

iBright™ Imaging Systems

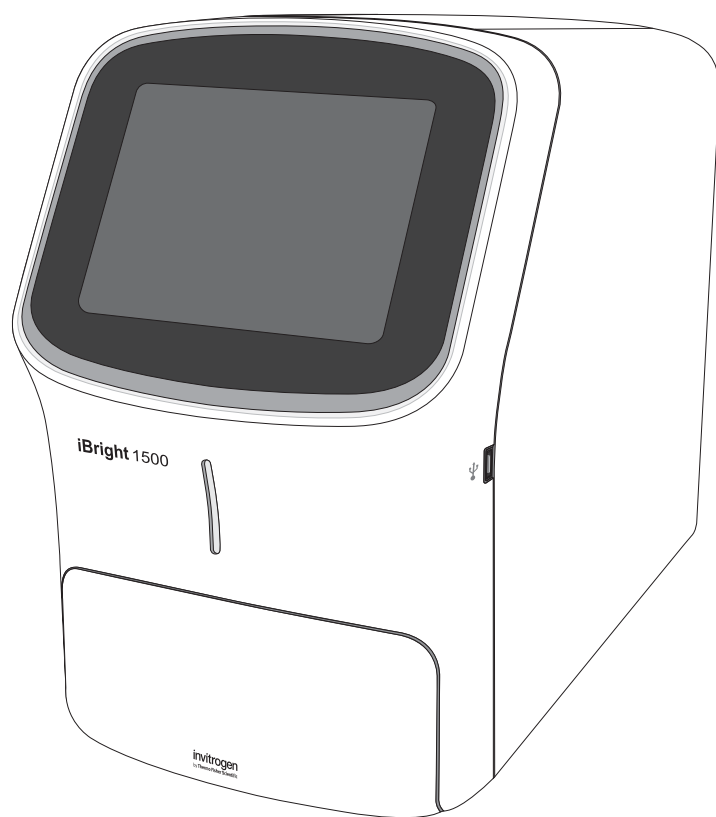
USER GUIDE

For use with the iBright™ CL1500 Imaging System and the iBright™ FL1500 Imaging System

Catalog Numbers A44114 (CL1500) and A44115 (FL1500)

Publication Number MAN0018592

Revision D.0



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



Revision history: Pub. No. MAN0018592

Revision	Date	Description
D.0	31 October 2022	<ul style="list-style-type: none">Updating workflows according to instrument firmware upgrades.Adding SmartRange™ HDR functionality and analysis of E-gel and colony images in Universal mode.
C.0	24 August 2020	Adding SAE chapter due to addition of SAE instrument functionality.
B.0	23 September 2019	<ul style="list-style-type: none">Updated the components list (see Table 1 on page 6).Updated the recommended excitation/emission wavelengths to image fluorescent Western blots (see "Fluorescent Blots (FL1500 model only)" on page 14).Updated the emission filters for the iBright™ FL1500 Instrument (see Table 4 on page 18).Added additional outputs for TPN-normalized data.Added additional information for resetting factory defaults, backing up, and restoring settings (see "Settings" on page 30).
A.0	7 June 2019	New document for iBright™ Imaging Systems.

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 IT.

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.

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

©2022 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■ CHAPTER 1	Product information	7
	Product description	7
	Contents	7
■ CHAPTER 2	Start, sign on, and configure the instrument	9
	Create an instrument profile	9
	Create an instrument account and link to Connect using a cloud account	9
	Mobile device	10
	PC using Connect account	10
	PC using instrument Connect app	10
	Link to Connect using an existing instrument account	11
	If the unit is off	11
	If the unit is on and in sleep mode	12
	If the unit is on and in active mode	12
■ CHAPTER 3	Operate the iBright™ imager	13
	Choose imaging mode and start image capture	13
	Chemi Blots	14
	Image using Chemi Blots	14
	Fluorescent Blots (FL1500 model only)	15
	Image using Fluorescent Blot	15
	Nucleic Acid Gels	17
	Image using Nucleic Acid Gel	17
	Protein Gels	18
	Image using Protein Gels	18
	Universal Mode	19
	Options available in modes	21
	Image Gallery	24
	View images	24

- **CHAPTER 4** Export and analyze images 25
 - Export images 25
 - Export new images 25
 - Export gallery images 25
 - Select an Export file type 26
 - Archive/Backup g2i image files 27
 - Analyze images 28
 - Analyze new images 28
 - Analyze gallery images 30
 - Analyze 96-well E-Gel™ images taken in Universal mode using the E-Gel™
96 application 30
 - Analyze colony images taken in Universal mode using the Colony
Count application 31

- **CHAPTER 5** Maintain the instrument 33
 - Settings 33
 - Manage accounts 35
 - Run instrument diagnostics 36
 - Clean the instrument 36
 - Replace the instrument fuses 37
 - Materials required 37
 - Replace the fuses 37
 - Move the instrument 38
 - Prepare for shipping 38

- **CHAPTER 6** Invitrogen™ iBright™ SAE Software Solution for 21 CFR
Part 11 Support 39
 - Security Audit e-Signature (SAE) for the iBright™ Imager 39
 - Activate your license key 40
 - Enable SAE mode 40
 - Disable SAE mode 41
 - Sign in to the instrument in SAE mode 41
 - Change your password in SAE mode 42
 - SAE user roles 42
 - SAE privileges for default user roles 43
 - Non-SAE user roles 43
 - Acquire images in SAE mode 44

Use the Gallery in SAE mode	44
File status	44
Filter Gallery by file status	44
Archive/Backup g2i image files in SAE mode	44
Commit changes	45
Sign data	46
Unlock files	46
Generate an audit history	47
View the e-Signature history	47
Export image files	47
Generate and export Analysis Report	48
■ APPENDIX A Instrument specifications	49
Instrument dimensions and specifications	49
Electrical requirements	49
Environmental requirements	49
Network requirements	50
Instrument and computer connections	51
■ APPENDIX B Safety	53
Symbols on this instrument	53
Location of safety labels on the instrument	55
Safety information for instruments not manufactured by Thermo Fisher Scientific	55
Instrument safety	55
General	55
Physical injury	56
LED (Light-Emitting Diode)	56
Electrical safety	57
Cleaning and decontamination	58
Instrument component and accessory disposal	58
Safety and electromagnetic compatibility (EMC) standards	59
Safety compliance	59
EMC	59
Environmental design	60
Chemical safety	62
Biological hazard safety	63

- **APPENDIX C** Documentation and support 64
 - Related documentation and references 64
 - Customer and technical support 65
 - Limited product warranty 65



Product information

IMPORTANT! Before using this product, read and understand the information in the “Safety” appendix in this document.

Product description

The iBright™ CL1500 Imaging System allows users to image chemiluminescent Western blots, DNA and RNA gels stained with fluorescent nucleic acid stains, and visible protein gels. The iBright™ FL1500 Imaging System adds the capability of imaging fluorescent Western blots using multiple dyes.

The iBright™ imagers use a simple design and intuitive workflows to deliver high-resolution images. They can be run directly from the touchscreen to start and create images, while also providing on-board software for image analysis. To further analyze data, the instrument can be integrated with the iBright™ Image Analysis Software available for desktop and Connect.

Contents

Table 1 Included imager system parts

Component	Quantity
iBright™ Imaging System	1
Power cord (region specific)	1
Quick Reference Guide	1
Contrast tray	1
iBright™ Imaging System Sample Blot	1
Sample stage/turntable	1
Product Information Sheet	1
Safe Imager™ Viewing Glasses	1


Table 2 Available accessories

Component	Part number
iBright™ System Contrast Tray	A56597
Sample stage	A33829
High-Power USB Wi-Fi Module	A26774
iBright™ E-Gel™ Tray Adapter	A56600 (11/22) A56599 (48/96)




Start, sign on, and configure the instrument

Create an instrument profile

1. Touch  to sign in to your instrument user account.
2. Touch **Get started**.
3. Touch **Create a Profile**.
4. Enter a **Screen name**.
5. Enter and confirm a four-digit pin.
6. Touch **Create profile**.
7. (Optional) If you want to link your instrument to your Connect account, see “Link to Connect using an existing instrument account” on page 11.



Create an instrument account and link to Connect using a cloud account

1. Touch  to sign in to your instrument user account.
2. Touch **Get Started**.
3. Click on **Connect**.
4. Link your account using one of these methods:
 - **Mobile device** — Connect using a QR generated by the instrument and scanned by the instrument Connect mobile app.
 - **PC** — Connect using a link code that is entered into the instrument Connect app or your online Connect account.
 - **Instrument** — Enter your Connect account information directly on the instrument.

Mobile device

1. Touch **Mobile Device**.
2. Launch instrument Connect mobile app.
3. Sign into your Connect account on the mobile app.
4. Click on **+**.
5. Click on **QR code**.
6. Scan QR code on the instrument.
Your Connect account should now be linked to your instrument account.

PC using Connect account


1. Touch **PC**.
2. Login to your Connect account online.
3. Click on .
4. Click on .
5. Select **iBright** from the drop-down menu.
6. Enter the linking code and click on **Send**.
Your Connect account should now be linked to your instrument account.

PC using instrument Connect app

1. Touch **PC**.
2. Launch the instrument Connect mobile app.
3. Sign into your Connect account on the mobile app.
4. Click on **+**.
5. Click on **Linking code**.
6. Enter the linking code and click **Send**.
Your Connect account should now be linked to your instrument account.

Link to Connect using an existing instrument account



Linking to Connect after account creation can only be done during the export process.

1. Click on .
2. Select an image.
3. Click on **Actions**.
4. Click on **Export**.
5. Click on **Choose destination**.
6. Click on **Cloud**.
7. Click on **Sign in**.
8. Connect using one of the methods outlined in “Create an instrument account and link to Connect using a cloud account” on page 9.


If the unit is off

1. Ensure the unit is plugged in and turn on using the switch on the back of instrument.
The instrument will go directly to the sign-in page.
2. Enter a **Screen name**.
3. Enter and confirm a 4-digit pin.


Note: For those not logged on as a user (guest), acquired images are saved to a guest gallery. The guest gallery is not private. Any logged in user can view and modify images residing in the guest gallery.

4. From the **Welcome** screen, choose the appropriate Mode from the drop-down menu, touch  and place sample in the center of the transilluminator glass.
5. Touch  to close drawer (do not physically push on the drawer). Instrument adjusts and displays a live-view sample image in the viewport.


If the unit is on and in sleep mode

1. Touch the screen to wake.
2. Touch  to sign in to use or set up a user profile. See “Create an instrument profile” on page 9 for instructions on how to create a user profile.



Note: For those not logged on as a user (guest), acquired images are saved to a guest gallery. The guest gallery is not private. Any logged-in user or other guest user can view and modify images residing in the guest gallery.

3. Enter a **Screen name**.
4. Enter and confirm a 4-digit pin.
5. From the Welcome screen, choose the appropriate Mode from the drop-down menu, touch Open Drawer, and place sample in the center of the transilluminator glass.
6. Touch  to close drawer (do not physically push the drawer). Instrument adjusts and displays a live-view sample image in the viewport.

If the unit is on and in active mode

1. If not already signed in, touch  to sign in or set up a new user profile. See “Create an instrument profile” on page 9 for instructions on how to create a user profile.

Note: For those not logged on as a user (guest), acquired images are saved to a guest gallery. The guest gallery is not private. Any logged-in user can view and modify images residing in the guest gallery.



2. From the **Welcome** screen, choose the appropriate Mode from the drop-down menu, touch  and place sample in the center of the turntable glass.
3. Touch  to close drawer (do not physically push the drawer). Instrument adjusts and displays a live-view sample image in the viewport.



Operate the iBright™ imager

Choose imaging mode and start image capture

On the **Welcome** screen:

1. Use the drop-down menu to select the desired mode.
 - **Chemi Blots** to image chemiluminescent substrates.
 - **Fluorescent Blots** to image fluorescent substrates.
 - **Nucleic Acid Gels** to image DNA and RNA gels.
 - **Protein Gels** to image visible and fluorescently stained protein gels.
 - **Universal** to image samples containing multiple readouts, such as chemiluminescence, fluorescence and/or colorimetric stains. Image display is similar to Fluorescent Blots and allows you to assign false color to any sample. Use Colony Count and E-Gel™ 96 applications to provide specialized workflows to ensure proper analysis for these applications.
2. Touch  to open drawer.
3. Place sample in the center of the viewing area.
4. Touch  to close drawer.

Note: In **Chemi**, **Fluorescent Blots**, and **Universal** modes the live view is illuminated using the white Epi-LED by default. To use the green transilluminator LED for illumination for live view and to capture of the membrane layer, touch the box **Use Trans Light** on the live-view screen.

Instrument automatically aligns, zooms, focuses, and acquires a live-view sample image.

Note: If any manual adjustments are required for zoom or focus, touch **More options ▶ Camera** and choose the appropriate menu item.

- To image in Chemi Blots mode, proceed to “Image using Chemi Blots” on page 14.
- To image in Fluorescent Blots mode, proceed to “Fluorescent Blots (FL1500 model only)” on page 15.
- To image in Nucleic Acid Gels mode, proceed to “Image using Nucleic Acid Gel” on page 17.
- To image in Protein Gels mode, proceed to “Image using Protein Gels” on page 18.
- To image in Universal Mode, proceed to “Universal Mode” on page 19.

Chemi Blots

Mode to image chemiluminescent Western blots.

Image using Chemi Blots

1. Touch Smart Exposure.

The imager acquires a series of short exposures, then renders a preview image and recommends an exposure time. This is not a real image.

2. (Optional) Touch More options ▶ Region of Interest. Region of Interest is for directing **Smart Exposure** to a specific region of the sample. Select the area to use as a **Smart Exposure** focal point.

3. (Optional) Adjust the exposure time using the following methods:


- Touch **+** or **-**.
- Finger swipe in the segmented dial.
- Touch the dial center box to select a preset exposure time or touch **Custom** to manually input a time.

The image preview updates in real-time.

4. Touch Enter (for a custom manual input).

5. (Optional) Touch Smart Range™ HDR to capture a high dynamic range image. HDR exposure time will display. Touch **Capture** to acquire the HDR image or uncheck **Smart Range™ HDR** to return to **Smart Exposure**.

Note: This feature is also available after image capture.

6. (Optional) Touch  to lock the camera settings (exposure time and zoom) to acquire additional images using the same settings.

Note: Once lock settings are applied, the sample auto-rotation feature will be disabled after the drawer has opened and closed. This feature is also available after image capture.

7. Touch Capture to acquire an image.

The captured image is displayed and automatically saves to the gallery.

8. If the image is acceptable, touch Export, Gallery, or Analyze. If the image is not acceptable, touch **More options** to optimize the image or touch **Trash** to remove the image, then return to step 2 to adjust exposure conditions.

Fluorescent Blots (FL1500 model only)

Mode to image fluorescent Western blots using a single channel or up to 4 channels using excitation wavelengths (FWHM) of approximately 455-485 nm, 515-545 nm, 608-632 nm, 610-660 nm, and 745-765 nm; approximate emission wavelengths (FWHM) 508-557 nm, 568-617 nm, 675-720 nm, 710-730 nm and 800-850 nm.

Image using Fluorescent Blot

1. Touch **Choose a dye** to select from one to four dyes from the pre-populated list.
Alternatively, to add a dye that is not in the pre-populated list, touch **Custom Dye** at the bottom left-hand corner of the screen. Select a custom dye that has previously been entered or enter the information for a new dye: dye name, emission wavelength, excitation wavelength, and false color.
 - a. To edit an existing dye, touch the dye name button. Touch **Dye** and change the assigned dye to any other pre-populated dye in the fluorescent dye list. Dyes in the pre-populated dye list are categorized by most recently used, name, and emission color.
 - b. Touch **False Color** and change the false color assigned to that dye.

Note: Each dye is assigned a false color to be used in the composite overlay.

 - c. To remove the assigned dye from the channel, touch **Remove Dye** at the bottom left-hand corner of the screen. Touch **Confirm**.
2. Touch **Done**.
3. Touch **Smart Exposure**. Alternatively, for manual exposure, touch **More options ▶ Manual Exposure**. For manual exposures, click on the dye to add a time for each dye selected. The imager will acquire a series of short exposures for each dye type selected. It will then render preview images and a recommended exposure time for each channel. These are not real images.
4. (Optional) Touch **More options ▶ Region of Interest**. **Region of Interest** is for directing **Smart Exposure** to a specific region of the sample. You can select different, same, or no region of interest for each channel.
5. To adjust or set a manual exposure time for any of the channels you are acquiring, select the channel (dye) that you want to change:
 - Touch **+** or **-**.
 - Finger swipe in the segmented dial.
 - Touch the dial center box to select a preset exposure time or touch **Custom** to manually input a time.

The image preview updates in real-time.

- 6.** Touch **Capture** to acquire an image.
The captured image is displayed and automatically saves to the gallery.

- 7.** If the image is acceptable, touch **Export**, **Gallery**, or **Analyze**. If the image is not acceptable, touch **More options** to optimize the image or touch **Trash** to remove the image, then return to step 2 to adjust exposure conditions.

Nucleic Acid Gels

Mode to image DNA and RNA gels stained with fluorescent nucleic acid stains such as Ethidium Bromide, SYBR™ Safe, SYBR™ Green (I or II), SYBR™ Gold and similar products. For imaging high-throughput E-Gel™ 96-well agarose gels, use the **E-Gel™ 96** specialty application in Universal mode.

Image using Nucleic Acid Gel

1. Touch **Smart Exposure**. Alternatively, to set a manual exposure time, touch the center of the dial. The imager acquires a series of short exposures and then renders a preview image and a recommended exposure time. This is not a real image.
2. (Optional) Touch **More options ▶ Region of Interest**. **Region of Interest** is for directing **Smart Exposure** to a specific region of the sample.
3. To adjust or set a manual exposure time, select a method:
 - Touch **+** or **–** within dial.
 - Finger swipe in the segmented dial.
 - Touch the dial center box to select a preset exposure time or touch **Custom** to manually input a time.

The image preview will update in real-time.

4. Touch **Capture** to acquire image with the indicated exposure time. Captured image appears on screen and automatically saves to the gallery.
5. If the image is acceptable, touch **Export**, **Gallery** or **Analyze**. If the image is not acceptable, touch **More options ▶ Image adjust** to optimize the image or touch **Trash** to remove the image, then return to step 2 to adjust exposure conditions.

Protein Gels

- **Protein Gels Visible:** USE THE WHITE SCREEN OR CONTRAST TRAY BETWEEN THE SAMPLE AND TRANSILLUMINATOR GLASS for samples including Coomassie-stained protein gels and silver-stained protein gels.
- **Protein Gels Fluorescent:** PLACE SAMPLE DIRECTLY ON THE TRANSILLUMINATOR GLASS for samples including No-Stain™-labeled protein gels and SYPRO™ dye-stained protein gels.

Image using Protein Gels

1. Select whether you are imaging a visible sample (**Protein Visible**) or fluorescent sample (**Protein Fluorescent**).
2. An auto-zoomed and focused **Live View** of the image is now displayed in the viewport.

IMPORTANT! For visual stained protein on nitrocellulose and PVDF membrane, use **Universal Mode ▶ Visible channel** without the white screen.

3. Touch **Smart Exposure**. Alternatively, to set a manual exposure time, touch the center of the dial. The imager acquires a series of short exposures and then renders a preview image and a recommended exposure time. This is not a real image.
4. (Optional) Touch **More options ▶ Region of Interest**. **Region of Interest** is for directing **Smart Exposure** to a specific region of the sample. You can select different, same, or no region of interest for each channel.
5. To adjust or set a manual exposure time, select a method:
 - Touch **+** or **-** within dial.
 - Finger swipe in the segmented dial.
 - Touch the dial center box to select a preset exposure time or touch **Custom** to manually input a time.

The image preview will update in real-time.

6. Touch **Capture** to acquire image with the indicated exposure time. Captured image appears on screen and automatically saves to the gallery.
7. If the image is acceptable, touch **Export**, **Gallery** or **Analyze**. If the image is not acceptable, touch **More options ▶ Image adjust** to optimize image or touch **Trash** to remove image, then return to step 2 to adjust exposure conditions.

Universal Mode

Mode to image samples containing one or more signal types (chemi, fluor, and/or visible).

Table 3 Illumination sources, excitation filters, and emission filters for the iBright™ CL1500 Instrument.

Illumination	Excitation Filters	Emission Filters
White Epi-LED	455-485 nm	400-700 nm 568-617 nm
ND filter ^[1]	400-700 nm	
Green trans	490-520 nm	

^[1] Neutral density filter decreases the intensity of the white Epi-LED for visible imaging.

Table 4 Illumination sources, excitation filters, and emission filters for the iBright™ FL1500 Instrument.

Illumination	Excitation Filters	Emission Filters
White Epi-LED	455-485 nm	400-700 nm 508-557 nm 568-617 nm 675-720 nm 710-730 nm 800-850 nm
	515-545 nm	
	608-632 nm	
	610-660 nm	
nIR Epi-LED	745-765 nm	
ND filter	400-700 nm	
Green trans	490-520 nm	

1. Touch **Choose Channel** to select from one to four different dyes or stains from the pre-populated list.

Alternatively, to add a dye that is not in the pre-populated list, touch **Custom Dye** at the bottom left hand corner of the screen. Select a custom dye that has previously been entered or enter the information for a new dye: dye name, emission wavelength, excitation wavelength, and false color.

- a. To edit an existing dye, touch the dye name button. Touch **Dye** and change the assigned dye to any other pre-populated dye in the fluorescent dye list. Dyes in the pre-populated dye list are categorized by most recently used, name, and emission color.
- b. Touch **False Color** and change the false color assigned to that dye.

Note: Each dye is assigned a false color to be used in the composite overlay.

- c. To remove this assigned dye from the channel, touch **Remove Dye** at the bottom left hand corner of the screen. Touch **Confirm**.

Channel refers to the type of signal you want to capture and the imaging conditions you want to use.

- **Chemi:** Channel for measuring chemiluminescent signals.
 - **Fluor [Epi-LED]:** Channel for measuring fluorescent signals that are excited using the appropriate Epi-LED/excitation filter combination and the appropriate emission filter.
 - **Fluor [Trans]:** Channel for measuring fluorescent signals that are excited by the green transilluminator and appropriate emission filter.
 - **Nucleic acid:** Channel for measuring fluorescent signals in gel that are excited by the green transilluminator and the 568-617 emission filter (stained DNA and RNA).
 - **Protein:** Channel for measuring fluorescent and visible signals in gel that are excited or illuminated by the green transilluminator and the 568-617 emission filter — stained or labeled protein.
 - **Visible:** Channel for measuring visible signals on an opaque media (nitrocellulose membrane, PVDF membrane, etc.) using the Epi-white LED with neutral density filter and appropriate emission filter.
 - **TPN:** Total protein normalization channel is for normalizing Western blot results to total protein load by lane.
 - **E-Gel™ 96:** Channel for measuring fluorescent signals in high throughput E-Gel™ 96 excited by the green transilluminator and the 568-617 nm emission filter. Choosing this channel also allows for an analysis workflow with deconvolution of image and data.
 - **Colony Count:** Channel for imaging bacterial petri dishes and cell culture plates using different fluorescent filters, trans illumination, and Epi-white LEDs. Choosing this channel also allows for a specialized colony analysis workflow.
2. Touch **Smart Exposure**. Alternatively, for manual exposure, touch **More options ▶ Manual Exposure**. The imager will acquire a series of short exposures for each dye type selected. It will then render preview images and a recommended exposure time for each channel. These are not real images.
 3. (Optional) Touch **More options ▶ Region of Interest**. **Region of Interest** is for directing **Smart Exposure** to a specific region of the sample.
 4. To adjust or set a manual exposure time for any of the channels that you are acquiring, select the channel (dye) that you want to change:
 - Touch **+** or **-** within dial.
 - Finger swipe in the segmented dial.
 - Touch the dial center box to select a preset exposure time or touch **Custom** to manually input a time.

Note: When recommended exposure time is adjusted, the image preview will update in real time.

5. Touch **Capture** to acquire all of the selected channels with the indicated exposure times. Captured image(s) appears in viewport and automatically saves to the gallery.
6. If the image is acceptable, touch **Export**, **Gallery**, or **Analyze**. If the image is not acceptable, touch **More options** ▶ **Image adjust** to optimize image or touch **Trash** to remove the image, then return to steps 3-5 to adjust exposure conditions.

Options available in modes

Each mode offers **More Options** on certain workflow screens to provide detailed camera and image adjustment.

Table 5 Camera

Sub option	Detail
Resolution/Sensitivity	Changes binning setting to increase resolution/decrease sensitivity or decrease resolution/increase sensitivity.
Zoom/Focus	Changes zoom level to increase or decrease image area. Focus is automatically adjusted based on zoom. Auto focus can be used to optimize sharpness. Note: Optical zoom 1X to 2X and digital zoom 2X to 8X.
Sample rotation	Imager mechanically rotates sample $\pm 10^\circ$ depending on orientation. Rotation is automatic, but also allows ability to adjust in More Options ▶ Camera Settings .

Table 6 Routines

Sub option	Detail
Multi-exposure	Series of 5 preset exposure times producing one image for each time period.
Signal accumulation	<p>Series of exposures where user defines the first capture (time), the last capture (cumulative capture time) and the total number of captures.</p> <p style="text-align: center;">Signal Accumulation (Example)</p> <div style="text-align: center;"> </div> <p>Note: Images acquired using Signal Accumulation are not recommended for publication or analysis. This feature is recommended for determining an optimal exposure time.</p>

Table 7 Image adjust

Sub option	Detail
Contrast	Grayscale differentiation between image features. Can use Auto contrast (high, medium, low) or view raw image with no contrast. User can manually adjust Gamma, White, Black using slider bars.
Channels and Layers	Allow user to navigate between individual channels associated with a multi-channel image; pertains only to Chemi Blot, Fluorescent Blot, and Universal modes (e.g., Channels: membrane, Alexa Fluor™ 488, Alexa Fluor™ 555, Composite). Layers allow toggling on and off different channels.
Zoom	For the CL1000, CL1500, FL1000, and FL1500 instruments, zoom increases or decreases the optical (2X) and digital zoom (4X) up to a total 8X zoom.
Invert	Produces a negative or positive (white to black or black to white) of the displayed image.
Saturation	False colors saturated white pixels (65,536) as red to differentiate from non-saturated pixels.

Table 8 Other options

Sub option	Detail
Region of Interest	Specify an area on the image to determine the optimal exposure time for that region using the Smart Exposure tool.
Manual exposure	Manually select an exposure time for sample.
Smart Exposure	Predicts optimal exposure for minimizing pixel saturation and maximizing dynamic range for a specific sample. Renders a preview of how the image will appear at the recommended exposure time.
Smart Range™ HDR	The Smart Range™ HDR algorithm mathematically combines a short and long exposure together into a single image that contains both high and low signal intensities that fall within the linear dynamic range. The new image maintains 16-bit resolution with lower noise and higher linear dynamic range.
Band excision (Nucleic Acid Gels mode)	Opens drawer and turns on green transilluminator to allow for safe band excision on nucleic acid gels.
Membrane overlay (Chemi blots, Fluorescent blots, and Universal Mode)	Allows visible membrane overlay to see any visible prestained markers, which aid in identifying molecular weights for unknown samples.

Table 8 Other options (continued)




Sub option	Detail
<p>2UP view (Fluorescent Blot mode)</p>	<p>After a Smart Exposure or Capture, images are displayed by default in a 2UP view.</p> <ul style="list-style-type: none"> • Top image displays individual channels in grayscale. Switch between individual channels by touching the channel you want to display and edit. After a dye is selected, that channel image is displayed in the top view and becomes editable so the user can increase or decrease the exposure time dial and see in real time the effect on the image preview. • Bottom image displays the color composite with channels displayed as different false colors overlaid on a black background. Individual channels can be toggled on and off in the composite image by touching  at the left of the appropriate channel name. • For the Visible channel color composite, the signal is false-colored overlaid on a white background. Example: This channel is used to display Coomassie-stained protein gels.
<p>1UP view (Fluorescent Blot mode)</p>	<p>After a Smart Exposure or Capture, images can also be displayed in a 1UP view.</p> <p>Image displays the grayscale and false-color composite images by toggling Gray or Color. User can switch between Individual channels by touching the dye name you want to display. Individual channels can be toggled on and off in the composite image by touching  on the left of the appropriate dye name.</p> <p>Note: Functionality for toggling between individual channels in grayscale view or toggling on and off individual channels in the false-color composite view is identical to the 2UP view.</p>
<p>Auto Enhance</p>	<p>Applies rolling-ball correction to the composite image of a multichannel image. It can be toggled on and off and only affects the visible image, not the raw data.</p>
<p>Camera lock settings</p>	<p>Allows user to lock the camera settings to take additional images with the same parameters.</p>

Image Gallery

The Image Gallery is an on-board storage location where your image files are kept for future viewing. In the gallery you can easily retrieve, analyze, and manage your image files. Any image acquired on the imager is automatically saved into a user's individual image gallery. For those not logged on as a user (guest), acquired images are saved to a common gallery that houses all guest image files.

View images

On the **Welcome** screen or mode drop-down menu:

1. Touch **Gallery** or .
2. Toggle between **Thumbnail view** and **List view** to see stored images as thumbnails or a table. Click the column headers to sort the images.
3. Touch an image to select it. Touch again to deselect images. To choose all stored images, touch **Select all**. Unselect using **Deselect all**.
4. To filter images, touch **Filter options** and enter the search criteria. Touch **Done**.
5. To select all of the images captured on a selected date, check the box on the left side for the appropriate date.







Export and analyze images

Export images

Export new images

From the "Welcome" screen:

1. Capture an image. See "Choose imaging mode and start image capture" on page 13.
2. Touch **Export**.
3. (Optional) Touch **Edit** to enter a new file name or comments. Touch **Save**.
4. Choose a **Destination**.
 -  Connect
 -  USB Drive
 -  Network Drive
 -  Network Printer

Note: The instrument requires a path for connection to a network drive or printer. Consult your IT department for questions regarding your network. The following information may be required to map a network drive:

- Drive location: Enter the IP address, server name and file pathway (use forward slashes).
Example: 192.168.1.100/shared/R&D Data/John_Smith/2016
 - Additional information: Enter the domain name, username, and password.
-

5. Choose a file type for export. See "Select an Export file type" on page 26.

Export gallery images

1. In the Gallery, select the desired image/images. See "View images" on page 24 .
2. Touch **Actions**.
3. Touch **Export**.
4. (Optional) Touch **Edit** to enter a new file name or comments. Touch **Save**.

5. Choose a Destination.

-  Connect
-  USB
-  Network Drive
-  Print

Note: The instrument requires a path for connection to a network drive or printer. Consult your IT department for questions regarding your network. The following information may be required to map a network drive:

- Drive location: Enter the IP address, server name and file pathway (use forward slashes).
Example: 192.168.1.100/shared/R&D Data/John_Smith/2016
 - Additional information: Enter the domain name, username, and password.
-

6. Choose a file type for export. See “Select an Export file type” on page 26.

Select an Export file type

When exporting images, select an appropriate file type for your application. From the **Export images** screen:

1. Touch **Choose File Type**.
2. Touch a file type.

File type	Description
Publication - Exports 24-bit color-adjusted images for data sharing and presentation	
TIFF	Best quality
JPEG	Medium quality
PNG	Medium quality
Analysis - Exports 16-bit black and white, non-adjusted images for external image analysis and processing	
TIFF	Best quality
G2i (iBright™ proprietary file type for iBright™ Analysis Software - available on desktop and Connect)	Best quality

Note: **Publication** file types allow you to export composite and all channels as separate image files. Touch the checkbox at the bottom of the screen to choose this option. If unchecked, then only the false-color composite image will be exported.

3. Touch **Next**.

Archive/Backup g2i image files

All iBright™ system users have an individual gallery for viewing and managing their own image files. iBright™ administrators can manage all user galleries simultaneously or individually.

1. Touch  **Gallery**.

Note: You can view the image files in the Gallery as thumbnails or a list. In list view, the user can sort by file name, imaging mode, file size (kB), or acquisition date.

2. Touch **Filter options**.


Note: Galleries can be filtered by keyword, date, imaging mode, and exposure mode (normal and HDR).

3. (Optional) Check the box **Include guest's images** to include the guest gallery.


4. Check the **Images** box for all users to view images from all user galleries.

5. (Optional) Check the **Images for all users** box, touch **Enter ID**, and choose the desired username of the individual gallery that you want to manage.

6. Touch **Done**. In the filtered Gallery, select the desired image file(s) or touch **Select All** to select all image files in the Gallery.

7. Touch **Export** to archive/backup the selected g2i file(s) or touch  to delete.

IMPORTANT! If an excessive number of files are selected, touching **Actions** is not recommended. This may result in long wait times or possibly a system crash. Wait time is directly proportional to the number of and total size of the selected image files.

Note: **Action** in the Gallery is intended for viewing image information, image adjustments, image analysis, file renaming, and file export with no more than 50 image files in the tray. Bulk export and delete should be done in the main Gallery using **Export** or  **Delete**, respectively.

Analyze images

Analyze new images

1. Capture an image. See “Choose imaging mode and start image capture” on page 13.
2. With the desired image open, touch **Analyze**.
The system automatically finds and displays analysis frames, lanes, and bands for the selected image file. Multi-channel images are displayed as individual channels that can be toggled or swiped between using the touchscreen. The system can identify 1-4 different analysis frames.
3. (Optional) Touch the **Sensitivity** dial to raise or lower the band display threshold to increase or decrease the number of bands displayed in each analysis frame.
4. (Optional) Touch **Edit analysis frame** to remove, add or edit existing analysis frames.
5. (Optional for fluorescent images) Touch the dye drop-down menu to select a particular dye or the membrane view.
6. (Optional) Touch **More options** to access image adjustments or to generate a report.

Image adjustments	Functions
Adjust lanes	Add, remove, or edit lanes.
Adjust bands	Add, remove, or edit bands.
Add markers	Identify a molecular weight lane for each analysis frame. A known standard can be used to approximate the molecular weight of unknown bands.
Adjust image	Images are by default displayed auto-contrasted (medium). User can change the auto-contrast to high or low setting or manually set black, white and gamma using the appropriate slider bar.
Normalization	Total Protein Normalization (TPN) uses a total protein label or stain to detect and quantify the total lane protein in a gel or on a membrane. Western blot targets are normalized to total protein to correct for loading and transfer anomalies. Alternatively, a housekeeping protein can be immuno-detected and quantified to approximate protein load for the same purpose.
Generate report	Prepare a pdf report displaying the images and analysis table associated with the image file. Prior to generation of the report, select the desired data to be included in the analysis table and edit the report name if needed. This file can be exported and/or printed.

After each image adjustment:

- a. Touch **Apply** to accept the changes.

- b. (Optional) Touch **Next** to display the analysis table. At the top of the table, touch ∇ to change the table output by:

Output	Units	Description (Lane and Band Table)
Volume	Intensity	Sum of pixel intensities contained in an identified band.
Weight	kDa	Molecular weight estimated for an unknown sample based on the sample's relative mobility in the gel versus the molecular weight and relative mobility of known standards.
Density	Intensity/Area	Average intensity for all pixels contained in an identified band (Volume divided by Area).
Local Background Corrected Volume	Intensity	Volume minus the local background intensity (2-pixel area surrounding the identified band region).
Local Background Corrected Density	Intensity/Area	Local Background Corrected Volume divided by the Area.
Migration Distance	R _f	Distance migrated divided by the gel length.
Band Purity	% percent	Divide the background-corrected band volume (intensity) by the sum of all background-corrected band volume (intensities) in the lane and multiply by 100.
Lane Purity ^[1]	% percent	Divide the background-corrected band volume (intensity) by the background-corrected total lane volume (intensity) and multiply by 100.

^[1] This output is only available in the report.


- c. For TPN-normalized data, the lane and band table has the following additional outputs:

Output	Units	Description (Lane and Band Table)
Total Lane Volume	Intensity	Sum of pixel intensities contained in an identified lane.
Background Volume (RB)	Intensity	Background determined using Rolling Ball Background Correction.
Adjusted Total Lane Volume	Intensity	Total Lane Volume minus background volume.
Normalization Factor	—	Normalization factors are proportional to the total lane protein signal with a reference lane nominally set to 1.000.
Area (band) ^[1]	Pixels	Total number of pixels contained within an identified band.
Normalized Corrected (band) Volume ^[1]	Intensity	Adjusted Total (band) Volume divided by the Normalization Factor of the lane.

^[1] This output is available in the report only.

Analyze gallery images

On the "Welcome" screen or capture screen Mode drop-down menu:

1. Touch **Gallery** or . Select one or multiple images.
2. Touch **Actions**.
3. If more than one image was initially selected, then choose one of the images within the scrollbar. If needed, edit, remove, or delete the selected image.
4. Touch **Analyze**.
The system automatically finds and displays analysis frames, lanes, and bands for the selected image file. Multi-channel images are displayed as individual channels that can be toggled or swiped between using the touchscreen. The system can identify 1-4 different analysis frames.
5. (Optional) Touch the **Sensitivity** dial to raise or lower the band display threshold to increase or decrease the number of bands displayed in each analysis frame.
6. (Optional) Touch **Edit analysis frame** to remove, add or edit existing analysis frames.
7. Touch **Apply** to accept any changes.
8. (Optional) Touch **More options** to adjust lanes or bands, add markers, access image adjustments, perform total protein normalization, or generate a report. For details, see "Analyze new images" on page 28.
For details on TPN normalized data outputs and TPN normalization channel lane outputs, see "Analyze new images" on page 28.
9. (Optional) To generate a report, touch **Generate report**. Select data to include in the report and edit the report name if desired. Press **Enter**. Touch **OK**.
10. Touch **Export**.
11. Select a destination and touch **Done**.
Analysis data is exported as a .csv file along with the report as a .pdf file.

Analyze 96-well E-Gel™ images taken in Universal mode using the E-Gel™ 96 application

1. Open the desired 96-well E-Gel™ from the gallery or viewport. Touch **Analyze**.
The system automatically finds and displays the analysis frame and well template for the selected image file. Lanes and bands are automatically detected after acquiring the frames and wells.
2. Touch **Edit analysis frame** to remove, add, or resize existing analysis frames.
3. Touch **Apply** to save changes. The system will run the well detection algorithm.
4. (Optional) If needed for improved analysis, touch **Reanalyze**.
5. Touch **Next** to analyze sample.

6. Touch **More options** to add, remove, or edit lanes and/or bands.

Image adjustments	Functions
Adjust lanes	Add, remove, or edit lanes.
Adjust bands	Add, remove, or edit bands.
Adjust image	Images are by default displayed auto-contrasted (medium). User can change the auto-contrast to high or low setting or manually set black, white and gamma using the appropriate slider bar.
Deconvolute	Simplify viewing of the image by separating individual samples into individual images.
Generate report	Prepare a pdf report displaying the images and analysis table associated with the image file. Prior to generation of the report, select the desired data to be included in the analysis table. This file can be exported and/or printed.

7. Touch **Apply** to save changes.
8. (Optional) To generate a report, touch **Generate report**. Select data to include in the report and edit the report name if desired. Press **Enter**. Touch **OK**.
9. Touch **Export**.
10. Select a destination and touch **Done**.
Analysis data is exported as a .csv file along with the report as a .pdf file.

Analyze colony images taken in Universal mode using the Colony Count application

1. With the desired image open, touch **Analyze**.
The system automatically finds and displays wells and colonies for the selected image file.
2. Touch **Edit analysis well** to add wells, remove wells, adjust well size, and adjust well shape.
3. Touch **Apply** to save changes.
4. Touch **Adjust gating** to change the colonies displayed based on area, brightness, and circularity.
5. Touch the well to be adjusted, then change the specific parameters.
6. Touch **Done** to save changes.
7. Select the proper channel and background based on the image.
8. Touch **Next** to analyze sample.

9. Touch **More options** to adjust colonies and image.

Image adjustments	Functions
Adjust colonies	Move and resize colonies.
Adjust image	Adjust contrast, channels, and layers.

10. (Optional) To generate a report, touch **Generate report**. Select data to include in the report and edit the report name if desired. Press **Enter**. Touch **OK**.
11. Touch **Export**.
12. Select a destination and touch **Done**.
Analysis data is exported as a .csv file along with the report as a .pdf file.



Maintain the instrument

Settings

Access Settings  from the **Home** screen.

Instrument Settings	
About Instrument	Provides all instrument details including serial number and firmware version.
Date and Time	Select a Time Zone, Date Format, and Time Format for your region.
Sleep Mode	Enable or disable a sleep mode by toggling the button.
Manage Users	Security mode Off = Guest users allowed. On = Login required for all users.
	Manage accounts See "Manage accounts" on page 35.
Instrument Name	Customize the instrument name.
Language	Select a display language (subject to availability).
Reset User Tips	Reset user tips for the logged-in guest.
Reset Factory Defaults	<p>Note: Selecting this reset erases all user data and gallery files.</p> <p>Resets instrument to factory settings and clears out existing user information. The instrument must be powered off, then powered on to apply the change.</p> <p>After resetting factory defaults, perform the new instrument set-up tasks. See Chapter 2, "Start, sign on, and configure the instrument".</p>
Backup and Restore	<p>Backup Systems Settings and restore settings.</p> <p>Note: This does not back up image files.</p> <p>To enable or disable automatic file backup, see "Network Drive" in the "Network Configuration" section (below).</p>
File Naming Convention	
User sets preference for file naming by selecting up to 3 of the following aspects: Mode , Binning , Custom Field User , Exposure Time , and Sequential Number . Date and Time are mandatory aspects. The aspects can be ordered by choice.	

Error Log
Displays instrument error history with sortable columns for Date , User , and Error Description . The Error Log is exportable.

Event Log
Displays instrument error history with sortable columns for Date , User , Event Type , and Cloud Status .

Software Update
Updates the instrument software package by downloading via USB dongle or connection to the cloud. Note: Cloud option requires a connection to your Connect account.


Network Configuration	
Wired	Searches for an ethernet connection.
Wireless	Searches for a wireless connection. Requires a USB wireless adapter (High-Power USB Wi-Fi Module, Cat. No. A26774).
Network Drive	<ul style="list-style-type: none"> • Enter a Drive location for connection. If needed, enter a Domain name, Username, and Password. • Toggle the button to enable or disable automatic g2i file export and backup upon logout.
Printer Configuration	Enter an IP address and choose between US Letter or A4 paper size.
Cloud Region	Select the cloud region associated with your location: China if you are located in China and Other if you are located anywhere other than China.

Capture History
Displays image capture history with sortable columns for Date , User , Capture Mode , Image Name , and Capture Status .


Service Tools	
Instrument Diagnostic	Runs diagnostic testing on the instrument hardware. Provides an end report showing the functional status of hardware items. Note: Running diagnostics is not required. Contact a service technician for any issues.
Ship Prep Mode	Allows removal of the turntable for shipping. Note: Instrument can only be powered off once turntable is removed.
Pixel Mapping	Pixel mapping is used to identify camera-specific pixel intensity differences that occur in long exposures and applies a correction for any pixel that deviates from the normal dark value.

Manage accounts

Manage accounts allows for editing of user types and PINs and deleting profiles.

1. Touch  and log in to your user account. Once you log in, you are automatically taken to the Welcome screen.

Note: You must be an Admin logged into your profile to manage accounts.


2. Touch , then click **All Accounts**.
 - All current user profiles will display.
 - Columns for **Users**, **Created (date)**, **User Type**, and **Account Type** are sortable.
3. Select the user whose account you want to manage.
4. Select **Manage** on the bottom left side of the screen. Change a standard user to an Admin user by sliding the button to **Admin**.

Note: You cannot change an Admin user to a standard user.

5. **Delete PIN** — If a standard user forgets their PIN number, then you can delete their PIN so they are prompted to create a new PIN the next time they log on.
6. **Delete Account** — If you delete a user's account, you will also delete all of the image files in their gallery.

Run instrument diagnostics

Instrument functions can be checked by a user to confirm proper hardware operating conditions.

1. Touch .
2. Touch **Service Tools**.
3. Touch **Instrument diagnostic**.
4. Touch **Start test**.

The instrument will check the instrument hardware for correct operation and settings. The test will take ~10 minutes to complete. The results of the test will display once complete. Detailed results can be exported for further review.

Note: Ensure the drawer is not blocked. The drawer will open as part of the testing. (Not applicable to iBright™ CL750 model.)

Clean the instrument

Repeated instrument use can cause spots and smudges on the sample stage, which can then decrease image quality. Clean the sample stage as needed.

Materials required

- Safety glasses
- Powder-free gloves
- Tissue, lint-free
- Deionized water
- Ethanol, 70% solution

Note: Avoid the use of detergents.

1. Open the drawer to expose the sample stage.
2. Lightly spray the glass surface with deionized water or a 70% ethanol solution.
3. Wipe the surface with a lint-free tissue until sufficiently clean.
4. Close the drawer and operate the instrument as normal.

Replace the instrument fuses

Materials required

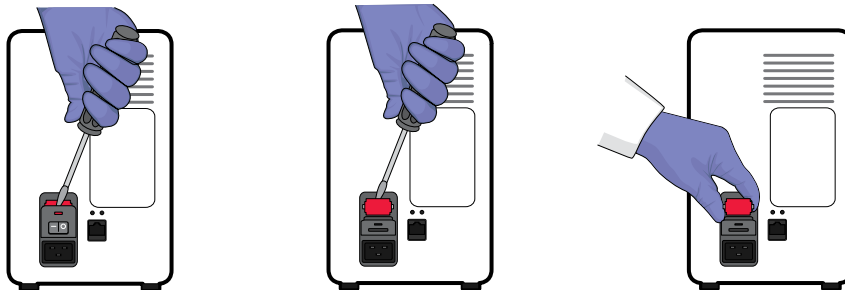
- Fuses, 1.6A, Time-Lag T, 250VAC, 5 × 20-mm (2)
- Safety glasses
- Powder-free gloves
- Screwdriver, flathead

Replace the fuses

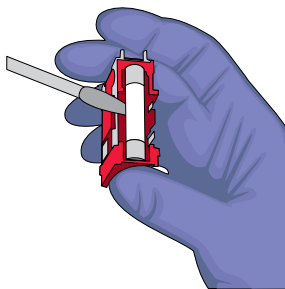


CAUTION! FIRE HAZARD. For continued protection against the risk of fire, replace fuses only with listed and certified fuses of the same type and rating as those currently in the iBright™ imager.

1. Power off, then unplug the iBright™ imager.
2. Using a flat-head screwdriver, pry open the fuse door, and remove the fuse holder.



3. Remove each fuse from its fuse holder and inspect it for damage. Carbon typically coats the inside of failed fuses.



4. Replace each failed fuse with a 1.6A, Time-Lag T, 250VAC, 5 × 20-mm Fuse.

Note: The voltage and amperage ratings are on the fuse holder.

5. Install the fuse holder.
6. Plug in, then power on the iBright™ imager. The installation is successful if the instrument powers on.

Note: Fuse failure can result from fluctuations in the supplied power to the iBright™ imager. To prevent further failures, consider installing an electrical protective device, such as a UPS or surge protector.

Move the instrument




CAUTION! PHYSICAL INJURY HAZARD. Do not attempt to lift the instrument or any other heavy objects unless you have received related training. Incorrect lifting can cause painful and sometimes permanent back injury. Use proper lifting techniques when lifting or moving the instrument. At least two people are required to lift it.

IMPORTANT! Moving your instrument can create subtle changes in the alignment of the instrument optics. Recalibrate the instrument if necessary.

- Ensure that the surface on which you place the instrument can support at least 50 kg (110 lbs).
- Ensure that the path to transport the instrument is clear of obstructions.
- At least two people are needed to lift and carry the instrument.
- Keep your spine in a good neutral position.
- Bend at the knees and lift with your legs.
- Do not lift an object and twist your torso at the same time.
- Coordinate your intentions with your assistant before lifting and carrying.

Prepare for shipping

1. Touch .
2. Touch **Service Tools**.
3. Touch **Ship prep mode**.
Drawer opens.
4. Remove turntable.
5. Touch **Next**.
6. Power off the instrument.



Invitrogen™ iBright™ SAE Software Solution for 21 CFR Part 11 Support

Security Audit e-Signature (SAE) for the iBright™ Imager

IMPORTANT! 21 CFR part 11 is a regulation that describes the criteria for acceptance by the U.S. Food and Drug Administration (FDA) for electronic records and electronic signatures. Part 11 is composed of procedural and technical requirements. Procedural requirements are the standard operating procedures instituted by the end user, and technical requirements are the functional characteristics of the compliance management software used.

This section is intended to provide instructions for using Security, Audit, and E-Signature (SAE) on board iBright™ Imaging Systems. Security, Audit, e-Signature (SAE) software provides the tools necessary for supporting 21 CFR Part 11 technical compliance including:

- Creating and maintaining user accounts
- Managing and enforcing password policies of all accounts
- Assigning, managing and enforcing access rights to all accounts
- Documenting and maintaining audit and e-Signature histories
- Permitting e-Signature approval for data verification

To implement the iBright™ SAE Software Solution for 21 CFR Part 11 support on iBright™ Imaging Systems, you need the following components installed, activated, and communicating:

- **SAE Administrator Console** – With the iBright™ Imager Analysis Application Profile, used to configure the SAE settings for iBright™ Imaging Systems
- **iBright™ SAE License** – Used to activate the SAE settings for the iBright™ instrument and desktop analysis software
- **iBright™ Instrument – SAE Mode** – Firmware connected with the SAE Administrator Console
- **iBright™ Analysis Software – Secure** – Desktop analysis software connected with the SAE Administrator Console


Note: For details on accessing iBright™ SAE Software Solution downloads and resources, see the "Related documentation and references" section.

Activate your license key

1. Log in as Local Admin on the instrument.

IMPORTANT! Only the Local Administrator (Local Admin) has the ability to activate the license key on the instrument. The license key only needs to be activated one time.

Note: Running an instrument qualification using Installation Qualification/Operational Qualification/Performance Qualification (IQ/OQ/PQ) is recommended before SAE mode is enabled on the iBright™ imager for the first time.

2. Touch  **Settings** on the right of the top toolbar.

3. Touch **SAE**.

4. Touch **SAE License**.

5. Enter the SAE License Activation Code and touch **Next**.

Note: If the license validation is successful, "Registration complete" appears.

Note: The license activation code is case sensitive.

6. Press **Done**. The activation code will be saved.


7. Enter the IP address and port of the SAE console location, then touch **Next**.

Note: IP address and port location (e.g., 10.X.X.X:8201/admin-console/index.html) are available in the address bar of the SAE Admin Console.

Enable SAE mode

1. Log in as Local Admin on the iBright™ instrument.

IMPORTANT! Only the Local Administrator (Local Admin) has the ability to enable SAE mode on the iBright™ imager.

2. Touch  **Settings** located on the right side of the top tool bar.

3. Touch **SAE**.

4. Touch **SAE mode**.


5. Move the slider to **Enable**, then touch **Next**.

6. If not already done during the license activation, enter the IP address and port of the SAE console location and then touch **Next**.

Note: IP address and port location (e.g., 10.X.X.X:8201/admin-console/index.html) are available in the address bar of the SAE Admin Console.

7. The system will require the SAE Administrator to enter their credentials to complete the process. Enter credentials and touch **Next** to complete the process.

Disable SAE mode

1. Log in as Local Admin or SAE Console Admin on the iBright™ instrument.
2. Touch  **Settings** located on the right side of the top tool bar.
3. Touch **SAE**.
4. Touch **SAE mode**.
5. Move the slider to **Disable**, then touch **Next**.
6. At this point the SAE Admin may need to enter their credentials and touch **Disable** to continue or **Cancel** to leave the SAE mode enabled.
7. If user touches **Disable**, then the following warning will pop up:

Note: SAE users will not be able to access their current images when SAE is disabled. You can backup or export those files before disabling SAE. Image files remain on the system and can be accessed when SAE mode is enabled again.

Note: It is recommended an SAE Admin perform this action. If a Local Admin performs this action then it will not be recorded in the audit history on the SAE Admin Console.

Sign in to the instrument in SAE mode

1. Enter your SAE username and password in the SAE login window.

Note: Your username and password must be created by the SAE Admin on the SAE Console prior to logging in.

2. Touch **Sign In**. After signing in, you may be prompted to change your SAE password.

IMPORTANT! The following symbols cannot be used in your password: + & % \

Note: The following characters can be used in addition to letters and numbers for passwords: - _ () [] {} / ~ ! @ # \$ * ^ = . , ' " ; : | < > ?

Change your password in SAE mode

1. Touch **User Profile** on the left side of the top toolbar.
2. Touch **Edit** and enter your old password, your new password, and then confirm your new password.

IMPORTANT! The following symbols cannot be used in your password: + & % \

Note: The following characters can be used in addition to letters and numbers for passwords: - _ () [] {} / ~ ! @ # \$ * ^ = . , ' " : ; | < > ?

Note: Your new password must meet the complexity requirements set by the SAE Console Admin.

Note: If you exceed the maximum number of sign-in attempts, you will be temporarily locked out of your account. See the SAE Console Admin.

Note: You may have limited access to iBright™ instrument functions based on your user role set by the SAE Console Admin. in the SAE Console.

SAE user roles

Role	Description
SAE Console Administrator (SAE Console Admin)	A default system role with full access privileges to the SAE Console. The SAE Console Admin creates new SAE user accounts, defines what privileges are granted to each SAE user role, and assigns user roles to all other SAE users.
iBright™ Administrator	A default SAE role with privileges to use and manage the iBright™ instrument and iBright™ Analysis Software – Secure.
iBright™ Scientist	A default SAE role with privileges to use the iBright™ instrument and iBright™ Analysis Software – Secure.

SAE privileges for default user roles

Category	Function	Default Roles		
		SAE Console Admin	iBright™ Admin	iBright™ Scientist
Image Management	Acquire, Analyze, Adjust, Export, and Delete	Yes	Yes	Yes
Security Configuration	Sign g2i File, Unlock g2i File	Yes	Yes	Yes
	Enable SAE on instrument [1]	Yes	No	No
	Disable SAE on instrument	Yes	Yes	No
Instrument Configuration	Administer Instrument, Update Instrument System Configuration	Yes	Yes	No
Service Tools	Instrument Diagnostics, Pixel Mapping, and Export Workspace	Yes	Yes	No

[1]

Enabling SAE on board iBright™ instruments requires the Local Admin (initiate) and the SAE Console Admin (approve).

Non-SAE user roles

Role	Description
Local Administrator (Admin)	Local Admin has full administrator access to the iBright™ instrument when SAE mode is disabled. When SAE mode is enabled, this non-SAE role is only allowed to enable and disable SAE mode and to change their own password.
Field Service Representative (FSE)	FSE must log into the iBright™ instrument when SAE mode is disabled to have full access to the Service Tools.

Acquire images in SAE mode

Image acquisition in SAE mode is the same as in non-SAE mode.

For iBright™ CL750 instruments, see the "Choose imaging mode and start image capture" section of the iBright™ CL750 Imaging System User Guide (MAN0018652).


For iBright™ CL1500/FL1500 instruments, see the "Choose imaging mode and start image capture" section of the iBright™ CL1500/FL1500 Imaging Systems User Guide (MAN0018592).

Note: In SAE mode, data is committed to the audit log as v1 at image acquisition.


Use the Gallery in SAE mode

File status

File status is depicted on the thumbnails in the Gallery using the following symbols:

- X – Uncommitted: Image file(s) with uncommitted changes
- ✓ – Committed: Image file(s) with committed changes (saved)
-  – Signed Data (locked): Image file(s) with committed changes and required e-Signatures

Filter Gallery by file status

1. To filter or sort image files in the iBright™ Gallery by file status, touch **Filter options**.
2. Check the appropriate boxes.
 - X – Uncommitted: Image file(s) with uncommitted changes
 - ✓ – Committed: Image file(s) with committed changes (saved)
 -  – Signed Data (locked): Image file(s) with committed changes and required e-Signatures

Archive/Backup g2i image files in SAE mode

All SAE users have an individual gallery for viewing and managing their own secured image files. SAE users with privileges to configure the iBright™ instrument can manage all user galleries simultaneously or individually.



1. Touch  **Gallery**.

Note: You can view the image files in the Gallery as thumbnails or a list. In list view, the user can sort by file name, imaging mode, file size (kB), or acquisition date.



2. Touch **Filter options**.

Note: Galleries can be filtered by acquisition date and imaging mode.

3. (Optional) Check the box **Include guest's images** to include the guest gallery.

4. Check the **Images** box for all users to view images from all user galleries.
5. (Optional) Check the **Images for all users** box, touch **Enter ID**, and choose the desired username of the individual gallery that you want to manage.
6. Touch **Done**. In the filtered Gallery, select the desired image file(s) or touch **Select All** to select all image files in the Gallery.
7. Touch  **Export** to archive/backup the selected g2i file(s) or touch  to delete.

IMPORTANT! If an excessive number of files are selected, touching **Actions** is not recommended. This may result in long wait times or possibly a system crash. Wait time is directly proportional to the number of and total size of the selected image files.

Note: Action in the Gallery is intended for viewing image information, image adjustments, image analysis, file renaming, and file export with no more than 50 image files in the tray. Bulk export and delete should be done in the main Gallery using  **Export** or  **Delete**, respectively.

Commit changes

In SAE mode, user changes (acquisition, analysis, adjustments, editing, renaming, and export) are recorded in the audit history for each image file. After image acquisition, the file is automatically committed, and the audit history is initiated. All additional changes are maintained as uncommitted changes and officially become part of the audit history when the user commits the change(s).

Note: The **Commit Changes** button is only available if the file(s) you are reviewing have uncommitted changes (X).

1. To commit changes for an image file(s) in the gallery, select the desired image file(s) and touch **Actions**.
2. Touch **Commit Changes**.

Note: The **Commit Changes** button is also available in the **More Options** menu in the acquisition, analyze, and adjust workflows.

3. Touch **Reason for Commit** and from the dropdown menu choose the appropriate reason.

Note: **Reason for Commit** will appear based on settings in the SAE Admin Console. If Required or Optional, users will have the ability to indicate the reason. If set to Silent, changes will be committed to the audit log, but no reason will be prompted.

4. Touch **Save**. If successful, a window will show **Changes committed**.

Note: Once complete, the file status changes from **Uncommitted** (X) to **Committed** (✓) and a new audit version is added to the audit history. The **Commit** button is dynamic and once changes are committed, it changes to **Sign Data** (see “Sign data” on page 46) to allow the user to perform that action.

Sign data

In SAE mode, committed image files can be e-signed by one or more users to verify data integrity. The SAE Console Admin can dictate what actions require signatures and the number of user roles that are required to e-sign for each action. Once an image file has been e-signed by all required signees, it becomes a signed file and is locked from future edits.

1. To sign a committed file, touch **Sign Data**.

Note: The **Sign Data** button is also available under **More options** in the acquisition, analyze, and adjust workflows.

2. Select the **Meaning** for signing the file from the drop-down menu.

IMPORTANT! If more than one e-Signature is required on the imager, all signatures must be collected before exiting the e-Signature workflow or the process must be repeated (i.e., all signers must be present).

3. Touch **Next**.

Note: A series of icons will be displayed for all required signatures. Role requirements will be listed below.

4. Select any **Press to Sign** button.

5. Enter your **Username** and **Password**.

6. Touch **Done**. If credentials are accepted the user returns to the **Sign Data** screen.

7. Repeat steps 4-6 until all required signatures are collected, the file status changes from Committed (✓) to Signed/Locked (🔒), and they are recorded in the e-Signature history with an active status.

Unlock files

In SAE mode, signed files are locked and must be unlocked by a user with unlock privileges before any further changes can be made to the file. When a file is unlocked, the file's e-Signature status changes to inactive.

1. In the Gallery, select the desired image file(s), touch **Actions** and then touch **Unlock**.


Note: The user will be warned that they are about to inactivate the signature.

Note: **Unlock** is also available under **More Options** in the acquisition, analyze, and adjust workflows.

2. The user must comment about why the file is being unlocked, then touch **Unlock**.

- A message informs you when the file has been successfully unlocked.
- e-Signature becomes inactive in the e-Signature history status.

Generate an audit history

1. Touch **Audit history** to review the audit history and the change table for the selected image file.
 - The audit history contains the following information: Version, Date & Time, Username, Role, Action, Reason and Comment.
 - **Audit history** is also available in the Adjust and Analyze workflows under **More options**.
2. Select one row in the audit history to view the image and change details for the selected and previous version or select any two rows to view the images and change details for the two selected versions.
3. Touch **Open** to view and compare the images from selected versions.
4. (Optional) Change the channel from the drop-down menu on top of the image to view a different channel image in a multichannel image.
5. (Optional) Select/deselect layer options on the top of image to show/hide the layers on image.
6. Touch **Audit Report** to generate the Audit History Report as a pdf file. The report contains an audit history table along with the image comparisons of the selected versions.
7. Touch  **Export** to export the Audit History Report to a USB drive, network drive, or network printer.


View the e-Signature history

Touch **e-Signature history** to review the e-Signature history.

Note: The **e-Signature History** contains the following information: Commit version, Signed Date (and time), Signed By, Signed Role, Meaning, and Status.

Note: **e-Signature History** is also available in the Adjust and Analyze workflows under **More options**.

Export image files

Touch  **Export** to export an image as a g2i file type.

Note: Other file extensions are available, but only g2i files can be securely exported from the iBright™ instrument and opened in the iBright™ Analysis Software – Secure.

Note: Uncommitted files cannot be exported.

Note: g2i files can be exported as signed or unsigned files based on the file status.

Generate and export Analysis Report

Touch **Generate Report** under the **More options** menu to generate a PDF report with image information, channel images, and densitometry data.

Note: When generating reports from an unsigned g2i file, you will be prompted to cancel the report generation and sign the file to generate the signed report; or proceed to generate unsigned report.

Note: Reports generated from signed g2i files can be exported only as signed reports.



Instrument specifications

Instrument dimensions and specifications

Specification	iBright™ Imager
Dimensions (L × W × H)	68.0 × 38.2 × 59.8 cm
Weight	52 kg (115 lbs)
Instrument clearance (Back)	15.25 cm (6 in.)

Electrical requirements



WARNING! For safety, the power outlet used for powering the instrument must be accessible at all times. In case of emergency, you must be able to immediately disconnect the main power supply to the instrument. Allow adequate space between the wall and the equipment so the power cord can be disconnected in case of emergency.

- Electric receptacle with grounding capability
- Maximum power dissipation: ~960 W (not including computer and monitor)
- Mains AC line voltage tolerances must be up to ±10 percent of nominal voltage

Device	Rated voltage	Circuit required	Rated frequency	Rated power
iBright™ Imager	100–240 ±10% VAC ^[1]	10 A	50/60 Hz	410 W

^[1] If the supplied power fluctuates beyond the rated voltage, a power line regulator may be required. High or low voltages can adversely affect the electronic components of the instrument.

Environmental requirements

Condition	Acceptable range
Installation site	Indoor use only
Electromagnetic interference	Do not use this device in close proximity to sources of strong electromagnetic radiation (for example, unshielded intentional RF sources). Strong electromagnetic radiation may interfere with the proper operation of the device.

(continued)

Condition	Acceptable range
Altitude	Between sea level and 2000 m (6500 ft.) above sea level
Operating conditions	<ul style="list-style-type: none"> Humidity: 15–80% relative humidity (noncondensing) Temperature: 15 to 30°C (59 to 86°F) <p>Note: For optimal performance, avoid rapid or extreme fluctuations in room temperature.</p>
Storage and transport conditions	<ul style="list-style-type: none"> Humidity: 20–80% relative humidity (noncondensing) Temperature: –30 to 60°C (–22 to 140°F)
Thermal output	During operation, the net thermal output, based on the actual current draw of the instrument, is expected to be approximately 960 W (3275 Btu/h).
Vibration	Ensure that the instrument is not adjacent to strong vibration sources, such as a centrifuge, pump, or compressor. Excessive vibration will affect instrument performance.
Pollution degree	<p>The instrument has a Pollution Degree rating of II. The instrument may only be installed in an environment that has nonconductive pollutants such as dust particles or wood chips. Typical environments with a Pollution Degree II rating are laboratories and sales and commercial areas.</p> <p>The noise output of the instrument is ≤ 55 dB(A) when running.</p>
Other conditions	<p>Ensure the instrument is located away from any vents that could expel particulate material onto the instrument components.</p> <p>Avoid placing the instrument and computer adjacent to heaters, cooling ducts, or in direct sunlight.</p>

Network requirements

The instrument:

- Is factory-configured for IPv4 TCP/IP communication and includes an Ethernet adapter (100/1,000 Mbps) with an RJ45-type connector for integrating the device into a local area network (LAN).
- Can alternatively be configured for wireless networking (High Power USB WiFi Module required, sold separately as an optional accessory).

The instrument can be configured for *either* wired or wireless networking, not both.

If a Thermo Fisher Scientific service representative is to install the instrument:

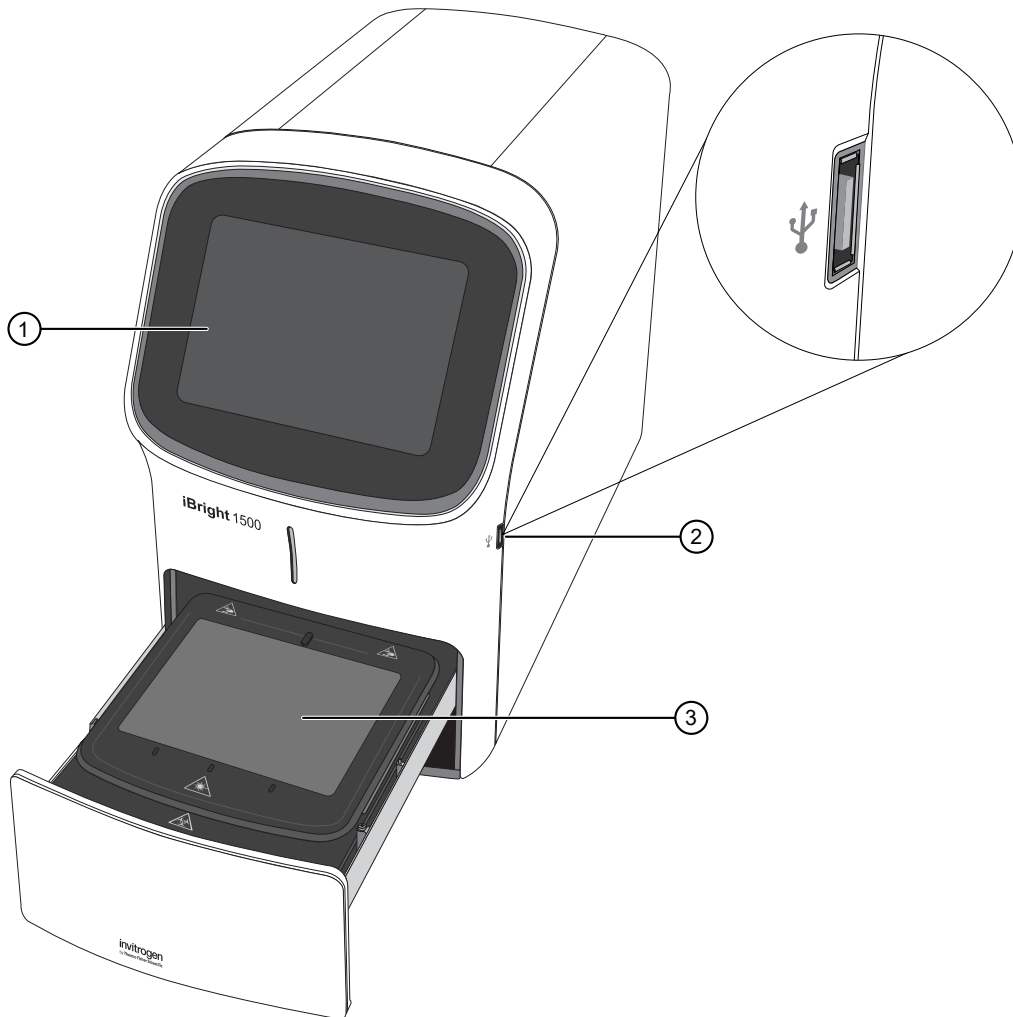
- If the instrument will be connected to a LAN, an active, tested network jack must be in place before the scheduled installation date.
- A representative from your information technologies department must be available during the installation to help connect the instrument to your network.

Required materials to network the instrument:

- Wired—Ethernet cable of sufficient length with RJ45 connectors
 - CAT5 cable for a 100 Mbps network connection
- or–
- Wireless—High Power USB WiFi Module (Cat. No. A26774, sold separately)

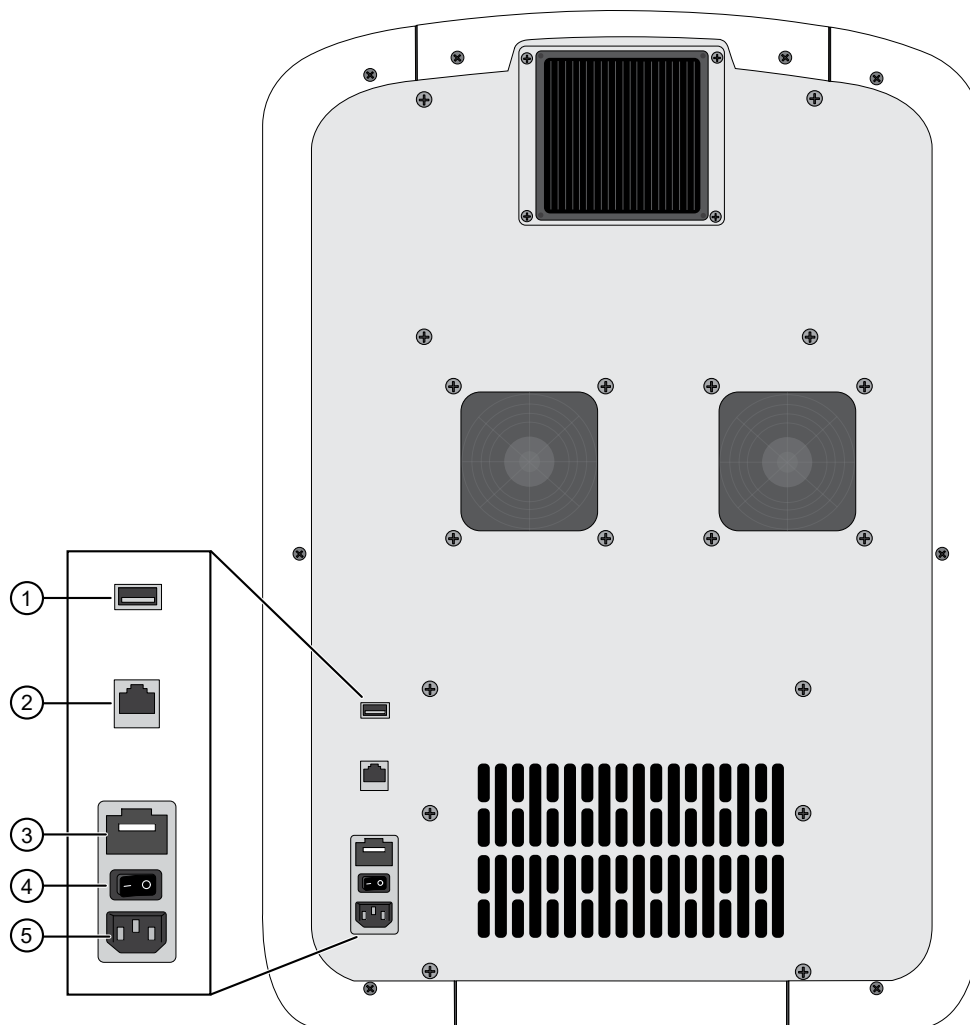
Instrument and computer connections

iBright™ Imager front view



- ① Touchscreen
- ② USB port
- ③ Sample stage/turntable

iBright™ Imager rear view



- ① USB port
- ② Ethernet port — RJ45 port for 10/100 Mbps ethernet communication with the instrument
- ③ Fuse cover
- ④ Power switch
- ⑤ Power port — 100-240 VAC



WARNING! GENERAL SAFETY. Using this product in a manner not specified in the user documentation may result in personal injury or damage to the instrument or device. Ensure that anyone using this product has received instructions in general safety practices for laboratories and the safety information provided in this document.

- Before using an instrument or device, read and understand the safety information provided in the user documentation provided by the manufacturer of the instrument or device.
- Before handling chemicals, read and understand all applicable Safety Data Sheets (SDSs) and use appropriate personal protective equipment (gloves, gowns, eye protection, and so on). To obtain SDSs, visit thermofisher.com/support.




Symbols on this instrument

Symbols may be found on the instrument to warn against potential hazards or convey important safety information. In this document, the hazard symbol is used along with one of the following user attention words:

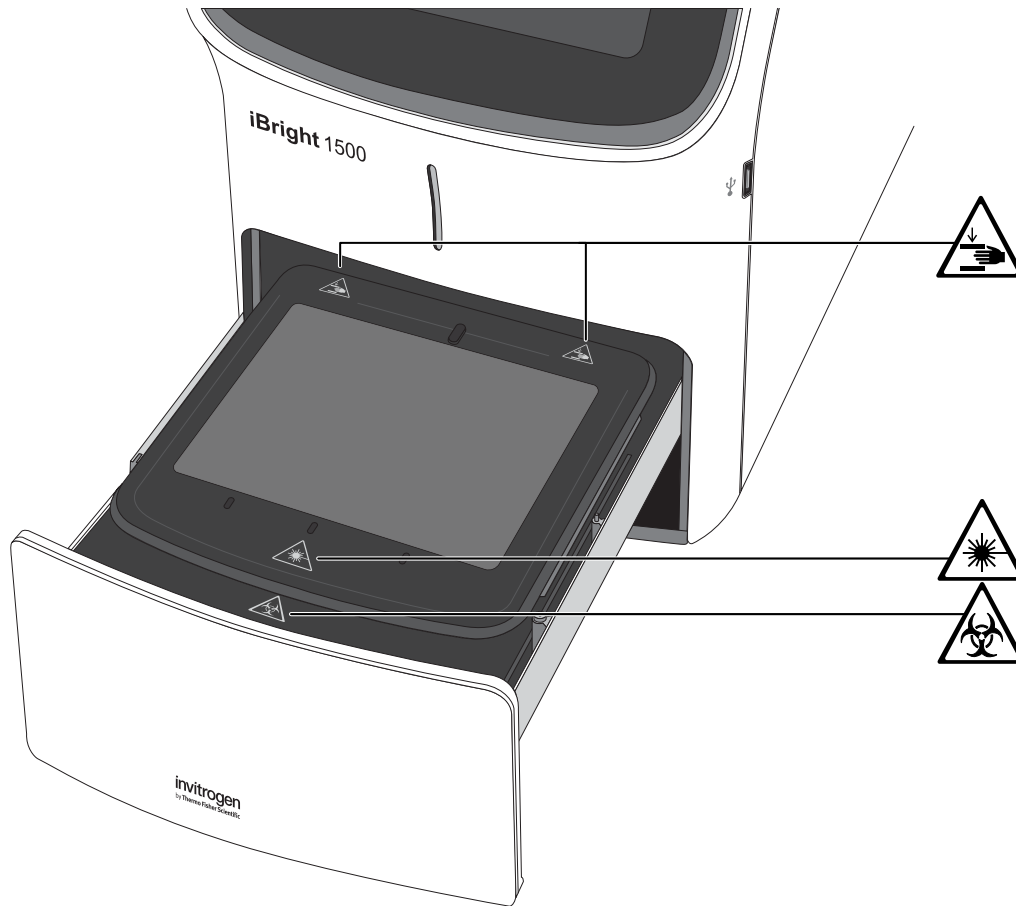
- **CAUTION!** – Indicates a potentially hazardous situation that, if not avoided, may result in minor or moderate injury. It may also be used to alert against unsafe practices.
- **WARNING!** – Indicates a potentially hazardous situation that, if not avoided, could result in death or serious injury.
- **DANGER!** – Indicates an imminently hazardous situation that, if not avoided, will result in death or serious injury.

Symbol	English	Français
	Caution, risk of danger Consult the manual for further safety information.	Attention, risque de danger Consulter le manuel pour d'autres renseignements de sécurité.
	Protective conductor terminal (main ground)	Borne de conducteur de protection (mise à la terre principale)

(continued)

Symbol	English	Français
	<p>Do not dispose of this product in unsorted municipal waste</p> <p> CAUTION! To minimize negative environmental impact from disposal of electronic waste, do not dispose of electronic waste in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provision and contact customer service for information about responsible disposal options.</p>	<p>Ne pas éliminer ce produit avec les déchets usuels non soumis au tri sélectif.</p> <p> MISE EN GARDE ! Pour minimiser les conséquences négatives sur l'environnement à la suite de l'élimination de déchets électroniques, ne pas éliminer ce déchet électronique avec les déchets usuels non soumis au tri sélectif. Se conformer aux ordonnances locales sur les déchets municipaux pour les dispositions d'élimination et communiquer avec le service à la clientèle pour des renseignements sur les options d'élimination responsable.</p>

Location of safety labels on the instrument



Safety information for instruments not manufactured by Thermo Fisher Scientific

Some of the accessories provided as part of the instrument system are not designed or built by Thermo Fisher Scientific. Consult the manufacturer's documentation for the information needed for the safe use of these products.

Instrument safety

General



CAUTION! Do not remove instrument protective covers. If you remove the protective instrument panels or disable interlock devices, you may be exposed to serious hazards including, but not limited to, severe electrical shock, laser exposure, crushing, or chemical exposure.

Physical injury



CAUTION! Moving Parts. Moving parts can crush, pinch and cut. Keep hands clear of moving parts while operating the instrument. Disconnect power before servicing.

LED (Light-Emitting Diode)



CAUTION! LED (light-emitting diode) HAZARD. Removing the protective covers and (when applicable) defeating the interlock(s) may result in exposure to the internal LED. LEDs can burn the retina, causing permanent blind spots. To ensure safe LED operation:

- Never look directly into the light beam.
- Wear proper eye protection and post a warning sign at the entrance to the laboratory if the LED protection is defeated for servicing
- Remove jewelry and other items that can reflect a light beam into your eyes or those of others
- Do not remove safety labels, instrument protective panels, or defeat safety interlocks.
- The system must be installed and maintained by a Thermo Fisher Scientific Technical Representative.

Electrical safety



WARNING! Ensure appropriate electrical supply. For safe operation of the instrument:

- Plug the system into a properly grounded receptacle with adequate current capacity.
- Ensure the electrical supply is of suitable voltage.
- Never operate the instrument with the ground disconnected. Grounding continuity is required for safe operation of the instrument.



AVERTISSEMENT ! Veiller à utiliser une alimentation électrique appropriée. Pour garantir le fonctionnement de l'instrument en toute sécurité :

- Brancher le système sur une prise électrique correctement mise à la terre et de puissance adéquate.
- S'assurer que la tension électrique est convenable.
- Ne jamais utiliser l'instrument alors que le dispositif de mise à la terre est déconnecté. La continuité de la mise à la terre est impérative pour le fonctionnement de l'instrument en toute sécurité.



WARNING! Power Supply Line Cords. Use properly configured and approved line cords for the power supply in your facility.



AVERTISSEMENT ! Cordons d'alimentation électrique. Utiliser des cordons d'alimentation adaptés et approuvés pour raccorder l'instrument au circuit électrique du site.



WARNING! Disconnecting Power. To fully disconnect power either detach or unplug the power cord, positioning the instrument such that the power cord is accessible.



AVERTISSEMENT ! Déconnecter l'alimentation. Pour déconnecter entièrement l'alimentation, détacher ou débrancher le cordon d'alimentation. Placer l'instrument de manière à ce que le cordon d'alimentation soit accessible.

Cleaning and decontamination



CAUTION! Cleaning and Decontamination. Use only the cleaning and decontamination methods specified in the manufacturer's user documentation. It is the responsibility of the operator (or other responsible person) to ensure the following requirements are met:

- No decontamination or cleaning agents are used that could cause a HAZARD as a result of a reaction with parts of the equipment or with material contained in the equipment.
- The instrument is properly decontaminated a) if hazardous material is spilled onto or into the equipment, and/or b) prior to having the instrument serviced at your facility or sending the instrument for repair, maintenance, trade-in, disposal, or termination of a loan (decontamination forms may be requested from customer service).
- Before using any cleaning or decontamination methods (except those recommended by the manufacturer), users should confirm with the manufacturer that the proposed method will not damage the equipment.



MISE EN GARDE ! Nettoyage et décontamination. Utiliser uniquement les méthodes de nettoyage et de décontamination indiquées dans la documentation du fabricant destinée aux utilisateurs. L'opérateur (ou toute autre personne responsable) est tenu d'assurer le respect des exigences suivantes:

- Ne pas utiliser d'agents de nettoyage ou de décontamination susceptibles de réagir avec certaines parties de l'appareil ou avec les matières qu'il contient et de constituer, de ce fait, un DANGER.
- L'instrument doit être correctement décontaminé a) si des substances dangereuses sont renversées sur ou à l'intérieur de l'équipement, et/ou b) avant de le faire réviser sur site ou de l'envoyer à des fins de réparation, de maintenance, de revente, d'élimination ou à l'expiration d'une période de prêt (des informations sur les formes de décontamination peuvent être demandées auprès du Service clientèle).
- Avant d'utiliser une méthode de nettoyage ou de décontamination (autre que celles recommandées par le fabricant), les utilisateurs doivent vérifier auprès de celui-ci qu'elle ne risque pas d'endommager l'appareil.

Instrument component and accessory disposal

To minimize negative environmental impact from disposal of electronic waste, do not dispose of electronic waste in unsorted municipal waste. Follow local municipal waste ordinances for proper disposal provision and contact customer service for information about responsible disposal options.

Safety and electromagnetic compatibility (EMC) standards

The instrument design and manufacture complies with the following standards and requirements for safety and electromagnetic compatibility.

Safety compliance

Reference	Description
EU Directive 2014/35/EU	European Union “Low Voltage Directive”
IEC 61010-1 EN 61010-1 UL 61010-1 CAN/CSA C22.2 No. 61010-1	<i>Safety requirements for electrical equipment for measurement, control, and laboratory use – Part 1: General requirements</i>
IEC 61010-2-081 EN 61010-2-081	<i>Safety requirements for electrical equipment for measurement, control and laboratory use – Part 2-081: Particular requirements for automatic and semi-automatic laboratory equipment for analysis and other purposes</i>
IEC 62471-1:2007 EN 62471-1:2007	Photobiological safety of lamps and lamp systems

EMC

Reference	Description
Directive 2014/30/EU	European Union “EMC Directive”
IEC 61326-1 EN 61326-1	<i>Electrical Equipment for Measurement, Control and Laboratory Use – EMC Requirements – Part 1: General Requirements</i>
AS/NZS 2064	<i>Limits and Methods of Measurement of Electromagnetic Disturbance Characteristics of Industrial, Scientific, and Medical (ISM) Radiofrequency Equipment</i>

(continued)

Reference	Description
ICES-001, Issue 4	Industrial, Scientific and Medical (ISM) Radio Frequency Generators
FCC Part 15 Subpart B (47 CFR)	U.S. Standard Radio Frequency Devices

This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to part 15 of the FCC Rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

Environmental design

Reference	Description
EU Directive 2012/19/EU	European Union “WEEE Directive” – Waste electrical and electronic equipment
EU Directive 2011/65/EU	European Union “RoHS Directive” – Restriction of hazardous substances in electrical and electronic equipment
EU Directive 2006/66/EC	European Union “Battery Directive”
GB/T 26572-2011	Requirements of concentration limits for certain restricted substances in electrical and electronic products
SJ/T 11364-2014	Marking for the restricted use of hazardous substances in electronic and electrical products

产品中有害物质的名称及含量 China EEP Hazardous Substances Information						
部件名称 Component Name	iBright™ CL1000 Instrument					
	iBright™ FL1000 Instrument					
	有害物质 Hazardous Substances					
	铅 (Pb)	汞 (Hg)	镉 (Cd)	六价铬 (Cr(VI))	多溴联苯 (PBB)	多溴二苯醚 (PBDE)
电路板 PCBA's	X	O	O	O	O	O
电源供应器 Power Supply	X	O	O	O	O	O

(continued)

产品中有害物质的名称及含量 China EEP Hazardous Substances Information						
电机组件 Electromechanical Assemblies	X	O	O	O	O	O
<p>本表格依据 SJ/T11364 的规定编制 This table is compiled according to SJ/T 11364 standard.</p> <p>O: 表示该有害物质在该部件所有均质材料中的含量均在 GB/T26572 规定的限量要求以下. Indicates that the concentration of the hazardous substance in all homogeneous materials for the part is below the relevant threshold of the GB/T 26572 standard.</p> <p>X: 表示该有害物质至少在该部件的某一均质材料中的含量超出 GB/T26572 规定的限量要. Indicates that the concentration of the hazardous substance in at least one homogenous material of the part is above the relevant threshold of the GB/T 26572 standard.</p>						

Chemical safety



WARNING! GENERAL CHEMICAL HANDLING. To minimize hazards, ensure laboratory personnel read and practice the general safety guidelines for chemical usage, storage, and waste provided below. Consult the relevant SDS for specific precautions and instructions:

- Read and understand the Safety Data Sheets (SDSs) provided by the chemical manufacturer before you store, handle, or work with any chemicals or hazardous materials. To obtain SDSs, see the “Documentation and Support” section in this document.
- Minimize contact with chemicals. Wear appropriate personal protective equipment when handling chemicals (for example, safety glasses, gloves, or protective clothing).
- Minimize the inhalation of chemicals. Do not leave chemical containers open. Use only with adequate ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer's cleanup procedures as recommended in the SDS.
- Handle chemical wastes in a fume hood.
- Ensure use of primary and secondary waste containers. (A primary waste container holds the immediate waste. A secondary container contains spills or leaks from the primary container. Both containers must be compatible with the waste material and meet federal, state, and local requirements for container storage.)
- After emptying a waste container, seal it with the cap provided.
- Characterize (by analysis if necessary) the waste generated by the particular applications, reagents, and substrates used in your laboratory.
- Ensure that the waste is stored, transferred, transported, and disposed of according to all local, state/provincial, and/or national regulations.
- **IMPORTANT!** Radioactive or biohazardous materials may require special handling, and disposal limitations may apply.



AVERTISSEMENT ! PRÉCAUTIONS GÉNÉRALES EN CAS DE MANIPULATION DE PRODUITS CHIMIQUES. Pour minimiser les risques, veiller à ce que le personnel du laboratoire lise attentivement et mette en œuvre les consignes de sécurité générales relatives à l'utilisation et au stockage des produits chimiques et à la gestion des déchets qui en découlent, décrites ci-dessous. Consulter également la FDS appropriée pour connaître les précautions et instructions particulières à respecter :

- Lire et comprendre les fiches de données de sécurité (FDS) fournies par le fabricant avant de stocker, de manipuler ou d'utiliser les matériaux dangereux ou les produits chimiques. Pour obtenir les FDS, se reporter à la section « Documentation et support » du présent document.
- Limiter les contacts avec les produits chimiques. Porter des équipements de protection appropriés lors de la manipulation des produits chimiques (par exemple : lunettes de sûreté, gants ou vêtements de protection).
- Limiter l'inhalation des produits chimiques. Ne pas laisser les récipients de produits chimiques ouverts. Ils ne doivent être utilisés qu'avec une ventilation adéquate (par exemple, sorbonne).
- Vérifier régulièrement l'absence de fuite ou d'écoulement des produits chimiques. En cas de fuite ou d'écoulement d'un produit, respecter les directives de nettoyage du fabricant recommandées dans la FDS.
- Manipuler les déchets chimiques dans une sorbonne.

- Veiller à utiliser des récipients à déchets primaire et secondaire. (Le récipient primaire contient les déchets immédiats, le récipient secondaire contient les fuites et les écoulements du récipient primaire. Les deux récipients doivent être compatibles avec les matériaux mis au rebut et conformes aux exigences locales, nationales et communautaires en matière de confinement des récipients.)
- Une fois le récipient à déchets vidé, il doit être refermé hermétiquement avec le couvercle fourni.
- Caractériser (par une analyse si nécessaire) les déchets générés par les applications, les réactifs et les substrats particuliers utilisés dans le laboratoire.
- Vérifier que les déchets sont convenablement stockés, transférés, transportés et éliminés en respectant toutes les réglementations locales, nationales et/ou communautaires en vigueur.
- **IMPORTANT !** Les matériaux représentant un danger biologique ou radioactif exigent parfois une manipulation spéciale, et des limitations peuvent s'appliquer à leur élimination.

Biological hazard safety



WARNING! BIOHAZARD. Biological samples such as tissues, body fluids, infectious agents, and blood of humans and other animals have the potential to transmit infectious diseases. Conduct all work in properly equipped facilities with the appropriate safety equipment (for example, physical containment devices). Safety equipment can also include items for personal protection, such as gloves, coats, gowns, shoe covers, boots, respirators, face shields, safety glasses, or goggles. Individuals should be trained according to applicable regulatory and company/ institution requirements before working with potentially biohazardous materials. Follow all applicable local, state/provincial, and/or national regulations. The following references provide general guidelines when handling biological samples in laboratory environment.

- U.S. Department of Health and Human Services, *Biosafety in Microbiological and Biomedical Laboratories (BMBL)*, 5th Edition, HHS Publication No. (CDC) 21-1112, Revised December 2009; found at:
<https://www.cdc.gov/labs/pdf/CDC-BiosafetymicrobiologicalBiomedicalLaboratories-2020-P.pdf>
- World Health Organization, *Laboratory Biosafety Manual*, 3rd Edition, WHO/CDS/CSR/LYO/2004.11; found at:
www.who.int/publications/i/item/9789240011311



Documentation and support

Related documentation and references

Product Details	Product Description	Webpage URL
iBright™ Analysis Software Connect	Cloud-based iBright™ Analysis Software Connect is supplied with one free terabyte of digital data storage.	http://thermofisher.com/Connect
iBright™ Imager Firmware	Firmware updates and patches for all iBright™ Instruments are available for download on the Thermo Fisher website.	http://thermofisher.com/iBrightAnalysis
iBright™ Analysis Software	iBright™ Analysis Software (desktop version) is for managing, adjusting, analyzing, and annotating G2i digital image files that were captured on an iBright™ Imager.	
iBright™ Analysis Software – Secure	iBright™ Analysis Software - Secure (desktop version) supports 21 CFR Part 11 compliance.	
SAE Administrator Console Software	SAE Administrator Console Software supports 21 CFR Part 11 compliance by providing security, audit, and e-Signature (SAE) for iBright™ Instruments and iBright™ Analysis Software – Secure.	
SAE License for iBright™	License for Security, Audit, and e-Signature (SAE) on iBright™ Instruments and iBright™ Analysis Software – Secure.	

Document	Publication number	Description
<i>iBright™ Imaging Systems Documentation</i>	100085109	Provides instructions for obtaining service and technical support for the iBright™ Imaging Systems.
<i>iBright™ Imaging Systems Quick Reference</i>	100085108	Provides basic instructions for performing imaging with the iBright™ Imaging Systems.



Customer and technical support

Visit thermofisher.com/support for the latest service and support information.

- Worldwide contact telephone numbers
- Product support information
 - Product FAQs
 - Software, patches, and updates
 - Training for many applications and instruments
- Order and web support
- Product documentation
 - User guides, manuals, and protocols
 - Certificates of Analysis
 - Safety Data Sheets (SDSs; also known as MSDSs)

Note: For SDSs for reagents and chemicals from other manufacturers, contact the manufacturer.

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 at 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.

