



CloneSelect Imager

OPERATOR MANUAL
SOFTWARE RELEASE 1.4



Contents

Introduction	5
CloneSelect Imager	5
System Features	5
Before Starting	5
Hardware Guide	6
Layout and Dimensions	6
Indicator Lights	6
Setup	8
Calibration and Alignment	9
Power-Up Procedure	9
Starting up CloneSelect Imager	10
CloneSelect Imager Software	10
Menu Options	11
File Menu	11
View Menu	11
Tools Menu	12
Help Menu	12
Main Navigation Screen	13
New Process Tab	13
Templates Tab	13
Recent Processes Tab	13
Open Process Button	13
CloneSelect Imager Processes	14
Image Microplate	14
Imaging Criteria Setup	14
Select Source	16
Load Plate	17
Focus	17
Alignment	19
Imaging	19
Results	19
Remove Plate	20
Finish	20
Review Results	21

Load Results	21
Analysing Results	25
Migration Application.....	38
Batch Microplate Imaging.....	46
E-Scope.....	46
How to Image Using E-Scope	46
Plate Alignment	47
Eject Plate Carrier	47
Data Analysis Examples	49
Monoclonality Workflow.....	49
Colony Forming Assay Workflow	52
Health and Safety	55
Transport and Storage	55
Lifting Points.....	55
External Covers.....	55
Electrical Safety	55
Drive Safety	55
Noise Levels	55
Maintenance.....	56
Cleaning CloneSelect Imager.....	56
Regular Maintenance	56
Daily.....	56
Weekly.....	56
Annually.....	56
General Precautions.....	56
Technical Specifications	57
Dimensions	57
Operating Environment	57
Electrical Supply	57
Symbols on Equipment.....	57
Appendix A	58
Compatible Plate Types	58
Appendix B	59
Well List File Format.....	59
Appendix C	60
Setting up Cell Number Estimation.....	60
Setting up a dilution series.....	60

Creating a cell number estimation formula	61
Appendix D	63
Requesting a License	63
Installing a License	66
Index	68
Obtaining Support	69

Introduction

CloneSelect Imager

CloneSelect™ Imager is an automated CCD camera-based system for imaging and confluence determination of cells growing in microplates.

Note: CloneSelect Imager is strictly for research use only and is not intended or recommended for the diagnosis of disease in humans or animals.

If the instrument is used in a manner not specified in this manual the protection provided by the equipment may be impaired.

For correct functioning of the system, users must have read and write access to the c:\ProgramData\Genetix\Fusion\CloneSelectImager config folder.

System Features

- Automated, rapid confluence determination of cells growing in microplates
- Automatically plots time course of cell growth
- Cell number estimation
- Monoclonality indication
- Colony forming assay
- Fast scan and full scan modes
- Standard and high resolution imaging
- Exports data and images for further analysis

Before Starting



Before using the instrument it is very important to read this manual and understand all the safety instructions. Then follow the procedures as set out in the Power Up section

Hardware Guide

Layout and Dimensions

Indicator Lights

The indicator lights show the status of the instrument.



On



Ready



Imaging



Vacuum On



Warning

Plate Carrier

The plate carrier consists of a glass plate surrounded by a vacuum bed seal. The microplate is placed on the plate carrier so that the skirt is over the vacuum bed seal. When the plate carrier retracts into the instrument for imaging, two pushers gently push the microplate into the back right corner of the plate carrier so that the plate is always in the correct position for imaging.

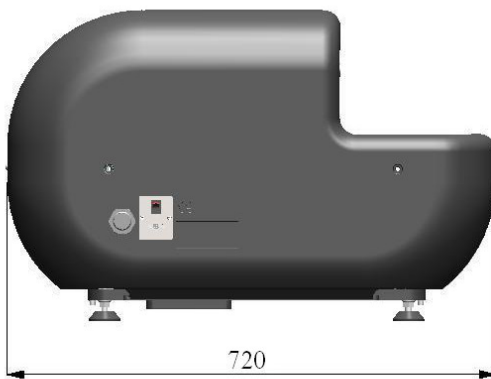
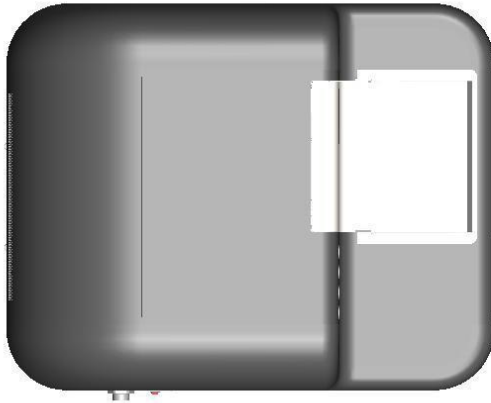
Before imaging, a vacuum is applied under the plate so that it is held down flat on the glass plate so that all the wells are in the same focal plane. This obviates the need to autofocus on every well.

Leveling Feet

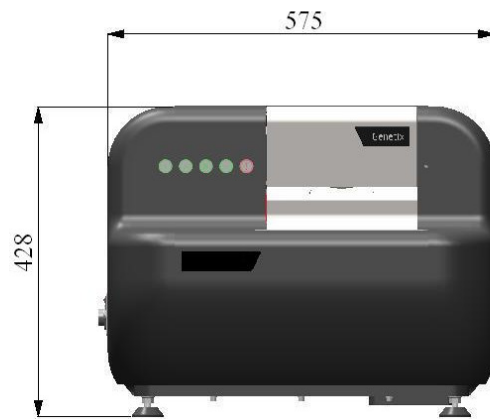
The leveling feet can be adjusted by loosening the locking nuts and turning the feet until the instrument is level and stable. Retighten the locking nuts once any adjustments have been made.

The overall layout of the instrument.

Top View



Side View



Front View

Note: All dimensions are in millimetres.

Setup

Locating CloneSelect Imager

CloneSelect Imager should be situated on a stable, level surface. Level the instrument by loosening the locking nuts and turning the leveling feet to ensure that it is stable and not able to move. Retighten the locking nuts after adjustment.

Workstation

CloneSelect Imager is supplied with a Dell T3500 Workstation with special hardware and software to support the imaging function. Do not attempt to use any other workstation to operate CloneSelect Imager.

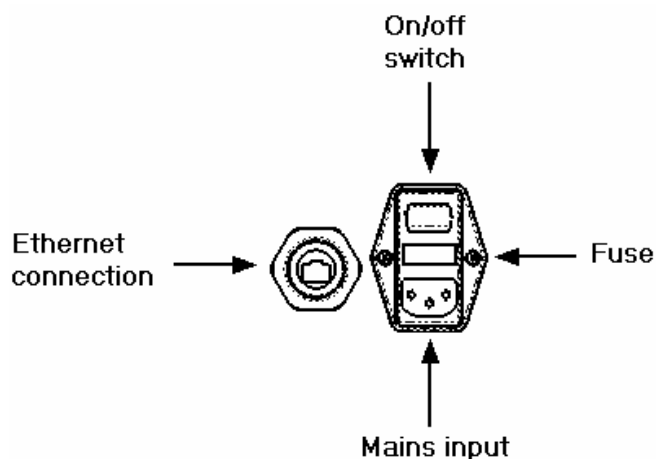
The workstation is supplied with Microsoft Windows 7™ and will require security configuration if it is to be connected to a network. If that is done, ensure that no changes are made to the configuration of the private network connection to the CloneSelect Imager since that will stop it operating.



Warning – CloneSelect Imager is optimized for use with Microsoft Windows 7™ and Microsoft Office™ with standard English settings. The performance of the system may be affected if these settings are changed.

Connecting CloneSelect Imager

The connections to **CloneSelect Imager** are on the connections panel on the left hand side of the instrument.



Connect the mains cables to **CloneSelect Imager**, the workstation and to the monitor before making the network connections.

The connection to CloneSelect Imager is a private Gigabit Ethernet link and the connection must be made to the workstation add-on card Gigabit Ethernet port.

Calibration and Alignment

Calibration and alignment are performed by a Molecular Devices engineer before CloneSelect Imager is shipped. If the instrument requires recalibration and alignment follow the procedure below:

- Select the E-Scope process on the Main Navigation screen.
- Select the PetriWell W1500 plate as the plate to be imaged.
- Start the run and insert the Fixed Cell plate (shipped with instrument) when prompted.
- Align on to well A1 of the Fixed Cell plate.
- On the image selection screen select the lower left hand quadrant image of well D1.
- Adjust the focus until the cells on the Fixed Cell plate are in focus.
- Image wells A12 and H12 and check that the focus is correct in each image (focus should not be changed at this point, just reviewed).

If the focus is not correct, contact Molecular Devices to arrange for the machine to be re-aligned.

Power-Up Procedure

The recommended power-up procedure is:

- Switch on CloneSelect Imager.
- Wait for the Ready light to illuminate.
- Leave CloneSelect Imager in the 'Ready' state for more than 5 minutes prior to use (in cases where the unit is cold to the touch, wait for 30 minutes prior to use).
- Switch on the workstation and start CloneSelect Imager software by clicking on the desktop shortcut. The instrument is now ready for use.

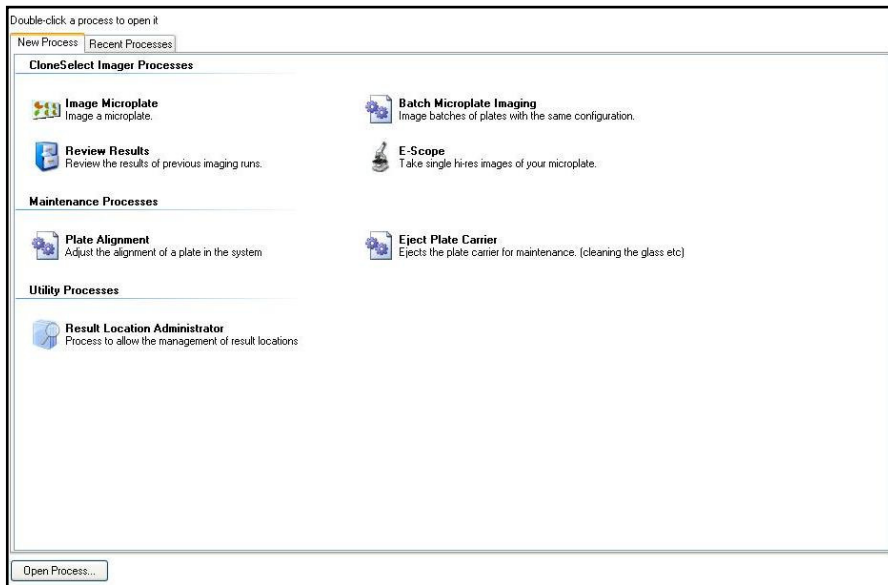
Rebooting CloneSelect Imager

If the instrument has to be rebooted, switch it off. Wait for 30 seconds then switch on again. Once the Ready light is illuminated the instrument is ready to run.

Starting up CloneSelect Imager

Before using CloneSelect Imager, please refer to the relevant sections of this manual for important setup, maintenance and safety information.

To start CloneSelect Imager software, double click the icon on Windows desktop. The software will connect to the instrument and display the Main Navigation Screen.



CloneSelect Imager Software

CloneSelect Imager software is based around the concept of the instrument running processes. The instrument has a default set of processes that can be modified and saved. Processes are managed from the **Main Navigation Screen**, which is the view displayed when the software starts.

Menu Options

The same menu is displayed in all the windows, but some item availability may vary between windows.

File Menu

Open Process

Allows previously saved processes to be opened.

Save Process

Saves the current process.

Save Process As

Allows the current process to be saved with a different name or location.

Close Process

Closes the current process and re-opens the Main Navigation Screen.

Save As Template

Saves the current process as a template.

Recent Processes

Allows a recent process to be selected and opened.

Switch User

Allows a user to log off and another user to sign on when User Authorization is licensed within the software.

Exit

Exits CloneSelect Imager software.

View Menu

Properties

The view of the process window when a process's properties are being set and prior to starting the process.

Progress

The view of the process window when a process is executing.

Administrative Properties

Can be used to alter the way properties are displayed. It is not recommended that users alter these properties.

Tools Menu

Configuration

Provides access to the configuration settings for the instrument. It is strongly recommended that only Molecular Devices staff alter these settings.

Hardware Model View

Accesses the datum positions for the instrument. It is strongly recommended that only Molecular Devices staff alter these settings.

Prepare Error Report

A wizard will create a data file containing the configuration and recent log files to help Molecular Devices Support diagnose recent problems.

Help Menu

About

Displays the version numbers of the software modules. Most of these have the same version number so, if asked to provide the software version number, give this number unless otherwise directed.

Online Support

Opens a web browser on the Online Support web page. This function only operates if the workstation has an internet connection, e.g. via a network.

CloneSelect Imager User Guide

An electronic summary of this operator manual is available; however this guide is the comprehensive overview of the CloneSelect Imager software.

Main Navigation Screen

The main screen enables navigation of CloneSelect Imager imaging and maintenance processes.

New Process Tab

The New Process tab provides **CloneSelect Imager Processes** for imaging microplates or reviewing results. It also contains **Maintenance Processes** for various setup features of the system.

CloneSelect Imager Processes

Image Microplate

Standard imaging of microplates.

Review Results

Review of results from previous runs including confluence, cell number estimation and monoclonality analysis.

Batch Microplate Imaging

Imaging of batches of plates with the same configuration.

E-Scope

High resolution imaging of microplates.

Maintenance Processes

Plate Alignment

Allows the alignment of a plate to be adjusted in the system.

Eject Plate Carrier

Ejects the plate carrier for cleaning and removal of microplates.

Utility Processes

Results Location Administrator

Allows the management of the results locations

Templates Tab

Saved templates can be retrieved by clicking on this tab. Template files can be saved as .gft files using the '**Save process as template...**' option found in **Image Microplate**.

Recent Processes Tab

Recent processes can be retrieved by clicking on this tab. Recent processes can be saved as .gfp files using the '**Save process**' option found in **Image Microplate**.

Open Process Button

Opens the same Open Process dialogue box accessed by the File > Open Process menu item.

CloneSelect Imager Processes

This section explains in detail how to run processes on CloneSelect Imager.

Image Microplate

This is the default process for imaging cells in microplates to determine confluence and carry out further analysis, such as cell number estimation and monoclonality.

In the **Main Navigation Screen** clicking on **Image Microplate**, or opening a previously saved Image Microplate-based process, will open the process. Once opened a summary window will be displayed in order for the process criteria to be set up.

Imaging Criteria Setup

Criteria for imaging a microplate are selected in this screen.



There are several options for selecting the way in which the process settings can be displayed for editing. The following options can be selected for editing criteria:

Summary

This is the default display option. The current settings for the process are displayed grouped in categories. Click on the category heading to call up a screen where the relevant settings can be altered via drop down lists.

Details

The current process settings are displayed in drop down lists from which they can be edited directly.

Guide

This option displays the process setup via a wizard style interface with Back and Next buttons to scroll through the settings.

If the options above are not used for editing the criteria the sections headers can be selected and drop down menus will be displayed for editing.

Source

Select the microplate to be imaged from the drop down list. The list contains a number of compatible microplates from a variety of manufacturers. If the desired plate does not appear in the drop down menu list, contact Molecular Devices.

Barcodes

Enable or disable barcode reading by ticking the box. If barcode reading is enabled and the instrument fails to read the barcode, or barcode reading is not enabled, a screen is displayed prompting for a barcode/plate name to be entered manually. The barcode/name can also be edited within **Review Results** at a later stage if required.

Setup

Review focus

If enabled the **Focus** window is displayed before the microplate is imaged to allow the focus and brightness settings to be checked before imaging a microplate. Individual images can also be saved from the **Focus** window.

Auto-focus before run

If enabled the instrument carries out an auto focus procedure before imaging the microplate.

Scan type

Select Full Scan to image the entire area of each well in the microplate. In Partial Scan an area from the centre of each well is imaged. This speeds up imaging and is suitable when the data from the centre of each well is representative of the well as a whole.

Enhance Images

Selecting this option will normalize / flatten the background of the images to enhance them for display.

Processing Types

The following processing types have been developed for cell and colony detection.

Cell Detection Method 1

Formally known as 'Typical Adherent Cell Confluence', this algorithm is ideal for settled suspension cells and adherent cells with good contrast.

Cell Detection Method 2

This algorithm helps to detect cell samples when contrast is low, cells are clumped or at the edge of the well.

Cell Detection Method 3

This algorithm is designed to detect colonies and distinguish them from any artefacts around the edges of the colonies.

Start

Once the process's properties have been set, click **Start** to enter the process.

Select Source

Wells selected to be imaged are shown in red, those not selected for imaging are shown in light blue. Left click on a well to select it and right click to deselect. Groups of wells can be selected or deselected by clicking the appropriate button and dragging over the wells.

The microplate type to be imaged can be altered via the drop down list.

Import

Displays an Open File dialogue box for importing a file containing a list of wells to be imaged. The format of the file is given in **Appendix B**.

Select All

Selects all the wells to be imaged.

Deselect All

Deselects all the wells.

Scan Options

Scan Type

There are two scan type options – Full and Partial Scan. A full scan images the whole well, so for a 96 well plate the imager takes four images per well to capture to entire well. A partial scan images the centre of the well only so for a 96 well plate the imager would take a single image. The partial scan results in a faster scan of all the wells to be imaged.

Review Focus

By selecting this tick box, the focus of the image will be reviewed before the scan begins.

Next

Once the wells to be imaged have been selected, click **Next** to move to the next stage. This button is present throughout the entire process but may be grayed out where its use is not applicable.

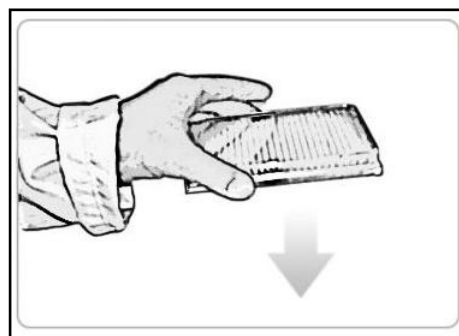
Cancel

Click **Cancel** to exit the process and return to the starting screen of the **Image Microplate** process. The button is present throughout the entire process but may be grayed out where its

use is not applicable.

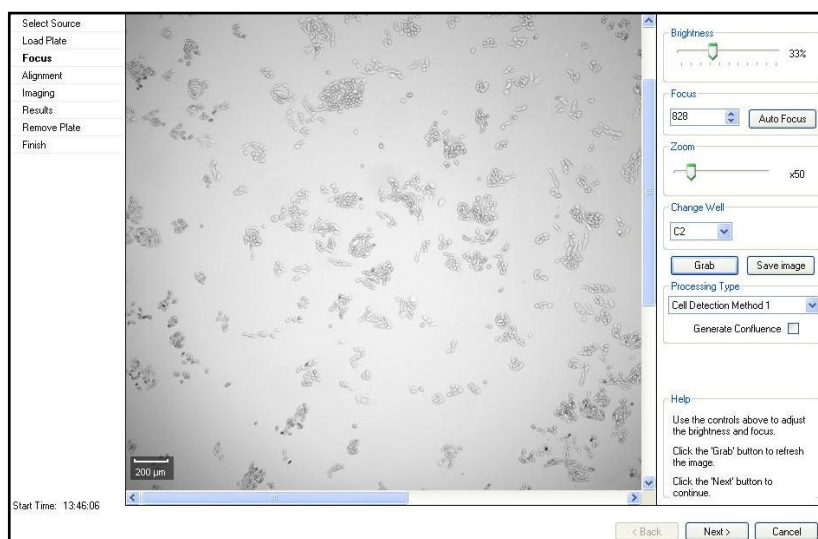
Load Plate

The plate carrier is ejected to enable the loading of the microplate. Ensure that well A1 is to the back right corner of the plate carrier.



Focus

The **Focus** window allows the image to be optimized before the scan takes place.



Brightness

The software automatically determines the optimum brightness setting for imaging. However, the setting can be adjusted via the **Brightness** slider. After adjusting the brightness the **Grab** button must be clicked to retake the image using the new settings. Saturated pixels will be displayed in red, so it is important not to make the image too bright.

Focus

The system stores the last focus setting used for each plate type. So after a plate type has been imaged for the first time the focus setting should not need to be altered significantly. After altering the focus setting the **Grab** button must be clicked to retake the image using the new setting. The altered focus setting now becomes the default focus setting for the plate type being used in this run.

Clicking the **Auto Focus** button will start the autofocus procedure. This may take a minute or two. If there is debris on the bottom of the plate or on the glass plate carrier the autofocus may focus on these instead of the cells. If this is the case, the focus will have to be adjusted manually.

Zoom

The **Zoom** slider can be used to zoom into the image. This can also be achieved by clicking on the image and using the mouse wheel to zoom into the image. While the zoom setting

figure (e.g. x 99) displayed to the right of the slider is black the system is zooming optically. When the figure turns red the system is zooming digitally and some pixilation of the image may be seen.

Change Well

By default focusing is performed on well A1 or the first well of the plate to be imaged if A1 is not selected to be imaged. Different wells can be selected to focus on from the drop down list. After selecting the desired well the **Grab** button must be clicked to display the newly selected well.

Grab

This **Grab** button must be clicked after altering any settings to retake the image.

Save Image

Clicking on **Save Image** will display a **Save File** dialogue that will allow the current image used for focusing to be saved. Images can be saved in .bmp, .jpg or .png format.

Processing Type

This drop down menu displays the three **Cell Detection Methods** to choose from. See page 15 for the description of the three **Cell Detection Methods**.

Generate Confluence

Selecting this box will identify and highlight in green all objects within the well.

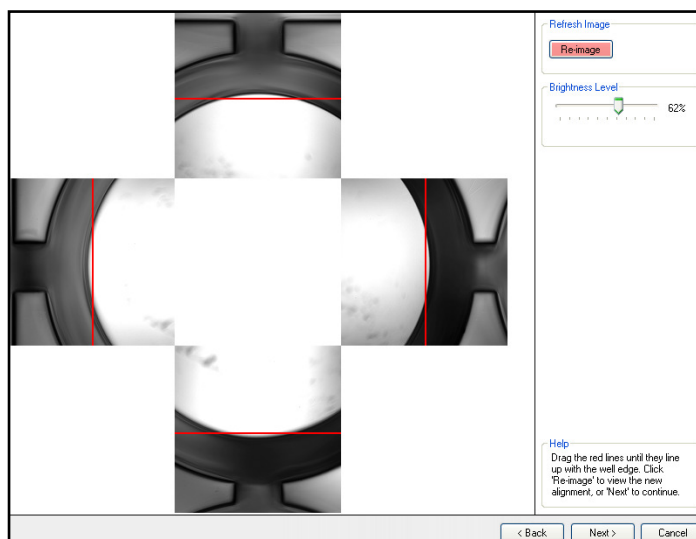
Scale Bar

A scale bar is displayed in the lower left of the window. This resizes automatically as the zoom level is altered.

Alignment

The **Alignment** window ensures the camera is properly aligned with the wells of the microplate.

Four images are taken over the north, south, east and west edges of well A1. The well edges should line up with the four red lines. If they do not line up, clicking and dragging the lines will correct the alignment.



Re-image

This button should be clicked if the **Re-image** button is highlighted in pink. This will recheck the alignment following adjustment.

Brightness Level

The brightness level of this image can be altered by moving the slider. **Note:** This will only change the brightness for this alignment image.

Imaging

During the imaging run, a schematic of the plate is displayed with the image frames superimposed in light blue. As images are captured and processed the frame changes color to light brown then green.

Once a frame has changed to green, it is possible to click on it to display the image in the View Image tab. Clicking on the **Confluence** tab will return the screen to the plate view.

Any unselected wells will not be imaged and are displayed as a blue square.



Results

The results of the **Image Microplate** process are displayed. The images can be reviewed here or they can be viewed via **Review Results** once the process has finished completely.

Remove Plate

When Imaging is complete the microplate will be ejected from the system. Within this screen there is the option to repeat the imaging with another plate or move onto the next step of ending the process.

Image Another Plate

Another plate can be imaged using the same configuration. This is achieved by clicking **Image Another Plate** instead of **Next**. The plate carrier remains in the ejected position and the **Select Source** screen is displayed ready to continue.

Finish

Clicking **Finish** will end the process.

The **'Image Microplate'** page will be displayed. Once returned to this page there are several options available.

'Save process' will save the current process to the **'Recent Processes'** tab for future access.

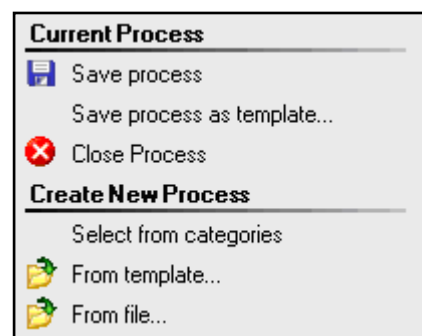
Note: A saved process can be changed.

'Save process as template...' will save the current process to the **'Template'** tab for future access.

Note: A saved template cannot be changed.

'Close Process' will close the process and return to the **Main Navigation Screen**.

Results can be viewed via **'Review Results'** after closing the process (see **Review Results** section).



There are also options to select and run a new process or a template from a saved location using the **'From template'** and **'From file'** locations.

Review Results

The **Review Results** process allows previous imaging runs to be reviewed and also to carry out further data processing, such as confluence, monitor cell growth, cell number estimation, monoclonality analysis and migration assay analysis.

When an **Image Microplate** process has completed the results are automatically displayed in a **Review Results** process.

Click on Review Results to start the process.

Load Results

The default location where imaging results are stored is the 'C:\Image Archive' folder. All the image and data files for each individual run are stored in a folder that is named with the date and time stamp for the run, e.g. 2005-11-01T11-52-10. Within the Image Archive folder, the run folders are automatically organized into separate folders for each month, e.g. 2006-10.

When opening the **Review Results** process on the CloneSelect Imager the user is directed straight to the results.

Plates imaged on the system are displayed as a list. If free space on the disk is running low, the bar at the top of the window appears orange and a warning is displayed in the top right corner. If disk space is getting critically low the bar is colored red.

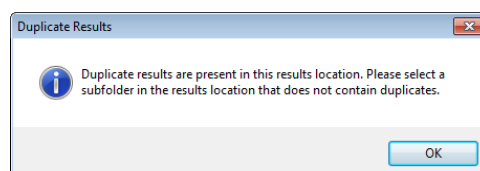
The screenshot shows the 'Load Results' window with a warning bar at the top: 'Warning: Only 21% Free Disk Space Available.' The main area contains a table with the following data:

Runs	Date	Barcode	Wellplate	Operator	Annotation	Disk Space
Archive	17/06/2011 13:53:28	006393	96 Well	Administrator	1 well has been imaged.	8.06 MB
Results	17/06/2011 13:50:37	003026	96 Well	Administrator	60 wells have been imaged.	423.59 MB
Finish	17/06/2011 09:31:55	003026	96 Well	Administrator	12 wells have been imaged.	84.45 MB

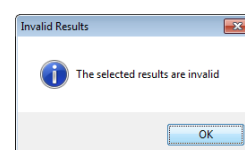
At the bottom of the window, there is a search bar with 'Find:' and 'In Field: All' dropdown, and a status bar showing 'Showing 29 of 29'. A context menu is open on the right side with options: View, Export, Archive, Delete, Rename, Results Selection, Results Location..., Other Applications, and Migration.

Right click on a selected run to display the context menu.

If there is more than one copy of the same data set present in the Image Archive folder, the following error message will be displayed when trying to open the data set. The duplicated data should be removed from the Image Archive folder.



If results have been deleted from the Image Archive folder, the data set will still appear in the Review Results list. If this data set is selected the following error message will appear.



Delete the data set from the Review Results screen to completely remove it.

Manage Results

VIEW RESULTS

To view results either double click on the data of interest or click to highlight the data and then click **View**.

Multiple data sets can be selected by holding down the **Shift** or **Control** keys while clicking the data of interest to select contiguous or non-contiguous sets of data, respectively.

EXPORT

Displays a wizard that allows data from the selected results to be exported. A variety of formats are available.

View results

Export results...

Archive results...

Reprocess results...

Delete...

Rename...

ARCHIVE

The results for runs stored in the default location (C:\Image Archive) can be archived to CD/DVD or the network/disk. Click **Archive** results to open the Archive window.

Note: Archive results is not available for results that are not stored in the default location. To move such results to other locations or burn to CD/DVD, use Windows Explorer.

Note: Archive to CD/DVD is only available with Windows XP.

Archive to CD/DVD

Select CD/DVD as the archive type.

Archive Name

Give the archive an appropriate name.

Media Type

Select the media type from the drop down list. An explanation of the media types is given at the right hand side of the screen.

Delete result images once written

If checked, the result files are deleted from the C:\Image Archive folder once the archiving has been performed. Select this if the archiving is being carried out to free space on the C Drive.

Archive

Click Archive to start burning the results to CD/DVD.

Archive to Network/Disk

Select Network/Disk as the archive type.

Select an archive type: CD / DVD Network / Disk

Archive to Network or Disk

Archive Location:

Remove local copy once move complete:

Help

Archiving to Network / Disk
 Select the location to which the data should be archived by clicking the 'Browse' button.
 To remove the data from the local archive after it has been copied to the new location, check the tick box.
 Click the 'Archive' button to begin archiving.

Archiving to CD / DVD
 Enter a name for the archive.
 Select a media type.
 CD-R for a writable CD.
 CD-RW for a re-writable CD.
 DVD5 for a single layer writable DVD.
 DVD9 for a dual layer writable DVD.
 To remove the stored image data from the computer when the disk has been written, check the tick box.
 Click the 'Archive' button to begin archiving.

Archive Location

Either type in the location for the archived files or click Browse to display a Save As Dialogue to select the location.

Remove local copy once move complete

If checked, the results files are deleted from the C:\Image Archive folder once the archiving has been performed. Select this if the archiving is being carried out to free space on the C Drive

Archive

Click Archive to send the results to the selected destination.

DELETE

Deletes all the images and results for the selected data set.

RENAME

Barcodes/plate names can be re-named in case of any misnamed plates which will subsequently be excluded from growth curve data. This can be achieved by clicking on the barcode/name to be changed and clicking **Rename**. A small dialogue box will appear in place of the barcode/name to enable a new barcode/name to be typed in its place.

Results Selection

RESULTS LOCATION

If results have been saved to a different location (not Image Archive) it is possible to browse to another location using this feature.

Other Applications

MIGRATION

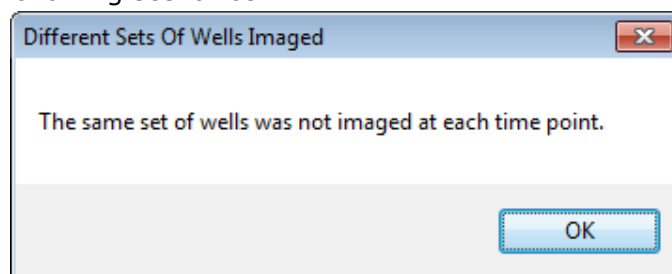
This option does not appear if the software has not been licensed for this application.

To view the data for a migration assay click on the relevant data set and click on the **Migration** link to open the software application.

The migration data analysis screen will open and the data is displayed within three separate tabs; Overview, Thumbnails and Well Data. When the data is opened the Thumbnails tab will be grayed out and display 'Loading...'. This allows extra time for the images to be loaded into this data tab. Meanwhile, data can be viewed in the other tabs.

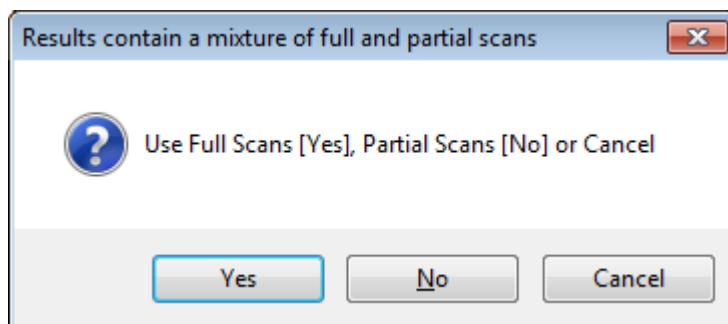
The Migration link is disabled and grayed out in the following scenarios:

- The selected results only have 1 time point.



The Migration link is enabled in the following scenarios:

- For a time point series, if different sets of wells have been imaged a message will appear informing that the migration application is unavailable and the Review Results screen will remain open to allow selection of other data.
- For a time point series, if there is a mix of Full and Partial scans with the same plate/barcode the following will happen:
 - If there is only one time point for each scan type, the migration option will not be applicable and will be grayed out.
 - If there is only one Full scan data set and more than one Partial data set, the migration will only be applicable to the Partial scan data and vice versa.
 - If there is more than one data set for both the Full and Partial scan a message will appear allowing the selection of the Full or Partial scans.



For more information on the Migration application see pages 38 - 44.

Analysing Results

Once an imaging process has completed the results are displayed in the **Results** window, which has several tabs: **Confluence**, **View Image**, **Cell Number**, **Loci Count**, **Plate Thumbnails**, **Growth Rate** and **Summary**. This window opens with the **Confluence** tab displayed by default.

The **Plate Type**, **Barcode** and **Processing Type** are all displayed above all the tabs to make it clear which data set is being viewed throughout the processing.

Confluence Tab

This tab displays the confluence level for each well, both graphically and as a list.

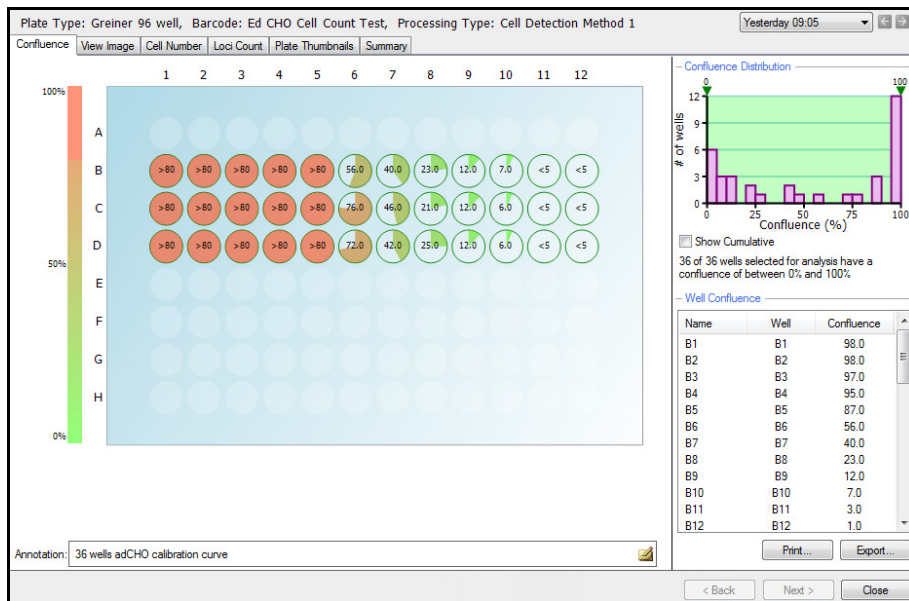
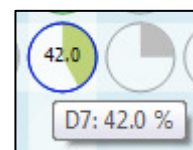


PLATE CONFLUENCE OVERVIEW

Each well is represented by a pie chart of the confluence level for that well. The pie charts are color-coded such that low confluence is green and high confluence is red with shades in between representing the intermediate levels.

Hovering the mouse over the pie chart area causes the percentage confluence for all the wells to be displayed. Confluence that is detected as lower than 5 percent is displayed as <5 and confluence above 80 percent is displayed as >80.



Hovering over a well causes a tool tip to appear displaying the well co-ordinate and the confluence for that well.

Clicking on a well will display the well image in the View Image tab. Right clicking on the overview will allow the overview to be copied to clipboard.

CONFLUENCE DISTRIBUTION

The bar chart at the top right of the window shows the number of wells with a given confluence level.

Dual Gates

The arrows pointing down at the top of the bar chart can be dragged across the chart. As both lower and upper arrows are dragged, the corresponding confluence data is displayed. Any wells that do not fall inside these lower and upper gates will be grayed out. The text displayed below the chart will update to give the number of wells that fall within the selected confluence (e.g. 10 out of 36 wells selected for analysis have a confluence of between xx % and xx %).

Show Cumulative

When checked the individual bars in the chart will merge to show the confluence as continuous data.

WELL CONFLUENCE

The confluence for each well is displayed in the list to the right. If desired the well name can be changed in this list by clicking twice on the well name of choice.

Print

This option will print a schematic presentation of the confluence overview in the plate format. The confluence values are displayed within the wells and the plate details are displayed at the top of the overview.

Export

This will enable the list of wells and the corresponding confluence to be exported as a .csv or .xml file. The Data Export wizard will guide the process of exporting the confluence data. Cell Number and Loci Count data can also be exported at this point (See the relevant sections for the description of exporting these data sets).

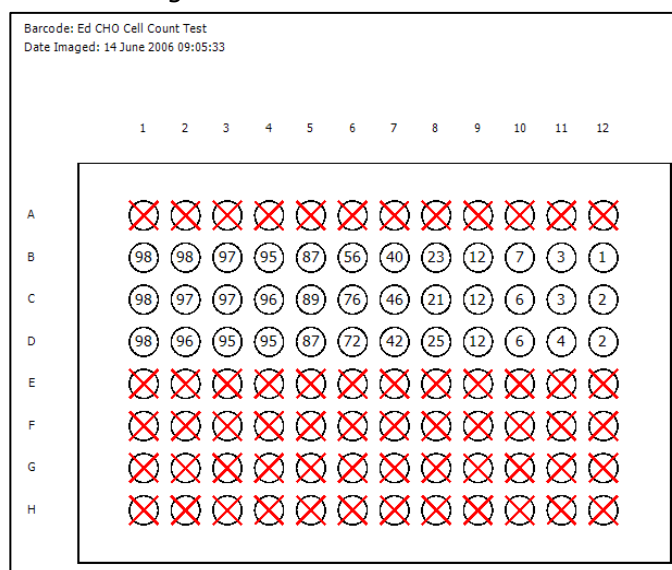
The wizard provides the option to export the data for the following time points:

Select which time points to export:

The currently selected time

The most recent time for each well plate

The complete time series



ANNOTATION

By default the **Annotation** text box displays the number of wells that have been imaged. This can be edited to more meaningful annotation. **Note:** This can not be changed when viewing the data through a remote data viewer.

View Image Tab

This tab displays the images for a selected well. The whole well is displayed.

WELL SCHEMATIC

By clicking on the image of the well currently being displayed, it is possible to navigate to different areas of the well and this is reflected on the image schematic to the right. The confluence levels of the currently selected whole well are displayed below the schematic.

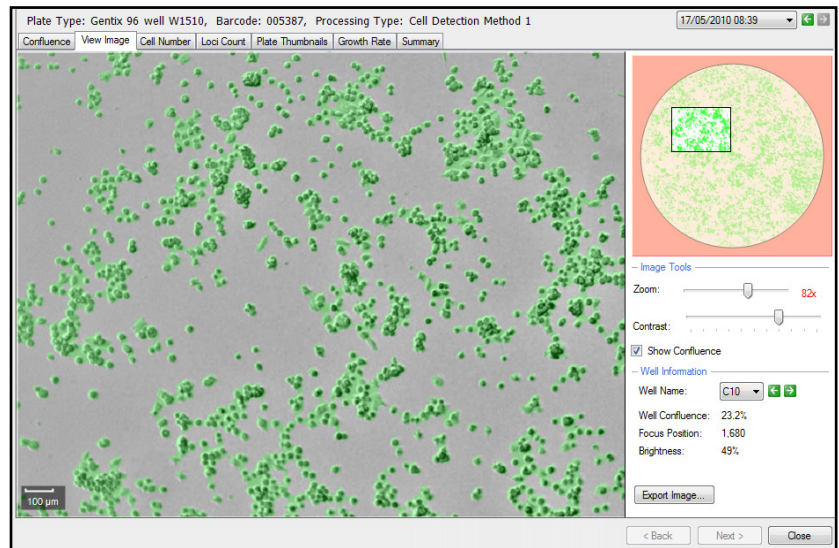


IMAGE TOOLS

Zoom

The slider can be moved either left or right to zoom out or in respectively. The lowest magnification of the image is 18x and the highest is 144x. When the figure turns red the system is zooming digitally and which may cause some pixilation of the image. When zoomed into an image, the zoomed area will be displayed on the image thumbnail to the right.

Contrast

The contrast of the image can be improved by moving the slider to the left or right.

Show Confluence

Checking this box will display the confluence detected within the well. Detected objects are displayed in green.

WELL INFORMATION

Well Name

Using the drop down menu or the arrows it is possible to toggle to the well of choice.

Well Confluence

The well confluence percentage is displayed here for the selected well.

Focus Position

The focus point of the image when captured is displayed here.

Brightness

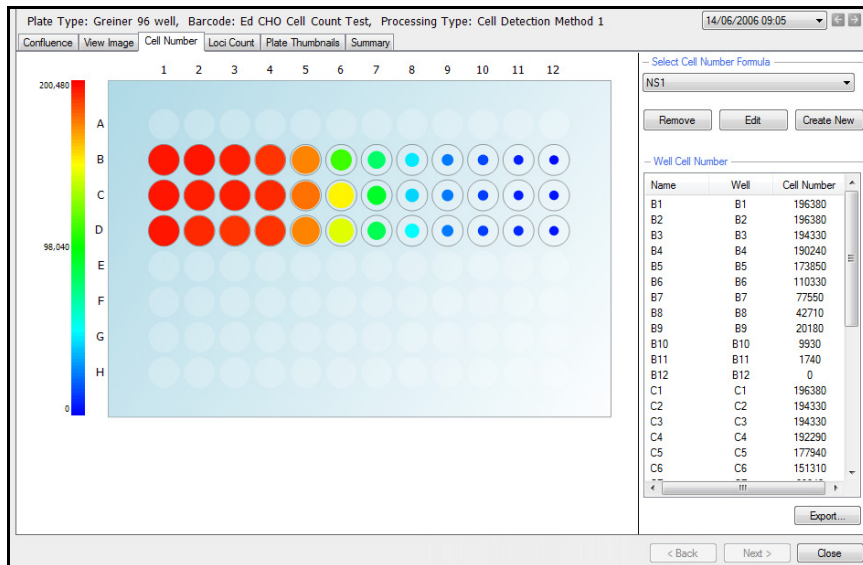
The brightness level used to capture the imaged is displayed as a percentage here.

EXPORT IMAGE

It is possible to the export the currently displayed image in .bmp, .jpg or .png format. If **Show Confluence** has been selected, the confluence overlay is displayed with the exported image. The zoomed position is saved within the image.

Cell Number Tab

This tab displays the estimated number of cells in each well, both graphically and as a list. The number of cells is estimated using a formula that can be created for each cell type from the confluence readings of a standard plate containing known numbers of cells. See **Appendix C** for a detailed procedure.



Each well contains a colored fill with the estimated cell number indicated by the color and size of the fill area, with a large red fill area representing a high cell number and a small blue fill area representing a low cell number. A color scale is shown to the left of the graphic.

Hovering the mouse over a well displays a tool tip giving the well co-ordinate and the estimated cell number for that well.

SELECT CELL NUMBER FORMULA

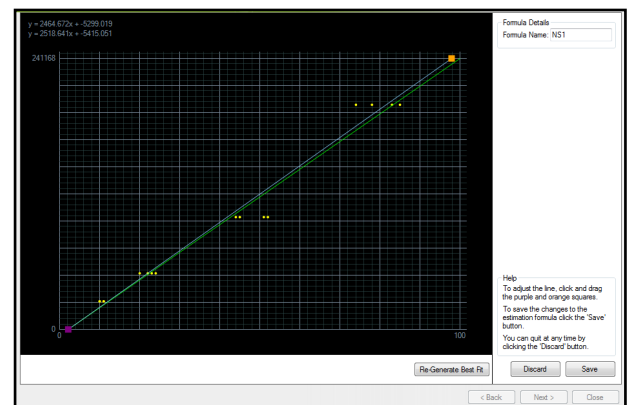
This drop down menu will list all the formulas created for estimating the number of cells per well.

Remove

Clicking this will remove the currently selected cell number formula from the drop down menu.

Edit

Clicking this will allow editing to be carried out on the currently selected cell number formula from the drop down menu. A graph of cell number against confluence for the formula is displayed. The handles at each end of the line can be dragged to edit the formula. The formula is displayed at the top left and the edited version is displayed immediately below it.



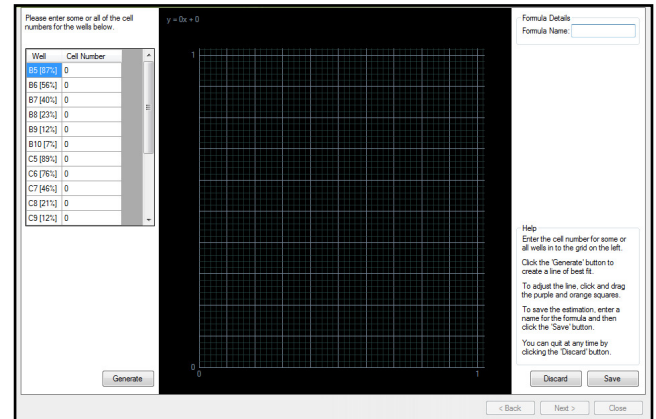
Discard can be clicked to abandon the edits and return to the **Cell Number** tab or **Save** can be clicked to save the changes made to the cell number estimation formula. The formula name can be changed in the top right hand box. There is also the option to **Regenerate Best Fit**. Clicking on this will automatically generate the line of best fit if required.

Create New

Clicking this will enable a new cell number estimation formula to be created.

Ensure that the confluence results displayed in the Confluence Tab are those for a standard plate containing known numbers of cells. The software will automatically select up to twelve wells with confluence values ranging between 10 % and 80 % and display them at the bottom of the graph.

The name for the new formula should be entered in the top right hand corner and the cell numbers for each well should be entered into the corresponding well number box to the left of the graph. Clicking Create after all relevant information is entered will create the new formula.

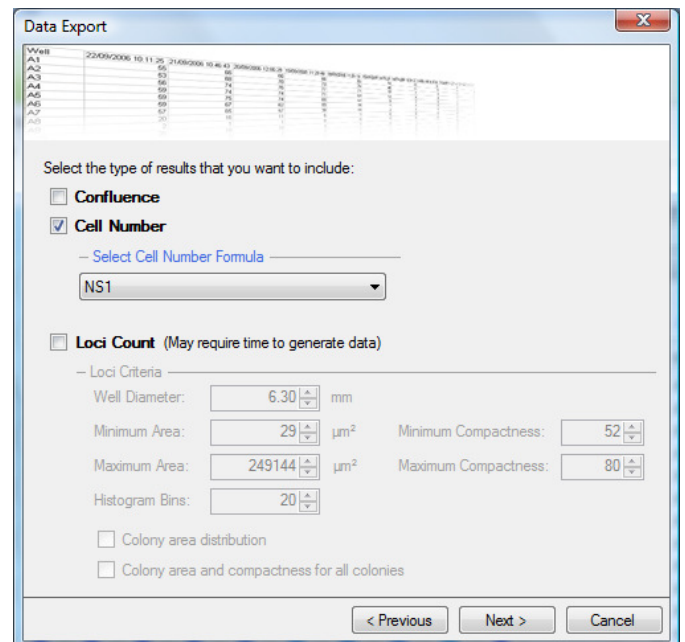


WELL CELL NUMBER

A list of estimated cell numbers is displayed to the right of the plate overview.

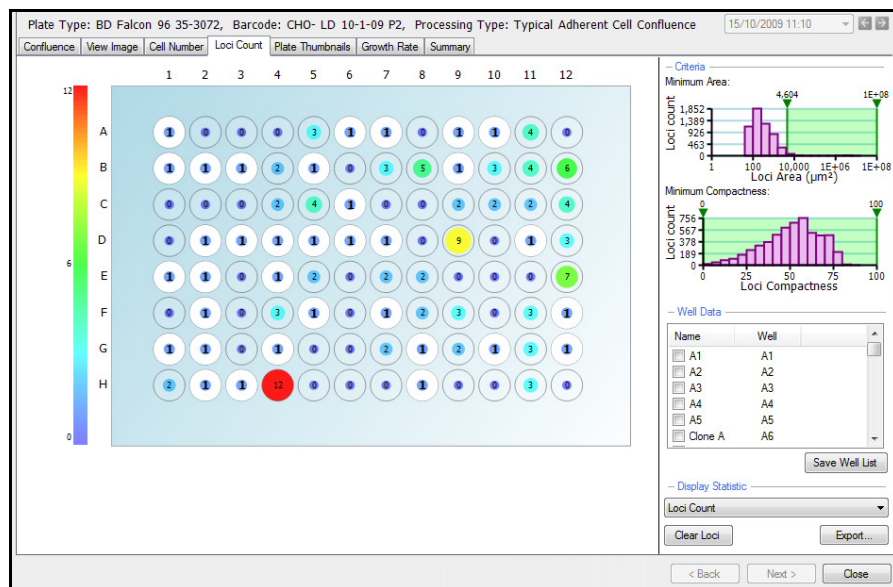
Export will enable the lists of wells and the corresponding Cell Number to be exported as a .csv or .xml file. The **Data Export** wizard will guide the process of exporting the Cell Number data.

Confluence and Loci Count data can also be exported at this point (See the relevant sections for this description of exporting these data sets).



Loci Count Tab

Generated images can be processed using the loci count feature. This detects and counts the number of loci of growth, i.e. cell colonies for applications including monoclonality verification and colony forming assays. If a plate is imaged multiple times during colony growth, the history of each well with one identified colony can be viewed for visual proof that the colony is derived from a single cell progenitor. For colony formation assays, the Loci Count feature can be used to count the number and size (area) of colonies in each well.



In the **Loci Count** tab, the plate view displays each well with a figure indicating the number of cell colonies found in the well. The default plate view displays the **Loci Count** (Under **Display Statistic**) however this can be changed to **Mean Loci Area** which will display this value within each of the wells.

To display the **Well View** for a well, click on the desired well in **Plate View** or in the **Well Data** list.

CRITERIA

Minimum Area

The loci count (frequency) is plotted against the loci area (μm^2) in a log scale. The bar chart has two, lower and upper, gates that can be moved accordingly to eliminate unwanted objects/debris within the well that artificially inflate the number of loci.

Minimum Compactness

The loci count (frequency) is plotted against the loci compactness. The bar chart has two, lower and upper, gates that can be moved accordingly to eliminate irregularly shaped objects from the loci count. Compactness is the relation between the area and the perimeter, expressed as a ratio of the actual area and that of a perfect circle with the same perimeter.

WELL DATA

This lists the wells. The well name can be changed if required by clicking twice on the well name and typing in the desired name.

Name	Well
<input checked="" type="checkbox"/> Clone A	A1
<input type="checkbox"/> A2	A2

DISPLAY STATISTIC

The drop down menu will display either the '**Loci Count**' or the '**Mean Loci Area**' on the overview.

The '**Clear Loci**' button allows the loci processing to be cleared. A prompt will appear warning that all the current loci data will be lost if continuing.

Export will enable the lists of wells and the corresponding Loci Count to be exported as a .csv or .xml file. The **Data Export** wizard will guide the process of exporting the Loci Count data.

Confluence and Cell Number data can also be exported at this point (See the relevant sections for this description of exporting these data sets).

The Loci Count criteria are automatically populated from the gated **Minimum Area** and **Minimum Compactness** graphs.

Histogram Bins: This option is used if the **Colony area distribution** box is selected. This value sets the number of equi-sized bins for exporting loci frequency counts. The default value is set to 20 bins.

Colony area distribution: When this box is selected the colony area distribution data is divided into the number of bins specified in the **Histogram Bins** and displayed with any other exported data.

Note: Only wells included/selected in the **Well Data** list are exported.

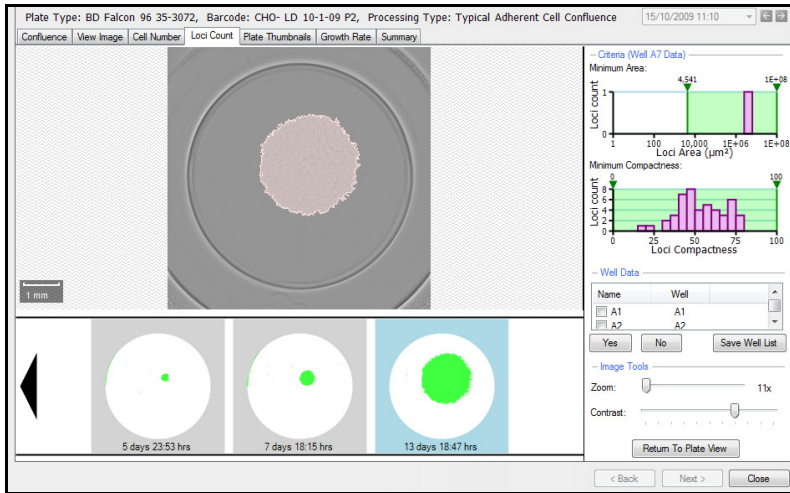
Colony area and compactness for all colonies: When this is selected a list of all colonies in all wells is exported with the area and compactness for each colony.

Note: There is no filtering so all the data is listed.

The screenshot shows the 'Data Export' dialog box. At the top, there is a table with columns for 'Well' and various data points. Below the table, the user is prompted to 'Select the type of results that you want to include:'. There are three checkboxes: 'Confluence', 'Cell Number', and 'Loci Count'. The 'Loci Count' checkbox is checked. Under 'Cell Number', there is a dropdown menu for 'Select Cell Number Formula' with 'NS1' selected. Under 'Loci Count', there is a section for 'Loci Criteria' with several input fields: 'Well Diameter' (5.90 mm), 'Minimum Area' (21210 μm²), 'Maximum Area' (100000000 μm²), and 'Histogram Bins' (20). There are also checkboxes for 'Colony area distribution' and 'Colony area and compactness for all colonies'. At the bottom, there are three buttons: '< Previous', 'Next >', and 'Cancel'.

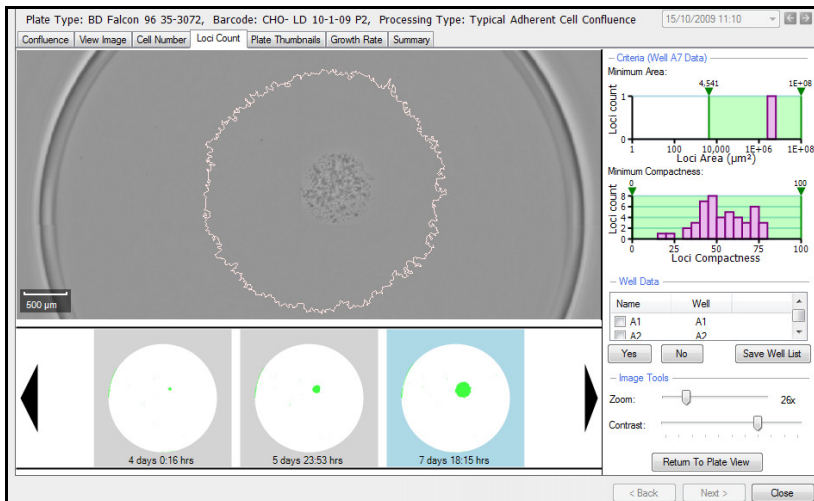
WELL VIEW

Clicking on the well of interest in the plate view of the **Loci Count Tab** will display the image for the well with an overlay over areas of growth detected according to the **Minimum Area** and **Minimum Compactness** criteria.



If the microplate has been imaged at several time points a film strip of schematics of the well at the various time points is displayed across the bottom of the screen with loci colored in green. Click on any of the schematics to display the well image at that time point with the loci from the latest image outlined.

This provides a means of checking if a colony has resulted from a single cell or from more than one cell. For an example of analyzing monoclonality data see the **Data Analysis Examples** section of this manual.



EXPORT IMAGE SEQUENCE

It is possible to export the sequence of images by clicking on the sequence with the right mouse button. The sequence can be saved as a .bmp, .jpg or .png file. The file will include the confluence percentage and time and day the image was acquired.

WELL DATA

It is possible to click on a well in the list to go to its corresponding **Well View** and check if the colony is monoclonal. There are three ways to move through the list and mark wells as monoclonal:

1. Click the Yes button to mark a well as monoclonal and scroll to the next well image. Click No to scroll to the next well without marking the well as monoclonal.
2. Click on a well to view its image and click on the check box to mark it as monoclonal.
3. Use the keyboard up/down arrow keys to scroll through the wells and the spacebar to mark the wells.

The selected wells in the list can be saved by clicking on the **Save Well List** beneath the list. This file can be saved as a .csv file. The file will display the plate barcode, run date, operator and the selected wells.

IMAGE TOOLS

Zoom

The slider can be moved either left or right to zoom out or in respectively. The lowest magnification of the image is 18x and the highest is 144x. When the figure turns red the system is zooming digitally and which may cause some pixilation of the image. When zoomed into an image, the zoomed area will be displayed on the image thumbnail to the right.

Contrast

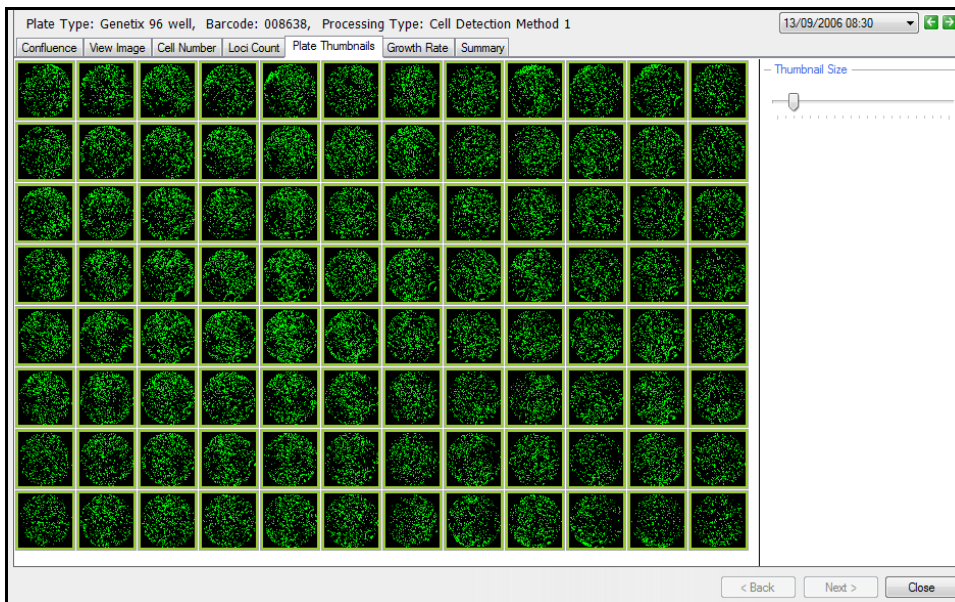
The contrast of the image can be improved by moving the slider to the left or right.

Return To Plate View

Clicking this button will return the screen to the plate view within the Loci Count Tab so other wells can be viewed.

Thumbnail Tab

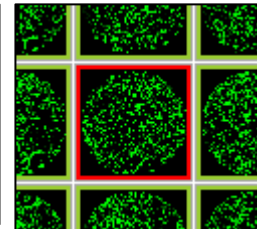
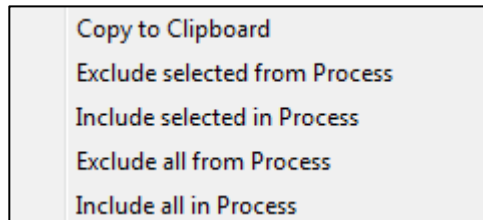
This tab displays the entire plate and the wells that have been imaged as thumbnails. The thumbnails display the confluence overlay of each of the imaged wells.



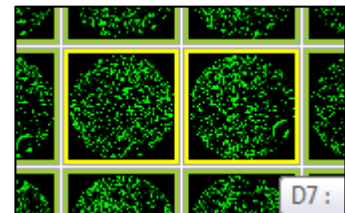
This tab allows wells to be included or excluded from analysis. This allows the analysis to be focused only on wells of interest. Any wells included or excluded from the **Thumbnails** tab will be updated in the **Confluence**, **Cell Number**, **Loci Count** and **Growth Rate** tabs as well.

This is achieved in the following ways.

- Right mouse clicking on the well of interest will display an options menu where a single or all wells can be excluded or included. Excluded wells are displayed with a red outline. Included wells are displayed with a green outline.



- Clicking the **Control** key on the keyboard and the left mouse button will allow several wells to be selected. These will be highlighted in yellow and the above menu will enable them to be included or excluded from the data set.



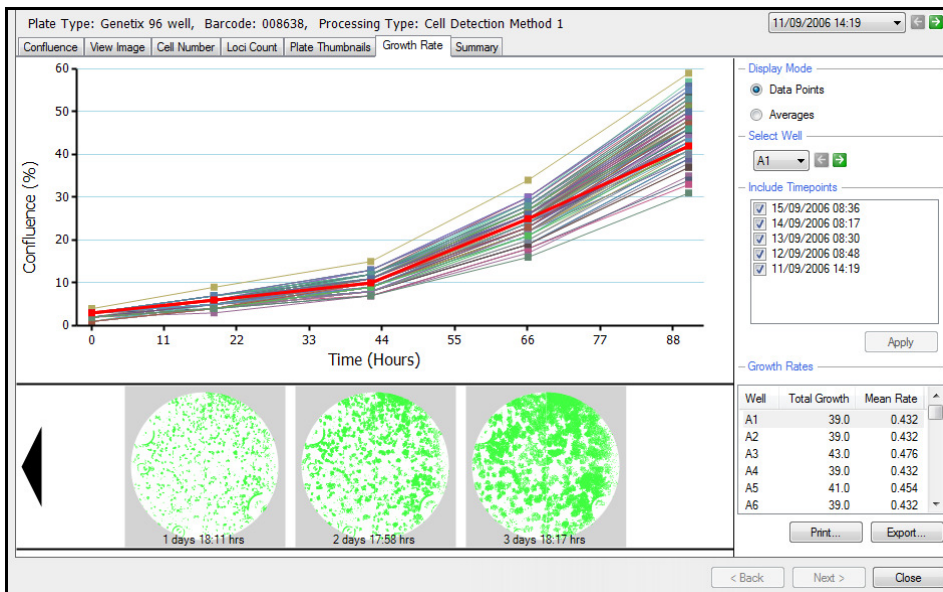
THUMBNAIL SIZE

This slider enables the size of the thumbnails to be increased or decreased.

Growth Rate Tab

If a microplate has been imaged previously, growth curves are displayed for all the wells in the plate. The system determines if the plate has been imaged by looking up its barcode in the stored results, so barcoded microplates should be used to enable this function. Named plates can also be used but the name must be typed in exactly the same each time the plate is imaged for the software to identify the series. The ability to rename plates also enables mis-named plates to be included in the growth curve data if required.

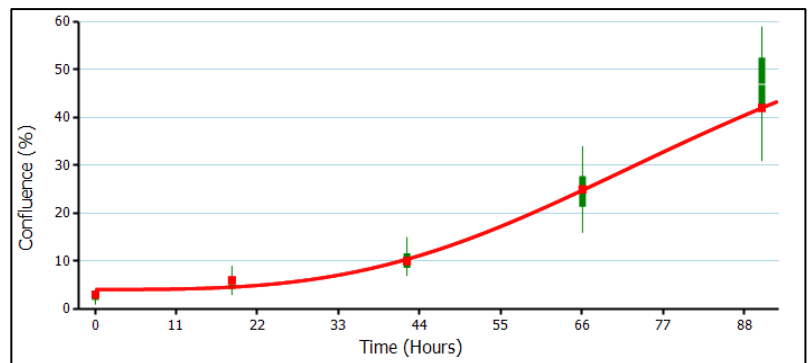
Hovering the mouse over each data point displays the well co-ordinate for that point.



DISPLAY MODE

The growth rate graph can be displayed with all the well data points by clicking the **Data Points** option or as averages by clicking the **Averages** option.

The thick red line within both graphs is displaying the growth rate for the selected well, e.g. A1. The well schematic below the graph is also displaying the captured images for the selected well.



SELECT WELL

This drop down menu allows the well to be changed.

INCLUDE TIMEPOINTS

This box allows various time points to be included or excluded from the growth rate output. This is achieved by selecting the time point to include or exclude and clicking **Apply**. The data will then either be included or excluded from the growth rate data. All plate scans with the same barcode will be automatically included in the growth rate data unless excluded at this point.

GROWTH RATES

This list displays the **Well**, **Total Growth** and **Mean Rate** for each well in the data set.

PRINT

This will print out a paper copy of the growth curves.

EXPORT

This will enable the list of wells and the corresponding confluence for each time point in the growth curve to be exported as a .csv or .xml file.

The **Data Export** wizard will guide the process of exporting the confluence data.

The wizard provides the option to export the data for the following time points:

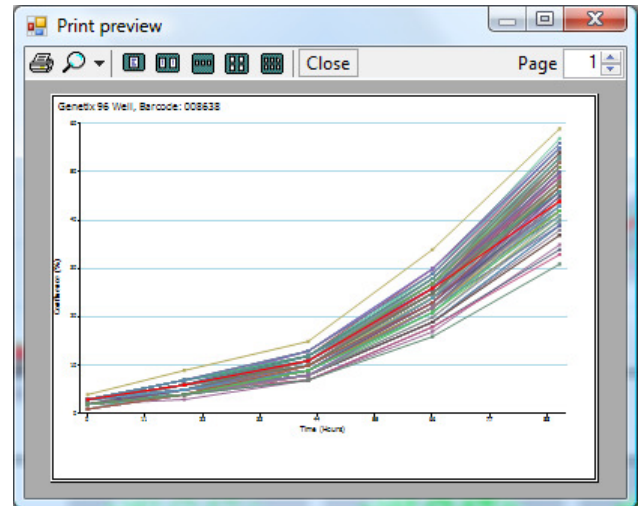
Select which time points to export:

The currently selected time

The most recent time for each well plate

The complete time series

Cell Number and Loci Count data can also be exported at this point (See the relevant sections for the description of exporting these data sets).



Data Export

Well	22/09/2006 10:11:25	21/09/2006 10:46:43	20/09/2006 12:48:28	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode	Barcode
A1	95	86	80																	
A2	95	86	80																	
A3	83	74	70																	
A4	89	74	70																	
A5	89	74	70																	
A6	87	74	70																	
A7	87	74	70																	
A8	87	74	70																	
A9	87	74	70																	
A10	87	74	70																	

Select the type of results that you want to include:

Confluence

Cell Number
 - Select Cell Number Formula:

Loci Count (May require time to generate data)

- Loci Criteria

Well Diameter: mm

Minimum Area: μm^2 Minimum Compactness:

Maximum Area: μm^2 Maximum Compactness:

Histogram Bins:

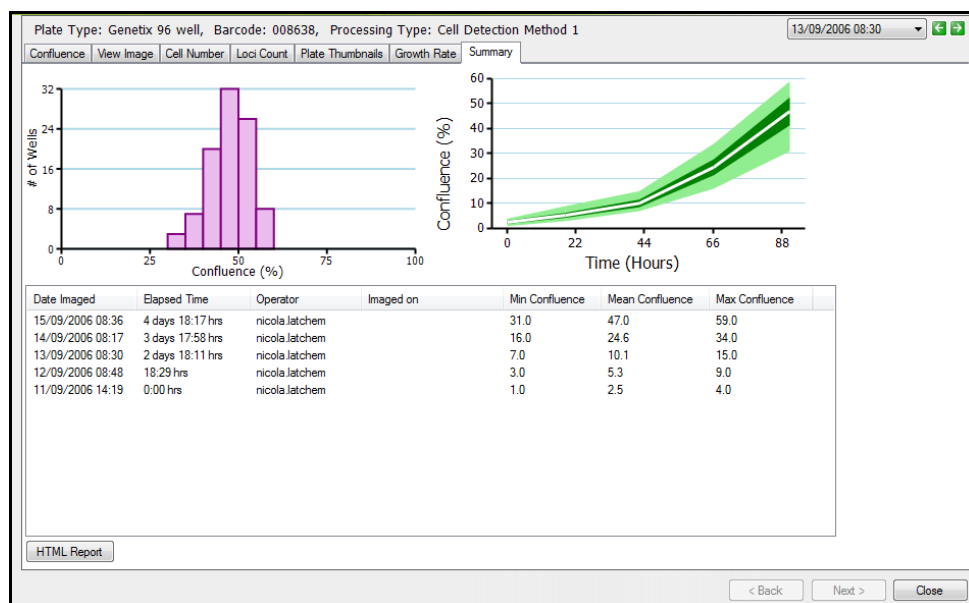
Colony area distribution

Colony area and compactness for all colonies

< Previous Next > Cancel

Summary Tab

This tab displays a summary of the confluence data, growth curve data and imaging credentials.



The table lists the Date Imaged, Elapsed Time, Operator, Imaged on, Min Confluence, Mean Confluence and Max Confluence for each time the plate was imaged.

HTML REPORT

This option enables a report to be generated containing the desired information in HTML format.

Various information, such as confluence, growth rate, cell number etc, relating to the data can be selected to be included in the report.

The loci data are automatically populated using the **Minimum** and **Maximum Area** and **Minimum** and **Maximum Compactness** data set in the **Loci Count Tab**.

Clicking **Generate** will require the report to be saved as a .html file before it is generated

HTML Report Wizard

Please select the sections that are to be included in the report:

- Confluence
- Growth Rate
- Cell Number

— Select Cell Number Formula —

NS1

- Plate Thumbnails
- Loci Count

— Loci Criteria —

Well Diameter: 6.90 mm

Minimum Area: 56 μm^2 Minimum Compactness: 0

Maximum Area: 161102 μm^2 Maximum Compactness: 100

Generate Cancel

EXIT

To exit Review Results click **Cancel** which will display the list of data in Review Results. Click **Cancel** again to exit this page, click **Finish** to close the process entirely and return to the Main Navigation Screen by clicking **Close Process**.

Migration Application

Migration analysis is carried out on data that has been imaged over a number of time points. For each data set analyzed the **Plate Type**, **Barcode** and **Processing Type** is displayed at the top of the screen.

Overview Tab

This tab provides an overview of the migration assay data. A heat map is generated and can be displayed using one of three parameters:

Migration Rate

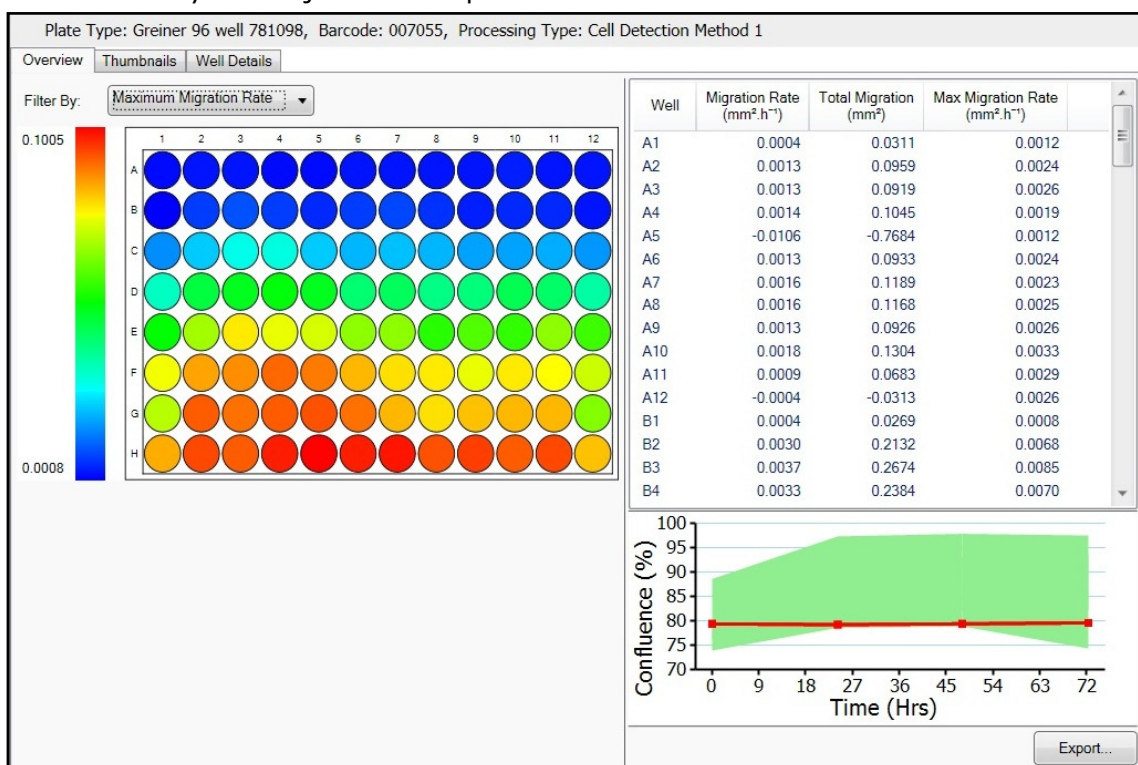
This is a measure of the rate of increase in cell movement (mm^2 per hour) over the entire time period.

Total Migration

This is a measure of the increase in area of cell movement (mm^2) over the entire time period.

Maximum Migration Rate

This is a measure of the largest rate of increase in cell movement (mm^2 per hour) between any two adjacent time points.



The table to the right of the heat map also displays these statistics for each well. Click on any of the table headers to reorder the data. It is possible to copy the heat map to the clipboard by right clicking anywhere on the map.

Click on any well of interest to highlight the well on the information table and on the Confluence (%) graph. Double click on any well of interest to hyperlink to the Well details tab.

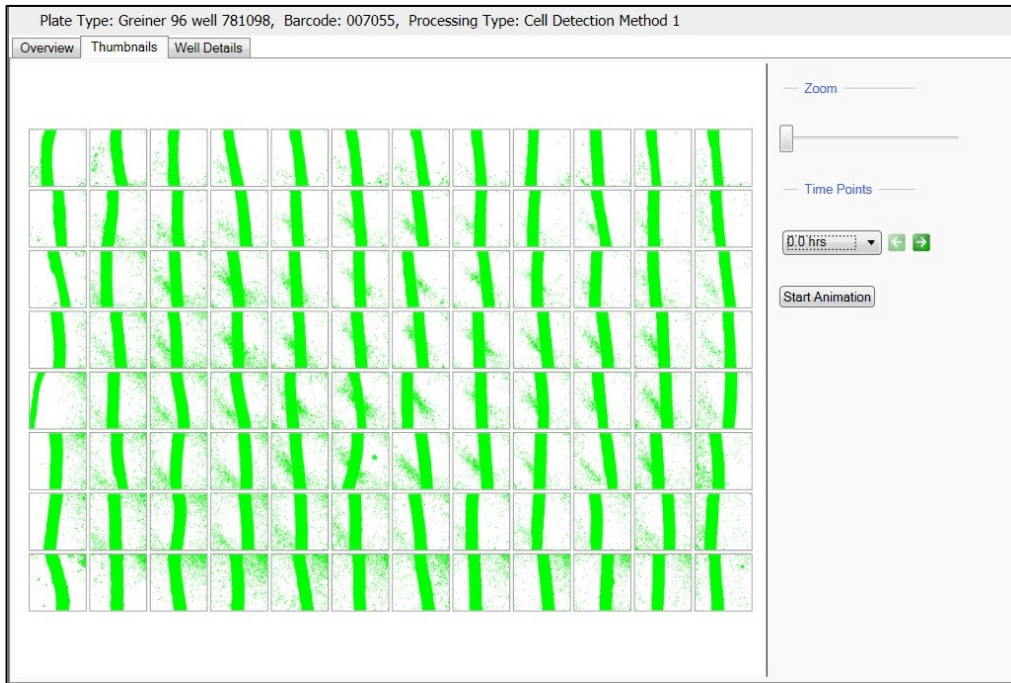
Export

The data for the well plate can be exported using this function. A wizard will guide the Data Export process. It is possible to export data for either Selected Time Points or for All Time points. It is also possible to export the confluence data if required. The data can be

exported as a CSV, Excel (XML) or XML file.

Thumbnails Tab

This tab provides an overview of each well image and the scratch/wound detection. The thumbnails can be copied to the clipboard by right clicking anywhere on the thumbnail map.



Zoom

Sliding the bar will enlarge/reduce the entire thumbnail map allowing navigation to any well of interest. It is possible to click and drag on the thumbnail map to scroll to the well of interest.

Time points

The drop down menu lists all of the time points the plate has been imaged. A time point can be selected from the drop down menu or by clicking on the arrows to the side of the menu. If the time points available are restricted by the End Time Point in the Graph display of the Well Details tab, only the selected time points will be displayed here.

Start Animation

Clicking this button will animate the thumbnails from the starting time point to the final time point in order to observe the migration taking place in the wells. If time points are restricted by the End Time Point option in the Graph display of the Well Details tab, animation will only occur using the time points available.

Well Details Tab

This tab provides all the details generated from individual wells. There are two displays available within this tab; Image Display and Graph Display.

IMAGE DISPLAY:

This display shows all the information gathered from the individual well images. A whole well image is displayed center screen with time point thumbnails below. Click on any of the thumbnails in the film strip to display the enlarged image for that time point.

The Display Graph header will change the view to the Graph view of the data.

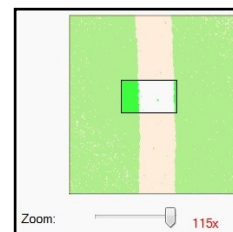
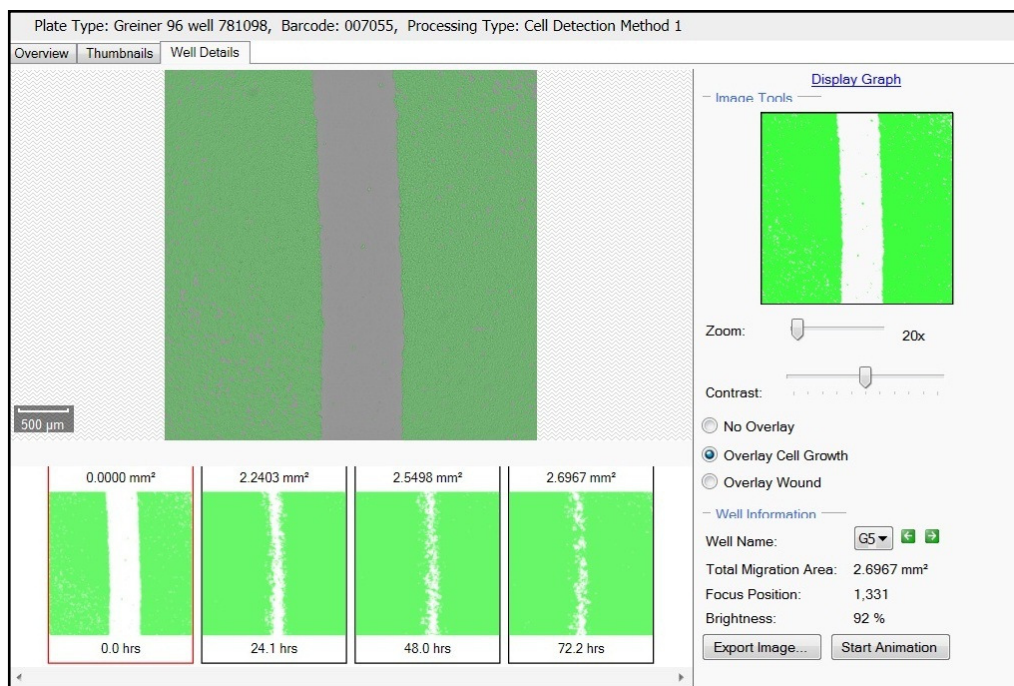


Image Tools

Zoom

Using the slide bar it is possible to zoom into any region of the well.

When zoomed into the image it is possible to click and drag to anywhere on the image. A small zoom box will appear on the navigation panel to identify the image region being viewed.

Contrast

It is possible to increase and decrease the contrast of the image using the slide bar, if required for better visualization of the cells.

No Overlay

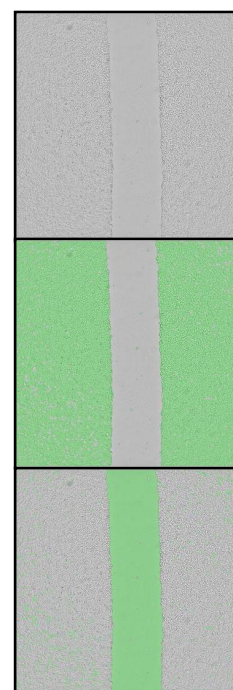
When selected no detection of any part of the image (cells or wound) is displayed.

Overlay Cell Growth

When selected only the detection of the region of cells is displayed in green.

Overlay Wound

When selected only detection of the regions without cells is displayed in green.



Well Information

Well Name

The displayed well can be selected from the drop down menu or by using the arrows to scroll through wells.

Total Migration Area

This value is a measure of the total cell growth in mm^2 over the entire time period for the selected well.

Focus Position

This is the focus position used for the microplate.

Brightness

This is the brightness value used for imaging.

Export Image

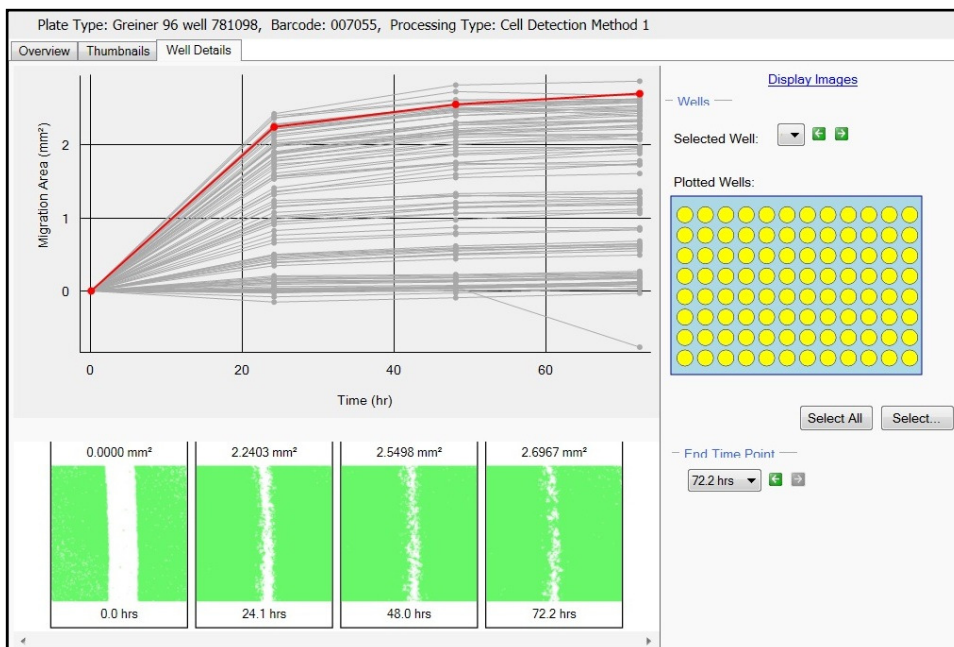
It is possible to export the well image as a .bmp, .jpeg or a .png.

Start Animation

If this option is selected each of the images in the time point series will be displayed as an animation to aid visualization of the migration of cells. The animation will always start from the first time point in the series. If time points are restricted by the End Time Point option in the Graph display of the Well Details tab, animation will only occur using the time points available.

GRAPH DISPLAY:

This display shows the information gathered from the individual wells and displays in graph format. The graph is displayed center screen with Migration Area (mm^2) plotted against Time (hours). A time point thumbnails of the selected well (denoted by a red line on the graph) is displayed below the graph.



Each thumbnail image in the film strip displays the total migration at that time point and time passed since the first time point in the series. When hovering over a data point on the graph for the selected well (denoted by a red line on the graph), the corresponding thumbnail will be highlighted in red.

The Display Images header will change the view to the Images view of the data.

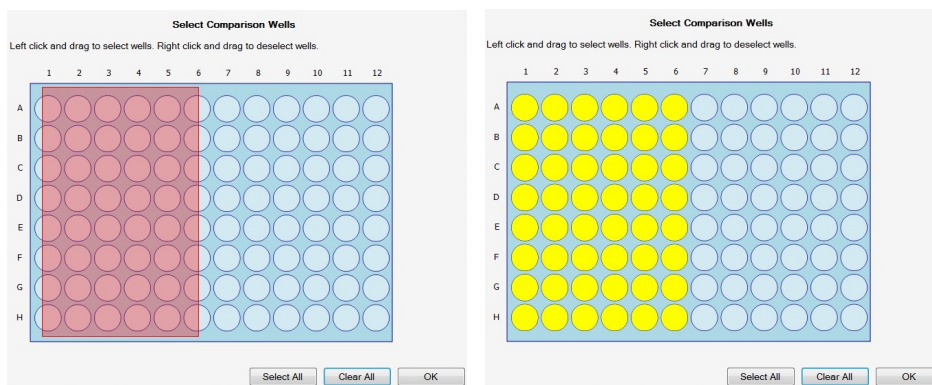
Wells

Selected Well

This is the well currently selected and displayed on the graph as a bold red line.

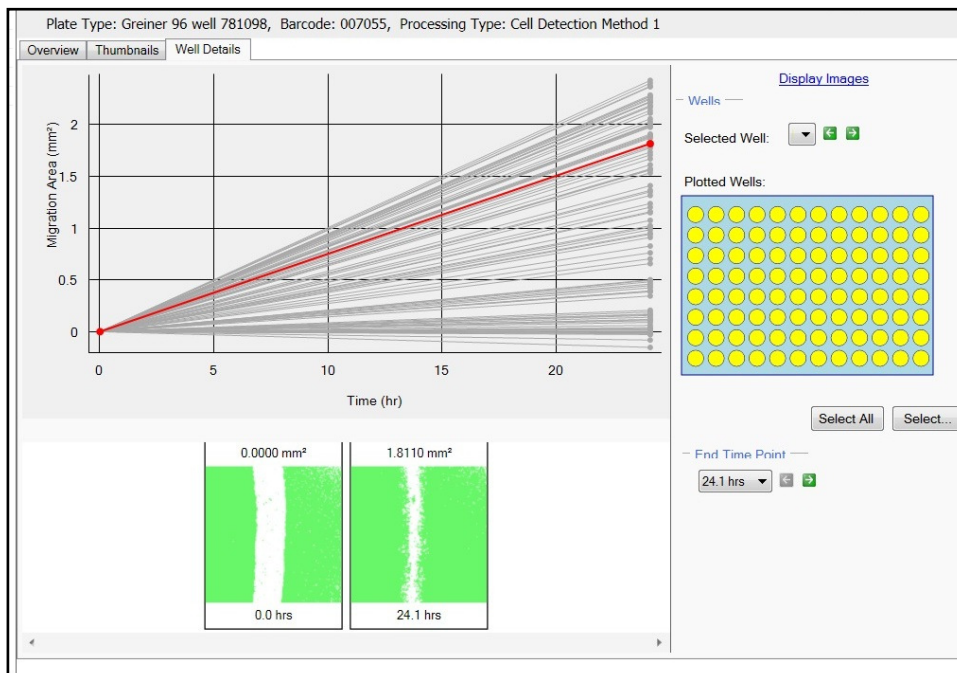
Plotted Wells

This plate map shows all of the wells that were imaged and are plotted on the graph. Any wells that are not imaged will be grayed out. Any wells can be included or excluded from this data by clicking on the **Select All** or **Select...** buttons. Wells can be selected/deselected using the left and right mouse buttons respectively. When wells are selected they are displayed in yellow. When wells are deselected they are displayed in blue. This feature enables the selection of wells to plot on the graph. This selection/deselection of wells only applies to the graph and not the overview.



End Time Point

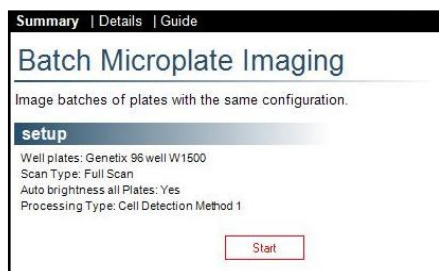
This drop down menu lists the time points of the imaged plate. It is possible to select the desired end time point directly from the menu or by scrolling through the time points using the arrows. When the end time point is selected the graph will display only the time points prior to and including the end point selected.



Note: When changing the End Time Point, only the selected time points will be displayed on the Thumbnails tab. Data on the Overview tab is also recalculated using the selected time points.

Batch Microplate Imaging

This process enables multiple plates to be imaged using the same configuration.



By clicking on the **Setup** header the appropriate plate type, scan type and auto brightness option can be selected.

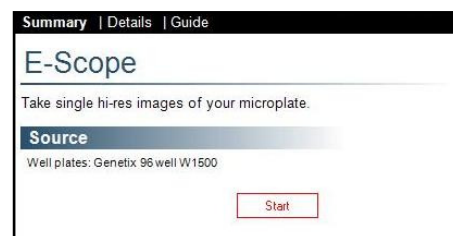
The Auto brightness all plates option automatically selects the appropriate brightness for the image. This can also be manually changed.

This process automatically looks for barcodes on plates but if these are not present a prompt will appear asking to manually insert a plate barcode/name. Once **Start** is clicked the process is the same as **Image Microplate**.

E-Scope

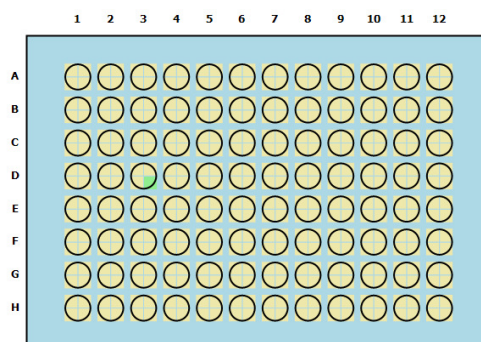
The **E-Scope** process provides a means of capturing higher resolution images of cells (1.85 μm) without performing any image analysis. Frames to be imaged can be freely selected by clicking with the mouse on the screen.

The setup of the **E-Scope** process is the same as **Image Microplate**.

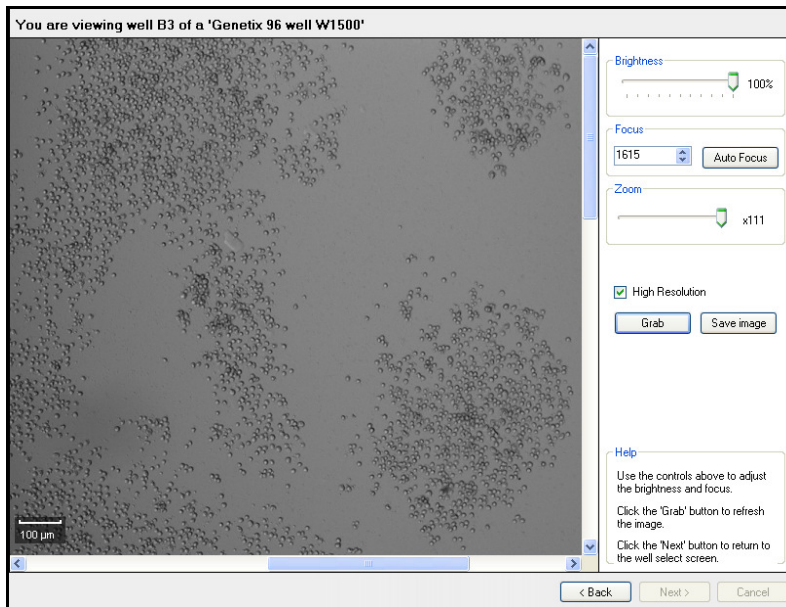


How to Image Using E-Scope

- Select the microplate type to be imaged by clicking on the Source Header and selecting the appropriate plate from the drop down list. Click **Apply** to set the changes.
- Click **Start** to commence the process.
- Load the plate making sure well A1 is placed to the back right corner of the plate carrier. Click **Next** to load the plate. **Cancel** can be clicked to exit the process at any point.
- When prompted, check that the camera is properly aligned with the wells of the microplate. The brightness levels of the image can be changed if required by moving the slider. Ensure the **Re-Image** button is clicked if any changes are made to recheck the alignment. Click **Next** once satisfied with the image and well alignment.
- The system displays a schematic of the microplate showing all the possible image frames.
- Clicking on a frame will prompt the system to capture and display a real time image.



- The brightness and focus of the image can be altered at this stage. **Grab** must be selected to set any changes made.



- The high resolution box can be selected to toggle between the high resolution (1.85 μm) and standard resolution (3.7 μm) imaging modes.
- If required the image can be exported as a .bmp, .jpg or .png file using the Export button.
- To select another image or well go back to the plate view screen by clicking the back button.
- To exit click the **Next** button on the plate view screen. This will eject the plate for it to be removed and the process is finished. Click **Finish** to exit the entire process and return to the E-Scope process start page.

Plate Alignment

This process enables the alignment of the plate within the system. This process is carried out in the same manner as **Alignment** within the **Image Microplate** process but as a stand alone process and no subsequent imaging.

Eject Plate Carrier

The **Eject Plate Carrier** process provides a means of ejecting the plate carrier for cleaning or removing a plate if it has been left in the instrument.

Clicking **Start** will eject the plate holder. Clicking **Next** will return the plate carrier into the instrument.

Click **Cancel** to exit the process and return to the starting page of the **Eject Plate Carrier** process.

Click **Finish** to return to the starting page of the **Eject Plate Carrier** process.

Data Analysis Examples

Here are some examples of how to use the features of the CloneSelect Imager software to analyze various data sets.

Monoclonality Workflow

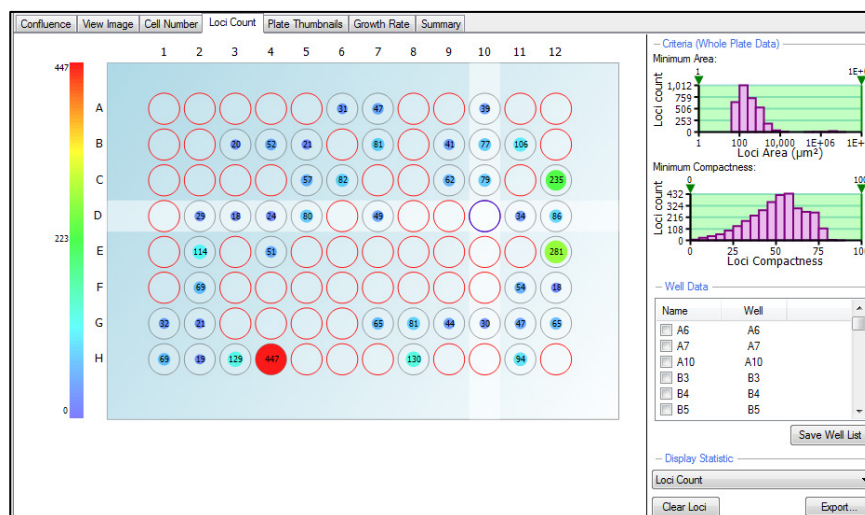
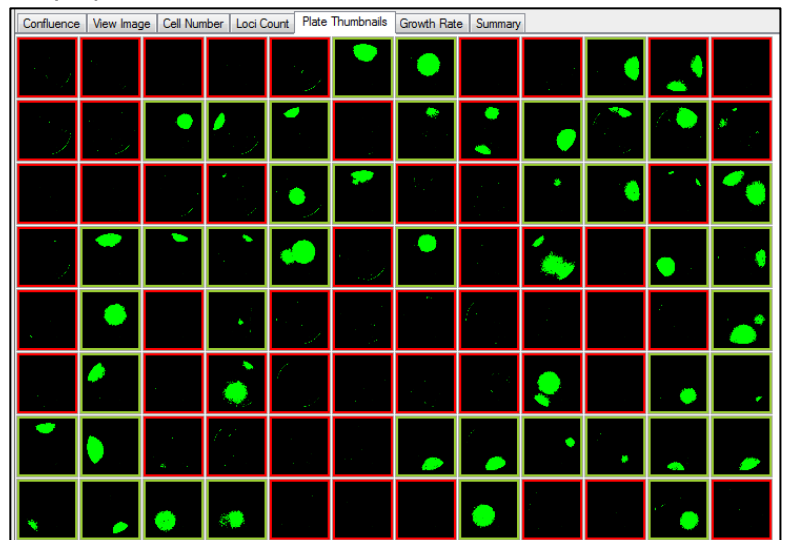
→ After imaging the plate of interest at several time points open the latest data point in Review Results. The software automatically opens the data in the first tab; Confluence.

→ Select the Thumbnails tab to display all the imaged wells. Exclude any wells that do not appear to have 1 colony within the well. (See the Plate Thumbnails tab section on how to exclude wells).

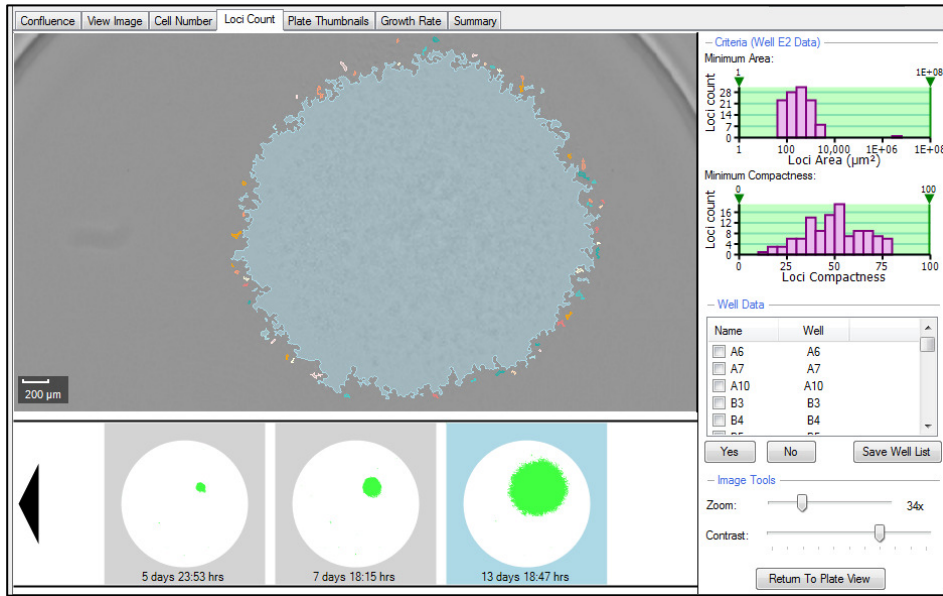
→ This will save time by not processing wells that do not contain 1 colony.

→ Open the Loci Count tab. The well diameter can be adjusted if required.

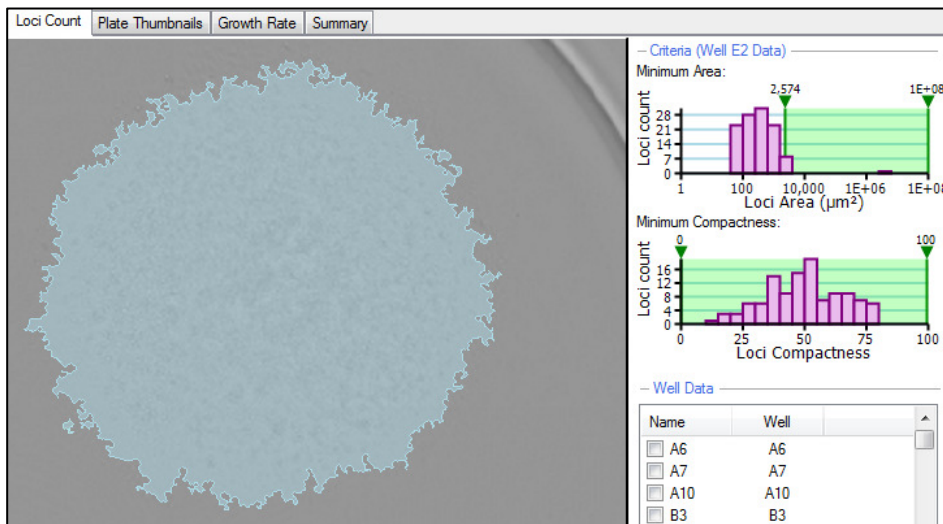
→ Click Process to produce the loci counts for the selected wells. The Loci Count tab screen will change to display an overview of all the selected wells and their corresponding loci count.



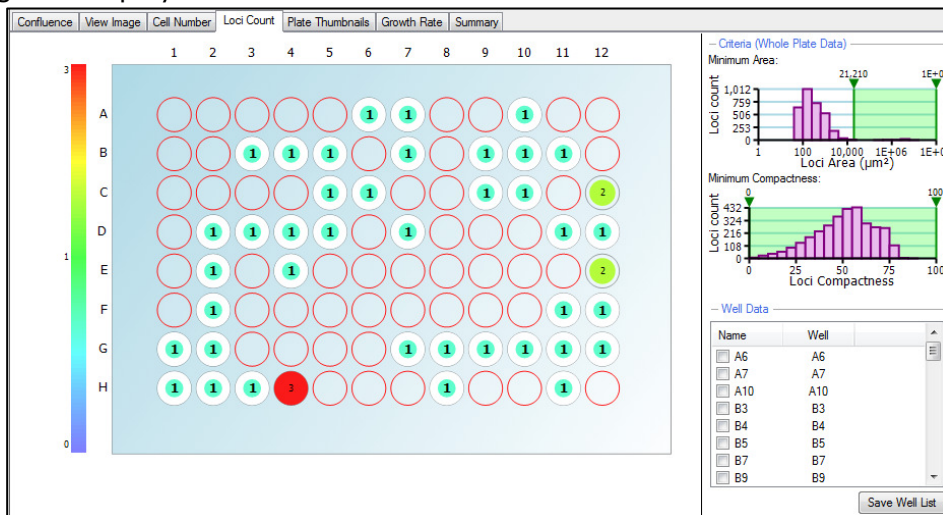
→ Click on one of the wells to display the well image. The image will show the loci processing.



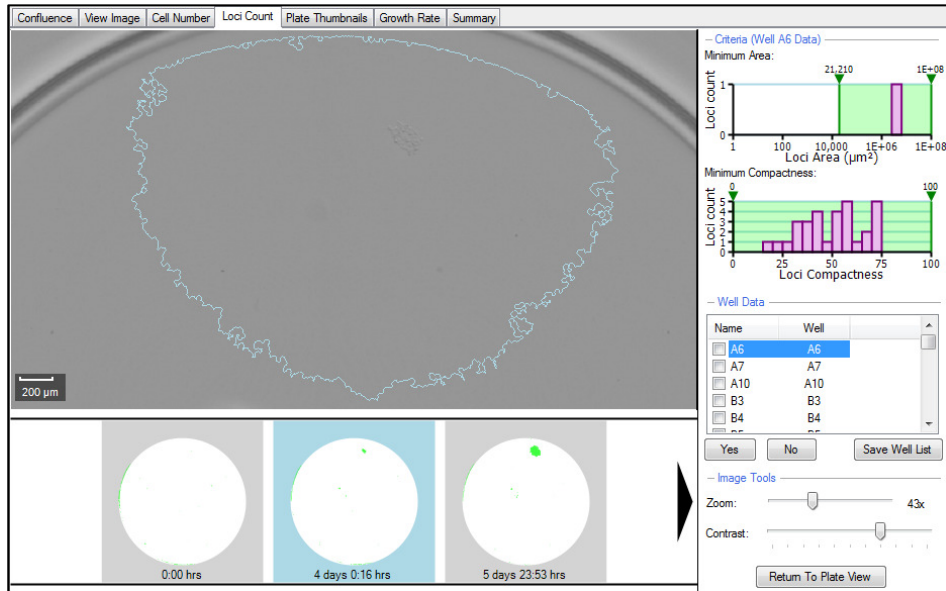
→ Move the Minimum Area and Minimum Compactness lower and uppers sliders on the graphs to gate out any unwanted debris and misshaped colonies within the wells.



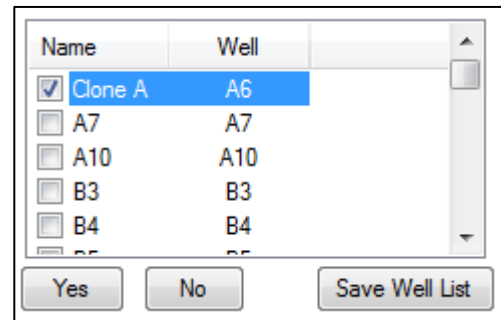
→ Click on Return To Plate View and check through some of the other wells to make sure all debris and misshapen colonies are excluded. The plate overview will reflect the processing and display the new loci count number in each of the wells.



→ Now by clicking through each well in the Well Data list, the well image will appear and mono-clonality can be determined by looking back through the series of well images and identifying if the colony originated from a single cell.



→ If the colony did originate from a single cell the well can be selected by either ticking the box or clicking the Yes button. The well name can also be changed to reflect the clone if required. Once all the wells have been checked, the list of selected wells containing a mono-clonal colony can be saved and printed for further use.



Colony Forming Assay Workflow

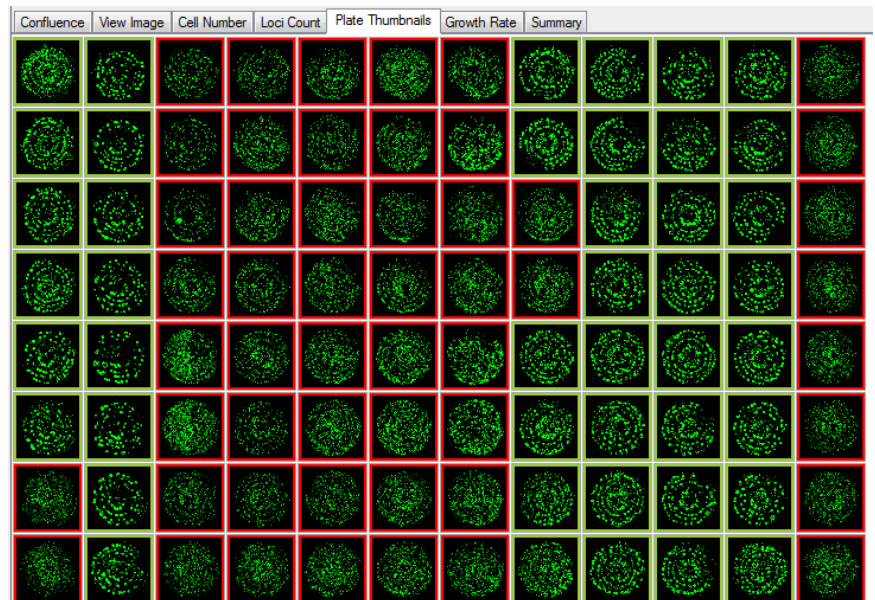
Analyzing colony forming assay data sets are very similar to gating monoclonality data, however the data export can be more detailed.

→ After imaging the plate of interest open the data point in Review Results. The software automatically opens the data in the first tab; Confluence.

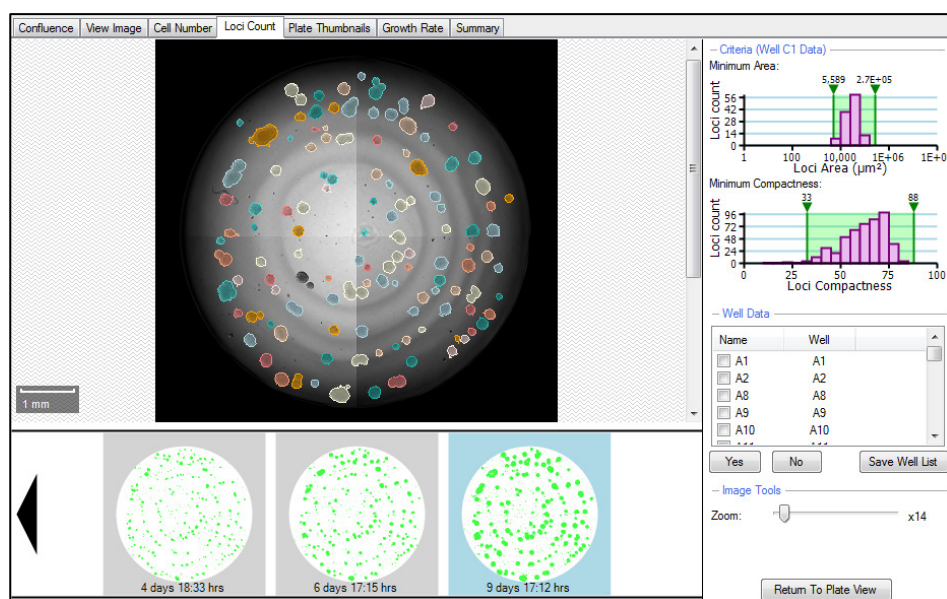
→ Select the Thumbnails tab to display all the imaged wells. Exclude any wells that have not formed colonies (See the Plate Thumbnails tab section on how to exclude wells).

→ Open the Loci Count tab. The well diameter can be adjusted if required.

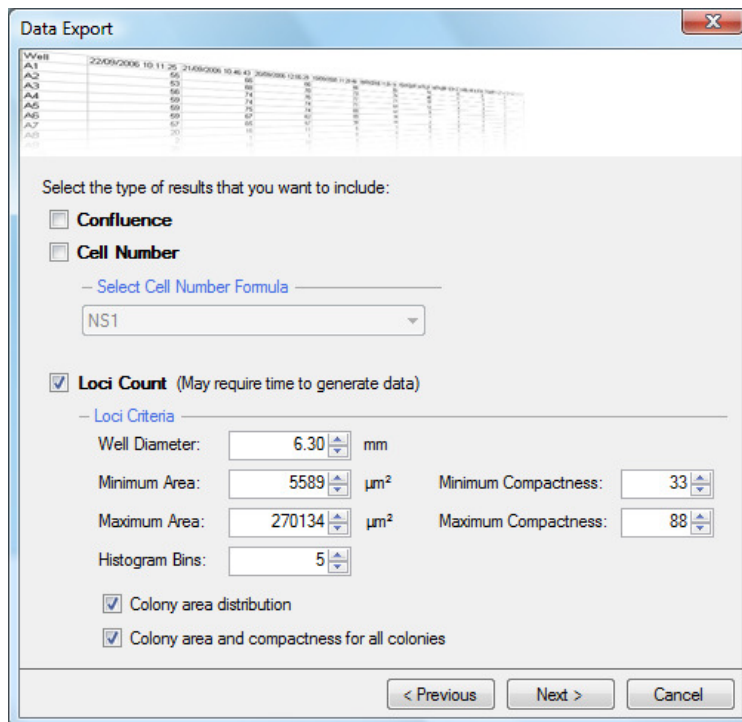
→ Click Process to produce the loci counts for the selected wells. The Loci Count tab screen will change to display an overview of all the selected wells and their corresponding loci count.



→ As above in the monoclonality work flow, click on a well to show the well image and gate out debris and misshapen colonies using the lower and upper sliders on the Minimum Area and Minimum Compactness histograms.



- Once satisfied with the gating of the data and the loci count subsequently produced, the data can be exported. The Well Data list is mainly used for accepting wells based on monoclonality so it is not necessary to check any of these wells. The data will be exported for all wells when exporting the data via the wizard.



- In the Data Export wizard the Loci Count option is selected and the criteria are automatically populated. The default number of bins is 20, but in the example it has been changed to 5 bins. The Colony area distribution and Colony area and compactness for all colonies are selected and Next is clicked to finish exporting the data.

Microsoft Excel - Colony Forming Assay Results.csv

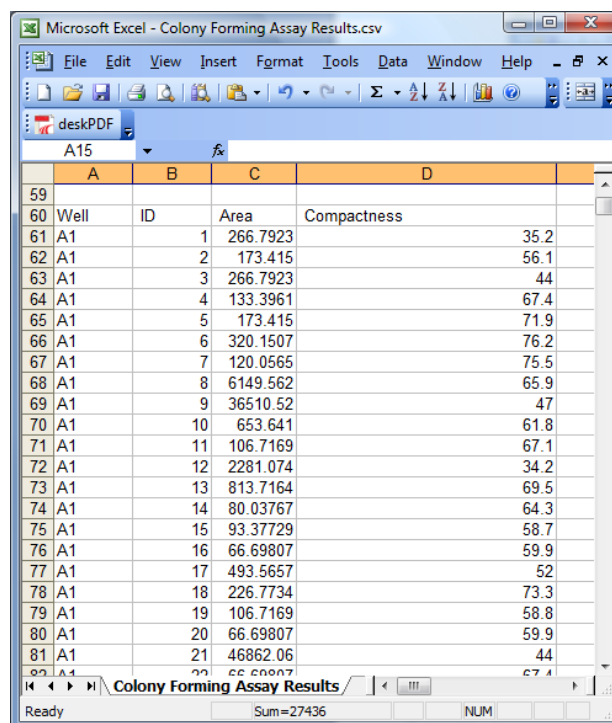
Colony Forming Assay Results									
Experiment Criteria									
1	Barcode=Exp 8								
2	PlateType=Grainer 96 Well								
3	WellRows=8								
4	WellColumns=12								
5	Operator=Andrea Gough								
6	RunDate=29/10/2007 09:12:47								
7	Annotation=96 wells have been imaged.								
8	LociWellDiameter=6.30								
9	LociMinimumArea=5589								
10	LociMaximumArea=270134								
11	LociMinimumCompactness=33								
12	LociMaximumCompactness=88								
13									
14	Name	Well	Loci Count	Mean Loci Area (microns squared)	Count 5589.00-58498.00	Count 58498.00-111407.00	Count 111407.00-164316.00	Count 164316.00-217225.00	Count 217225.00-270134.00
15	A1		129	27178	119	10	0	0	0
16	A2		85	41978	67	15	3	0	0
17	A8		105	41712	83	19	3	0	0
18	A9		121	34479	105	16	0	0	0
19	A10		71	45914	54	15	1	1	0
20	A11		99	38229	82	13	3	1	0
21	B1		118	32922	103	13	2	0	0
22	B2		65	49658	45	17	2	1	0
23	B8		92	44808	71	15	6	0	0
24	B9		108	39249	91	16	1	0	0
25	B10		61	45797	42	18	1	0	0
26	B11		79	45865	56	21	1	1	0
27	C1		116	34832	103	11	2	0	0
28	C2		89	50474	49	13	4	2	0

- Within the exported results produced, a summary of the experiment criteria is displayed in the first rows the spreadsheet (top left).

When the Colony area distribution is selected, the following information is produced for each well:

- Name (if changed within the software)
- Well number (eg A1)
- Loci Count (this is the total)
- Mean Loci Area
- Bins equally splitting the loci area range into equal groups so that each identified loci will fall into one of the corresponding groups based on its size. For example, out of the 129 identified loci within well A1, 119 of these fall into the first bin with a loci area range of $5589\mu\text{m}^2$ to $58498\mu\text{m}^2$.

When selecting the **Colony area and compactness for all colonies**, the area and compactness of each identified loci is displayed:



	A	B	C	D
59				
60	Well	ID	Area	Compactness
61	A1	1	266.7923	35.2
62	A1	2	173.415	56.1
63	A1	3	266.7923	44
64	A1	4	133.3961	67.4
65	A1	5	173.415	71.9
66	A1	6	320.1507	76.2
67	A1	7	120.0565	75.5
68	A1	8	6149.562	65.9
69	A1	9	36510.52	47
70	A1	10	653.641	61.8
71	A1	11	106.7169	67.1
72	A1	12	2281.074	34.2
73	A1	13	813.7164	69.5
74	A1	14	80.03767	64.3
75	A1	15	93.37729	58.7
76	A1	16	66.69807	59.9
77	A1	17	493.5657	52
78	A1	18	226.7734	73.3
79	A1	19	106.7169	58.8
80	A1	20	66.69807	59.9
81	A1	21	46862.06	44

Health and Safety

Transport and Storage

CloneSelect Imager must be stored and transported in temperatures within the range -25°C to $+55^{\circ}\text{C}$.

Lifting Points

CloneSelect Imager is intended for bench top operation. Care should be taken when moving CloneSelect Imager. It is recommended that 2 people lift it.

External Covers



Warning – If any of the external covers on CloneSelect Imager are removed the power supply is not automatically interrupted. If it is necessary to remove any of the external covers you must ensure that the power is switched off first and do not attempt to use the robot until the covers are replaced.

Electrical Safety



CloneSelect Imager must be connected to a properly earthed power outlet to protect users from the risk of electric shock. The main chassis of the machine is earthed together with all associated electrical components. Do not remove any of the fixed covers, as there are no user serviceable parts inside. All internal work should be referred to Molecular Devices approved service personnel.

Drive Safety



There is a potential pinch hazard with the drive mechanism. Please ensure that you do not attempt to load the plate until the drive has fully extended.

Noise Levels

During normal operation the level of airborne noise emitted by the system will not exceed 70db measured at 1 meter.

Maintenance

Cleaning CloneSelect Imager

The machine should be cleaned each day.

Clean the machine using 80% ethanol or dilute detergent and a lint-free cloth. Organic solvents and abrasive cleaners should not be used, as they will damage the cover.

Do not pour cleaning solution directly on to the machine or other objects; apply using a suitable lint-free cloth.

Regular Maintenance

Daily

- Ensure that CloneSelect Imager is free from dirt and dust (see Cleaning CloneSelect Imager - above).

Weekly

- Check operation of equipment
- Check calibration
- Check for damage

Annually

- Maintenance by manufacturer

We would strongly recommend that maintenance be carried out regularly and by a Molecular Devices-approved service engineer.

Maintenance contracts can be obtained from Molecular Devices, see Contact Details on page 69.

General Precautions

- All waste must be disposed of according to local regulation.
- Do not use in explosive environments.

Technical Specifications

Manufactured by Molecular Devices (New Milton) Ltd.

Dimensions:

Total assembled:

Size: 575mm (width) x 720mm (depth) x 438mm
(height)

Excluding ancillary equipment

Weight: 45 kg

Operating Environment:

Indoor use only

Temperature: 10°C to 40°C

Humidity: 20 to 80% non-condensing

Altitude: Up to 2000M

Mains supply: +/- 10% Rated Voltage

Transient overvoltage: Installation Category (Overvoltage category) II

Rated pollution: Pollution degree 2

Electrical Supply:




Voltage: 100 - 240V AC 50-60 Hz single phase

Power: 100VA

Connections: IEC Input

Fuses: F1 - F3A

Symbols on Equipment

Symbol	Meaning
	Beware moving parts. Refer to user manual for operating instructions
	The year of manufacture is given on the serial number label which is located on the rear of the unit
	European Economic Area (EEA) This mark on the product indicates compliance with the following EEC Directives: <ul style="list-style-type: none"> - 98/37/EEC, relating to The Supply of Machinery (Safety) Regulations 1992 - 89/336/EEC with amendments 92/31/EEC, relating to Electromagnetic Compatibility - 73/23/EEC, the Low Voltage Directive

Appendix A

Compatible Plate Types

The standard CloneSelect Imager is configured to use the following plate types:

Manufacturer	Catalogue number	Well format
Molecular Devices	W1500, W1505, W1510, W1515, W1550, W1555, W1560, W1565	96
	W1300, W1305, W1310, W1315, W1350, W1355, W1360, W1365	24
	W1100, W1105, W1110, W1115, W1150, W1155, W1160, W1165	6
	W1000, W1005, W1007, W1009, W1050, W1055, W1057, W1059	1
Corning / Costar	3595, 3596	96
	3524	24
	3513	12
	3506	6
Greiner	655180	96
	662160	24
	657160	6
Nunc	140675	6

Note: A manufacturer often supplies several plates of the same well format with different catalogue numbers. In many cases these plates have the same dimensions, so if your plate type does not appear in the above table, contact Molecular Devices (New Milton) Ltd.

Appendix B

Well List File Format

This appendix describes the format of a **Well List** file to be imported into an **Image Microplate** process via the **Select Source** window.

The file containing a list of wells to image should be a text file (.txt) with the plate type as the first line. Each well to be imaged is then entered on a new line. Note that each file describes a single plate only.

PetriWell 96 well

A6

B5

E8

etc.

Appendix C

Setting up Cell Number Estimation

The confluence values measured by CloneSelect Imager can be used to non-invasively estimate the number of cells in each well. This is achieved by a cell number estimation formula. Because of the variation in size or properties of different cell types (e.g. adherent versus suspension cells) it is essential to generate a cell number estimation formula for each cell type.

Generation of the formula on CloneSelect Imager requires preparation of a dilution series of your cell type with known cell number. The following section provides guidelines for setting up a dilution series:

Setting up a dilution series

- Check viability of your cell line (e.g. using Trypan Blue). It is essential that viability is high (95-100%).
- Prepare a cell suspension in complete medium with the following density:
 - 1 000 000 (10⁶) cells/ml for suspension cells (e.g. hybridomas)
 - 250 000 cells/ml for adherent cells
- Create a dilution series:

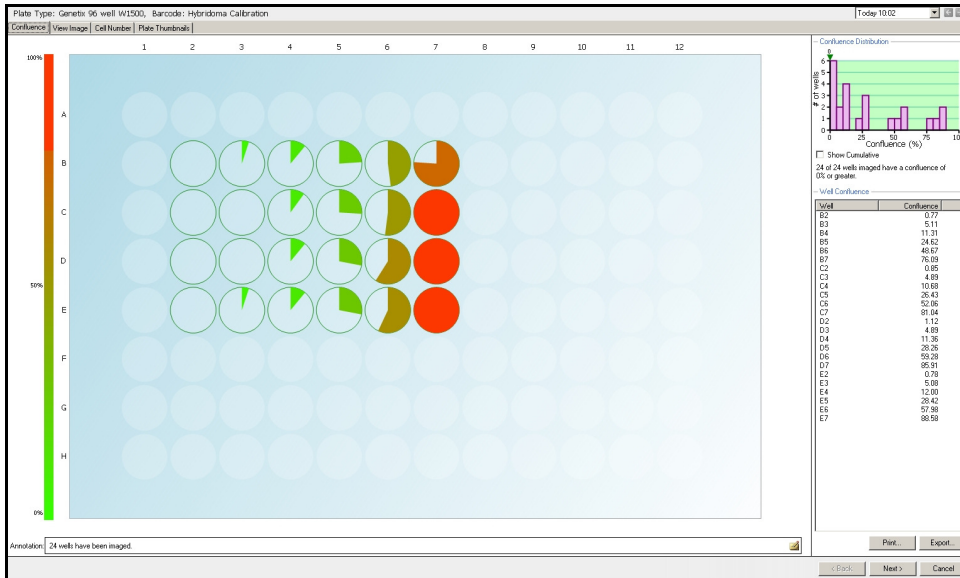
% dilution	Cell suspension (µl)	Medium (µl)	Number of cells in 200µl	
			Suspension cells	Adherent cells
100	2000	0	200 000	50 000
75	1500	500	150 000	37 500
50	1000	1000	100 000	25 000
25	500	1500	50 000	12 500
12.5	250	1750	25 000	6 250
6.25	125	1875	12 500	3 125
3.125	62.5	1938	6 250	1562

- Dispense 200µl aliquots of each dilution into a 96-well plate, plating out at least 4 replicates for each dilution. It is recommended to dispense the cell samples vertically to ensure an even spread of cells.
- Incubate the plate to allow the cells to settle. For suspension cells it is recommended to leave for 30 minutes, and for adherent cells leave at least 4 hours to attach.

Creating a cell number estimation formula

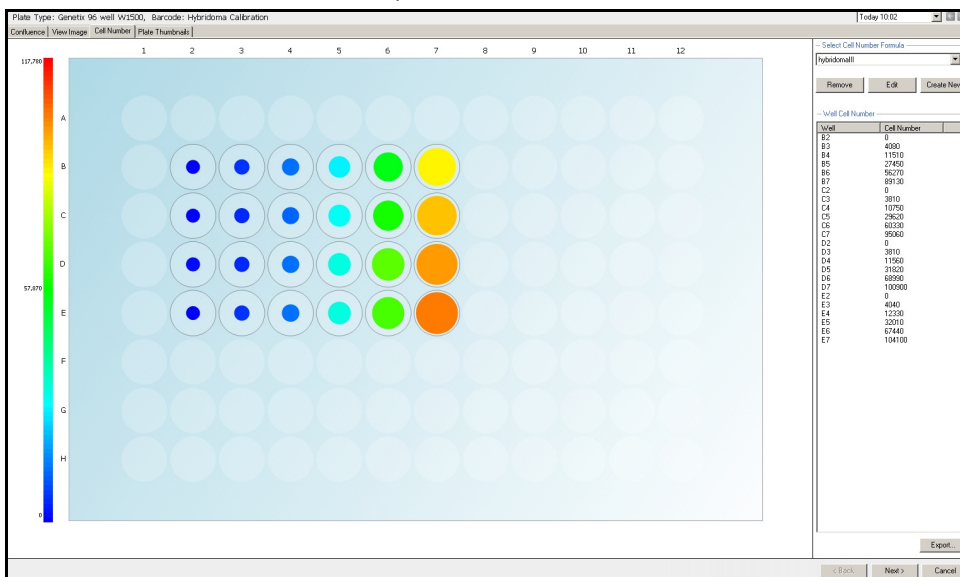
→ Image the dilution series on CloneSelect Imager.

→ In the Confluence tab, check that the dilution points with confluence between 5% and 80% have acceptable replicates.



Confluence tab with expected dilution point / confluence percentage

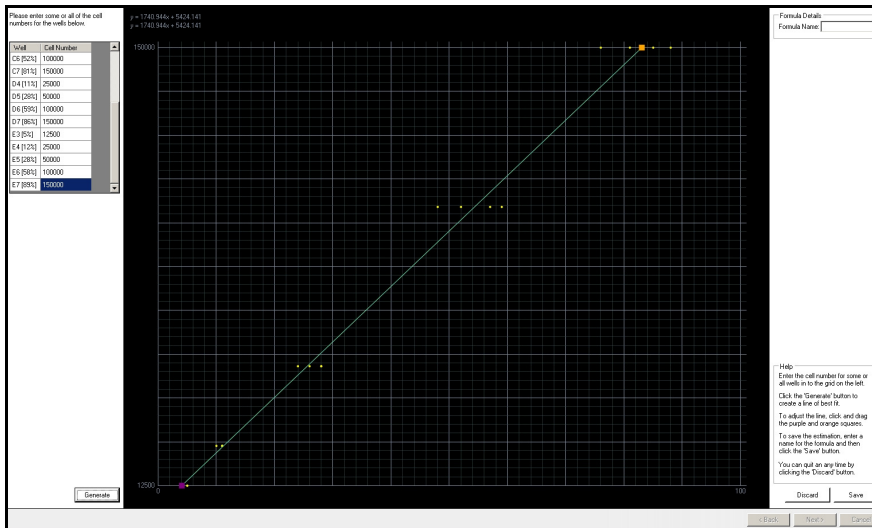
→ In the **Cell Number** tab, click on the **Create New** button.



Cell Number tab. Viewed with a previously defined cell number formula

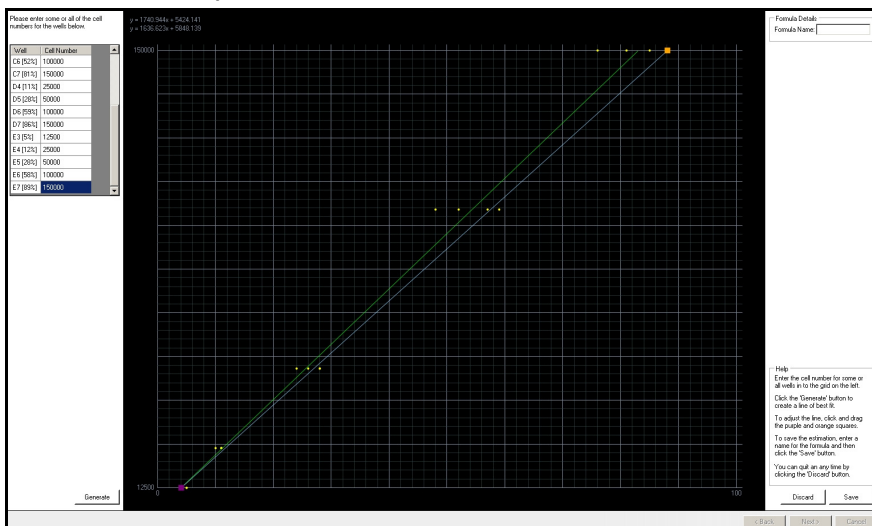
→ This opens the Cell Number Graphical Display window.

- For wells with confluence between 5% and 80%, input the cell number associated with the wells into the table in the left panel.
- Click on the **Generate** button to create a graphical plot of the data and the statistical equation shown in the top left corner. If data has been inputted incorrectly it can easily be re-entered and corrected using the **Generate** button.



Cell Number Graphical Display window showing plotted data.

- The line of best fit can be manually adjusted by dragging the handles at the end of the line. The formula is displayed at the top left and the adjusted version is displayed immediately below it. All cell estimate results are based on the adjusted version.



Two lines visible, showing the manually moved line in grey

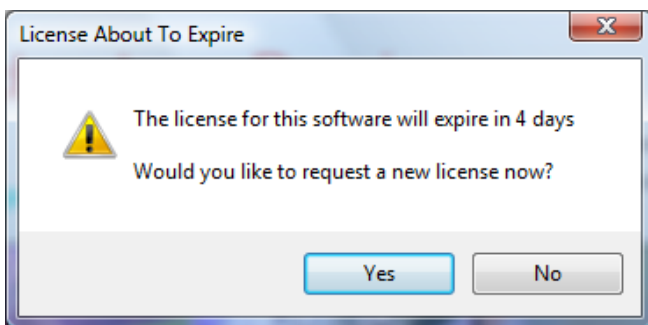
- Assign a name to the data set in the **Formula Name** box.
- Click **Save** to save the cell number estimation formula and return to the **Cell Number** tab. Click **Discard** to abandon the changes.

→ The new formula can then be used to estimate cell numbers in plates containing the same cell type.

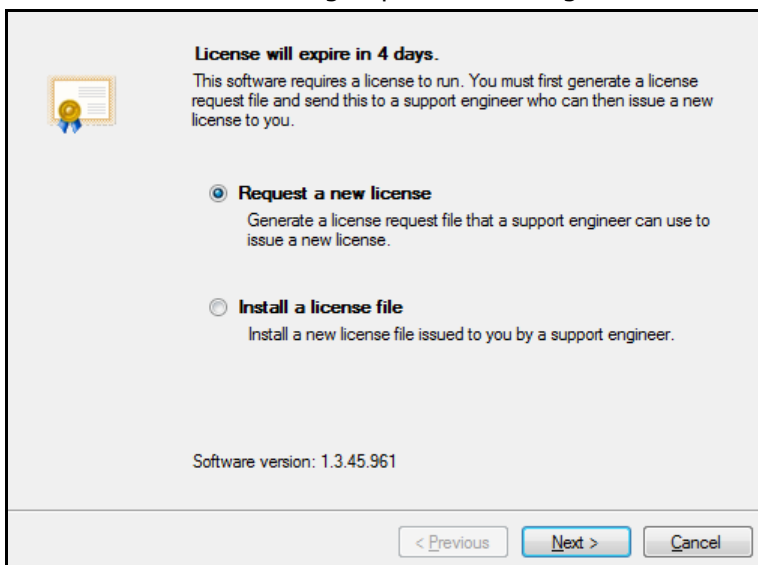
Appendix D

Requesting a License

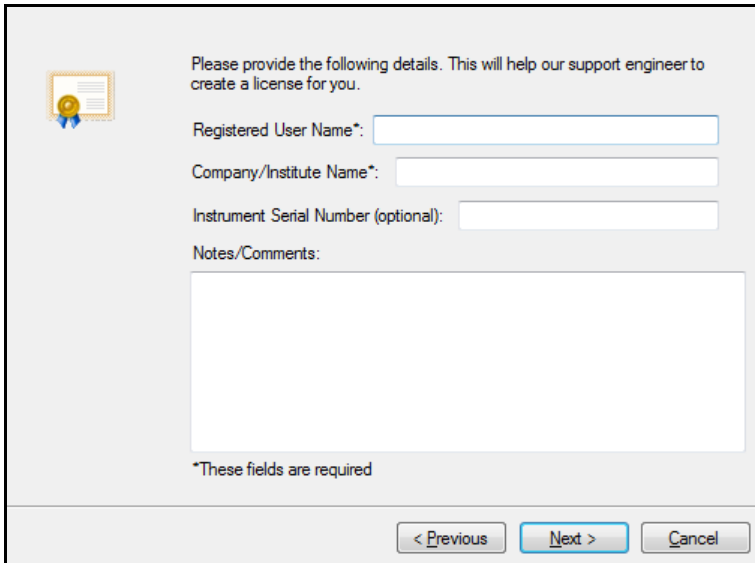
If the following message is displayed after starting the CloneSelect Imager software, a new license will need to be requested. Create a license request and then e-mail it to Customer Support.



→ Click **Yes**. This brings up the following window:



→ Select "Request a new license" and click Next.



Please provide the following details. This will help our support engineer to create a license for you.

Registered User Name*:

Company/Institute Name*:

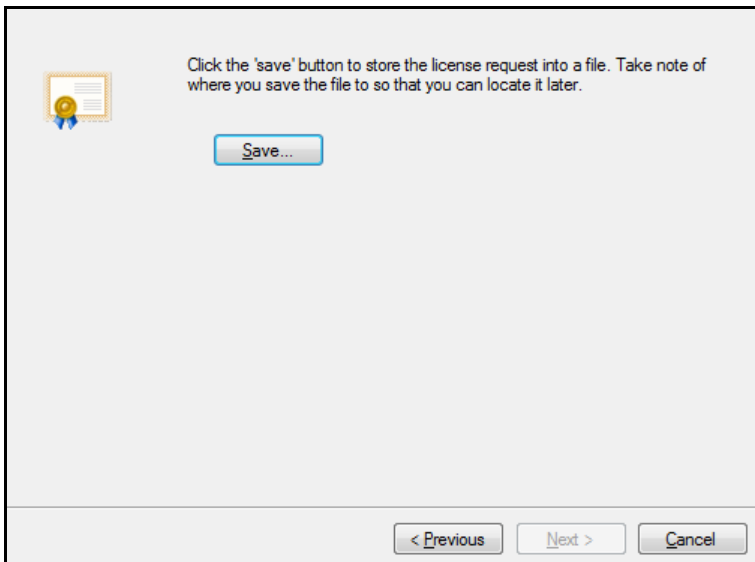
Instrument Serial Number (optional):

Notes/Comments:

*These fields are required

< Previous Next > Cancel

→ Enter the relevant details. Click **Next**.

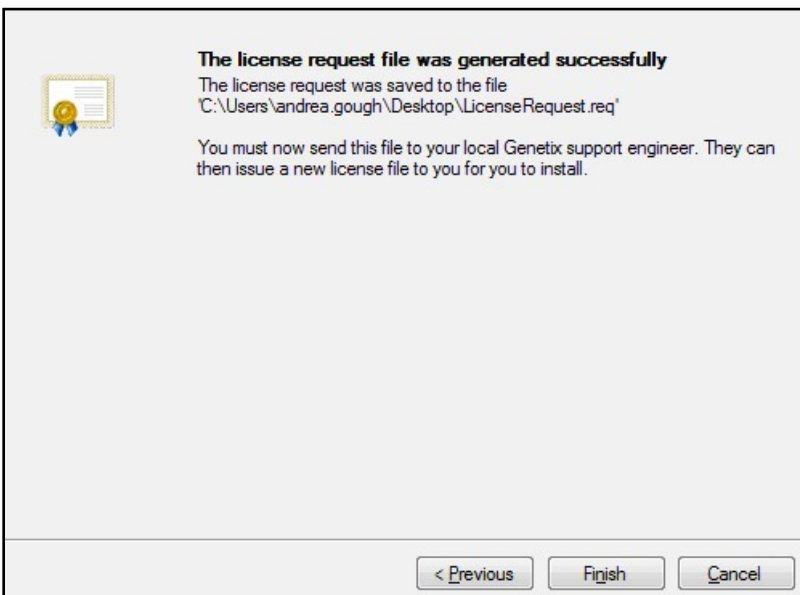
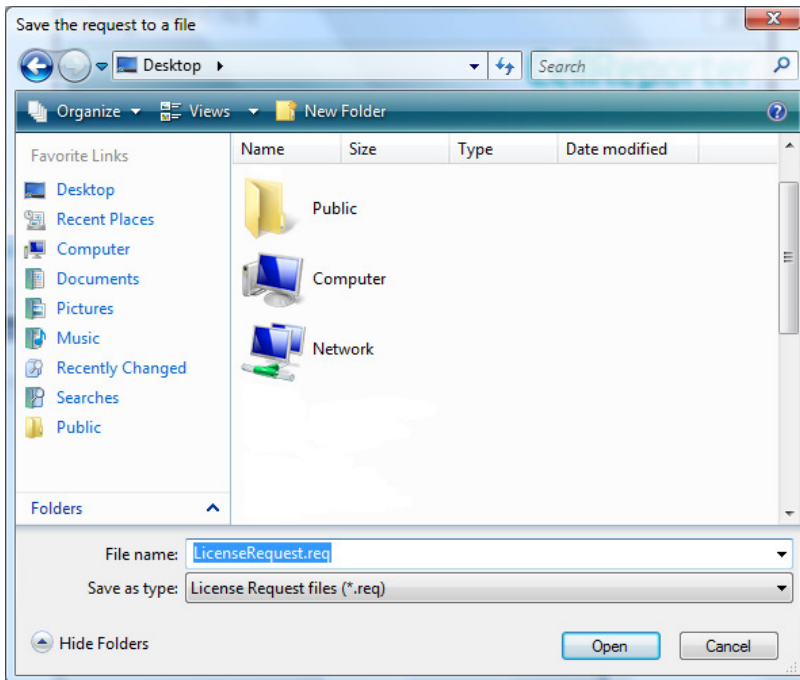


Click the 'save' button to store the license request into a file. Take note of where you save the file to so that you can locate it later.

Save...

< Previous Next > Cancel

→ Save file to a location where it can be attached to an e-mail.

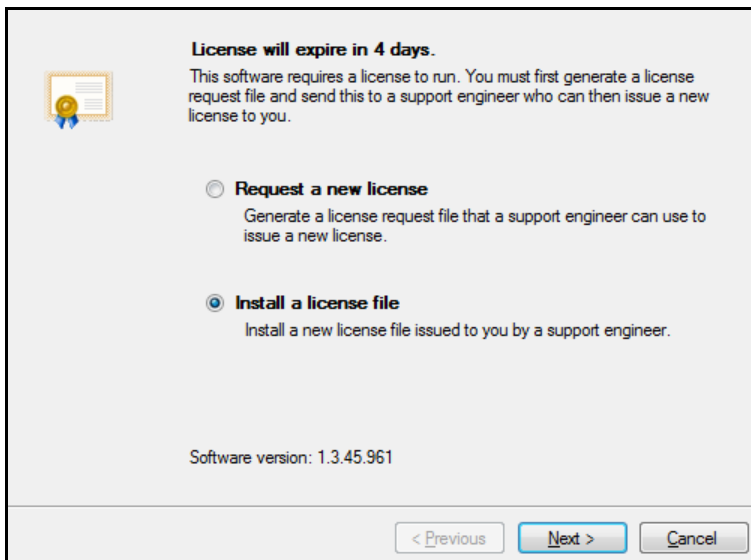


→ E-mail this **.req** file to a Molecular Devices representative.

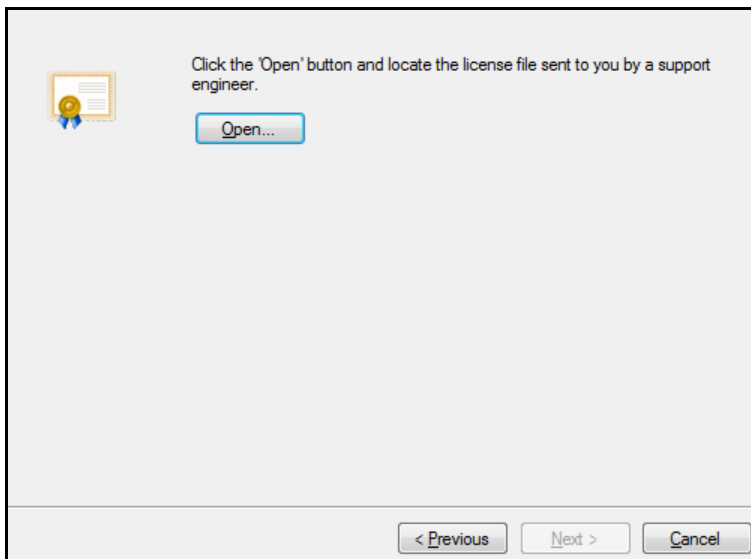
→ Click **Finish**.

Installing a License

- When a license file is received from a Customer Support representative, save the file to a place accessible on the computer of your instrument.
- Re-open the CloneSelect Imager software.
- Select "Install a license file":

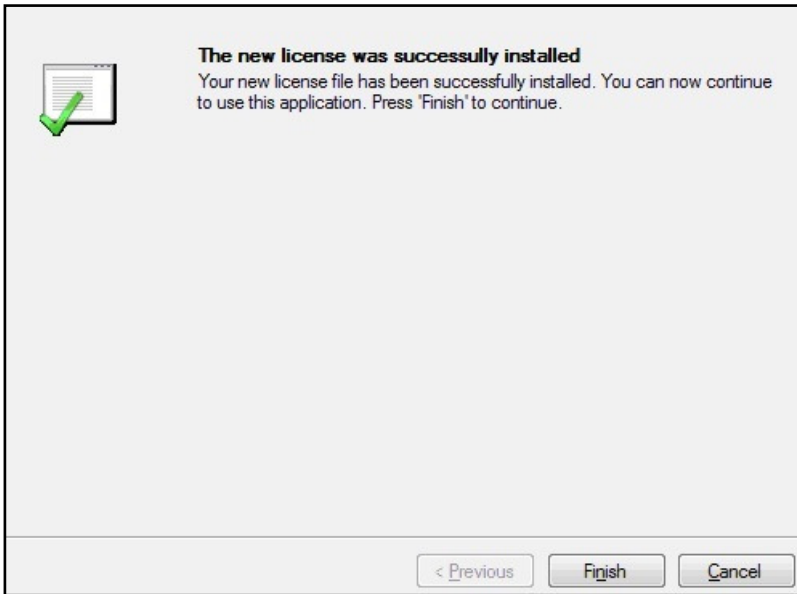


- Click **Next**.



- Click **Open**, and browse to the location where the license (*.License) file was saved.
- Open the file.

→ A window should appear to indicate that the license file was successfully installed:



→ Click **Finish**. The software will now start.

If for any reason the license does not appear to work as expected please contact Customer Support for assistance.

Index

A	
Alignment	9
Annotation	26
Archive	22
Auto focus before run	15
Averages	35
B	
Barcodes	15
C	
Calibration	9
Cancel	16
Cell Detection Method 1	15
Cell Detection Method 2	15
Cell Detection Method 3	15
Close Process'	20
colony forming assay	48
Contrast	27, 33
D	
Deselect All	16
Details	14
Display Statistic	31
Dual Gates	26
E	
Enhance Images	15
Export Image Sequence	32
F	
Focus	17
Full Scan	15
G	
Grab	18
Guide	14
H	
Histogram Bins	31
HTML Report	37
I	
Image Another Plate	20
Image Microplate	14
Import	16
L	
Leveling Feet	6
Load Results	21
M	
Main Navigation Screen	10
Mean Loci Area	30
Menu Options	11
Minimum Area	30
Minimum Compactness	30
N	
Next	16
P	
Partial Scan	15
Plate Carrier	6
Power-Up Procedure	9
R	
Rebooting	9
Re-image	19
Remove Plate	20
Rename	23
Return To Plate View	46
Review focus	15
Review Focus	16
S	
Save process as template...	20
Save process'	20
Scale Bar	18
Scan Options	16
Scan type	15
Select All	16
Show Cumulative	26
Source	15
Summary	14
W	
Well Data	30
Well Name	27
Workstation	8

Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest possible level of technical service.

Our support web site, www.moleculardevices.com/support.html, has a link to the Knowledge Base with technical notes, software upgrades, safety data sheets, and other resources. If you do not find the answers you are seeking, follow the links to the Technical Support Service Request Form to send an email message to a pool of technical support representatives.

You can contact your local representative or contact Molecular Devices Technical Support by telephone at 800-635-5577 (U.S. only) or +1 408-747-1700.

In Europe call +44 (0) 118 944 8000.

Please have your instrument serial number or Work Order number, and your software version number available when you call.

For research use only. Not for use in diagnostic procedures.

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/productpatents>

Product manufactured by Molecular Devices, LLC.
1311 Orleans Drive, Sunnyvale, California, United States of America 94089.

Molecular Devices, LLC is ISO 9001 registered.

© 2015 Molecular Devices, LLC.
All rights reserved.1