

ClonePix 2

CHANGE HISTORY



Contents

- Changes **3**
 - Build 1.2.89.1155 **3**
 - Build 1.2.81.1091 **3**
 - Build 1.2.80.1071 **3**
- Contact Details **5**

Changes

Build 1.2.89.1155

Changes and fixes made since version 1.2.81.1091, up to and including build 1.2.89.1155:

1. Fixed a problem where older results no longer reprocessed correctly.
2. Implemented the whole plate view for images.
3. Increased the speed of navigating results that are located across a network.

Build 1.2.81.1091

Changes and fixes made since version 1.2.80.1071, up to and including build 1.2.81.1091:

1. Replaced Image Resolution with Manual Image Resolution.
2. Started removing leading and trailing spaces from optical configuration names on save.
3. Altered the dispersal function in the Pick Run process to aspirate odd, dispense odd, aspirate even, dispense even.

Build 1.2.80.1071

Changes and fixes made since version 1.2.67.1030, up to and including build 1.2.80.1071:

1. Removed Leica Microsystems branding.
2. Fixed the issue where trans illumination could be erroneously used instead of epi illumination in commissioning processes.
3. Extended the Image Rotation process to take four images and removed the image overlaps.
4. Added a data validation step to the Image Resolution process.
5. Altered the Image Resolution process so that it involves a plate rather than a pin.
6. Removed an extra pin fire from the Drive Alignment process.
7. Fixed an issue in Coldfire Interpreter which caused booleans not to be processed correctly.
8. Enabled the software to be installed on Windows 7 Service Pack 1.
9. Fixed the issue that caused picked colonies not to be auto logged (when the option is selected) in the Pick Run process.
10. Added Molecular Devices branding.
11. Fixed the issue with the derived minimum and maximum colony diameter calculations and saved processes.

12. Addressed the issue that can occur with Drive Alignment and Drive Resolution processes when Genetix plates are not present in the consumables list.
13. Resolved the issue in Image Resolution that can cause a crash when less than two points are checked in the data visualisation stage.

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*®, *MRich*®, SlidePath™, Data Arena™ and Image Arena™ are trademarks of Molecular Devices (New Milton) Ltd.

Copyright © 2012 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 March 2012