

User Bulletin

Applied Biosystems 3130/3130x/ Genetic Analyzers and
Applied Biosystems 3730/3730x/ DNA Analyzers

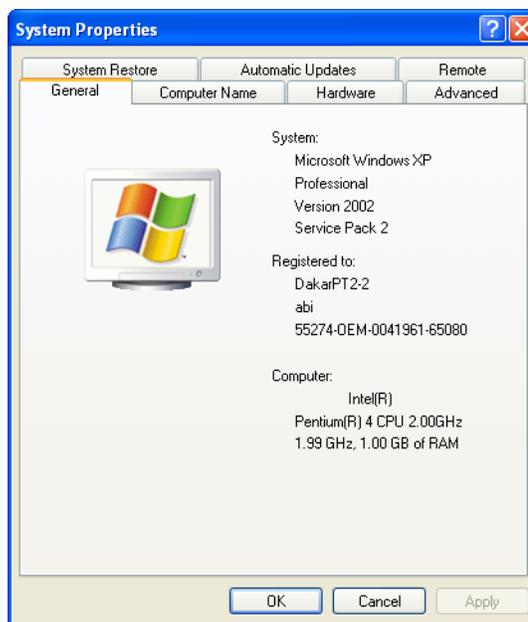
February 16, 2005

SUBJECT: Data Collection Software v3.0 and Windows® XP Service Pack 2

In This User Bulletin	This user bulletin includes the following topics:
	Configuring the Windows Security 2 Monitoring Window Security Alerts 6

Purpose After you have upgraded your Windows® XP operating system to Service Pack 2, you must configure Windows security to accommodate Applied Biosystems Data Collection Software v3.0. The purpose of this user bulletin is to provide instructions for configuring security settings to ensure that Data Collection Software v3.0 functions properly with your Windows XP Service Pack 2 operating system.

Assumptions This user bulletin assumes that you have successfully upgraded your operating system to Service Pack 2. To verify that the upgrade was successful, select **My Computer > Properties** to open the System Properties window.



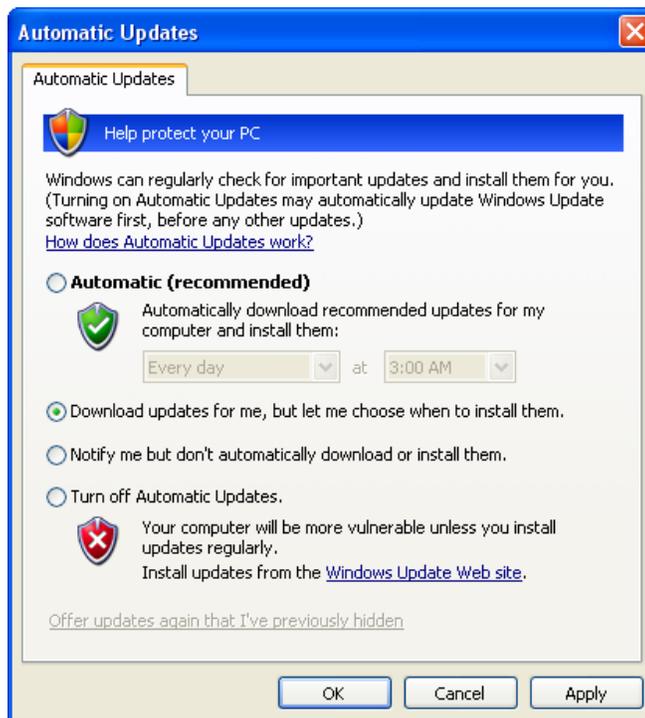
Configuring Windows Security

Configuring the Windows Security Center

1. Select **Start > Settings > Control Panel** and click **Security Center** to open the Windows Security Center window.

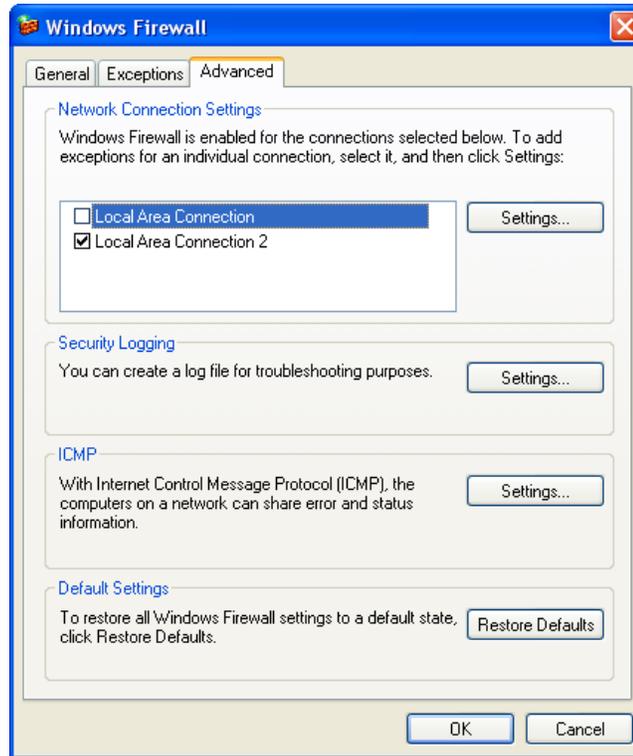


2. In the Security Center window, click **Automatic Updates** to set security for updates.
3. Select **Download updates for me, but let me choose when to install them**, then click **OK** to return to the Security Center.



Note: Selecting Automatic (recommended) may cause the computer to unexpectedly reboot when updates are automatically downloaded and installed. Run data could be lost when the computer reboots.

4. In the Security Center window, click **Windows Firewall** and select the **Advanced** tab.



5. In the **Network Connection Settings** pane, deselect the network connection that communicates between the instrument and computer.

Note: To determine which network connection is dedicated for private communication between the instrument and computer, type `ipconfig` from a command prompt. The network connection with a static IP Address of **192.168.0.1** is the one which should be disabled. If the firewall is enabled for the private network connection, the instrument will fail to boot and the status lights are solid yellow. Check that the external network connection is selected to ensure the firewall is enabled for the external network.

```

C:\WINDOWS\System32\cmd.exe
Microsoft Windows XP [Version 5.1.2600.1]
(C) Copyright 1985-2001 Microsoft Corp.

C:\Documents and Settings\Administrator>ipconfig

Windows IP Configuration

Ethernet adapter Local Area Connection:

    Connection-specific DNS Suffix  . : 
    IP Address . . . . . : 192.168.0.1
    Subnet Mask . . . . . : 255.255.255.0
    Default Gateway . . . . . : 

Ethernet adapter Local Area Connection 2:

    Connection-specific DNS Suffix  . : pe-c.com
    IP Address . . . . . : 167.116.114.98
    Subnet Mask . . . . . : 255.255.255.0
    Default Gateway . . . . . : 167.116.114.1

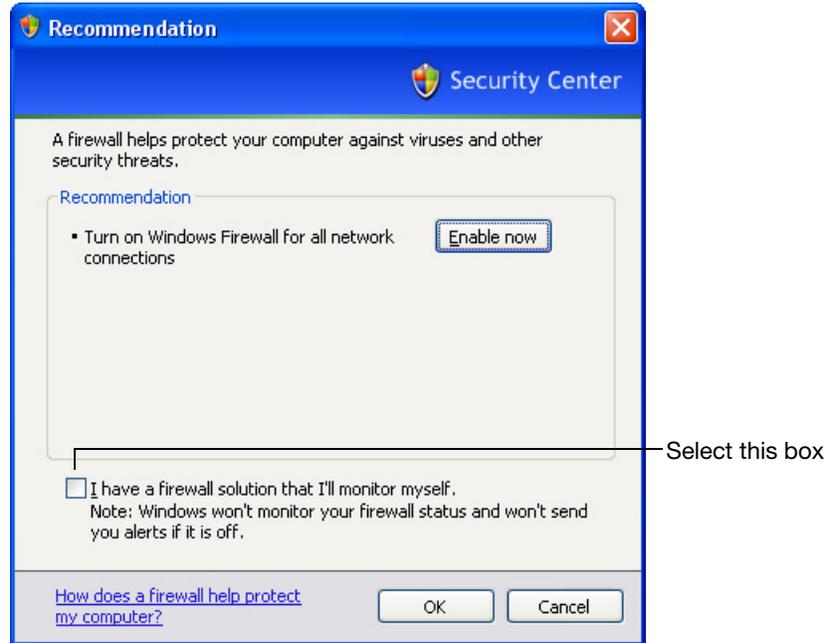
C:\Documents and Settings\Administrator>
    
```

Click **OK** to return to the Security Center window.

6. (Optional) You may disable the warning **Your computer might be at risk** when the Windows Firewall private network connection is disabled. To disable the warning:
 - a. In the Security Center under Firewall, click **Recommendations**.



- b. Select **I have a firewall solution that I'll monitor myself** then click **OK** to return to the Security Center.



7. Click **OK** to close the Windows Security Center.

Windows XP Service Pack 2 has been successfully configured for Data Collection Software v3.0.

Monitoring Window Security Alerts

After you have installed the Windows XP service pack 2 (SP2) upgrade and you launch any of the applications below, the Windows operating system reports Windows Security Alerts such as the following:

Data Collection Software v3.0



Click **Unblock** to unblock the following Data Collection Software v3.0 alerts:

- RunJavaW
- FMQServer
- Java
- Daemon3730
- Viewer3730

Note: The name for the Daemon and Viewer alert messages may vary depending on your instrument model.

Autoanalysis Manager v3.0



GeneMapper® Software v3.7 and GeneMapper® ID Software v3.2



Sequencing Analysis Software v5.2 and SeqScape® Software v2.5



Click **Unblock** to dismiss these alerts and continue using the software.

Note: Go to **Windows Security Center > Windows Firewall > Exceptions** tab to locate the unblocked security alerts.

© Copyright 2005, Applied Biosystems. All rights reserved.

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

Information in this document is subject to change without notice. Applied Biosystems assumes no responsibility for any errors that may appear in this document. This document is believed to be complete and accurate at the time of publication. In no event shall Applied Biosystems be liable for incidental, special, multiple, or consequential damages in connection with or arising from the use of this document.

Notice to Purchaser: License Disclaimer

Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under patent claims owned or otherwise controlled by Applied Biosystems, either expressly or by estoppel.

TRADEMARKS:

Applied Biosystems, GeneMapper, and SeqScape are registered trademarks and AB (Design), and Applied Biosystems are trademarks of Applied Biosystems Corporation or its subsidiaries in the U.S. and/or certain other countries.

Microsoft, Windows, and Windows XP are registered trademarks of the Microsoft Corporation in the United States and other countries.

All other trademarks are the sole property of their respective owners.



iScience. To better understand the complex interaction of biological systems, life scientists are developing revolutionary approaches to discovery that unite technology, informatics, and traditional laboratory research. In partnership with our customers, Applied Biosystems provides the innovative products, services, and knowledge resources that make this new, **Integrated Science** possible.

Worldwide Sales and Support

Applied Biosystems vast distribution and service network, composed of highly trained support and applications personnel, reaches 150 countries on six continents. For sales office locations and technical support, please call our local office or refer to our Web site at www.appliedbiosystems.com.

Applied Biosystems is committed to providing the world's leading technology and information for life scientists. Applied Biosystems Corporation consists of the Applied Biosystems and Celera Genomics businesses.

Headquarters

850 Lincoln Centre Drive
Foster City, CA 94404 USA
Phone: +1 650.638.5800
Toll Free (In North America): +1 800.345.5224
Fax: +1 650.638.5884

02/2005