



CloneSelect Imager FL Software Release Notes

This Release Notes document is distributed as an addendum for the initial release of the CloneSelect® Imager FL software, version 1.8.21.

This version of the CloneSelect Imager FL Software is installed by your Molecular Devices representative on a computer supplied by Molecular Devices. The operating system of this computer is currently Windows 10.

Known Issues

The following is a known limitation of the CloneSelect Imager FL software.

When you run the Image Microplate Multi Channel process for a full plate with wells that contain more than 2000 objects, this version of the software cannot handle the volume of data collected and an error will result.

We are currently working to resolve this issue for a future version of the software.

Get Help

For more information about how to use the CloneSelect Imager FL software, see the CloneSelect Imager FL User Guide or the application help.

The most up-to-date user guides are available on the Molecular Devices Knowledge Base at: support.moleculardevices.com.

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.

Please have the instrument serial number (on the rear of the instrument), or reagent kit lot number, and any related sample data files available when you call.

