

## New Installation of Sentinel System Driver 7.3.0

**Note:** This procedure assumes that you do not have a version of the sentinel drivers currently installed.

### Files required

The files below are Zip archives. Please download them and unpack each. The result should be an application named *SSDCleanup11* and an application named *Sentinel Protection Installer 7.3.0*.

- *Installer for Sentinel System Driver 7.3.0:*  
[ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel%20Drivers/Sentinel Protection Installer 7.3.0.zip](ftp://ftp.meta.moleculardevices.com/pub/uic/software/Sentinel%20Drivers/Sentinel%20Protection%20Installer%207.3.0.zip)

**ARTICLE #**  
T20079

**PRODUCTS**  
MetaMorph®  
MetaFluor®  
MetaVue®

**CREATED**  
29-April-2008

**LAST UPDATED**  
29-November-2009

### Procedure

1. If using a USB Memory Key, unplug the Memory Key from the USB port.  
Steps 2 -3 deleted for a new installation
4. Install SSD 7.3.0 using the *Sentinel Protection Installer 7.3.0.exe* application.
5. Start the application *C:\Program Files\Common Files\SafeNet Sentinel\Sentinel System Driver\SetupSysDriver*.
6. 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.
7. If the computer has a parallel port, skip steps 7a through 7e. Otherwise, custom-configure the Sentinel driver to act as though one were available, even though it ought not be used for a key.
  - a. Click the *Add* button. This opens the *Add Port* dialog.
  - b. The dialog's *Use this port?* field is set to *Yes*. Change it to *No*.
  - c. The dialog's *Bus Address* field is set to *0*. Change it to *378*.
  - d. The dialog's *Bus Type* field is set to *ISA*. Change it to *PCIBus*.
  - e. Choose *OK*. When asked, confirm that you wish to change the settings of this port. This closes the *Configure Port* dialog.
8. The Sentinel System Driver will converse with all parallel ports for which the entry in the *Use?* Column is *Yes*. The Driver need not converse with ports to

which the key is not attached. For each port to which the key is not attached, follow steps 8a-8c.

- a. Highlight the line for the parallel port and click on the *Edit* button.
- b. This displays the *Configure Port* dialog.
- c. The dialog's *Use this port?* field is set to *Yes*. Change it to *No*, and choose *OK*. When asked, confirm that you wish to change the settings of this port. This closes the *Configure Port* dialog.

9. Choose *OK*. This closes the *Sentinel Driver* dialog.

10. Exit the *Sentinel System Driver* application.

11. Attach the memory key.