

MultiClamp™ 700B

Commander Software

Software Release Notes

MultiClamp 700B Commander Software 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.

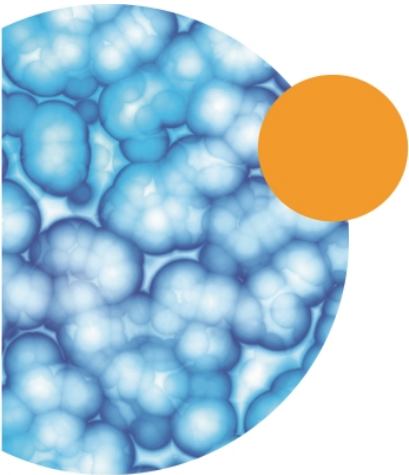
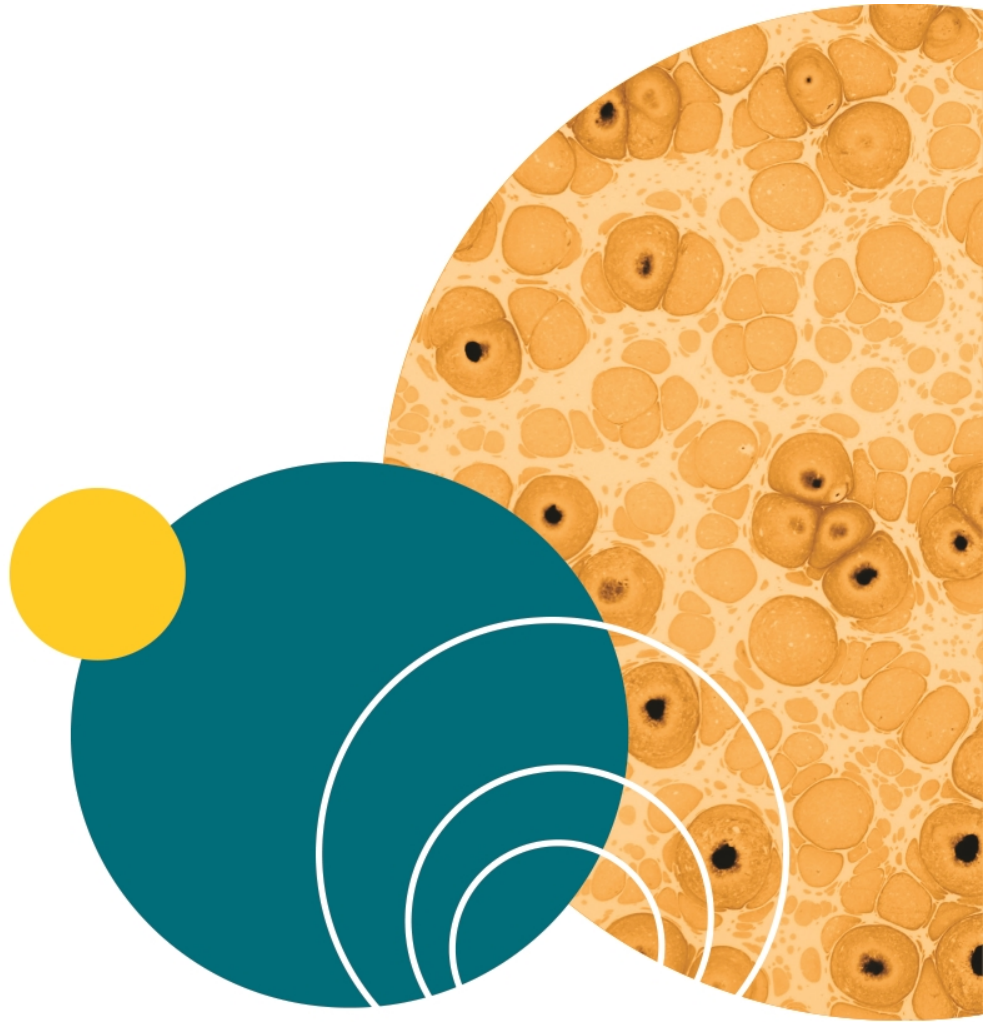
The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: <http://www.moleculardevices.com/productpatents>

Product manufactured by Molecular Devices, LLC.
1311 Orleans Drive, Sunnyvale, California, United States of America 94089.
Molecular Devices, LLC is ISO 9001 registered.
©2016 Molecular Devices, LLC.
All rights reserved.

Contents

- Chapter 1: MultiClamp Commander Software General Information 5**
 - Computer System Requirements 5
 - Software Installation 6
 - Obtaining Support 6
- Chapter 2: MultiClamp Commander Software Version 2.2: Software Release Notes .. 7**
 - New in MultiClamp Commander Software Version 2.2 7



Chapter 1: MultiClamp Commander Software General Information

MultiClamp Commander Software is the amplifier interface. In addition to the values that are generally telegraphed by hardware connections on conventional amplifiers (cell capacitance, filter cutoff frequency, and output gain), MultiClamp Commander Software provides five additional signal settings: command sensitivity, operating mode (voltage/current clamp), scaled output signal, scale factors, and scaling units of the output signal. These additional settings allow the data acquisition software, such as pCLAMP™ Software Version 10, to automatically configure stimulus and recording signals based on the commander software settings. The MultiClamp Commander Software interface runs independently of data acquisition software and therefore can be used with most data acquisition systems in stand-alone mode.

The following topics are included in this chapter:

- [Computer System Requirements on page 5](#)
- [Software Installation on page 6](#)
- [Obtaining Support on page 6](#)

Computer System Requirements

MultiClamp Commander Software version 2.2 can be installed on a computer with the following minimum system specifications:

Operating System

- Windows 10, 64-bit
- Windows 7, 64-bit

Hardware

- 1.5 GHz or higher processor
- 2 GB RAM or more
- 1 USB 2.0 port

Software Installation

The following procedure explains how to install the MultiClamp Commander Software version 2.2.

1. Double-click the **MultiClamp_2_2_x.exe** MultiClamp 700B Commander Software installation file. A Welcome to the MultiClamp 700B Commander Software Install Wizard dialog is displayed.
2. Click **Next**.
3. In the **License Agreement** dialog, select **I accept the terms of the license agreement**, and click **Next**.
4. In the **Destination Folder** dialog, the **Install MultiClamp 700B Commander Software to** field displays the default installation directory. To change the installation directory, click **Change**, navigate to the desired directory, then click **OK**.
5. Click **Next**.
6. In the **Select Program Folder** dialog, leave the displayed default **Program Folder** settings, and click **Next**.
7. Click **Next**.
8. When the installation is complete, click **Finish** to exit the wizard.

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 possible level of technical service.

Our support web site, www.moleculardevices.com/support, has a link to the Knowledge Base with technical notes, software upgrades, safety data sheets, and other resources. If you do not find the answers you are seeking, follow the links to the Technical Support Service Request Form to send an email message to a pool of technical support representatives.

You can contact your local representative or contact Molecular Devices Technical Support by telephone at 800-635-5577 (North America only) or +1 408-747-1700. In Europe call +44 (0) 118 944 8000.

To find regional support contact information, visit www.moleculardevices.com/contact.

Please have your instrument serial number or Work Order number and your software version number available when you call.

Chapter 2: MultiClamp Commander Software Version 2.2: Software Release Notes

The MultiClamp Commander Software version 2.2 update is a minor release. The following is a summary of the changes incorporated in this revision as compared to version 2.1.0.16, the last general release of the MultiClamp Commander Software.

- [New in MultiClamp Commander Software Version 2.2 on page 7](#)

New in MultiClamp Commander Software Version 2.2

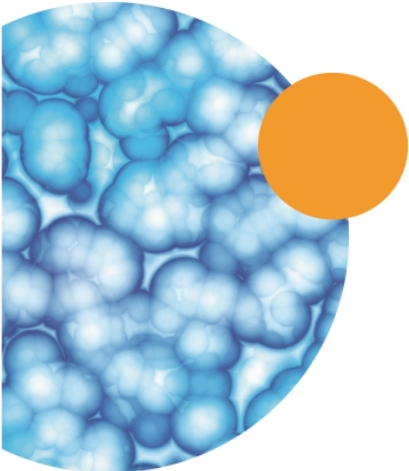
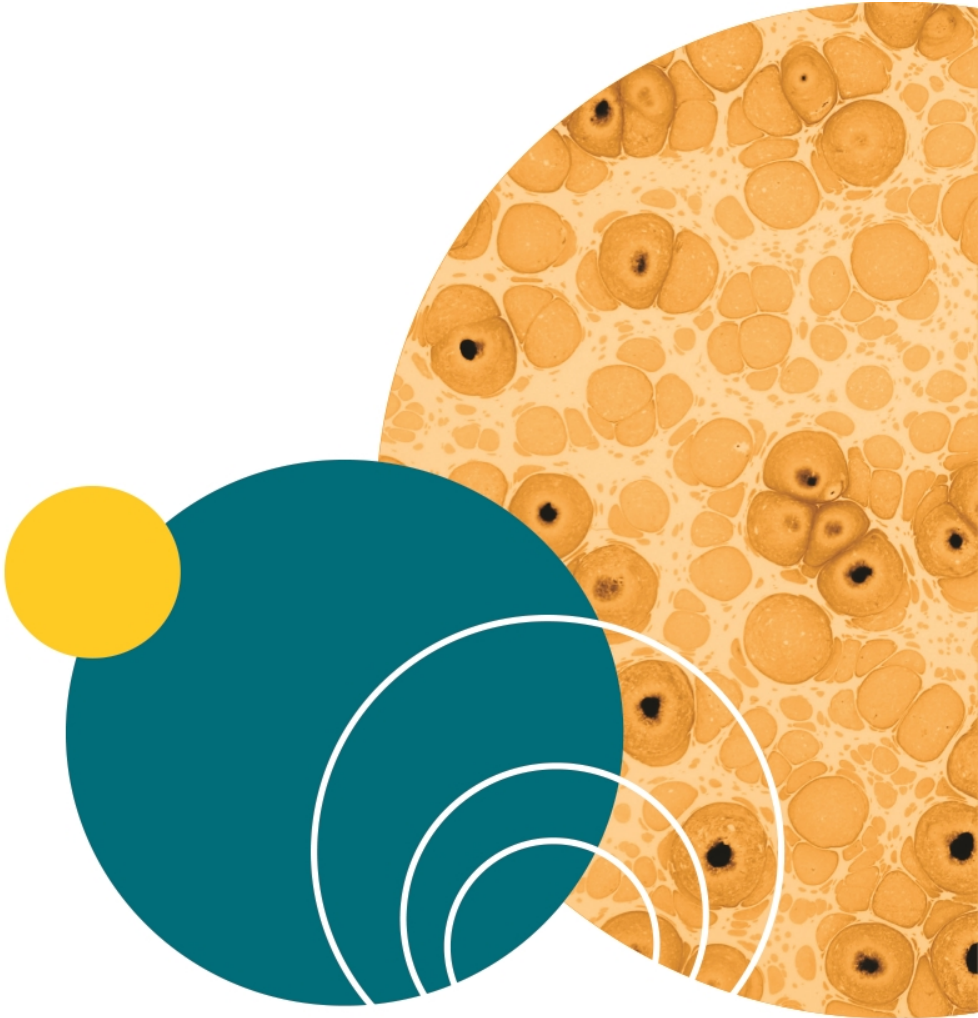
The following new features are included in MultiClamp Commander Software version 2.2.

Native 64-Bit Software

MultiClamp Commander Software version 2.2 runs on 64-bit Windows 7 or Windows 10 operating systems only. Includes the AxMultiClampMsg.dll file for use with 64-bit third-party software.



Note: MultiClamp Commander Software version 2.1.0.16 is available with the same functionality to run on 32-bit Windows 7 or Windows 10 operating systems.



Contact Us

Phone: [+1-800-635-5577](tel:+18006355577)
Web: moleculardevices.com
Email: info@moldev.com

Visit our website for a current listing of worldwide distributors.