## invitrogen



# **EVOS<sup>™</sup> Onstage Incubator**

Catalog Number AMC1000

Doc. Part No. ZP-PKGA-0633 Pub. No. MAN0008920 Rev. A.0

> Thermo Fisher S C I E N T I F I C

For Research Use Only. Not for use in diagnostic procedures.

## Product description

The EVOS<sup>™</sup> Onstage Incubator (Cat. No. AMC1000) is an optional accessory for the EVOS<sup>™</sup> FL Auto and EVOS<sup>™</sup> FL Auto 2 Imaging Systems. It allows you to incubate your cells on the automatic X-Y axis stage to capture images from the same sample over long periods of time and to record time lapse movies.

The EVOS<sup>TM</sup> Onstage Incubator consists of a Environmental Chamber that is placed on the automatic X-Y stage of the imaging system and a separate Control Unit that supplies the power and gas (air or air-CO<sub>2</sub> premix, CO<sub>2</sub>-only, and N<sub>2</sub>-only). The EVOS<sup>TM</sup> Onstage Incubator is controlled by the same software that controls the imaging system.

## Standard items included

- · Stagetop environmental chamber
- Control unit
- Master stage plate
- · Vessel holder for multi-well plates
- Light shield retainer (included only when required)
- Cable with 6-pin connector
- Cable, USB A-to-B
- Heated hose with temperature control
- Gas line, 1/8" ID, 1/4" OD
- Push-to-connect gas line adaptor (3 each)
- Standard-head open-end wrench
- Hex screw driver
- Power Cord, Type A (North America)

Note: For more information on the EVOS<sup>™</sup> Onstage Incubator, refer to the EVOS<sup>™</sup> FL Auto or the EVOS<sup>™</sup> FL Auto 2 Imaging System User Guide, available for download at thermofisher.com.

## Install the master plate

 Remove the X-Y stage base plate 1 from the X-Y stage 2 by unscrewing the four 3.0-mm screws on the base plate (indicated by red arrows).



 Secure the onstage incubator master plate (2) to the X-Y stage (2) using the four thumb screws (indicated by red arrows).



### Assemble the environmental chamber

- Place the incubator chamber ④ on the onstage incubator master plate ⑤ and secure it in place using the four 2.0-mm hex screws (indicated by red arrows).
- Attach the vessel holder/adaptor (5) to the incubator chamber (4) using the four thumb screws.
- Place the heated glass lid () with the no-fog glass window on the incubator chamber (). The heated glass lid is guided and held secure in its place by the two magnets on its rim.
- Place the light shield 
   with tinted plastic window on top of the heated glass lid. Use of the light shield is required for fluorescence imaging applications.
- If desired, place the light shield cover (3) on the light shield for fluorescence imaging applications.
- 6. If required, secure the Stagetop Environmental Chamber to the X-Y stage using the light shield retainer (not shown).





## Setup for operation

- Plug power cord into the power input jack 

   on the back of the Control Unit and the wall outlet.
- 2. Plug USB cable into the USB control cable jack (10) on the Control Unit and the USB port on the computer.
- **3.** Connect each gas line to the appropriate gas tank by pushing the tubing into the PTC (push-to-click) connector until it clicks into place. Pull on tubing slightly to ensure a tight connection.
- **4.** Attach the gas lines to the Control Unit via the PTC connectors for the appropriate gas intake port.
  - If using pre-mixed air, attach to Port 1: Air In ①
  - If using compressed air and CO<sub>2</sub>, attach to Port 1: Air In ① and Port 3: CO<sub>2</sub> In ③
  - For oxygen displacement, attach to Port 1: Air In 1 and Port 2: N<sub>2</sub> In 12
- 5. Plug the 6-pin sensor data cable <sup>(2)</sup> to the Stagetop Environmental Chamber and the appropriate input jack on the Control Unit.
- 6. Assemble the water reservoir and add warm water to the max fill line through the fill hole (3). Do not overfill the water reservoir.
- 7. Place the water reservoir into the Control Unit with the fill holes to the front and close the lid <sup>(1)</sup>.
- 8. Attach the heated hose 😗 between the Stagetop Environmental Chamber and the Control Unit.
- 9. Plug the hose heater cable (1) to the connector on the heated hose.

Note: For detailed instructions on using the EVOS<sup>™</sup> Onstage Incubator, including configuring gas connections and calibrating the oxygen sensor, see the EVOS<sup>™</sup> FL Auto or the EVOS<sup>™</sup> FL Auto 2 Imaging System User Guide.









#### 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-andconditions.html. If you have any questions, please contact Life Technologies at 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. MAN0008920

Revision	Date	Description
A.0	23 June 2017	Rebrand document, replace images with EVOS <sup>™</sup> FL Auto 2.
1.0	17 July 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 | 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 | thermofisher.com/askaquestion

thermofisher.com

