



# EVOS™ XL Core

Imaging System for Transmitted  
Light Applications

Catalog Number AMEX1000

Doc. Part No. ZP-PKGA-0583

Pub. No. MAN0007718

Rev. A.0

# Overview

The EVOS™ XL Core Imaging System has two types of controls: manual and onscreen. Manual controls include the illumination wheel, focus knobs, Freeze and Save buttons, objective turret, and the optional mechanical stage X-Y axis knobs. Onscreen controls appear at the bottom of the screen when the mouse rolls over the controls area. Using a mouse is optional; you can operate the EVOS™ XL Core Imaging System using only manual controls.

**Note:** The EVOS™ XL Core Imaging System User Guide (Pub. No. MAN0007989) is provided on the USB flash drive. It is also available for download from the EVOS™ XL Core Imaging System product page at [thermofisher.com](http://thermofisher.com).

## Operating environment

- ▶ Place the EVOS™ XL Core Imaging System on a level surface away from vibrations from other pieces of equipment.
- ▶ Allow at least 5 cm (2 in) of free space at the back of the LCD housing to allow for proper ventilation and prevent overheating of electronic components.
- ▶ Set up the instrument away from direct light sources, such as windows. Ambient room lighting can enter the imaging path and affect the image.
- ▶ Operating temperature range: 4°–32°C (40°–90°F).
- ▶ Relative humidity range: 30–90%.



**Note:** Onscreen controls are available only when mouse is installed. Roll over bottom of screen with mouse to activate controls.

## Dimensions

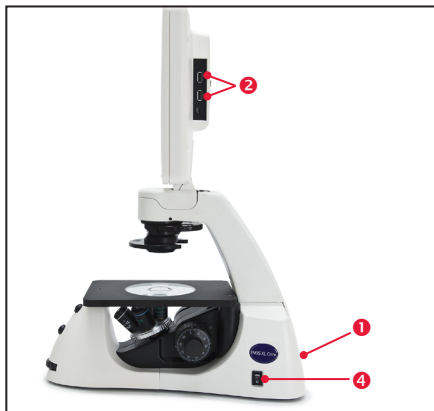
The EVOS™ XL Core Imaging System will fit in cell culture hoods that are at least 18 inches (457 mm) deep. If your cell culture hood is smaller, it may be necessary to turn the instrument at a slight angle to fit.

|                   |                   |
|-------------------|-------------------|
| DEPTH             | 16.0 in (40.6 cm) |
| WIDTH             | 12.5 in (31.8 cm) |
| HEIGHT, TRANSPORT | 13.5 in (34.3 cm) |
| HEIGHT, DISPLAY   | 21.0 in (55.3 cm) |

# Setup

1. Connect the **power adapter** to the **power jack 1** on the back of the instrument base and plug the cord into an outlet.
2. Plug the **wireless mouse receiver** into one of the **USB ports 2** located on the right side of the LCD screen. Install the battery and switch ON the mouse.
3. *Optional:* If the instrument is equipped with a mechanical stage, install the mechanical stage following the instructions provided in the user guide.
4. Unpack the **glass stage plate 3** and insert it into the stage opening (if a mechanical stage is installed, move it all the way back to expose the stage opening).
5. Switch the instrument ON using the **power switch 4** on the right side of the instrument. Startup takes approximately one minute. If the display remains black after startup, adjust the illumination.

**Note:** For detailed setup instructions, refer to the *EVOS™ XL Core Imaging System User Guide*.



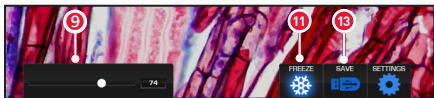
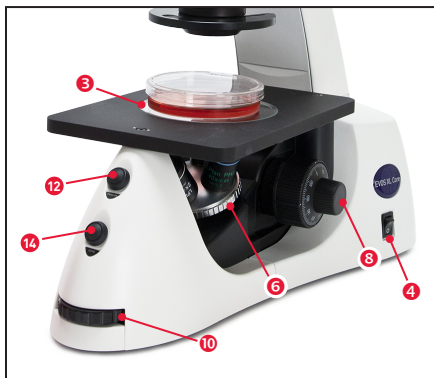
## Basic operation

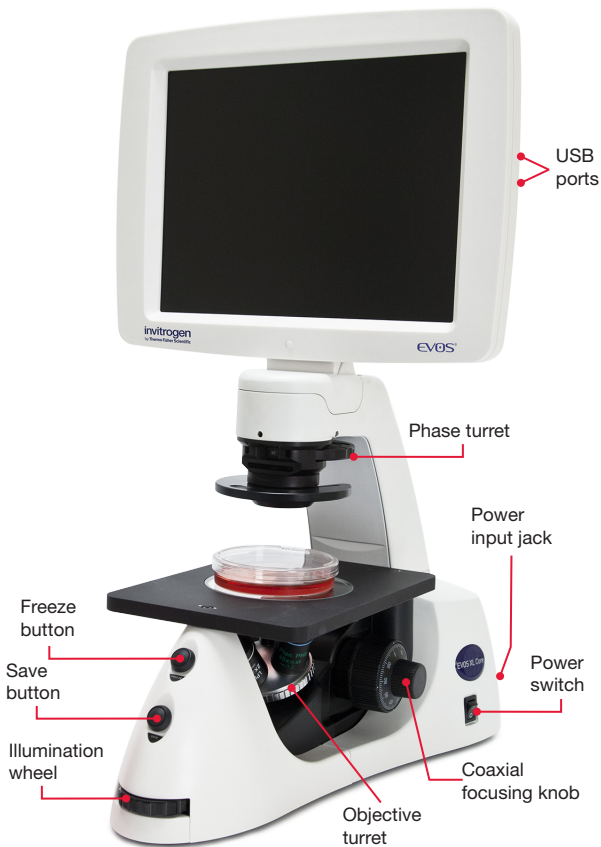
1. Turn on the instrument using the **power switch 4** on the right side of the base.
2. Plug a **USB flash drive** into one of the **USB ports 2** on the right side of the instrument.
3. Place the sample on the **glass stage plate 3**. If the instrument is equipped with a mechanical stage, secure a vessel with the **stage clip 5** to hold the sample over the stage plate.

**Note:** Place slides with coverslips face up.

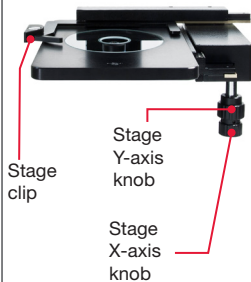
4. Set the **objective turret 6** to the desired magnification.
5. Turn the **phase turret 7** to the position that corresponds to the selected objective and contrast method.
6. Bring the sample into focus using the **coaxial focus knob 8**.
7. Adjust the illumination intensity if necessary, using the **onscreen illumination slider 9** or the **illumination wheel 10**.
8. Use the onscreen **Freeze/Live button 11** or the manual **Freeze button 12** to acquire the image.
9. Use the onscreen **Save button 13** or the manual **Save button 14** to save the image.
10. Click the onscreen **Freeze/Live button 11** or the manual **Freeze button 12** to return to live viewing.

**Note:** The Freeze/Live button toggles between Freeze and Live.





#### Mechanical Stage Controls\*



\* The attachable Mechanical Stage (Cat. No. AMEP4712) is an optional feature. See the *EVOS™ XL Core Imaging System User Guide* for installation and removal instructions.

## Limited Product Warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at [www.thermofisher.com/us/en/home/global/terms-and-conditions.html](http://www.thermofisher.com/us/en/home/global/terms-and-conditions.html). If you have any questions, please contact Life Technologies at [www.thermofisher.com/support](http://www.thermofisher.com/support).

Information in this document is subject to change without notice.

## DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Revision history Pub. No. MAN0007718

| Revision | Date            | Description              |
|----------|-----------------|--------------------------|
| A.0      | 02 October 2017 | Rebrand document.        |
| 1.0      | 18 March 2013   | Basis for this revision. |

## Important Licensing Information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.



Manufacturer: Life Technologies Corporation | 22025 20th Ave SE St #100 | Bothell, WA 98021

## Trademarks

All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

© 2017 Thermo Fisher Scientific Inc. All rights reserved.

[thermofisher.com/support](http://thermofisher.com/support) | [thermofisher.com/askaquestion](http://thermofisher.com/askaquestion)

[thermofisher.com](http://thermofisher.com)

02 October 2017

**ThermoFisher**  
S C I E N T I F I C