

Xcalibur[®]

SEQUEST[®] Cluster

Installation and Operation Manual

XCALI-97160 Revision B

July 2007

© 2007 Thermo Fisher Scientific Inc. All rights reserved.

BioWorks™ and Finnigan™ are trademarks and Xcalibur® is a registered trademark of Thermo Fisher Scientific. Microsoft® and Microsoft Windows® are registered trademarks of Microsoft Corporation. SEQUEST®/SEQUEST® and ICAT® are registered trademarks of the University of Washington. PVM™ is a trademark of Oak Ridge National Laboratory.

Thermo Fisher Scientific Inc. provides this document to its customers with a product purchase to use in the product operation. This document is copyright protected and any reproduction of the whole or any part of this document is strictly prohibited, except with the written authorization of Thermo Fisher Scientific Inc.

The contents of this document are subject to change without notice. All technical information in this document is for reference purposes only. System configurations and specifications in this document supersede all previous information received by the purchaser.

Thermo Fisher Scientific Inc. makes no representations that this document is complete, accurate or error-free and assumes no responsibility and will not be liable for any errors, omissions, damage or loss that might result from any use of this document, even if the information in the document is followed properly.

This document is not part of any sales contract between Thermo Fisher Scientific Inc. and a purchaser. This document shall in no way govern or modify any Terms and Conditions of Sale, which Terms and Conditions of Sale shall govern all conflicting information between the two documents.

Release history:

Software requirements: BioWorks 3.3.1 SP1 and Xcalibur 2.0.6.

For Research Use Only. Not regulated for medical or veterinary diagnostic use by U.S. Federal Drug Administration or other competent authorities.

WEEE Compliance

This product is required to comply with the European Union's Waste Electrical & Electronic Equipment (WEEE) Directive 2002/96/EC. It is marked with the following symbol:



Thermo Fisher Scientific has contracted with one or more recycling or disposal companies in each European Union (EU) Member State, and these companies should dispose of or recycle this product. See www.thermo.com/WEEERoHS for further information on Thermo Fisher Scientific's compliance with these Directives and the recyclers in your country.

WEEE Konformität

Dieses Produkt muss die EU Waste Electrical & Electronic Equipment (WEEE) Richtlinie 2002/96/EC erfüllen. Das Produkt ist durch folgendes Symbol gekennzeichnet:



Thermo Fisher Scientific hat Vereinbarungen mit Verwertungs-/Entsorgungsfirmen in allen EU-Mitgliedsstaaten getroffen, damit dieses Produkt durch diese Firmen wiederverwertet oder entsorgt werden kann. Mehr Information über die Einhaltung dieser Anweisungen durch Thermo Fisher Scientific, über die Verwerter, und weitere Hinweise, die nützlich sind, um die Produkte zu identifizieren, die unter diese RoHS Anweisung fallen, finden sie unter www.thermo.com/WEEERoHS.

Conformité DEEE

Ce produit doit être conforme à la directive européenne (2002/96/EC) des Déchets d'Equipements Electriques et Electroniques (DEEE). Il est marqué par le symbole suivant:



Thermo Fisher Scientific s'est associé avec une ou plusieurs compagnies de recyclage dans chaque état membre de l'union européenne et ce produit devrait être collecté ou recyclé par celles-ci. Davantage d'informations sur la conformité de Thermo Fisher Scientific à ces directives, les recycleurs dans votre pays et les informations sur les produits Thermo Electron qui peuvent aider la détection des substances sujettes à la directive RoHS sont disponibles sur www.thermo.com/WEEERoHS.

Contents

Preface	xi
About This Guide.	xi
Related Documentation	xi
Safety and Special Notices	xi
Contacting Us	xii
Assistance	xii
Changes to the Manual and Online Help	xii
 Chapter 1 Introduction to Clusters	 1
What are Clusters?.	1
How SEQUEST Cluster Works	2
 Chapter 2 Setting Up the SEQUEST Cluster	 5
Required Software	6
Customer Supplied Software	6
Thermo Scientific Supplied Software.	6
Setting Up the Master Node	6
Installing and Configuring Microsoft Windows 2000 Server.	7
Installing and Configuring Microsoft Windows 2003 Server.	12
Installing Master Node Software	16
Setting Up the Slave Nodes: Microsoft Windows 2000 Professional	20
Installing and Configuring Microsoft Windows 2000 Professional	20
Installing Slave Node Software	25
Updating the Microsoft Windows Registry	27
Setting Up the Slave Nodes: Microsoft Windows XP Professional.	28
Installing and Configuring Microsoft Windows XP Professional.	28
Installing Slave Node Software	32
Setting Up the Slave Nodes: Red Hat Linux (Version 8.0)	35
Installing Red Hat Linux 8.0.	35
Configuring Red Hat Linux 8.0	37
Setting Up the Slave Nodes: Red Hat Linux (Version 9.0)	40
Installing Red Hat Linux 9.0.	40
Configuring Red Hat Linux 9.0	42
Setting Up the Slave Nodes: Red Hat Enterprise Linux 3.0.	45
Installing Red Hat Enterprise Linux 3.0 ES.	45
Configuring Red Hat Enterprise Linux 3.0 ES	47

Setting Up the Slave Nodes: Red Hat Enterprise Linux 4.0	49
Installing Red Hat Enterprise Linux 4.0 ES	49
Configuring Red Hat Enterprise Linux 4.0 ES	51
Setting Up the Slave Nodes: Fedora 3.0	53
Installing Fedora 3.0	54
Configuring Fedora 3.0	56
Setting Up the Slave Nodes: Fedora 4.0	58
Installing Fedora 4.0	58
Configuring Fedora 4.0	60
Setting Up the Slave Nodes: Fedora 5.0	62
Installing Fedora 5.0	62
Configuring Fedora 5.0	64
Setting Up the Slave Nodes: Fedora 6.0	67
Installing Fedora 6.0	67
Configuring Fedora 6.0	69
Updating from SEQUEST Cluster 3.1 or SEQUEST Cluster 3.1 SR1 to SEQUEST Cluster 3.3.1	71
Final Configuration	72
Configuration Using the Administrator Login	72
Configuration Using the Proteomics Login	78
After the Final Configuration	80
Files Modified During the Installation	80
 Chapter 3 Using the SEQUEST Cluster	81
SEQUEST Cluster Manager	81
Starting and Stopping the PVM Software	82
Managing Databases	83
Submitting SRF Files to the Cluster for Processing	84
Configuring BioWorks for Use with SEQUEST Cluster	85
Configuring BioWorks on the Cluster Master Node	86
Configuring BioWorks on a Remote Client Computer	88
Running a SEQUEST Search on the Cluster	90
Starting a Search Using the SEQUEST Search Dialog Box	90
Starting a Search Using the SEQUEST Batch Search Dialog Box	93
Starting a Search Using the Search Selected Samples Dialog Box	94
PVM Console	96
 Chapter 4 Troubleshooting	101
Computer Name and Login Account Name Are the Same	102
Account Name and WRSHD Login Are Not the Same	102
AutoLogon Feature of Microsoft Windows 2000 Is Not Set	102
Need Additional Linux Subsets	103
Extra Configuration Entries in WRSHD	104
WRSHD Cannot Create Service	104
WRSHD Will Not Start on Microsoft Windows 2000 / Windows XP	105

RSH Does Not Work Between Microsoft Windows 2000 Server / Windows 2003 Server and Linux Nodes	105
RSH Is Not Working Correctly.....	106
PVM Will Not Start	108
Extra Files From an Aborted PVM Session	109
Poor Cluster Performance	110
Cluster Searches Will Not Run Through SEQUEST Queue.....	110
Appendix A Installation Checklists	111
Master Computer Checklist: Windows 2000 Server	112
Master Computer Checklist: Windows 2003 Server	113
Slave Computer: Microsoft Windows 2000 Checklist.....	114
Slave Computer: Microsoft Windows XP Checklist.....	115
Slave Computer: Red Hat Linux / Fedora Checklist	116

Preface

About This Guide

Welcome to Thermo Fisher Scientific, Xcalibur SEQUEST® Cluster. This SEQUEST Cluster Installation and Operation manual provides information on how to set up and use your SEQUEST Cluster with BioWorks 3.3.1 SP1.

Related Documentation

In addition to this guide, Thermo Fisher Scientific provides the following documents for SEQUEST Cluster:

- BioWorks User Guide

Safety and Special Notices

Make sure you follow the precautionary statements presented in this guide. The safety and other special notices appear in boxes.

Safety and special notices include the following:



CAUTION Highlights hazards to humans, property, or the environment. Each CAUTION notice is accompanied by an appropriate CAUTION symbol.

IMPORTANT Highlights information necessary to prevent damage to software, loss of data, or invalid test results; or may contain information that is critical for optimal performance of the system.

Note Highlights information of general interest.

Tip Helpful information that can make a task easier.

Contacting Us

There are several ways to contact Thermo Fisher Scientific.

Assistance

For new product updates, technical support, and ordering information, contact us in one of the following ways:

Visit Thermo Scientific on the Web

www.thermo.com/finnigan

Contact Technical Support

Phone: 1-800-685-9535
Fax: 1-561-688-8736
E-mail: TechSupport.C+MS@thermofisher.com
Knowledge base: <http://www.thermokb.com>

Find software updates and utilities to download at <http://sjsupport.thermo.com>.

Contact Customer Service

For ordering information in the U.S. and Canada:

Phone: 1-800-532-4752
Fax: 1-561-688-8731
Web site: www.thermo.com/finnigan

Changes to the Manual and Online Help

To suggest changes to this guide or to the Help, use one of the following:

- Fill out a reader survey online at www.thermo.com/lcms-techpubs.
- Send an e-mail message to the Technical Publications Editor at techpubs.finnigan-lcms@thermofisher.com.

Introduction to Clusters

The SEQUEST® Cluster is an optional utility that is accessible from BioWorks™, a Thermo Scientific post-processing software application. SEQUEST Cluster is licensed separately from the main BioWorks application. This chapter provides a basic introduction to cluster technology.

Contents

- [What are Clusters?](#)
- [How SEQUEST Cluster Works](#)

What are Clusters?

Cluster-based supercomputers have been around since the mid 1990s. These systems have provided a more economical path to extreme computing power for larger groups of users than was possible with the earlier proprietary supercomputers.

The basic setup of a cluster involves multiple computers that are connected together using networking technology and specialized software such as Parallel Virtual Machine (PVM™) or Message Passing Interface (MPI).

The design of the SEQUEST Cluster follows the “Beowulf” design model. Beowulf was the outgrowth of the research work of Thomas Sterling and Donald Becker at NASA. The Beowulf design model is a design for high-performance parallel computing clusters on inexpensive personal computer hardware.

Clusters have been built by using nearly every computer operating system that is currently available. Thermo Scientific supports several operating systems with the SEQUEST Cluster product: Microsoft® Windows® 2000, Microsoft Windows XP, Red Hat Linux, Red Hat Enterprise Linux, and Fedora.

How SEQUEST Cluster Works

SEQUEST Cluster software uses PVM software to pass data from one node in the cluster (the *master*, or management node) to another node or nodes in the cluster (the *slave*, or compute nodes).

The master node of the SEQUEST Cluster contains all of the data files from mass spectrometers and FASTA databases, plus the DTA and OUT or SRF files that are generated by SEQUEST. The slave nodes contain only the FASTA databases. All of the search data that is needed by the slave nodes is passed between the master and slave nodes through software connections known as sockets. Thus, this data is never stored on the hard disk of a slave nodes.

When you start SEQUEST search on the master node of the cluster, the program *sequest_master* is launched. This program identifies all of the slave nodes that have been configured and licensed to run with the SEQUEST Cluster and then launches at least one copy of the *sequest_slave* program on these nodes.

The actual number of copies of *sequest_slave* that are launched on each slave node depends upon how many CPUs are present and on how many CPUs are licensed through Thermo Scientific. If these systems are dual CPU computers, then two copies of *sequest_slave* are launched on each slave node. The master node of the cluster also starts a single copy of the *sequest_slave* program, even if it has two CPUs. For example, in a 16-node cluster where each node contains a dual CPU computer, there would be 33 instances of *sequest_slave* running.

The *sequest_master* program reads the SEQUEST search parameters file (*sequest.params* file), and then passes the search parameters to the slave nodes. When all of the slave nodes are ready, the *sequest_master* program starts reading the DTA files and passes this information to the slave nodes. Because the master computer waits to be requested for information by the slave nodes before sending the information, the *sequest_master* program is said to be acting in a passive mode.

As each slave node receives the DTA data from the master node, the master node keeps a record of which node has what data. When a *sequest_slave* program is finished processing the data, it sends a message back to the master node that it has finished processing that DTA data, and then it asks for more data. The *sequest_master* program updates the internal record for the previous data to indicate that the data has been processed, and then sends the information from another DTA file to that slave node. This process is repeated until all of the DTA files have been processed.

At this point, the *sequest_master* program checks the internal list of processed DTA files to ensure that all of the files are completely processed. If any of the files are not completely processed, this indicates a problem with one of the slave nodes. The *sequest_master* then resubmits these DTA files to other nodes in the cluster to be processed again.

The design of the *sequest_master* program helps the overall processing of the DTA files in the event that any nodes in the cluster became unavailable. If a node becomes unavailable, the *sequest_master* program does not receive a message requesting more data from the node, and

therefore does not try to send any more DTA data to this node. While this occurrence does result in a longer search time (because there are fewer nodes processing the data), it does not cause the entire search to fail.

In addition, the *sequest_master* programs receives the contents of the OUT file from the slave nodes through PVM and then writes the output to the hard disk of the master node. The slave nodes do not directly write the output to the hard disk of the master node because writing files across a network connection can be extremely slow, which would slow down DTA file processing.

Thermo Scientific recommends that the FASTA databases be stored on each node of the cluster instead of using a shared directory or folder on the master node of the cluster. If you use a shared directory or folder for the FASTA databases, you might experience nodes dropping out of the search. This increases the overall time for the search.

Setting Up the SEQUEST Cluster

This chapter guides you through the process of setting up the SEQUEST Cluster.

Contents

- Required Software
- Setting Up the Master Node
- Setting Up the Slave Nodes: Microsoft Windows 2000 Professional
- Setting Up the Slave Nodes: Microsoft Windows XP Professional
- Setting Up the Slave Nodes: Red Hat Linux (Version 8.0)
- Setting Up the Slave Nodes: Red Hat Linux (Version 9.0)
- Setting Up the Slave Nodes: Red Hat Enterprise Linux 3.0
- Setting Up the Slave Nodes: Red Hat Enterprise Linux 4.0
- Setting Up the Slave Nodes: Fedora 3.0
- Setting Up the Slave Nodes: Fedora 4.0
- Setting Up the Slave Nodes: Fedora 5.0
- Setting Up the Slave Nodes: Fedora 4.0
- Final Configuration
- After the Final Configuration
- Files Modified During the Installation

IMPORTANT When you load the operating system and configure the master and slave nodes, consistency is extremely important. Set up each slave node in an identical fashion. This will make subsequent maintenance of the system much easier.

Required Software

To set up the SEQUEST cluster, you need the software listed in this section.

Customer Supplied Software

Some or all of this software may be available from your computer hardware vendor:

- Master node:

Microsoft Windows 2000 Server with Service Pack 4 or Microsoft Windows 2003 Server with Service Pack 1

- Slave nodes:

Microsoft Windows 2000 Professional with Service Pack 4 or Microsoft Windows XP Professional with Service Pack 2

or

Red Hat Linux 8.0 or 9.0, or Red Hat Enterprise Server 3.0 or 4.0, or Fedora 3.0 or 4.0

Thermo Scientific Supplied Software

The following software is supplied by Thermo Fisher Scientific:

- Parallel Virtual Machine (PVM™) 3.4.3 for Microsoft Windows 2000/XP
- WRSHD RSH for Microsoft Windows NT/2000/XP
- Xcalibur® Bioworks™ 3.3.1 SP1
- SEQUEST Cluster 3.3.1

Note Xcalibur 1.4 SR1 or 2.0.6 is required to install Bioworks 3.3.1 SP1, but it is not needed to run the SEQUEST Cluster.

Setting Up the Master Node

The master node must be configured with the Microsoft Windows 2000 Server or Microsoft Windows 2003 Server operating system.

This section assumes that the master node has not been configured with the Microsoft Windows 2000 Server or Microsoft Windows 2003 Server operating system. If the master node has already been configured, then skip the corresponding steps in the procedure below.

This section also assumes that you are familiar with installing Microsoft Windows 2000 Server or Microsoft Windows 2003 Server.

This section contains the following topics:

- [Installing and Configuring Microsoft Windows 2000 Server](#)
- [Installing and Configuring Microsoft Windows 2003 Server](#)
- [Installing Master Node Software](#)

Installing and Configuring Microsoft Windows 2000 Server

❖ To install and configure the Microsoft Windows 2000 Server operating system

1. Install Microsoft Windows 2000 Server on the master computer.

More information about installing Microsoft Windows 2000 Server can be found in Chapters 3 and 4 of the *Getting Started* manual that comes with Microsoft Windows 2000 Server.

2. Install additional drivers for Microsoft Windows 2000 Server to recognize the hard disk(s) in your system, if required. If any special drivers are required, then consult the manual that comes with the computer hardware.
3. Define disk partitions for the operating system: Create a small 4 GB partition for the operating system, and then commit the rest of the space on the hard disk to a separate partition that will store the FASTA databases, and DTA and OUT or SRF files. You will set up this unused portion of the drive later.
4. Format the drive using the NTFS file system.

After you install the server software, the computer restarts and opens the Microsoft Windows 2000 Server Setup program.

5. Use the Microsoft Windows 2000 Server Setup program to set up the server software.

During the setup, you are asked to enter the base operating system license, license mode and number of licenses that were purchased. Additional information on choosing the license mode can be found on pages 46-47 and page 76 of the *Microsoft Windows 2000 Server Getting Started* manual.

Note When you choose a name for the system during configuration, do not use your user account name, or the PVM software will not work correctly.

Do not configure this server to be a network server (that is, DHCP, DNS, WINS, Print Server, Application Server, and so on) for the other computers on your network. If you want to run Proteomics Browser on the master computer, then you need to install the Internet Information Server (IIS) software. Refer to the *BioWorks Installation Guide* for more information on installing IIS and Proteomics Browser.

The Microsoft Windows 2000 Server Setup program restarts the computer after all of the basic setup has been completed.

6. Configure the Microsoft Windows 2000 Server software:

- a. Log into the master computer.
- b. Configure the network cards.

Configure the primary network card to work with your corporate computer network, and configure the secondary network card to work with the dedicated network setup between the master and slave computers.

Note When you are configuring the network, do not disable the Client for Microsoft Networks option. This option is required to share drives between individual computers that are running Microsoft Windows 2000.

Use a private address space number for the network containing the slave computers (that is, a secondary network card). For example, use 192.168.66.250 with the master computer. Then, use the IP addresses 192.168.66.1 through 192.168.66.xxx with each second slave computer that is part of the cluster.

- c. Because the slave computers will not be a part of a domain, edit the `C:\WINNT\system32\drivers\etc\hosts` file and add the name and IP address for each slave node in your cluster as shown in [Figure 1](#).

7. Install Service Pack 4 for Microsoft Windows 2000 Server. After the installation is complete, the computer restarts.

8. Log in as *administrator*.

Figure 1. Master node hosts file

```
# Copyright (c) 1993-1999 Microsoft Corp.
#
# This is a sample HOSTS file used by Microsoft TCP/IP for Microsoft Windows.
#
# This file contains the mappings of IP addresses to host names. Each
# entry should be kept on an individual line. The IP address should
# be placed in the first column followed by the corresponding host name.
#
# The IP address and the host name should be separated by at least one
# space.
#
# Additionally, comments (such as these) may be inserted on individual
# lines or following the machine name denoted by a '#' symbol.
#
127.0.0.1 localhost
192.168.66.250master
192.168.66.1node1
192.168.66.2node2
192.168.66.3node3
192.168.66.4node4
192.168.66.5node5
192.168.66.6node6
192.168.66.7node7
192.168.66.8node8
```


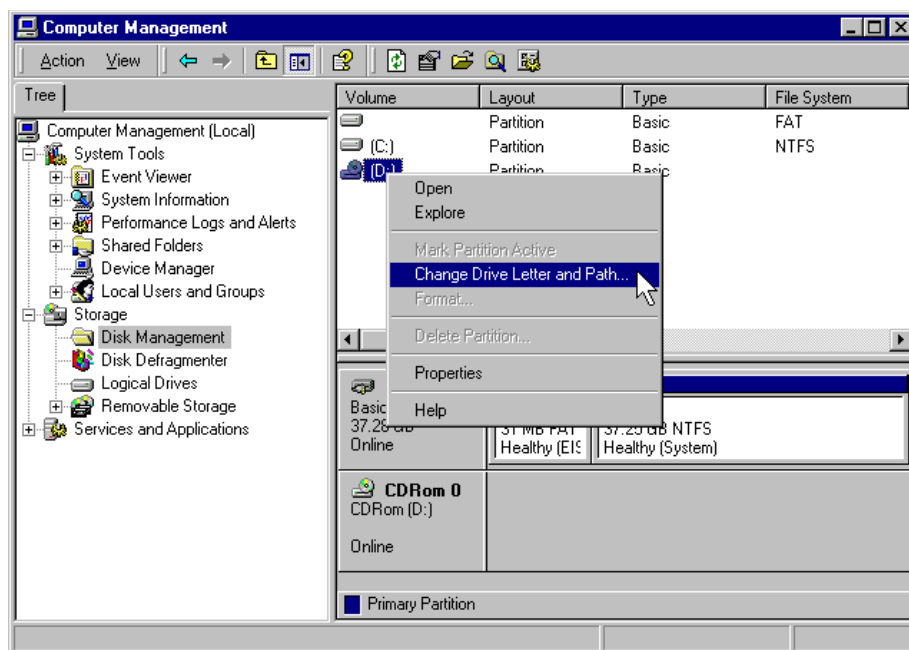
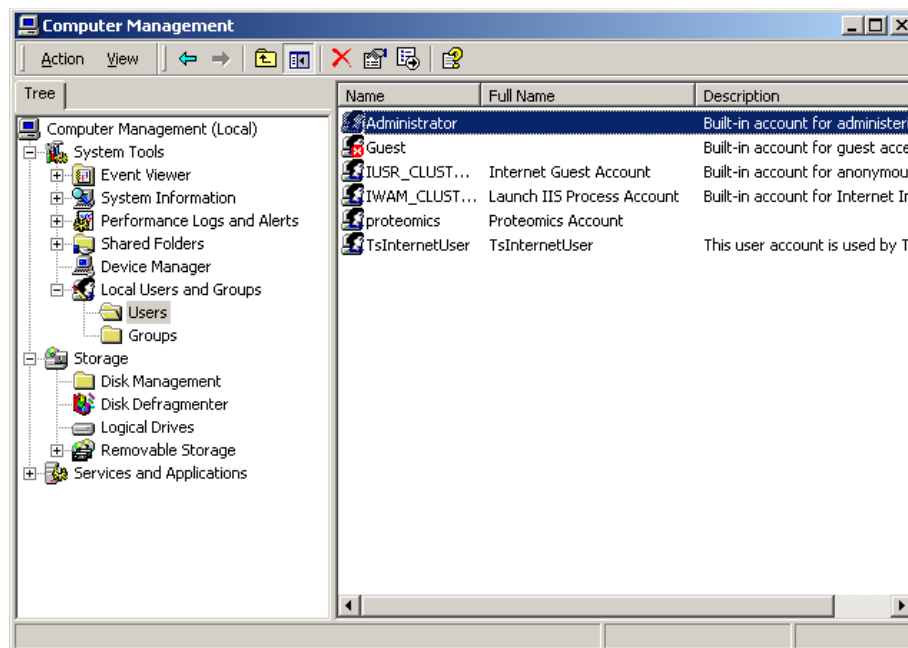
9. Format the unused portion of the hard disk and set the drive letter for this partition to D:\ as follows:
 - a. Choose  **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open the Administrative Tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 2](#).
 - d. Format the drive using the NTFS file system.
 - e. In the right pane, right-click on the CD-ROM (D:) entry and choose **Change Drive Letter and Path** from the shortcut menu.
 - f. Select an open/unused drive letter for the CD-ROM. You can now format the unused portion of the hard drive as D:\.

Figure 2. Computer Management tool



10. After you format the D:\ drive, create a directory named *Database* on that drive. You will use this directory to store the FASTA databases that are needed by SEQUEST.
11. Configure the graphics card.
12. Create a new user named *proteomics*:
 - a. Choose **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open the Administrative Tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 3](#).
 - d. In the left pane, double-click **Local Users and Groups**, and then select **Users**.

Figure 3. Computer Management tool



- e. Choose **Action > New User** and fill in the information in the dialog box. Enter “proteomics” as a user name. Click **Create**.
13. Add the new user proteomics to the Power Users and Administrators groups.
 - a. Click **proteomics** tab.
 - b. Choose **Action > Properties** to open the Proteomics Properties dialog box.
 - c. Click **Member Of** tab and click **Add** to open the Select Groups dialog box.
 - d. In the upper pane, click on the Power Users group to select it, and then click **Add**.
 - e. Click **OK** to close the Select Groups dialog box.
 - f. Click **OK** to close the Proteomics Properties dialog box.
 - g. Close the Computer Management tool.
 14. Install the Thermo Fisher Scientific-supplied software for the master node, use the procedure described in [“Installing Master Node Software”](#) on [page 16](#).

Installing and Configuring Microsoft Windows 2003 Server

❖ To install and configure the Microsoft Windows 2003 Server operating system

1. Install Microsoft Windows 2003 Server on the master computer.

More information about installing Microsoft Windows 2003 Server can be found in Chapters 3 and 4 of the *Getting Started* manual that comes with Microsoft Windows 2003 Server.

2. Install additional drivers for Microsoft Windows 2003 Server to recognize the hard disk(s) in your system, if required. If any special drivers are required, then consult the manual that comes with the computer hardware.
3. Define disk partitions for the operating system: Create a small 6 GB partition for the operating system, and then commit the rest of the space on the hard disk to a separate partition that will store the FASTA databases, and DTA and OUT or SRF files. You will set up this unused portion of the drive later.
4. Format the drive using the NTFS file system.

After you install the server software, the computer restarts and opens the Microsoft Windows 2003 Server Setup program.

5. Use the Microsoft Windows 2003 Server Setup program to set up the server software.

During the setup, you are asked to enter the base operating system license, license mode and number of licenses that were purchased. Additional information on choosing the license mode can be found on pages 46-47 and page 76 of the *Microsoft Windows 2003 Server Getting Started* manual.

Note When you choose a name for the system during configuration, do not use your user account name, or the PVM software will not work correctly.

Do not configure this server to be a network server (that is, DHCP, DNS, WINS, Print Server, Application Server, and so on) for the other computers on your network. If you want to run Proteomics Browser on the master computer, then you need to install the Internet Information Server (IIS) software. Refer to the *BioWorks Installation Guide* for more information on installing IIS and Proteomics Browser.

The Microsoft Windows 2003 Server Setup program restarts the computer after all of the basic setup has been completed.

6. Configure the Microsoft Windows 2003 Server software.
 - a. Log into the master computer.
 - b. Configure the network cards.

Configure the primary network card to work with your corporate computer network, and configure the secondary network card to work with the dedicated network setup between the master and slave computers.


Note When you are configuring the network, do not disable the Client for Microsoft Networks option. This option is required to share drives between individual computers that are running Microsoft Windows 2003.

Use a private address space number for the network containing the slave computers (that is, a secondary network card). For example, use 192.168.66.250 with the master computer. Then, use the IP addresses 192.168.66.1 through 192.168.66.xxx with each second slave computer that is part of the cluster.

- c. Because the slave computers will not be a part of a domain, edit the `C:\WINNT\system32\drivers\etc\hosts` file and add the name and IP address for each slave node in your cluster as shown in [Figure 4](#).
7. Install Service Pack 1 for Microsoft Windows 2003 Server. After the installation is complete, the computer restarts.
8. Log in as *administrator*.

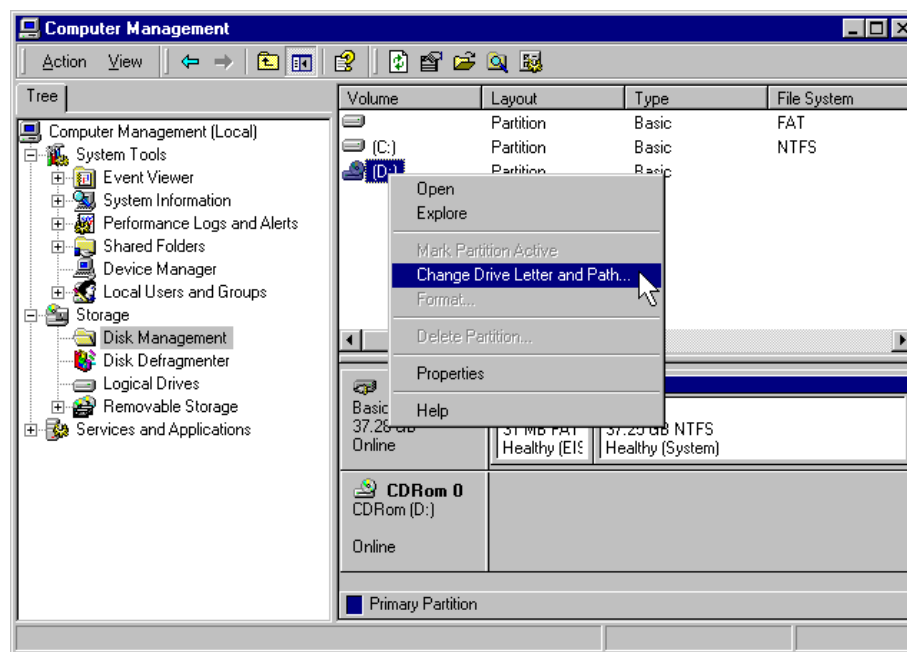
Figure 4. Master node hosts file

```
# Copyright (c) 1993-1999 Microsoft Corp.
#
# This is a sample HOSTS file used by Microsoft TCP/IP for Microsoft Windows.
#
# This file contains the mappings of IP addresses to host names. Each
# entry should be kept on an individual line. The IP address should
# be placed in the first column followed by the corresponding host name.
#
# The IP address and the host name should be separated by at least one
# space.
#
# Additionally, comments (such as these) may be inserted on individual
# lines or following the machine name denoted by a '#' symbol.
#
127.0.0.1 localhost
192.168.66.250master
192.168.66.1node1
192.168.66.2node2
192.168.66.3node3
192.168.66.4node4
192.168.66.5node5
192.168.66.6node6
192.168.66.7node7
192.168.66.8node8
```

9. Format the unused portion of the hard disk and set the drive letter for this partition to D:\.
 - a. Choose  **Start > Settings > Control Panel** to open the Control Panel.

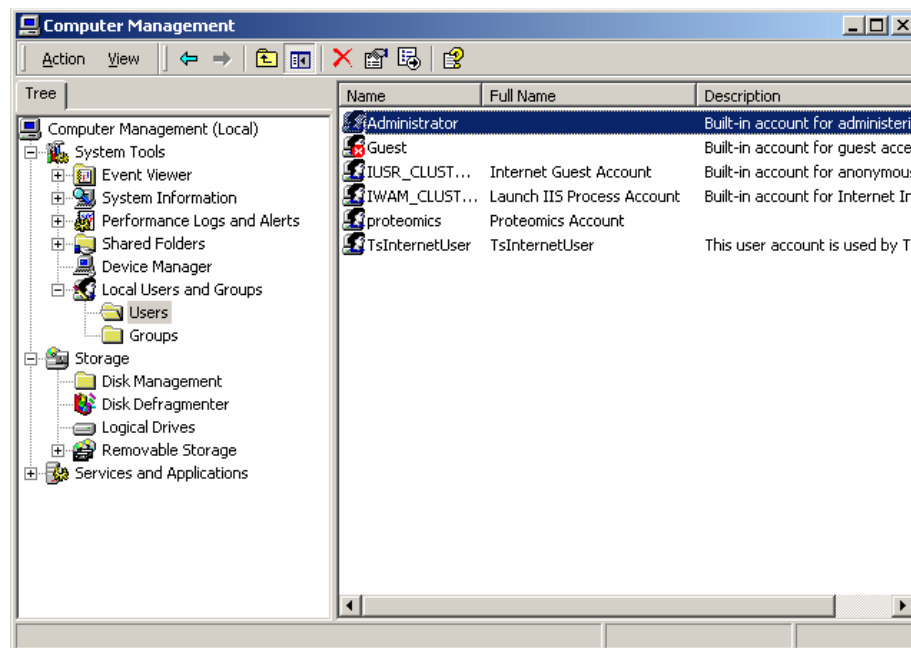
- b. Double-click **Administrative Tools** to open the Administrative Tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 5](#).
 - d. Format the drive using the NTFS file system.
 - e. In the right pane, right-click on the CD-ROM (D:) entry and choose **Change Drive Letter and Path** from the shortcut menu.
 - f. Select an open/unused drive letter for the CD-ROM. You can now format the unused portion of the hard drive as D:\.
10. After you format the D:\ drive, create a directory named *Database* on that drive. Use this directory to store the FASTA databases that are needed by SEQUEST.
 11. Configure the graphics card.

Figure 5. Computer Management tool



12. Create a user named *proteomics*.
 - a. Choose **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open the Administrative Tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 6](#).
 - d. In the left pane, double-click **Local Users and Groups**, and then select **Users**.

Figure 6. Computer Management tool



- e. Choose **Action > New User** and fill in the information in the dialog box. Enter “proteomics” as a user name. Click **Create**.
13. Add the new user proteomics to the Power Users and Administrators groups.
 - a. Right-click on **proteomics** and choose **Properties** from the shortcut menu to open the Proteomics Properties dialog box.
 - b. Click **Member Of** tab and click **Add** to open the Select Groups dialog box.
 - c. In the upper pane, click on the Power Users group to select it, and click **Add**.
 - d. Click **OK** to close the Select Groups dialog box.
 - e. Click **OK** to close the Proteomics Properties dialog box.
 - f. Close the Computer Management tool.
 14. Install the Thermo Fisher Scientific-supplied software for the master node, use the procedure described in “[Installing Master Node Software](#)” on page 16.

Installing Master Node Software

Install the following necessary software onto the master node as described in the procedures in this topic:

- [Installing Xcalibur, BioWorks, and the FASTA Databases](#)
- [Installing PVM](#)
- [Installing RSH](#)
- [Installing SEQUEST Cluster Software](#)

Installing Xcalibur, BioWorks, and the FASTA Databases

❖ To install Xcalibur, BioWorks and the FASTA databases on the master node

1. Insert the Xcalibur 1.4 SR1 or 2.0.6 CD into the computer and install Xcalibur onto the C: drive. Remove the Xcalibur CD.
2. Insert the BioWorks 3.3.1 SP1 CD into the computer and install Bioworks 3.3.1 SP1 onto the C: drive.

When you are prompted for the location of the FASTA database and SEQUEST results, if you created a D: drive on the system, then select a directory in that drive. For example:

```
D:\Database (Fasta Database Location)
D:\Sequest (SEQUEST Results)
```

3. Install the FASTA databases.
4. Remove the BioWorks 3.3.1 SP1 CD.

Note For more detailed instructions for installing BioWorks and the FASTA databases, refer to the *BioWorks Installation Guide*.

Installing PVM

❖ To install PVM on the master node


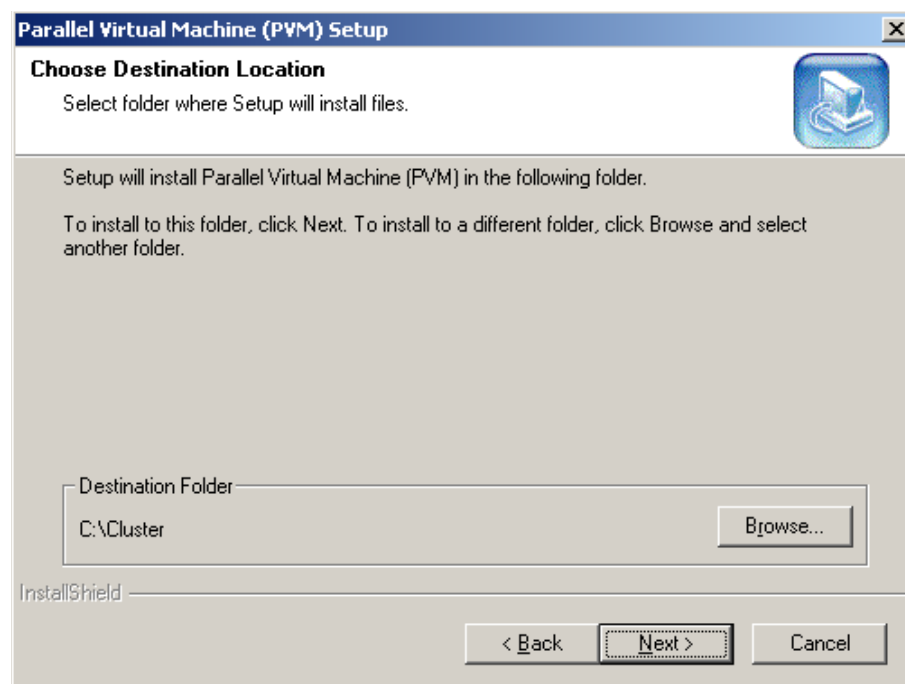
1. Insert the SEQUEST Cluster CD into the computer. This CD contains the SEQUEST Cluster, PVM, and RSH software.
2. In Microsoft Windows Explorer ( **Start > Programs > Accessories > Windows Explorer**) open the *PVM* folder on the CD.
3. Double-click **Setup.exe** to open the Parallel Virtual Machine Setup dialog box. See [Figure 7](#).
4. On the Choose Destination Location page, choose **C:\Cluster** and click **Next**. The installer creates the *C:\Cluster* folder and the correct sub-folders.
5. On the Start Copying Files page, click **Next** to begin the installation.

Figure 7. PVM Setup dialog box



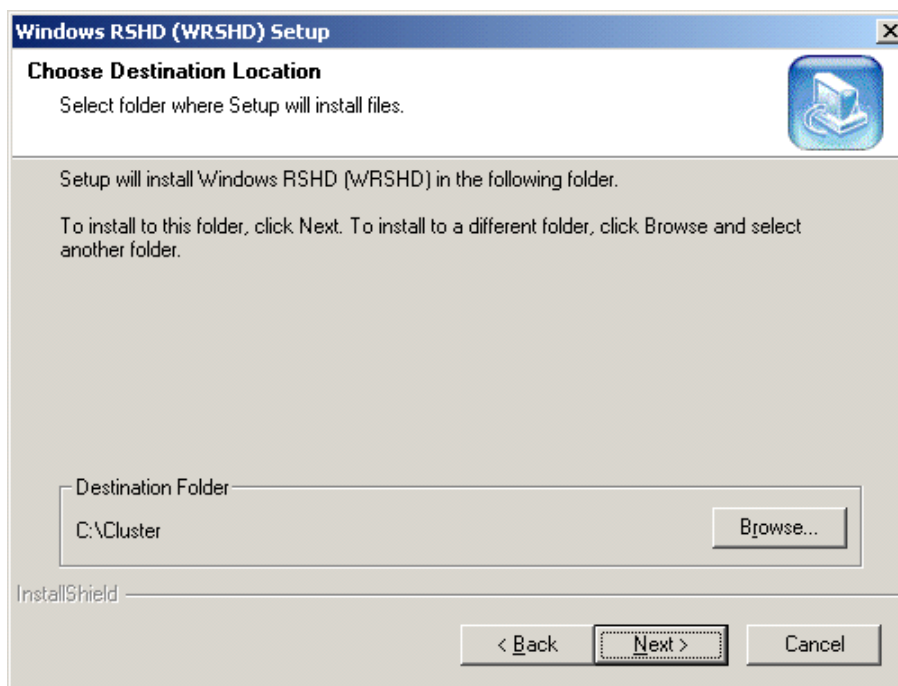
6. In Microsoft Windows Explorer, locate C:/Cluster/temp folder.
7. Select the folder and then, right-click to access the shortcut menu. Click **Properties** from the shortcut menu.
8. Click **Security** tab.
9. Set the Users Group to Full Control, select the **Allow** checkbox.
10. Click **OK** to accept the changes.

Installing RSH

❖ To install RSH on the master node

1. In Microsoft Windows Explorer, open the *RSH* folder on the SEQUEST Cluster CD.
2. Double-click **setup.exe** to open the RSH Installer. See [Figure 8](#).
3. On the Choose Destination Location page, select the **C:\Cluster** folder.
4. Click **Next** to begin the installation.
5. When the installation is complete, the Setup Complete page is displayed. You do not need to select either of the check boxes on this page. Click **Finish**.

Figure 8. Windows RSHD Setup dialog box



6. Open an MS-DOS Console window (**Start > Programs > Accessories > Command Prompt**).
7. Change to the directory `c:\cluster\wrshd` and type the command:

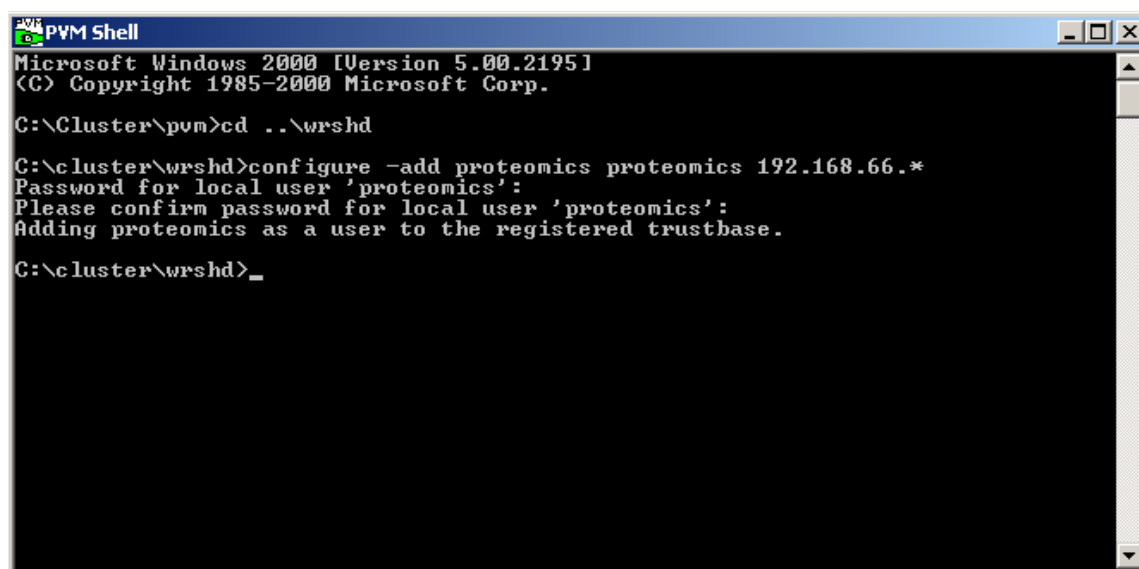
```
rshd.exe -install
```

This command installs the rshd daemon as a service.
8. Type the command:

```
configure -add proteomics proteomics 192.168.66.*
```

You are prompted for the password for the proteomics account.
9. Enter the password twice. See [Figure 9](#).
10. Close the MS-DOS console window.

Figure 9. Output from configure --add command



```

PVM Shell
Microsoft Windows 2000 [Version 5.00.2195]
(C) Copyright 1985-2000 Microsoft Corp.

C:\Cluster\pvm>cd ..\wrshd

C:\cluster\wrshd>configure --add proteomics proteomics 192.168.66.*
Password for local user 'proteomics':
Please confirm password for local user 'proteomics':
Adding proteomics as a user to the registered trustbase.

C:\cluster\wrshd>_

```

Installing SEQUEST Cluster Software

❖ To install SEQUEST Cluster software on the master node

1. In Microsoft Windows Explorer, open the *Cluster* folder on the SEQUEST Cluster CD.
2. Double-click **setup.exe** to open the SEQUEST Cluster Installer.
3. Accept all of the default settings.
4. Remove the SEQUEST Cluster CD.

The configuration of the master node is now complete.

5. Restart your computer.

You can now configure the slave nodes as described in the next sections:

- To configure the slave nodes with Microsoft Windows 2000 Professional, go to [“Setting Up the Slave Nodes: Microsoft Windows 2000 Professional”](#) on page 20.
- To configure the slave nodes with Microsoft Windows XP Professional, go to [“Setting Up the Slave Nodes: Microsoft Windows XP Professional”](#) on page 28.
- To configure the slave nodes with Red Hat Linux 8.0, go to [“Setting Up the Slave Nodes: Red Hat Linux \(Version 8.0\)”](#) on page 35.
- To configure the slave nodes with Red Hat Linux 9.0, go to [“Setting Up the Slave Nodes: Red Hat Linux \(Version 9.0\)”](#) on page 40.
- To configure the slave nodes with Red Hat Enterprise Linux 3.0 ES, go to [“Setting Up the Slave Nodes: Red Hat Enterprise Linux 3.0”](#) on page 45.

- To configure the slave nodes with Red Hat Enterprise Linux 4.0 ES, go to [“Setting Up the Slave Nodes: Red Hat Enterprise Linux 4.0”](#) on page 49.
- To configure the slave nodes with Fedora 3.0, go to [“Setting Up the Slave Nodes: Fedora 3.0”](#) on page 53.
- To configure the slave nodes with Fedora 4.0, go to [“Setting Up the Slave Nodes: Fedora 4.0”](#) on page 58.
- To configure the slave nodes with Fedora 5.0, go to [“Setting Up the Slave Nodes: Fedora 5.0”](#) on page 62.
- To configure the slave nodes with Fedora 6.0, go to [“Setting Up the Slave Nodes: Fedora 6.0”](#) on page 67.

Setting Up the Slave Nodes: Microsoft Windows 2000 Professional

This procedure assumes that the slave nodes have not been configured with the Microsoft Windows 2000 Professional operating system. If the slave nodes have already been configured, then you can skip the corresponding steps in the procedure below.

This section contains the following topics:

- [Installing and Configuring Microsoft Windows 2000 Professional](#)
- [Installing Slave Node Software](#)
- [Updating the Microsoft Windows Registry](#)

Note The following procedures must be repeated for **each** slave node in the cluster. Configure each slave node **identically** to the other nodes.

Installing and Configuring Microsoft Windows 2000 Professional

❖ To install and configure the Microsoft Windows 2000 Professional operating system

1. Install Microsoft Windows 2000 Professional on each slave computer.

More information about installing Microsoft Windows 2000 Professional can be found in Chapters 3 and 4 of the *Getting Started* manual that comes with Microsoft Windows 2000.

2. Install additional drivers for Microsoft Windows 2000 to recognize the hard disk(s) in your system, if required. If any special drivers are required, then consult the manual that comes with the computer hardware.

3. Define the disk partitions for the operating system. Create a small 4GB partition for the operating system and then commit the rest of the space on the hard disk to a separate partition that will store the fasta databases, dta and out files. You will set up this unused portion of the drive later.

4. Format the drive using the NTFS file system.

After you install the server software, the computer restarts and opens the Microsoft Windows 2000 Professional Setup program.

5. Use the Microsoft Windows 2000 Professional Setup program to set up the software.

During the setup, you are asked to enter the base operating system license. Enter the license that was provided with the Microsoft Windows 2000 software.

6. When you are asked for a name for this computer, name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.

When you are finished with the setup, the Microsoft Windows 2000 Professional Setup program restarts the computer. You can now configure the Microsoft Windows 2000 software.

7. When the Microsoft Windows 2000 Professional software restarts, the Network Identification Wizard dialog box appears. The dialog box prompts:

“Who can log on to this computer”.

Select the option titled **Microsoft Windows always assumes the following user has logged onto this computer** and enter the user account “Administrator” and the password for the administrator account. These steps activate the AutoLogon feature. Later, you will change the account used with the AutoLogon feature.

8. Configure the network card. The slave computers are connected to the secondary network card in the master computer through the Ethernet hub, and therefore they should have IP address in the same group as the secondary network card in the master computer.

Use a private address space number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Note When you are configuring the network, do not disable the Client for Microsoft Networks option. This option is required in order to share drives between individual computers running Microsoft Windows 2000.

9. Because the slave computers are not be a part of a domain, edit the C:\WINNT\system32\drivers\etc\hosts file and add the name and IP address for the master computer and each slave node in your cluster (see [Figure 10](#)).

Figure 10. Slave node hosts file

```
# Copyright (c) 1993-1999 Microsoft Corp.

#

# This is a sample HOSTS file used by Microsoft TCP/IP for
# Microsoft Windows.

#

# This file contains the mappings of IP addresses to host names. Each
# entry should be kept on an individual line. The IP address should
# be placed in the first column followed by the corresponding host
# name.

# The IP address and the host name should be separated by at least
# one space.

#

# Additionally, comments (such as these) may be inserted on
# individual lines or following the machine name denoted by
# a '#' symbol.

#
127.0.0.1          localhost
192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```


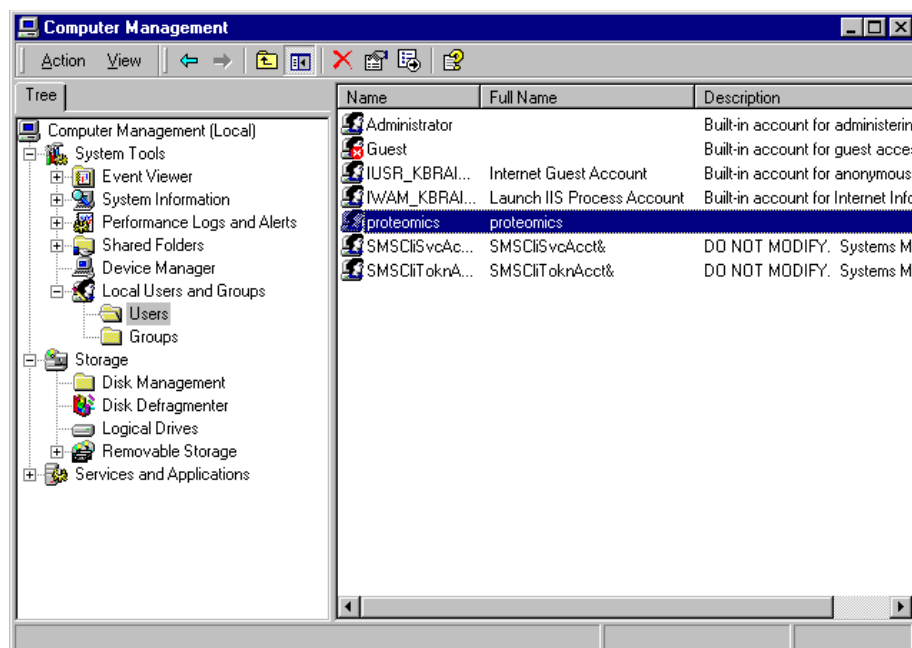
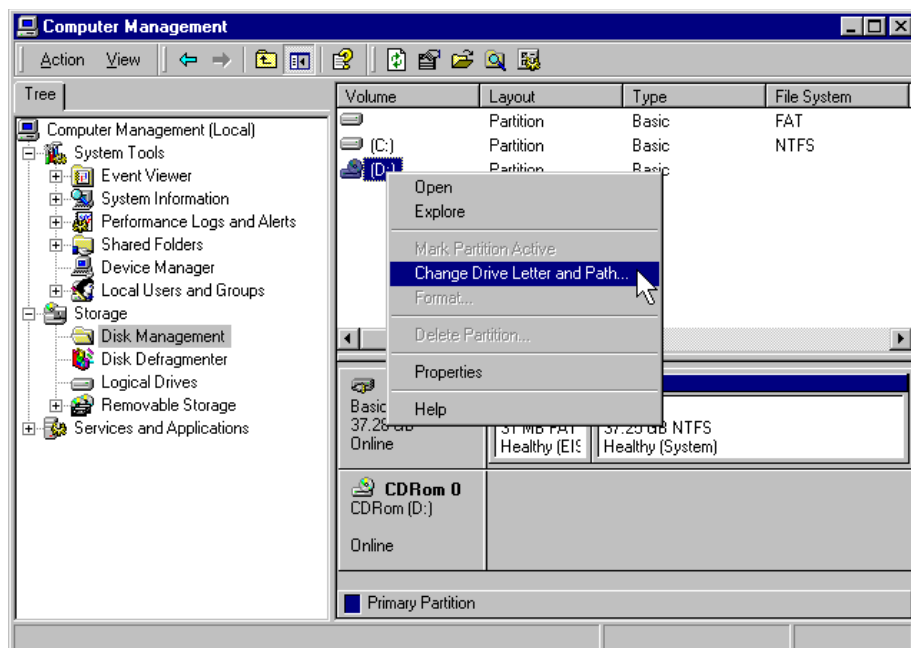
10. Install Service Pack 4 for Microsoft Windows 2000 onto the slave computer. When the installation of Service Pack 4 is finished, the computer will restart.
11. Log in as **administrator**.
12. Configure the graphics card.
13. Create a user named proteomics:
 - a. Choose  **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open administrative tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 11](#).
 - d. In the left pane, double-click **Local Users and Groups**, and then select **Users**.

Figure 11. Computer Management tool

- e. Choose **Action > New User** and fill in the information in the dialog box. Enter “proteomics” as a user name. Ensure that you select the **Password Never Expires** check box. Click **Create**.
14. Add the new user proteomics to the Power Users group.
 - a. Click **proteomics**.
 - b. Choose **Action > Properties** to open the Proteomics Properties dialog box.
 - c. Click **Member Of** tab and click **Add** to open the Select Groups dialog box.
 - d. In the upper pane, click on the Power Users group to select it, and then click **Add**.
 - e. Click **OK** to close the Select Groups dialog box.
 - f. Click **OK** to close the Proteomics Properties dialog box.
 - g. Close the Computer Management tool.
 15. Format the unused portion of the hard disk and set the drive letter for this partition to D:\ as follows:
 - a. Choose **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open administrative tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 12](#).

Figure 12. Computer Management tool



- d. Format the drive using the NTFS file system.
 - e. In the right pane, right-click on the CD-ROM (D:) entry and choose **Change Drive Letter and Path** from the shortcut menu.
 - f. Select an open/unused drive letter for the CD-ROM. You can now format the unused portion of the hard drive as D:\.
16. After you format the D:\ drive, create a directory named *Database* on that drive. You will use this directory to store the FASTA databases that are needed by SEQUEST.
 17. Create a directory on the C:\ drive named *Cluster*. This will be used when loading the PVM software. Share the C:\Cluster directory as *Cluster*.
 18. The *Database* and *Cluster* directories must be shared for the Cluster Manager program to work correctly. When you share these directories, ensure that the proteomics account has permission to access these share points.

Installing Slave Node Software

Install the following necessary software onto the slave node as described in the procedures in this topic:

- [Installing PVM](#)
- [Installing RSH](#)

Note The Xcalibur, Bioworks Browser, SEQUEST Cluster, and IIS web server software programs do not need to be installed on the slave nodes.

Installing PVM

❖ To install PVM on each of the slave nodes


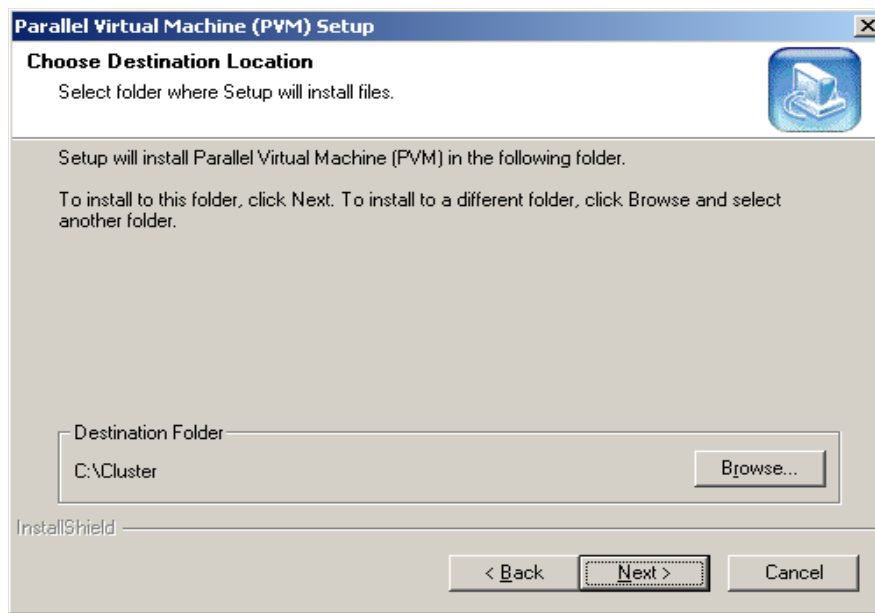
1. Insert the SEQUEST Cluster CD into the computer. This CD contains the SEQUEST Cluster, PVM, and RSH software.
2. In Microsoft Windows Explorer ( **Start > Programs > Accessories > Windows Explorer**) open the *PVM* folder on the CD.
3. Double-click **Setup.exe** to open the Parallel Virtual Machine Setup dialog box. See [Figure 13](#).
4. On the Choose Destination Location page, choose **C:\Cluster** and click **Next**. The installer creates the *C:\Cluster* folder and the correct sub-folders.
5. On the Start Copying Files page, click **Next** to begin the installation.

Figure 13. PVM Setup dialog box

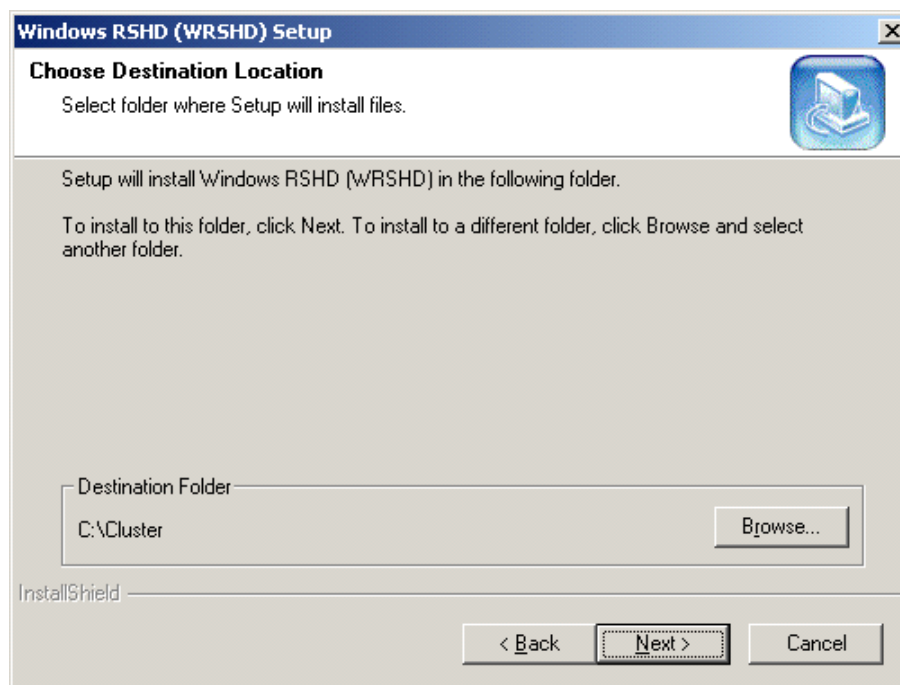


Installing RSH

❖ To install RSH on each of the slave nodes

1. In Microsoft Windows Explorer, select the *RSH* folder on the SEQUEST Cluster CD.
2. Double-click **setup.exe** to open the RSH Installer. See [Figure 14](#).
3. On the Choose Destination Location page, select the **C:\Cluster** folder.
4. Click **Next** to begin the installation.
5. When the installation is complete, the Setup Complete page is displayed. Do not select either of the check boxes on this page. Click **Finish**.

Figure 14. Windows RSHD Setup Dialog Box.



6. Open an MS-DOS Console window (**Start > Programs > Accessories > Command Prompt**).
7. Change to the directory *C:\Cluster\wrshd* and type the command:

```
rszd.exe -install
```

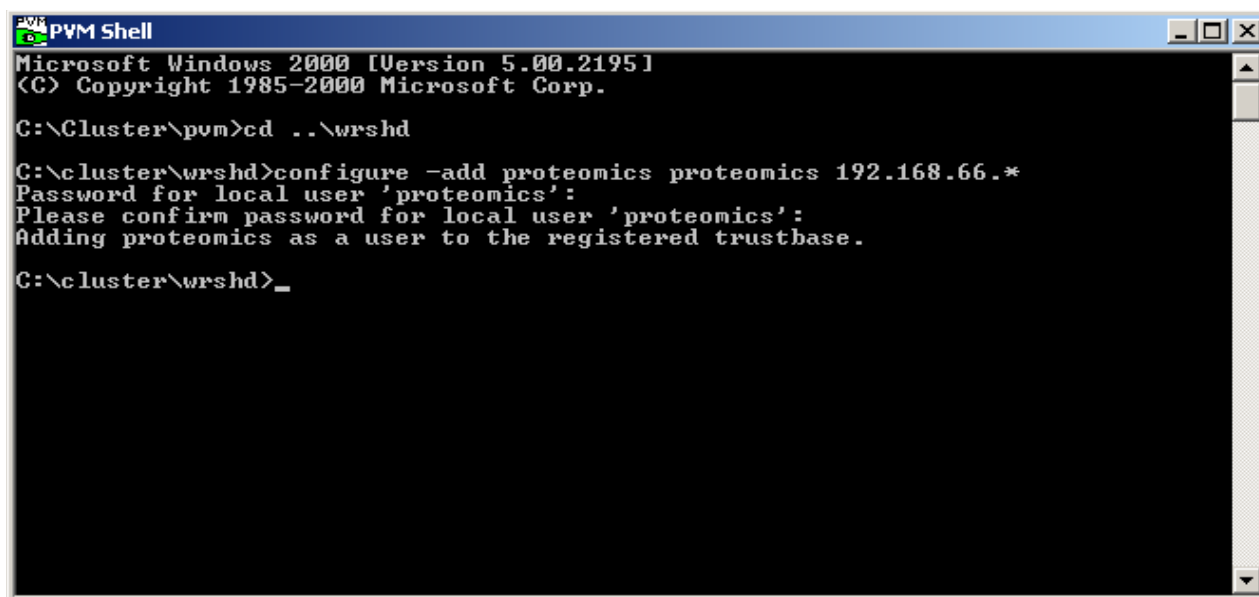
This command installs the rszd daemon as a service.
8. Type the command:

```
configure -add proteomics proteomics 192.168.66.*
```

You are prompted for the password for the proteomics account.
9. Enter the password twice. See [Figure 15](#).

10. Close the MS-DOS console window.

Figure 15. Output from `configure --add` command



```

PVM Shell
Microsoft Windows 2000 [Version 5.00.2195]
(C) Copyright 1985-2000 Microsoft Corp.

C:\Cluster\pvm>cd ../wrshd

C:\cluster\wrshd>configure --add proteomics proteomics 192.168.66.*
Password for local user 'proteomics':
Please confirm password for local user 'proteomics':
Adding proteomics as a user to the registered trustbase.

C:\cluster\wrshd>_

```

Updating the Microsoft Windows Registry

The user name that is used with the AutoLogin feature of Microsoft Windows needs to be updated to the user *proteomics*.

❖ To update the AutoLogin user name

1. Choose **Start > Run** to open the Run dialog box.
2. Type “regedit” and then click **OK** to open the Registry Editor.
3. Browse to the folder:

HKEY_LOCAL_MACHINE\SOFTWARE\Microsoft\Windows NT\CurrentVersion\Winlogon

4. Choose **Edit > New String Value** to add a new string value.
5. Name the new string value **DefaultUserName**.
6. Click **DefaultUserName** to select it and choose **Edit > Modify**.
7. Type the user name that you want to use to automatically log on. For example: *proteomics*.
8. Add a new string value and name it **DefaultPassword**.
9. Modify **DefaultPassword**: set it to the password for the user entered in step 7. For example, the password for *proteomics*.
10. Add a new string value and name it **AutoAdminLogon**.

11. Modify AutoAdminLogon: set it to 1 to enable auto logon.
12. Close the Registry Editor.

The configuration of the slave node is now complete.

13. Restart the computer.

When all of the slave nodes have been configured, go to [“Final Configuration”](#) on [page 72](#).

Setting Up the Slave Nodes: Microsoft Windows XP Professional

This procedure assumes that the slave nodes have not been configured with the Microsoft Windows XP Professional operating system. If the slave nodes have already been configured then you can skip the corresponding steps in the procedure below.

This procedure assumes that you are familiar with installing Microsoft Windows XP Professional.

This section contains the following topics:

- [Installing and Configuring Microsoft Windows XP Professional](#)
- [Installing Slave Node Software](#)

Note The following procedures must be repeated for **each** slave node in the cluster. Configure each slave node **identically** to the other nodes.

Installing and Configuring Microsoft Windows XP Professional

❖ To install and configure the Microsoft Windows XP Professional operating system

1. Install Microsoft Windows XP Professional on each slave computer.

More information about installing Microsoft Windows XP Professional can be found in Chapters 3 and 4 of the *Getting Started* manual that comes with Microsoft Windows XP.

2. Install additional drivers for Microsoft Windows XP to recognize the hard disk(s) in your system, if required. If any special drivers are required, then consult the manual that comes with the computer hardware.
3. Define the disk partitions for the operating system. Create a small 6GB partition for the operating system and then commit the rest of the space on the hard disk to a separate partition that will store the FASTA databases, and DTA and OUT or SRF files. You will set up this unused portion of the drive later.
4. Format the drive using the NTFS file system.

After you install the server software, the computer restarts and opens the Microsoft Windows XP Professional Setup program.

5. Use the Microsoft Windows XP Professional Setup program to set up the software.

During the setup, you are asked to enter the base operating system license. Enter the license that was provided with the Microsoft Windows XP software.

6. When you are asked for a name for this computer, name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.

When you are finished the setup, the Microsoft Windows XP Professional Setup program restarts the computer. You can now proceed configuring the Microsoft Windows XP software.

7. When the Microsoft Windows XP Professional software restarts, the Network Identification Wizard dialog box appears. The Network Identification Wizard dialog box prompts:

“Who can log on to this computer”.

Select the option titled **Microsoft Windows always assumes the following user has logged onto this computer** and enter the user account Administrator and the password for the administrator account. These steps activate the AutoLogon feature. Later the account used with the AutoLogon feature will be changed.

8. Configure the network card. The slave computers are connected to the secondary network card in the master computer through the Ethernet hub, and therefore they should have IP address in the same group as the secondary network card in the master computer.

Use a private address space number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Note When you are configuring the network, do not disable the Client for Microsoft Networks option. This option is required in order to share drives between individual computers running Microsoft Windows XP.

9. Because the slave computers are not part of a domain, edit the C:\WINNT\system32\drivers\etc\hosts file and add the name and IP address for the master computer and each slave node in your cluster (see [Figure 16](#)).

Figure 16. Slave node hosts file

```
# Copyright (c) 1993-1999 Microsoft Corp.

#

# This is a sample HOSTS file used by Microsoft TCP/IP for
# Microsoft Windows.

#

# This file contains the mappings of IP addresses to host names. Each
# entry should be kept on an individual line. The IP address should
# be placed in the first column followed by the corresponding host
# name.

# The IP address and the host name should be separated by at least one
# space.

#

# Additionally, comments (such as these) may be inserted on
# individual lines or following the machine name denoted by
# a '#' symbol.

#

127.0.0.1          localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```


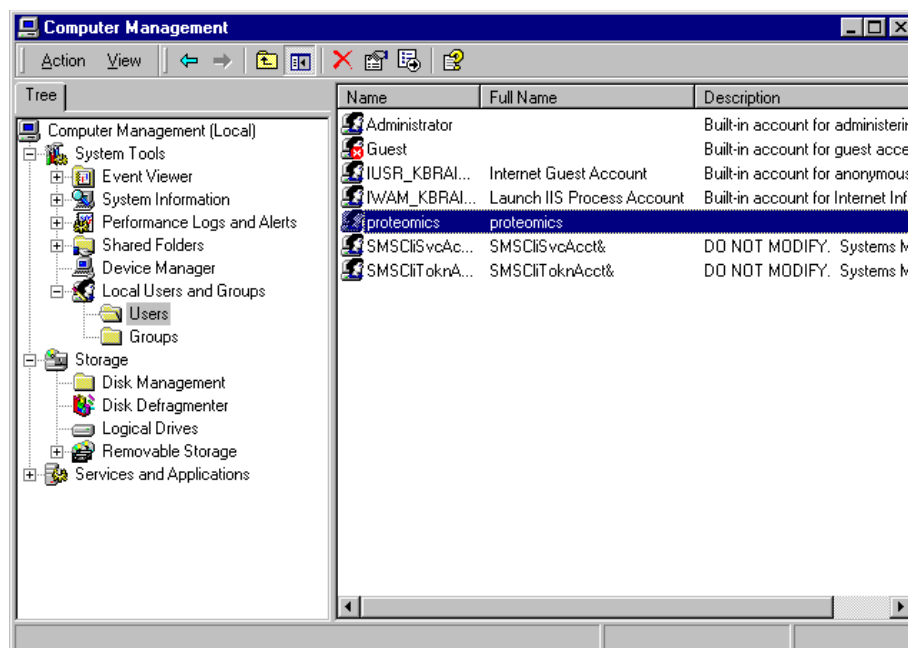
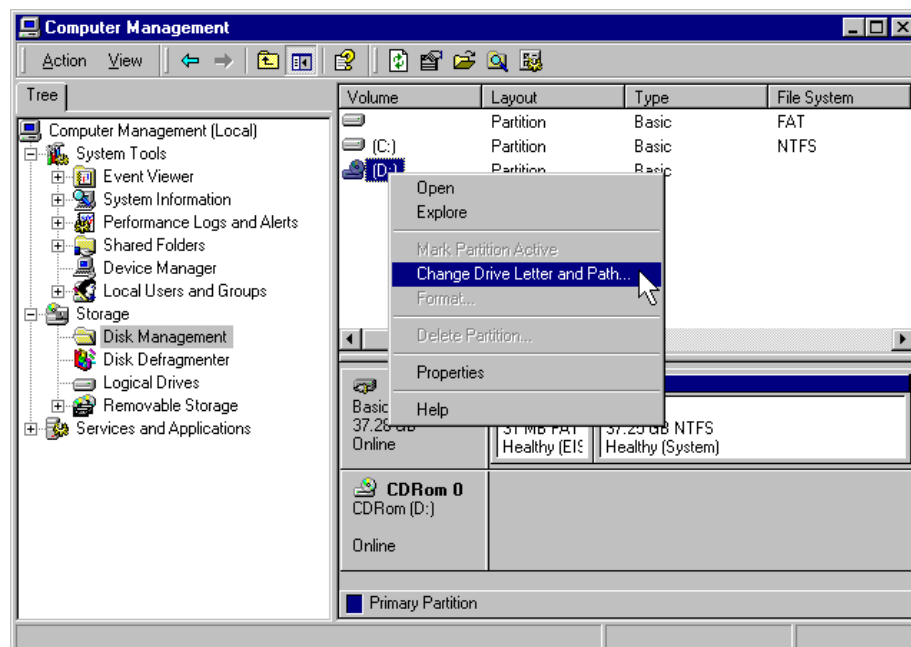
10. Install Service Pack 2 for Microsoft Windows XP onto the slave computer. When the installation of Service Pack 2 is finished, the computer will restart.
11. Log in as **administrator**.
12. Configure the graphics card.
13. Create a user named proteomics:
 - a. Choose  **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open administrative tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 17](#).
 - d. In the left pane, double-click **Local Users and Groups**, and then select **Users**.

Figure 17. Computer Management tool



- e. Choose **Action > New User** and fill in the information in the dialog box. Enter “proteomics” as a user name. Ensure that you select the **Password Never Expires** check box. Click **Create**.
14. Add this proteomics user to the Power Users group:
- a. Click **proteomics**.
 - b. Choose **Actions > Properties** to open the Proteomics Properties dialog box.
 - c. Click **Member Of** tab and click **Add** to open the Select Groups dialog box.
 - d. In the upper pane, click the **Power Users** group to select it, and then click **Add**.
 - e. Click **OK** to close the Select Groups dialog box.
 - f. Click **OK** to close the Proteomics Properties dialog box.
 - g. Close the Computer Management tool.
15. Format the unused portion of the hard disk and set the drive letter for this partition to D:\ as follows:
- a. Choose **Start > Settings > Control Panel** to open the Control Panel.
 - b. Double-click **Administrative Tools** to open administrative tools.
 - c. Double-click **Computer Management** to open the Computer Management tool. See [Figure 18](#).

Figure 18. Computer Management tool



- d. Format the drive using the NTFS file system.
 - e. In the right pane, right-click on the CD-ROM (D:) entry and choose **Change Drive Letter and Path** from the shortcut menu.
 - f. Select an open/unused drive letter for the CD-ROM. You can now format the unused portion of the hard drive as D:\.
16. After you format the D:\ drive, create a directory named *Database* on that drive. You will use this directory to store the FASTA databases that are needed by SEQUEST.
 17. Create a directory on the C:\ drive named *Cluster*. This will be used when loading the PVM software. Share the C:\Cluster directory as *Cluster*.
 18. The *Database* and *Cluster* directories must be shared for the Cluster Manager program to work correctly. When you share these directories, ensure that the proteomics account has permission to access these share points.

Installing Slave Node Software

Install the following necessary software onto the slave node as described in the procedures in this topic:

- [Installing PVM](#)
- [Installing RSH](#)

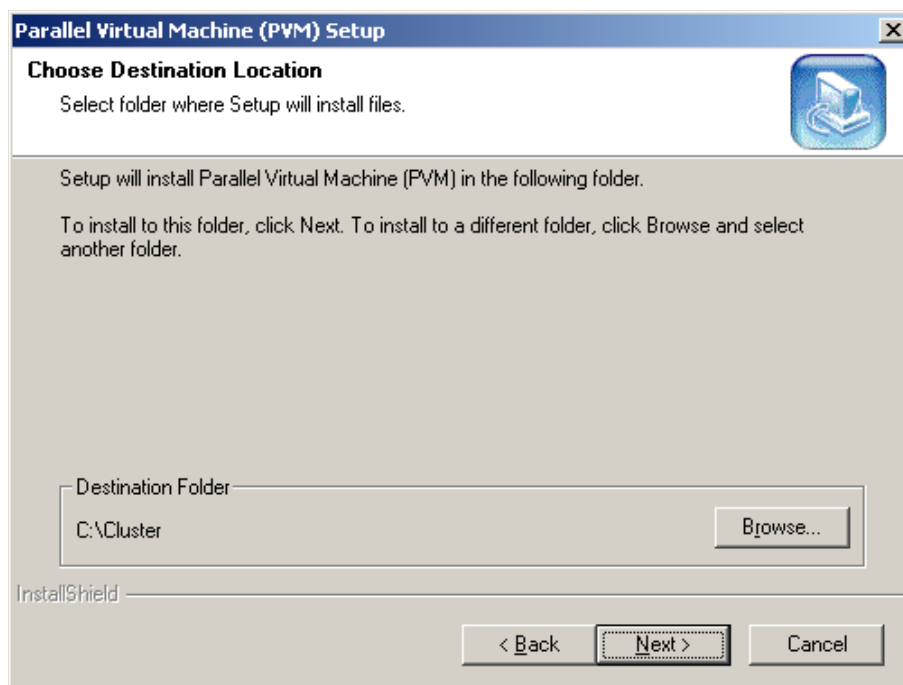
Note The Xcalibur, Bioworks Browser, SEQUEST Cluster, and IIS web server software programs do not need to be installed on the slave nodes.

Installing PVM

❖ To install PVM on each of the slave nodes

1. Insert the SEQUEST Cluster CD into the computer. This CD contains the SEQUEST Cluster, PVM, and RSH software.
2. In Microsoft Windows Explorer (Start > Programs > Accessories > Windows Explorer) open the *PVM* folder on the CD.
3. Double-click **Setup.exe** to open the Parallel Virtual Machine Setup dialog box. See Figure 19.
4. On the Choose Destination Location page, choose *C:\Cluster* and click **Next**. The installer creates the *C:\Cluster* folder and the correct sub-folders.
5. On the Start Copying Files page, click **Next** to begin the installation.

Figure 19. PVM Setup dialog box



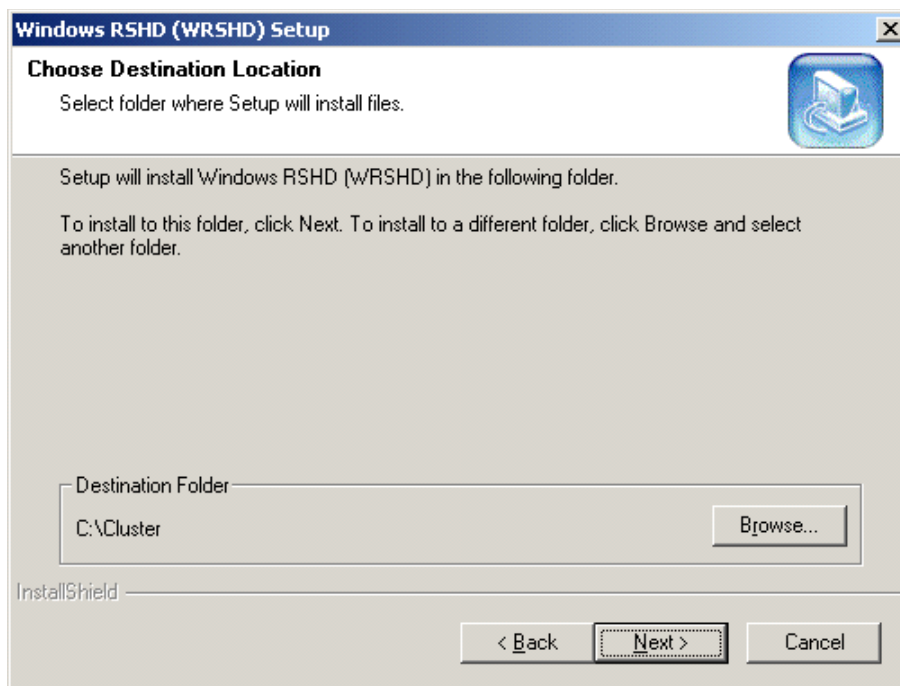
Installing RSH

❖ To install RSH on each of the slave nodes

1. In Microsoft Windows Explorer, select the *RSH* folder on the SEQUEST Cluster CD.
2. Double-click **setup.exe** to open the RSH Installer. See Figure 20.
3. On the Choose Destination Location page, select the **C:\Cluster** folder.
4. Click **Next** to begin the installation.

- When the installation is complete, the Setup Complete page is displayed. Do not select either of the check boxes on this page. Click **Finish**.

Figure 20. Windows RSHD Setup Dialog Box.



- Open an MS-DOS Console window (**Start > Programs > Accessories > Command Prompt**).
- Change to the directory C:\Cluster\wrshd and type the command:

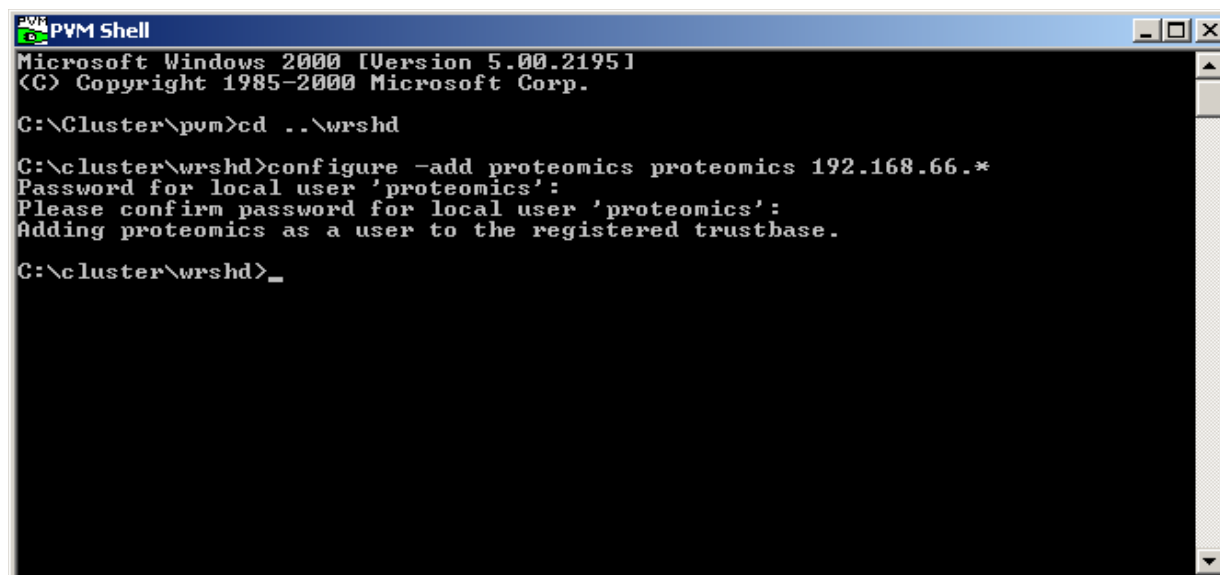
```
rshd.exe -install
```

This command installs the rshd daemon as a service.
- Type the command:

```
configure -add proteomics proteomics 192.168.66.*
```

You are prompted for the password for the proteomics account.
- Enter the password twice. See [Figure 21](#).
- Close the MS-DOS console window.

Figure 21. Output from configure --add command



```
PVM Shell
Microsoft Windows 2000 [Version 5.00.2195]
(C) Copyright 1985-2000 Microsoft Corp.

C:\Cluster\pvm>cd ../wrshd

C:\cluster\wrshd>configure -add proteomics proteomics 192.168.66.*
Password for local user 'proteomics':
Please confirm password for local user 'proteomics':
Adding proteomics as a user to the registered trustbase.

C:\cluster\wrshd>_
```

Setting Up the Slave Nodes: Red Hat Linux (Version 8.0)

This procedure assumes that you are familiar with installing Red Hat Linux 8.0. Additional information regarding Red Hat Linux can be found in *Red Hat Linux 8 UNLEASHED* by Bill Ball (ISBN #067232458X).

This section contains the following topics:

- [Installing Red Hat Linux 8.0](#)
- [Configuring Red Hat Linux 8.0](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node **identically** to the other nodes.

Installing Red Hat Linux 8.0

This procedure assumes that the Red Hat Linux operating system has not already been installed on the slave nodes. If Red Hat Linux has already been installed then skip to [“Configuring Red Hat Linux 8.0”](#) on [page 37](#).

❖ To install Red Hat Linux

1. Insert the Red Hat Linux 8.0 Installation 1 of 3 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing ENTER or by waiting for a few seconds.

3. In the Welcome to Red Hat Linux 8.0 page, click **Next**.
4. In the Language Selection page, select the language you want to use and click **Next**.
5. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
6. In the Mouse Configuration page, select what type of mouse is connected to the computer and click **Next**.
7. In the Installation Type page, select **Custom** from the list and click **Next**.
8. In the Disk Partitioning Setup page, unless you are familiar with installing Red Hat Linux, select the **Automatically partition** option and click **Next**.
9. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option and click **Next**.
10. In the Partitioning page, click **Next**.
11. In the Boot Loader Configuration page, select the **GRUB boot loader** and click **Next**.
12. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Deselect the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up. The Netmask is automatically entered.
 - c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.

Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
13. In the Firewall Configuration page, select the **No Firewall** option and then click **Next**.
14. In the Additional Language Support page, click **Next**.
15. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.

16. In the Account Configuration page, enter the root password for the system and click **Next**.

The Account Configuration page allows a user account to be created but it limits the length of the account to 8 characters. Therefore, the name *proteomics* cannot be used. Because of this limitation, do not create an account at this point.

17. In the Authentication Configuration page, click **Next**.

The installation program takes a few moments to build the list of packages that make up Red Hat Linux.

18. In the Package Group Selection page, deselect the options:

- X Window, GNOME Desktop Environment
- Graphical Internet, Text Based Internet
Office/Productivity, Sound and Video
- Graphics
- Printing Support

19. Enable the **FTP Server** option.

20. Enable the **Scientific and Engineering** option. Then display the sub-packages dialog box. Enable the **PVM** option, and de-select all of the other options.

21. Enable the **Networking Servers** option, and then display the sub-packages dialog box. Enable the **RSH Server** option.

22. Click **Next** to install the RedHat Linux software.

23. Skip the Boot Diskette Creation and go to the Congratulations screen. Follow the instructions to complete the installation.

Configuring Red Hat Linux 8.0

❖ To configure Red Hat Linux

1. Log into your computer as *root*.
2. Run the service configuration program:

```
/usr/sbin/ntsysv
```
3. Scroll down the list and make sure the *rexec*, *rsh*, and *vsftpd* options are active. You can toggle an option by pressing **SPACE** on the keyboard.
4. Disable the sendmail option.
5. Press **TAB** to select the **OK** option and press **ENTER**.

6. While the previous operation enabled the FTP software, it will not allow inbound FTP until the */etc/xinetd.d/vsftpd* file is modified. Open the existing file with this command:

```
vi /etc/xinetd.d/vsftpd
```

7. Change *yes* to *no* in the line that begins with *disable* (see [Figure 22](#)).

Figure 22. vsftpd configuration file for Red Hat Linux 8.0

```
# default: off

# description: The vsftpd FTP server serves FTP connections. It \
# uses normal, unencrypted usernames and passwords for \
# authentication.

service ftp
{
    disable          = no
    socket_type      = stream
    wait             = no
    user             = root
    server           = /usr/sbin/vsftpd
    nice             = 10
}
```

8. Save the file.
9. Create the proteomics user account with this command:

```
adduser proteomics
```

10. Set the password for this account with this command:

```
passwd proteomics
```

11. Create a directory for the FASTA databases and change the permissions on that directory:

```
mkdir /usr/database
chmod 777 /usr/database
```

12. Change the permissions of the */usr/share/pvm3/bin/LINUX* directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUX
```

13. Create a configuration file (sequest.ini) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```

14. Add the following line to the configuration file and then save the file:

```
DatabasePath: /usr/database
```

15. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

16. Add the following lines to the file and then save the file:

```
Master_computer_name proteomics
```


Where *Master_computer_name* is replaced with the name you used for the master computer.

17. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

18. Add the following line to the file and then save the file:

```
+ proteomics
```

19. Update the permissions for the *.rhosts* file:

```
chmod 600 /.rhosts
```

20. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 23](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 23. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

21. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 24](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 24. *.bashrc* file for Red Hat Linux 8.0

```
# .bashrc

# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUX
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

22. Save the changes to the file. The slave node is now configured.

23. Restart the computer with this command:

```
shutdown -r now
```

24. Repeat steps 1 through 23 for each of the remaining nodes.

When all of the slave nodes have been configured, go to [“Final Configuration”](#) on [page 72](#).

Setting Up the Slave Nodes: Red Hat Linux (Version 9.0)

This procedure assumes that you are familiar with installing Red Hat Linux 9.0. Additional information regarding Red Hat Linux can be found in *Red Hat Linux 9 UNLEASHED* by Bill Ball (ISBN #0672325888).

This section contains the following topics:

- [Installing Red Hat Linux 9.0](#)
- [Configuring Red Hat Linux 9.0](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node identically to the other nodes.

Installing Red Hat Linux 9.0

This procedure assumes that the Red Hat Linux operating system has not already been installed on the slave nodes. If Red Hat Linux has already been installed then skip to [“Configuring Red Hat Linux 9.0”](#) on [page 42](#).

❖ To install Red Hat Linux

1. Insert the Red Hat Linux 9.0 Installation 1 of 3 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing **ENTER** or by waiting for a few seconds.
3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Red Hat Linux 9.0 page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Mouse Configuration page, select what type of mouse is connected to the computer and click **Next**.
8. In the Installation Type page, select **Custom** from the list and click **Next**.

9. In the Disk Partitioning Setup page, unless you are familiar with installing Red Hat Linux, select the **Automatically partition** option and click **Next**.
10. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option and click **Next**.
11. In the Partitioning page, click **Next**.
12. In the Boot Loader Configuration page, select the **GRUB boot loader** and click **Next**.
13. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Deselect the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up. The Netmask will automatically be entered.
 - c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.

Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
14. In the Firewall Configuration page, select the **No Firewall** option and then click **Next**.
15. In the Additional Language Support page, click **Next**.
16. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
17. In the Account Configuration page, enter the root password for the system and click **Next**.

The Account Configuration page allows a user account to be created but it limits the length of the account to 8 characters. Therefore, the name *proteomics* cannot be used. Because of this limitation, do not create an account at this point.
18. In the Authentication Configuration page, click **Next**.

The installation program will take a few moments to build the list of packages that make up Red Hat Linux.

19. In the Package Group Selection page, deselect the options:
 - X Window, GNOME Desktop Environment
 - Graphical Internet, Text Based Internet Office/Productivity, Sound and Video
 - Graphics
 - Printing Support
20. Enable the **FTP Server** option.
21. Enable the **Scientific and Engineering** option. Then display the sub-packages dialog box. Enable the **PVM** option and deselect all of the other options
22. Enable the **Networking Servers** option and then display the sub-packages dialog. Enable the **RSH Server** option.
23. Click **Next** to install the Red Hat Linux software.
24. Skip the Boot Diskette Creation and go to the Congratulations screen. Follow the instructions to complete the installation.

Configuring Red Hat Linux 9.0

❖ To configure Red Hat Linux

1. Log into your computer as *root*.
2. Run the service configuration program:

```
/usr/sbin/ntsysv
```
3. Scroll down the list and make sure the *rexec*, *rsh*, and *vsftpd* options are active. You can toggle an option by pressing **SPACE** on the keyboard.
4. Disable the sendmail option.
5. Press **TAB** to select the **OK** option and press **ENTER**.
6. While the previous operation enabled the FTP software, it will not allow inbound FTP until the necessary file is modified. Open the */etc/vsftpd/vsftpd.conf* file, type:

```
vi /etc/vsftpd/vsftpd.conf
```
7. In the *vsftpd.conf* file, change *yes* to *no* in the line that begins with *anonymous enable* (see [Figure 25](#)).

Figure 25. vsftpd configuration file for Red Hat Linux 9.0

```
# default: off
# description: The vsftpd FTP server serves FTP connections. It \
# uses normal, unencrypted usernames and passwords for \
# authentication.
service ftp

{
    anonymous_enable = no
    socket_type      = stream
    wait            = no
    user            = root
    server          = /usr/sbin/vsftpd
    nice            = 10
}
```

8. Save the file.

9. Create the proteomics user account with this command:

```
adduser proteomics
```

10. Set the password for this account with this command:

```
passwd proteomics
```

11. Create a directory for the FASTA databases and change the permissions on that directory:

```
mkdir /usr/database
chmod 777 /usr/database
```

12. Change the permissions of the */usr/share/pvm3/bin/LINUXI386* directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUXI386
```

13. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```

14. Add the following line to the configuration file and then save the file:

```
DatabasePath: /usr/database
```

15. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

16. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.

17. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

2 Setting Up the SEQUEST Cluster

Setting Up the Slave Nodes: Red Hat Linux (Version 9.0)

18. Add the following line to the file and then save the file:

```
+ proteomics
```

19. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

20. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 26](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 26. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

21. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 27](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 27. *.bashrc* file for Red Hat Linux 9.0

```
# .bashrc
# User specific aliases and functions

# Source global definitions
if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUXI386
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

22. Save the changes to the file. The slave node is now configured.

23. Restart the computer with this command:

```
shutdown -r now
```

24. Repeat steps 1 through 23 for each of the remaining nodes.

When all of the slave nodes have been configured, go to “[Final Configuration](#)” on [page 72](#).

Setting Up the Slave Nodes: Red Hat Enterprise Linux 3.0

This procedure assumes that you are familiar with installing Red Hat Enterprise Linux 3.0 ES. This section contains the following topics:

- [Installing Red Hat Enterprise Linux 3.0 ES](#)
- [Configuring Red Hat Enterprise Linux 3.0 ES](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node identically to the other nodes.

Installing Red Hat Enterprise Linux 3.0 ES

This procedure assumes that the Red Hat Enterprise Linux operating system has not already been installed on the slave nodes. If Red Hat Enterprise Linux has already been installed then skip to “[Configuring Red Hat Enterprise Linux 3.0 ES](#)” on [page 47](#).

❖ To install Red Hat Enterprise Linux

1. Insert the Red Hat Enterprise Linux Installation 1 of 4 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing **ENTER** or by waiting for a few seconds.
3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Red Hat Enterprise Linux page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Mouse Configuration page, select what type of mouse is connected to the computer and click **Next**.
8. In the Disk Partitioning Setup page, unless you are familiar with installing Red Hat Enterprise Linux, select the **Automatically partition** option and click **Next**.
9. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option and click **Next**.
10. In the Partitioning page, click **Next**.
11. In the Boot Loader Configuration page, select the **GRUB boot loader** and click **Next**.
12. In the Network Configuration page, set up the network configuration parameters:

- a. Click **Edit** in the Network Devices box.
- b. Deselect the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Enter the netmask if it is not entered automatically.

- c. Click **OK** to save the changes.
- d. Select the **Manually Set the Hostname** option and enter the hostname.
Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
- e. Enter a gateway address for this network and then click **Next** to save these changes.
- f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.

13. In the Firewall Configuration page, select the **No Firewall** option and then click **Next**.
14. In the Additional Language Support page, click **Next**.
15. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
16. In the Set Root Password page, enter the root password for the system and click **Next**.
17. In the Package Defaults page, select **Customize the set of packages to be installed** and click **Next**.
18. Make the following selections and deselections:
 - a. In the Package Group Selection page, deselect the following options if they are selected:
 - Desktops - X Window, GNOME Desktop Environment
 - Applications - Editors, Graphical Internet, Text Based Internet, Office/Productivity, Graphics, Games and Entertainment
 - Servers - Server Configuration Tools, Web Server, Windows File Server
 - Development - Development Tools
 - System - Printing Support, Administration Tools
 - b. Select the following options in the Servers section:

- Applications - Scientific and Engineering Programs
 - Servers - FTP Server, Legacy Network Server.
- c. After selecting the **Legacy Network Server** option, click **Details** and select **rsh-server**.
 - d. After selecting the **Scientific and Engineering Programs** option, click **Details** and select **PVM**. Deselect all other options.
 - e. Click **Next**.
19. Click **Next** to install the Red Hat Linux software.
 20. In the Congratulations screen, follow the instructions to finish the installation.

Configuring Red Hat Enterprise Linux 3.0 ES

❖ To configure Red Hat Enterprise Linux

1. Log into your computer as *root*.
2. Run the service configuration program:

```
/usr/sbin/ntsysv
```
3. Scroll down the list and make sure the *rexec*, *rsh*, and *vsftpd* options are active. You can toggle an option by pressing **SPACE** on the keyboard.
4. Disable the sendmail option.
5. Press **TAB** to select the **OK** option and press **ENTER**.
6. Create the proteomics user account use this command:

```
adduser proteomics
```
7. Set the password for this account use this command:

```
passwd proteomics
```
8. Create a directory for the fasta databases and change the permissions on that directory:

```
mkdir /usr/database  
chmod 777 /usr/database
```
9. Change the permissions of the */usr/share/pvm3/bin/LINUXI386* directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUXI386
```
10. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```
11. Add the following line to the configuration file and then save the file:

```
DatabasePath: /usr/database
```

12. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

13. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.

14. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

15. Add the following line to the file and then save the file:

```
+ proteomics
```

16. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

17. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 28](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 28. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

18. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 29](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 29. `.bashrc` file for Red Hat Enterprise Linux 3.0 ES

```
# .bashrc
# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUXI386
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

19. Save the changes to the file. The slave node is now configured.

20. Restart the computer with this command:

```
shutdown -r now
```

21. Repeat steps 1 through 20 for each of the remaining nodes.

When all of the slave nodes have been configured, go to [“Final Configuration”](#) on [page 72](#).

Setting Up the Slave Nodes: Red Hat Enterprise Linux 4.0

This procedure assumes that you are familiar with installing Red Hat Enterprise Linux 4.0 ES. This section contains the following topics:

- [Installing Red Hat Enterprise Linux 4.0 ES](#)
- [Configuring Red Hat Enterprise Linux 4.0 ES](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node **identically** to the other nodes.

Installing Red Hat Enterprise Linux 4.0 ES

This procedure assumes that the Red Hat Enterprise Linux 4.0 ES operating system has not already been installed on the slave nodes. If Red Hat Enterprise Linux 4.0 ES has already been installed then skip to [“Configuring Red Hat Enterprise Linux 4.0 ES”](#) on [page 51](#).

❖ To install Red Hat Enterprise Linux

1. Insert the Red Hat Enterprise Linux Installation 1 of 4 CD into your computer and then restart your computer.

2. When you are prompted for an installation mode, select the graphical mode by pressing **ENTER** or by waiting for a few seconds.
3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Red Hat Enterprise Linux page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Disk Partitioning Setup page, unless you are familiar with installing Red Hat Enterprise Linux, select the **Automatically partition** option and click **Next**.
8. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option and click **Next**.
9. In the Partitioning page, click **Next**.
10. In the Boot Loader Configuration page, select the **GRUB boot loader** and click **Next**.
11. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Disable the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Enter the netmask if it is not entered automatically.
 - c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.

Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
12. In the Firewall Configuration page, select the **No Firewall** option and change the Enable SELinux option to **Disabled**. Click **Next**.
13. In the Additional Language Support page, click **Next**.

14. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
15. In the Set Root Password page, enter the root password for the system and click **Next**.
16. In the Package Defaults page, select **Customize the set of packages to be installed** and click **Next**.
17. Make the following selections and deselections:
 - a. In the Package Group Selection page, deselect the following options if they are selected:
 - Desktops - X Window, GNOME Desktop Environment
 - Applications - Editors, Graphical Internet, Text Based Internet, Office/Productivity, Graphics, Games and Entertainment, Sound and Video
 - Servers - Server Configuration Tools, Web Server, Windows File Server
 - Development - Development Tools
 - System - Printing Support, Administration Tools
 - b. Select the following options in the Servers section:
 - Applications - Scientific and Engineering Programs
 - Servers - FTP Server, Legacy Network Server.
 - c. After selecting the **Legacy Network Server** option, click **Details** and select **rsh-server**.
 - d. After selecting the **Scientific and Engineering Programs** option, click **Details** and select **PVM**. Deselect all other options.
 - e. Click **Next**.
18. Click **Next** to install the Red Hat Linux software. An installation progress dialog box appears. The dialog box prompts which CDs are required during the software loading process.
19. In the Congratulations screen, follow the instructions to finish the installation.

Configuring Red Hat Enterprise Linux 4.0 ES

❖ To configure Red Hat Enterprise Linux

1. Log into your computer as *root*.
2. Run the service configuration program:

```
/usr/sbin/ntsysv
```

2 Setting Up the SEQUEST Cluster

Setting Up the Slave Nodes: Red Hat Enterprise Linux 4.0

3. Scroll down the list and make sure the *rexec*, *rsh*, and *vsftpd* options are active. You can toggle an option by pressing **SPACE** on the keyboard.

4. Disable the sendmail option.

5. Press **TAB** to select the **OK** option and press **ENTER**.

6. Create the proteomics user account with this command:

```
adduser proteomics
```

7. Set the password for this account with this command:

```
passwd proteomics
```

8. Create a directory for the FASTA databases and change the permissions on that directory:

```
mkdir /usr/database  
chmod 777 /usr/database
```

9. Change the permissions of the */usr/share/pvm3/bin/LINUXI386* directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUXI386
```

10. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```

11. Add the following line to the configuration file and then save the file:

```
DatabasePath: /usr/database
```

12. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

13. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.

14. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

15. Add the following line to the file and then save the file:

```
+ proteomics
```

16. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

17. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 30](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 30. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

18. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 30](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 31. *.bashrc* file for Red Hat Enterprise Linux 4.0 ES

```
# .bashrc
# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUXI386
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

19. Save the changes to the file. The slave node is now configured.
20. Restart the computer with this command:


```
shutdown -r now
```
21. Repeat steps 1 through 20 for each of the remaining nodes. When all of the slave nodes have been configured, go to [“Final Configuration”](#) on [page 72](#).

Setting Up the Slave Nodes: Fedora 3.0

This procedure assumes that you are familiar with installing Fedora 3.0. This section contains the following topics:

- [Installing Fedora 3.0](#)
- [Configuring Fedora 3.0](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node **identically** to the other nodes.

Installing Fedora 3.0

This procedure assumes that the Fedora Linux operating system has not already been installed on the slave nodes. If Fedora Linux has already been installed then go to “[Configuring Fedora 3.0](#)” on [page 56](#).

❖ To install Fedora 3.0

1. Insert the Fedora Linux Installation 1 of 4 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing **ENTER** or by waiting for a few seconds.
3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Fedora Linux page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Installation Type page, select **Custom** and click **Next**.
8. In the Disk Partitioning Setup page, unless you are familiar with installing Fedora, select the **Automatically partition** option and click **Next**.
9. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option and click **Next**.
10. In the Partitioning page, click **Next**.
11. In the Boot Loader Configuration page, select the **GRUB boot loader** and click **Next**.
12. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Deselect the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Enter the netmask if it is not entered automatically.

- c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.
Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
13. In the Firewall Configuration page, select the **No Firewall** option and change the Enable SELinux option to **Disabled**. Click **Next**.
 14. In the Additional Language Support page, click **Next**.
 15. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
 16. In the Set Root Password page, enter the root password for the system and click **Next**.
 17. Make the following selections and deselections:
 - a. In the Package Group Selection page, deselect the following options if they are selected:
 - Desktops - X Window, GNOME Desktop Environment
 - Applications - Editors, Graphical Internet, Text Based Internet, Office/Productivity, Graphics, Games and Entertainment, Sound and Video
 - Servers - Server Configuration Tools, Web Server, Windows File Server
 - Development - Development Tools
 - System - Printing support, Language Support
 - b. Select the following options in the Servers section:
 - Applications - Science and Engineering Programs
 - Servers - FTP Server, Legacy Network Server.
 - c. After selecting the **Legacy Network Server** option, click **Details** and select **rsh-server**.

- d. After selecting the **Scientific and Engineering Programs** option, click **Details** and select **PVM**. Deselect all other options.
 - e. Click **Next**.
18. Click **Next** to install the Red Hat Linux software. An installation progress dialog box appears. The dialog box will prompt which CDs are required during the software loading process.
19. In the Congratulations screen, follow the instructions to finish the installation.

Configuring Fedora 3.0

❖ To configure Fedora 3.0

1. Log into your computer as *root*.
2. Run the service configuration program:

```
/usr/sbin/ntsysv
```
3. Scroll down the list and make sure the *rexec*, *rsh*, and *vsftpd* options are active. You can toggle an option by pressing **SPACE** on the keyboard.
4. Disable the sendmail option.
5. Press **TAB** to select the **OK** option and press **ENTER**.
6. Create the proteomics user account with this command:

```
adduser proteomics
```
7. Set the password for this account with this command:

```
passwd proteomics
```
8. Create a directory for the fasta databases and change the permissions on that directory:

```
mkdir /usr/database  
chmod 777 /usr/database
```
9. Change the permissions of the */usr/share/pvm3/bin/LINUXI386* directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUXI386
```
10. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```
11. Add the following line to the configuration file and then save the file:

```
DatabasePath: /usr/database
```
12. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

13. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.

14. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

15. Add the following line to the file and then save the file:

```
+ proteomics
```

16. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

17. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 32](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 32. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

18. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 33](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 33. *.bashrc* file for Fedora 3.0

```
# .bashrc
# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
. /etc/bashrc
fi

PVM_ARCH=LINUXI386
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

19. Save the changes to the file.

The slave node is now configured.

20. Restart the computer with this command:

```
shutdown -r now
```

21. Repeat steps 1 through 20 for each of the remaining nodes.

When all of the slave nodes have been configured, go to “[Final Configuration](#)” on [page 72](#).

Setting Up the Slave Nodes: Fedora 4.0

This procedure assumes that you are familiar with installing Fedora 4.0. This section contains the following topics:

- [Installing Fedora 4.0](#)
- [Configuring Fedora 4.0](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node identically to the other nodes.

Installing Fedora 4.0

This procedure assumes that the Fedora 4.0 operating system has not already been installed on the slave nodes. If Fedora 4.0 has already been installed then skip to “[Configuring Fedora 4.0](#)” on [page 60](#).

❖ To install Fedora 4.0

1. Insert the Fedora Installation 1 of 4 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing **ENTER** or by waiting for a few seconds.
3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Fedora Linux page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Installation Type page, select **Custom** and click **Next**.
8. In the Disk Partitioning Setup page, unless you are familiar with installing Fedora, select the **Automatically partition** option and click **Next**.

9. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option and click **Next**.
10. In the Partitioning page, click **Next**.
11. In the Boot Loader Configuration page, select the **GRUB boot loader** and click **Next**.
12. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Disable the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. Therefore, they should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Enter the netmask if it is not entered automatically.

- c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.

Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
13. In the Firewall Configuration page, select the **No Firewall** option and change the **Enable SELinux** option to **Disabled**. Click **Next**.
14. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
15. In the Set Root Password page, enter the root password for the system and click **Next**.
16. In the Package Defaults page, select Customize the set of packages to be installed and click **Next**.
17. Make the following selections and deselections:
 - a. In the Package Group Selection page, deselect the following options if they are selected:
 - Desktops - X Window, GNOME Desktop Environment
 - Applications - Editors, Graphical Internet, Text Based Internet, Office/Productivity, Graphics, Games and Entertainment, Sound and Video

- Servers - Server Configuration Tools, Web Server, Windows File Server
 - Development - Development Tools
 - System - Printing Support, Language Support
- b. Select the following options in the Servers section:
 - Applications - Scientific and Engineering Programs
 - Servers - FTP Server, Legacy Network Server.
 - c. After selecting the **Legacy Network Server** option, click **Details** and select **rsh-server**.
 - d. After selecting the **Scientific and Engineering Programs** option, click **Details** and select the **PVM** option. Deselect all other options.
 - e. Click **Next**.
18. Click **Next** to install the Fedora software. An installation progress dialog box appears. The dialog box will prompt which CDs are required during the software loading process.
 19. In the Congratulations screen, follow the instructions to finish the installation.

Configuring Fedora 4.0

❖ To configure Fedora 4.0

1. Log into your computer as *root*.
2. Run the service configuration program:

```
/usr/sbin/ntsysv
```
3. Scroll down the list and make sure the **rexec**, **rsh**, and **vsftpd** options are active. You can toggle an option by pressing **SPACE** on the keyboard.
4. Deselect the **sendmail** option.
5. Press **TAB** to select the **OK** option and press **ENTER**.
6. Create the proteomics user account with this command:

```
adduser proteomics
```
7. Set the password for this account with this command:

```
passwd proteomics
```
8. Create a directory for the FASTA databases and change the permissions on that directory:

```
mkdir /usr/database  
chmod 777 /usr/database
```

9. Change the permissions of the `/usr/share/pvm3/bin/LINUXI386` directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUXI386
```

10. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```

11. Add the following line to the configuration file and then save the file:

```
DatabasePath: /usr/database
```

12. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

13. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.

14. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

15. Add the following line to the file and then save the file:

```
+ proteomics
```

16. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

17. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 34](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 34. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

18. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 35](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 35. *.bashrc* file for Fedora 4.0

```
# .bashrc
# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUXI386
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

19. Save the changes to the file. The slave node is now configured.

20. Restart the computer with this command:

```
shutdown -r now
```

21. Repeat steps 1 through 20 for each of the remaining nodes.

When all of the slave nodes have been configured, go to “[Final Configuration](#)” on [page 72](#).

Setting Up the Slave Nodes: Fedora 5.0

This procedure assumes that you are familiar with installing Fedora 5.0. This section contains the following topics:

- [Installing Fedora 5.0](#)
- [Configuring Fedora 5.0](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node identically to the other nodes.

Installing Fedora 5.0

This procedure assumes that the Fedora 5.0 operating system has not already been installed on the slave nodes. If Fedora 5.0 has already been installed then go to “[Configuring Fedora 5.0](#)” on [page 64](#).

❖ To install Fedora 5.0

1. Insert the Fedora Installation 1 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing ENTER or by waiting for a few seconds.

3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Fedora Linux page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Disk Partitioning Setup page, unless you are familiar with installing Fedora, select the **Automatically partition** option. Click **Next**.
8. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option. Click **Next**.
9. In the Partitioning page, click **Next**.
10. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Disable the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Enter the netmask if it is not entered automatically.
 - c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.

Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
11. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
12. In the Set Root Password page, enter the root password for the system and click **Next**.
13. In the Package Defaults page, disable the **Office Package** option and select **Customize Now** option. Click **Next**.

14. In the Package Group Selection page, deselect the following options:
 - Desktop Environments - All items in this section should be deselected.
 - Applications - All items in this section should be deselected except for Engineering and Scientific.
 - Development - All items in this section should be deselected.
 - Servers - All items in this section should be deselected except FTP Servers and Legacy Network Servers.
 - Base - All items in this section should be deselected except for Base and Legacy Software Support.
 - Language - All items in this section should be deselected.
15. Select the **Scientific and Engineering Programs** option. Click **Details** and select the **PVM** option. Deselect all other options.
16. Select the **Legacy Network Server** option. Click **Details** and select **telnet** and **rsh-server**.
17. Select the **Legacy Software Server** option. Click **Details** and select **capability library** options.
18. Click **Next**.
19. Click **Next** to install the Fedora software. An installation progress dialog box appears. The dialog box will prompt which CDs are required during the software loading process.
20. In the Congratulations screen, follow the instructions to finish the installation.

Configuring Fedora 5.0

❖ To configure Fedora 5.0

1. After the computer reboots and if the Setup Agent displays, complete the following:
 - a. Select the Firewall option and disable the firewall.
 - b. Select the Systems Services option and enable the **rexec**, **rsh**, and **vsftpd** options.
 - c. Disable the sendmail option.
 - d. Select **OK** and press ENTER.
2. Log into your computer as *root*. If you ran the Setup Agent in [step 1](#) then go to [step 9](#). [Step 9](#) modifies the selinux configuration file.
3. Run the service configuration program:

```
/usr/sbin/ntsysv
```
4. Scroll down the list and make sure the **rexec**, **rsh**, and **vsftpd** options are active. You can toggle an option by pressing **SPACE** on the keyboard.

5. Disable the **sendmail** option.
6. Press **TAB** to select **OK** and press **ENTER** to save the changes.
7. Disable the firewall with this command:

```
system-config-securitylevel-tui
```
8. Select the disable firewall option and press **TAB** to select **OK** and press **ENTER** to save the changes.
9. Change the selinux (Security Enhanced Linux) configuration file to enable FTP uploads:
 - a. Access the configuration file with this command:

```
vi /etc/selinux/config
```
 - b. Change the line that begins with SELINUX to:

```
SELINUX=permissive
```
 - c. Save the file and the changes.
10. Create the proteomics user account with this command:

```
adduser proteomics
```
11. Set the password for this account with this command:

```
passwd proteomics
```
12. Create a directory for the FASTA databases and change the permissions on that directory:

```
mkdir /usr/database  
chmod 777 /usr/database
```
13. Change the permissions of the `/usr/share/pvm3/bin/LINUX` directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUX
```
14. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```
15. Add the following line to the configuration file (*sequest.ini*) and then save the file:

```
DatabasePath: /usr/database
```
16. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```
17. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.
18. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```
19. Add the following line to the file and then save the file:

```
+ proteomics
```

20. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

21. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 36](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 36. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```

22. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 37](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 37. *.bashrc* file for Fedora 5.0

```
# .bashrc
# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUX
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

23. Save the changes to the file. The slave node is now configured.

24. Restart the computer with this command:

```
shutdown -r now
```

25. Repeat steps 1 through 20 for each of the remaining nodes.

When all of the slave nodes have been configured, go to [“Final Configuration”](#) on [page 72](#).

Setting Up the Slave Nodes: Fedora 6.0

This procedure assumes that you are familiar with installing Fedora 6.0. This section contains the following topics:

- [Installing Fedora 6.0](#)
- [Configuring Fedora 6.0](#)

Note Repeat the following procedures for **each** slave node in the cluster. Configure each slave node identically to the other nodes.

Installing Fedora 6.0

This procedure assumes that the Fedora 6.0 operating system has not already been installed on the slave nodes. If Fedora 6.0 has already been installed then go to “[Configuring Fedora 6.0](#)” on [page 69](#).

❖ To install Fedora 6.0

1. Insert the Fedora Installation 1 CD into your computer and then restart your computer.
2. When you are prompted for an installation mode, select the graphical mode by pressing **ENTER** or by waiting for a few seconds.
3. If you are prompted to test the CD-ROM, select **Skip**.
4. In the Welcome to Fedora Linux page, click **Next**.
5. In the Language Selection page, select the language you want to use and click **Next**.
6. In the Keyboard Configuration page, select what type of keyboard is connected to the computer and click **Next**.
7. In the Disk Partitioning Setup page, unless you are familiar with installing Fedora, select the **Automatically partition** option. Click **Next**.
8. In the Automatic Partitioning page, select the **Remove All Partitions On This System** option. Click **Next**.
9. In the Partitioning page, click **Next**.
10. In the Network Configuration page, set up the network configuration parameters:
 - a. Click **Edit** in the Network Devices box.
 - b. Disable the **Configure using DHCP** option and enter the IP address and netmask for this computer.

The slave computers will be connected to the secondary network card in the master computer through the Ethernet hub. They should have IP addresses in the same group as the secondary network card in the master computer.

Use a private address space (AS) number for the network containing the slave computers. For example: use 192.168.66.1 for the first slave computer, 192.168.66.2 for the second slave computer, and so on until all of the slave computers have been set up.

Enter the netmask if it is not entered automatically.

Disable the IP6 address field.

- c. Click **OK** to save the changes.
 - d. Select the **Manually Set the Hostname** option and enter the hostname.
Name the first slave node in your cluster *node1*, then use name *node2* for the second node in the cluster, and so on for the remaining nodes.
 - e. Enter a gateway address for this network and then click **Next** to save these changes.
 - f. A warning box appears, but because the slave computers do not need DNS, click **Continue** in the warning box.
11. In the Time Zone Selection page, select the correct time zone for your location and click **Next**.
 12. In the Set Root Password page, enter the root password for the system and click **Next**.
 13. In the Package Defaults page, disable the **Office Package** option and select **Customize Now** option. Click **Next**.
 14. In the Package Group Selection page, deselect the following options:
 - Desktop Environments - All items in this section should be deselected.
 - Applications - All items in this section should be deselected except for Engineering and Scientific.
 - Development - All items in this section should be deselected.
 - Servers - All items in this section should be deselected except FTP Servers and Legacy Network Servers.
 - Base - All items in this section should be deselected except for Base and Legacy Software Support.
 - Language - All items in this section should be deselected.
 15. Select the **Scientific and Engineering Programs** option. Click **Details** and select the **PVM** option. Deselect all other options.
 16. Select the **Legacy Network Server** option. Click **Details** and select **telnet** and **rsh-server**.
 17. Select the **Legacy Software Server** option. Click **Details** and select **capability library** options.
 18. Click **Next**.

19. Click **Next** to install the Fedora software. An installation progress dialog box appears. The dialog box will prompt which CDs are required during the software loading process.
20. In the Congratulations screen, follow the instructions to finish the installation.

Configuring Fedora 6.0

❖ To configure Fedora 6.0

1. After the computer reboots and if the Setup Agent displays, complete the following:
 - a. Select the Firewall option and disable the firewall.
 - b. Select the Systems Services option and enable the **rexec**, **rsh**, and **vsftpd** options.
 - c. Disable the **sendmail** option.
 - d. Select **OK** and press **ENTER**.
2. Log into your computer as *root*. If you ran the Setup Agent in [step 1](#) then go to [step 9](#). [Step 9](#) modifies the selinux configuration file.
3. Run the service configuration program:


```
/usr/sbin/ntsysv
```
4. Scroll down the list and make sure the **rexec**, **rsh**, and **vsftpd** options are active. You can toggle an option by pressing **SPACE** on the keyboard.
5. Disable the **sendmail** option.
6. Press **TAB** to select **OK** and press **ENTER** to save the changes.
7. Disable the firewall with this command:


```
system-config-securitylevel-tui
```
8. Select the disable firewall option and press **TAB** to select **OK** and press **ENTER** to save the changes.
9. Change the selinux (Security Enhanced Linux) configuration file to enable FTP uploads:
 - a. Access the configuration file with this command:


```
vi /etc/selinux/config
```
 - b. Change the line that begins with SELINUX to:


```
SELINUX=permissive
```
 - c. Save the file and the changes.
10. Create the proteomics user account with this command:


```
adduser proteomics
```
11. Set the password for this account with this command:


```
passwd proteomics
```

12. Create a directory for the FASTA databases and change the permissions on that directory:

```
mkdir /usr/database
chmod 777 /usr/database
```

13. Change the permissions of the */usr/share/pvm3/bin/LINUX* directory with this command:

```
chmod 777 /usr/share/pvm3/bin/LINUX
```

14. Create a configuration file (*sequest.ini*) for the SEQUEST executable program:

```
vi /usr/share/pvm3/sequest.ini
```

15. Add the following line to the configuration file (*sequest.ini*) and then save the file:

```
DatabasePath: /usr/database
```

16. Create the *hosts.equiv* file:

```
vi /etc/hosts.equiv
```

17. Add the following line to the file and then save the file:

```
Master_computer_name proteomics
```

Where *Master_computer_name* is replaced with the name you used for the master computer.

18. Create the *.rhosts* file:

```
vi ~proteomics/.rhosts
```

19. Add the following line to the file and then save the file:

```
+ proteomics
```

20. Update the permissions for the *.rhosts* file:

```
chmod 600 ~proteomics/.rhosts
```

21. Because the slave computers will not be set up in a domain, modify the existing */etc/hosts* file (see [Figure 38](#)) to include the host names of each slave node:

```
vi /etc/hosts
```

Figure 38. */etc/hosts* file

```
# Do not remove the following line, or various programs
# that require network functionality will fail.

127.0.0.1          localhost.localdomain localhost

192.168.66.250     master
192.168.66.1       node1
192.168.66.2       node2
192.168.66.3       node3
192.168.66.4       node4
192.168.66.5       node5
```


22. Add the PVM environment variables to the existing *.bashrc* file that is in the proteomics account with this command (see [Figure 39](#) for edits):

```
vi ~proteomics/.bashrc
```

Figure 39. *.bashrc* file for Fedora 6.0

```
# .bashrc
# User specific aliases and functions
# Source global definitions

if [ -f /etc/bashrc ]; then
    . /etc/bashrc
fi

PVM_ARCH=LINUX
PVM_ROOT=/usr/share/pvm3
PVM_RSH=/usr/bin/rsh
PVM_TMP=/tmp
PVM_VERSION=3.4.4

export PVM_ARCH PVM_ROOT PVM_RSH PVM_TMP PVM_VERSION
```

23. Save the changes to the file. The slave node is now configured.

24. Restart the computer with this command:

```
shutdown -r now
```

25. Repeat steps 1 through 20 for each of the remaining nodes.

When all of the slave nodes have been configured, go to [“Final Configuration”](#) on [page 72](#).

Updating from SEQUEST Cluster 3.1 or SEQUEST Cluster 3.1 SR1 to SEQUEST Cluster 3.3.1


If you have not yet installed Xcalibur 1.4 SR1 or 2.0.6 and BioWorks 3.3.1 SP1 on the master node of the cluster, do that before proceeding. Refer to the *BioWorks Installation Guide* for more information on installing and licensing BioWorks 3.3.1 SP1.

Once the SEQUEST Cluster 3.3.1 software has been loaded onto the master node of the cluster, you will need to update the configuration of the clusters.

Before starting this process, make sure that SEQUEST and PVM are not running on the system. Open the PVM console and then type **halt** at the PVM prompt to stop the PVM application. For more information on the PVM console, refer to [“PVM Console”](#) on [page 21](#).

❖ To update to SEQUEST Cluster 3.3.1

1. Make sure you are logged on to the cluster master node using an account with Administrator privileges.
2. Start Cluster Manager:

- a. Choose  **Start > Programs > Xcalibur > BioWorks Browser** to Start BioWorks.
- b. Choose **Tools > Cluster Manager** to open the Cluster Manager application.
3. Click the Configuration tab to display the Configuration page.
4. Select the operating system that is installed on each node of your cluster. Once you have selected the OS that is running on each node, click **Save Configuration** to update the configuration.
5. Click **Copy SEQUEST** to copy the SEQUEST executable to each node of the cluster.
6. Reboot the entire cluster.

Note The Copy SEQUEST and Save Configuration buttons are available only when you are logged into an account with Administrator privileges.

Final Configuration

This section has the following topics:

- [Configuration Using the Administrator Login](#)
- [Configuration Using the Proteomics Login](#)

Note You do not need to complete the procedures described in this section if you are upgrading from a previous SEQUEST Cluster installation.

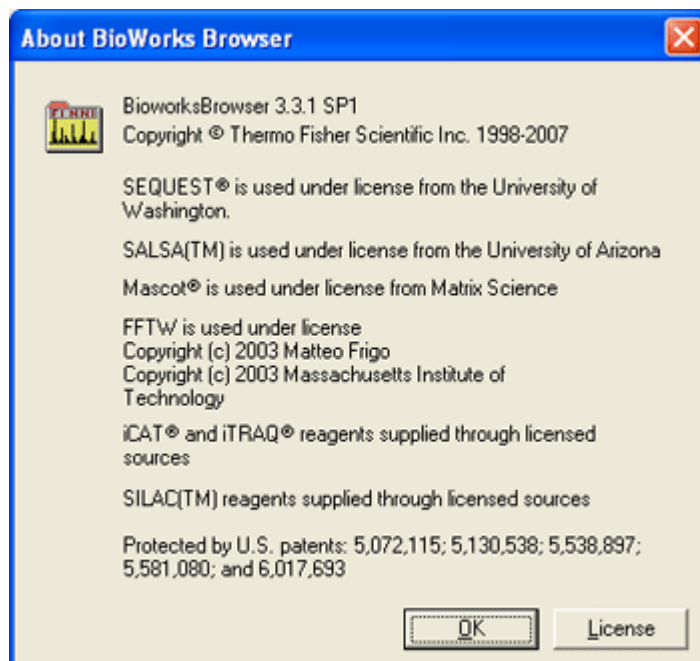
Configuration Using the *Administrator* Login

❖ To complete the final configuration using the administrator login

1. Log into the master node as *administrator*.
2. Start BioWorks Browser.

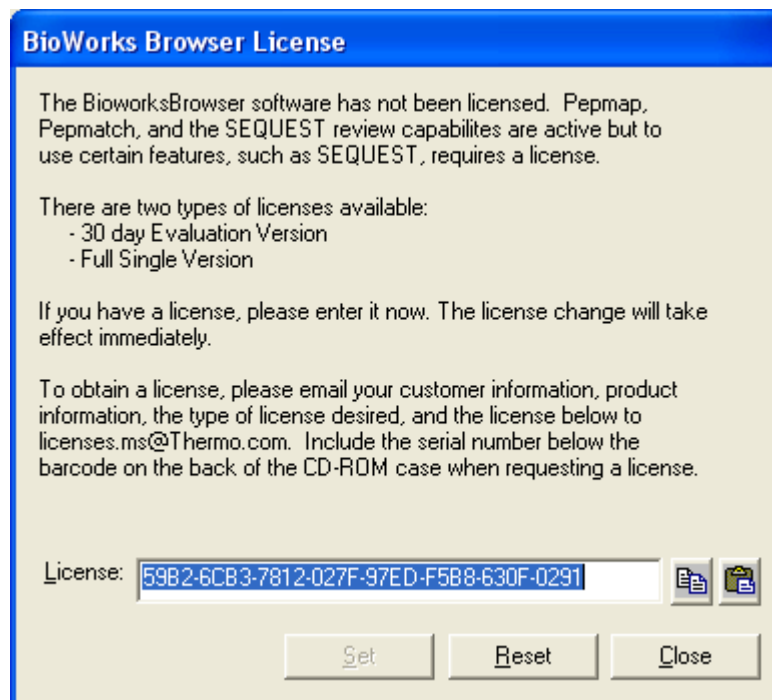
BioWorks Browser starts, but it only allows you to review data. To use the SEQUEST and Cluster software, you need a license key.
3. Choose **Help > About Bioworks Browser** to open the About BioWorks Browser dialog box (see [Figure 40](#)).
4. Click **License**.

Figure 40. About Bioworks Browser dialog box



5. Copy the license key from the License dialog box (see [Figure 41](#)) and e-mail it to licenses.ms@thermo.com. Request that Thermo Scientific enable the SEQUEST/SEQUEST and Cluster features in the updated license. Include the serial number found on the BioWorks 3.3.1 SP1 CD case with your request.
6. When you receive the updated license from Thermo Scientific, enter it into the BioWorks Browser License dialog box. Then, click **Set**.

Figure 41. Bioworks Browser License dialog box



7. Start BioWorks Browser and then open the Cluster Manager program by choosing **Tools > Cluster Manager**. Like BioWorks Browser, the Cluster Manager (see Figure 42) will ask for a license the first time it is started.
8. Copy the license key from the License Manager and e-mail it to licenses.ms@thermo.com. In the e-mail for this license, include the Purchase Order number for your SEQUEST Cluster, the number of CPUs that are licensed, and the serial number from the SEQUEST Cluster CD case.

Figure 42. Cluster Manager License dialog box

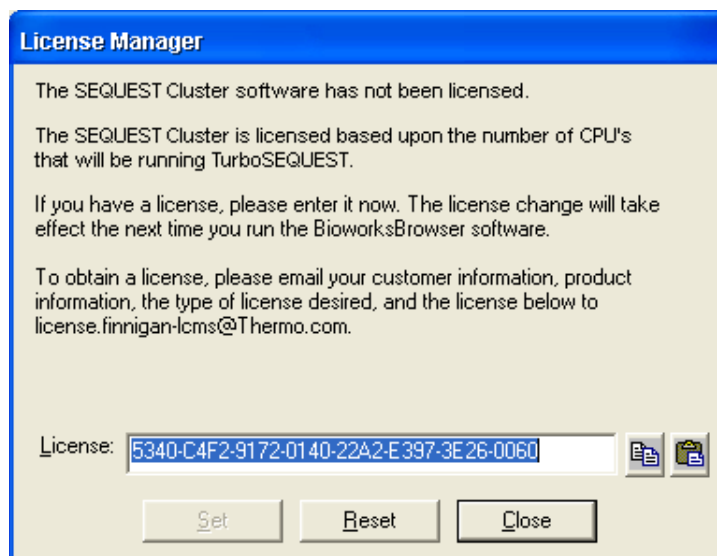


Figure 43. Cluster Manager dialog box

Cluster Manager

Status | Databases | Cluster Daemon | **Configuration**

General Info

Windows

Cluster Share Point:

Database Share Point:

Linux

Cluster Directory:

Database Directory:

User Login Name:

User Password:

Nodes

Total Number of CPU's Licensed:

	Computer	# CPU	OS Type	Status
1	node1	2	RedHat 8.0	
2	node2	2	RedHat 8.0	
3	node3	2	RedHat 8.0	
4	node4	2	RedHat 8.0	
5	node5	2		
6	node6	2	Windows 2000	
7	node7	2	RedHat 7.2	
8			RedHat 8.0	
9			RedHat 9.0	
10			RedHat Enterprise 3.0	

9. When you receive the updated license from Thermo Fisher Scientific, enter it into the License Manager dialog box. Then, click **Set**.
10. Start the Cluster Manager again. The default display for the Cluster Manager is shown in [Figure 43](#). Select the Configuration page (see [Figure 44](#)).
11. If you are using Linux, you must enter your User Login Name and User Password.
12. Do one of the following:
 - If your slave nodes are running Microsoft Windows 2000 or Microsoft Windows XP, then in the Windows group box, enter the share points that were created during the Slave Node Configuration in the **Cluster Share Point** and **Database Share Point** text boxes. Normally, these values are:

Cluster (sharepoint for the PVM software)

Database (sharepoint for the FASTA databases)

- If your slave nodes are running Red Hat Linux, Red Hat Enterprise Linux, or Fedora, then in the Linux group box, enter the directory location of the PVM software and FASTA databases into the **Cluster Directory** and **Database Directory** text boxes. Normally, these values are:

/usr/share/pvm3 (location of PVM software)

/usr/database (location of FASTA databases)

Because the account *proteomics* was created during the installation of the Linux software, enter that login name and its password in the User Login Name and User Password text boxes.

Figure 44. Cluster Manager dialog box - Configuration page, showing node information

Cluster Manager

Status | Databases | Cluster Daemon | **Configuration**

General Info

Windows:

Cluster Share Point: Cluster

Database Share Point: Database

Linux:

Cluster Directory: /usr/share/pvm3

Database Directory: /usr/database

User Login Name:

User Password:

Nodes

Total Number of CPU's Licensed: 16

	Computer	# CPU	OS Type	Status
1	node1	2	RedHat 8.0	
2	node2	2	RedHat 8.0	
3	node3	2	RedHat 8.0	
4	node4	2	RedHat 8.0	
5	node5	2	RedHat 8.0	
6	node6	2	RedHat 8.0	
7	node7	2	RedHat 8.0	
8				
9				
10				

About ClusterManager... License... Save Configuration Copy SEQUEST

Exit Help

IMPORTANT Do not enter your computer's full qualified domain name. Enter your computer name such as, my_laptop.

13. For each slave node of your cluster, enter the computer name and number of CPUs that are present in that computer. Then, select the type of OS that is installed on that slave node.

14. Click **Save Configuration** to create the PVM configuration file.

Note The PVM Configuration is found in the directory *C:\Cluster* and is named *pvm_hosts.txt*. Example PVM hosts files are shown in [Figure 45](#) and [Figure 46](#).

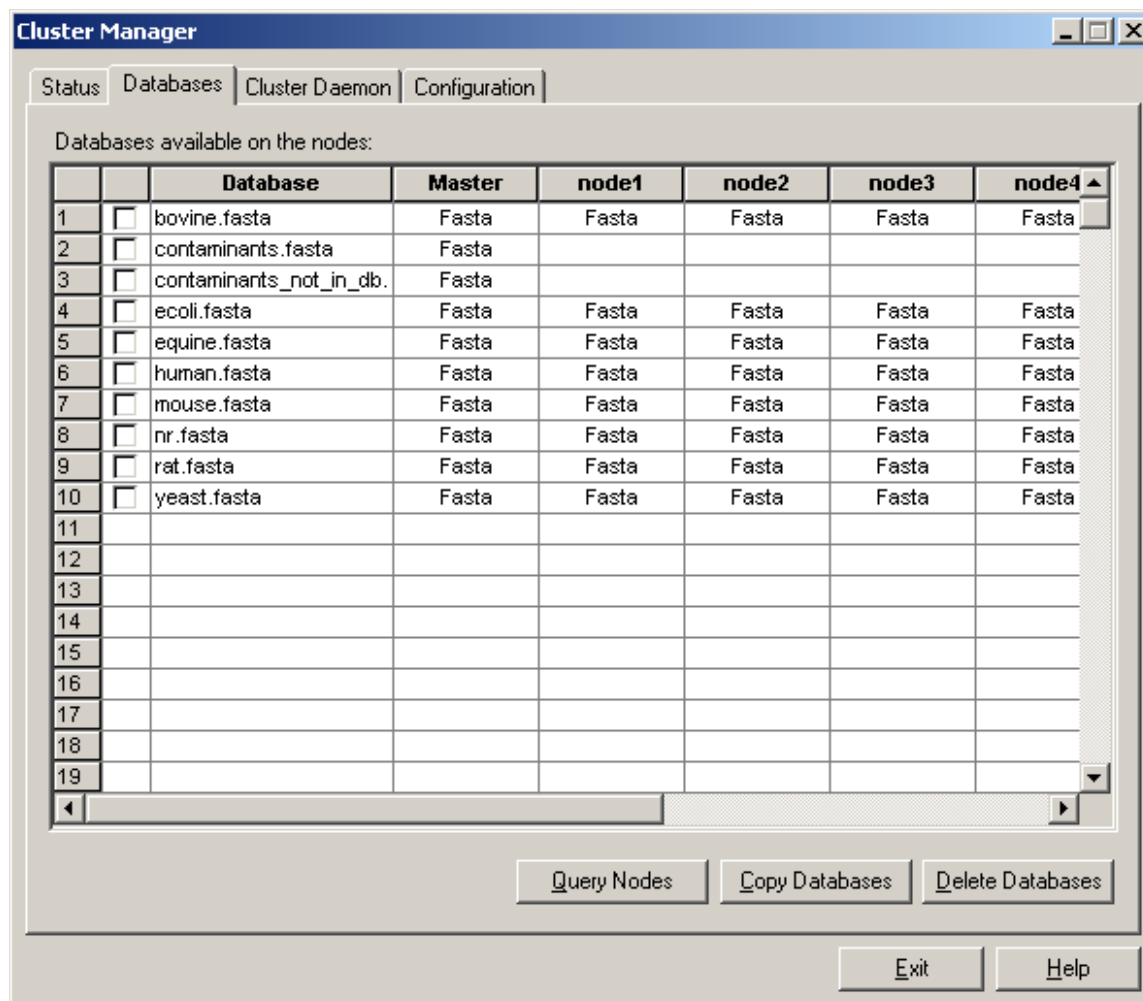
[Figure 46](#) is a Red Hat Linux 7.2 or 8.0 example. If you are using, a later version of Red Hat Linux, replace LINUX with i386 for each node path.

```
node1 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node2 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node3 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node4 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node5 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node6 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node7 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
node8 dx=c:\cluster\pvm\lib\win32\pvmd3.exe sp=2
```

```
node1 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node2 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node3 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node4 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node5 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node6 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node7 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
node8 dx=/usr/share/pvm3/lib/LINUX/pvmd3 sp=2
```

15. With the Configuration page set up, you can also copy the Sequest task to each slave/compute node of the cluster by clicking **Copy SEQUEST**.
16. Now that the Cluster Manager is configured, you can copy the FASTA databases that were installed onto the master computer to the slave/compute nodes. Do this through the Databases page of the Cluster Manager (see [Figure 45](#)).
17. To copy the FASTA databases to the slave nodes, select each database that you want to copy (by clicking in the second column of the table) and then click **Copy Database**. When the databases have been copied, the program automatically updates the display, by performing the same action as **Query Nodes**.

Figure 45. Cluster Manager dialog box - Databases page



Configuration Using the *Proteomics* Login

❖ To complete the final configuration using the proteomics login

1. Log out of the computer (as *administrator*) and log in to the computer as *proteomics*.
2. Open a MS/DOS console window and change to the directory *C:\Cluster\wrshd*.
3. Type one of the following:

- If the slave node is running Microsoft Windows 2000 or Microsoft Windows XP:

```
rsh node1 c:\cluster\wrshd\id.exe
```

This command displays the output shown in [Figure 46](#).

Figure 46. Output from id.exe command

```

Select PVM Shell
Microsoft Windows 2000 [Version 5.00.2195]
(C) Copyright 1985-2000 Microsoft Corp.

c:\cluster\pvm>cd ..\wrshd

C:\Cluster\wrshd>rsh cluster1 c:\cluster\wrshd\id.exe
(Using stderr:) You are logged on at cluster1 as proteomics.
(Using stdout:) You are logged on at cluster1 as proteomics.
Please enter a string:
Thanks, I got: '' on Standart Input. Bye!
One more time:
Thanks again. Bye

C:\Cluster\wrshd>_

```

- If the slave node is running Red Hat Linux, Red Hat Enterprise Linux, or Fedora:

```
rsh node1 "cat /etc/hosts"
```

This command displays the contents of the hosts file on the slave node, as shown in [Figure 47](#).

Figure 47. Output from "cat /etc/hosts" command

```

PVM Shell
(C) Copyright 1985-2000 Microsoft Corp.

c:\cluster\pvm>cd ..\wrshd

C:\Cluster\wrshd>rsh node1 cat /etc/hosts
127.0.0.1      localhost.localdomain  localhost
172.20.3.25   clustertwo node0
172.20.3.1    node1
172.20.3.2    node2
172.20.3.3    node3
172.20.3.4    node4
172.20.3.5    node5
172.20.3.6    node6
172.20.3.7    node7
172.20.3.8    node8
172.20.3.9    node9
172.20.3.10   node10
172.20.3.11   node11
172.20.3.12   node12
172.20.3.13   node13
172.20.3.14   node14
172.20.3.15   node15
172.20.3.16   node16

C:\Cluster\wrshd>_

```

After the Final Configuration

If the master and slave nodes are configured as described in this chapter, then the following locations will contain the SEQUEST program files:

Microsoft Windows 2000 Server / Windows 2003 Server (master node)

```
c:\cluster\pvm\bin\WIN32\sequest_master.exe  
c:\cluster\pvm\bin\WIN32\sequest_slave.exe
```

Microsoft Windows 2000 Professional / Windows XP Professional (slave node)

```
c:\cluster\pvm\bin\WIN32\sequest_slave.exe
```

Red Hat Linux / Red Hat Enterprise Linux / Fedora (slave node)

```
/usr/share/pvm3/bin/LINUX/sequest_slave
```

Files Modified During the Installation

The files that are modified during installation include the following:

Microsoft Windows 2000 Server / Windows 2003 Server (master node)

```
C:\cluster\pvm  
C:\cluster\wrshd  
C:\cluster\temp
```

Microsoft Windows 2000 Professional / Windows XP Professional (slave node)

```
C:\cluster\pvm  
C:\cluster\wrshd  
C:\cluster\temp
```

Red Hat Linux / Red Hat Enterprise Linux / Fedora (slave node)

```
/etc/hosts  
/etc/hosts.equiv  
/home/proteomics/.rhosts  
/home/proteomics/.bashrc  
/usr/share/pvm3/sequest.ini  
/usr/database
```

Using the SEQUEST Cluster

This chapter describes how to use the SEQUEST cluster.

Contents

- [SEQUEST Cluster Manager](#)
- [Configuring BioWorks for Use with SEQUEST Cluster](#)
- [Running a SEQUEST Search on the Cluster](#)
- [PVM Console](#)

SEQUEST Cluster Manager

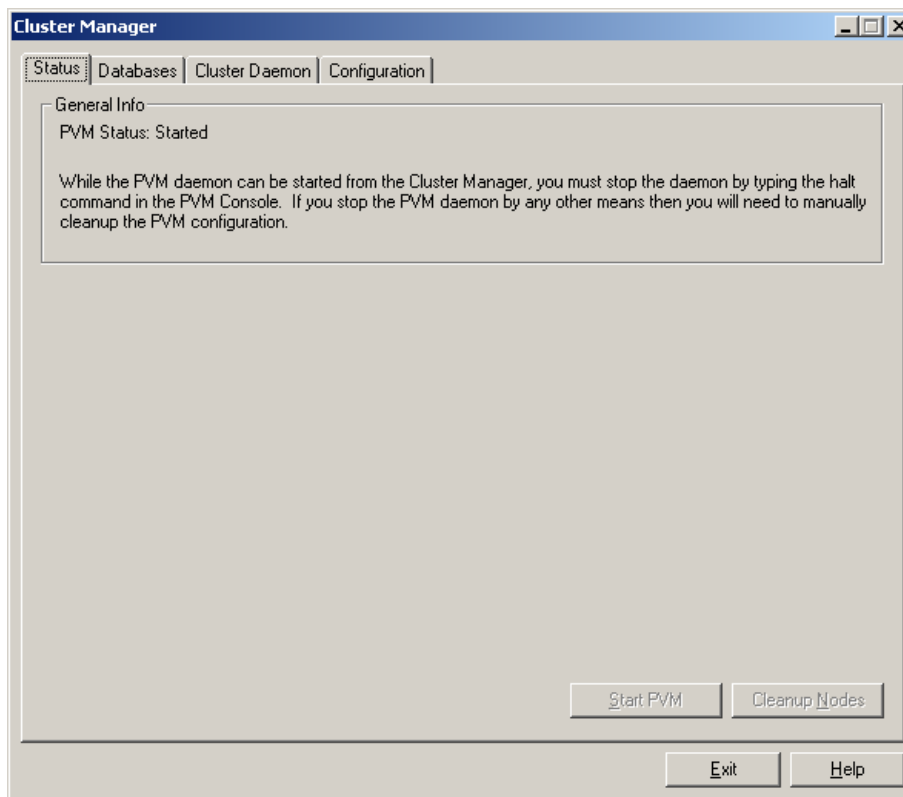
The Cluster Manager program provides tools to do the following:

- Install the SEQUEST Cluster license
- Configure the PVM nodes
- Distribute the SEQUEST application to the slave nodes
- Start and stop the PVM software
- Maintain the FASTA databases on each slave node
- Submit SRF files to the Cluster for processing

In “[Final Configuration](#)” on [page 72](#), the Cluster Manager was used to install the SEQUEST Cluster license, to configure the PVM nodes, and to distribute the SEQUEST application to the slave nodes. Refer to that section for information on these tasks. The remaining tasks are discussed in this section.

When you start the Cluster Manager program (in the BioWorks window, choose **Tools > Cluster Manager**), the Status page is displayed. See [Figure 48](#). The purpose of the Status page is to start PVM and to clean up the PVM operation files on each node of the cluster.

Figure 48. Cluster Manager dialog box - Status page



Starting and Stopping the PVM Software

When the Cluster Manager is open, it periodically checks to see if PVM is running, and, if it is, the Cluster Manager displays a status message. When PVM is running the Start PVM and Cleanup Nodes buttons are unavailable.

To start the PVM software, click **Start PVM**. This starts the pvmd3 program, which acts as the main console for PVM. The PVM console runs in a separate console window. This window must be running for the PVM software to function correctly. [Figure 48](#) displays the Status page when PVM is running.

To stop the PVM software, you must display the MS/DOS Console window that was started by the Cluster Manager and type “halt”, and then press **ENTER**. Do not close the PVM console window using the X button in the upper-right corner of the window. Doing so does not stop PVM on each of the slave nodes. If this occurs, you must manually terminate the PVM files on each node of the cluster by clicking **Cleanup Nodes**.

When PVM is running, it creates some operation files (*.pvmd) in the PVM temporary directory. If you stop PVM using the halt command, these files are automatically removed. However, if you do not stop PVM by using the halt command, these files are not removed. PVM will not start if old versions of these operation files are found in the temporary directory. Click **Start PVM** to automatically remove the operation files.

Once PVM has been started, the Cluster Manager can be closed without it affecting the running PVM.

More information about the PVM Console can be found in “[PVM Console](#)” on [page 96](#).

Managing Databases

You use the Databases page of the Cluster Manager (see [Figure 49](#)) to copy and remove FASTA and indexed databases from the slave nodes of the cluster.

Note Some of the functionality of this page of the Cluster Manager was already demonstrated in “[Final Configuration](#)” on [page 72](#).

For the Databases page to work properly, it needs to know where the databases are stored on the slave nodes. This information was specified in “[Final Configuration](#)” on [page 72](#). For slave nodes that are using:

- Windows 2000 or Windows XP, a shared directory named Database was created on each slave node. This share point is specified on the Configuration page.
- Linux or Fedora, the actual directory location of the databases must be specified on the Configuration page because an ftp connection is used to copy the databases to the slave nodes.

When Cluster Manager is started, only the local databases will be displayed. To find out what databases are present on the other nodes, click **Query Nodes**.

❖ To copy a database

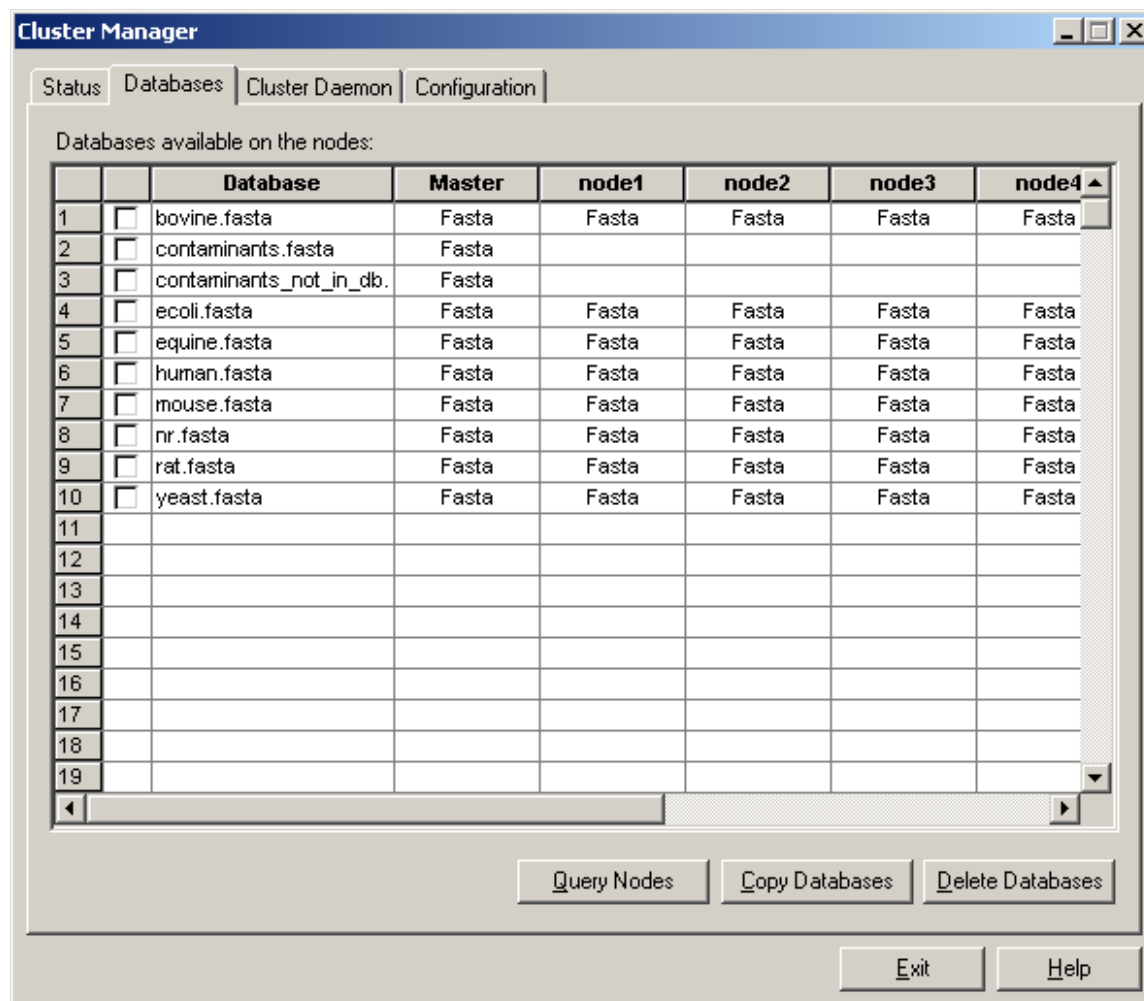
1. To copy a database from the master node, select the check box in the first column of the table that corresponds to the database that you want to copy.
2. Click **Copy Databases**.

The application copies the database to each node of the cluster. While the databases are being copied, a status dialog box appears displaying which database is being copied to which node. The program copies the FASTA database plus the associated files for that database (*.idx*, *.hdr*, and *.dgt*).

❖ To delete a database

1. To delete databases from the nodes, select a database.
2. Click **Delete Databases**. While the databases are being deleted, a status dialog box appears displaying which database are being deleted from which node.

Figure 49. Cluster Manager dialog box - Databases page



Submitting SRF Files to the Cluster for Processing

The Cluster Daemon allows a remote instrument to submit Unified Search (SRF) files to the SEQUEST Cluster for processing. The Cluster Daemon continuously watches a specified directory for the presence of SRF files. When it finds the files, it submits a SEQUEST search to the SEQUEST Queue.

The Cluster Daemon requires a shared directory on the cluster. This directory is specified in the Inbound Folder text box on the Cluster Daemon page. See [Figure 50](#).

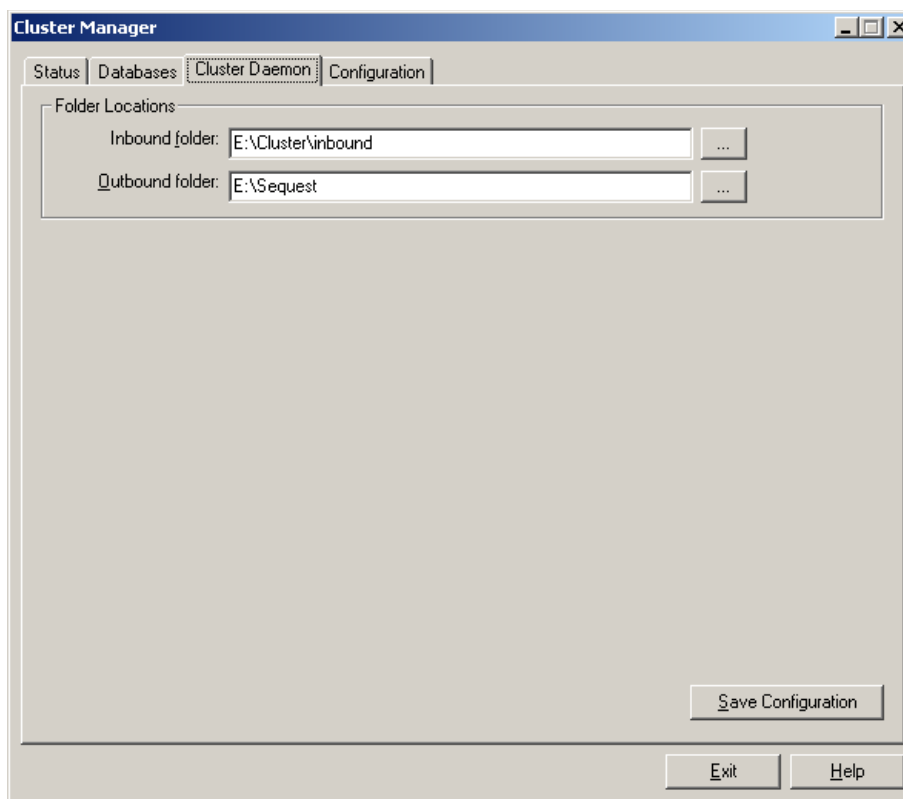


Figure 50. Cluster Manager dialog box - Cluster Daemon page

Share this folder as you did with the PVM and Database folders during the configuration of the slave nodes.

When searches are complete, the SRF files containing the results can be found in the folder specified in the Outbound Folder text box.

Configuring BioWorks for Use with SEQUEST Cluster

SEQUEST Cluster 3.3.1 allows you to start searches on the cluster in two ways. Searches can be started using BioWorks

- Running on the master node of the cluster.
- Running on a properly configured remote client computer.

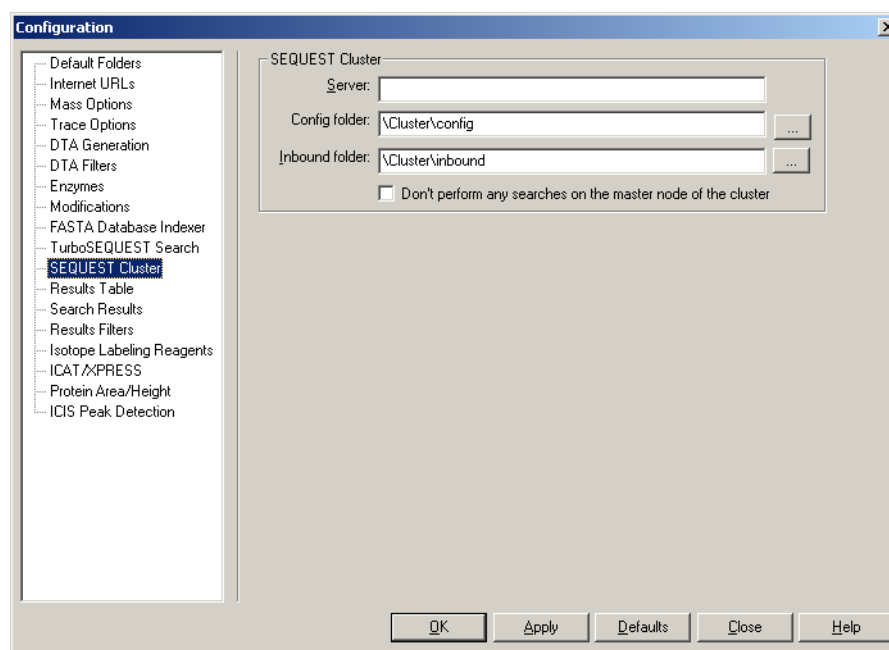
To use either of these types of searches you must first configure BioWorks 3.3.1 SP1 to work with the cluster, as described in the following sections.

Configuring BioWorks on the Cluster Master Node

❖ To configure BioWorks on the master node of SEQUEST Cluster

1. Choose **Start > Programs > Xcalibur > BioWorks Browser** to launch BioWorks.
2. Choose **Options > Configuration**. The Configuration dialog box appears.
3. Select **SEQUEST Cluster** from the tree control to display the SEQUEST Cluster page.
See [Figure 51](#).

Figure 51. Configuration dialog box - SEQUEST Cluster page



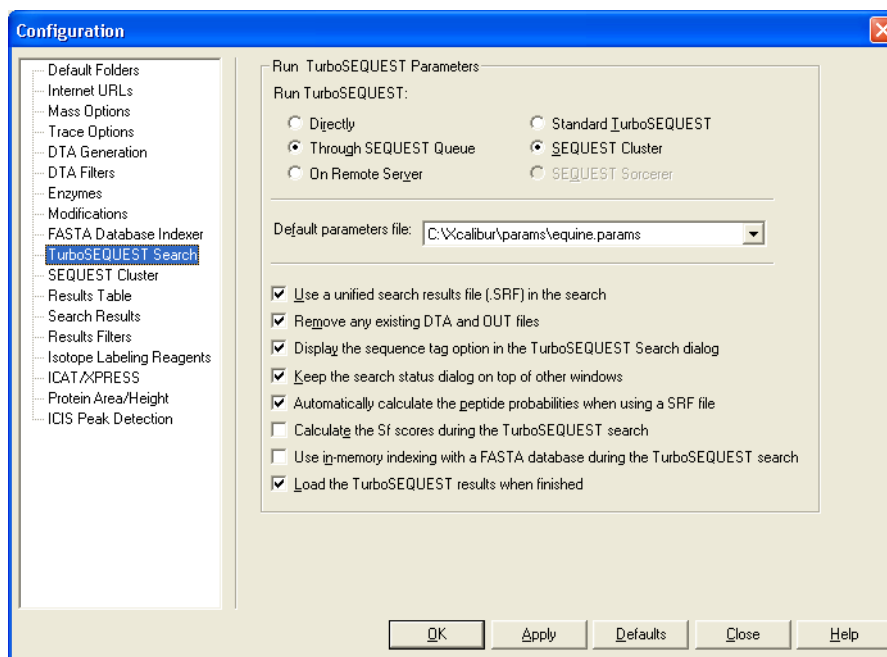
4. Set the parameters on the SEQUEST Cluster page:
 - a. Leave the **Server** text box blank.
 - b. Use the default values in the Config Folder and Inbound Folder text boxes.
 - c. If you do not want to perform searches on the master node of the cluster, select the **Don't Perform Any Searches on the Master Node of the Cluster** check box.
5. Click **OK** to save the settings and close the Configuration dialog box.

After you have set the SEQUEST Cluster configuration options, you can also set the SEQUEST Search options to run searches on the cluster by default. These configuration settings are optional, and can be changed when the search is launched.

❖ **To set SEQUEST Search options on the master node**

1. Choose **Start > Programs > Xcalibur > BioWorks Browser** to launch BioWorks.
2. Choose **Options > Configuration**. The Configuration dialog box appears.
3. Select **SEQUEST Search** from the tree control to display the SEQUEST Search page. See [Figure 52](#).
4. Set the parameters in the Run SEQUEST Parameters group box:
 - a. Select the **SEQUEST Cluster** option.
 - b. Select how you want to run the search:
 - Select **Directly** to run the search directly on the cluster.
 - Select **Through SEQUEST Queue** to run the search through the SEQUEST Queue. This option is recommended if multiple users can submit searches to SEQUEST Cluster at the same time. (For example, if users can submit searches from both the master node and remote client computers.)
5. Set the other parameters on the SEQUEST Search page. For more information on these parameters, refer to Chapter 2: “Configuration” in the *BioWorks User Guide*.
6. Click **OK** to save the settings and close the dialog box.

Figure 52. Configuration dialog box - SEQUEST Search page, showing settings for searches from the master node of the cluster



Configuring BioWorks on a Remote Client Computer

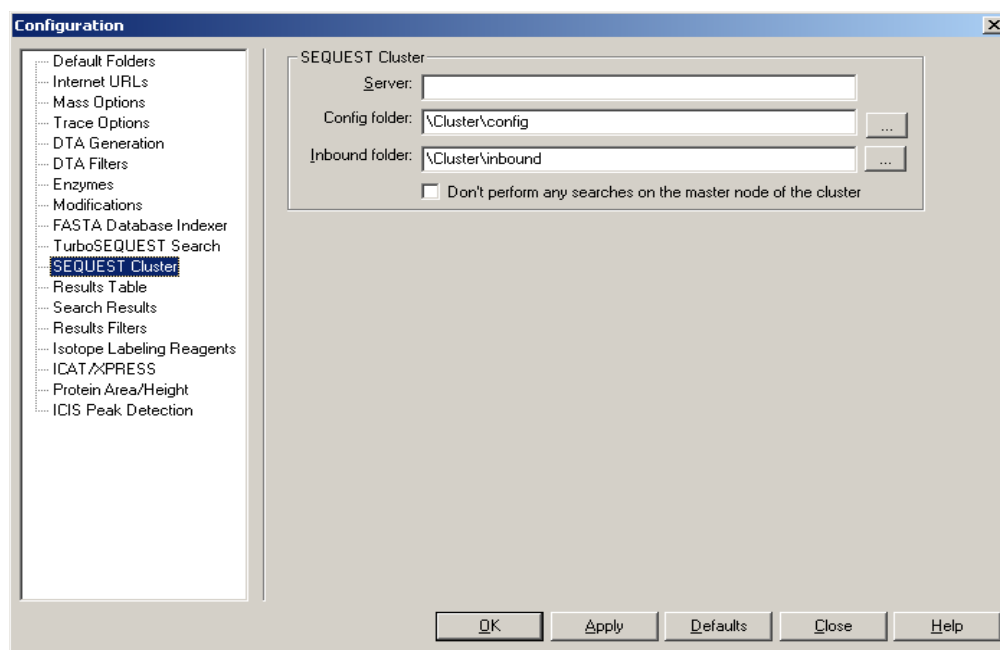
To launch searches on the cluster, the remote client computer must meet the following requirements:

- The client computer must be capable of accessing the cluster over the network. Consult your IT department if you need assistance in network configuration.
- The client computer must have BioWorks 3.3.1 SP1 installed and licensed for use with SEQUEST Cluster. Refer to the *BioWorks Installation Guide* for information on installing BioWorks and obtaining a license.
- BioWorks on the client computer must be properly configured to work with the cluster. This section describes how to configure BioWorks 3.3.1 SP1 for use with the cluster.

❖ To configure BioWorks for remote searches on SEQUEST Cluster

1. Choose **Start > All Programs > Xcalibur > BioWorks Browser** to launch BioWorks.
2. Choose **Options > Configuration**. The Configuration dialog box appears.
3. Select **SEQUEST Cluster** from the tree control to display the SEQUEST Cluster page. See [Figure 53](#).

Figure 53. Configuration dialog box - SEQUEST Cluster page



4. Set the parameters on the SEQUEST Cluster page:
 - a. In the **Server** text box, enter the server name of the SEQUEST Cluster master node.
 - b. Leave the default values in the Config Folder and Inbound Folder text boxes.
5. Click **OK** to save the settings and close the Configuration dialog box.

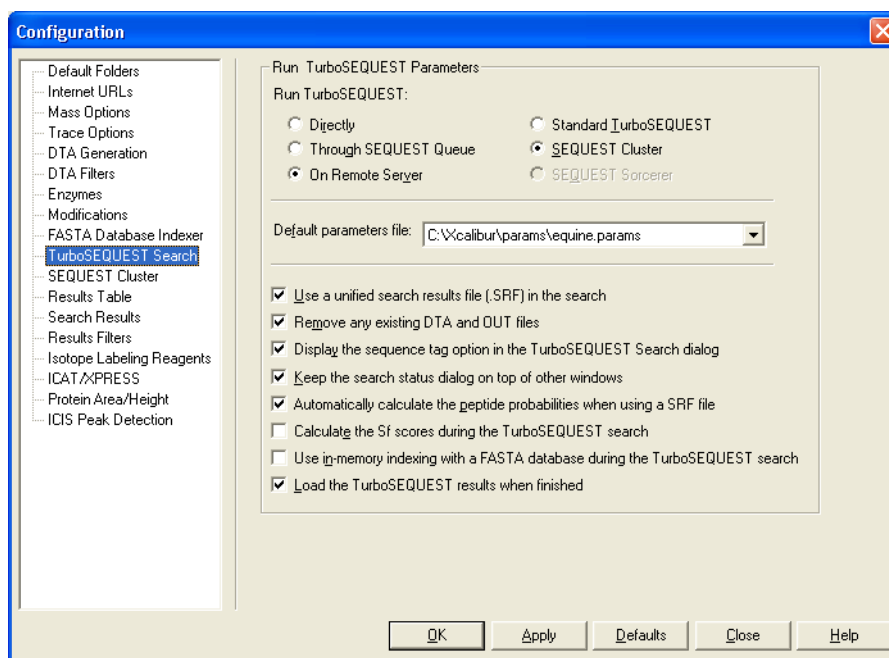
Note The **Don't Perform Any Searches on the Master Node of the Cluster** check box has no effect if it is selected in the configuration of a remote client computer. This option is only valid when running a search from the master node of the cluster.

After you have set the SEQUEST Cluster configuration options, you can also set the SEQUEST Search options to run searches on the cluster by default. These configuration settings are optional, and can be changed when the search is launched.

❖ **To set SEQUEST Search options on a remote client computer**

1. Choose **Start > Programs > Xcalibur > BioWorks Browser** to launch BioWorks.
2. Choose **Options > Configuration**. The Configuration dialog box appears.
3. Select **SEQUEST Search** from the tree control to display the SEQUEST Search page. See [Figure 54](#).
4. Set the parameters in the Run SEQUEST Parameters group box ([Figure 54](#)):
 - a. Select the **SEQUEST Cluster** option button.
 - b. Select the **On Remote Server** option button.
5. Set the other parameters on the SEQUEST Search page. For more information on these parameters, refer to Chapter 2: "Configuration" in the *BioWorks User Guide*.
6. Click **OK** to save the settings and close the dialog box.

Figure 54. Configuration dialog box - SEQUEST Search page, showing settings for a remote client computer



Running a SEQUEST Search on the Cluster

The SEQUEST Cluster can be run in BioWorks from the following locations:

- SEQUEST Search dialog box
- SEQUEST Batch search dialog box
- Search Selected Samples dialog box

The procedure for starting a search is similar whether you are starting the search from the master node of the cluster or from a remote client computer. Where there are differences, these differences are noted in the procedures that follow.

Starting a Search Using the SEQUEST Search Dialog Box

This topic describes only the basic steps necessary for running a search on the SEQUEST Cluster. For detailed instructions on setting SEQUEST search parameters, refer to Chapter 3: “Performing Search” in the *BioWorks User Guide*.

❖ To start a SEQUEST Search on the cluster

1. Start with the RAW data file you want to search open in BioWorks.
2. Choose **Actions > SEQUEST Search**. The SEQUEST Search dialog box appears.
3. Click the **SEQUEST Queue** tab to display the SEQUEST Queue page. See [Figure 55](#).

Figure 55. SEQUEST Search dialog box - SEQUEST Queue page



4. Select where to run the search:
 - If you are running the search from the master node of the SEQUEST Cluster, select either the **Directly** option button or the **Through SEQUEST Queue** option button.
 - If you are running the search from a remote client computer, select the **On Remote Server** option button.
5. Select the **SEQUEST Cluster** option.
6. Click the **Basic** tab to return to the Basic page.
7. In the **Search Parameters** list, select a parameters file to use in the search.
8. Click **Edit** to display the SEQUEST Search Parameters dialog box.
9. In the **Database** list, select the database you want to use in the search.

You must select a database that has been copied to the cluster master node and slave nodes. If you are running a search from a remote client computer, databases that have been copied to the master node are marked “Cluster” in the list.

To check which databases have been copied to the master node and the slave nodes, use the Database page of the SEQUEST Cluster Manager as described in [“Managing Databases”](#) on [page 83](#).

Figure 56. SEQUEST Search Parameters dialog box, showing a Cluster database selected on a remote client computer

SEQUEST Search Parameters

Basic | Modifications

General Parameters

Parameters file: C:\Xcalibur\params\AvePrecFrag-PartialEnz-equine.param ...

Database: equine (FASTA) ...

Type: ☒ Protein ☐ Nucleotide

Mass type: Average precursor and fragments

Enzyme: Trypsin(KR)

Enzyme limits: Partially enzymatic - cleaves at either end

Missed cleavage sites: 2

Peptide tolerance: 2.0000 Units: ☒ AMU ☐ MMU ☐ PPM

Fragment ions tolerance: 1.0 Units: ☒ AMU ☐ MMU

Number results scored: 250

☐ Remove precursor ion from spectra

Ions and Ion Series Calculated

☐ A Ions ☒ B Ions ☐ C Ions ☒ Y Ions ☐ Z Ions

Output Options

Peptide matches reported: 10

☐ Report duplicate peptide matches 10

Save Defaults Close Help

10. Set the other parameters in the SEQUEST Search Parameters dialog box. Refer to “Setting the SEQUEST Search Parameters” in the *BioWorks User Guide* for more information.
11. When you have finished settings parameters, click **Save** to save your changes and close the SEQUEST Search Parameters dialog box.
12. Set the other parameters in the SEQUEST Search dialog box. For more information, refer to “Setting and Saving the Search Parameters” of the *BioWorks User Guide*.

Note When you run a search on the cluster from a remote client computer, only SRF files are supported. Searches in DTA and OUT format are not supported.

If you are searching from a remote client computer, you must copy the SRF file from the cluster back to the remote client computer when the search is finished in order to view your results on the remote client computer.

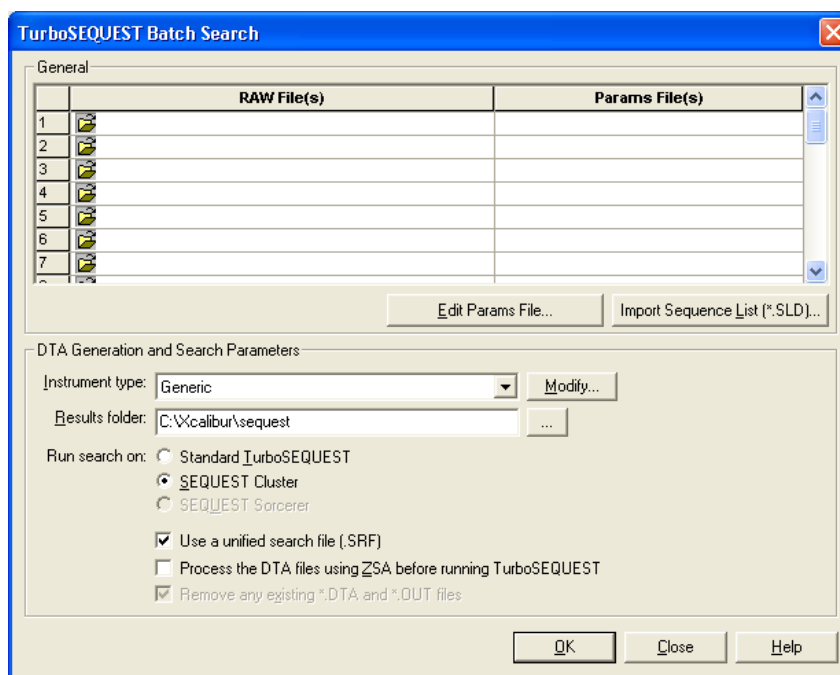
Starting a Search Using the SEQUEST Batch Search Dialog Box

This section describes only the basic steps required to do a search on the SEQUEST Cluster. For detailed instructions on setting SEQUEST search parameters, refer to Chapter 3: “SEQUEST Search” in the *BioWorks User Guide*.

❖ To start a SEQUEST Batch Search on the cluster

1. From the BioWorks Browser window, open the SEQUEST Batch Search dialog box by choosing **Actions > SEQUEST Batch Search**. See [Figure 57](#).

Figure 57. SEQUEST Batch Search dialog box



2. Select the **Run Search On: SEQUEST Cluster** option button.
3. Set the other parameters in the SEQUEST Batch Search dialog box. Refer to “Performing a SEQUEST Batch Search” on page 3-17 of the *BioWorks User Guide* for more information.
4. For each .params file you have selected to use in the search, edit the .params file to ensure that the correct database is selected:
 - a. Click **Edit Params File** to open the SEQUEST Search Parameters dialog box.
 - b. In the **Parameters File** text box, enter the name of the .params file you want to edit.
 - c. In the **Database** list, select the database you want to use in the search.

You must select a database that has been copied to the cluster master node and slave nodes. If you are running a search from a remote client computer, databases that have been copied to the master node are marked “Cluster” in the list. You can check which databases have been copied to the master node and the slave nodes using the Database page of the SEQUEST Cluster Manager as described in “[Managing Databases](#)” on [page 83](#).

- d. Set the other parameters in the SEQUEST Search Parameters dialog box. Refer to “Setting the SEQUEST Search Parameters” of the *BioWorks User Guide* for more information.
 - e. Click **Save** to save the changes and close the SEQUEST Search Parameters dialog box.
 - f. Repeat steps a through e for each .params file you want to edit.
5. To Start the search click **OK**.

Note When you run a search on the cluster from a remote client computer, only SRF files are supported. Searches in DTA and OUT format are not supported.

If you are searching from a remote client computer, you must copy the SRF file(s) from the cluster back to the remote client computer when the search is finished in order to view your results on the remote client computer.

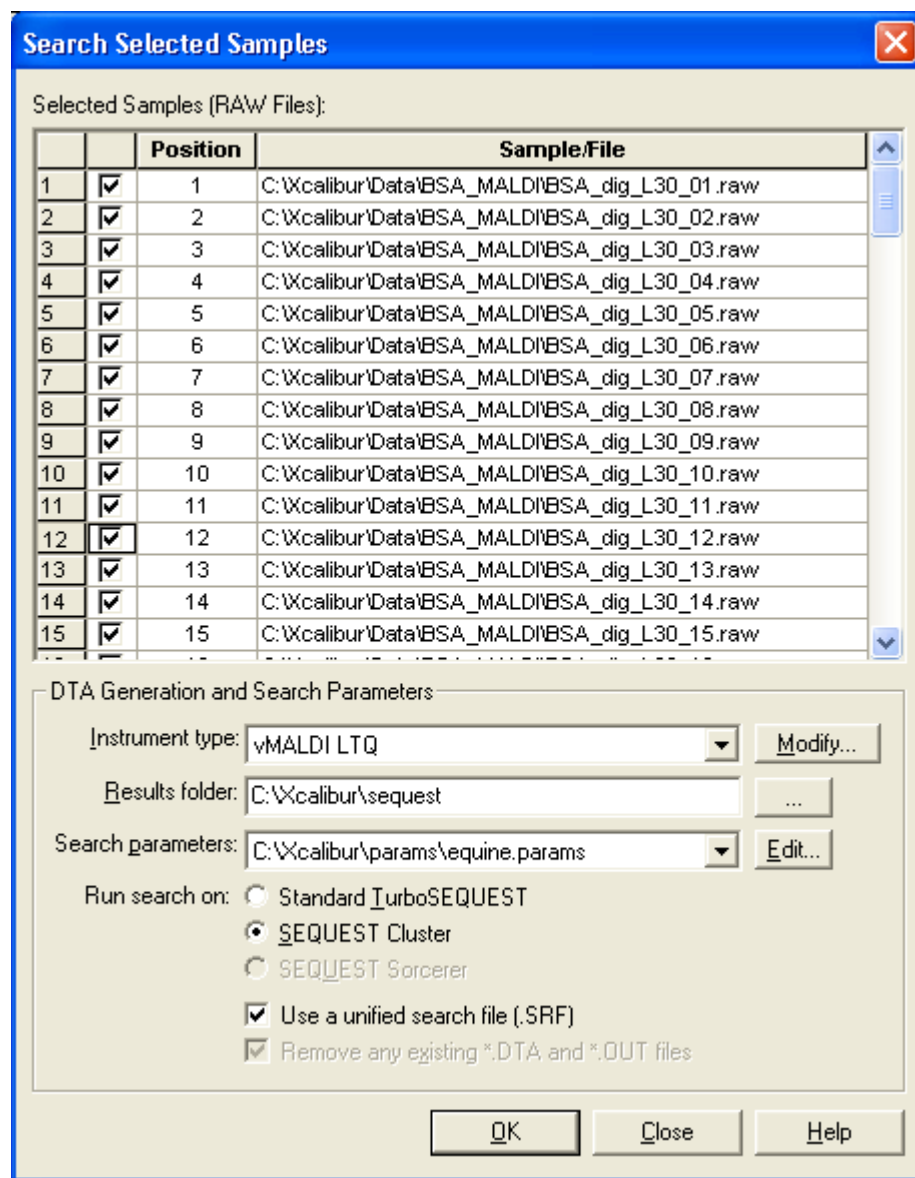
Starting a Search Using the Search Selected Samples Dialog Box

This section describes only the basic steps required to do a search on the SEQUEST Cluster. For detailed instructions on setting SEQUEST search parameters, refer to Chapter 3: “SEQUEST Search” in the *BioWorks User Guide*.

❖ To start a SEQUEST Search from the Search Selected Samples dialog box

1. From the BioWorks Browser window, load a set of MALDI Samples into the MALDI Plate view as described in “Searching vMALDI Samples in BioWorks” of the *BioWorks User Guide*.
2. Select the spots in the plate view corresponding to the samples you want to search, and click **Search Selected Samples** to display the Search Selected Samples dialog box.

Figure 58. Search Selected Samples dialog box



3. Select the **Run Search On: SEQUEST Cluster** option.
4. Set the other parameters in the Search Selected Samples dialog box. Refer to “Searching vMALDI Samples in BioWorks” in the *BioWorks User Guide* for more information.
5. For the .params file you have selected to use in the search, edit the .params file to ensure that the correct database is selected:
 - a. Click **Edit** to open the SEQUEST Search Parameters dialog box.
 - b. In the **Database** list, select the database you want to use in the search.

You must select a database that has been copied to the cluster master node and slave nodes. If you are running a search from a remote client computer, databases that have been copied to the master node are marked “Cluster” in the list.

To check which databases have been copied to the master node and the slave nodes use the Database page of the SEQUEST Cluster Manager as described in [“Managing Databases”](#) on [page 83](#).

- c. Set the other parameters in the SEQUEST Search Parameters dialog box. Refer to “Setting the SEQUEST Search Parameters” of the *BioWorks User Guide* for more information.
 - d. Click **Save** to save the changes and close the SEQUEST Search Parameters dialog box.
6. Start the search by clicking **OK**.

Note When you run a search on the cluster from a remote client computer, only SRF files are supported. Searches in DTA and OUT format are not supported.

If you are searching from a remote client computer, you must copy the SRF file(s) from the cluster back to the remote client computer when the search is finished in order to view your results on the remote client computer.

PVM Console

When the PVM daemon application is running, it maintains a console mode that allows you to interact with the daemon. The PVM Console contains its own set of commands that are documented in the PVM Manual (*PVM: Parallel Virtual Machine: A Users' Guide and Tutorial for Network Parallel Computing (Scientific and Engineering Computation)* by Al Geist, Adam Bequelin, and Jack Dongarra (ISBN# 0262571080)). This section highlights only the commands that a SEQUEST Cluster user may need.

When PVM is running on the SEQUEST Cluster, an MS-DOS console window will be present. If this window is maximized, the following display is shown (see [Figure 59](#)).

Figure 59. PVM Console

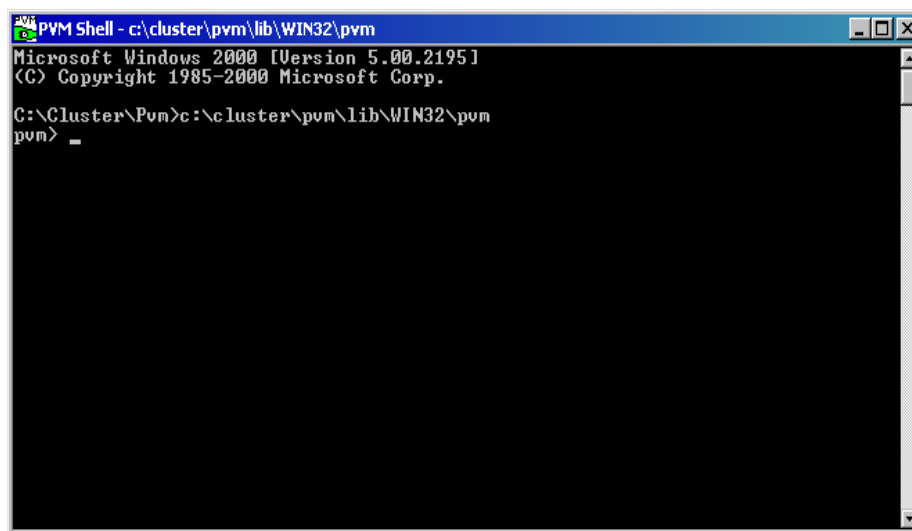


Table 1 contains those PVM commands that are used with the SEQUEST Cluster.

Table 1. PVM commands and descriptions

Command	Description
conf	Display the virtual machine configuration
halt	Stops the PVM daemon
help	Display the PVM console help
jobs	Displays a list of running jobs
kill	Terminates a PVM task
ps	Lists the PVM tasks
reset	Kills a running tasks
version	Lists the PVM version

The *conf* command (see Figure 60) lists the computers that are running with the current PVM session. It is helpful in diagnosing if all of the nodes have been started. The Speed column indicates how many CPUs are being used on each node of the cluster. By default, the Speed value is used to provide a weighting as to how fast each machine is on the cluster. The default value is 1000. Values less than 1000, indicate how many CPUs are being used on that node. So, for the nodes cluster1 to cluster8, 2 CPUs are being used in this configuration.

Figure 60. Example conf command

```

c:\cluster\pvm\lib\WIN32\pvm.exe
pvm> conf
conf
9 hosts, 1 data format
      HOST      DTID      ARCH      SPEED      DSIG
clusterone    40000    WIN32      1000  0x00408841
cluster1      80000    WIN32        2  0x00408841
cluster2      c0000    WIN32        2  0x00408841
cluster3     100000    WIN32        2  0x00408841
cluster4     140000    WIN32        2  0x00408841
cluster5     180000    WIN32        2  0x00408841
cluster6     1c0000    WIN32        2  0x00408841
cluster7     200000    WIN32        2  0x00408841
cluster8     240000    WIN32        2  0x00408841
pvm>

```

The *ps* and *kill* commands (see Figure 61) are very similar to their UNIX counterparts. The *ps* command is used to list the active PVM processes on each node of the cluster. The *ps ax* command lists all the process and gives additional information about the processes. The TID value is process ID number and is used with the *kill* command to remove a process. The command *kill 1c0046* removes the process associated with task 1c0046. The *reset* command removes all processes from the PVM process list.

Figure 61. Example of ps and kill commands

```

c:\cluster\pvm\lib\WIN32\pvm.exe
pvm> ps ax
ps ax
      HOST      TID      FLAG 0x COMMAND
clusterone    40001    4/c c:\cluster\pvm\bin\WIN32\hoster.exe
clusterone    <cons>    4/c -
clusterone    40065    4/c sequest27_output.exe
clusterone    40066    4/c sequest27_slave.exe
clusterone    4006b    4/c sequest27_output.exe
clusterone    4006c    4/c sequest27_slave.exe
cluster1      80041    4/c sequest27_slave.exe
cluster1      80042    4/c sequest27_slave.exe
cluster3     100001    4/c pvmgs.exe
cluster3     100042    4/c sequest27_slave.exe
cluster3     100043    4/c sequest27_slave.exe
cluster3     100046    4/c sequest27_slave.exe
cluster3     100047    4/c sequest27_slave.exe
cluster5     180041    4/c sequest27_slave.exe
cluster5     180042    4/c sequest27_slave.exe
cluster5     180045    4/c sequest27_slave.exe
cluster5     180046    4/c sequest27_slave.exe
cluster6     1c0041    4/c sequest27_slave.exe
cluster6     1c0042    4/c sequest27_slave.exe
cluster6     1c0045    4/c sequest27_slave.exe
cluster6     1c0046    4/c sequest27_slave.exe
pvm> kill 1c0046

```

The *version* command (see Figure 62) provides information about the PVM version.

The *halt* command stops the PVM daemon on each node of the cluster and removes the pvmd and pvml session files that are created on each node of the cluster. This command should be run before any maintenance is done to the cluster if the PVM program is running.

```

c:\cluster\pvm\lib\WIN32\pvm.exe

pvm>
pvm> version
version
3.4.3
pvm> conf
conf
9 hosts, 1 data format
      HOST      DTID      ARCH      SPEED      DSIG
clusterone  40000  WIN32    1000  0x00408841
cluster1    80000  WIN32     2  0x00408841
cluster2    c0000  WIN32     2  0x00408841
cluster3   100000  WIN32     2  0x00408841
cluster4   140000  WIN32     2  0x00408841
cluster5   180000  WIN32     2  0x00408841
cluster6   1c0000  WIN32     2  0x00408841
cluster7   200000  WIN32     2  0x00408841
cluster8   240000  WIN32     2  0x00408841
pvm>
pvm> version
version
3.4.3
pvm>

```

Figure 62. Output from version command

PVM depends on five environmental variables to work correctly. [Table 2](#) lists and describes these variables.

Table 2. Environmental variables used by PVM

Variable	Microsoft Windows 2000	Red Hat LINUX
PVM_ARCH	WIN32	LINUX or LINUXI386
PVM_ROOT	C:\Cluster\Pvm	/usr/share/pvm3
PVM_TMP	C:\Cluster\Temp	/tmp
PVM_RSH		/usr/bin/rsh
PVM_VERSION	3.4.3	3.4.4

Note Depending upon your configuration, some of these values may vary from what is shown above.

Troubleshooting

If you are having trouble using your cluster, refer to the following common problems and solutions as given in this chapter.

Contents

- Computer Name and Login Account Name Are the Same
- Account Name and WRSHD Login Are Not the Same
- AutoLogon Feature of Microsoft Windows 2000 Is Not Set
- Need Additional Linux Subsets
- PVM Will Not Start
- Extra Configuration Entries in WRSHD
- WRSHD Cannot Create Service
- WRSHD Will Not Start on Microsoft Windows 2000 / Windows XP
- RSH Does Not Work Between Microsoft Windows 2000 Server / Windows 2003 Server and Linux Nodes
- RSH Is Not Working Correctly
- Extra Files From an Aborted PVM Session
- Poor Cluster Performance
- Cluster Searches Will Not Run Through SEQUEST Queue

Computer Name and Login Account Name Are the Same

PVM cannot work on a system that uses the same name for a computer and the login account under which PVM is run. For example, if you use *proteomics* as the login account name then you do not want to name any of the computers *proteomics*.

Account Name and WRSHD Login Are Not the Same

Because the RSH is set up with the account name *proteomics*, you must be logged in as *proteomics* for the PVM to work correctly.

Note If you decided to use a different account name (instead of *proteomics*) as the login name then you use that account name when configuring the RSH software.

AutoLogon Feature of Microsoft Windows 2000 Is Not Set

If during the initial setup of the slave nodes, the automatic logon feature has not been set, then you can manually configure the software using the following procedure.

Microsoft Windows includes a feature that enables you to configure the computer to automatically logon to the network, bypassing the Winlogon dialog box. To enable this function you need to add several new values to the registry key.

❖ To configure the AutoLogon feature

1. Choose **Start > Run** to open the Run dialog box.
2. Type **regedit** and then click **OK** to open the Registry Editor.
3. Locate the following entry in the registry:

```
HKEY_LOCAL_MACHINE\SOFTWARE\Microsoft\Microsoft Windows  
NT\CurrentVersion\Winlogon.
```

4. Add a new string value named *DefaultUserName* and set it to the username you want to use to automatically logon. For example: **proteomics**.
5. Add a new string value named *DefaultPassword* and set this to the password for the user entered above. For example, the password for **proteomics**.
6. Add a new string value named *AutoAdminLogon* and set it to 1 to enable auto logon.
7. Exit the Registry Editor and restart your computer.

Once you have configured AutoLogon, Microsoft Windows does not stop at the login box and wait for you to enter the login information; instead, it automatically proceeds to the desktop corresponding to the default username.

Note The password is stored in registry, which means anyone who has access to the machine has access to the password. You can bypass this function by holding down <Shift> during the boot or logoff process. It is also important to note that if the DontDisplayLastUserName value is enabled, the auto logon feature does not function.

Need Additional Linux Subsets

This procedure shows how to load the RSH Server, PVM and VSFTPD subsets. Other subsets may also be installed using this procedure but you will have to use the name of the specific subset instead of the RSH Server, PVM or VSFTPD names. Additionally, these other subsets may be on a different CD-ROM than is shown in these procedures—so you might have to search each of the CD-ROMs that comes with the Red Hat software.

Note The locations and package names used in this example are based upon Red Hat Linux 8.0.

❖ To load a Linux subset after the operating system has been installed

1. Search for any RSH related subsets that are installed on the system:

```
rpm -qa | grep -i rsh
```

2. Search for the PVM or VSFTPD subsets that you would replace rsh with either pvm or ftp.

3. Install the RSH-Server software, insert the CD-ROM labeled *Installation CD 2 of 3* into the CD-ROM drive, and type:

```
mount /dev/cdrom
```

4. Change to the RedHat/RPMS directory:

```
cd /mnt/cdrom/RedHat/RPMS
```

5. Install the RSH-Server package:

```
rpm -Uhv rsh-server-0.17-10.i386.rpm
cd /
umount /dev/cdrom
```

6. Install the VSFTPD and PVM software, insert the CD-ROM labeled *Installation CD 3 of 3* into the CD-ROM drive and type:

```
mount /dev/cdrom
```

7. Change to the RedHat/RPMS directory:

```
cd /mnt/cdrom/RedHat/RPMS
```

8. Install the VSFTPD package:

```
rpm -Uhv wu-ftpd-2.6.2.-8.i386.rpm
```

9. Install the PVM software:

```
rpm -Uhv pvm-3.4.4-6.i386.rpm
```

10. Remove the CD-ROM:

```
umount /dev/cdrom
```

Extra Configuration Entries in WRSHD

If you need to remove a configuration entry from the WRSHD program, use this command:

```
configure -remove username
```

Where you replace *username* with the name that you want to remove.

Note To run this command you must be logged in as *administrator*.

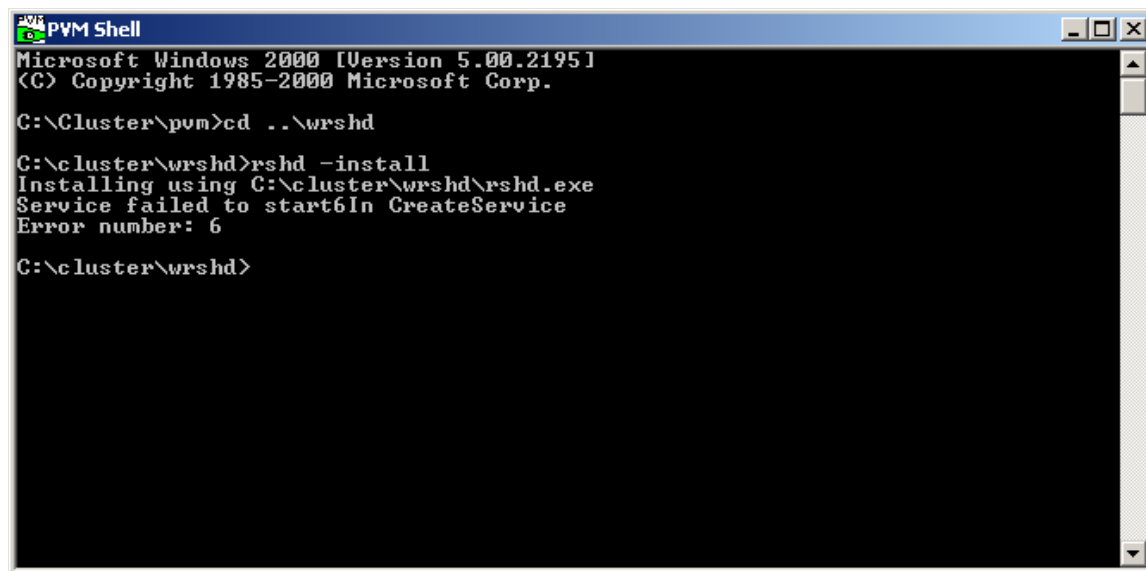
WRSHD Cannot Create Service

If WRSHD was already installed on the system, removed using the Uninstall Microsoft Windows command, and then re-installed on the system, then use this command:

```
rshd -install
```


might generate an error (see [Figure 63](#)).

Figure 63. WRSHD CreateService error



In this case, do the following:


1. Uninstall the software.

2. Choose  **Start** > **Run** to open the Run dialog box.
3. Type “regedit” and then click **OK** to open the Registry Editor.
4. Locate the following entry in the registry:

```
HKEY_LOCAL_MACHINE\SYSTEM\CurrentControlSet\Services\BsdRshd.
```

5. Delete the entry and close the Registry Editor.

WRSHD Will Not Start on Microsoft Windows 2000 / Windows XP

The WRSHD on Microsoft Windows 2000/ Windows XP is run as a Microsoft Windows Service ( **Start** > **Settings** > **Control Panel** > **Administrative Tools** > **Services**) under the name *BsdRshd*.

You can stop and start the daemon from this dialog box by double-clicking on the *BsdRshd* name in the list of services and clicking on the Start or Stop button in the *BsdRshd Properties* dialog box.

To manually start the daemon, stop the BsdRshd daemon, and then, from a console window, use these commands:

```
c:
cd \cluster\wrshd
rshd -d
```

When the BsdRshd daemon is started in debug mode the program prints information to the console window as it receives commands. You can use this information to debug the RSH problems.

Additional information about the WRSHD program can be found in the help file *wrshd.hlp* in the directory *c:\cluster\wrshd*.

RSH Does Not Work Between Microsoft Windows 2000 Server / Windows 2003 Server and Linux Nodes

If the RSH does not work between the Windows 2000 Server / Windows 2003 Server master node and a Linux slave node, then check the log file */var/log/messages* on the Linux node to see what is wrong and correct the problem.

To view the contents of this value use this command:

```
cat /var/log/messages
```

RSH Is Not Working Correctly

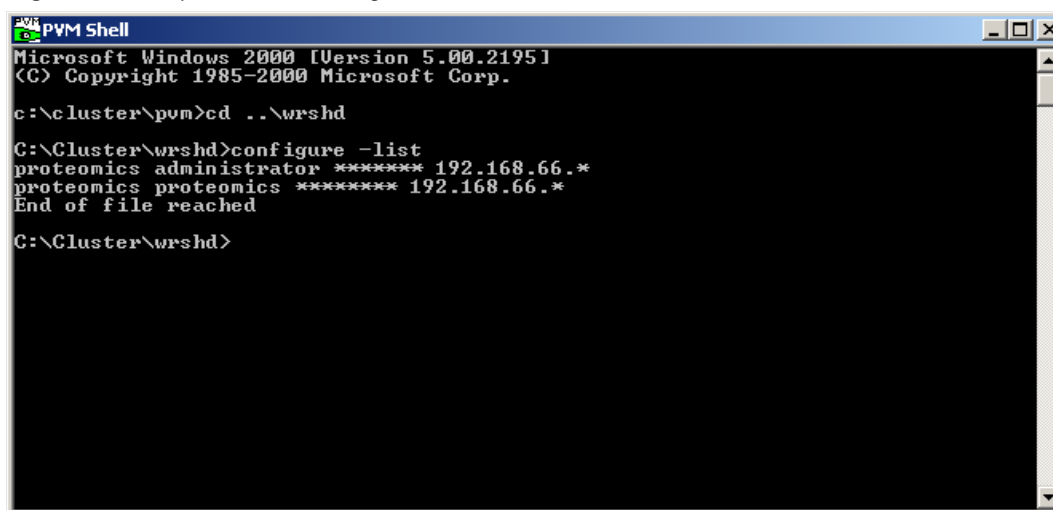
If it appears that the RSH isn't working correctly, you can test it with the following steps:

1. Log in to the master computer as *administrator*.
2. Open a MS/DOS console window and change to the directory *C:\Cluster\wrshd*.
3. Check the RSH configuration, use this command:

```
configure -list
```

This command displays the output shown in [Figure 64](#).

Figure 64. Output from the configure command



```
PVM Shell
Microsoft Windows 2000 [Version 5.00.2195]
(C) Copyright 1985-2000 Microsoft Corp.

c:\cluster\pvm>cd ..\wrshd

C:\Cluster\wrshd>configure -list
proteomics administrator ***** 192.168.66.*
proteomics proteomics ***** 192.168.66.*
End of file reached

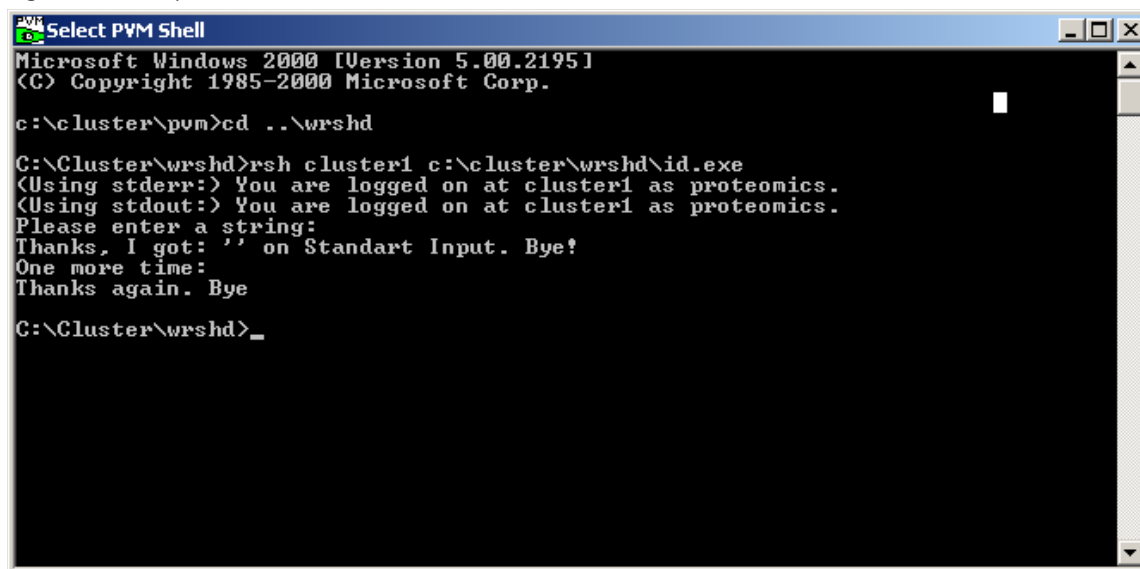
C:\Cluster\wrshd>
```

4. Type this command to test the RSH command if the slave node is running Microsoft Windows 2000 / Windows XP:

```
rsh node1 -l proteomics c:\cluster\wrshd\id.exe
```

This command displays the output shown in [Figure 65](#).

Figure 65. Output from id.exe command



```

Select PVM Shell
Microsoft Windows 2000 [Version 5.00.2195]
(C) Copyright 1985-2000 Microsoft Corp.

c:\cluster\pvm>cd ..\wrshd

C:\Cluster\wrshd>rsh cluster1 c:\cluster\wrshd\id.exe
(Using stderr:) You are logged on at cluster1 as proteomics.
(Using stdout:) You are logged on at cluster1 as proteomics.
Please enter a string:
Thanks, I got: '' on Standart Input. Bye!
One more time:
Thanks again. Bye

C:\Cluster\wrshd>_

```

5. If the slave node is running Red Hat Linux, use this command:

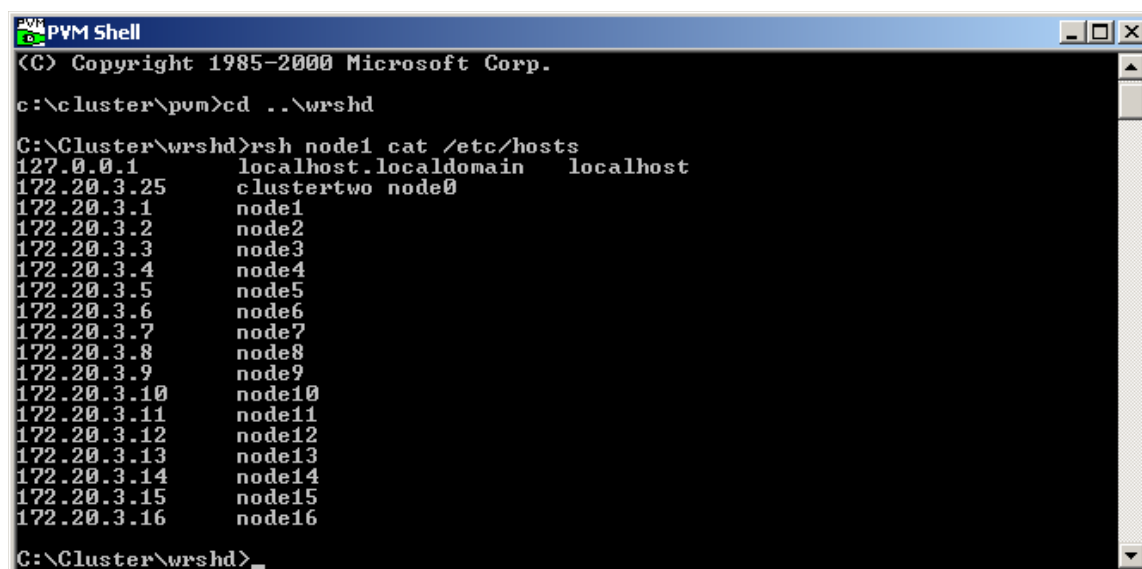
```
rsh node1 -l proteomics "cat /etc/hosts"
```

This command displays the contents of the hosts file on the slave node as shown in [Figure 66](#).

6. You can repeat these commands between the slave nodes and master computer by logging into one of the slave nodes and use this command:

```
rsh master "c:\cluster\wrshd\id.exe"
```

Figure 66. Output from "cat /etc/hosts" command



```

PVM Shell
(C) Copyright 1985-2000 Microsoft Corp.

c:\cluster\pvm>cd ..\wrshd

C:\Cluster\wrshd>rsh node1 cat /etc/hosts
127.0.0.1      localhost.localdomain  localhost
172.20.3.25   clustertwo node0
172.20.3.1    node1
172.20.3.2    node2
172.20.3.3    node3
172.20.3.4    node4
172.20.3.5    node5
172.20.3.6    node6
172.20.3.7    node7
172.20.3.8    node8
172.20.3.9    node9
172.20.3.10   node10
172.20.3.11   node11
172.20.3.12   node12
172.20.3.13   node13
172.20.3.14   node14
172.20.3.15   node15
172.20.3.16   node16

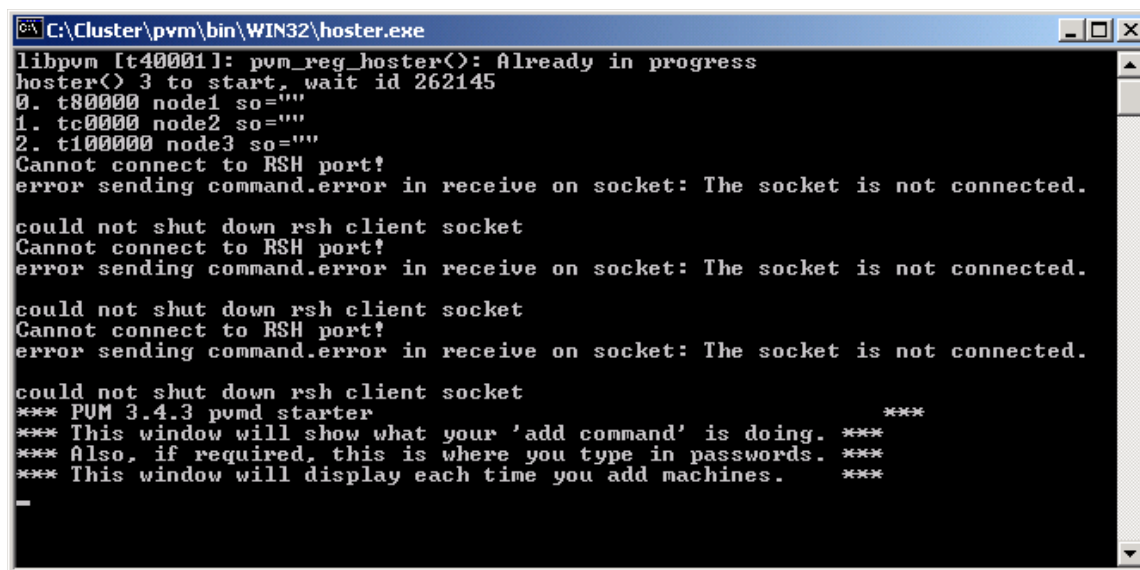
C:\Cluster\wrshd>_

```

PVM Will Not Start

If, during the startup of PVM, any error messages occur (see [Figure 67](#)), the messages are written to the *pvm.proteomics* file in the directory *c:\cluster\temp*. This file is a text file that can be opened in Microsoft Notepad. Additional information is included in this file that is not shown in the console window. Check this file first if PVM does not start.

Figure 67. PVM Startup with errors



```
C:\Cluster\pvm\bin\WIN32\hoster.exe
libpvm [t400011]: pvm_reg_hostor(): Already in progress
hoster(>) 3 to start, wait id 262145
0. t80000 node1 so=""
1. tc0000 node2 so=""
2. t100000 node3 so=""
Cannot connect to RSH port!
error sending command.error in receive on socket: The socket is not connected.

could not shut down rsh client socket
Cannot connect to RSH port!
error sending command.error in receive on socket: The socket is not connected.

could not shut down rsh client socket
Cannot connect to RSH port!
error sending command.error in receive on socket: The socket is not connected.

could not shut down rsh client socket
*** PVM 3.4.3 pvm starter ***
*** This window will show what your 'add command' is doing. ***
*** Also, if required, this is where you type in passwords. ***
*** This window will display each time you add machines. ***
```

In most cases, if PVM was running correctly in the past, you can fix this type of problem by restarting the master node of the cluster. If, after restarting the master computer, the problem still exists, then you need to reboot the entire cluster.

If the error messages in the PVM console state that the PVM nodes cannot be found, then the binding order for the network cards might not be correct. This message is also found in the *pvm* file in the PVM temp directory (normally, *c:\cluster\temp*).

Windows 2000 Server and Windows 2003 Server associate the computer name with the networking protocol that is considered the initial protocol in the loading order for the network cards. In Windows 2000 Server or Windows 2003 Server, this order can be changed with the following steps:

1. From the Windows Taskbar, choose **Start > Settings > Control Panel**.
2. Double-click on the Network and Dial-up Connections icon.
3. Choose **Advanced > Advanced Settings** from the menu.
4. In the Connections box, select the adapter that you want to change the binding order on. After you select the adapter, the Bindings for Local Area Connection box displays the current binding order.

5. Use the arrows on the right side of the box to move the protocol you choose to the top of the list and then click **OK**.
6. If you are prompted to restart the computer, click **Yes**.

If PVM starts on the master node but it does not start on the slave nodes when you are using Linux, then verify that the PVM_RSH variable is set. If it has been set then verify that it is pointing to the correct RSH version.

Extra Files From an Aborted PVM Session

If you stop PVM without entering the *halt* command in the PVM console, then you might have to manually cleanup the PVM session files. This can happen if you close the PVM Console window or if the computer crashes.

If the system is running and you close the PVM console window, then you need to manually stop the pvmd daemon. On Microsoft Windows 2000 or XP, this is done in the Task Manager, whereas on Red Hat Linux or Fedora, you need to use the *ps* and *kill* commands. If you are unfamiliar with the operation of either of these tools, then consult with your System Administrator.

In the PVM temp directory (normally *c:\cluster\temp* on Microsoft Windows 2000 / XP or */tmp* on Red Hat Linux / Fedora), some files begin with *pvmd* and *pvml*. The *pvmd* file contains the process id for the active PVM session. The *pvml* file is a log file that contains status information for the current PVM session.

Note The extension for these files might vary based upon the OS that you are running. On Microsoft Windows, these files will use an extension that is based upon the account name, whereas on Linux the files will use the UID value for the account name.

❖ To delete these files (pvmd.* and pvml.*) on a Windows 2000 or XP computer

1. Open an MS/DOS console window.
2. Enter these commands:

```
c:
cd \cluster\wrshd
del pvm*.*
```

❖ To delete these files (pvmd.* and pvml.*) on Red Hat Linux or Fedora computer

Enter these commands:

```
cd /tmp
rm pvm*.*
```

Poor Cluster Performance

If it appears that the cluster is taking longer to perform searches than it was previously requiring, then check the process list in the PVM console window. Wait until no SEQUEST searches are running on the cluster. Then, in the PVM Console window enter the following command:

```
ps -ax
```

This command lists the PVM processes that are running on the cluster. If there are processes remaining for the *sequest_slave*, then you will want to delete these processes. This is done with either the *kill* or *reset* command in the pvm console window.

Cluster Searches Will Not Run Through SEQUEST Queue

If a SEQUEST Cluster search is set up to run on the SEQUEST Queue and the search is not running, then you will want to check 2 items on the master node.

The SEQUEST Queue works with the service SEQUEST Daemon. Windows services run under the SYSTEM account. The PVM software is (normally) run under the proteomics account. When a task connects to PVM, on Windows, it looks for a pvmd file where the file extension matches the account the program is being run under. Therefore, if you run a search directly from Bioworks Browser, PVM looks for the pvmd.proteomics file, whereas a search run from the SEQUEST Daemon service is looking for a pvmd.system file. If PVM can't find the appropriate file, it will exit from the task that is trying to connect to PVM. If PVM is started through the Cluster Manager program, then the pvmd.system file will be created. If you manually start PVM (from a MS-DOS console window or through some other means), you will need to manually copy the pvmd.proteomics file to pvmd.system.

If normal SEQUEST searches can be run through the SEQUEST Queue but the SEQUEST Cluster searches will not run on the SEQUEST Queue, then verify that the PVM environment variables (PVM_ROOT, PVM_ARCH, PVM_TMP, PVM_VERSION) are set as SYSTEM variables in the **Advanced > Environment Variables** page of the System Properties dialog. If the PVM values are only present in the User section, then add them to the SYSTEM section and click **OK** when you are finished. Then stop the PVM console and reboot the master node.

If a SEQUEST Cluster search is set up to run on the SEQUEST Queue and the search is not running, then you will need to check the permissions of the Cluster Temp folder. Locate the C:/Cluster/temp folder and set the security level for USERS to Full Control.

Installation Checklists

Contents

- Master Computer Checklist: Windows 2000 Server
- Master Computer Checklist: Windows 2003 Server
- Slave Computer: Microsoft Windows 2000 Checklist
- Slave Computer: Microsoft Windows XP Checklist
- Slave Computer: Red Hat Linux / Fedora Checklist

Master Computer Checklist: Windows 2000 Server

Task	Status
Install and configure Microsoft Windows 2000 Server	
Install Microsoft Windows 2000 Service Pack 4	
Configure network cards	
Install PVM	
Install and configure of RSH	
Install of Xcalibur 1.4 or 2.0	
Install of Bioworks 3.3.1	
Install and configure of SEQUEST Cluster 3.3.1	
Install of FASTA Databases	
Configure of PVM Nodes in Cluster Manager	
Distribute SEQUEST program to slave nodes	
Distribute FASTA databases to slave nodes	

Master Computer Checklist: Windows 2003 Server

Task	Status
Install and configure Microsoft Windows 2003 Server	
Install Microsoft Windows 2003 Server Service Pack 1	
Configure network cards	
Install PVM	
Install and configure of RSH	
Install of Xcalibur 1.4 or 2.0	
Install of Bioworks 3.3.1	
Install and configure of SEQUEST Cluster 3.3.1	
Install of FASTA Databases	
Configure of PVM Nodes in Cluster Manager	
Distribute SEQUEST program to slave nodes	
Distribute FASTA databases to slave nodes	

Slave Computer: Microsoft Windows 2000 Checklist

[illegible]

Slave Computer: Microsoft Windows XP Checklist

[illegible]

Slave Computer: Red Hat Linux / Fedora Checklist

[illegible]

Index

A

account name and WRSHD login are not the same [102](#)
 administrator login, configuration using [72](#)
 autologin feature is not set [102](#)

B

background information [1](#)
 batch searching [93](#)
 BioWorks, installing [16](#)

C

cluster
 background information [1](#)
 how it works [2](#)
 setting up [5](#)
 Cluster Manager [81](#)
 computer name same as login account name [102](#)
 configuring
 BioWorks
 on master node [86](#)
 on remote client computer [88](#)
 Fedora 3.0 [56](#)
 Fedora 4.0 [60](#)
 Fedora 5.0 [64](#)
 Fedora 6.0 [69](#)
 final configuration [72](#)
 Red Hat Enterprise Linux 3.0 [47](#)
 Red Hat Enterprise Linux 4.0 [51](#)
 Red Hat Linux version 8.0 [37](#)
 Red Hat Linux version 9.0 [42](#)
 using the administrator login [72](#)
 using the proteomics login [78](#)

D

databases, managing [83](#)

E

extra configuration entries in WRSHD [104](#)

extra files from an aborted PVM session [109](#)

F

FASTA Databases, installing [16](#)
 final configuration [72](#)

I

installation
 checklists [111](#)
 installing
 BioWorks [16](#)
 FASTA Databases [16](#)
 Fedora 3.0 [54](#)
 Fedora 4.0 [58](#)
 Fedora 5.0 [62](#)
 Fedora 6.0 [67](#)
 master node software [16](#)
 PVM, master node [16](#)
 PVM, slave node [25, 33](#)
 Red Hat Enterprise Linux 3.0 [45](#)
 Red Hat Enterprise Linux 4.0 [49](#)
 Red Hat Linux 8.0 [35](#)
 Red Hat Linux 9.0 [40](#)
 RSH [17](#)
 RSH, slave node [26, 33](#)
 SEQUEST Cluster software [19](#)
 slave node software [25, 32](#)
 Xcalibur [16](#)
 installing and configuring
 Windows 2000 [20](#)
 Windows 2000 server [7](#)
 Windows 2003 server [12](#)
 Windows XP [28](#)

L

Linux
 Fedora 3.0
 configuring [56](#)
 installing [54](#)

- Fedora 4.0
 - configuring 60
 - installing 58
- Fedora 5.0
 - configuring 64
 - installing 62
- Fedora 6.0
 - configuring 69
 - installing 67
- Red Hat Enterprise Linux 3.0
 - configuring 47
 - installing 45
- Red Hat Enterprise Linux 4.0
 - configuring 51
 - installing 49
- Red Hat Linux
 - setting up the slave nodes with version 9.0 58, 62, 67
- Red Hat Linux 8.0
 - configuring 37
 - installing 35
- Red Hat Linux 9.0
 - configuring 42
 - installing 40
- subsets, need additional 103

M

- managing databases 83
- master node
 - installing software 16
 - setting up 6

N

- need additional Linux subsets 103

P

- poor cluster performance 110
- processing SRF files 84
- proteomics login, configuration using 78
- PVM
 - commands 97
 - console 96
 - installing, master node 16
 - installing, slave node 25, 33
 - starting and stopping 82
 - will not start 108

R

- registry, updating 27, 35
- remote client
 - configuring BioWorks on 88

- required software 6
- RSH
 - does not work between nodes 105
 - installing, master node 17
 - installing, slave node 26, 33
 - is not working correctly 106

S

- SEQUEST
 - batch search, using 93
 - running 90
 - running from the Search Selected Samples dialog box 94
- SEQUEST Cluster Manager 81
- SEQUEST Cluster software, installing 19
- SEQUEST cluster, how it works 2
- setting up
 - cluster 5
 - master node 6
 - slave nodes with Windows 2000 20, 28
 - slave nodes, Red Hat Linux 58, 62, 67
- slave nodes
 - setting up with Windows 2000 20, 28
 - software, installing 25, 32
- SRF files, processing 84
- starting and stopping PVM software 82
- submitting SRF files for processing 84

U

- updating the Microsoft Windows registry 27, 35
- using
 - SEQUEST batch search 93
 - SEQUEST Cluster 81

W

- Windows 2000
 - Professional, installing and configuring 20
 - Server, installing and configuring 7
- Windows 2003
 - Server, installing and configuring 12
- Windows XP
 - Professional, installing and configuring 28
- WRSHD
 - cannot create service 104
 - will not start 105

X

- Xcalibur, installing 16