

E-Gel® Imager Gel Documentation System

Cat. nos. 44666611, 44666612, 44666613

Rev. Date: 6 August 2011

Part no. 25-1071 MAN0004953

Directions for installation of the E-Gel® Imager system are described below. For additional details on installation, image capture, and image analysis, refer to the **E-Gel® Imager Gel Documentation System** manual included in the software CD, or contact Technical Support.

E-Gel® Imager Camera Hood



E-Gel® Imager Light Base



E-Gel® Imager Adaptor Base



Kit contents

Each E-Gel® Imager system kit contains an E-Gel® Imager Camera Hood, an E-Gel® Imager Base, an E-Gel® Imager Universal Filter, an E-Gel® iBase™ Light Diffuser, a USB Software Activation Dongle for quantitation software, and two CDs containing the software and manuals for use with the system. See the **E-Gel® Imager Gel Documentation System** manual for more information, descriptions of parts, and unpacking instructions.

Installation overview

The installation procedure for the E-Gel® Imager Gel Documentation System can be completed in five steps:

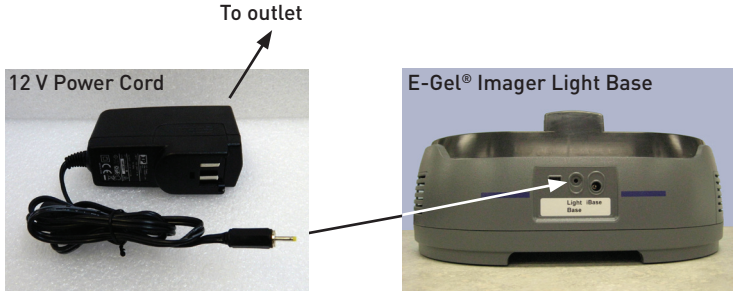
1. Hardware installation
2. GelCapture software installation
3. Camera driver installation
4. Camera calibration
5. GelQuant Express software installation

Intended Use

For research use only. Not intended for any animal or human therapeutic or diagnostic use.

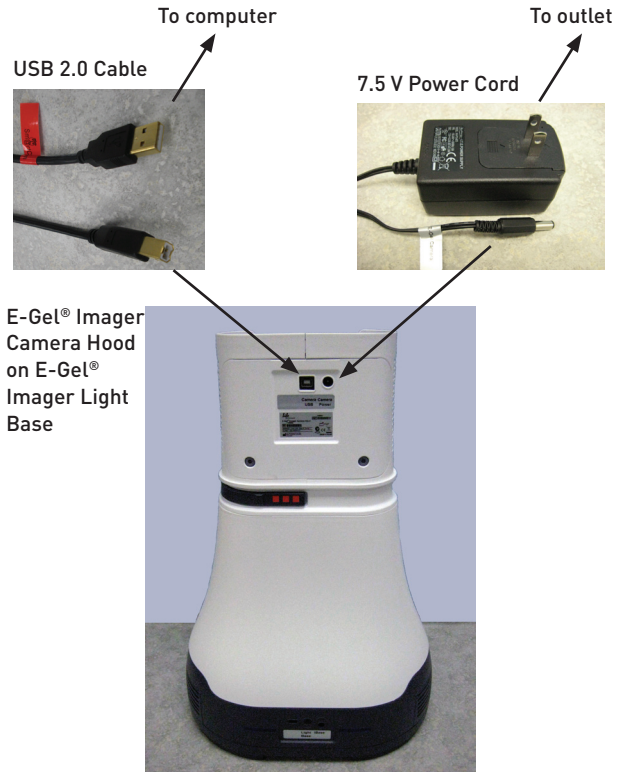
E-Gel® Imager Light Base installation

1. Place the E-Gel® Imager Light Base on a flat surface.
2. Plug the light base power cable into the light base power port.



E-Gel® Imager Camera Hood installation

1. Place the E-Gel® Imager Camera Hood on top of the E-Gel® Imager Light Base.
2. Plug the E-Gel® Imager Camera Hood power supply cable into the camera hood power port.
3. Plug the USB 2.0 cable into the camera USB port of the camera hood and connect the other end into the USB 2.0 port of your computer.



Minimum PC requirements	<p>To install and operate the E-Gel® Imager system and software, your PC must meet the following minimum requirements:</p> <ul style="list-style-type: none">• Intel® Core™ 2 Duo processor, 1.8 GHz.• Minimum 1 GB RAM of memory (2 GB recommended).• 32 bit Windows XP Pro (English version, SP 3) OR Windows 7 Professional (English version) operating system.• Minimum monitor resolution of 1024 × 768 pixels.• A free USB 2.0 port (Not compatible with USB 1.0 or 1.1).
Software installation guidelines	<ul style="list-style-type: none">• Log in as user with Administrator privileges (for installation only). If you do not have Administrator privileges, contact your IT department for help.• Change PC power setting to “Never Hibernate”.• Software installation requires .NET Framework 3.5 for GelCapture, and .NET Framework 4 for GelQuant Express. The .NET Framework installation process is automatically launched if not already present on your computer.
GelCapture software installation	<ul style="list-style-type: none">• Install GelCapture software before turning on the E-Gel® Imager and installing camera drivers.• Follow the instructions provided by the installation wizard to install the software for Windows XP Professional.• For manual installation instructions to install the software for Windows 7 Professional, refer to the E-Gel® Imager Gel Documentation System manual provided on the software CD.
Camera driver installation	<ul style="list-style-type: none">• Before installing the camera drivers, make sure that the GelCapture software is installed on the computer.• Follow the instructions provided by the installation wizard to install the drivers. Refer to the E-Gel® Imager Gel Documentation System manual for details and manual installation instructions.
Camera calibration	<p>Calibrate the camera in the E-Gel® Imager immediately after hardware set up and software installation by launching the GelCapture software. Follow the installation wizard to perform the camera calibration process.</p>
GelQuant Express software installation	<ul style="list-style-type: none">• Follow the instructions provided by the installation wizard to install the software for Windows XP Professional.• For manual installation instructions to install the software for Windows 7 Professional, refer to the GelQuant Express Analysis Software manual provided on the software CD.• The GelQuant Express Software Activation Dongle must be inserted into the computer to use GelQuant Express. If the software activation dongle is not detected, a warning message will appear.

Observation	Cause	Solution
E-Gel® Imager does not turn on after pressing either camera hood or E-Gel® Imager Base ON/OFF button	Loose power supply cable	Verify that the camera hood power supply and light base power supply connections are secure and properly connected to an electrical outlet.
	Loose internal power supply button cables	Contact technical support.
The camera is not recognized	Camera not turned on	Make sure the camera is turned on.
	Cables not properly connected	Verify that the camera USB 2.0 and power supply cables are connected properly.
	Camera drivers did not install successfully	Uninstall GelCapture, then reinstall GelCapture and the camera drivers.
No image on screen	E-Gel® Imager Light Base not turned on	Make sure the E-Gel® Imager Light Base is turned on.
	Possible time delay each time images are changed	Wait a few seconds for images to be displayed.
A black image is seen on screen	Exposure time too short	Increase the exposure time.
	The iris is closed	Turn the Manual Iris Dial to increase the aperture.
Images out of focus	Incorrect focus settings	Adjust the focus settings using the Manual Focus.
Improper illumination intensity	Button is receiving incorrect voltage	Contact technical support.
	Low incoming illumination source voltage	Contact technical support.
Duration of E-Gel® Safe Imager™ transillumination too short	E-Gel® Safe Imager™ transilluminator turned on with short press (30 second transillumination)	Press the button on the E-Gel® Safe Imager™ for 2 seconds. The transilluminator turns on for 5 minutes before automatic shutdown.

Limited Use Label License

Limited Use Label License: Research Use Only: The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial services of any kind, including, without limitation, reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For technical support, email tech_support@invitrogen.com.
 For country-specific contact information, visit www.invitrogen.com.
 To reorder visit www.invitrogen.com

