

ClonePix FL

Installing the Data Viewer Software

Overview

The ClonePix FL Data Viewer software enables you to view results generated by the ClonePix FL on a PC remote from the instrument. You can either access a copy of the results files, for example on a DVD, or you can access the folder on the instrument PC using Windows folder sharing if the two PCs are networked together.

Requirements

To install the data viewer software, the PC must meet the following minimum requirements:

Operating System:	Microsoft Windows XP with Service Pack 3 or: Microsoft Windows Vista with Service Pack 1 Note: currently only 32-bit operating systems are supported.
Memory:	2GB minimum.
Processor:	1GHz Pentium compatible processor.
Display:	1024x768 32-bit minimum

Note: You must have local administrator privileges to install the software onto your PC. If your PC is attached to a corporate IT network, you may need assistance from your IT support department in order to install the software.



Installation

Prerequisites

The ClonePix FL Data View software requires the Microsoft .NET Framework 3.5 Service Pack 1 to be installed. Some PCs may already have this, so it is important to check whether it is installed before proceeding.

To check if the .NET Framework 3.5 SP1 is installed, open up Control Panel and click the 'Add/Remove Programs' icon. Look through the list of installed programs for 'Microsoft .NET Framework 3.5 SP1'. You might see versions 1.0, 1.1 or 2.0, however these are not sufficient to run the software. You must have version 3.5 SP1 installed.

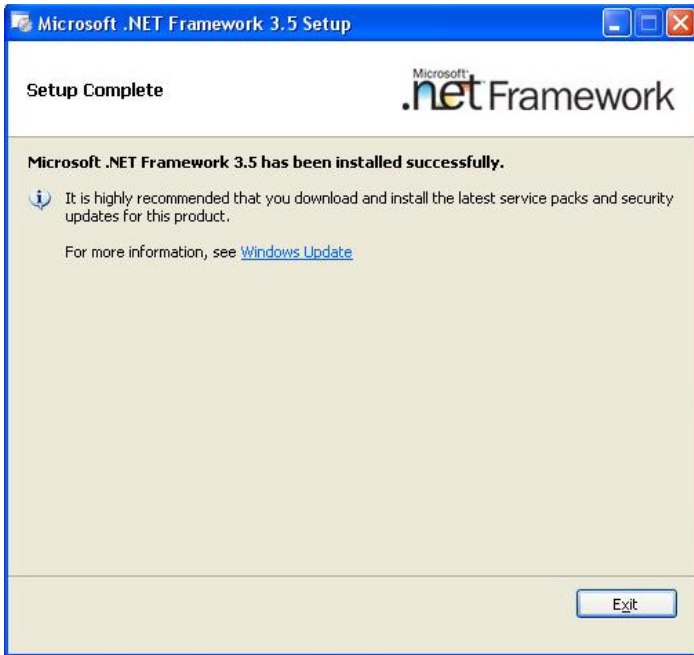
If the .NET Framework 3.5 SP1 is installed, move on to the next section, "Installing Microsoft DirectX (Nov 2007)", otherwise follow the instruction below on how to install .NET Framework 3.5 SP1.

Installing Microsoft .Net 3.5 SP1

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.



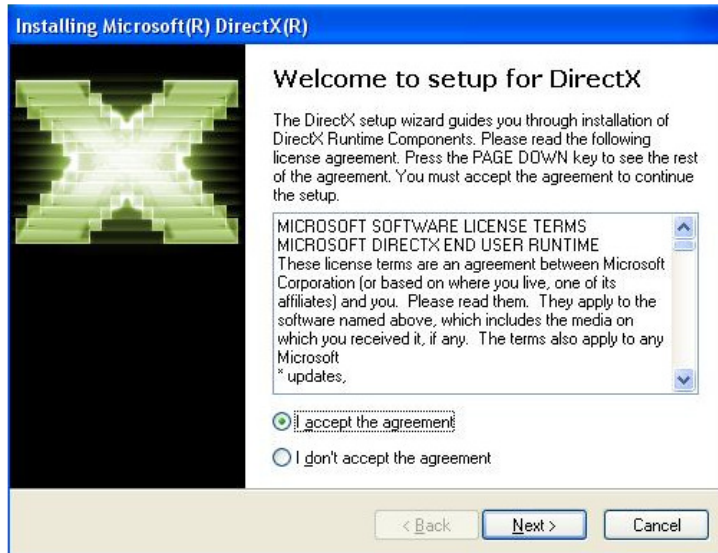
4. The install will now start. **WARNING:** This may take over an hour on some PCs.
5. When shown the 'Setup Complete' screen, click the 'Exit' button.



6. If prompted, restart the computer.

Installing Microsoft DirectX (Nov 2007)

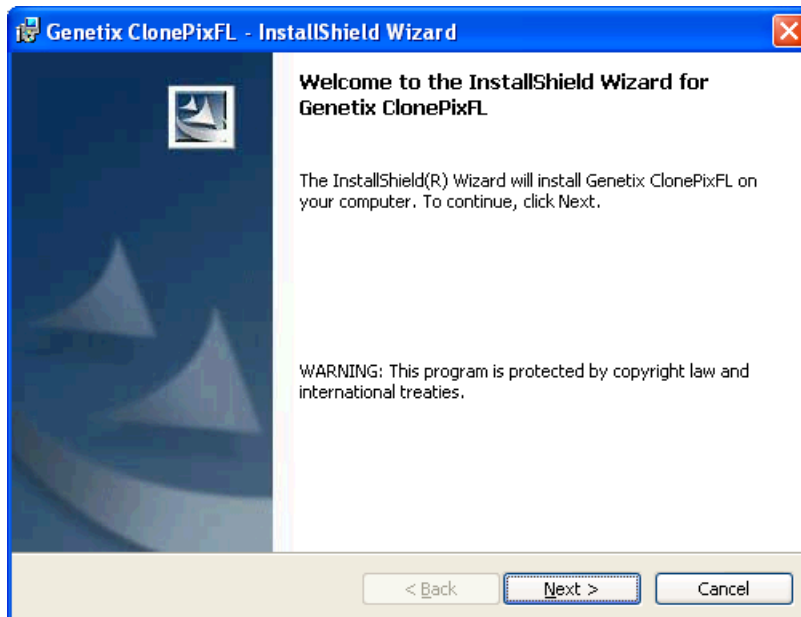
1. Insert the recovery DVD and browse to "DVD:\Prerequisites\DirectX 9.0c (November 2007)"
2. Run the 'DXSETUP.exe' file found in that folder.
3. Accept the licence agreement and then click the 'Next' button.



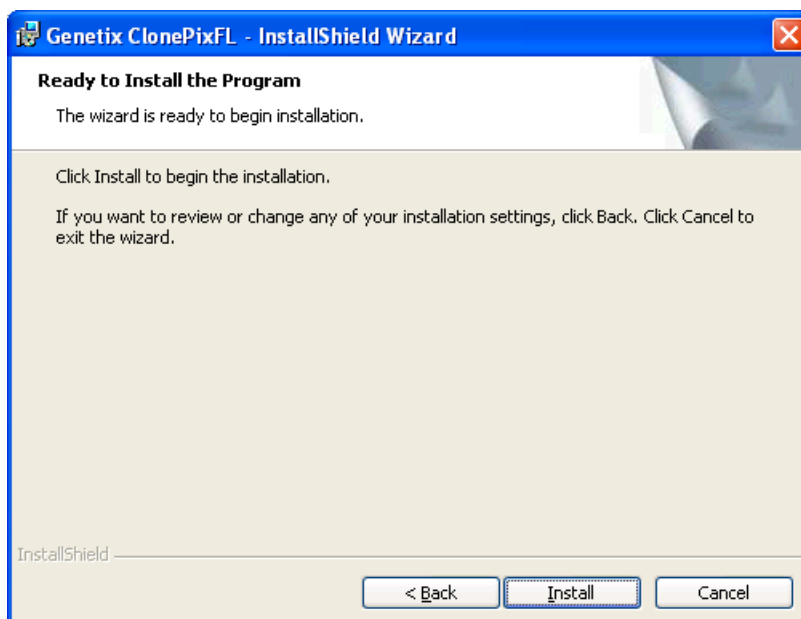
4. Click the 'Next' button.
5. On the 'Installation Complete' screen, click the 'Finish' Button.

Installing the ClonePix FL Data Viewer Software

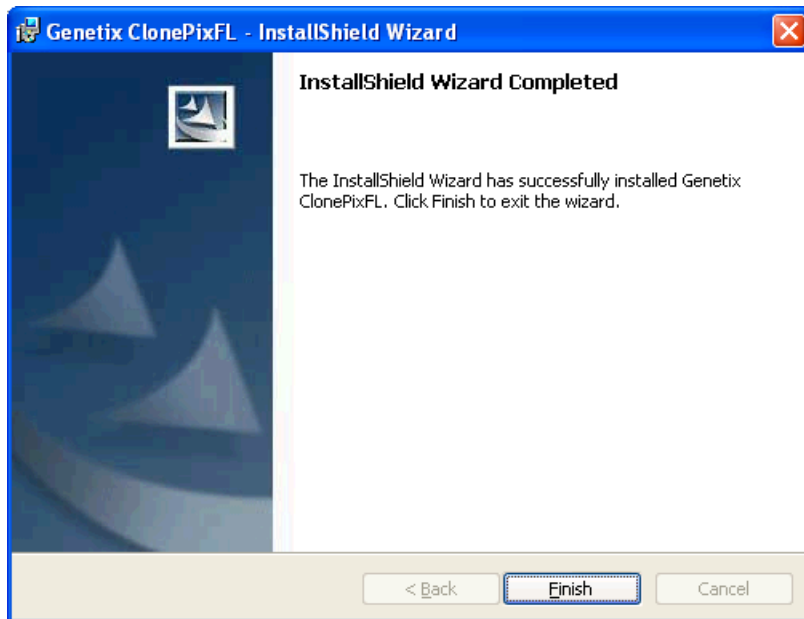
1. Navigate to the "Setup\ClonePix FL Data Viewer x.x.xx.xxxx" folder on the Recovery DVD.
2. Run the file "Setup ClonePixFL Viewer.exe".



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.

Licensing the ClonePix FL Data Viewer

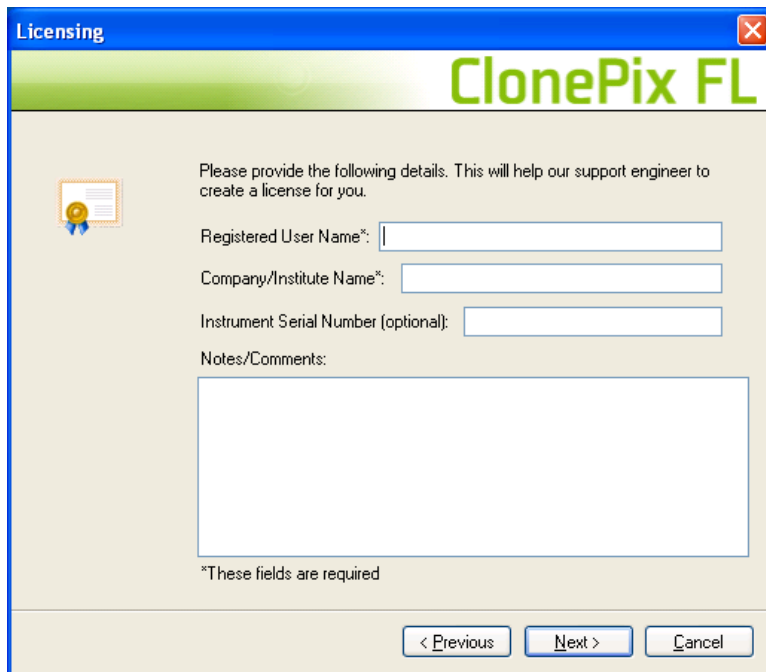
Once the ClonePix FL Data Viewer is installed, you must then license it for use. To do this, run the software from the 'Genetix' folder in the start menu. When the software first runs, it will detect that a license is required. It will prompt you to generate a license request file. You must send this file to Genetix, where we can then generate a license file for you and send it back. You must then import this file into the software, which will then be licensed and ready to use. These steps are outlined below.

1. Run the software. It will detect that a license is required and prompt you:



Ensure that 'Request a new license' is selected, then click the 'Next' button.

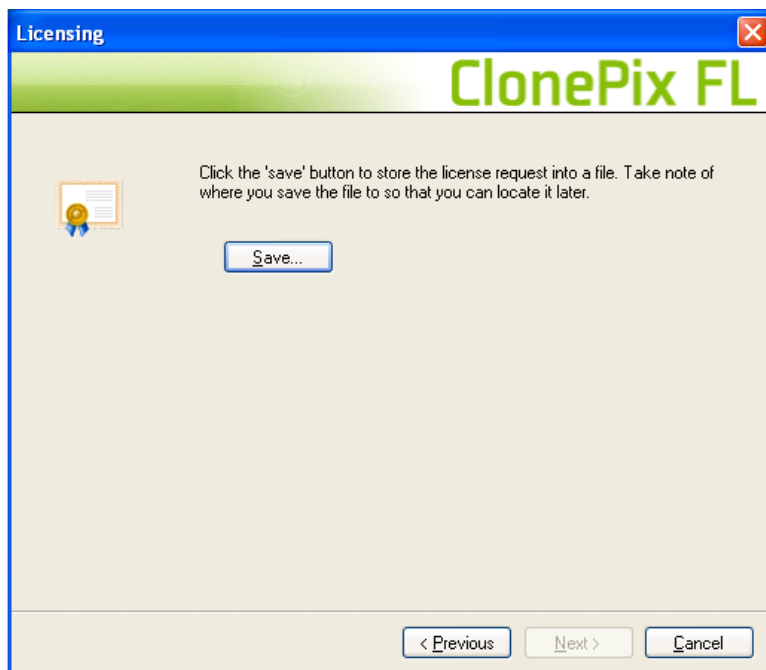
2. The software will prompt you for your details:



The image shows a software dialog box titled "Licensing" for "ClonePix FL". The window has a blue title bar with a close button. Below the title bar is a green header with the "ClonePix FL" logo. The main area contains a document icon with a ribbon, followed by the text: "Please provide the following details. This will help our support engineer to create a license for you." There are three input fields: "Registered User Name*", "Company/Institute Name*", and "Instrument Serial Number (optional)". Below these is a larger text area for "Notes/Comments:". A note at the bottom states "*These fields are required". At the bottom of the dialog are three buttons: "< Previous", "Next >", and "Cancel".

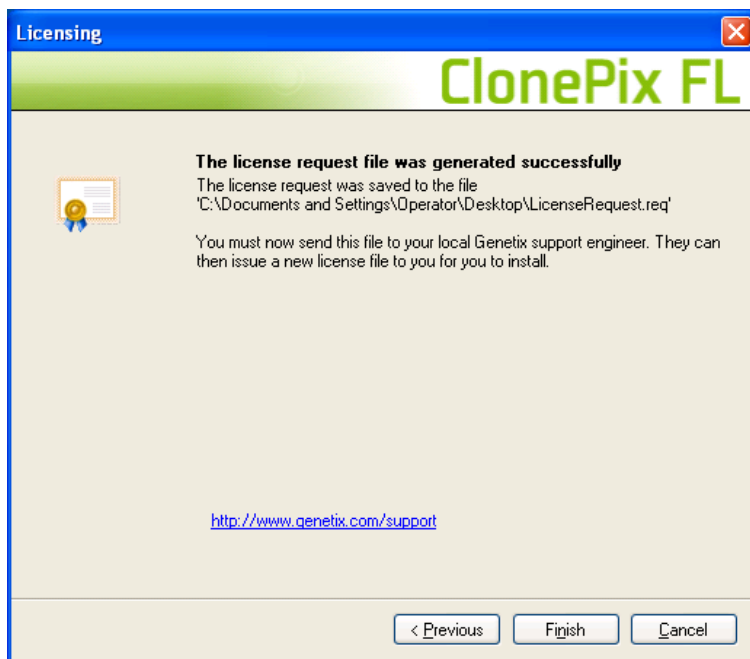
Fill these out and then click the 'Next' button.

3. The software prompts you to save the license request file. Click the 'Save' button and browse to a location that you can access later (e.g. the computer desktop).



The image shows the same "Licensing" dialog box for "ClonePix FL". The main area now contains a document icon with a ribbon and the text: "Click the 'save' button to store the license request into a file. Take note of where you save the file to so that you can locate it later." A "Save..." button is centered in the dialog. At the bottom are three buttons: "< Previous", "Next >", and "Cancel".

4. A message will be displayed when the file is saved successfully:



You must now email this file to Genetix so that we can issue a license file to you. You can use the website link in the dialog to access our Customer Support contact details, or you can email the request file to your local sales representative who will deal with it for you.

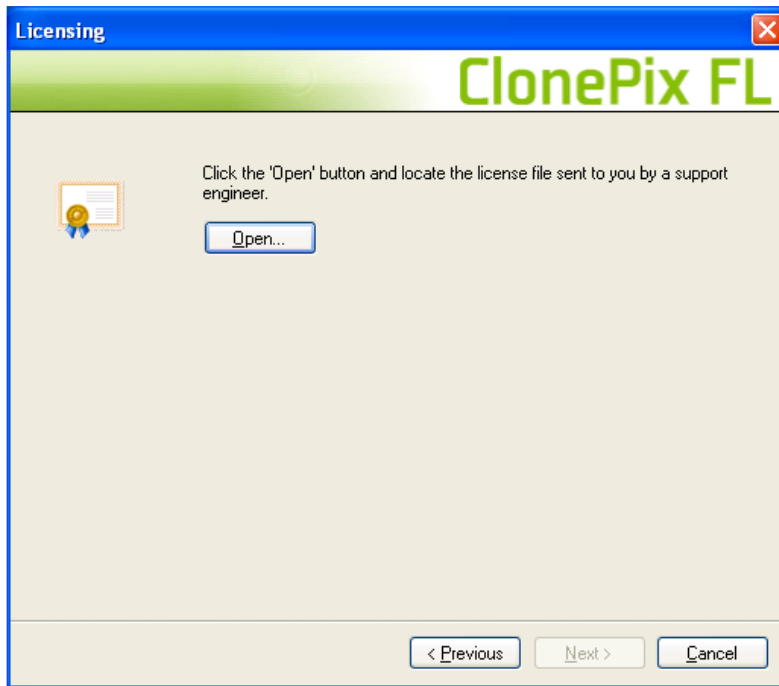
Click the 'Finish' button to return to the start of the licensing process. If you wish, you can cancel the dialog to close the software. You can return to the licensing process by starting the software again.

5. Once you have received your license file from Genetix, you can then install it. First save it out of the email to a location that you can access later (such as the computer desktop). Then return to the licensing process (by starting the software again if necessary).

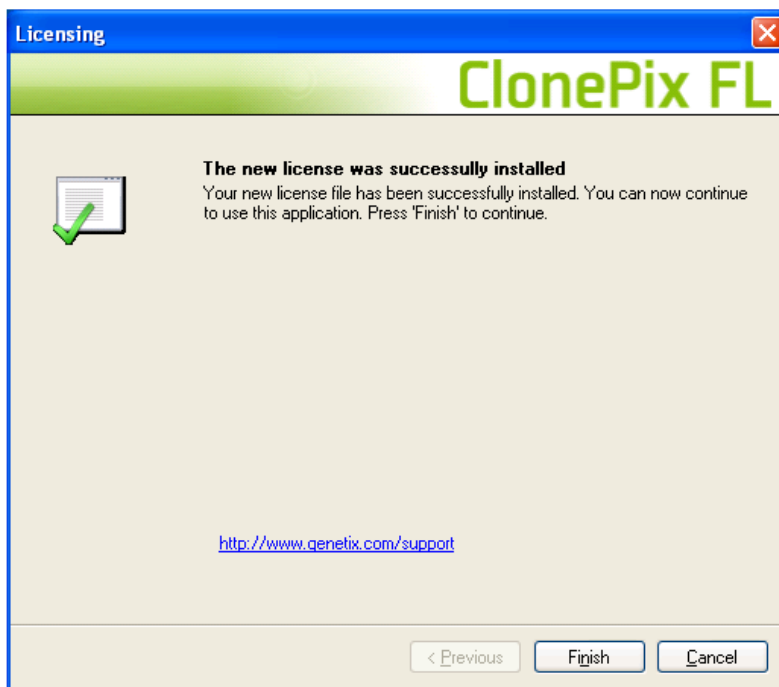


Ensure that 'Install a license file' is selected, then click the 'Next' button.

6. To install the license file, click the 'Open' button and browse to the location you have saved the license file to.



7. Once the license file is installed, a message will be displayed.

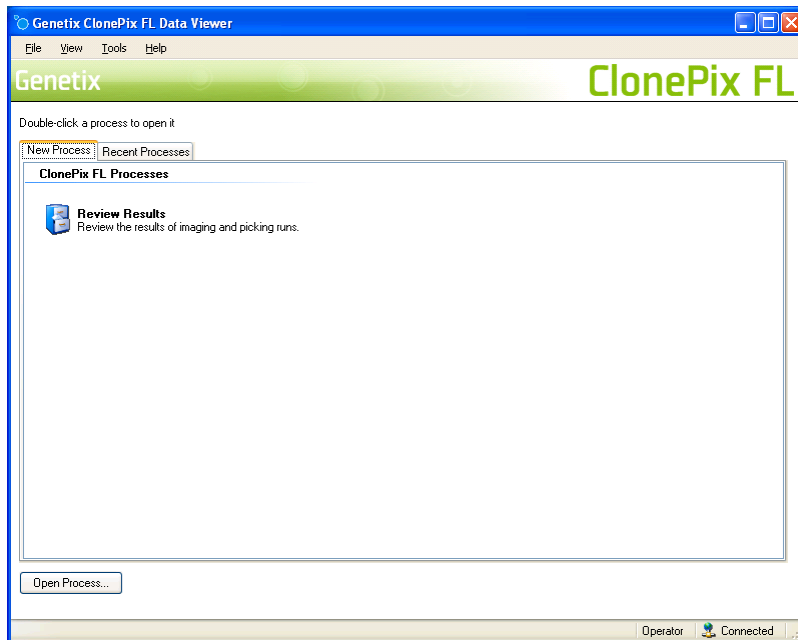


If the license failed to install for some reason, a message giving details will be displayed. Contact Genetix Customer Support for assistance if required.

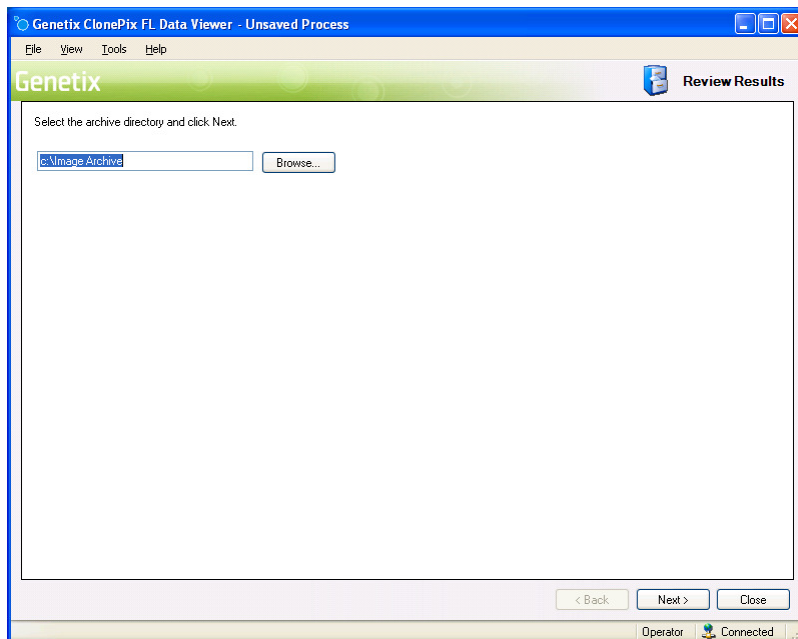
If the license installs successfully, click 'Finish' to start the software.

Using the ClonePix FL Imager Data Viewer

When the software starts, the main process screen is displayed:



Double-click the 'Review Results' icon to start the process:



To view the results, you must browse to the location where they are stored. Click the 'Browse...' button, and browse for the location of the results. This could be a local folder on the PC, a CD-ROM or DVD, or it could be a shared folder on the instrument PC if the two computers are networked together.

Once you have selected the location, the software will catalogue all of the results and display a list of them to you. You can then view the results in the same way as you can on the instrument PC.