

Arcturus^{XT}™ Laser Capture Microdissection System Operating Software Installation Instructions

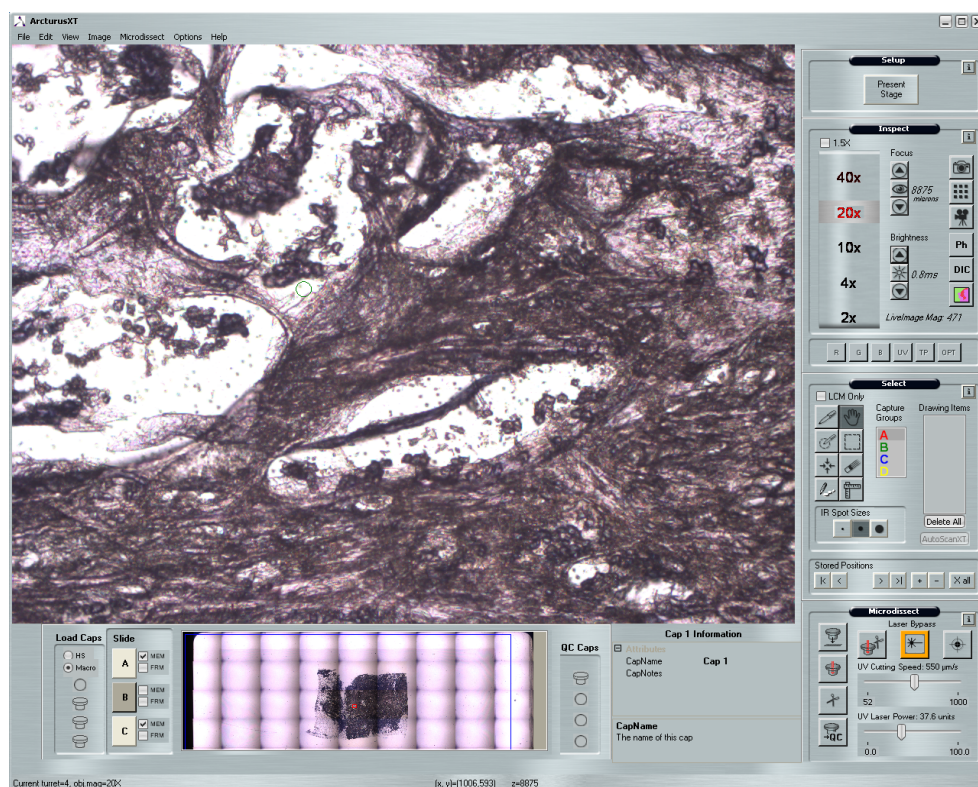
To download updated software, follow these instructions. For best results, do not deviate from any of these steps. If you have any questions, contact technical support at 1-800-831-6844, option 5.

1	Delete previous versions	Before downloading the new version of Arcturus ^{XT} operating software from the website, delete any previous versions of the operating software from your flash drive.
2	Download latest version	Go to the Applied Biosystems website to download and save the Arcturus ^{XT} version 3.2.exe (or newest) file to the flash drive: http://www.appliedbiosystems.com .
3	Locate folder	On the C-drive of your computer, locate the "ArcturusXT" folder in the following location: C:\Program Files\LifeTechnologies. Note: If you have version 3.1 (or older) installed on your system, the file path will be C:\ProgramFiles\MolecularDevices.
4	Copy folder	Copy the "ArcturusXT" folder by right-clicking on the folder and selecting Copy from the pop up window. Do <i>not</i> create a copy of the folder by simply re-naming it.
5	Paste new folder	Right-click anywhere on the window and select Paste . A new folder appears titled "Copy of ArcturusXT".
6	Rename folder	Rename this new folder "Copy of ArcturusXT 3.1.0.0" (use the old version number).
7	Select Set Program Access and Defaults	Open the Start menu located in the lower-left corner of your desktop, and click on Set Program Access and Defaults .
8	Select Change or Remove Programs	In the Add or Remove Programs window, click Change or Remove Programs in the column on the left side of the window. The right side populates.
9	Click Remove button	Click on "ArcturusXT", and then click the Remove button. Do <i>not</i> delete the "ArcturusXT" folder on the C-drive by right-clicking on the folder and selecting Delete . The desktop shortcuts for the Arcturus ^{XT} software and for the Hardware (HW) Configurator will be missing, but the "ArcturusXT" folder will still be on the C-drive. Do not delete this folder or any of its content. The existing files in this folder contain current instrument configurations and calibrations.
10	Download new version	Download the new version.exe file from your flash drive and follow the prompts.
11	Select Everyone	When a dialog box appears asking who this installation is for, select Everyone .
12	Follow prompts	Follow the prompts to complete the software installation. When you have finished, the shortcuts for the Arcturus ^{XT} operating software and the Arcturus ^{XT} Hardware Configurator will again appear on the desktop.

13 Initialize software

Initialize the Arcturus^{XT} operating software and click on **Help**, and then **About** in the menu bar to verify that the new software version has been installed. The primary screen should look like the figure below.

This is the primary screen for the Arcturus^{XT}™ LCM System version 3.2, issued November of 2010.



For Research Use Only. Not intended for any animal or human therapeutic or diagnostic use.

Limited Use Label License: Product Disclaimer

The products in this Bulletin may be covered by one or more Limited Use Label License(s). Please refer to the respective product documentation or the Applied Biosystems website under www.appliedbiosystems.com for the comprehensive license information. By use of these products, the purchaser accepts the terms and conditions of all applicable Limited Use Label Licenses. These products are sold for research use only, and are not intended for human or animal diagnostic or therapeutic uses unless otherwise specifically indicated in the applicable product documentation or the respective Limited Use Label License(s). For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

Purchase of this software product alone does not imply any license under any process, instrument or other apparatus, system, composition, reagent or kit rights under owned or otherwise controlled by Life Technologies, either expressly, or by estoppel.

©2010 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

Part Number 4458766 Rev. B 11/2010

