

Sentinel Protection Server v7.4.2 Installation

Overview

A clean Install of MIS v7.5 from the MIS v7.5 CD will install the Sentinel Protection Server v7.4.0 for the memory key dongle. There is a known security issue with this version of the Sentinel driver. SafeNet's website describes the issue as follows:

ARTICLE #
T20069

PRODUCTS
MetaMorph®
MetaFluor®
MetaVue®

CREATED
02-May-2008

LAST UPDATED
06-Nov-2009

Description: Sentinel Protection Installer 7.4.0 [Security Patch]

Content: Web Vulnerability Scanner scans indicate that our Sentinel Server is prone to Directory Traversal Attack. If you paste the link below, you will be able to download the files from any system that is running the Sentinel Keys Server or Sentinel Protection Server. Below is the link for Sentinel Keys Server.

<http://192.168.0.1:7002/../../../../../../../../Windows/notepad.exe>

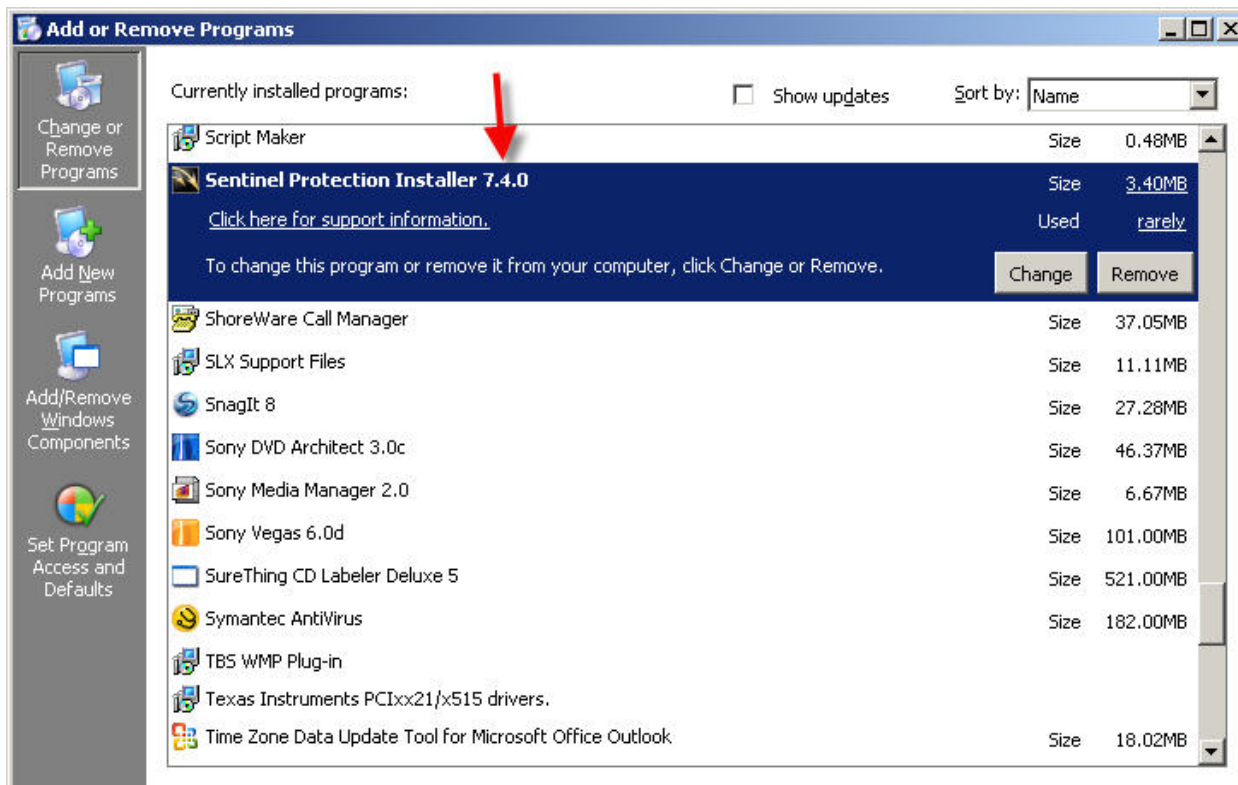
The link below gives access to the root. Now, if anyone knows where the file is located, will be able to get the file, even if he is not authorized.

<http://192.168.0.1:7002/../../../../../../../../>

This security issue is for **v7.4.0 only**, as earlier versions of the driver are not effected by this vulnerability, and the latest version of the Sentinel Protection Server (v7.4.2) fixes this security issue. If you upgraded from an earlier MIS v7.x, chances are the installed Sentinel Protection Server from that version was not changed.

How to check which Sentinel Protection Server version you have installed

1. Open the Add Remove Programs and scroll down until you see the Sentinel Protection Server listed



2. If the version is listed as v7.4.0, you will need to update the Sentinel Protection Installer to fix the security issue.

Clean Installation of Sentinel Protection Server 7.4.2

Files required

The files below are Zip archives. Please download them and unpack each. The result should be an application named *SSDCleanup* and an application named *Sentinel Protection Installer 7.4.2*.

32-Bit OS (Windows 2000, XP, Vista)

- *SSDCleanup Zip archive:*
<ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel%20Drivers/SSDCleanup-Win32/SSDCleanup.exe>
- *Installer for Sentinel System Driver 7.4.2:*
<ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel Drivers/Sentinel Protection Installer 7.4.2.zip>

64-Bit OS (XP x64, Vista 64-bit)

- *SSDCleanup Zip archive:*

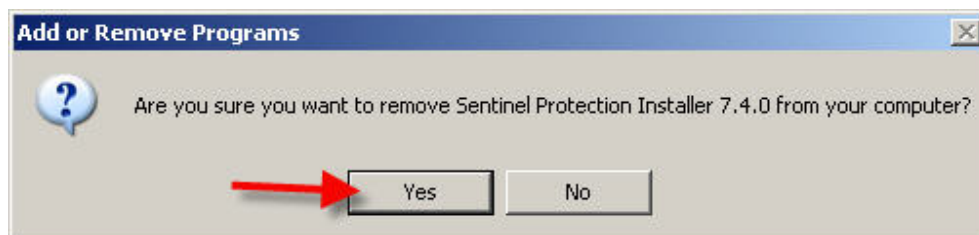
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

[ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel Drivers/SSDCleanup-Win64/SSDCleanupx64.exe](ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel%20Drivers/SSDCleanup-Win64/SSDCleanupx64.exe)

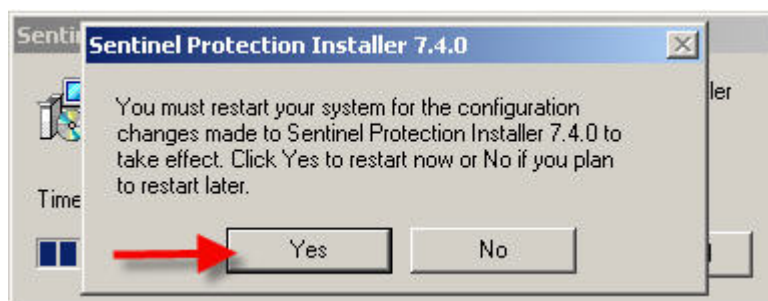
- o *Installer for Sentinel System Driver 7.4.2:*
[ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel Drivers/Sentinel Protection Installer_7.4.2.zip](ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel%20Drivers/Sentinel%20Protection%20Installer_7.4.2.zip)

Procedure

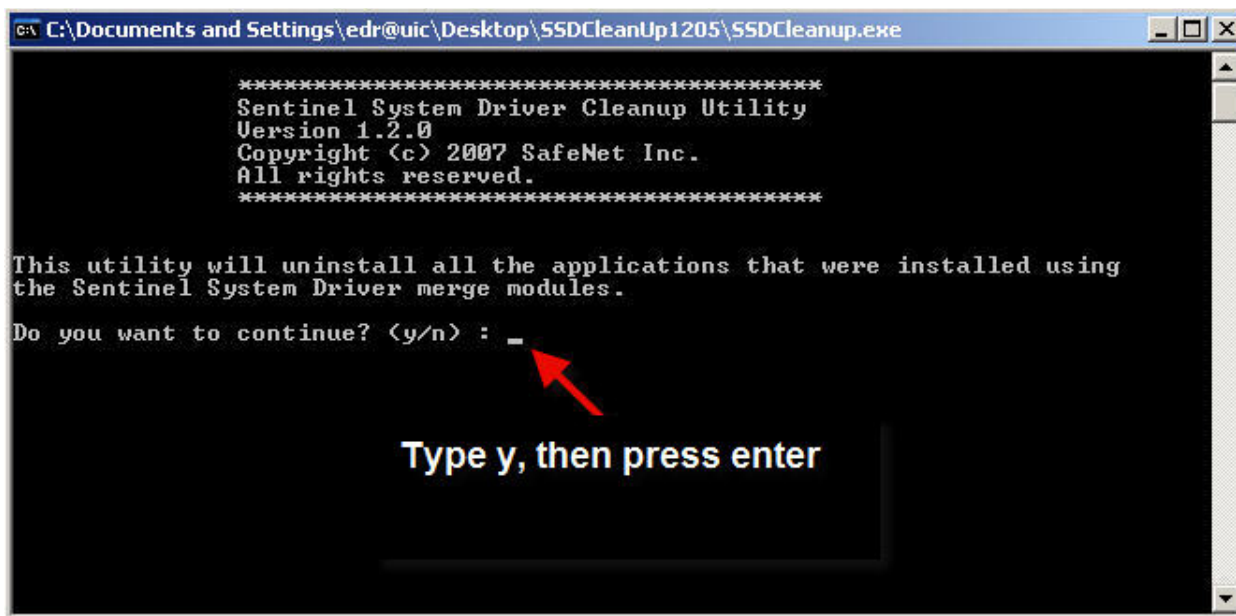
1. If using a USB Memory Key, unplug the Memory Key from the USB port.
2. Uninstall the existing driver using the *Add or Remove Programs* control panel by following steps 2a-2d.
 - a. If the control panel's list of currently installed programs includes a *Sentinel System Driver*, the computer has Version 7.3.0 or earlier of the driver is installed; You do not need up update the driver.
 - b. If the control panel's list of currently installed programs includes a *Sentinel Protection Installer*, the computer has Version 7.4.0 of the driver installed; remove it.
 - c. Click Yes when asked to Remove Sentinel Protection Installer 7.4.0



- d. Click Yes to Reboot system.



3. Run *SSDCleanup*. Restart the computer if told to do so.



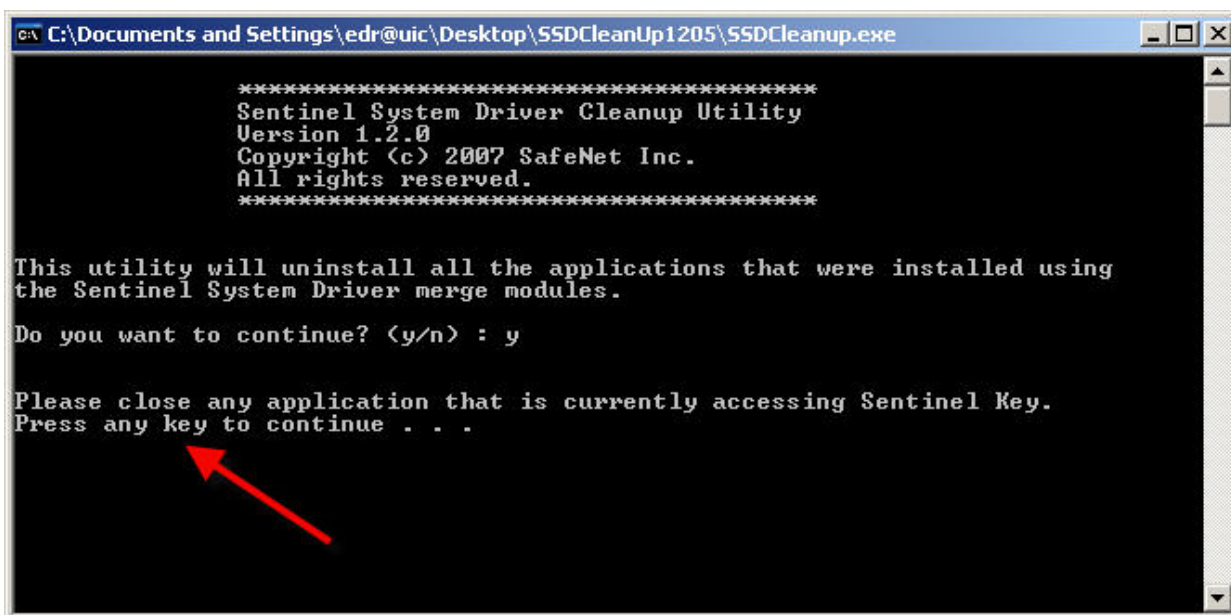
```
C:\Documents and Settings\edr@uic\Desktop\SSDCleanUp1205\SSDCleanup.exe

*****
Sentinel System Driver Cleanup Utility
Version 1.2.0
Copyright (c) 2007 SafeNet Inc.
All rights reserved.
*****

This utility will uninstall all the applications that were installed using
the Sentinel System Driver merge modules.

Do you want to continue? (y/n) : _

Type y, then press enter
```



```
C:\Documents and Settings\edr@uic\Desktop\SSDCleanUp1205\SSDCleanup.exe

*****
Sentinel System Driver Cleanup Utility
Version 1.2.0
Copyright (c) 2007 SafeNet Inc.
All rights reserved.
*****

This utility will uninstall all the applications that were installed using
the Sentinel System Driver merge modules.

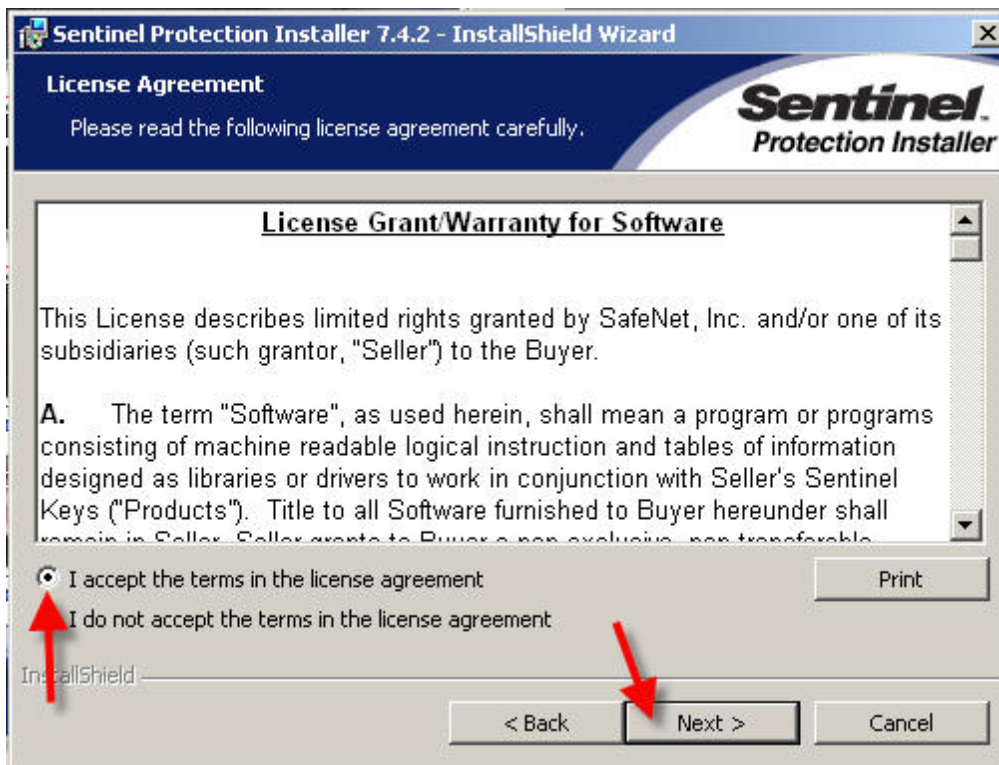
Do you want to continue? (y/n) : y

Please close any application that is currently accessing Sentinel Key.
Press any key to continue . . .
```

4. Install Sentinel Protection Server 7.4.2 using the *Sentinel Protection Installer 7.4.2* application.



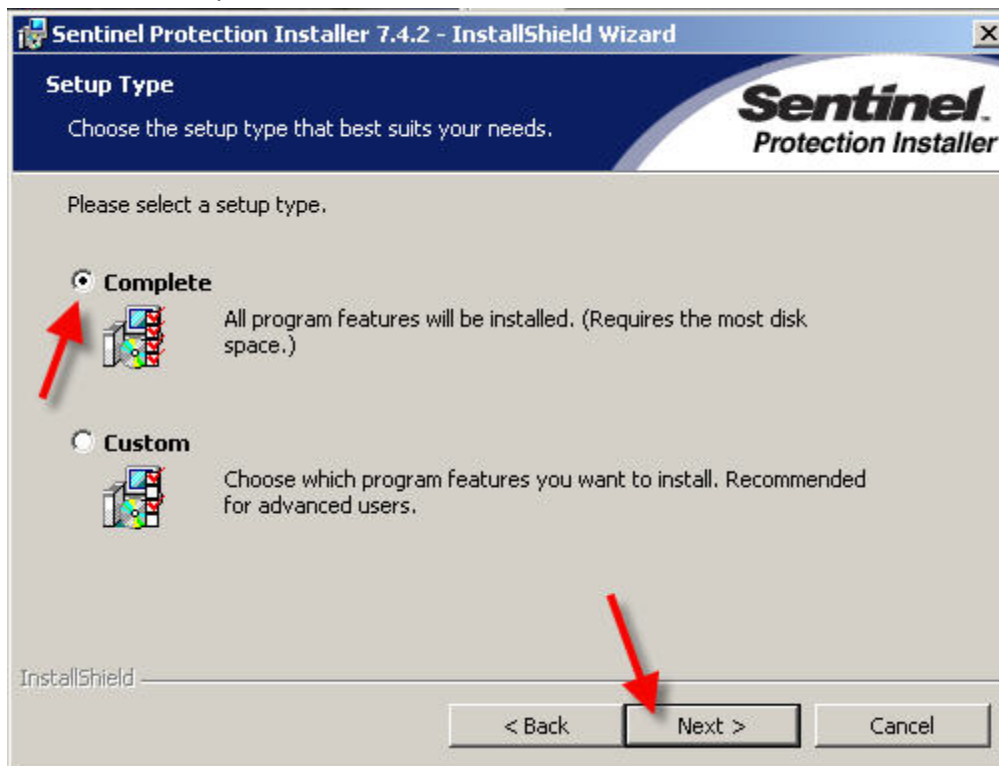
a. Click Next



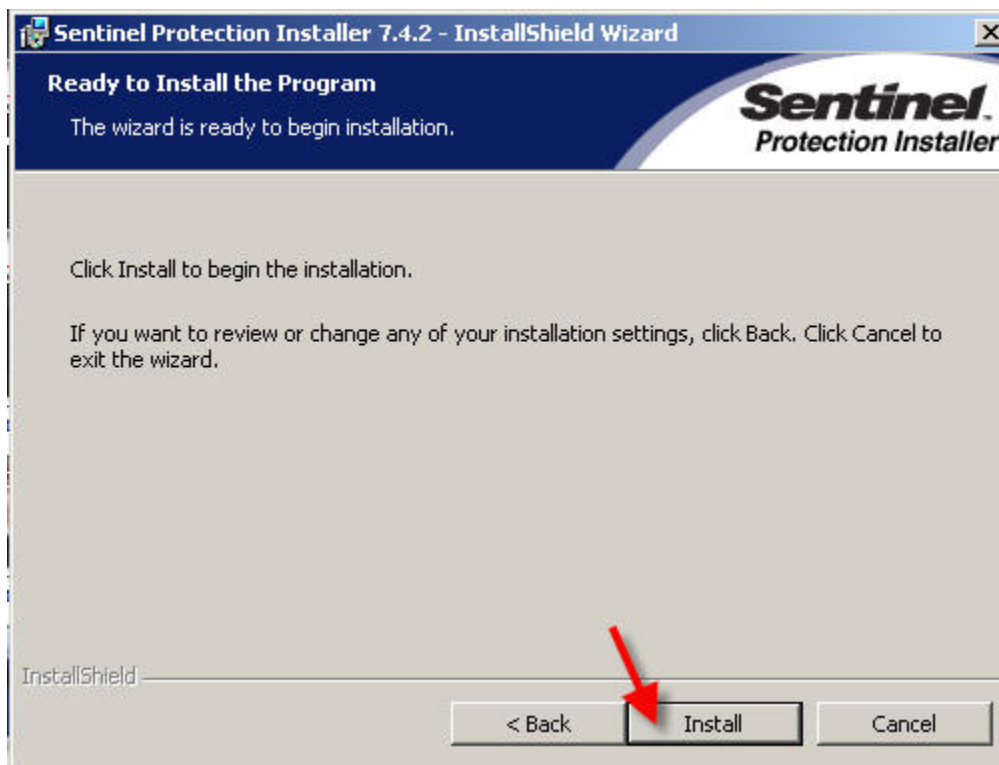
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES.

MOLECULAR DEVICES, the Molecular Devices logo, METAMORPH, META IMAGING SERIES, METAFUOR and the MetaMorph logo are trademarks of Molecular Devices, Inc. ©2010 Molecular Devices, Inc. Printed in the U.S.

b. Select "I accept..." then click Next



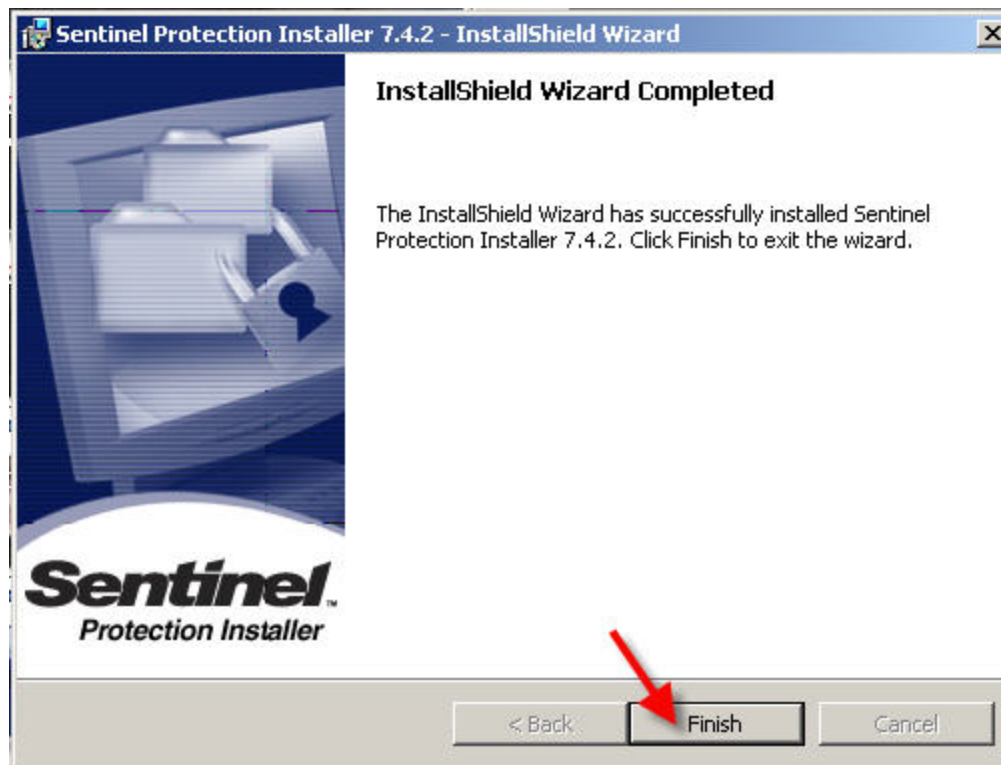
c. Select Complete, then click Next.



d. Click Install



- e. If installing a Network License memory key that other offline seats will be connecting to this system, click Yes. If installing a local single license memory key, click No.



- f. Click Finish to close the installer

Verify installation

1. Start the application *C:\Program Files\Common Files\SafeNet Sentinel\Sentinel System Driver\SetupSysDriver*.
2. Click on the *Configure Driver* button. This opens the *Sentinel Driver* dialog, with one line for the USB ports and one line for each LPT port that the Sentinel Driver recognizes.
3. Choose *OK*. This closes the *Sentinel Driver* dialog.
4. Exit the *Sentinel System Driver* application.
5. Attach the memory key and Launch MetaMorph, MetaFluor, or MetaVue program.

If there are any further questions, please contact Technical support at 800-635-5577 x1820 or support.dtn@moldev.com.