

ClonePix FL Upgrading the Software

- Overview
- PC Requirements
- Backing up Previous Version Configuration
- Removing the Previous Version
- Installing Microsoft .Net 3.5 SP1
- Installing Microsoft DirectX (Nov 2007)
- Installing the ClonePix FL Software
- Enabling Changeable Process Property Defaults
- Re-Installing Previous Version of Software

Overview

This document describes how to upgrade the ClonePix FL software. To set up a PC from scratch, follow the 'Installing ClonePix FL Software' instructions instead.

PC Requirements

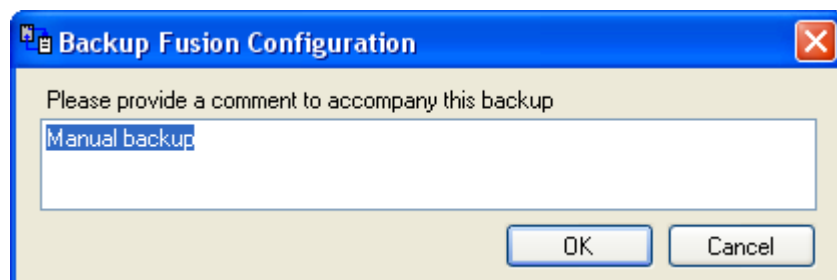
Versions 1.2.3.842 and later of the software require 2GB of memory in order to operate.

The software requires Windows XP with Service Pack 3 applied. If you see an error message referring to an incorrect version of Windows Installer, apply Windows XP Service Pack 3 to resolve the issue.

Backing up Previous Version Configuration

It is important to backup the previous software configuration in case you need to reinstall it for any reason.

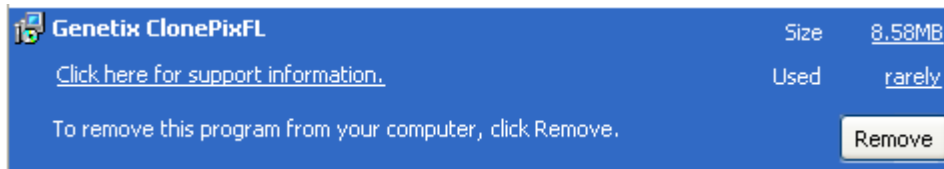
1. In Windows, open All Programs > Genetix > ClonePix FL > Configuration Manager.
2. In the Configuration Manager, click 'Backup to file' and enter the current software version number and date.



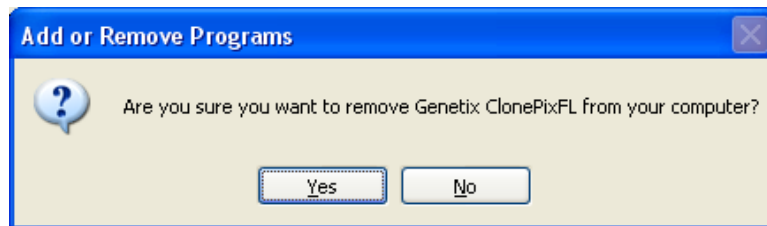
3. Click 'OK' and you will be prompted to save the backup file. Save the file to your desktop.

Removing the Previous Version

1. In Windows, open the Control Panel.
2. Double click the 'Add or Remove Programs' icon.
3. Select the 'Genetix ClonePix FL' item from the list and then click the 'Remove' button.



4. When asked if you are sure that you want to remove 'Genetix ClonePix FL', click the 'Yes' button.



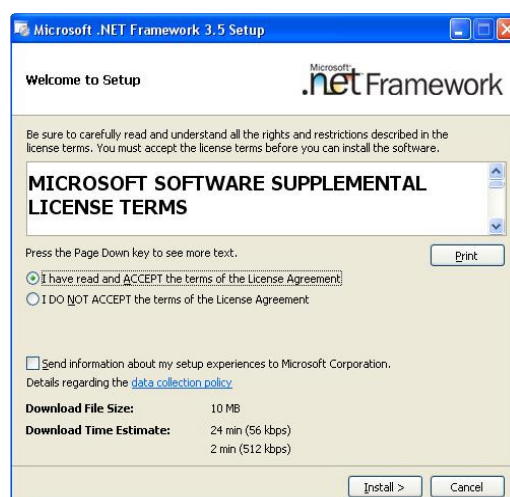
5. Once complete, check that all software is removed by opening All Programs. If the software is fully removed you should not see the Genetix folder.
6. If the software is not fully removed, right click on the folder and delete.

If you are upgrading from version 1.2.3.842 and later, skip forward to the section 'Installing the ClonePix FL software'.

Installing Microsoft .Net 3.5 SP1

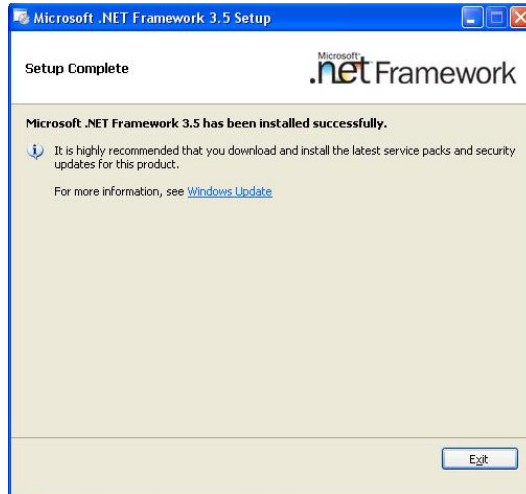
Microsoft .NET Framework 3.5 SP1 is required for the software to run. You can check whether it is currently installed by looking in the 'Add or Remove Programs' section of the Windows Control Panel. Be sure to check that version 3.5 SP1 is installed, rather than the older 3.5 without SP1. If it is not installed, follow these steps to install it.

1. Insert the recovery DVD and browse to `DVD:\Prerequisites\NET Framework 3.5 SP1`
2. Run the `'dotnetfx35.exe'` file found in that folder.
3. Accept the licence agreement and then click the 'Install >' button.



4. The install will now start. **WARNING:** This may take quite some time.
5. When shown the 'Setup Complete' screen, click the 'Exit' button.





6. If prompted, restart the computer.

Installing Microsoft DirectX (Nov 2007)

1. Insert the recovery DVD and browse to “DVD:\Prerequisites\DirectX 9.0c (November 2007)”
2. Run the ‘DXSETUP.exe’ file found in that folder.
3. Accept the licence agreement and then click the ‘Next’ button.

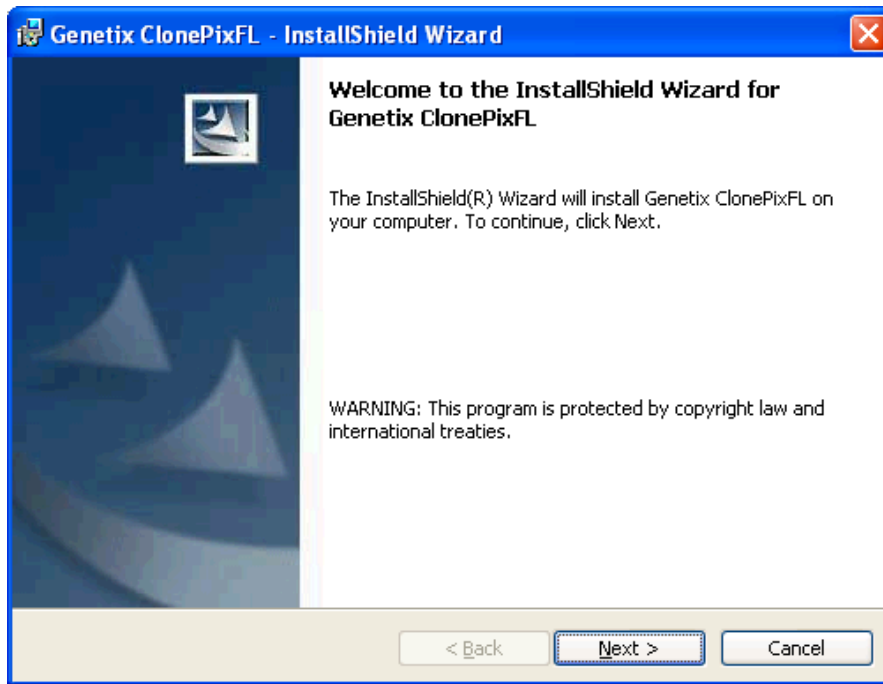


4. Click the ‘Next’ button.
5. On the ‘Installation Complete’ screen, click the ‘Finish’ Button.

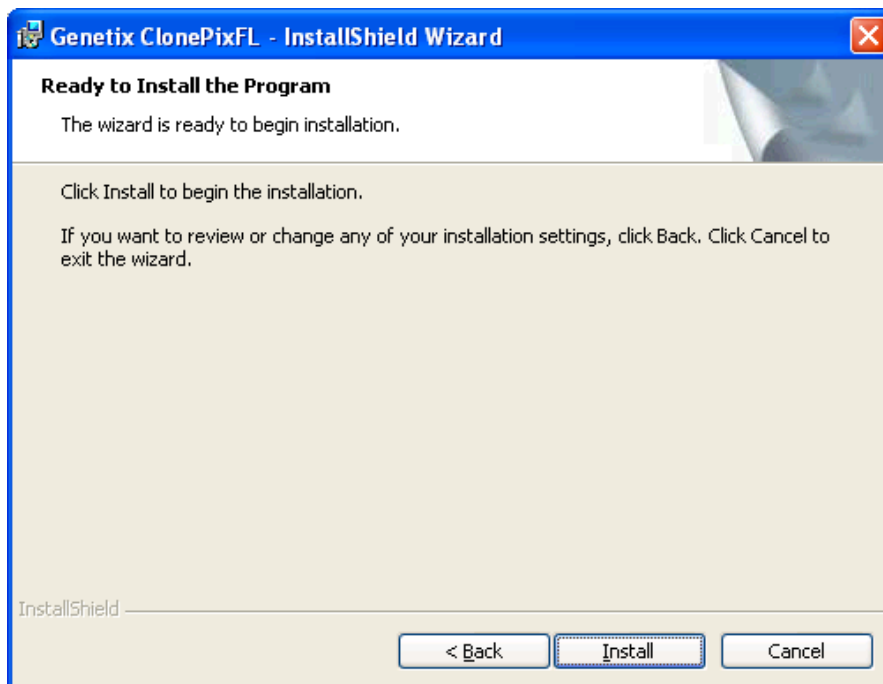
Installing the ClonePix FL Software

1. Navigate to the “Setup\ClonePix FL x.x.xx.xxxx” folder on the Recovery DVD.
2. Run the file “Setup ClonePixFL.exe”.



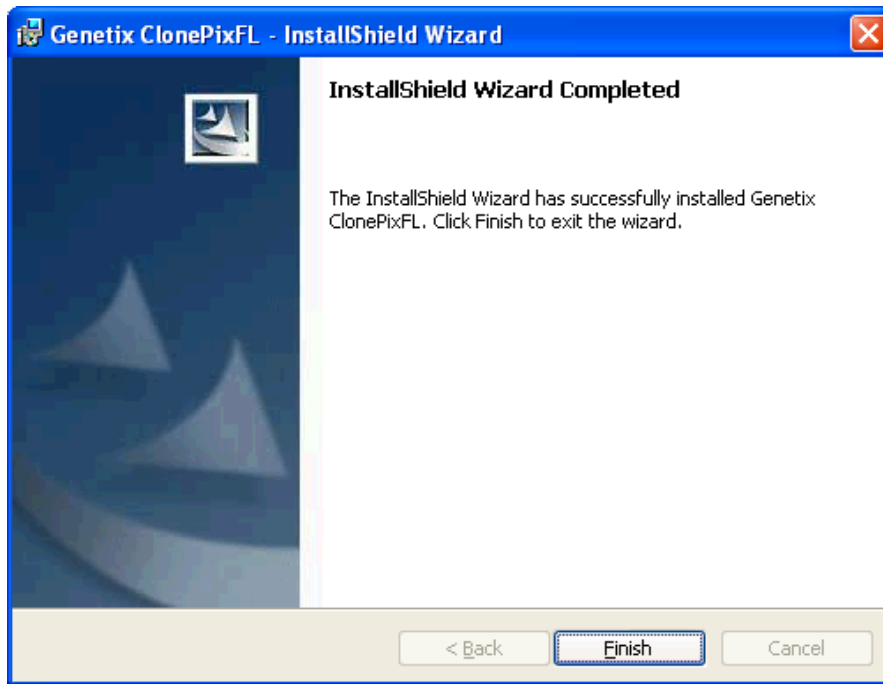


3. Click "Next" to continue.



4. Click "Install" to install the software.





5. Click “Finish” to complete the installation.
6. Run the “ClonePix FL” software from the “Genetix” folder on the Start menu. The software will prompt for a license – follow the on-screen instructions and acquire a valid license file from a Genetix Support representative.

Note: When updating software it is important to delete and recreate your templates. Templates created in previous version software may produce errors when opened in later software versions.

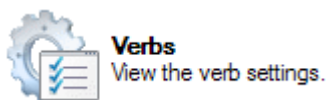
Enabling Changeable Process Property Defaults

A new feature available in this version of the software is the ability to change the default settings for new processes. You can use this to change the default well plate type, for example.

If you are upgrading from version 1.2.3.842 and later, this feature should have already been set up so you can skip these steps.

To enable this feature, follow these steps.

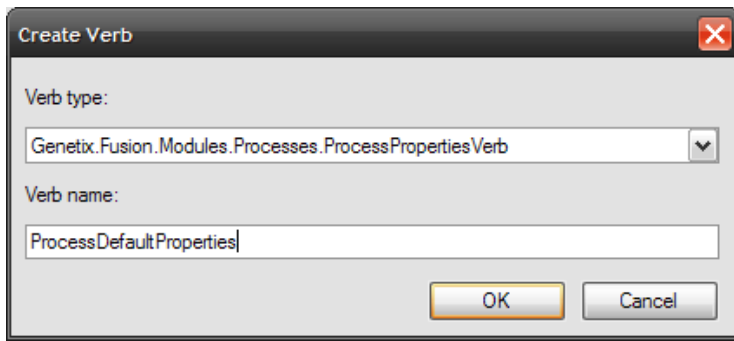
1. Open the ClonePix FL application.
2. Select ‘Tools’ and then ‘Configuration..’ from the menu bar.
3. Click the ‘Verbs’ icon.



4. Click the ‘Add...’ button.
5. From the ‘Verb type:’ dropdown box, select ‘Genetix.Fusion.Modules.Processes.ProcessPropertyVerb’.



- In the 'Verb name:' text box, type 'ProcessDefaultProperties'.



- Click the 'OK' button.
- Click the 'Close' button.
- Close the software.

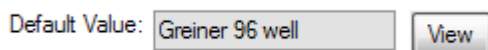
Changing the Default Properties for a Process

In order to change the default values for new processes, follow these optional steps:

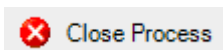
- Open the ClonePix FL application.
- Open the process for which the default properties are to be changed.
- Select 'View' and then 'Administrate Properties' from the menu bar.
- Select the property that you would like to change from the list.

Group	Name	Caption	Type	Default Value
Source	WellPlates	Well plates	Genetix.Fusion....	Greiner 96 well
Barcodes	readbarcode	Read barcode	System.Boolean	Yes
Setup	reviewfocus	Review focus	System.Boolean	Yes
Setup	autofocus	Auto focus before run	System.Boolean	No
Setup	Scan Type	Scan Type	Genetix.Fusion....	Full Scan
Setup	EnhanceImages	Enhance Images	System.Boolean	Yes
Setup	PipeLine	Processing Type	Genetix.Fusion....	Typical Adherent Cell Co...

- Click the 'View' button next to the 'Default Value:' option.



- Select your new default value, then click the 'OK' button.
- Repeat steps 4 to 6 for each property that you would like to alter.
- Select 'View' and then 'Properties' from the menu bar.
- Click the 'Close Process' item from the menu on the right of the window.



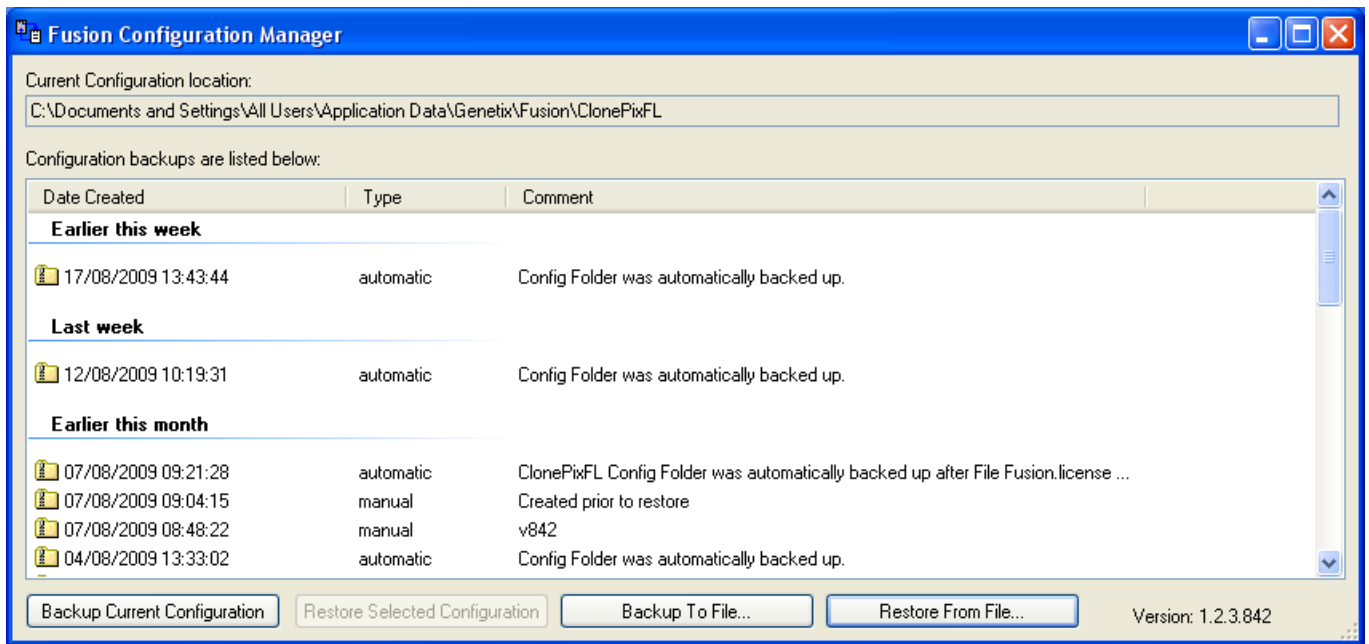
- Open the process again. Your new default properties should now be shown.



Re-Installing Previous Version of Software

If for any reason the newly installed version of software is not functioning correctly you can reinstall the previous version.

1. Remove the newly installed software as described above in the **'Removing Previous Version'** section. You do not have to backup the configurations for this version.
2. Insert the recovery DVD for the previous version of software and follow the instructions as in the **'Installing the ClonePix FL Software'** section above. The older version of software will function correctly using the newly installed prerequisites so there is no requirement to adjust these.
3. Once the software is installed, open All Programs > Genetix > ClonePix FL > Configuration Manager.
4. In Configuration Manager, click on 'Restore From File'.



5. Open the previously saved configuration from the desktop.
6. Once the configuration is restored, exit the Configuration Manager.
7. Run the "ClonePix FL" software from the "Genetix" folder on the Start menu. The software will prompt for a license – follow the on-screen instructions and acquire a valid license file from a Genetix Support representative.

