



## Action Required for ImageXpress® Micro XL and XLS System Owners!

A device driver must be updated before proceeding with the MetaXpress® Software installation if your instrument is one of the following:

- ImageXpress Micro XLS System
- ImageXpress Micro XL System with serial number 139000 and higher
- ImageXpress Micro XL System with a replacement camera

### To update the driver:

1. Visit the Molecular Devices Support site: [www.moleculardevices.com/support](http://www.moleculardevices.com/support)
2. Click **Visit the Knowledge Base**.
3. Click **Resource Center**.  
You are not required to log in to the knowledge base.
4. Search for article number **20092**.
5. Download the driver: **3.11.10015.0.zip**
6. Unzip the downloaded file.
7. Run **AndorSDK3Setup-3.11.10015.0.exe**.

### 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>



Required Driver Update

1



## Action Required for ImageXpress® Micro XL and XLS System Owners!

A device driver must be updated before proceeding with the MetaXpress® Software installation if your instrument is one of the following:

- ImageXpress Micro XLS System
- ImageXpress Micro XL System with serial number 139000 and higher
- ImageXpress Micro XL System with a replacement camera

### To update the driver:

1. Visit the Molecular Devices Support site: [www.moleculardevices.com/support](http://www.moleculardevices.com/support)
2. Click **Visit the Knowledge Base**.
3. Click **Resource Center**.  
You are not required to log in to the knowledge base.
4. Search for article number **20092**.
5. Download the driver: **3.11.10015.0.zip**
6. Unzip the downloaded file.
7. Run **AndorSDK3Setup-3.11.10015.0.exe**.

### 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>



Required Driver Update

1

