Affymetrix® Laboratory Information Management System (LIMS)



LIMS Manager User's Guide

Version 3.0

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



700339 Rev. 3

Trademarks

Affymetrix[®], GeneChip[®], EASITM, ATM, ATM

Microsoft[®] is a registered trademark of Microsoft Corporation.

Oracle[®] is a registered trademark of Oracle Corporation.

GeneArray® is a registered U.S. trademark of Agilent Technologies, Inc.

Limited License

PROBE ARRAYS, INSTRUMENTS, SOFTWARE AND REAGENTS ARE LICENSED FOR RESEARCH USE ONLY AND NOT FOR USE IN DIAGNOSTIC PROCEDURES. NO RIGHT TO MAKE, HAVE MADE, OFFER TO SELL, SELL, OR IMPORT OLIGONUCLEOTIDE PROBE ARRAYS OR ANY OTHER PRODUCT IN WHICH AFFYMETRIX HAS PATENT RIGHTS IS CONVEYED BY THE SALE OF PROBE ARRAYS, INSTRUMENTS, SOFTWARE OR REAGENTS HEREUNDER. THIS LIMITED LICENSE PERMITS ONLY THE USE OF THE PARTICULAR PRODUCT(S) THAT THE USER HAS PURCHASED FROM AFFYMETRIX.

Patents

Software products may be covered by one or more of the following patents: U.S. Patent Nos. 5,733,729; 5,795,716; 5,974,164; 6,066,454; 6,090,555; 6,185,561 and 6,188,783; and other U.S. or foreign patents.

Copyright

©1999 - 2001 Affymetrix, Inc. All rights reserved.





CHAPTER 1 Getting Started	3
System Requirements	3
Installing LIMS Manager	4
LIMS Manager Overview	11
Features in Version 3.0	11
Workflow-Based Tracking	12
Data Security	12
LIMS and Affymetrix® Software Applications	13
Summary	15
User Interface Orientation	16
Icons Used in LIMS	17
Tool/Toolbar Items	18
Glossary	18
File and Object Types Across All Tabs	18
Tab-Specific Terminology	19
Documentation	22
Conventions Used	22
Your Feedback is Welcome	24
Technical Support	24

CHAPTER 2 Process Tab	27
Process Tab Window	27
Process Tab Functions	29
Data Archive	30
Data Unarchive	31
Assume Ownership	32
Delete Process Data	32
Delete Probe Array Type	34
Export	35
Views in Process Tab Window	36
Print an Experiment	38
Register a LIMS Server	39
Unregister a LIMS Server	40
Server Events	40
Server Tasks	40
Task List Commands	42
Find Function	45
Sample Filter	47

CHAPTER 3 Publish Tab

Publish Tab Window	53
Publish Databases	55
Deleting	56
Probe Array Type	56
Publish Data	57
Creating Publish Databases	60
Database Size Considerations	60
Microsoft® SQL Server Publish Database	61
Oracle [®] Publish Database	63

53

CHAPTER 4 Import Tab	67
Import Tab Window	67
Entering and Updating Data	69
Importing Data	70
Selecting an Import File Path	70
Identifying a Sample	71
Associating Files	73
Start Import	75
Edit Sample	75
Remove and Remove All	76
Sample Filter	76

CHAPTER 5 Userset Tab

Userset Tab Window	81
Userset Definition	83
Creating a Userset	84
Modifying a Userset	85
Define Parameter Values	85
Delete a Probe Array Type from a Userset	88
Delete a Userset	89

CHAPTER 6 Template Tab	93
Template Tab Window	93
Templates	95
Defining or Editing Attributes and Values	97
Deactivating and Activating	101
Renaming	103



81

W	Vorking with Attributes	103
	Multi selecting	104
	Inserting	104
	Deactivating	104
	Activating	105
	Copying	105
	Cutting	106
	Pasting	106

CHAPTER 7 Roles Tab	109
Roles Tab Window	109

Creating a Role	111
Defining Membership	115
Deleting Roles	116

APPENDIX A	Print Samples	119
Diagram Only		119

Diagram & Properties	120

APPENDIX B	FAQs	123
LIMS FAQs		123

Ind	ex	127



Chapter 1





Welcome to Affymetrix[®] Laboratory Information Management System (LIMS). LIMS Manager v 3.0 is a stand-alone database application designed to assist the system administrator in managing the data created by Affymetrix[®] Microarray Suite (formerly GeneChip[®] Analysis Suite). LIMS Manager imports and archives probe array data, provides security and allows the system administrator to assign users and set user privileges.

This chapter provides you with the information you need to get started with Affymetrix[®] LIMS Manager, including system requirements and installation procedures. Additionally, an overview of LIMS Manager and other Affymetrix[®] software applications is presented.

System Requirements

To use Affymetrix[®] LIMS Manager your system must meet the following hardware and software minimum requirements:

- 500 MHz Processor
- 128 MB RAM
- Microsoft Windows NT[®] or Windows 2000 Workstation
- Available space on hard drive = 1 GB
- Microsoft[®] Internet Explorer 5.5 with SP2
- Microsoft[®] Data Access Components (MDAC) 2.5 SP1

Installing LIMS Manager

During the installation of Affymetrix[®] LIMS Manager 3.0 previous version(s) are not removed. If installed in the same directory, the installation program overwrites the previous version's files.

After completing the installation of LIMS Manager 3.0, the program asks you to reboot the system. Restart the computer. Upon logging in, the system registers information into the registry and displays a message on screen informing you that the installation has finished.

Note

You must be logged in as administrator to install the LIMS Manager 3.0 software.

LIMS Manager 3.0 is not required to run MAS 5.0 in LIMS mode.

The screen captures depicted in this manual may not exactly match the windows displayed on your screen.

▲ WARNING

If at any time during the installation a dialog box with a red 'X' appears, this indicates that you logged on with a username that does not have the local administrative privileges required to install the software. Log out and log back in under a user name with proper privileges.

- Insert the Affymetrix[®] LIMS 3.0 Client CD-ROM. If the autorun feature does not start the program:
 - a. Click Start \rightarrow Run.
 - **b.** Type *<cd drive letter*>:\setup.exe.
 - c. Click OK.
 - \Rightarrow A splash screen is displayed, followed by the Welcome window (Figure 1.1).



Figure 1.1 Welcome window

2. Click Next.

 \Rightarrow Several consecutive Software License Agreement windows appear.



Figure 1.2 Software License Agreement w

Software License Agreement window



- **3.** Review the contents and click **Yes** in each window to accept the terms of the licensing agreement.
 - \Rightarrow A Question window appears regarding the path location in reference to an upgrade (Figure 1.3).



✓ Note

For upgrades, we recommended that you install LIMS Manager v 3.0 in the same location as the previous version and overwrite the existing files.

- 4. If this is an upgrade you are asked if you want the latest version of LIMS Manager in the same location as the previous version.
 - Click **Yes** to install in defined location.
 - \Rightarrow The files from the previous version are overwritten.
 - Click No to install in new location. In the dialog box that appears, define the new path location desired.
 - \Rightarrow The Choose Destination Location window appears (Figure 1.4).



Figure 1.4 Choose Destination Location window

- Select the destination to install LIMS Manager 3.0. Click Next. C:\GeneChip is the default location.
 - \Rightarrow Files are copied.
- 6. Click Next.
 - \Rightarrow The Select Components window appears (Figure 1.5).



Figure 1.5 Select Components window - configuration disk



- 7. Select Yes if you have the LIMS configuration disk, No if you do not.
- 8. If Yes, go to step 9. If No, go to step 11.
- 9. Browse to the location of LIMS configuration disk.

The file needed is called **GeneChipDB.ini** and resides on the LIMS server. The file is located at <server>\gclims\library. Browse to the file location.

Setup needs the GeneChip LIMS configuration disk			
÷	Please insert the GeneChip LIMS configuration disk. Use the browse button to select the path.		
Path:	Browse	J	
\mathbb{R}	DK Cancel		
Figure 1.6			

Browse to File Location

- 10. Click OK. Go to step 14.
- **11.** Enter the name of the LIMS server and click **Next** (Figure 1.7).



Figure 1.7 Enter Information window - enter LIMS server name

 Enter the name of the domain the server resides on with the global group name that contains the LIMS users. (Domain Name\Global Group) (Figure 1.8).



Figure 1.8

Enter Information window - enter global group name

- 13. Click Next.
 - \Rightarrow The Setup Not Complete window appears.
- **14**. Select **Yes**, **I** want to restart my computer now to complete the installation.





Reboot to continue installation

- **15.** Log in as an administrator to complete the installation.
- **16.** Click **Finish** in the Setup dialog box.

LIMS Manager Overview

Affymetrix[®] LIMS Manager v 3.0 is the software tool that provides you with a convenient way to manage your GeneChip[®] expression data. Improved performance in process database, improved access security throughout and improved query performance are just some of the features new in version 3.0 designed to enhance the management capabilities of LIMS Manager.

Features in Version 3.0

- Improved query performance for better database access
- Improved database management and data security features
- Improved access security throughout, including administrator defined user access level for added data security
- Easily track experiment status
- Ability to import historical data from a local hard drive or an external storage device (ZIP, JAZ, CD-ROM) and store on LIMS server
- Supports experiment and sample attribution
- Support for new data analysis algorithms
- Expanded Oracle[®] database support to include Oracle 8.1.7 both publish and process database
- Expanded Microsoft[®] SQL database support to include SQL 2000 both publish and process database
- Support for Microsoft Windows NT[®] and Microsoft[®] Windows 2000



Workflow-Based Tracking

Affymetrix[®] LIMS Manager software tracks experiments and processes. Its database provides storage and management for expression data.

You can access experimental data from any workstation on the network through the Affymetrix[®] LIMS server. Each user with access to the LIMS system can set up samples, experiments and publish data.

The LIMS Manager software processes experimental data through a workflow-based tracking system. The software tracks the stages of sample registration, experiment setup, hybridization, scan, grid alignment, analysis processes and publishing. This tracking allows you to follow the status of an experiment from sample preparation through analysis.

Data Security

The Affymetrix[®] LIMS software incorporates data security from the Windows 2000 computer. The logon name verifies that the user has access to the Affymetrix[®] LIMS database.

Users can access the Affymetrix[®] LIMS database only if they have been granted authorization. Access is provided to the LIMS server by requesting privileges from the system administrator who is responsible for assigning users' rights and permissions.

LIMS protects users' data from unauthorized changes. A user who creates, generates and analyzes an experiment is the original owner of the data. With security settings defined in the Roles tab, a user who is not the owner of the data will only have read access but is allowed to create new analyses without overwriting the original data. Security settings can also specify that only the owner of the data can overwrite the original data files.

LIMS and Affymetrix® Software Applications

Affymetrix[®] LIMS Manager is the second of three software tools from Affymetrix for gene expression analysis (Figure 1.10).





Use Affymetrix[®] Microarray Suite to create, analyze and publish expression analysis data files. Next, import these files to a server database using LIMS Manager. Finally, access the publish database using Affymetrix[®] Data Mining Tool (DMT) to further analyze the data by filtering, pivoting and graphing. All three products are separate Windows applications available from Affymetrix.





Workstations can contain any of the following software combinations:

- Affymetrix Microarray Suite: instrument control; File Mode: data acquisition and intensity analysis, LIMS mode: Data publishing, retrieval and sample tracking.
- ◆ LIMS: batch processing Access to sample IDs
- ◆ Data Mining Tool: Mining of published data Data visualization using graphs, tables



Summary

The Affymetrix[®] LIMS and Data Mining Tool applications ensure that multiple users have simultaneous access to the database from any workstation. When used together, Affymetrix[®] Microarray Suite, Affymetrix[®] LIMS and Affymetrix[®] Data Mining Tool software applications generate experimental and research data for scientific analysis. Using these powerful tools in conjunction, users are able to review and analyze data in a robust and dynamic manner.



User Interface Orientation

Active Server			Me Tool	enu Bar bar	
LIMS Manager - [LIMS Server Manager] Die View Iools Process Publich Import Userset Template Boles Window INSTALL	Help				_ & ×
Process Publich () Import () Userset () Template () Roles IMS Servers INSTALL Expression Probe Arey Type Expression Process Database					
	Local Task + Export + Import Delete Archive * Create	Task Description	Started 08:02:42 08:02:42 08:02:42 08:02:42 08:02:42 08:02:42	Status Available Available Available Available Available	
For Help, press F1					NUM

LIMS Manager interface

Icons Used in LIMS

Table 1.1

LIMS Manager Icons

A	Sample	Ĩ	Template	r	Properties Information
6 2	Project	۲	Global	2	Gray bullet indicates processes that have been completed.
8	Experiment Information / .exp file	-8-	Server	۲	Green bullet indicates processes waiting to commence.
88	Image of Scanned Probe Array / .dat file		Process Database (yellow)	a /	Red check mark on .exp, .dat, .cel or .chp icon(s) indicates that the file already exists in the Publish Database.
	Averaged Intensities / .cel file	T	Publish Database (gray)	6+	Export Status
	Analysis Output / .chp file	0+	Import to Database	6+	Import Status
	Probe Array type	Þ	Domain	N	Delete Status
	Chip	Q	User Group	H	Archive Status
Ú	Vessel	d,	User	6*	Create Publish Database Status
ĒĒ	Userset	2 2	Role	T	Filter



Tool/Toolbar Items

Most toolbar or **Tool** menu items are available across all tabs. Detailed descriptions of these commands are presented in subsequent sections of this manual.



Glossary

The following terms are used frequently throughout this manual. For your convenience the terms are listed under the tabs where they are most frequently found.

File and Object Types Across All Tabs

Automate	Automate is a monitored task whose status, either on or off, is viewed in the Local Task pane. Automate On indicates LIMS is automatically tracking an experiment through data analysis after a scan of the probe array has successfully completed.
.cel	The .cel averaged intensity file. Also known as "cell" file, this file is automatically created after grid alignment. The cell averaging analysis calculates the average intensities of each cell and assigns it an x and y coordinate, and can be used to re-analyze data with different algorithm patterns. The .cel file is one of two file types you can migrate to a publish database using Affymetrix [®] Microarray Suite.
.chp	Also known as a "chip" file, this analysis file is the output generated by the analysis of a .dat and .cel file. The .chp is the second of two file types you can migrate to a publish database using Affymetrix [®] Microarray Suite.

.dat	The .dat file is the image of the scanned probe array and is created by the analysis.
.exp	The .exp file is created by the user and contains information defined by the experiment template used. The experiment name becomes the file name for subsequent files generated in the analysis. The Import function presents the concept of moving files and on this tab there is an experiment (.exp) file. The process database contains files and data, but does not contain experiment files; only experiment and sample data.
Sample	A set of experiments.

Tab-Specific Terminology





📔 Process 🖺	Publish 🗇+ Import 🔛 Userset 🛛 🝸 Template 🗍 🕵 Roles 🗌			
Affymetrix® Analysis Data Model	The Affymetrix [®] Analysis Data Model is the relational database schema Affymetrix uses to store experiment data so that results may be filtered and mined by analysis tools. Previously referred to as the GATC data model, the Affymetrix [®] Analysis Data Model includes additional tables to support mapping, spotted arrays and new expression results ¹ .			
	The Affymetrix [®] Analysis Data Model is publicly available to support open access to experiment information generated and managed by Affymetrix [®] software.			
Intensities	See .cel file on page 18.			
Publish Database	An Affymetrix [®] Analysis Data Model database residing on a server containing select process data that have been published. You may have several publish databases on a single server.			
Publish	Publishing is the act of migrating sample data to a Microsoft [®] SQL or Oracle [®] publish database and is a function of Microarray Suite v 4.0 or above.			
×				



1 Affymetrix® LIMS Manager 3.0 does not support mapping or spotted array data.

👔 Process 👔	Publish 🗇 Import 🧮 Userset 🖷 Template 🕼 Roles		
Userset	A userset is a predefined prarameter set used when running an expression analysis.		
Independent Parameter	Defined in the userset, independent parameters are static settings not influenced by probe array selection.		
Dependent Parameter	Defined in the userset, dependent parameters are settings that are specific to the particular probe array type selected.		



	process. The registration process is a function of Affymetrix [®] Microarray Suite.
Deactivate	On the Template tab, the Deactivate command does not permanently remove an existing template or attribute. The deactivate action only

renders the item inactive. The item does not appear in Affymetrix[®] Microarray Suite until the **Activate** command is used.





Documentation

This manual provides a detailed outline for all tasks associated with Affymetrix[®] LIMS Manager.

Additionally, the LIMS Manager installation CD includes an electronic version of this user's guide. The on-line documentation is presented in Adobe Acrobat[®] format (a .pdf file) and is readable with the Adobe Acrobat[®] Reader software, available free from Adobe at http://www.adobe.com. The electronic user's guide is printable, searchable and fully indexed.

Conventions Used

Various conventions are used throughout the manual to help illustrate the procedures described. Explanations of these conventions are provided below.

Steps

Instructions for procedures are written in a step format. Immediately following the step number is the action to be performed. On the line below the step there may be the following symbol: \Rightarrow . This symbol defines the system response or consequence as a result of user action; what you see and what has happened that you may not see.

Following the response, additional information pertaining to the step may be found and is presented in paragraph format. For example:

- 9. Click Yes to continue.
 - \Rightarrow The **Delete** task proceeds.

In the lower right pane the status is displayed.

To view more information pertaining to the delete task, right-click **Delete** and select **View Task Log** from the shortcut menu.

Font Styles

Bold fonts indicate names of commands, buttons or options within a dialog box. When asked to enter specific information, such input appears in italics within the procedure being outlined. For example:

- 1. To select another server, enter the server name in the Oracle Alias box.
- Enter DMT_2_Tutorial in the Publish Database box, then click Register.
 - \Rightarrow The tutorial database is available to DMT.

Screen Captures

The steps outlining procedures are frequently supplemented with screen captures to further illustrate the instructions given.

V Note The screen captures depicted in this manual may not exactly match the windows displayed on your screen.

Additional Comments

Throughout the manual, text and procedures are occasionally accompanied by special notes. These additional comment types and their meanings are described below.

TIP	Information presented as tips provide helpful advice or shortcuts for completing a task.
✔ Note	The Note format presents information pertaining to the text or procedure being outlined.
! CAUTION	Caution notes advise you that the consequence(s) of an action may be irreversible and/or result in lost data.
▲ WARNING	Warnings alert you to situations where physical harm to person or damage to hardware is possible.



Your Feedback is Welcome

Affymetrix Technical Publications is dedicated to continually improving the quality of our documentation and helping you get the information that you need. We welcome any comments or suggestions you may have regarding this manual. Please contact us at:

technical_publications@affymetrix.com

Technical Support

Affymetrix provides technical support to all licensed users via phone or e-mail. To contact Affymetrix Technical Support:

Affymetrix Inc.

3380 Central Expressway Santa Clara, CA 95051 USA Tel: 1-888-362-2447 (1-888-DNA-CHIP) Fax: 1-408-731-5441 E-mail: support@affymetrix.com

Affymetrix UK Ltd.,

Voyager, Mercury Park, Wycombe Lane, Wooburn Green, High Wycombe HP10 0HH United Kingdom **Tel:** +44 (0) 1628 552550 **Fax:** +44 (0) 1628 552585 **E-mail:** supporteurope@affymetrix.com

www.affymetrix.com



Chapter 2





The initial view presented upon opening Affymetrix[®] LIMS Manager is the **Process** tab. The **Process** tab shows you the data in the system, where it is found and allows you to view environmental property information regarding the data.

This chapter orients you to the layout of the **Process** tab window and describes in detail the functions and tasks accomplished here.

Process Tab Window

The **Process** tab window is divided into three panes (Figure 2.1). The first pane on the left side displays a tree that identifies the active server, process database and data types (**Samples**, **Projects** and **Probe Arrays**). The data type subfolders contain the **Sample Name**, or **Probe Array Name** files for expression data. The tree stops at this level.

The second pane in the upper right corner takes the **Sample Name** or the **Project Name**, selected in the first pane, copies it as the top node and lists the associated experiment(s) underneath. The first time you click an experiment in this pane, the information is read and a '+' is displayed next to the experiment name. Click the '+' to display processes completed with inputs and outputs, or associated .dat files. Alternatively, double-click an experiment in this pane and the information is read and processes or associated files are displayed. Experiment information in this pane is displayed in either a **Process view** or a **File view**.

The third pane in the lower right corner displays either a Property Information or Local Task status view, depending on the item selected. Property Information lists environment details pertaining to the selected item. This information is defined by the template used during the **Sample Registration** process completed in Affymetrix[®] Microarray Suite. The Local Task status view displays the status of processes that may be occurring in the database, such as **Export**, **Import**, **Delete**, **Archive** and **Create**. This view allows you to monitor these jobs while concurrently working on additional tasks within the LIMS system.



Figure 2.1

Process tab: Process View (2) of selected sample (1) with environmental property information displayed(3)

Process Tab Functions

The primary functions of the **Process** tab include archiving, unarchiving, deleting, exporting, assuming ownership and printing data. These commands are found on the menu bar under **Process**.

The primary data types in the **Process** tab are sample, project, experiment and file. Each data type has associated functions whose commands are found under **Process** on the menu bar, or accessed by right-clicking the data type and selecting from the shortcut menu that appears (Figure 2.2).





Data Archive

Archiving data moves specified files and stores them in an external location. Although archived files no longer exist in the process database, their representative icon remains in the tree and appears gray to alert you that this data is available through the **Unarchive** process.

It is not possible in the current version of LIMS to archive experiments. You may only archive specific data files (.dat, .cel and .chp).

Archiving Data Files

- 1. On the menu bar select **Process** \rightarrow **File View**, if necessary.
- In the upper right pane, click to select the file to be archived. Selecting a parent file also includes all related children files in the archive process.
- **3**. On the menu bar select **Process** \rightarrow **Archive**.

Alternatively, right-click the file to archive and select **Archive** from the shortcut menu that appears.

 \Rightarrow The Select Archive File Path window appears (Figure 2.3).

Select Archive File Path	
Current path:	
c:\program files\affymetrix	
Directories:	ОК
c:\ 🔺	Cancel
🗁 affymetrix	
MAS 🗾	
Dri <u>v</u> es:	
E C:	Network

Figure 2.3 Select Archive File Path window

4. Select a **Drive** from the drop-down list and browse to find desired directory.
5. Click OK.

 \Rightarrow The **Archive** Local Task status is displayed in the lower right pane.

Double-click **Archive** at any time to view additional information pertaining to the archive status, or right-click **Archive** and select **View Task Log** from the shortcut menu that appears.



Data Unarchive

TIP

The icon(s) for data that have been archived lose their color and appear gray. The **Unarchive** command recalls the archived data and restores it to the process database.

Unarchiving Data Files

1. Click to select the file to unarchive.

Selecting a parent file also includes all related children files in the unarchive process.

If the archived file cannot be located a **File Location** dialog appears. Use this dialog box to define the location of the desired files.

On the menu bar select Process → Unarchive, or right-click the file and select Unarchive from the shortcut menu.

✓ Note

The Process window does not automatically refresh upon completion of the **Archive** or **Unarchive** task. Therefore, if you unarchive a data file the icon remains gray even though the task was completed successfully. To refresh the view select **View** \rightarrow **Refresh** from the menu bar.



Assume Ownership

Change the **User** attribute of data to enable the reassignment of data when necessary for security or management purposes. Ownership of samples, experiments or analyses may be assumed when the **Take Ownership Process Data** option is defined for the user in the **Roles** tab.

Assuming Ownership

- 1. Click to select the appropriate sample, experiment or analyses.
- **2.** From the menu bar select **Process** \rightarrow **Assume Ownership**.
 - \Rightarrow Your user name appears as **User** in the properties window in the lower right pane.

OR

- 1. Right-click the appropriate sample, experiment or analyses.
- 2. Select Assume Ownership from the shortcut menu.
 - \Rightarrow Your user name appears as **User** in the properties window in the lower right pane.

Delete Process Data

You may delete samples, experiments, .cel, .chp analyses and usersets. Deleting samples deletes all associated experiments and analyses. Deleting experiments deletes all associated analyses.

! CAUTION

The **Delete** option permanently removes data from the database. There is no restore function to recover deleted data.

Deleting Process Data

- **1**. Click to select the item to delete.
- **2.** From the menu bar select **Process** \rightarrow **Delete**.
 - \Rightarrow A warning / confirmation window appears (Figure 2.4).





- 3. Click Yes to continue.
 - \Rightarrow The **Delete** Local Task status is displayed in the lower right pane.

Double-click **Delete** at any time to view additional information pertaining to the delete status, or right-click **Delete** and select **View Task Log** from the shortcut menu that appears.

OR

- **1**. Right-click an item to delete.
- 2. Select **Delete** from the shortcut menu that appears.
 - \Rightarrow A warning / confirmation window appears (Figure 2.4).
- **3.** Click **Yes** to continue.
 - \Rightarrow The **Delete** Local Task status is displayed in the lower right pane.

Double-click **Delete** at any time to view additional information pertaining to the delete status, or right-click **Delete** and select **View Task Log** from the shortcut menu that appears.

✓ Note

Deleting process data is different from deleting publish data (see Publish Data, on page 57). Deleting published data removes only the data that has been published into a publish database. However, the data source in the process database remains available.



Delete Probe Array Type

Remove an unused or unwanted probe array type from the process database from the left hand pane in the **Process** tab window.

Deleting a Probe Array Type from the Process Database

- 1. Under the **Probe Array Type** folder in the left hand pane, click to expand the **Expression** folder.
- 2. Click to select the probe array type of interest.
- 3. On the menu bar select **Process** \rightarrow **Delete**, or right-click the probe array type of interest and select **Delete** from the shortcut menu that appears.
 - \Rightarrow A warning / confirmation window appears (Figure 2.5).

LIMS Manager 🛛 🔀			
⚠	Delete probe array Hu6800?		
	Yes No		
Figure 2.5			

Delete warning / confirmation window

- 4. Click Yes to delete the probe array type from the process database.
 - \Rightarrow The probe array type is removed and no longer appears listed in the left hand pane.



When a probe array type is deleted the associated usersets are also removed. Data created with these usersets remains, but is not accessible.

Export

Note

The **Export** command allows you to copy all data including sample, experiment, .dat, .cel and .chp files and send them to an external location for storage.

Exporting Data

- 1. In the left pane, click to select the sample of interest.
- **2**. On the menu bar select **Process** \rightarrow **File View**, if necessary.

You may use **Process View** if exporting the entire sample or experiment and related files.

3. In the upper right pane, click to select the data to export.

Selecting an object takes that object and all associated data with it. For example, selecting a sample takes the sample, experiment and any related .dat, .cel and .chp files.

The sample and experiment data are exported to an Affymetrix[®] Microarray Suite v4.x .exp file.

- 4. On the menu bar, select **Tools** \rightarrow **File Path**, or click the \swarrow button on the toolbar.
 - \Rightarrow The Select Export File Path window appears (Figure 2.6).

Select Export File Path	
Current Path:: h:\global_gene 700292	
Directories:	OK Cancel
Drives:	Network

Figure 2.6 Select Export File Path window



- **5.** Select a **Drive** from the drop-down list and browse to find desired export destination directory, then click **OK**.
- On the menu bar select Process → Export, or right-click the file and select Export from the shortcut menu that appears.
- 7. Click OK.

 \Rightarrow The **Export** Local Task status is displayed in the lower right pane. Double-click **Export** at any time to view additional information pertaining to the **Export** status, or right-click **Export** and select **View Task Log** from the shortcut menu that appears.

Views in Process Tab Window

The upper right pane of the **Process** tab window displays either a **File View** or a **Process View** of the selected sample or data file. The **Process View** identifies the stages completed by the experiment displayed as gray bullets, and stages that are pending are displayed as green bullets. The **File View** displays the experiment and all related files. For samples of both views, see **Figure 2.7**.

Changing Views

- **1**. Click to select a sample in the left pane.
- 2. From the menu bar select **Process** \rightarrow **Process View** (or **File View**, as appropriate).

OR

- 1. Right-click a sample in the pane on the left side of the window.
- 2. From the shortcut menu that appears, select **Process View** or **File View** as appropriate.









Print an Experiment

The **Print Experiment** command allows you to obtain a printout of experimental information. Two print options are available (described below), and samples of each are provided in *Appendix A* on page 119.

Printing an Experiment

- 1. From the upper right pane in the **Process** tab window, click to select an experiment to print.
- 2. From the menu bar click **Process**.
- 3. Select Print Experiment \rightarrow Diagram Only or Diagram & Properties.
 - Diagram Only option displays the experiment process and lists the inputs and outputs for each stage.
 - Diagram & Properties option displays the experiment processes, input and outputs and provides property information relevant to each stage.
 - \Rightarrow The experiment information is printed on the default printer.

OR

- 1. From the upper right pane in the **Process** tab window, right-click the experiment to print.
- 2. Select **Print** from the shortcut menu.
- **3.** Select Print Experiment \rightarrow Diagram Only or Diagram & Properties.
 - Diagram Only option displays the experiment process and lists the inputs and outputs for each stage.
 - Diagram & Properties option displays the experiment processes, input and outputs and provides property information relevant to each stage.
 - \Rightarrow The experiment information is printed on the default printer.

Register a LIMS Server

Registering a LIMS Server means to add a server and make it available to LIMS Manager.

Registering a Server

- 1. From the menu bar select **Tools** \rightarrow **Register LIMS Server**, or click the button on the toolbar.
 - \Rightarrow The Register LIMS Server window appears (Figure 2.8).

Register LIMS Server
Domain / Sql Server List AFFYMETRIX CISLAB SWDEVELOPMENT SWDEVELOPMENT SWTEST01 SWTEST02 SWTEST02 SWTESTSERVER
OK Cancel

Figure 2.8 Register LIMS Server dialog box

- 2. Search the tree and select the desired server to register.
- 3. Click OK.
 - \Rightarrow The selected server is added to the drop-down list on the toolbar and appears listed in the left pane of the **Process** tab (Figure 2.1).

Note

The newly registered server becomes the active server.

 \Rightarrow If the server selected is not LIMS aware, an error message is displayed (Figure 2.9).





Register LIMS Server error message

🗸 Note

An error message indicates that there is no LIMS system installed, or that there is an installation error. For further information, please refer to the 'Affymetrix[®] LIMS Installation and Administration Guide.'

Unregister a LIMS Server

Right-click to select the server to remove and select **Unregister LIMS Server** from the shortcut menu that appears.

Server Events

The **Server Events** option displays a log allowing you to view events occurring on the active LIMS server.

To view the Server Events log: Select Tools \rightarrow Server Events, or click the (button on the toolbar.

Server Tasks

Use the **Server Tasks** option to view a task window that displays the publishing queue. Review the task list at any time to monitor all publishing occurring on the active server (Figure 2.10).

To view Server Tasks: Select Tools \rightarrow Server Tasks or click the \clubsuit button on the toolbar.

Click icon for menu.	Select a column to define the primary sort order. Click the column heading once to sort the task list in an ascending manner, click again to sort descending.					
<u>H</u> estore	User 🗖	Workstation	Description V	Туре 🌂	Started 🌂	Updated 🛰 🔺
Move	AFFYMETRIX\jcampm	NTWKQJQDU	Items From GeneC	Publish	2001-04-11 16:03:46	2001-04-11 16:07:18
<u>S</u> ize	AFFYMETRIX\rbjorg	NTWKDRWKO	(1) Items From GeneC	Publish	2001-04-11 15:52:47	2001-04-11 15:56:09
Minimize	AFFYMETRIX\rbjorg	NTWKDRWKU	(1) Items From GeneC	Publish	2001-04-11 15:43:46	2001-04-11 15:47:08
Maximize	AFFYME I RIX\msavoy	NTWK1X6E6	(1) Items From GeneU	Publish	2001-04-11 15:35:12	2001-04-11 15:38:39
Cancel	AFFYME I RIX\msavoy	NTWKIX6E6	(1) Items From Genet	Publish	2001-04-06 14:20:51	2001-04-06 14:27:28
Bestart	AFFTMETRIXMIJO	NTWKUMHUL	(10) Items From Gene	Publish	2001-04-06 12:34:02	2001-04-06 14:04:51
- Hostan	AFFTMETRIX/rbjorg	NTWKDRWKU	(2) Items From GeneU	Publish	2001-04-06 11:08:50	2001-04-06 12:31:58
View <u>T</u> asks		NTWKDHWKU	(2) Items From GeneC	Publish	2001-04-06 11:08:33	2001-04-06 14:04:46
View Task <u>I</u> tems	AFFTMETHIA MISAVOY	NTWRIADED	(2) Items From GeneC	Publish	2001-04-05 17:35:52	2001-04-05 17:44:46
View <u>A</u> ll Items	AFFTMETHIA INSAVOY	NTWRIADED	(2) Items From Genet	Publish	2001-04-05 14:45:36	2001-04-05 14:52:39
Save As	AFFTMETHIA INSAVOY	NTWRIADED	(2) Items From Genet	Publish	2001-04-04 13:21:50	2001-04-04 13:23:21
Delete Task Items		NTWRIADED	(2) Items From GeneC	Publish	2001-04-04 11:49:12	2001-04-04 11:49:30
	AFFTMETDIXGamor	NTWKIAGED	(2) Items From GeneC	F ublish Dublish	2001-04-04 03:38:00	2001-04-04 11:13:37
<u>C</u> lose Alt+F4	AFEYMETRIX/campin	NTWK0JQDU	(3) Items From GeneC	Publish	2001-03-10 10.44.12	2001-03-13 10.46.10
		NTWK0JQDU	(3) Items From GeneC	Publish	2001-03-13 11.47.00	2001-03-13 11.47.41
	AFFYMETRIX\icampin	NTWK00000	(3) Items From GeneC	Publish	2001-03-13 11:41:22	2001-03-13 11:32:43
	Art The This (campin	NT WROUGDO	joj tems From Genee	T GDIGTT	2001 00 10 11.02.01	
Publish Service Active						
Identifies a new item in the queue, pending processing.						
	a plata d taol				serv	ice status
✓ Indicates a completed task.						
Indicates that an error occurred and the task was not completed.						

Figure 2.10

Server Tasks window



Task List Commands

Further action may be taken on items in the Task List window. Access the command menu for the task list either by clicking the 📑 icon in the upper left corner of the window (Figure 2.10), or by right-clicking a task (Figure 2.11).

Viewing Tasks

The Task List window displays the status of each task on the server when **View Tasks** is selected (Figure 2.11).

🔣 SWTES	iTO1 Task List							×
Task ID	Status	User	Workstation	Description	Туре	Started	Updated	
75	COMPLETE	AFFYMETRIX\	NTWK68GPN	(9) Items From	Publish	2001-04-11 16:03:46	2001-04-11 16:07:18	
🔀 74	ERROR	AFFYMETRIX\	NTWK68GPN	(43) Items From	Publish	2001-04-11 15:52:47	2001-04-11 15:56:09	
✓ 73	COMPLETE	AFFYMETRIX\	Cancel	ems From	Publish	2001-04-11 15:43:46	2001-04-11 15:47:08	
🔀 72	ERROR	AFFYMETRIX\	Bestart	ms From	Publish	2001-04-11 15:35:12	2001-04-11 15:38:39	
71	COMPLETE	AFFYMETRIX\	- ngotan	ms From	Publish	2001-04-06 14:20:51	2001-04-06 14:27:28	
70	COMPLETE	AFFYMETRIX\	View <u>T</u> asks	ms From	Publish	2001-04-06 12:34:02	2001-04-06 14:04:51	
🔀 69	ERROR	AFFYMETRIX\	View Task <u>I</u> tems	ns From	Publish	2001-04-06 11:08:50	2001-04-06 12:31:58	
🔀 68	ERROR	AFFYMETRIX\	View <u>All Items</u>	ms From	Publish	2001-04-06 11:08:33	2001-04-06 14:04:46	
🔀 67	ERROR	AFFYMETRIX\		ms From	Publish	2001-04-05 17:35:52	2001-04-05 17:44:46	
🔀 66	ERROR	AFFYMETRIX\	Sa <u>v</u> e As	ms From	Publish	2001-04-05 14:45:36	2001-04-05 14:52:39	
🔀 65	ERROR	AFFYMETRIX\	<u>D</u> elete Task Iter	ns ms From	Publish	2001-04-04 13:21:50	2001-04-04 13:23:21	
✓ 64	COMPLETE	AFFYMETRIX\	NTWKQIQDU	(2) Items From	Publish	2001-04-04 11:49:12	2001-04-04 11:49:30	
✓ 63	COMPLETE	AFFYMETRIX\	NTWK68GPN	(3) Items From	Publish	2001-04-04 09:38:00	2001-04-04 11:13:37	
✓ 62	COMPLETE	AFFYMETRIX\	NTWK68GPN	(3) Items From	Publish	2001-03-15 10:44:12	2001-03-15 10:46:10	
✓ 61	COMPLETE	AFFYMETRIX\	NTWKQJQDU	(6) Items From	Publish	2001-03-13 11:47:05	2001-03-13 11:47:41	
✓ 60	COMPLETE	AFFYMETRIX\	NTWKQJQDU	(6) Items From	Publish	2001-03-13 11:41:22	2001-03-13 11:41:59	
✓ 59	COMPLETE	AFFYMETRIX\	NTWKQUQDU	(6) Items From	Publish	2001-03-13 11:32:07	2001-03-13 11:32:43	⊒
1							•	
						Publish	Service Active	

Figure 2.11

Right-click task for shortcut menu

Viewing Task Items

The Task List window displays all items for a specified task when **View Task Items** is selected (Figure 2.12).

🔣 Items For T	ask #73						
Task ID	Publish Type	Item Name	Database	Overwrite Type	Intensity Flag	Item Status	Updal
73	Experiment	venu1071	LIMS_MGR	1	1	COMPLETE	10-11
73	CEL File	venu1073.CEL	LIMS_MGR	1	1	COMPLETE	10-11
73	CHP File	venu1073.CHP	LIMS_MGR	1	1	COMPLETE	10-11
73	Experiment	Hu6800subD	LIMS_MGR	1	1	COMPLETE	10-11
73	CEL File	Hu6800subD	LIMS_MGR	1	1	COMPLETE	10-11
73	CHP File	Hu6800subD	LIMS_MGR	1	1	COMPLETE	10-11
73	Experiment	venutest10111	LIMS_MGR	1	1	COMPLETE	10-11
73	CEL File	venutest10111	LIMS_MGR	1	1	COMPLETE	10-11
73	CHP File	venutest10111	LIMS_MGR	1	1	COMPLETE	10-11
73	Experiment	exp1	LIMS_MGR	1	1	COMPLETE	10-11
73	CEL File	probimg.CEL	LIMS_MGR	1	1	COMPLETE	10-11
73	CHP File	probimg.CHP	LIMS MGR	1	1	COMPLETE	10-11
						Publish Servic	e Active

Figure 2.12

View Task Items window

Viewing All Items

All items for all tasks are displayed in the Task List window when **View All Items** is selected.

Saving the Task List

The **Save As** command allows you to save displayed items from the Task List window as a .txt file.

- Click the sicon in the upper left corner of the Task List window and select Save As from the drop-down menu. Alternatively, right-click a task in the list and select Save As from the shortcut menu that appears.
 - \Rightarrow The Save As dialog box appears (Figure 2.13).







- 2. From the Save in drop-down list, browse to the desired destination.
- **3.** Name the file and click **Save**.
 - \Rightarrow The task list is saved as a text file in the defined destination.

Deleting Task Items

The Delete Task Items command removes all listed task items.

- **1**. Click to select a task.
- 2. Right-click the task and select **Delete Task Items** from the shortcut menu that appears.

 \Rightarrow A warning / confirmation window appears (Figure 2.14).



Task Delete warning / confirmation window

- 3. Click Yes to delete all tasks on and before the date and time indicated.
 - \Rightarrow Items are deleted and Task List window is empty.

Cancelling an Active Publish Task

Right-click a pending task and select **Cancel** from the shortcut menu to stop the publish process.

Restarting a Cancelled Publish Task

Right-click a task that has been cancelled or stopped due to error and select **Restart** to reactivate the publish process.

Find Function

Use the **Find** feature to query **Samples**, **Experiments** or analyses within the process database for **Export**, **Archive**, **Assume Ownership** or **Delete** functions. When preforming one of these functions after a **Find** command, only the selected objects are affected by the action and do not include any related children objects.

V Note The **Find** function only searches the process database.

Finding a Sample, Experiment or File

- 1. From the menu bar select **Tools** \rightarrow **Find**, or click the **M** button on the Toolbar.
 - \Rightarrow The Find window appears (Figure 2.15).



📴 LIMS Manage	<mark>r - [Find1</mark> Tools Find	l] d Window Help					_ 문 ×
DBSERVER04		- <u>Tungen</u> , <u>T</u> obb	🍫 🖻 🗛 🝸	?			
Select Object Ad	Ivanced Filte	r Date Filter		I			
Sample names:		Sample type:					Find Now
[All Samples] Experiment name		[All Types] Probe array type:					Stop
[All Experiments]		[All Types]	Select a	column h	eading to define	the	<u>C</u> lear
Sample project:	/ ا ا	Assays:	primary list in an	sort order ascendin	. Click once to so g manner, click a	rt the gain to	Q
User:		File type:	sort des	cending.	•	-	
[All Users]	-	[All Types]					
Comments		Description		$/ \setminus $	$\langle \rangle$		
<u>'</u>				- <u></u>			
Sample Name (17)	Assay	Sample Project	Sample Lype	User	Uate	Description	Lomments A
	Expression	NEW2 Meuse	(null) (mull)	student03	Oct 20 1333 01:41FM		N
Terrest	Expression	mutest	(null) mutest	alew	Mar 24 2000 03:47PM		2
	Expression	Wisc	total BNA	lwang	Mar 30 2000 10:06AM		
🗒 TBHQ	Expression	Wisc	total RNA	lwang	Mar 30 2000 10:11AM		
🗒 Uof W	Expression	Uof W	total RNA	lwang	Mar 30 2000 10:21AM		
🗒 TBHQ1	Expression	Uof W	total RNA	lwang	Mar 30 2000 10:25AM		
🗐 polyA_spikes	Expression	polyA_spikes	(null)	lwang	Apr 04 2000 05:58PM		
🧱 ctest	Expression	mytest	mytest	ctruon	Feb 26 2001 06:04PM		
al_temp	Expression	al_temp	al_temp	alew	May 08 2001 04:10PM		_
 ∙							•
For Help, press F1							NUM

Figure 2.15

Find window with Advanced Filter tab displayed



All three tabs work in union. For Example, if <u>sample</u> is selected as the **Object Type** in the **Select Object** tab, when you switch to the **Advanced Filter** tab, only those boxes that pertain to finding a <u>sample</u> will be active.

Multiple Find windows are allowed.

2. Click either the Select Object, Advanced Filter or Data Filter tab and make selections from the drop-down lists, or type in requested information.

Further refine your search by selecting additional tabs.

3. Click Find Now.

- \Rightarrow Matching results are displayed at the bottom of the window.
- 4. Click to select the sample, experiment or file of interest.

The commands available for the selected object are found under **Find** on the menu bar, or right-click the object to access these commands from the shortcut menu (Figure 2.17).

To select multiple contiguous objects to act upon: click an object in the list, then SHIFT-click another item anywhere either above or below the first selected. All objects between the two selections are selected.

To select multiple non-contiguous objects to act upon: click an object in the list, then CTRL-click another item anywhere in the list. Repeat as necessary for additional items.

Follow the procedure(s) for the **Export**, **Archive/Unarchive**, **Delete** and **Assume Ownership** commands outlined earlier in this chapter. There are two differences when performing these procedures after a **Find** task; **①** the commands are found under **Find** on the menu bar, and **②** the **Export**, **Assume Ownership** and **Archive/Unarchive** procedures only affect the specific file selected and do not include any associated data.

Sample Filter

The **Filter** function is used to reduce the number of samples displayed in a tree of samples. The **Filter** icon becomes active in the **Process** tab and the **Import** tab when you click a node under a **Samples** folder and remains active from session to session until the parameters are changed.

Using the Sample Filter

- 1. Click the Samples folder.
 - \Rightarrow The tree expands displaying samples and the **Filter** icon on the toolbar becomes active.
- **2.** Click the **Filter** icon $\overline{\mathbf{Y}}$.
 - \Rightarrow The Sample Filter dialog box appears (Figure 2.16).



Sample Filters	×
Sample name:	
[All Samples]	_
Sample project:	
[All Projects]	▼
Sample type:	
[All Types]	•
Role:	User:
[All Roles]	[All Users]
Filter by sample crea	ate date range
From : Tuesday ,	May 15, 2001 💌
To: Tuesday ,	May 15, 2001 💌
	Cancel

Figure 2.16 Sample Filter dialog box

3. From the drop-down boxes, define filter parameters.

The **Role** and **User** filter options are mutually exclusive. You may select one or the other, but not both.

- 4. Click Apply.
 - \Rightarrow The Samples are filtered and the sample list is adjusted accordingly.

✔ Note

The filter parameters remain active from session to session. You must reset the parameters when a different display is desired.



Figure 2.17 Right-click objects for shortcut menus after a Find procedure





Chapter 3





The **Publish** tab allows you to create new publish databases and delete published data from a database. All publish databases are password-protected and can only be accessed using the password defined during the creation of the database.

This chapter orients you to the layout of the **Publish** tab and describes in detail the functions and tasks accomplished here.

Publish Tab Window

There are three panes in the **Publish** tab window (Figure 3.1). The first pane on the left side displays a tree showing the active LIMS server, all publish databases residing on the server, all supported probe array types and their contents.

The second pane becomes active and is filled with a create database dialog when a **Create Database** command is initiated.

The third pane in the lower right corner is the Local Task status view that displays the status of processes that may be occurring in the database, such as **Export**, **Import**, **Delete**, **Archive** and **Create**. This view allows you to monitor these jobs while concurrently working on additional tasks within the LIMS system.



Figure 3.1

The three panes of the Publish tab window

Publish Databases

A new security feature in LIMS Manager v3.0 is password protection for publish databases. Passwords are defined during the creation of the database and can be changed on the server using the LIMS Sync utility. For more information see *Affymetrix LIMS Installation and Administration Guide*.

Accessing Publish Databases

- 1. From the **Publish** tab, click the database you want to access.
 - \Rightarrow A Publish Database Login window appears.
- 2. Enter the database **Password** and click **Login**.

🗸 Note

Passwords are case-sensitive.

 \Rightarrow The publish database tree expands, displaying the **Probe Array Type** folder (Figure 3.2).





Deleting

The **Delete** function allows you to delete data or a probe array type from a publish database.

Probe Array Type

Probe array types that are no longer used may be removed from the publish database. When a probe array type is deleted all published data associated with it, from the parent down, are also deleted.

Deleting a Probe Array Type

- 1. Click the **Probe Array Type** folder under the publish database of interest.
- 2. Click to select the probe array type to delete.
- **3**. From the menu bar select **Publish** \rightarrow **Delete**
 - \Rightarrow A warning window appears (Figure 3.3).

LIMS Manager 🛛 🔀		
⚠	Delete probe array HG_U95A?	
	Yes <u>N</u> o	



- 4. Click Yes to delete the selected probe array type.
 - \Rightarrow The probe array type is removed from the publish database and all associated data is deleted.

OR

- **1**. Right-click the probe array type to delete.
- 2. Select **Delete** from the shortcut menu that appears.
 - \Rightarrow A warning / confirmation window appears (Figure 3.3).

- **3**. Click **Yes** to delete the selected probe array type.
 - \Rightarrow The probe array type is removed from the publish database and all associated data is deleted.

Publish Data

Deleting publish data removes the specified experiments or analyses from the publish database. When deleting a published data file, all associated files below are also removed.

I CAUTION

Deleting publish data permanently removes data from the database. Deleted data can be restored by repeating the publish procedure in Affymetrix [®] Microarray Suite only if the information still resides on a process database.

Deleting Published Data

- **1.** From the drop-down list on the toolbar, select the server where the publish database resides.
- **2.** In the first pane on the left side of the window, select the database of interest and expand the tree.
- **3.** Click to select the data to delete.

You may either select a single analysis file or an experiment.

- 4. On the menu bar select **Publish** \rightarrow **Delete**.
 - \Rightarrow A warning / confirmation window appears (Figure 3.4).



Figure 3.4 Analysis delete warning / confirmation window



5. Click **Yes** to continue.

⇒ The Delete task proceeds. An information window appears asking you to please wait while published file(s) are deleted. In the lower right pane the status is displayed.

To view more information pertaining to the delete task, double-click **Delete** in the lower right pane, or right-click **Delete** in the lower right pane and select **View Task Log** from the shortcut menu (Figure 3.6).

OR

- 1. From the drop-down list on the toolbar, select the server where the publish database resides.
- **2.** In first pane on the left side of the window, select the database of interest and expand the tree.
- **3.** Right-click the data to delete and select **Delete** from the shortcut menu (Figure 3.5).





- \Rightarrow A warning window appears (Figure 3.4).
- 4. Click Yes to continue.
 - \Rightarrow The **Delete** task proceeds.

An information window appears asking you to please wait while published file(s) are deleted. In the lower right pane the status is displayed.

To view more information pertaining to the delete task, double-click **Delete** in the lower right pane, or right-click **Delete** in the lower right pane and select **View Task Log** from the shortcut menu (Figure 3.6).



Active Delete publish data task



Creating Publish Databases

Create new publish databases at any time from the **Publish** tab of Affymetrix[®] LIMS Manager. In addition to creating the new database, the system also creates the necessary ODBC DSNs for publishing data. Several key pieces of information are required to create a publish database. This information includes:

- Server on which to create the database
- Username and password with appropriate permissions for creating databases on the selected server
- Size of database to create
- Proper drive and path for the database files

The server name selected on the **Create Publish Database** pages is represented by an ODBC data source name description. This ODBC DSN specifies the location of an administration database that is used to create the actual publish database. This DSN is normally created during the installation process of the Affymetrix[®] LIMS system.

Database Size Considerations

When creating a Microsoft[®] SQL Server publish database, the initial size of the database can be between 128 MB and 1024 MB. This size corresponds to approximately the following storage capabilities (based on a probe array with 270 x 270 cells, 1800 probe sets per probe array):

- 8 MB per experiment (includes experiment information, cell file and chip file)
- 8 MB per library file
- 128 MB database = 4 library files and 12 experiments
- 1024 MB database = 16 library files and 110 experiments

Creating an Oracle[®] publish database has only 2 size options, **Small** and **Large**. The small database takes approximately 6 GB of space and the large is almost 12 GB. These publish databases have enough room to publish approximately 500 and 1000 experiments respectively.

The size of a Microsoft[®] SQL Server database can easily be increased once the database is created. For information on increasing the size of a SQL Server publish database, see *Affymetrix*[®] *LIMS Installation and Administration Guide*.

Microsoft[®] SQL Server Publish Database

Affymetrix[®] LIMS Manager 3.0 features increased security with a password-protected publish database. Publish databases can only be accessed with the password defined at the time the publish database was created.

Creating a Microsoft® SQL Server Database

- 1. From the drop-down list on the toolbar, select the server where the publish database will reside.
- 2. Click to select the server icon in the left pane of the **Publish** tab.
- 3. On the menu bar select **Publish** → **Create SQL Database**, or rightclick the server icon in the left pane and select **Create SQL Database**.
 - \Rightarrow The Create SQL Database dialog box appears in the upper right pane (Figure 3.7).

Atabase name: //yPublish /atabase password: infirm database password: i	
MyPublish Intracesses ionfirm database password: ionfirm database password: intracesses ata device ath: Size (MI D:\GeneChip\Data float intracesses intracesses Size (MI D:\GeneChip\Data float intracesses Size (MI D:\GeneChip\Data intracesses intracesses intracesses intracesses Size (MI D:\GeneChip\Data intracesses intracesses </th <th></th>	
atabase password: infiim database password: in	
ata device	
ionfim database password: ata device ata device ath: Size (MI D:\GeneChip\Data 1024 dex device ath: Size (MI D:\GeneChip\Data 256 bg device	
ata device Tath: Size (MI D:\GeneChip\Data 1024 dex device ath: Size (MI D:\GeneChip\Index 256 bg device	
ata device	
ata device	
tath: Size (M) D:\GeneChip\Data 1024 idex device	
D:\GeneChip\Data 1024 idex device):
idex deviceSize (Mi zahr:Size (Mi):\GeneChip\Index256 ug device	
ah: Size (Mi D:\GeneChip\Index 256 g device	
256 256 256 256 256 256 256 256 256 256	I):
og device	
ath: Size (Mi):
D:\GeneChip\Log 256	
Install Default Clear Close	

Figure 3.7

Create SQL Database dialog box



- 4. Enter the following information in the **Database** section:
 - Database name: <name for the new publish database>
 - Database password: <password of your choice>
 - Confirm database password: <re-enter password to confirm>

The publish database name is limited to 22 characters in length.

- 5. Enter the following information in the **Data device**, **Index device** and **Log device** sections:
 - Path: Use the complete drive description, including drive letter, colon and backslash (E:\Affymetrix DBDevices).

```
✓ Note
```

Note

The drive letter specifies the drive on the actual publish database server. If the publish database is being created on a remote server, make sure the selected drive letter is available on that server.

- Size (MB): The size depends on the number of experiments that will be published to the database and the amount of available space on the drive selected. Microsoft[®] SQL Server databases are relatively easy to expand after creation; therefore, creating a smaller database and expanding it as necessary is recommended.
- 6. Click Install.
 - \Rightarrow It takes several minutes to create the database and allocate the amount of drive space requested (2 -15 minutes or more, depending on the speed of the server and drive).

Once created, the database is available for publishing without further configuration. If the newly created database does not appear in the **Select Publish Database**(s) drop-down list, verify that the database and the ODBC DSN were created properly.

Oracle[®] Publish Database

✔ Note

The Oracle® publish databases are much larger than the default Microsoft® SQL Server databases. Verify that there is adequate space on the server before creating an Oracle database. The **Small** Oracle publish database requires 6 GB and the **Large** database requires 12 GB. These sizes correspond to publishing approximately 500 and 1000 experiments respectively.

Creating an Oracle® Publish Database

- 1. From the drop-down list on the toolbar, select the server where the publish database will reside.
- 2. Click to select the server icon in the left pane of the **Publish** tab.
- On the menu bar select Publish → Create Oracle Database, or rightclick the server icon and select Create Oracle Database from the shortcut menu.
 - \Rightarrow The Create Oracle Database dialog box appears in the upper right pane (Figure 3.8).

Publish Oracle alias name: SIG03-	
Database	Data device
Database name:	Server device path:
Database password:	Database size:
	Small (500)
Confirm database password:	
<u>I</u> nstall <u>D</u> efault	<u>Clear</u> Clo <u>s</u> e





- 4. In the **Database** section enter the following information:
 - Database name: <name for the new publish database>
 - Database password: password of your choice>
 - Confirm database password: <re-enter password to confirm>

The publish database name is limited to 22 characters in length.

- 5. In the **Data Device** section enter the following information:
 - Server Device Path: This is the complete path for the database files. Use the complete drive description for Windows 2000. For example: E:\Affymetrix DBdevices\ Be sure to include the trailing slash.

🗸 Note

Note

The datafile path directory used specifies the file location on the actual publish database server. If the publish database is being created on a remote server, make sure the path specified is available on that server.

- Database Size: from the drop-down list. The Small database requires approximately 6 GB of drive space and the Large requires approximately 12 GB.
- 6. Click Install.
 - \Rightarrow It takes several minutes to create the database and allocate the amount of drive space requested (15 30 minutes or more, depending on the speed of the server and drive).

Once created, the database is available for publishing without further configuration. If the newly created database does not appear in the **Select Publish Database**(s) drop-down list, verify that the database and the ODBC DSN were created properly.



Chapter 4




Data created in Affymetrix[®] Microarray Suite and stored on a local hard drive is moved, or copied, to a process database on a LIMS server with LIMS Manager via the **Import** tab. The additional tasks of editing and updating data are also accomplished from the **Import** tab.

This chapter orients you to the layout of the **Import** tab and describes in detail the tasks accomplished here.

Import Tab Window

The window of the **Import** tab is divided into four panes (Figure 4.1). The first pane displays all registered servers and process databases and lists the samples associated with expression assays.

The second pane in the upper right corner is tabbed with a separate file type on each of the three tabs. The tabs display the corresponding files found at the defined import file path.

In the upper center of the **Import** tab window is the **Data to Import** pane, a staging area where you identify the sample and associate the files of interest with the sample.

The fourth pane found in the lower right corner displays either a Property Information or Local Task status view. Property Information lists environment details pertaining to the selected item. This information is primarily obtained from the **Sample Registration** process and includes information such as **Sample Project**, **Sample Type**, **Initial Stage**, **Description**, **Comments**, **Date**, **User** and **Template**.

The Local Task status view displays the status of processes that may be occurring in the database, such as **Export**, **Import**, **Delete**, **Archive** and **Create**. This view allows you to monitor these jobs while concurrently working on additional tasks within the LIMS system.



Figure 4.1

The four panes of the Import tab window

Entering and Updating Data

Only the **Import** tab requires your input to enter or update data. The following rules apply for all user input data fields:

Table 4.1Rules for Data Input

Data Input	Rule
Character length	≤ 64 characters
Invalid Characters	\/:*?"<> '~´,{}[]
Required Data fields	Required fields are indicated by blue titles and may not be left blank
Data fields	Names may not duplicate existing names

Table 4.2

Valid Field Lengths

Editing Samples Attributes	Length
Sample Project	64
Sample Type	64
Description	64
Comments	64
Sample Name	64



Importing Data

The **Import** function imports or copies data files created in Affymetrix[®] Microarray Suite stored on a local hard drive and moves them to a process database on a LIMS server. It is possible to import data to multiple LIMS Servers. Importing data is a three-step process consisting of selecting a file path, identifying a sample and associating files with the sample.

Selecting an Import File Path

- On the menu bar select Tools → File Path, or click the button on the toolbar.
- **2.** Select a **Drive** from the drop-down list and browse to find the desired files.
- 3. Click OK.
 - \Rightarrow The file path defined appears above the upper right pane, and the data files at that location are listed under the **EXP**, **CEL** and **CHP** tabs (Figure 4.2).

File Path: h:\global_gene 700292\huheart and lung file	s-
EXP CEL CHP	
heart1.EXP	
heart2vheart1.EXP	
Lung1.EXP	
Lung2vheart1.EXP	
Lung2vLung1.EXP	
S0498a4x.EXP	
S0498aA4x.EXP	
S0498bA4x.EXP	
S0598aA4.EXP	
S0598bA4.EXP	

Figure 4.2 File Path location and data type tabs



The default file path is the same path assigned to Affymetrix[®] Microarray Suite experimental data.

Identifying a Sample

LIMS offers the ability of tracking both the sample and projects associated with an array (experiment). Affymetrix[®] Microarray Suite starts with the array but does not make project or sample associations. Thus, after selecting a file path, the next step requires you to rebuild this association. This is accomplished by providing a project and sample for each imported array. You do this by creating a new sample or importing data into an existing sample.

Creating a New Sample

The **Create Sample** option, found on the **Import** tab, captures sample information and is available for all assay types. **Create Sample** differs from the **Register Sample** option, available in Affymetrix[®] Microarray Suite that captures the sample information defined by the template used.

- 1. In the left pane, click the **Expression** folder.
- 2. On the menu bar select **Import** \rightarrow **Create Sample**, or right-click the folder and select **Create Sample** from the shortcut menu.
 - \Rightarrow The Create Sample dialog box appears (Figure 4.3).

	required fields
eate Expression Sample	
Sample information Sample name:	
Sample project:	
Sample type:	
Description:	×
Comments:	
OK	Cancel

Figure 4.3 Create Sample dialog box



- 3. Enter the Sample name
- 4. Enter the **Sample project** and **Sample type**, or select from the dropdown lists.
- 5. Enter **Descriptions** and **Comments**, if desired.
- 6. Click OK.
 - \Rightarrow The sample created appears in the center, **Data to Import** pane.

VNote Once a sample is created, all sample information may be edited except the **Sample Name**. You may only edit new samples and not existing ones. See "Edit Sample" on page 75.

Selecting an Existing Sample

The second method of identifying a sample is to select an existing sample from the process database tree.

- 1. In the first pane, expand the **Expression** folder and locate the sample of interest.
- Click and drag the sample from the first pane and drop into the Data to Import pane.
 - ⇒ The sample and all associated files appear in the staging area with red check marks indicating that these files already exist in the database (Figure 4.4).



Associating Files

The third and final step of the import data task is to associate files with a sample. Drag and drop an .exp file, and the .cel and .chp files with the same prefix are automatically associated.

Additional files with different prefixes may also be added. This is done in a hierarchical manner. First match the .exp file(s) to the sample. Then match the .cel file(s) to the .exp, and finally the .chp to the .cel.

V Note The filenames of imported files need to conform to the naming conventions within LIMS (see page 69). You may need to rename files to ensure that the filenames conform.

Associating .exp with Sample

The **EXP** tab displays all files located at the designated path that contain the .exp extension. The import process also migrates the .dat file associated with the .exp file via a common filename. Importing an .exp file that does not have an associated .dat file within the designated folder results in an error message. These .dat files without associated .exp files are not available for import.

- 1. Click and drag the .exp file from the upper right pane and drop on top of the sample you wish to associate it with in the **Data to Import** pane.
 - \Rightarrow The .exp file appears on the branch below the sample name with .cel and .chp files containing the same prefix (Figure 4.5).



2. Repeat step 1 to associate additional .exp files with the sample.



Note

3. If other .cel or .chp files with different prefixes are to be included, continue with the following steps.

You can only associate files from probe array types installed in the LIMS database. If a particular probe array type is not installed, an error message appears when you try to associate files from that probe array type.

Associating .cel with .exp

- **1**. Select the **CEL** tab.
- 2. Click and drag the .cel file from the upper right pane and drop on top of the .exp file you wish to associate it with in the **Data to Import** pane.
 - \Rightarrow The .cel file appears in the branch below the .exp file in the **Data to Import** pane (Figure 4.6).



Figure 4.6

Associating a .cel file with an .exp

3. Repeat step 2 for additional .cel files.



To avoid having to associate a .cel and .chp file with each .exp file imported into the process database, simply import only the .exp file. Then recreate the .cel and .chp files using Affymetrix® Microarray Suite in LIMS mode. This minimizes potential error from mistakenly dragging a .chp file onto the wrong .cel file. Contact Affymetrix Technical Support for help on batch analysis methods in Affymetrix® Microarray Suite, or refer to Affymetrix® Microarray Suite User's Guide.

Associating .chp with .cel

- **1**. Select the **CHP** tab.
- 2. Click and drag the .chp file from the upper right pane and drop on top of the .cel file you wish to associate it with in the **Data to Import** pane.
 - \Rightarrow The .chp file appears in the branch below the .cel file in the **Data to Import** pane.
- **3.** Repeat step 2 for additional files to associate.

Start Import

From the menu bar select **Import** \rightarrow **Start Import**, or right-click the sample and select **Start Import** from the shortcut menu.

 \Rightarrow The import task proceeds and the status is displayed in the lower right pane.

Double-click **Import** to view more information pertaining to the import task, or right-click **Import** in the lower right pane and select **View Task Log** from the shortcut menu.

Edit Sample

A sample can only be edited in LIMS Manager before it has been imported into a process database. After a sample has been imported, or created via Microarray Suite, it can no longer be edited.

- 1. Click to select a sample to edit in either the **Data to Import** pane that has not been imported.
- 2. On the menu bar select **Import** \rightarrow **Edit Sample**, or right-click the sample and select **Edit Sample** from the shortcut menu.
- 3. Make changes as necessary. You may not alter the Sample Name.
- 4. Click OK.



Remove and Remove All

To correct errors that occur as you build the sample to import, use the **Remove** or **Remove** All command.

- In the Data to Import pane, click to select the file to remove. Select a node without children and the Remove command is active. Select a parent node and the Remove All command is active and erases the parent file and all associated children files from the Data to Import pane.
- On the menu bar select Import → Remove or Remove All as appropriate, or right-click the file to remove and select from the shortcut menu.
 - \Rightarrow The file(s) disappear from the **Data to Import** pane.

Sample Filter

The **Filter** function is used to reduce the number of samples displayed in a tree of samples. The **Filter** icon becomes active in the **Process** tab and the **Import** tab when you click a node under a **Samples** folder and remains active from session to session until the parameters are changed.

Using the Sample Filter

- 1. Click the **Samples** folder.
 - \Rightarrow The tree expands displaying samples and the **Filter** icon on the toolbar becomes active.
- **2.** Click the **Filter** icon $\overline{\mathbf{Y}}$.
 - \Rightarrow The Sample Filter dialog box appears (Figure 4.7).

ample Filters			×
Sample name:			
[All Samples]			•
Sample project:			
[All Projects]			•
Sample type:			
[All Types]			•
Role:		User:	
[All Roles]	•	[All Users]	•
Filte	r by sample crea	ate date range	
From :	Tuesday ,	May 15, 2001	7
To:	Tuesday ,	May 15, 2001	Y
L.	Apply	Cancel	
aure 4.7			

- Sample Filter dialog box
- **3.** From the drop-down boxes, define filter parameters.

The **Role** and **User** filter options are mutually exclusive. You may select one or the other, but not both.

- 4. Click Apply.
 - \Rightarrow The Samples are filtered and the sample list is adjusted accordingly.

🗸 Note

The filter parameters remain active from session to session. You must reset the parameters when a different display is desired.





Chapter 5





The userset is a predefined set of parameters used when running an expression analysis in Affymetrix[®] Microarray Suite. Creating and modifying a userset is accomplished from this tab in Affymetrix[®] LIMS Manager.

This chapter orients you to the layout of the **Userset** tab window and describes in detail the functions and tasks accomplished here.

Userset Tab Window

The window of the **Userset** tab is divided into three panes (Figure 5.1). The first pane displays the active servers and process databases, and lists defined usersets and probe array types associated with the expression assay.

The second pane in the upper right corner is the userset definition area where existing usersets are reviewed or edited and where new usersets are created.

The third pane found in the lower right corner displays the Local Task status view that displays the status of processes that may be occurring in the database, such as **Export**, **Import**, **Delete**, **Archive** and **Create**. This view allows you to monitor these jobs while concurrently working on additional tasks within the LIMS system.

You only have access to this tab if the appropriate privileges are defined for you in the **Roles** tab.



Figure 5.1

The three panes of the Userset tab window

Userset Definition

The algorithm involved in expression probe array analysis has approximately 18 user-defined parameters that are set before an analysis is run. Rather than requiring you to adjust these parameters for each analysis, a predefined userset, created by an administrator or other authorized user, frees you to simply select a userset and run the analysis. The function of the userset is to create defined parameters to be used when running an expression analysis.



Creating a Userset

The userset is created in Affymetrix[®] LIMS Manager and used by Microarray Suite when analyzing a probe array. The procedure outlined below creates the basic shell of a new userset, giving it a name and defining probe array types.

Creating a Userset

- 1. Type a unique userset name in the **Create set name** box.
- From the Probe Array Types list box, select a probe array type and click >> to move to the Probe array types used list box (Figure 5.3). Repeat this step to add as many probe array types as desired.

Current server: SWTEST01		
Probe Array Types : HG_U95B HG_U95C HG_U95D HG_U95E Hu6800subA Hu6800subB Hu6800subB Hu6800subC	Probe array types used: Hu6800 Hu6800subA Hu6800subB	Existing userset name(s): JC1 MeetaUserset Userset_0720_001 VetUserSet0718_1 VetUserSet0718_2 VetUserSet0718_2 Create set name: Name Set
Scaling Normalization Probe Mask • User Defined Scale F • All Probe Sets 1 • Selected Probe Sets 1 • Belected Probe Sets Hu6800subB_ClassA.MSK Hu6800subB_ClassA.MSK Hu6800subB_ClassA.MSK Hu6800subB_ClassA.MSK Hu6800subB_ClassA.MSK	Baseline Parame actor Alpha1 Tau = Gamm. Gamm. Gamm. Perturb	ter = 0.040000 2 = 0.060000 0.015000 alt = 0.002500 alt = 0.002500 a2H = 0.003000 a2H = 0.003000 a2L = 0.003000 pation = 1.220000
Create Set	Clear D	efaults

Figure 5.3 Creating a new userset

- 3. Click Create Set.
 - \Rightarrow The name of the new userset appears in the **Existing userset** name(s) list box in the upper right corner and in the left hand pane.
- 4. Proceed to *Modifying a Userset*.

Modifying a Userset

The 18 user-defined parameter settings are used by the algorithm when running an expression analysis. An existing userset is modified by defining these parameter values.

A userset may also be modified by removing a probe array type. See page 88 for more information.

Define Parameter Values

An independent parameter is a static setting that is not influenced by probe array selection. A dependent parameter is a setting that is specific to the particular probe array type selected. This section outlines how to modify these settings.



For a detailed explanation of parameters, refer to the Affymetrix[®] Microarray Suite User's Guide.



Defining Independent Parameters

- **1.** From the independent parameters list box (Figure 5.4) click to select a parameter of interest.
 - \Rightarrow The selected parameter name and current value are displayed in the box above.

Scaling Normalization Probe Ma	isk Baseline	Gamma1H	
User Defined Sc. All Probe Sets Selected Probe Sets Hu6800subB_ClassA_MSK Hu6800subB_ClassAB_MSK Hu6800subB_ClassAB_MSK Hu6800subB_ClassC_MSK HU6800suB_CLASS HU680SUB HU680SUBB_ClasSC_MSK HU680SUBB_CLASS HU680SUBB_CLASS HU680	ale Factor	0.002500 Alpha1 = 0.040000 Alpha2 = 0.050000 Tau = 0.015000 Gamma1H = 0.002500 Gamma2H = 0.002500 Gamma2H = 0.003000 Gamma2L = 0.003000 Petturbation = 1.220000	Modify
Create S	et Clear	Defaults	



Selecting Independent Parameters

- 2. Type a value in the box above to edit.
- 3. Click Modify.

New values do not take effect until Update Set or Create Set is clicked.

4. Repeat for additional values.

Click **Defaults** to reset independent and dependent parameter values.

5. Click Update Set.

Defining Dependent Parameter Values

- 1. From the **Probe array types used** list box, click to select a probe array type (Figure 5.5).
 - \Rightarrow The selected probe array information appears in the dependent parameter box.
- 2. Select a method for how the data will be analyzed by clicking the appropriate tab; Scaling, Normalization, Probe Mask or Baseline.

Current server: SWTEST01 Probe Array Types : HG_U95B HG_U95D HG_U95D Hu6800subA Hu6800subA Hu6800subD Hu6800subD	Probe array types used: Hu68003ubA Hu6800subA Hu6800subC Hu6800subC Hu6800subD	Existing userset name(s): MeetaUserset Userset_0720_001 VctUserSet0718_1 VctUserSet0718_2 VctUserSet12345678901234 Create set name: New_Set
Scaling Normalization Probe Mask Image: Constraint of the set of the s	Baseline Para actor 0.00 Alpi Tau Gar Gar Gar Perf	Modify 100000 Modify 1a1 = 0.040000 100000 1a2 = 0.050000 100000 1a1 = 0.002500 100000 1mma1L = 0.002500 100000 1mma2L = 0.003000 1000000 1urbation = 1.220000 1000000
Create Set	Clear	Defaults

Figure 5.5

Defining parameter values

3. Click Update Set.

4. If necessary, repeat for additional probe arrays listed.



Delete a Probe Array Type from a Userset

You may remove a probe array type from an existing userset in one of two ways.

Removing a Probe Array Type Via the Userset Definition Pane

- 1. Click to select the probe array type to delete in the **Probe array type used** list box in the userset definition pane.
- **2**. Click <<.
 - \Rightarrow The probe array type is removed from the **Probe array type used** list box and no longer appears under the userset in the left hand pane.

Removing a Probe Array Type Via the Left Hand Pane

 In the left hand pane click (+) to expand the userset of interest (Figure 5.6).



2. Click to select the probe array type to delete.

- **3.** From the menu bar select Userset \rightarrow Delete, or right-click the probe array type to delete and select Delete from the shortcut menu (Figure 5.6).
 - \Rightarrow A warning / confirmation window appears.



- 4. Click **Yes** to delete.
 - \Rightarrow The probe array type is removed from the userset.

Delete a Userset

Use the delete command to remove a userset that is no longer needed. Deleted usersets are removed in LIMS Manager and are unavailable to Microarray Suite users.

Deleting a Userset

1. In the left hand pane, click to select the userset to delete.





- 2. From the menu bar select Userset \rightarrow Delete, or right-click the userset to delete and select Delete from the shortcut menu (Figure 5.8).
 - \Rightarrow A warning / confirmation window appears (Figure 5.9).

LIMS Ma	LIMS Manager 🛛 📈		
⚠	Delete the userset New Set?		
<u>Yes</u> <u>N</u> o			
Figure 5.	9		
Delete w	arning / confirmation v	vindow	

- 3. Click Yes to delete userset.
 - ⇒ The userset is deleted and removed from both the userset list in the left hand pane and the Existing userset name(s) list box in the upper right corner.



Chapter 6





A template is a form created by the administrator, or other authorized user in Affymetrix[®] LIMS Manager, and is used to define the attributes and values for both the experiment and sample registration process. The registration process is a function of Affymetrix[®] Microarray Suite.

This chapter orients you to the layout of the **Template** tab window and describes in detail the functions and tasks accomplished here.

Template Tab Window

The window of the **Template** tab is divided into three panes (Figure 6.1). The first pane displays the active server, the process database and the **Experiment** and **Sample** templates that reside there.

The second pane in the upper right corner is the template attributes area. This pane displays the attributes and values of the template selected in the first pane.

The third pane found in the lower right corner displays the Local Task status view that displays the status of processes that may be occurring in the database, such as **Export**, **Import**, **Delete**, **Archive** and **Create**. This view allows you to monitor these jobs while concurrently working on additional tasks within the LIMS system.



The three panes of the template tab window

Templates

Creating a New Template

- **1.** Click to select either the **Experiment Templates** or **Sample Templates** folder.
- From the menu bar select Templates → Create, or right-click the templates folder and select Create from the shortcut menu that appears (Figure 6.2).



 \Rightarrow The Create Template name dialog box appears (Figure 6.3).

Create Template	
E. coli Project	
ОК	Cancel
W	

Figure 6.3 Create Template name dialog box



- 3. Type a unique name for the template type in the box and click **OK**.
 - \Rightarrow The new name appears under the appropriate template folder in the left hand pane, and the upper right hand pane displays the blank template (Figure 6.4).

Process 💼 Publish 🕞 Import 🔢 Userset 📱 Template	💱 Roles	E. coli	Project: experime	ent template attril	butes	
·		Name	Туре	Required	Values	
Cimdaffy Figure 1 Templates						_
Exp template1						_
TW Lab						_
E Sample Templates				_		_
- T Sample Temp1						
						_
reg3						
				_		_
						_
				_		_
			Apply	<u>U</u> ndo		
	Local	Task Task Desc	ription	Started	Status	
	C+Exp C+Imp	ort		10:56:07 10:56:07	Available	
	De	ete		10:56:07	Available	
	Arc	hive		10:56:07	Available	
	Cre	ate		10:56:07	Available	

4. Proceed to Defining or Editing Attributes and Values.

Defining or Editing Attributes and Values

The process of defining attributes for a new template, or editing an attribute(s) for an existing template is virtually identical. The procedure below is written as if working with a blank template, however it can also be used for the purpose of editing attributes.

Defining Attributes and Values

- 1. Click to select a template name in the first pane.
 - \Rightarrow The template appears in the second pane.
- 2. Double-click a cell under the Name column.
 - \Rightarrow The mouse pointer becomes an I-beam, and the other cells in the row are blocked (Figure 6.5).

E. coli Project: experiment template attributes				_	
Name Type Required Values					
Scientist					

Figure 6.5

Creating an attribute: Naming

- **3.** Type an attribute name and press ENTER.
 - \Rightarrow The attribute name is inserted, and a number appears in the left column.
- **4.** Click the cell under the **Type** column, then click the **•** button that appears.
- **5.** Define the data type for the attribute by selecting from the drop-down list.
 - Integer: whole number
 - Float: floating point number
 - String: limited to 255 characters
 - Date: date
 - Time: time
 - Controlled: limits user response to options listed



E. coli Project: experiment template attributes					
	Name	Туре		Required	Values
1	Scientist		•		
		Integer			
		Float			
		String			
		Controlled			

Creating an attribute: defining data type

- \Rightarrow The data type is defined and appears in the cell (Figure 6.6).
- 6. Click the cell under the **Required** column, then click the **▼** button that appears (Figure 6.7).

	E. coli Project: experiment template attributes			
	Name	Туре	Required	Values
1	Scientist	Controlled	.	
			No No	
-			- <u> </u>	



Creating an attribute: defining if required

7. Select Yes if the information is required or No if it is not.

A **Yes** selection requires the Microarray Suite user to enter a value before the sample or experiment can be registered.

- \Rightarrow The response entered appears in the cell.
- **8.** If the data type you selected was controlled, double-click the cell under the **Values** column, otherwise go to step 11.
 - \Rightarrow A Controlled Values Edit List dialog box appears (Figure 6.8).



Creating an attribute: double-click for Controlled Values Edit List dialog box

9. Define a controlled value by typing the desired value options in the box. Press ENTER after each option (Figure 6.9).

4
Cancel

Figure 6.9

Controlled Values Edit List dialog box

- **10.** Click **OK** when finished.
 - \Rightarrow Values appear in drop-down list box under the Values column (Figure 6.10).



	E. coli Project: experiment template attributes					
	Name	Туре	Required	Values		
1	Scientist	Controlled	Yes	N		
				Jokerst V		
				Kerr		
-				Meyer		

Defined controlled values

11. Click Apply.

The defined attributes do not take effect until you click Apply.

 \Rightarrow A warning message appears if this is an update to an existing template (Figure 6.11).

LIMS Manager 🛛 🗙		
There may be existing experiments that use template 'expt_temp07' Applying changes may obsolete existing attribute values. Do you wish to continue?		
	Cancel	
Figure 6	. 11 warning message	

- **12**. Click **OK** to continue.
- **13.** Repeat above steps to add additional attributes to the template.
- 14. Click Apply.
 - \Rightarrow The attributes are applied to the template.



Save time creating and defining attributes in a new template by copying the desired attribute(s) from existing templates and pasting them into a new template. See page 105.

Deactivating and Activating

Templates are used in Affymetrix[®] Microarray Suite during the experiment or sample registration process. Deactivating a template in Affymetrix[®] LIMS Manager removes it from use in Microarray Suite.

Deactivating a Template

- 1. Click to select the template to deactivate from the left hand pane.
- From the menu bar, select Template → Deactivate, or right-click the template to deactivate and select Deactivate from the shortcut menu that appears.
 - \Rightarrow A warning window appears (Figure 6.12).

LIMS Ma	anager 🔀
8	All samples using this template will be updated
Figure 6	12 te warning window

- 3. Click **OK** to deactivate.
 - \Rightarrow The template icon remains in the tree, but appears gray to indicate an inactive template (Figure 6.13).



Deactivated templates do not appear in Microarray Suite.

Activating a Template

- 1. Click to select the template to activate from the left hand pane.
- From the menu bar, select Template → Activate, or right-click the template to activate and select Activate from the shortcut menu that appears.
 - \Rightarrow The template is activated and the icon is restored to full color (Figure 6.13).




Renaming

Renaming a Template

- 1. Click to select the template to rename from the left hand pane.
- From the menu bar, select Template → Rename, or right-click the template to rename and select Rename from the shortcut menu that appears.
- 3. Type a unique template name in the box, and press ENTER.

🗸 Note

You may not duplicate an existing template name for a particular template type.

Working with Attributes

🕎 LIMS Manager - [LIMS Server Manager]						_ 8 ×
Eile View Iools Process Publish Import Userset Template	<u>Boles</u> <u>Window</u> <u>H</u> elp					_ 8 ×
INSTALL Seate DeActi	vate					
Process Publish 🕂 Import 📰 Userset		E. coli F	Project: experim	ent template attrib	outes	1
EIMS Servers	e ▶ Insert	Name	Туре	Required	Values	
🗄 😭 Cimdaffy	De-Activate 1	Scientist	Controlled	Yes		•
Experiment Templates	Copy 2	Date	Time	Yes		
T Exploring action	Cut Basto 3	Experiment Name	String	Yes		
E. coli Project	Activate 4	Lot Number	Integer	No		
🖻 🧰 Sample Templates	5	Description	String	No		
Bartell #12	¹ 6	Comments	String	No		
Bartell #15	7	Database	Controlled	Yes		
reg1						
reg2						
reg3						
1						

Figure 6.14 Templates menus and commands



Multi selecting

To select multiple contiguous attributes to act upon: click an attribute number in the left column, then SHIFT-click another number anywhere either above or below the first selected. All attributes between the two selections are selected.

To select multiple non-contiguous attributes to act upon: click an attribute number in the left column, then CTRL-click another number anywhere in the list. Repeat as necessary for additional attributes.

Inserting

Inserting an Attribute

- 1. Select the attribute below the insertion point and click the attribute number in the left column.
- From the menu bar, select Template → Attribute → Insert, or click the attribute number to select the row then right-click and select Insert from the shortcut menu that appears.
 - \Rightarrow A blank row is inserted in the template.
- **3**. Define the attribute.

For further information, refer to the section *Defining or Editing Attributes and Values* on page 97.

4. Click Apply.

Deactivating

Deactivating an Attribute

- 1. Click the attribute number in the left column.
- From the menu bar, select Template → Attribute → Deactivate, or click the attribute number to select the row then right-click and select Deactivate from the shortcut menu that appears.
 - \Rightarrow The attribute appears grayed-out, but remains in the attribute list.
- 3. Click Apply.
 - \Rightarrow A warning message appears (Figure 6.15).



Figure 6.15 Warning message

- 4. Click **OK** to deactivate the attribute.
 - \Rightarrow The attribute becomes inactive.

✓ Note

In Microarray Suite, deactivated attributes appear for the samples and experiments that have already been saved using the attribute, but cannot be modified. Deactivated attributes are not shown for new samples and experiments.

Activating

Activating an Attribute

- 1. Click the attribute number in the left column.
- From the menu bar, select Templates → Attribute → Activate, or click the attribute number to select the row then right-click and select Activate from the shortcut menu that appears.
 - \Rightarrow The attribute no longer appears gray.
- 3. Click Apply.
 - \Rightarrow The attribute becomes active.

Copying

Copying an Attribute

- 1. Click the attribute number in the left column.
- From the menu bar, select Template → Attribute → Copy, or click the attribute number to select the row then right-click and select Copy from the shortcut menu that appears.
 - \Rightarrow The information is copied to the clipboard and is available for a paste action.



Cutting

Cutting an Attribute

- 1. Click the attribute number in the left column.
- 2. From the menu bar, select **Template** \rightarrow **Attribute** \rightarrow **Cut**, or click the attribute number to select the row then right-click and select **Cut** from the shortcut menu that appears.
 - \Rightarrow The information is removed from the template, copied to the clipboard and is available for a paste action.

Pasting

Pasting an Attribute

- **1**. Click the attribute number in the left column.
- From the menu bar, select Template → Attribute → Paste, or click the attribute number to select the row then right-click and select Paste from the shortcut menu that appears.
 - \Rightarrow The information on the clipboard is pasted on the template as the last item in the list.
- 3. Click Apply.



Chapter 7





A role is a defined set of access privileges created by the administrator to control access to data by users. Creating and defining roles provides the system administrator with a security feature that allows granting various levels of access to different users, or different groups of users.

This chapter orients you to the layout of the **Roles** tab window and describes in detail the functions and tasks accomplished here.

Roles Tab Window

The window of the **Roles** tab is divided into three panes (Figure 7.1). The first pane on the left side identifies the server(s) and all the roles that have been created on that server. The branch immediately under the server lists the role names, and is known as the Role Definitions branch (Figure 7.3). LIMS Administrator and the LIMS User are default roles and cannot be deleted. The LIMS Administrator has the ability to add additional roles.

The process database name for the server is listed below the Role Definitions branch in the first pane. The branch below the process database name lists all role names that have been added to the process database. This is the User Definitions branch and is where users are added and membership to roles is made (Figure 7.3). Double-click a role name to expand the branch and view individual user names listed.

The default view for the second pane in the upper right corner lists the domains that reside on the network. Click a role name in the Role Definitions branch of the first pane and the corresponding template defining the rights and privileges of the role replaces the default view in this pane (Figure 7.2).

The third pane in the lower right corner is the Local Task status view that displays the status of processes that may be occurring in the database, such as **Export**, **Import**, **Delete**, **Archive** and **Create**. This view allows you to monitor these jobs while concurrently working on additional tasks within the LIMS system.

📴 LIMS Manager - [LIMS Server Manager]				_ 8 ×
Install	<u>Ereate Role</u> Delete Role			
🖻 Process 🔐 Publish 🔂 Import 📰 Userset 🕅 Template		unction	Other User Data Access	
Role Definitions NSTALL Mit Submitstrator Mit Mit Mit Submitstrator Mit Mit Mit Submitstrator Mit Mit Mit Mit Mit Submitstrator Mit Mit Mit Mit Mit Mit Mit Mit Mit Mit	<u>B</u> emove User	READ PROCESS DATABASE IPDATE PROCESS DATABASE UPDATE PROCESS DATABASE ARCHIVE PROCESS DATA DELETE PROLESS DATA DELETE PUBLISH DATA IMPORT PROCESS DATA EXPORT PROCESS DATA MAINTAIN ROLES DELETE SERVER TASKS MAINTAIN USERSETS CREATE PUBLISH DATABASE Local Task Task Description Feport Popot Popot Popot Popot Create	Yes Yes Yes Yes Yes Yes Yes Yes Yes Yes	2
				NUM

Figure 7.1

The three panes in the Roles tab window

🖃 🛞 Domains	Function	Other User Data Access
	READ PROCESS DATABASE	Yes
🚊 🖽 CISLAB	UPDATE PROCESS DATABASE	No
- 🔐 Domain Admins	UPDATE PUBLISH DATABASE	Yes
📿 📿 Domain Guests	ARCHIVE PROCESS DATA	No
📿 📿 Domain SDLOCAL	DELETE PROCESS DATA	Yes
📿 Domain SDREMOTE	DELETE PUBLISH DATA	Yes
💭 Domain Users	IMPORT PROCESS DATA	Yes
	EXPORT PROCESS DATA	Yes
C Seagate Info	TAKE OWNERSHIP PROCESS DATA	Yes
	MAINTAIN ROLES	Yes
	DELETE SERVER TASKS	Yes
	MAINTAIN USERSETS	Yes
	CREATE PUBLISH DATABASE	No
Figure 7.2		

Default View and Template View of the second pane of the Roles tab window

Creating a Role

A role provides a user with a set of access privileges. All LIMS users are assigned to a role. Click a role in the Role Definitions branch of the first pane (Figure 7.3) and see a template of information about that role appear in the second pane listing role functions. Every role has identical functions associated with it that are static. It is the privileges given to the user to access that function that are modified.

Creating a role is a three-fold process: naming the new role, defining the functions associated with the role, i.e., what the user can and cannot do, and defining role membership.





Creating a New Role

- 1. Click an existing role in the Role Definitions branch in the first pane.
- 2. On the menu bar select Roles \rightarrow Create Role, or right-click an existing role and select Create Role from the shortcut menu that appears.
 - \Rightarrow The Create Role dialog box appears (Figure 7.4).

Figure 7.4

The Create Role dialog box

- **3.** Type a role name in box and click **OK**.
 - \Rightarrow The new role appears in the Role Definitions branch in the first pane.

🖌 Note

A role name cannot be left blank or duplicate an existing name.

- 4. Click and drag a role name from the Role Definitions branch and drop it on top of the process database name in the User Definitions branch below.
 - \Rightarrow The new role appears in the User Definitions branch.

Defining Privileges

- 1. Click the role name in the Role Definitions branch in the first pane.
 - \Rightarrow The corresponding template defining the rights and privileges of the role appear in the second pane (Figure 7.5).

	Function	Other User Data Access
	READ PROCESS DATABASE	No
	UPDATE PROCESS DATABASE	No
	UPDATE PUBLISH DATABASE	No
	ARCHIVE PROCESS DATA	No
	DELETE PROCESS DATA	No
	DELETE PUBLISH DATA	Yes
	IMPORT PROCESS DATA	Yes
	EXPORT PROCESS DATA	No
	TAKE OWNERSHIP PROCESS DATA	Yes
	MAINTAIN ROLES	Yes
	DELETE SERVER TASKS	Yes
	MAINTAIN TEMPLATES	Yes
	MAINTAIN USERSETS	Yes
	CREATE PUBLISH DATABASE	No
J	<u> </u>	
F	igure 7.5	

Default privilege settings

The first column of boxes refers to the rights a role member is given with regards to their own data. The second column is the **Function** or privilege. The third column refers to the rights the role member is given with regards to **Other User Data Access**. Table 7.1 defines the various privilege functions.

- 2. Click to select the functions users will have with their own data.
 - \Rightarrow The selected functions appear with a check mark in the box.
- **3.** In the third column, click to toggle **Yes/No** to define the functions a user will have with **Other User Data Access**.



Other User Data Access is always Yes and cannot be toggled for the following functions: Take Ownership, Maintain Roles, Import Process Data and Delete Publish Data.

🗸 Note

The LIMS Administrator role cannot be modified.

Table 7.1

Role Privileges

Function	Description
Read Process Database	Allows a user to browse sample, experiment and analyses in formation in the process database.
Update Process Database	Allows a user to modify sample and experiment properties information in the process database.
Update Publish Database	Allows a user to overwrite experiment and analyses information in the publish database.
Archive Process Data	Allows a user to move and restore data files from and to the LIMS server.
Delete Process Data	Allows a user to delete sample, experiment, analyses, userset and probe array type information from the process database.
Delete Publish Data	Allows a user to delete experiment and analyses information from any publish database.
Import Process Data	Allows a user to migrate $\text{Affymetrix}^{\circledast}$ Microarray Suite datafiles into the process database.
Export Process Data	Allows a user to copy sample, experiment, and analyses files from the LIMS server to a defined location.
Take Ownership of Process Data	Allows a user to assume ownership of a sample, experiment and analyses information with the ability to overwrite a file with new contents (i.e., reanalyzing with different parameters).
Maintain Roles	Allows a user to manage role, role membership and set privileges.
Delete Server Tasks	Allows a user to delete publish tasks on or before a given date.
Maintain Templates	Allows a user to manage templates and their attributes.
Maintain Usersets	Allows a user to manage Expression Analysis parameter sets.
Create Publish Database	Allows a user to create a publish database.

Defining Membership

Adding a User

- 1. Click to expand the domain name tree in the upper right pane.
 - \Rightarrow The tree expands displaying the user groups residing on the domain.
- 2. Traverse the tree to find the desired user name.
- **3.** Click and drag the user name and drop it on top of the role in the User Definitions branch in the first pane.
 - \Rightarrow The user is added to the role, and the role name expands to display membership (Figure 7.6).





Click and drag the icon to collectively add an entire user group to a role.

Removing a User

- 1. Click to expand the role in the User Definitions branch of the first pane.
- 2. Click to select the user to delete.
- On the menu bar select Roles → Remove User, or right-click the user name and select Remove User from the shortcut menu.
 - \Rightarrow The user name is removed from role membership.



Deleting Roles

Note

LIMS User and **LIMS Administrator** are default roles and cannot be deleted.

Deleting a Role

- **1**. Click to select the role to delete.
- On the menu bar select Roles → Delete Role, or right-click the role to delete and select Delete Role from the shortcut menu.
 - \Rightarrow A warning window appears (Figure 7.7).

LIMS Manager 🛛 🔀
Delete role TechSupport?
Yes <u>N</u> o
Figure 7.7
LIMS warning window

- 3. Click Yes to delete.
 - \Rightarrow The role is deleted and disappears from both the Role Definitions and the User Definitions branches in the left pane.



Appendix A





The following are samples of two types of print output available in LIMS Manager.

Diagram Only

AFFYMETRIX													
\mathcal{N}			ProcessView	Report of	fI	Exp	eriment	:Wisc	Control	(Diagram	Only)	5/15/01	12:33:37
Sample Name	· T	IOF W		Date			Mar 30 2	000 10	:21AM	Server		DBSERVER04	
Sample Project	t: U	Jof W		Stage		:	Hybridiz	ation		Database		LIMS3	
Sample Type	: t	otal	RNA	Template	:	:				User	:	lwang	
Comments	:												
Description													



Figure A.1 Print Diagram Only

Diagram & Properties

		ProcessView	Report	of	Experiment:Wisc	Control		5/15/01 12:33:56
Sample Name : Sample Project:	UOF W Uof W	Date Stage	:	Mar Hyb:	30 2000 10:21AM ridization	Server Database	:	DBSERVER04 LIMS3
Sample Type : Comments : Description :	total RNA	Templat	e :	-		User	:	lwang

ß	Wisc Control <experiment></experiment>
	Probe Array Type : Hu6800
	Comment : (null)
	User : lwang
	Date : Mar 30 2000 10:21AM
	Auto : OFF
	Template :
	Sample Registration
	(O) 🗐 Uof W <sample></sample>
	Uof W:GE <vessel></vessel>
	Comment :
	Barcode : NEW2
	Type : Tube
	🕘 Experiment Setup
	(I) 🗐 Uof W <sample></sample>
	(0) 🛛 Wisc Control <experiment></experiment>
	Hybridization
	Operator : lwang
	Hyb Date : Jan 25 2000 11:35
	Hyb Comments :
	Hyb Description :
	Hyb Station ID : 0
	Hyb Station # : 1
	Hyb Protocol : EukGE-WS2
	Hyb Module : 1
	Hyb Stages : Completed
	Hyb Reagents :
	Hyb LotNumber :
	(I) 🏢 Uof W:GE <vessel></vessel>
	Wisc_Control <chip></chip>
	Probe Array Lot# : 97957
	Barcode : Wisc_Control
	Probe Array Type : Hu6800
	0 Scan
	Operator : iwang
	Scan Date : Jan 25 2000 10:44AM
	Scan Comments :
	Scaller Type : HP
	Scannel ID : Scan Filter : 570
	Scan Temperature . 0
	Scan Pixel Size : 3.000000
	Scan Line Length : 0.00000
	Scan # of Lines : 0
	Scan Horizontal Offset : 0.000000
	Scan Vertical Offset : 0.000000
	# of Scans : 2
	(I) 🝙 Wisc Control <chip></chip>
	(0) Wisc Control.dat
	Path : \\NTLIMS04\GCLims\Data\Wisc_Control.dat
	Barcode : Wisc_Control
	Created By : lwang
	Edited By : lwang
Ц	

Figure A.2

Print Diagram and Properties (page 1 of 4, only)



Appendix B





This appendix lists frequently asked questions (FAQs) regarding Affymetrix[®] LIMS Manager. If after reviewing these questions you still need assistance, please contact Affymetrix Technical Support.

LIMS FAQs

- **Q:** I want to view information about other expression samples. I try to select another sample by clicking its icon but the software refuses to switch to a different sample. What should I do?
- **A:** Throughout the program, click the text representing the sample or experiment name rather than the icon.
- Q: How can I change the "export to" or "import from" location?
- A: On the menubar, click the icon that depicts a folder opening (see page 18). If you are in the **Process**, **Publish**, or **Find** tabs, this will select a target location for exported files. If you are in the **Import** tab, this will select a location to import files.
- **Q:** How do I use the **Find** command to search the Publish database?
- A: The Find command only works with the Process database.
- **Q:** Using the **Find** command, I know some of my samples (projects or experiments) are in the process database. However, the results continually produce an empty set. Why does this happen?
- A: In the Find window, the Select Object tab, the Advanced Filter tab, and the Data Filter tab are additive. Thus if you select experiments with the text string *"xxxAssay"* using the Select Object tab and then limit the search to *Project X* using

the **Advanced Filter** tab, then the **Find** command is looking for experiments with the text string "xxxx" that are members of Project X.

- **Q:** I want to delete (assume ownership, export, or archive) multiple files (experiments or samples) at one time. However I cannot select multiple files at one time within the process database. How can I accomplish this?
- A: You can SHIFT-select or control-select multiple files, samples, or experiments at one time within the Find window. The majority of the functions available within the Process tab are also available within the Find window.
- **Q:** I want to archive a sample but the software won't let me choose that function. How do I archive a sample?
- A: Currently, archiving can be performed on experiments and files, but not samples. A list of available functions within the **Find** window is listed below:

Files: archive, assume ownership, delete, and export

Experiments: assume ownership, delete, export and rename

Samples: assume ownership, delete, and rename

- Q: How do I get out of the Find window and back to the process database?
- A: Simply close **Find** window or go to the Windows menu and select "LIMS server manager."
- Q: Can I view property information about files in the publish database?
- A: No, this information is not available. But you are able to view property information about files in the process database from either the **Publish** or the **Process** tabs.
- **Q:** I can't find my .dat file under the LIMS **Import** tab. I have confirmed that the .dat file exists on the local drive and that I have the import location mapped correctly. Why doesn't LIMS let me import my .dat file?
- A: LIMS will import .exp files and the associated .dat files with the same filename. Make certain that an .exp file that contains the exact same filename is contained within the same folder as your .dat file. As a last resort, open your .dat file in

Affymetrix[®] Microarray Suite. If an .exp file does not exist for that .dat, Affymetrix[®] Microarray Suite will automatically create one.

- **Q:** When I try to open a Sample folder, only a few samples are displayed and I know that there should be more in the database. What happened to the samples?
- A: The Sample filter remains active during successive sessions. Reset Sample Filter parameters. See *Sample Filter*, on page 47.





Α

activate attribute 105 template 101 active server 16, 39 Affymetrix software applications 15 data flow 14 Data Mining Tool 15 Technical Publications 24 technical support 24 archive data 30

assume ownership 32

attribute

activate 105 controlled values 98 copy 105 deactivate 104 define 97 edit 97 insert 104 shortcut tip 100 types 97

С

controlled value 98 create sample 71 vs. register sample 71

D

data archive 30 assume ownership 32 export 35 import 70 security 12 unarchive 31 database accessing 55 create 53 create Oracle publish 63

create publish 60 create SQL publish 61 security 55 size considerations 60 deactivate attribute 104 template 101 delete probe array type 34 probe array type from publish database 56 probe array type from userset 88 process data 32 publish data 57-59 userset 89 dependent parameter 87 documentation conventions used 22 on-line 22

E

edit sample 75 export process data 35

F

file path archive 30 default 70 export 35 icon 18 import 70 file type 18 file view 36 filter find function 46 samples import tab 76 process tab 47 find 45-49 shortcut commands 49 window 46

G

glossary 18

L

icons 17 import data 70-75 associating files 73 creating a sample 71 identifying a sample 71 select file path 70 selecting a sample 72 steps 70 import tab 67 appearance of 68 remove 76 remove all 76 terms used 20 independent parameter 85 installing LIMS Manager 4

L

LIMS Manager and Affymetrix software 13 FAQs 123 features 11 installing 4 interface orientation 16 requirements 3 terms 18 LIMS Sync 55

М

manual on-line 22

0

on-line documentation 22 Oracle creating database 63

Ρ

print diagram & properties 38 diagram only 38 samples 119

privilege settings default 113 functions defined 114

probe array type delete from process database 34 delete from userset 88

process tab 27

appearance of 28 change views 36 data types 29 file view 36 functions 29 process view 36 shortcut menus 29 terms used 19 views 27

process view 36

publish database 60-64 create 60-64 create a SQL 61 create an Oracle 63 create Oracle 63 create SQL 61 delete probe array type 56 size considerations 60

publish tab 53 appearance of 54 delete publish data 57

R

remove all 76 requirements 3 role add a user 115 create role 111 define membership 115 define privileges 113 definition of 111 delete a role 116 delete a user 115 roles tab 109 appearance of 110 privilege functions defined 114 terms used 21

S

sample edit 75 samples filter 47, 76 server

register 39 unregister 40

server events 40

server tasks 40-45 cancel an active task 45 restart a cancelled task 45 view all items 43 view task 42 view task items 43

SQL database

create 61

т

technical support 24 template activate 101 creating new 95–96 deactivate 101 define attribute 97–100 template tab 93

appearance of 94 terms used 21

terminology 18 toolbar 18

υ

unarchive data 31 userset creating 84 defined 83 defining dependent parameter 87 defining independent parameter 85 delete 89 modify 85 usersets tab 81 appearance of 82 definition pane 83 terms used 21

W

window general orientation 16 icons 17 import tab 67 process tab 27 publish tab 53 roles tab 109 template tab 93 toolbar 18 userset tab 81

workflow-based tracking 12