



FlashGel™ Camera and Capture Software

Software Install Instructions and User's Guide for version 3.1

Compatible with FlashGel™ Camera serial numbers LNZ/2012 and higher

For support on units with serial numbers smaller than LNZ/2012, please contact Scientific Support.

Introduction

Capture data from the FlashGel™ System right at your bench top. Say good bye to darkrooms and UV light. Complete separation and documentation safely, at your bench, in minutes. This simple digital camera in an enclosed hood connects directly to your Windows® laptop, PC, or PDA via USB cable.

Real-time separation and documentation

- Complete your gel run and image capture in just 5 minutes.
- Watch DNA migrate – on the FlashGel™ Dock or on your computer screen.
- Photograph gels at the bench – without UV light

The FlashGel™ Camera offers

- Sharp, clear, high-resolution images
- Simple user interface
- Small, compact design

Quick Start Guide

Note: Do not plug camera into computer until after software is installed.

1. Insert install CD into drive.

Note If AutoRun is disabled on your machine, open CD folder and execute the Setup.exe file.

2. Walk through setup wizard (Figure 1).
 - a. Select destination location.
 - b. Select start menu folder.
 - c. Click Install.

Note: Three drivers will be loaded during the install process. Some operating systems may have an additional prompt to install these drivers. You'll want to install these.

Figure 1: Setup Wizard



3. ****Important:** After install is complete, also run update.exe from the CD by opening the CD folder and doubleclicking the update.exe file.

Using FlashGel™ Capture

Starting the application

Please ensure the camera is connected and the camera hood is placed securely over the FlashGel™ Dock.



Figure 2: FlashGel™ Capture Desktop Icon

Launch the application from the desktop icon or Start Menu. (Figure 2).

Viewing a FlashGel™ Image

As soon as the application starts, you will see the software interface. (Figure 3).

Figure 3



Note: The camera is setup for the low-light conditions expected when viewing a FlashGel™ Cassette.

Note: If you remove the hood, the image will appear bright.

Turn on the FlashGel™ Dock light. The camera will automatically adjust its exposure settings to show an optimized image. You will then be able to monitor the separation of fragments during the course of the experiment.

Figure 4



Capturing and saving an image

To capture a still image at any time, simply click the Capture button. (Figure 5).



Figure 5: Capture Button

This action will show a Save As dialog and give you the opportunity to enter a filename and choose between three image types; BMP, JPG, or TIF.

On subsequent occasions (and even after restarting the application), FlashGel™ Capture will remember the name and location of the last file you saved. Furthermore, the program will append a sequential number to the last filename you saved, meaning that you simply have to click Save to continue. e.g. if your initial filename was “Test”, subsequent filenames would be “Test1”, “Test2”, “Test3”....

Printing an image

To print an image at any time, simply click the Print button. (Figure 6).



Figure 6: Print Button

This action will show a print dialog that will enable you to select the printer of choice and subsequently print the image.

Camera Specifications

Hood Dimensions:	10 cm [W] x 11 cm [L] x 16 cm [H]
Camera Type:	Digital/CMOS
Image File Type:	.JPG, .TIF, .BMP
Image Resolution:	1280 x 960
Bit Depth:	8 bit

Software/Hardware Requirements

Windows® XP (SP2 or higher), Windows Vista® or Windows® 7 (32-bit or 64-bit)
1 GHz
512MB RAM
USB 1.1 port (2.0 recommended)

Ordering Information

Cat. No.	Product	Size/Format
57040	FlashGel™ Camera	One size
57065	FlashGel™ Camera Pack	Includes FlashGel™ Dock and FlashGel™ Camera.
57062	FlashGel™ Device Pack	Includes FlashGel™ Dock, FlashGel™ Power Supply, and FlashGel™ Camera.

Related Products

FlashGel™ Kits

Cat. No.	Product	Size/Format
57025	FlashGel™ Dock	One size
57063	FlashGel™ DNA Kit	Includes FlashGel™ DNA Cassettes 1.2% 12+1 well single tier 9pk, FlashGel™ Loading Dye, and FlashGel™ Marker 100 bp – 4 kb.
57024	FlashGel™ RNA Kit	Includes FlashGel™ RNA Cassettes 1.2% 12+1 well single tier 9pk, RNA Marker, Sample Buffer, and Molecular Biology Water.
57064	FlashGel™ Recovery Kit	Includes FlashGel™ Recovery Cassettes 1.2% 8+1 well double tier 9pk, FlashGel™ Recovery Buffer, FlashGel™ Loading Dye, FlashGel™ QuantLadder, and Visualization Glasses.

Warranty and Liability

This product was produced utilizing the highest practical standards of materials, workmanship, and design. Lonza Rockland, Inc. warrants that the product has been tested and will meet or exceed published specifications. This warranty is valid only if the product has been operated and maintained according to the instructions provided.

Lonza Rockland, Inc. warrants this product to be free from defects in materials and workmanship under normal service for one year from date of shipment. If the product proves defective during this period, Lonza Rockland, Inc., will repair or replace it at Lonza Rockland, Inc.'s option, free of charge, if returned to Lonza

Rockland, Inc. postage prepaid. This warranty does not cover: damage in transit, damage caused by carelessness, misuse or neglect, normal wear through frequent use, damage caused by solvent corrosion, damage caused by improper handling or user alteration, nor unsatisfactory performance as a result of conditions beyond Lonza Rockland, Inc.'s control. Lonza Rockland, Inc. shall not be liable for any damage or injury to persons or property of any kind arising out of the purchase or use of the product. Any claim made by the buyer against Lonza Rockland, Inc. with respect to the product shall be waived by buyer unless made within thirty days of delivery. In no event shall Lonza Rockland, Inc.'s liability with respect to any such claim exceed the purchase price actually paid by buyer. Lonza Rockland, Inc., shall in no event be liable for incidental nor consequential damages, including without limitation, lost profits, loss of income, loss of business opportunities, loss of use and other related damages, however caused, nor any damage arising from the incorrect use of the product. Each party hereto shall be excused from liability for failure to perform due to any cause beyond its reasonable control, except that this clause shall not operate to relieve buyer of the obligation to pay the purchase price for merchandise delivered.

License and Trademark Information

Some components and technology of the FlashGel™ System are sold under licensing agreements. The nucleic acid stain in this product is manufactured and sold under license from Molecular Probes, Inc. and the FlashGel™ Cassette is sold under license from Invitrogen IP Holdings, Inc. and is for use only in research applications or quality control, and is covered by pending and issued patents. Dark Reader is a trademark of Clare Chemical Research, Inc. The FlashGel™ Dock technology contains Clare Chemical Research, Inc. Dark Reader® Transilluminator technology and is covered under US Patents 6,198,107; 6,512,236; and 6,914,250. The electrophoresis technology is licensed from Temple University and is covered under US Patent 6,905,585.

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

Manufactured for Lonza Rockland, Inc.

Microsoft and Vista are trademarks of Microsoft Corporation. All other trademarks herein are marks of the Lonza Group or its affiliates.

© 2012 Lonza Rockland, Inc.
All rights reserved.