

ClonePix FL Change History

Changes and fixes made since version 1.2.13.971, up to and including build 1.2.15.1032:

1. Ensured during the Pick Run process that source well, source barcode or any statistic involved in the current groups cannot be deselected.
2. Fixed the issue where pressing the back button at the results stage of the Pick Run process caused the results to be saved in a subfolder.
3. Fixed the issue that caused picked colonies not to be auto logged (when the option is selected) in the Pick Run process.
4. Enabled the software to be installed on Windows 7 Service Pack 1.

Changes and fixes made since version 1.2.4.846, up to and including build 1.2.13.971:

1. Fixed an issue in the Imaging and Pick Run process start screens where the start button can be drawn on top of the process property summary.
2. Removed the option to export statistics of picked colonies into Excel XML from an imaging run.
3. Ensured that the illumination cover is closed before asking for destination plates to be loaded.
4. Fixed the issue that allowed a de-selected group to be attempted to be deleted.
5. Stopped repeating group information in both image settings and picking summary sections of the HTML report.
6. Fixed an anomaly in how a plate being left out on the bed is dealt with at the start of imaging.
7. Added a well index property to the graph screens.
8. Fixed the cancel button implementation in Prepare for Pick Run process.
9. Added a back button implementation to Prepare for Pick Run process.
10. Added a plate map for each destination plate in the picking summary section of the HTML reports.
11. Fixed the cancel button implementation in Replace Picking Head process.
12. Added a UV sanitise step to the Replace Picking Head process.
13. Added a warning when pressing the back button at the Select Groups stage of the Pick Run process.
14. Added Match Destination to Source as an option in the Pick Run process.
15. Corrected the definition of the Normalized Intensity property.
16. Fixed the issue where changes to certain Image Settings were not being saved during the Pick Run process.
17. Added measures to combat the apparent memory leak in high volumes of large area plates in the Pick Run process.
18. Removed unnecessary reconnections to the Anorad motion controller.
19. Corrected the description of the Feature ID property.
20. Replaced the message box asking the user if they want to continue using the loaded plate with a question within Fusion.
21. Tailored the load plate messages to the situation they are being displayed in.
22. Started sending back the current destination plate before asking for source plates to be loaded.
23. Altered the instructions on some screens in source well datuming.
24. Removed unnecessary line breaks from the Picking Summary in Pick Run process.
25. Removed the summary panel from the Rank Plot screen.
26. Fixed an issue on the Imaging and Pick Run process start screens where the incorrect prime configuration could be preselected.
27. Fixed an issue where running the same Utility Process twice would cause an error.
28. Fixed an issue where running the same Picking Head Management process twice would cause an error.
29. Standardised the default dry time throughout the software.
30. Ensured that "Use each colony size when calculating exterior statistics" is off by default in Imaging and Pick Run processes.
31. Ensured that the pick results are saved even when the Pick Run is cancelled during the picking stage.
32. Fixed an issue in the HTML report where the incorrect detection parameters could be displayed.
33. Made process templates read only on saving them and added measures to prevent them being overwritten.
34. Added colony information for the "All Features" node in the Groups tab of Preview and Results.

Changes and fixes made since version 1.2.3.842, up to and including build 1.2.4.846:

1. Fixed an issue that prevented 'organise by well' from correctly organising the destination plate by source well.
2. Fixed an issue that prevented 'order by' during picking from working when a statistic that has units was specified.
3. Prevented an issue that can occur during results analysis from terminating the application.
4. Added drive tuning parameters.
5. Added new branding to the splash screen.

Changes and fixes made since version 1.0.1.791, up to and including build 1.2.3.842:

1. Made the 3D image snap to the last selected feature when the mouse cursor leaves the image.
2. Removed the Hardware model view from general display.
3. Changed the Rank Plot histogram to a stacked histogram, with group colours so that group distribution can be seen.
4. Fixed an issue where the plate row and column pitches were not used correctly.
5. Fixed an issue where dependant devices cannot connect if they are created in a specific order.
6. Fixed an issue where the incorrect emission filter was used when only some of the filters were licensed.
7. Improved the handling of plates in stackers.
8. Fixed an issue where the 'Reprocess' button did not reprocess all images, even when the user asked to do so.
9. Fixed issue where defaults verb would not store default process values.
10. Fixed an issue where the picking process could hang after the pause button was pressed.
11. Removed unnecessary icons from the process selection screen.
12. Fixed incorrect destination plate name displayed in the summary screen during a pick process.
13. Destination plate templates added. Allows selected destination wells to be left empty.
14. Added destination wells display to Output Results screen and reports.
15. Retired first deposit well option (superseded by destination plate templates).
16. Added ability to load images from broken optical configuration names that were truncated at full-stops in version 1.0.1.753 or earlier.
17. Added check to ensure destination plate definition is valid at beginning of picking run
18. Altered Output Results screen layout.
19. Colour coded the deposit wells on the results screen and enhanced the user interaction with that control.
20. Corrected issues with versioning in *Imaging*, *Prepare For Picking* and *Picking* saved processes.
21. Altered question and error text when getting a plate.
22. Added process property validation for Optical Configurations.
23. Improved the way administrate properties deals with 'hidden' process properties.
24. Fixed back button problems in Picking and Imaging processes.
25. Added units to statistic caption in scatter plot and histogram selection lists.
26. Added ability to make external area diameter specific to each individual colony diameter, rather than average colony diameter.
27. Added ability to allow the user to set the size of the external statistics diameter.
28. Added the new external statistics parameters to the imaging summary and HTML report.
29. Adjusted the order of the groups listed in the summary screen and HTML report to reflect correct priority.
30. Adjusted scatter plot and rank plot tool tips to include feature ID and units for axis statistics.
31. Adjusted review results process to reduce time taken to collate results.
32. Switched to .NET Framework 3.5 SP1.
33. Fixed error when deleting groups, expressions and rules.
34. Made scatter and rank plot not reset zoom when changing groups.
35. Ensured that the illumination cover is closed during UV sterilisation.
36. Modified the HTML report generation so that the report title defaults to the 'report name', which itself has been renamed to 'folder name' to more accurately reflect its purpose.
37. Display the exterior statistics diameter in mm for the mean colony diameter on the detection page, to enable the user to better understand the affect of changing the multiplier value.
38. Fixed issue when reprocessing previous results stored at a read-only location.
39. Modified dispersal to always disperse all eight pins, irrespective of how many pins were picked, to minimise time taken.
40. Retained 3D view position when switching between fluorophores.
41. Enabled 'Goto pick results' menu item in statistics table when colony has been picked.
42. Changed name of prepare for picking process to "Prepare for Pick Run".

Changes and fixes made since version 1.0.1.753, up to and including build 1.0.1.791:

1. Hourglass/Wait Display/Progress bars added at processing times, robotic movements etc.
2. HTML reports, the reporting facility has encompassed a number of requests to output information, including:
 - Custom Fields
 - Image Property Settings
 - Pick Property Settings
 - Colony Statistics
 - Scatter Plot
 - Rank Plot & Frequency Histogram
 - Pick Summary
3. Back button on Imaging & Picking processes.
4. Automated logging of statistics during a pick run. The feature is activated from the PickVerb properties and has the same functionality as the data export wizard.
5. The Output Results tab has a new layout that displays the statistic the colonies were 'Ordered By'.
6. The pre-run summary has been enhanced to include all the imaging/picking run property settings.
7. There is now a barcode option to automatically assign a barcode if a barcode fails to be read.
8. Extra column in Review Results to signify if the data was from an imaging or picking run. Also the operator column is now working and the name of the machine the run was performed has been added as an extra column.
9. Right clicking a colony now has an option to 'Go to Results Output'.
10. There is now a right click option on the Image, Scatter Plot and Rank Plot that allows the user to copy to the clipboard and then paste the bitmap into Paint, Word etc.
11. The focus bar has indicators to signify which way is up or down.
12. On the Groups tab the number of colonies found is displayed alongside each Group name.
13. The Group property histogram slide adjusters now inherit the same colour as its Group.
14. Excel xml export is activated in export statistics. This option will export each source plates colonies to a separate Excel worksheet.
15. The real time picking screen has a modification to the way the pick information is displayed. It is in a grid format and 0000 is no longer used to align the colony counts.
16. A request for confirmation is now given to the operator asking to confirm if they wish to cancel when cancelling on swapping cassettes in batch processing.
17. Select & Export statistic buttons are now located on the statistics tab.
18. The Import button on the well plate selector is now disabled.
19. The well plate image selector now only displays one tracked image, previously it was displaying four tracked images depending on the mouse position.
20. Wash defaults changed to 3 purge cycles, 3 wash cycles and 10 seconds in the dryer.
21. There is now an edge excluded group that rejects colonies close to the edge of the plate. You will notice the well mask is drawn at the edge of the well. This was added to counteract a problem that was found in detecting proximity of neighbouring colonies.
22. Override purge cycles now defaults to 30.
23. Barcode is now referred to as Source Barcode and Destination Barcode.
24. There was an issue when configuration names included a period(.).
25. Select stats window has a select/deselect all option.
26. Pick all colonies option has been reorganised with an option to limit the number of colonies.
27. 'Override purge cycles' renamed to 'Change wash bottle', head moves clear before the dialog box is presented, and the dialog has a cancel option
28. Starting a run without pressing the robot reset button gives a better error message
29. Starting the run without air gives an error message when opening/closing the illumination cover
30. Well alignment screen displays rectangle images now, not squares
31. Report wizard button moved to data table, next to data export button

32. Image ruler buttons added to image panel
33. Tool tabs simplified (group boxes etc. removed)
34. Added Composite image generation
35. Display equation of line of best fit on the scatter plot
36. Indicate group membership on pop-up menus for selected colonies.
37. Optimised many aspects of the user interface to make it more responsive.