



CellXpress.ai

Automated Cell Culture System

Software Version 1.4.1

Release Notes



This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

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.

Copyright notice for AutoQuant 2D RealTime Deconvolution: © Media Cybernetics, Inc.

Patents: www.moleculardevices.com/patents

Product manufactured by Molecular Devices, LLC., a part of Danaher Corporation.
3860 N. First Street, San Jose, California, 95134, United States of America.
Molecular Devices, LLC is ISO 9001 registered.
©2025 Molecular Devices, LLC.
All rights reserved.



Contents

Chapter 1: CellXpress.ai Automated Cell Culture System	4
About This Guide	4
Product Documentation	4
Obtaining Support	5
Chapter 2: Issues Addressed	6
Droplet Formation Reduced After Dispensing to Liquid Waste (MON-3717)	6
Email Server Authentication No Longer Prevents System Notifications (MON-3707)	6
System Successfully Acquires a Large Number of Images (MON-3833)	6
Software Prevents Unsupported Image Tiling Settings (MON-3903)	6

Chapter 1: CellXpress.ai Automated Cell Culture System

1

The CellXpress.ai® Automated Cell Culture System is an AI-driven cell culture innovation hub that gives your team total control and oversight on demanding cell culture feeding and passaging schedules. The CellXpress.ai system minimizes time in the lab while maintaining a 24/7 schedule for growing and scaling multiple stem cell lines, spheroids, or organoids.

About This Guide

This guide is intended for the scientist using the CellXpress.ai system. It describes the issues addressed in version 1.4.1 of the CellXpress.ai Automated Cell Culture System Software.

The information in this guide is subject to change without notice. We recommend that you review the guide on the Molecular Devices Knowledge Base at support.moleculardevices.com for the most up-to-date information.

Product Documentation

The following guides are available on the Molecular Devices Knowledge Base at support.moleculardevices.com:

- *CellXpress.ai Pre-Installation Guide*
- *CellXpress.ai Safety Guide*
- *CellXpress.ai System User Guide*
- *CellXpress.ai Back Port Table User Guide*
- *CellXpress.ai Tumor Spheroid Application Guide*
- *CellXpress.ai Release Notes*



A hard copy of the *CellXpress.ai Safety Guide* is also included in the package with a new system.

In addition, the CellXpress.ai software includes context-sensitive Help that you can access from within the software. Just click the  **Help** icon to view Help for the current page.

 **Tip:** We recommend that you review the documentation before using the CellXpress.ai system.

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 level of technical service.

Our Support website—www.moleculardevices.com/service-support—describes the support options offered by Molecular Devices, including service plans and professional services. It also has a link to the Molecular Devices Knowledge Base, which contains documentation, technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance, you can submit a request to Molecular Devices Technical Support.

Technical Support

To contact Molecular Devices Technical Support, submit a support request through the Molecular Devices Knowledge Base at support.moleculardevices.com.

You can also submit a support request by phone. For regional support contact information, go to www.moleculardevices.com/contact.

Documentation

Review the product documentation on the Molecular Devices Knowledge Base at support.moleculardevices.com. In addition, online Help is available within the CellXpress.ai software.

Training

Molecular Devices provides training on the general operation of the CellXpress.ai system at the time of installation. Contact Molecular Devices Technical Support for details on training after installation.

This section describes the issues addressed in CellXpress.ai software version 1.4.1.

Droplet Formation Reduced After Dispensing to Liquid Waste (MON-3717)

Improvements to liquid class definitions reduce droplet formation below the tips after dispensing to liquid waste. This lowers the chance of falling droplets, which can contaminate the liquid handler worktable.

Email Server Authentication No Longer Prevents System Notifications (MON-3707)

Email notifications from the CellXpress.ai system are now delivered as expected when using an email server that requires username / password authentication.

System Successfully Acquires a Large Number of Images (MON-3833)

With sufficient storage space, the system can now acquire many images per plate. For example, an acquisition that includes many sites, wells, channels, and Z-stacks can produce up to 30,000 images, and the experiment runs as expected.

Software Prevents Unsupported Image Tiling Settings (MON-3903)

The software now prevents image imaging tiling settings that will result in tiled images that exceed the maximum allowed size.



Visit us on Spectranet.

Contact Us

Phone: +1-800-635-5577
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.