

CloneSelect Imager

CHANGE HISTORY



Contents

Changes	3
Build 1.3.73.1073.....	3
Build 1.3.56.1043.....	3
Build 1.3.51.990.....	3
Build 1.3.45.961.....	3
Build 1.2.4.871.....	4
Build 1.2.2.848.....	4
Build 1.1.0.818.....	4
Build 1.0.1.807.....	4
Build 1.0.1.795.....	5
Contact Details	6

Changes

Build 1.3.73.1073

Changes and fixes made since version 1.3.56.1043 up to and including build 1.3.73.1073:

1. Added new licensable feature: Cell Migration to enable interpretation of cell migration.
2. A new pipeline named 'Very Low Contrast' has been added to improve confluence calculation when using very low contrast cell lines.
3. Fixed bug. If the results folder for a plate is deleted or moved a meaningful message is now displayed instead of an exception alert.

Build 1.3.56.1043

Changes and fixes made since version 1.3.51.990, up to and including build 1.3.56.1043:

1. Fixed a number of issues with the export of data.
2. Option to support multiple results locations has been added.
3. Two extra functions have been added to the automation interface.

Build 1.3.51.990

Changes and fixes made since version 1.3.45.961, up to and including build 1.3.51.990:

1. Fixed Batch Plate Imaging process so that user selected cell detection methods are used correctly.
2. Fixed data export of loci information to the Excel xml file format.
3. Fixed system hang during imaging when enhance images is disabled and images are being stored in collections.
4. Fixed crash in whole well image display when contrast is altered.
5. Moving between time points on the loci tab now does not cause the zoom to reset.
6. The pick list on the loci tab is now not reset if the user moves to a different tab.
7. The pick list on the loci tab now tracks the user selected item.

Build 1.3.45.961

Changes and fixes made since version 1.2.4.871, up to and including build 1.3.45.961:

1. A new cell detection method has been added to improve detection of low contrast cell lines.
2. A new cell detection method has been added to improve detection of colonies.
3. Barcode or plate name can be changed in results review.
4. Time sequence images in bmp, png and jpg formats are exportable from the loci tab of image review.
5. Area histogram on the loci tab now uses μm^2 instead of mm^2 .
6. Area and compactness histograms on the loci tab allow gating of large as well as small features.
7. Number of histogram bins can be set when exporting loci data.

8. Users can now give a descriptive name to a well on the confluence, cell number and loci tabs.
9. Wells can be excluded from loci processing on the thumbnail view tab.
10. Images can be stored in a collection on disk instead of individual files. This decreases the time taken to move, backup and delete results.

Build 1.2.4.871

Changes and fixes made since version 1.2.2.848, up to and including build 1.2.4.871:

1. Fixed a problem when loci count analysis occurred outside of mask.
2. Added export of compactness and area data for every found feature to csv, xml and excel xml files.

Build 1.2.2.848

Changes and fixes made since version 1.1.0.818, up to and including build 1.2.2.848:

1. Changed automated imaging process to support image compression
2. Added 'Are you sure' prompt when pausing an ICCMS process
3. Enforced requirement for images to be saved to local disk
4. Added error report generation facility

Build 1.1.0.818

Changes and fixes made since version 1.0.1.807, up to and including build 1.1.0.818:

1. Resolved memory leak when loading image files
2. Added more decimal places for loci area
3. Minor performance improvements

Build 1.0.1.807

Changes and fixes made since version 1.0.1.795, up to and including build 1.0.1.807:

1. Is now compatible with the new OEM integration kit (WCF-based)
2. Fixes an issue where the plate row and column pitches were not used correctly
3. Ensures that the magnification caption is visible on the image display when plates are being imaged
4. Removes the black line drawn around the red mask area to make it less intrusive
5. Removes the Hardware model view from general display
6. Renames the ColdFire upload/download operations to make it more obvious what is happening
7. Fixes an issue where manual barcoding was not working correctly
8. Fixes a problem where dependant devices cannot connect if they are created in a specific order

Build 1.0.1.795

Changes and fixes made since version 0.4.0.428, up to and including 1.0.1.795:

1. Full well view for all microplate formats
2. Improved auto focus with feedback
3. Improved growth curves display with selectable wells, well thumbnails and fitted curves
4. Summary page showing when the plate was imaged and growth trends
5. Faster loci counting with improved criteria display using histograms
6. Added 'back' button during plate focusing etc.
7. Store name of PC images are taken on to facilitate tracking data from multiple instruments
8. Selectable confluence generation algorithms
9. Added ability to remove cell estimation formulae
10. Ability to toggle confluence overlay display in focus screen
11. Added ability to export graph images
12. Added HTML plate report
13. Export clone size classes when loci counting
14. Added support for Genetix plate loader
15. Numerous usability and performance improvements

Contact Details

Molecular Devices

Queensway, New Milton
Hampshire BH25 5NN, UK

Tel: +44 (0) 1425 624 600
Fax: +44 (0) 1425 624 700
Web: www.moleculardevices.com/genetix

For all technical queries please contact your nearest Customer Support group. Visit www.moleculardevices.com/genetix for latest contact details.

Trademarks

ClonePix™, CloneSelect™, CellReporter™, HalfBD®, 'WRich®, SlidePath™, Data Arena™ and Image Arena™ are trademarks of Molecular Devices (New Milton) Ltd.

Copyright © 2011 by Molecular Devices (New Milton) Ltd

All rights reserved. No part of this publication may be reproduced, stored in a retrieval system, or transmitted, in any form by any means, electronic, mechanical, by photocopying, recording, or otherwise, without the prior written permission of Molecular Devices (New Milton) Ltd.

Information furnished by Molecular Devices (New Milton) Ltd is believed to be accurate and reliable; however, no responsibility is assumed by Molecular Devices (New Milton) Ltd, for its use; nor for any infringements of patents or other rights of third parties which may result from its use. No license is granted by implication or otherwise under any patent rights of Molecular Devices (New Milton) Ltd.

Revised June 2011