

Windows 7 Professional (64-bit) for MassHunter Workstation

Setup Guide

Step 1. Set up Windows 7 2
Step 2. Rename the network adapters 7
Step 3. Set the network adapter IP addresses 10
Step 4. Change firewall settings for the LC-MS instrument network adapter 12
Step 5. Confirm that Telnet Client and TFTP Client are enabled 16
Step 6. Disable the HP Support Assistant 18
Step 7. Set up Microsoft Update 22
Step 8. Install MassHunter programs 23

Before you can install MassHunter software on Windows 7 Professional, you need to set up the operating system.

Note that installations on computers not supplied by Agilent are supported on a best-effort basis only. Additional support can be purchased from the Agilent field support team.

Minimum Requirements

The minimum computer specifications to run MassHunter programs on Windows 7 Professional (64-bit) are:

- HP Z400 Xeon W3520 Quad Core Processor
- 8 GB of 1333 MHz ECC RAM
- + 4 \times 500 GB SATA HDD in a RAID 1+0 configuration (mirrored and striped)
- 80 GB C partition and a 851 GB D partition
- For MassHunter Acquisition, additional network adapter



Step 1. Set up Windows 7

The first time that Windows 7 Professional starts up after installation, you are prompted for setup information.

🕑 eji Set	Up Windows		
	Sountry or region:	ows ⁻⁷ Professional	-
	Time and currency:	inglish (United States)	-
	Keyboard layout:	15	
ø		Copyright © 2009 Microsoft Corporation	All rights reserved.

1 When prompted, select your **Country** or region, Time and currency, and **Keyboard layout**, then click **Next**.

Only these operating systems are supported:

- English (United States)
- Japanese (for MassHunter Qualitative Analysis and Quantitative Analysis programs)
- Simplified Chinese (for MassHunter Qualitative Analysis and Quantitative Analysis programs)

	Ninde	OWS' / Profession	nal
Choose a use	er name for your <u>account</u> an	d name your computer to distinguish	it on the network.
	Type a user nan	ne (for example, John):	
	user	name	
	user-PC	Lindin.	
3		Copyright © 2009 Microsoft Corpo	oration. All rights reserve
			Next
🔵 😝 Set Up V	Vindows		Next
ei set Up V	Vindows sword for your accour	A	Next
Set a pass	Vindows sword for your accour	it	Next
Set a pass Creating a p unwanted u	Vindows sword for your accour assword is a smart security p sets. Be sure to remember yo	it recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass	Vindows sword for your accour assword is a smart security p sers. Be sure to remember yo word (recommended):	It recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass	Vindows sword for your accour assword is a smart security p sets. Be sure to remember yo word (recommended):	it recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass Retype your	Vindows sword for your accour assword is a smart security p sers. Be sure to remember yo word (recommended): password:	It recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass Retype your	Vindows sword for your accour assword is a smart security p sers. Be sure to remember yo word (recommended): password:	it recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass Retype your Type a pass	Vindows sword for your accour assword is a smart security p sers. Be sure to remember yo word (recommended): password: word hint:	it recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass Retype your Type a pass	Vindows sword for your accour assword is a smart security p sers. Be sure to remember yo word (recommended): password: word hint:	it recaution that helps protect your use ur password or keep it in a safe place	r account from
Set a pass Creating a p unwanted u Type a pass Retype your Type a pass Choose a wn B you forgel	Vindows sword for your accour assword is a smart security p sers. Be sure to remember yo word (recommended): password: word hint: ord or phrase that helps you your password, Windows w	It recaution that helps protect your use ur password or keep it in a safe place	r account from

2 Create a user name and password for a user that will be a member of the Administrators group.

For Agilent bundle computers, create a new user other than "admin." The user name "admin" with password "3000hanover" is already created when the Emergency Recovery Disk is used.

Step 1. Set up Windows 7



 Read Microsoft's license terms for Windows 7 Professional. Select
 I accept the license terms and click Next.

4 Click **Ask me later** when prompted to turn on Microsoft Updates, then click **Next**.

If you choose to turn on Microsoft Updates, be aware that an update can automatically restart your computer, which could cause MassHunter programs to fail.



Agilent Technologies DataSystem End User License Agreement	
Welcome to your Agilent Technologies DataSystem You must agree to the end user license agreement before using the software on this computer	
Your HP Workstation contains preinstalled software programs. Please read the Software License Agreement before proceeding.	î
CAREFULLY READ THIS LICENSE AGREEMENT AND LIMITED WARRANTY STATEMENT BEFORE PROCEEDING TO OPERAIE THIS EQUIPMENT. RIGHTS IN THE SOFTWARE ARE OFFERED ONLY ON THE CONDITION THAT THE CUSTOMER AGREES TO ALL TERMS AND CONDITIONS OF THE LICENSE AGREEMENT. PROCEEDING TO OPERATE THE EQUIPMENT INDICATES YOUR ACCEPTANCE OF THESE TERMS AND CONDITIONS. IF YOU DO NOT AGREE WITH THE TERMS OF THE LICENSE AGREEMENT, YOU MUST NOW EITHER REMOVE THE SOFTWARE FROM YOUR HARD DISK DRIVE AND DESTROY THE MASTER DISKETTES, OR RETURN THE COMPLETE COMPUTER AND SOFTWARE FOR A FULL REFUND.	
* announcements and *	-
LAgreeCancel	

5 Set up the time and time zone for the computer.

- 6 If you are prompted to accept the End User License Agreement, click I Agree.
- 7 If you prompted to reboot the computer, ignore this message. Do not restart the computer.

Step 1. Set up Windows 7



- 8 Right-click the **Computer** icon on the Desktop and select **Properties**.
- **9** Confirm that your computer meets minimum requirements:
 - Windows 7 Professional
 - 8 GB of RAM
 - 64-bit system type

If the incorrect operating system is installed, get media and a license for Windows 7 Professional (64-bit) and reinstall the operating system.

- **10** Restart the computer.
- **11** After you restart the computer, if the screen resolution is not properly reset, then:
 - a Right-click the Desktop and click **Screen resolution**.
 - **b** Change the resolution to the correct resolution for the monitor.

The MassHunter Data Acquisition, Qualitative Analysis, and Quantitative programs require a minimum screen resolution of 1280 × 1024.

For efficient data processing, set the resolution to 1680×1050 or higher. Agilent LC/MS instruments ship with a minimum monitor resolution of 1680×1050 .

For maximum picture clarity, operate LCD monitors at their native resolution. Lower than native resolution causes blurred images.

Step 2. Rename the network adapters

- 1 Select Start > Control Panel.
- 2 From the Control Panel, select Network and Internet > Network and Sharing Center > View network status and tasks.
- **3** From the left menu, select **Change adapter settings**.

S S S S S S S S S S S S S S S S S S S	ernet > Network and Sharing Center - 4 Search Control Panel
Control Panel Home Change adapter settings	View your basic network information and set up connections •••••••••••••••••••••••••••••
settings	user-PC Internet (This computer)
	View your active networks Connect to a network You are currently not connected to any networks.
	Change your networking settings
	Set up a new connection or network Set up a wireless, broadband, dial-up, ad hoc, or VPN connection; or set up a router or access point.
	Connect to a network Connect or reconnect to a wireless, wired, dial-up, or VPN network connection.
	Choose homegroup and sharing options Access files and printers located on other network computers, or change sharing settings.
	Troubleshoot problems Diagnose and repair network problems, or get troubleshooting information.
See also	
Internet Options	
Windows Firewall	

You will see two network connections, Local Area Connection and Local Area Connection 2. Neither are connected at this time.

Step 2. Rename the network adapters



- **4** Identify the onboard network adapter:
 - **a** Connect a network cable to the onboard network adapter on the back of the computer.
 - **b** Note which Local Area Connection is enabled.

Typically, the onboard network adapter is Local Area Connection.

- **5** Right-click the Local Area Connection for the onboard network adapter and select **Rename**. Change the name to HOUSE connection.
- 6 Right-click the other Local Area Connection icon and select **Rename**. Change the name to LC-MS Instrument.

Step 2. Rename the network adapters

00-	(a) // Natural and Internal. A Matural Connections. A		
Organize	Wetwork and internet * Network connections * Search Network connections *	-	0
	HOUSE Connection Network Broadcom NetXtreme Gigabit Eth KC-MS Instrument Network cable unplugged Broadcom NetXtreme Gigabit Eth		

aworking	Sharing		
Connect us	sing:		
🔮 Broa	dcom NetXtren	ne Gigabit Etherne	t #2
			Configure
his c <u>o</u> nne	ction uses the	following items:	
✓ Image: Fill ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → ✓ → Insta	le and Printer S roadcom Advar ternet Protocol ternet Protocol nk-Layer Topol nk-Layer Topol all	haring for Microso need Server Progriv Version 6 (TCP/If Version 4 (TCP/If ogy Discovery Ma ogy Discovery Re	ft Networks am Driver 2v6) 2v4) pper I/O Driver sponder P <u>r</u> operties
Descriptio Transmis wide are across d	on ssion Control Pr a network prote iverse intercon	rotocol/Internet Pr ocol that provides nected networks.	otocol. The default communication
		_	

Step 3. Set the network adapter IP addresses

- General Alternate Configuration You can get IP settings assigned automatically if your network supports this capability. Otherwise, you need to ask your network administrator for the appropriate IP settings. Obtain an IP address automatically Use the following IP address: — IP address: Subnet mask: Default gateway: Obtain DNS server address automatically Use the following DNS server addresses: Preferred DNS server: Alternate DNS server: Validate settings upon exit Advanced... OK Cancel
- **3** In the General tab:

Properties.

(TCP/IPv4).

a Make sure **Obtain an IP address** automatically is selected.

1 Right-click **HOUSE connection** and select

2 Double-click Internet Protocol Version 4

b Click **OK** twice to return to the Network Connections dialog box.

LC-MS1	Instrument Properties
Networking	g Sharing
Connect	
Connect	
E Br	oadcom NetXtreme Gigabit Ethemet
	Configure
This con	nection uses the following items:
	Client for Microsoft Networks
🗹 🔒	QoS Packet Scheduler
🗹 📙	File and Printer Sharing for Microsoft Networks
M	Broadcom Advanced Server Program Driver
	Internet Protocol Version 6 (TCP/IPv6)
	Internet Protocol Version 4 (TCP/IPv4)
	Link-Layer Topology Discovery Mapper I/O Driver
	Link-Layer Topology Discovery Nesponder
l <u>n</u> :	stall Uninstall Properties
Descrip	ption
Transr	mission Control Protocol/Internet Protocol. The default
wide a	area network protocol that provides communication
001000	
	OK Cancel
tarm at Day	
ternet Pro	stocol version 4 (TCP/IPv4) Properties
General	
Veuene	act ID actions preisted submatically if your patwork support
this capa	ability. Otherwise, you need to ask your network administrato
for the a	ppropriate IP settings.
-	
Obt Obt	tain an IP address automatically

Use the following IP address:

Obtain DNS server address automatically
 Use the following DNS server addresses:

IP address:

Subnet mask:

Default gateway:

Preferred DNS server:

Alternate DNS server:

Validate settings upon exit

- 4 Right-click LC-MS Instrument and select Properties.
- 5 Double-click Internet Protocol Version 4 (TCP/IPv4).

- 6 In the General tab:
 - a Select Use the following IP address.
 - **b** For **IP address**, type 192.168.254.1.

The Subnet mask automatically populates as 255.255.255.0. Leave the Default Gateway and DNS Server lines empty.

c Click **OK** twice to return to the Network Connections panel.

192.168.254.1

255.255.255.0

.

OK

Advanced...

Cancel

. . .

. .

Step 4. Change firewall settings for the LC-MS instrument network adapter

- **1** Return to the Control Panel view.
- 2 Click System and Security > Windows Firewall.
- 3 From the left menu, click Advanced settings

😋 🕞 🗢 🕍 🕨 Control Panel 🕨	System and Security Windows Firewal	I 👻 🍫 Sea	rch Control Panel	Q
Control Panel Home Allow a program or feature through Windows Firewall Change notification settings Turn Windows Firewall on or off Restore defaults Advanced settings Troubleshoot my network	Help protect your computer v Windows Firewall can help prevent has through the Internet or a network. How does a firewall help protect my co What are network locations? What are network locations? Dependence of the protect my co What are network locations? Dependence of the protect my co Windows Firewall state: Incoming connections: Active public networks: Notification state:	vith Windows Firewa ckers or malicious software omputer? te) networks oorts or coffee shops On Block all comm the list of allo None Notify me wh program	II from gaining access Not C C nections to programs wed programs	to your computer
See also Action Center				
Network and Sharing Center				



4 Under Overview, click Windows Firewall Properties.

5 In the Domain Profile tab, next to Protected network connections, click Customize.

Windows Firewa	all with Advanced Sec	urity on Loca	Computer Pr
Domain Profile	Private Profile Public	Profile IPse	c Settings
Specify beh domain. State	avior for when a comput	er is connecte	d to its corporate
	rewall state: On (reco		nded) 🔻
	Inbound connections:	Bloc	k (default) ▼
	Outbound connection	s: Allo	w (default) 🔹
	Protected network co	nnections:	Customize
Settings	Specify settings that con Firewall behavior.	rol Windows	Customize
Logging t	Specify logging settings f roubleshooting.	or	Customize
Learn more a	about these settings		ancel Anoly

6 Clear the LC-MS Instrument check box, and then click OK.

Protected Network Connections for the Do	omain Profile 📃 🗖 💌
Select the check box for each connection ye protect.	ou want Windows Firewall to help
Network connections:	
HOUSE Connection	
Learn more about these settings	
	OK Cancel

7 Repeat the step 5 and step 6 for the **Private Profile** and **Public Profile** tabs.

Protected Network Connections for the Private Profile
Select the check box for each connection you want Windows Firewall to help protect.
Network connections:
HOUSE Connection LC-MS Instrument
Leam more about these settings OK Cancel
Protected Network Connections for the Public Profile
Select the check box for each connection you want Windows Firewall to help protect.
Network connections:
HOUSE Connection LC-MS Instrument
Leam more about these settings OK Cancel

- 8 Click **OK** to close the Windows Firewall with Advanced Security on Local Computer Properties dialog box.
- **9** Close the Windows Firewall with Advanced Security window.

Step 5. Confirm that Telnet Client and TFTP Client are enabled

Step 5. Confirm that Telnet Client and TFTP Client are enabled

1 Return to the Control Panel view and click **Programs.**



- 2 Under Programs and Features, click Turn Windows features on or off.
- **3** Verify that the **Telnet Client** and **TFTP Client** check boxes are selected.

These clients are required for the LC/MS Firmware Update Tool and for GC/MS instruments.

Windows Features	x
Turn Windows features on or off	0
To turn a feature on, select its check box. To turn a feature off, clear it: check box. A filled box means that only part of the feature is turned or	5 1.
🗉 💷 Microsoft Message Queue (MSMQ) Server	-
🕀 🔳 🕒 Print and Document Services	
Remote Differential Compression	
RIP Listener	
🗉 💷 🔐 Simple Network Management Protocol (SNMP)	
Simple TCPIP services (i.e. echo, daytime etc)	
Tablet PC Components	=
V Telnet Client	
Telnet Server	
TFTP Client	
Windows Gadget Platform	
🕞 🥅 🖟 Windows Process Activation Service	Ŧ
OK Cance	!

4 Press OK and close the Control Panel.

Step 6. Disable the HP Support Assistant

Step 6. Disable the HP Support Assistant

If your computer is a Hewlett-Packard computer, you must disable the HP Support Assistant so that it does not run scans while MassHunter programs are running.

- 1 Click Start > All Programs > HP > HP Support Assistant.
- 2 Click Cancel.





3 Click the **Settings** icon in the lower left corner. Make sure you click the icon and *not* the **Settings** label.

Back			Settings		
Health Analysis Tune up	General				
Select an analysis schedule:					
Frequency:	Day of the week:	Approxima	te time:		
Never					
Please select how you woul	d like to receive upd	ates from HP:			
Install important updates aut	o <i>matically</i> d updates				
Check for updates but let me					
Never check for updates					
HP Privacy Statement	view information coll	ected by HP			
				Close	

4 In the Health Analysis tab, change Frequency to Never.

= Documentaria - word	Pau		
	HP Support Assistant		- 🗆 X
Deals	Catting		
Dack	Jettings		
Health Analysis Iune	up General		
Select tacks performed (during a tune-up		
Set Restore Point			
Empty Recycle bin			
Clear Internet Explo	rer's temporary Internet files		
Clear Temporary fol	lder -		
V Microsoft Windows	Disk Defragment		
Select a tune-up schedu	ile:		
Frequency:	Day of the month: Approximate time:		
Never			
		Close	Save
Product Nam	e Loading Product Number Loading		
Settings Serial Number	er Loading Warranty		

5 In the **Tune up** tab, change **Frequency** to **Never**.

6 Click Save and then close the program.

Step 7. Set up Microsoft Update

Agilent recommends that Microsoft Update is turned on if possible. However, an update can cause your computer to restart unexpectedly during data acquisition and cause you to lose data.

Do this step to set up Windows 7 so that it updates your computer only with user permission.

- 1 In the Control Panel, click Systems and Security > Windows Update > Change settings.
- 2 Under Important updates, select Download updates but let me choose whether to install them.
- 3 Click OK and close the Control Panel.

When the Control Panel is closed, if the computer is connected to the Internet, Windows immediately checks for updates. The list of suggested updates can be large.

	of sectors.	- • x
Control Panel > System and Security > Windows Update > Change settings	Search Control Panel	٩
Choose how Windows can install updates When your computer is online, Windows can automatically check for important updates and install them using these settings. When new updates are available, you can also install them before shutting down the computer. How does automatic updating help me? Important updates Download updates but let me choose whether to install them Install new updates: Every day at 3:00 AM Choose me recommended updates the same way I receive important updates Who can install updates Who can install updates Mode all updates to install updates on this computer Note Windows Update might update itself automatically first when checking for other updates. Read our privacy statement online.		
🛞 OK Cancel		

Step 8. Install MassHunter programs

• Install MassHunter programs.

Follow the appropriate installation or upgrade guide. Note that on Windows 7, some MassHunter programs require that you run the installation program as administrator. The installation guide tells you which steps require that you right-click the installation command and select **Run as administrator**.

www.agilent.com

In This Book

This guide describes how to set up Windows 7 Professional (64-bit) for MassHunter Workstation, to prepare for MassHunter software installation.

Windows[®] is a registered trademark of Microsoft Corporation in the United States and other countries.

 ${}^{\textcircled{C}}$ Agilent Technologies, Inc. 2011

Revision A, April 2011

G3336-90016

