

SoftMax® Pro Software Version 5.4.1 (and later) Installation Addendum

This addendum provides information about installing SoftMax® Pro Software version 5.4.1 or later on a system with a Windows 7 operating system.

Support for the Windows 7 Operating System

SoftMax Pro Software v5.4.1 and later support installation on Windows 7, 32-bit and 64-bit systems.

Installing SoftMax Pro Software

To install SoftMax Pro Software, follow these steps:

1. Insert the SoftMax Pro CD into the CD-ROM drive.
2. **Windows:** Open My Computer and then double-click the CD-ROM drive icon to launch the installer.
Mac: Double-click the CD icon to launch the installer.
3. Double-click the SoftMax Pro Software Installer icon and follow the screen prompts.

During the installation, Windows 7 User Account Control (UAC) requires user confirmation in two places in order to complete the installation process. In the event that the installation is performed by a non-administrator, Windows 7 will also require administrator's authorization.



Note: The User Account Control warning window often appears behind the SoftMax Pro Software's installation window and you need to acknowledge it in order to complete the installation.

Using MDC FileServer

Due to increased security features present in Windows 7, to enable TCP/IP access to the user accounts file for SoftMax Pro GxP software, the following configuration steps are required.

Creating a New Inbound Rule

In the **New Inbound Rule Wizard** dialog box:

1. Select **Protocols and Ports**.
2. Click **ICP**.
3. Click **Specify local ports**.
4. Type **9001** in the entry field.
5. Click **Next** to continue the configuration.

Creating a New Outbound Rule

In the **New Outbound Rule Wizard** dialog box:

1. Select **Protocols and Ports**.
2. Click **ICP**.
3. Click **Specify local ports**.
4. Type **9001** in the entry field.
5. Click **Next** to continue the configuration.

Driver Installation

Due to elevated security of a Windows 7 64-bit operating system, the USB instrument drivers are not installed automatically. The instruments affected are the SpectraMax® L Microplate Reader and the StakMax® Microplate Handling System.

When connecting for the first time to either the SpectraMax L reader or StakMax system. Perform the following steps to install the drivers:

1. To access the **Device Manager** select **Start > Control Panel > System**.
2. Double-click the unknown device with the yellow warning icon.
3. Select the **Driver** tab and click **Update Driver**.
4. Select **Browse my computer**.
5. Select the SoftMax Pro Software folder's installation path.
6. On the **Windows Security warning** screen, select **Install this driver software anyway**.