

CloneSelect Imager

SOFTWARE UPGRADE GUIDE



Contents

| | |
|--|----------|
| Overview | 3 |
| Upgrading | 3 |
| Removing Previous Version | 3 |
| Installing New Version..... | 4 |
| Contact Details | 6 |

Overview

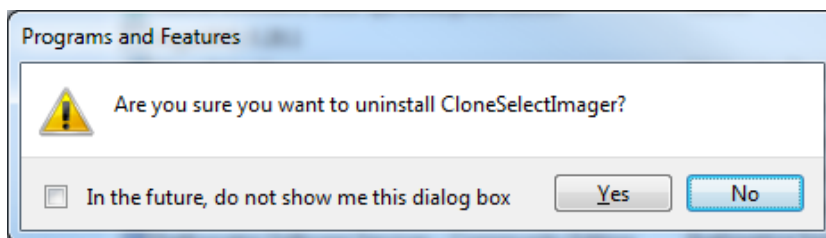
This document describes how to upgrade the CloneSelect Imager software from version 1.2.4.871 or higher. Please note that instruments that are licensed to use the OEM Automation kit and are integrated into a robot workcell must not be upgraded. The new version of software is not compatible with the old version, and the links between the third party software and CloneSelect Imager would have to be rewritten.

NOTE: Take a backup of your configuration prior to this upgrade.

Upgrading

Removing Previous Version

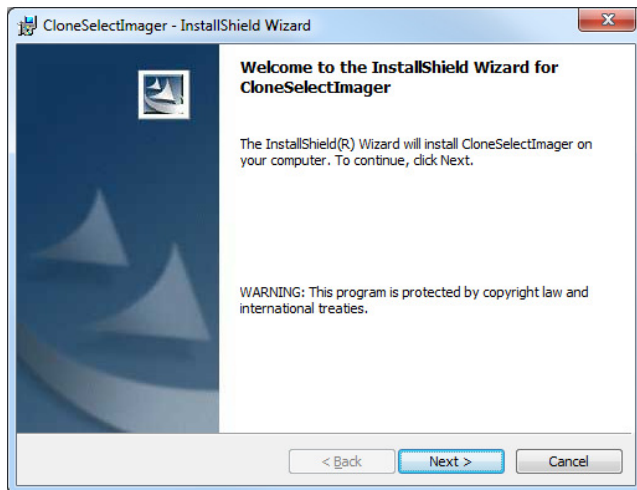
1. In Windows, open the Control Panel.
2. Double click the 'Add or Remove Programs' icon.
3. Right click the 'CloneSelectImager' item on the list and select the 'Uninstall' menu item.
4. When asked if you are sure that you want to remove 'CloneSelectImager', click the 'Yes' button.



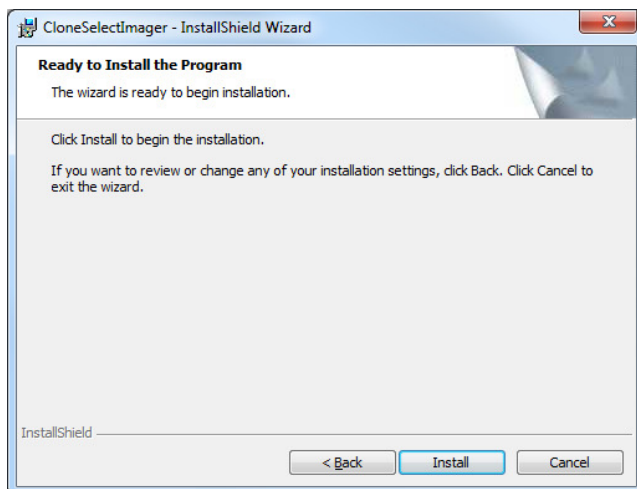
Installing New Version

This step is required when upgrading from any previous version.

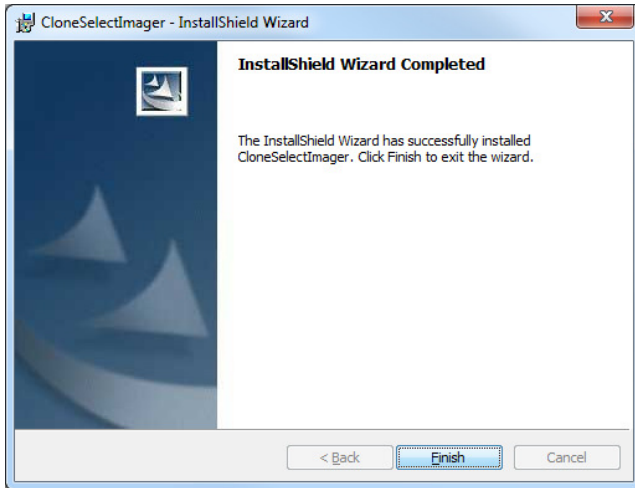
1. Insert the recovery DVD and browse to "DVD:\Setup\CloneSelect Imager x.x.xx.xxx" (Please Note: 'x.x.xx.xxx' will be the software version number.).
2. Run the 'Setup CloneSelectImager.exe' file found in that folder.
3. On the welcome screen, click the 'Next' button.



4. On the next screen, click the 'Install' button.



5. On the completion screen, click the 'Finish' button.



6. The installation is complete.

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