

iGlow™ Gel Documentation System

Catalog Numbers IGLOW and IGLOW220

Pub. No. MAN1002055 Rev. A



WARNING! Read the User Guide and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves. The User Guide is available at thermofisher.com/support.

Product description

The Invitrogen™ iGlow™ Gel Documentation System enables fast, user-friendly imaging of nucleic acid and protein gels.

Product Information

Table 1 iGlow™ Gel Documentation System – Shipment Contents

Category	Components list	Quantity	Catalog Number
Instrument	iGlow™ Gel Documentation System, 110V	1	IG0001
	iGlow™ Gel Documentation System, 220V		IG0002
Accessories	UV Tray iGlow™ ^[1]	1	IGA011
	Blue Tray iGlow™ ^[1]		IGA012
	White Tray iGlow™ ^[1]		IGA013
	Tray Holder iGlow™ ^[1]		IGA014
	Safety Guard (UV Shield iGlow™) ^[1]		IGA015
	Power cord (region specific)		–
Documents	Packaging Checklist	1	–
	Quick Reference Guide		–
(Optional) Accessories	Thermal Printer ^[1,2]	N/A	A45235
	USB Wi-Fi Module ^[1,2]		A26774

^[1] Can be purchased separately.

^[2] Not included in the shipment.

Note: For more detailed information and safety symbols on this instrument, refer to iGlow™ User Guide (Pub. No. [MAN1001999](#)).

Getting Started

1. Remove the instrument from the packaging.
2. Place the system on a clean, stable surface.
3. Plug in the power cord provided.
4. Turn on the imager using the **power switch on the top-right side** of the instrument.

User Accounts & Initial Setup

Default Accounts

The system includes two default user accounts:

Default Accounts	User ID	Password
Administrator	Administrator	AA
Standard User	Public	aa

IMPORTANT! Security Requirement: Both default accounts must reset their passwords at first login.

First-time setup (Administrator)

1. Sign-in using the Administrator account.
2. Update the default password.
3. Set the password policy (applies to all users).

Note:

- We recommend configuring the password policy during installation, as it applies to all user accounts. Changing the policy later may require existing users to reset their passwords.
- Other administrative responsibilities include configuring system settings, such as time zone, date format, and time format.

Creating Additional Users

1. From the welcome page, select **Create**.
2. Enter a **Username**.
3. Enter a **Password** then re-enter the password to confirm.
4. Select **Create** to complete account creation.
5. On the Welcome screen, select the username from the dropdown list.
6. Enter the newly created password to sign in and verify the account.

Note: For details on password requirements, refer to the iGlow™ User Guide (Pub. No. [MAN1001999](#)).

Imaging

Table 2

Illumination	Description	Recommended sample tray
UV	Use UV light to image or excise bands from nucleic acid or protein gels stained with UV excitable stains.	UV Tray
Blue	Use Blue light to image or excise bands from nucleic acid or protein gels stained with Blue excitable stains.	Blue Tray
White	Use white light to image or excise bands from protein gels stained with colorimetric stains.	White Tray

Acquiring an Image

1. From the **Welcome** screen, select your user account and sign-in to access the **Home screen**.
2. Open the drawer and place the desired **sample tray** on the transilluminator.
 - Use the alignment guides to ensure correct positioning.

Note: The system automatically detects the tray and enables the correct illumination source.

3. Place the samples in the **center of the tray**.
4. Close the drawer.
5. Tap the **highlighted illumination mode** to capture the image.
 - The system automatically calculates the optimal exposure time and captures an image.

- The image appears in the **Image Preview** screen.
6. (Optional) Retake images using a manual exposure, or by adjusting the field of view (FOV) or light intensity.
 7. (Optional) Select a predefined **Method** to take a new image using saved imaging conditions.
 8. When satisfied:
 - (Optional) Rename the file.
 - Save the image to a USB, or a Network Drive location.

Gallery

The on-board Gallery contains all image files captured under the logged-in user profile. Image files appear as thumbnails, and users can filter, search, delete, rename, and export files.

Export

The iGlow™ Gel Documentation System supports export of image files to a network drive or USB drive in 16-bit iBA format (a proprietary analysis file format compatible with iBright™ Analysis Software) as well as in publication-quality (TIF and JPG) formats.

Refer to the iGlow™ Gel Documentation System User Guide (Pub. No. [MAN1001999](#)) for additional information on system use, data handling, and maintenance.

Resources

Table 3

iGlow™ Gel Documentation System Docs	User Guide (Pub. No. MAN1001999) Quick Reference Guide (Pub. No. MAN1002055)
iBright™ Analysis Software: iBright™ Analysis Software, Desktop Version	https://www.thermofisher.com/us/en/home/technical-resources/software-downloads/ibright-western-imager.html

Limited product warranty

Life Technologies Corporation and its affiliates warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale at www.thermofisher.com/us/en/home/global/terms-and-conditions.html. If you have questions, contact Life Technologies at www.thermofisher.com/support.



Thermo Fisher Scientific | 3747 N. Meridian Road | Rockford, Illinois 61101 USA
 For descriptions of symbols on product labels or product documents, go to thermofisher.com/symbols-definition.

Revision history: Pub. No. MAN1002055A

Revision	Date	Description
A	13 May 2026	New document for iGlow™ Gel Documentation System.

The information in this guide is subject to change without notice.

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. 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 THE DOCUMENT OR THE PRODUCT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2026 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.