 (Five Additional fields-manually entered as freeform text)

3. When you have completed adding the information, click **Save and Close** or just **Close** to exit without saving.
 - When you click **Close** only, the following events may occur.
 - If you made no edits, the screen will close.
 - If you made edits, a pop-up asking **Would you like to save changes?** (with Yes, No, and Cancel) appears.
 - Click **Yes** in the dialog box and the software collects an e-signature, and then saves and closes the screen.
 - Click **No** and the software discards changes and closes the screen.
 - Click **Cancel** and the software returns to the Specimen ID additional information screen with edits preserved.

The Assay Information Screen—Adding Reagent Information

The Assay Information screen, or the Assay Home screen, provides you with a summary of all the specimen information, test request logs and pertinent reagent information for particular assay type.

The Assay Home screen has the following tabs/sub-screens:

- Reagent Information
- Specimen Report
- Test Request Log

In any worklist window, in the Assay Name field, click on the Assay Name.

- The Assay Home (aka Assay Information) window appears (Figure 1.13). Reagent Information screen is the default tab.

Reagent Information Tab

Click the **Reagent Information** tab (Figure 1.13) to view current reagent information.



NOTE: If you select specimen IDs from the list on the left, then scan the reagent kit barcode, the software parses the reagent information and enters the lot number and expiration date into the correct fields for all selected test requests without manual intervention.

Figure 1.13 The Reagent Information tab

Assay Name: miRNA 2.0 ASM v1.0

GeneChip® miRNA 2.0 Assay For Research Use Only.

Bolded Reagent Kit Name indicates multiple old values exist. **Bolded Data** indicates new value to be saved. Enter Expiration Date as YYYY-MM-DD.

Specimen Report | Test Request Log | **Reagent Information**

Test Requests		Reagent Information				
Specimen ID	Registration Date ▲	Reagent Kit Name	Part Number	Manufacturer	Lot Number	Expiration Date
H919800		FlashTag™ Biotin HSR RNA Labeling Kit (10 rxn)	HSR10FTA	Genisphere	7008471	2022-01-01
H919801		FlashTag™ Biotin HSR RNA Labeling Kit (30 rxn)	HSR30FTA	Genisphere	5906242	2024-02-03
H919802		GeneChip® Hybridization, Wash, and Stain Kit	900720	Affymetrix	3981867	2024-12-29
H919804		GeneChip® Eukaryotic Hybridization Control Kit	900454	Affymetrix	1016926	2022-09-17
H919805						
H919806						
H919808						
H919809						
H919809						

About... Save Close



NOTE: Select one or more test requests on the Test Requests list on the left pane. Data is entered only for the test requests that you have selected.

In the miRNA 2.0 ASM v1.0 (array P/N 901752), you can view the following reagent kit information:

- Reagent kit name
 - FlashTag™ Biotin HSR RNA Labeling Kit (10 rxn)
 - FlashTag™ Biotin HSR RNA Labeling Kit (30 rxn)
 - GeneChip™ Hybridization, Wash, and Stain Kit - P/N: 900720
 - GeneChip™ Eukaryotic Hybridization Control Kit - P/N: 900454
- Reagent kit part number information
- Reagent kit manufacturer information

- Reagent lot number
- Reagent expiration date

The Reagent Kit Name, Manufacturer, Part Number, Lot No. and Expiration Date are not required to run the assay.

AMDS provides the ability to transfer the Reagent Kit Name, Manufacturer, Part Number, Lot No. and Expiration Date to the server, along with other test request information.

1. In the Reagent Information tab, select the test requests in the test request list on the left side of the screen. Scan the barcodes from the reagents you used.

The software will enter the corresponding lot number and expiration date into the appropriate fields. Remember you cannot make any edits to a test request after the array associated with that particular test request has started scanning on the GCS3000Dx v.2 scanner.

2. Click the **Save** button.
3. When you click **Close** only, the following events may occur.
 - If you made no edits, the screen will close.
 - If you made edits, a pop-up **Would you like to save changes?** with Yes, No, and Cancel buttons appears.
 - Click **Yes** in the dialog box and the software saves and closes the screen.
 - Click **No** and the software discards the edits and returns to the Reagent Information screen.
 - Click **Cancel** and the software returns to the Reagent Information screen with edits preserved.



NOTE: You can also enter the reagent kit information manually by selecting test requests in the left side of the screen, then placing the cursor in to the lot number and expiration date fields for the appropriate kits and typing in the correct information.

The miRNA 2.0 ASM v1.0 remembers the association between a reagent kit and expiration date. Scenarios are as follows:

1. Enter a lot number and expiration date for the first time, click **Save**.
 - AMDS saves the data.
2. Enter a lot number already associated with another test request. AMDS auto-populates the expiration date that you had previously used.
 - Click **Save** and the data is saved for all associated test requests.
3. Enter a lot number already associated with another test request. AMDS auto-populates the expiration date that you had previously used. Modify expiration date and click **Save**.
 - The software will display a pop-up with the following message (or words to this effect): “Do you want to save this expiration date for all other test requests with this lot number?”
 - Click **Yes**, **No** or **Cancel**.
 - Yes saves this information for all test requests,
 - No discards all changes.
 - Cancel returns to the screen with edits preserved.

Special Note Regarding the Changing of Expiration Dates

If you change the expiration date for a previously entered lot number (see step 3 above), and choose to save this change for all the test requests with this lot number, the application displays the following message:

“You have changed the expiration date for <reagent kit name> Lot <lot number> from <old date> to <new date>.”

“This affects <count> Test Requests.”

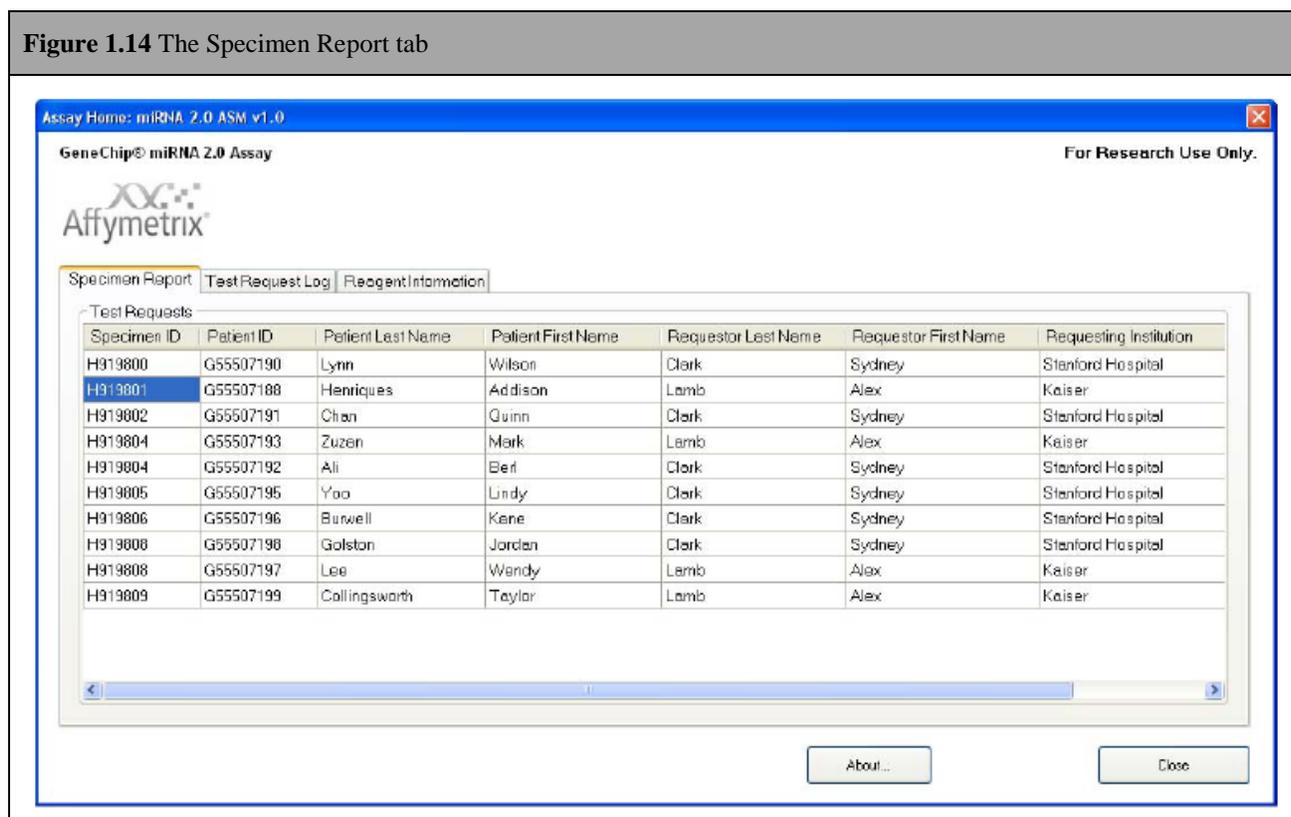
The application remembers the new expiration date and updates all the test requests with those lot number(s), which have not yet started scanning, with the correct expiration date(s).

If you click **Cancel**, the software returns to the Reagent Information Screen with all edits preserved.

Specimen Report Tab

Click the **Specimen Report** tab (Figure 1.14) to view all the specimen IDs associated with that assay type (in this case the GeneChip miRNA 2.0 Assay Software Module v1.0).

Figure 1.14 The Specimen Report tab



Click the **Specimen Report** tab to view all specimen IDs associated with that assay type. You can view the:

- Specimen ID
- Patient ID
- Patient Last Name
- Patient First Name
- Requestor Last Name
- Requestor First Name
- Requesting Institution
- Request Date

Test Request Log Tab

Click the **Test Request Log** tab (Figure 1.15) to view all the test requests associated with that assay type. Each test request will include information regarding the following:

- Date (of the log entry)
- Time (of the log entry)
- User (user logged in when the log entry was created)
- Type (of log entry)
- Subsystem (associated with log entry)
- Short Message (associated with the log entry)
- Long Message (related to the selected short message)

Figure 1.15 The Test Request Log tab

Assay Home: miRNA 2.0 ASM v1.0

GeneChip® miRNA 2.0 Assay For Research Use Only.

Specimen Report Test Request Log Reagent Information

Specimen ID	Date	Time	User	Type	Subsystem	Short Message
H919800	8/25/2011	5:00 PM	System	Workflow	WebService	State Change to Create dState
H919801	8/25/2011	5:06 PM	3173dd24-f38b-...	Workflow	GUI	Successfully saved test request info
H919802	8/25/2011	5:06 PM	3173dd24-f38b-...	Workflow	GUI	State Changed to ReadyForAnalysis
H919804	8/25/2011	5:14 PM	3173dd24-f38b-...	Workflow	GUI	Successfully saved test request info
H919804	8/25/2011	5:14 PM	3173dd24-f38b-...	Workflow	GUI	State Changed to ReadyForAnalysis
H919805						
H919806						
H919808						
H919808						
H919809						

Long Message

Saved these [Name|Value] pairs: [Species] [Patient ID] [G55507188] [Patient Last Name] [Patient First Name] [Sex] [Date of Birth] [Specimen Type] [Specimen Size Unit] [Specimen Size Value] [Method of Collection] [Collection Date] [Collection Time] [Requestor Last Name] [Requestor First Name] [Requesting Institution] [Request Date] [Additional info 1] [Additional info 2] [Additional info 3] [Additional info 4] [Additional info 5]

About... Close

Batching Additional Test Request Information

If you want to enter or edit the same information for multiple test requests at the same time, you can use the Batch Edit feature.



NOTE: You cannot associate any additional information with test requests after scanning has started. You cannot save your edits, and, if you attempt to save, the software displays an error message.

1. In the Active Worklist or any of the worklist windows, select your test requests.
2. Click the **Batch Edit** button.
 - The Assay Batch Information dialog box appears (Figure 1.16)

The left column lists the selected Specimen IDs

3. Enter or change in the right dialog box:
 - Sex (of the test request subject)
 - Specimen Type
 - Requestor Last Name
 - Requestor First Name
 - Requesting Institution
 - Request Date
 - Additional Info 1-5 (five user-defined fields)
4. Click the **Save and Close** button to save your changes and close the window.
5. Click the **Close** button. If you did not make any edits, the Assay Batch screen closes.

If you made any edits, a Yes, No, Cancel window appears.

- Click **Yes**, **No** or **Cancel**.
 - Yes saves this information.
 - No discards all changes.
 - Cancel returns you to the Assay Batch screen with edits preserved.



NOTE: You must save this information before you begin the scanning step. This is to ensure the safety of the data. Do not attempt to add or edit this information after scanning.

Figure 1.16 The Assay Batch Information dialog box

Assay Batch Information: miRNA 2.0 ASM v1.0

GeneChip® miRNA 2.0 Assay For Research Use Only.

Affymetrix

Test Requests in Assay Batch

Specimen ID
H919801
H919802
H919805

Bolded Field Name indicates multiple old values exist. **Bolded Value** indicates new value to be saved.
Enter Date as YYYY-MM-DD. Time as HH:MM:SS am/pm.

Field Name	Field Value
Sex	Female
Specimen Type	Blood
Requestor Last Name	
Requestor First Name	
Requesting Institution	
Request Date	
Additional info 1	
Additional info 2	
Additional info 3	
Additional info 4	
Additional info 5	

Save and Close Close

If the selection contains one or more test requests that have either started or completed scanning, when you click the **Batch Edit** button, the application displays the Assay Batch Information screen with the statement **Test Requests have passed scanning step. Data cannot be changed.**

Gridding Manually

You can manually grid your results if an error arises in the automatic gridding function.

Before you attempt to grid manually, you must be familiar with the gridding function referenced in the *GeneChip™ System 3000Dx v.2 User Guide* (P/N 08-0261), the *GeneChip™ Command Console™ 1.0 User's Guide* (P/N 702569), or the *GeneChip™ System 3000Dx (for EU) User's Guide* (P/N 08-0136) or other documentation. The AMDS manual gridding procedure is based on the gridding procedure outlined in these manuals.

A gridding failure for a test request will trigger a manual grid alignment alert. To remedy this gridding failure:

1. Click on the alert that is visible on the right side of the page in the Alerts panel.
2. View the alert, then click the **Resolve** button.
3. Enter your user ID and password. Click **OK**. The *DatImageViewer* window appears.
4. Manually adjust the grid.
5. Click the **Save** button.
6. Close the *DatImageViewer*.

Transferring Data

AMDS automatically transfers the completed assay data to the URL that you set up when you originally installed your assay. The descriptor labels and data that AMDS transfers to the non-AMDS system include:

- The following file types:

- .ARR file
- .AUDIT
- .CEL file
- .DAT file
- .log file
- .MD5 file

The .MD5 file is a checksum file. You can use a utility such as FastSum (Windows™) or md5sum (Linux) to confirm that AMDS has transferred all the files correctly.

- All Specimen Information descriptor labels and their corresponding data fields which may be filled in or not
- All tracked Reagent Information descriptor labels and their corresponding data fields which may be filled in or not

The AMDS confirms that the data has completed transfer successfully by the following two criteria:

- the assay module detects no exceptions upon transferring
- the assay module confirms that every file listed as transferred indeed exists on the server

Once AMDS transfers the files to the server, AMDS will place all the associated files in a folder named miRNA_2.0_ASM_v1.0. This folder resides on the server. All data files are intermingled. Upon transfer, the software assigns file names: test request creation date and time + "_" + 3 digits + "_" + specimenID + original extension (.DAT, .CEL, etc.). This makes it easy to determine which files are associated with a particular test request.



NOTE: Upon successful data transfer, AMDS deletes, from the local AMDS hard drive, the .dat and .cel files for the transferred test requests.

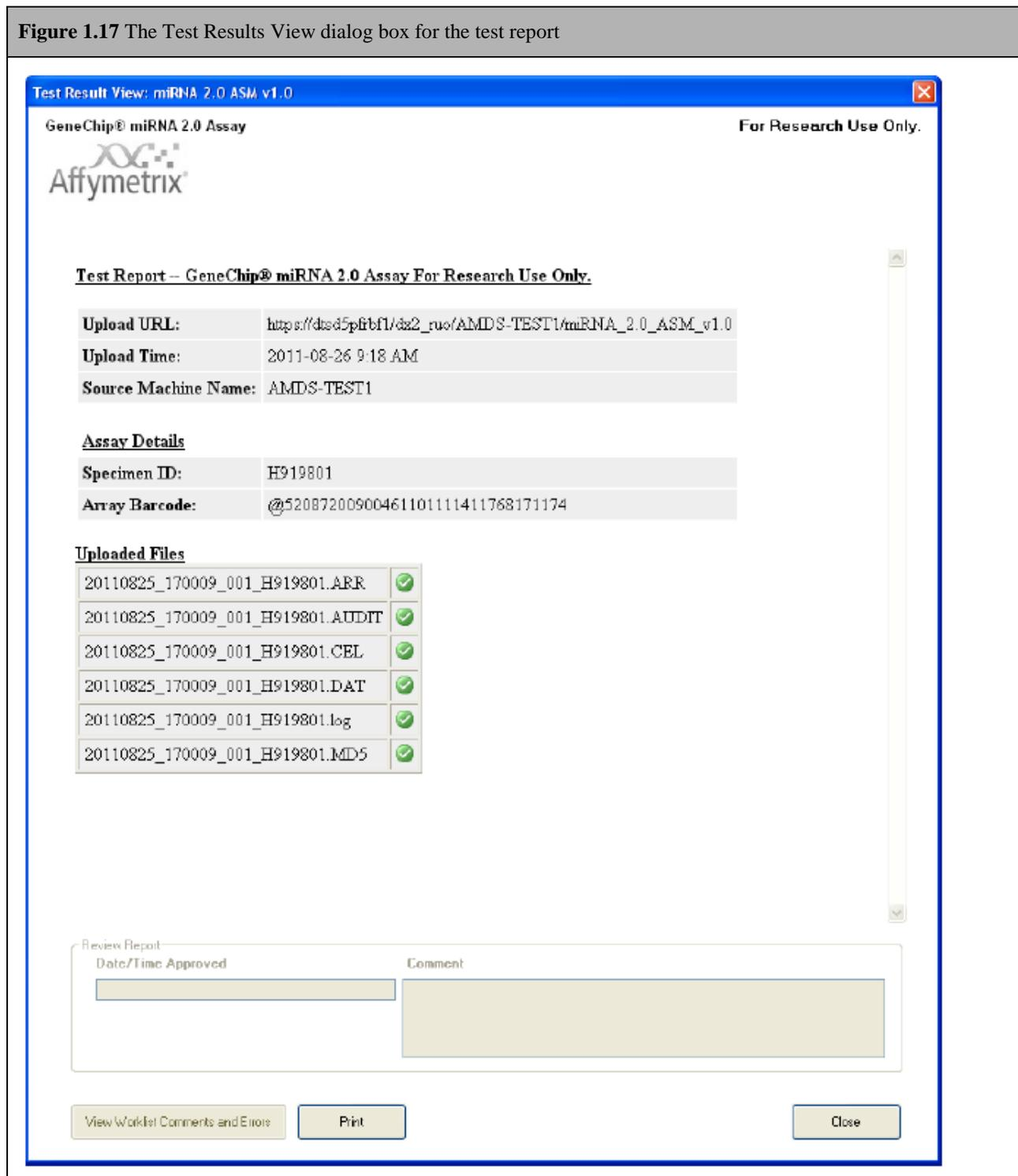
Reviewing the Test Report

When you have completed all the remaining assay steps (Registration, Hybridization, Wash/Stain, Scanning, etc.), the test request moves to the Non-Active Worklist with a hyperlinked date/time stamp in the Review Results field of the Test Request record.

To view the test result report, you must be in the Non-Active Worklist window.

1. Click the date/time hyperlink for the desired test request record. The Test Results dialog box appears (Figure 1.17).

Figure 1.17 The Test Results View dialog box for the test report



The reports displays:

- Upload details
 - Upload URL (where the files went)
 - Upload Time
 - Source Machine Name
- Assay Details
 - Specimen ID
 - Array Barcode
- Uploaded files (the list of the files that were transferred)
 - .ARR file
 - .AUDIT
 - .CEL file
 - .DAT file
 - .log file
 - .MD5 file
- Click the **Close** button to close the results screen without doing anything.
- Click the **Print** button to print the results to the default printer.
- Click the View Worklist Comments and Errors button.
 - The Audit Log Comments and Errors window opens (Figure 1.18). This window displays any comments or errors that the user or AMDS application associated with that test request.

Figure 1.18 The Audit Log Comments and Errors window

Audit Log				
Date/Time	Workflow Step	Type	User	Message
2008-07-14 16:36:23	Other	Error	dz2absuper	The algorithm upload attempt was unsuccessful. Id: 232ada63-707c-4c2a-b50b-ccb5560c025c
2008-07-14 16:25:04	Gridding	Error	dz2absuper	Gridding - Alignment failed. Id: 12842ebd-1430-4ed3-9f17-ds3d0e2e164a

Close

Ordering Information

Table 1.1 lists the miRNA 2.0 ASM v1.0 reagents and associated part number.

Table 1.1

Name	P/N	Supplier
FlashTag™ Biotin HSR RNA Labeling Kit (10 rxn)	HSR10FTA	Genisphere
FlashTag™ Biotin HSR RNA Labeling Kit (30 rxn)	HSR30FTA	Genisphere
GeneChip™ Hybridization, Wash, and Stain Kit	900720	Thermo Fisher
GeneChip™ Eukaryotic Hybridization Control Kit	900454	Thermo Fisher

Documentation and support

Obtaining support

Technical support

For the latest services and support information for all locations, visit **www.thermofisher.com**.

At the website, you can:

- Access worldwide telephone and fax numbers to contact Technical Support and Sales facilities
- Search through frequently asked questions (FAQs)
- Submit a question directly to Technical Support (**thermofisher.com/support**)
- Search for user documents, SDSs, vector maps and sequences, application notes, formulations, handbooks, certificates of analysis, citations, and other product support documents
- Obtain information about customer training
- Download software updates and patches

Safety Data Sheets (SDS)

Safety Data Sheets (SDSs) are available at **thermofisher.com/support**.

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at **www.thermofisher.com/us/en/home/global/terms-and-conditions.html**. If you have any questions, please contact Life Technologies at **www.thermofisher.com/support**.

For support visit thermofisher.com/support or emailtechsupport@lifetech.com

thermofisher.com

23 January 2017

ThermoFisher
SCIENTIFIC