

Countess™ 3 FL Automated Cell Counter

USER GUIDE

Catalog Number AMQAF2000

Publication Number MAN0019567

Revision F



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



Revision history: MAN0019567 F (English)

Revision	Date	Description
F	28 January 2026	<ul style="list-style-type: none">• New section, "Preset Protocols" was created under Chapter 3 - Protocols.• Chapter 5 - Cell count and cell viability assays was changed to Cell counting in brightfield.• Chapter 6 - Fluorescence assays was changed to Cell counting with fluorescent cells.• New Over-view section added to Chapter 5 and 6.• Content, images and tables were edited and formatted in the manual.
E.0	5 September 2023	Update latest draft of product manuals with latest screens and user workflow for SW#3
D.0	16 September 2022	Adding network-drive saving, Operational Qualification (OQ) functionalities, and scatter plot functionality.
C.0	9 June 2021	Changing "Templates" functions to "Protocols". Adding additional settings in protocol editing/creation to include Setup, Calculators, and Save.
B.0	28 January 2021	Add component to Contents table. Update technical specifications.
A.0	9 November 2020	New user guide for Countess™ 3 FL Automated Cell Counter.

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

DISCLAIMER: TO THE EXTENT ALLOWED BY LAW, THERMO FISHER SCIENTIFIC INC. AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF THE DOCUMENT OR THE PRODUCT.

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

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

©2020-2026 Thermo Fisher Scientific Inc. All rights reserved.

Contents

■	CHAPTER 1	Product information	7
		Contents	7
		Product description	8
		Upon receiving the instrument	8
		Register your instrument	8
		Exterior instrument parts	9
■	CHAPTER 2	Installation	10
		Operating environment	10
		Install the instrument	10
		Turn ON the instrument	10
		Sign in to the instrument with a new Thermo Fisher™ Connect Platform account	11
		Sign in to the instrument with an existing Thermo Fisher™ Connect Platform account ...	13
■	CHAPTER 3	Protocols	14
		Protocols screen	14
		Protocol functions	15
		Preset protocols	15
		Count parameters	16
		Load a protocol	17
		Create a protocol	18
		Edit a protocol	18
		Import a protocol	19
		Export a protocol	20
		Delete a protocol	22
■	CHAPTER 4	Slide preparation and operation	23
		Recommendations	23
		Load Countess™ Cell Counting Chamber Slide	24
		Load the Countess™ Reusable Slide	25

■	CHAPTER 5	Cell counting in brightfield	27
	Overview		27
	Count cells in brightfield		27
	Capture and count		27
	Next steps		30
	View results		30
	Results screen for colorimetric viability assays		30
	Identify objects counted		31
	Results screen		31
	Identify cells counted in cell count and cell viability assays		31
	Histogram count results		33
	View histogram		33
	View scatter plot		34
	Gate count results		34
	Gating screen		34
	Gate count results		35
	Save as new protocol		35
	Edit and save a new protocol		35
■	CHAPTER 6	Cell counting with fluorescent cells	37
	Count cell fluorescence		37
	Overview		37
	Capture and count		37
	Rapid capture		37
	Customize capture		38
	Next steps		40
	View results		41
	Results screen for BF or FL-based counts		41
	Results screen for FL-based counts		41
	Results screen for BF-based counts		42
	Identify objects counted		43
	Identify cells counted in fluorescence assays		43
	Histogram count results		43
	View histogram for cell fluorescence assays		43
	View scatter plot		45
	Gate count results		45
	Gating screen		45
	Gate count results		46
	Save as new protocol		46
	Edit and save a new protocol		46

■	CHAPTER 7	Calculators	48
		Pre-Dilution Calculator	48
		Cell Splitting Calculator	50
■	CHAPTER 8	Save results	52
		Save screen	52
		Save procedure	52
		Report	55
		Report file	55
■	CHAPTER 9	Settings overview	57
		Settings screen	57
		Connect to a wireless network	57
		Connect to a network drive	57
		Software update	58
		Guidelines for software update	58
		Update the Countess™ 3/3 FL Cell Counter software	58
		Set the date and time	59
		Change your cloud region	62
■	CHAPTER 10	Change EVOS™ light cubes	64
		Install or change an EVOS™ light cube	64
■	CHAPTER 11	Maintenance	66
		Instrument care	66
		General guidelines for care	66
		Power supply	66
		Clean the cell counter	67
		Introduction	67
		Clean the touchscreen	67
		Clean the instrument case	67
		Decontaminate the instrument	68
■	CHAPTER 12	Installation/Operational Qualification	69
		Generate and activate a license	69
		Execute the Installation Qualification (IQ)	70
		Execute the Operational Qualification (OQ)	75

■	APPENDIX A	Troubleshooting	82
■	APPENDIX B	Technical specifications	85
		Physical characteristics	85
		Technical specifications	85
		Optics	85
		Analysis slide	86
		Networking requirements	86
■	APPENDIX C	EVOS light cubes	87
		LED illumination	87
		EVOS™ light cubes	87
■	APPENDIX D	Countess™ 3 FL Automated Cell Counter and accessories ...	89
		Accessory products	89
■	APPENDIX E	CSV file format definition	91
		Overview	91
■	APPENDIX F	Safety	95
		Chemical safety	96
		Biological hazard safety	97
■	APPENDIX G	Documentation and support	98
		Customer and technical support	98
		Limited product warranty	98



Product information

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

Contents

The Countess™ 3 FL Automated Cell Counter is shipped with the components listed below.

Component	Quantity
Countess™ 3 FL Automated Cell Counter	1 each
Power cord with 4 adapters (for U.S./Canada/Taiwan/Japan, Europe, or UK)	1 each
Countess™ Cell Counting Chamber Slides (50 slides/box)	1 box
Countess™ 3 FL Disposable Slide Holder	1 each
Countess™ 3 FL Reusable Slide Holder	1 each
Countess™ 3 FL Light Cube Removal Tool	1 each
Countess™ 3 USB Drive	1 each
Countess™ 3 FL Automated Cell Counter Quick Reference Card	1 each
Wi-Fi Dongle	1 each
Trypan Blue Stain	1 each
SafeCount™	1 each

Product description

The Countess™ 3 FL Automated Cell Counter is a fully automated, 3-channel cell counter and assay platform that uses EVOS™ light cube technology, exceptional optics, and image analysis algorithms to analyze fluorescently labeled cells or colorimetric viability stained samples (for example, Trypan Blue) in suspension.

- The Countess™ 3 FL Automated Cell Counter offers an intuitive user interface with the option to save data and generate a report, which can then be transferred to a PC using the USB drive supplied with the instrument or available separately.
- The cells to be counted are loaded into the instrument either in disposable Countess™ Cell Counting Chamber Slides or in glass Countess™ 3 FL Reusable Slides (page 24). Each chamber slide contains two enclosed chambers to hold the sample to allow you to measure two different samples or perform replicates of the same sample.
- The instrument takes 10–20 seconds per sample for a typical cell count in the brightfield channel and is compatible with a wide variety of eukaryotic cells. In addition to cell count and viability, the Countess™ 3 FL Automated Cell Counter also provides information on cell size.
- In addition to the brightfield channel, the Countess™ 3 FL Automated Cell Counter can accommodate two interchangeable EVOS™ fluorescent light cubes (page 87), enabling it to be used for multiple-fluorescence research applications.
- When equipped with EVOS™ light cubes, the Countess™ 3 FL Automated Cell Counter can be used to perform fluorescence assays for cells in suspension, including simultaneous counts of cells stained with two different fluorescent dyes, GFP and TxRed expression, apoptosis, and cell viability (live, dead, and total cells).
- For laboratories operating under 21 CFR Part 11 requirements, the Countess™ 3 and Countess™ 3 FL instruments can be configured with the SAE console. This console enables controlled user access, comprehensive electronic record auditing, and the use of electronic signatures. The user interaction with the SAE Admin Console is described in the SAE Admin Console user guide for use with Countess™ 3 and 3 FL Automated Cell Counters (Pub. No. [MAN0025371](#))

Upon receiving the instrument

Examine the instrument carefully for damage incurred during transit. Ensure that all parts of the instrument, including accessories listed above, are included with the product. Damage claims must be filed with the carrier; the warranty does not cover in-transit damage.

See “Install the instrument” on page 10 for instructions on installing the instrument.

Register your instrument

Visit www.thermofisher.com/registercountess to register your instrument. You will be asked to supply the serial number, your name, and your contact details. Registering your instrument ensures that you will receive notifications of software upgrades and information on new assays for use with the Countess™ cell counter.

Exterior instrument parts



- ① **Touchscreen display:** The 7-inch capacitive display is the main user interface of the Countess™ 3 FL Automated Cell Counter. It contains the buttons for all instrument functions and displays data from the cell count.
- ② **Front USB port:** Allows you to transfer and save the cell count data and image to an external computer for record keeping and printing purposes. You can use the USB drive supplied with the instrument or any other standard, FAT32-, exFAT- or NTFS-formatted USB drive for data transfer. If desired, you can plug in a USB mouse into the rear USB port for instrument control.
- ③ **Slide port:** Used to insert the analysis slide containing the sample into the counter.
The Countess™ 3 FL Automated Cell Counter accepts both the disposable Countess™ Cell Counting Chamber Slides and the glass Countess™ 3 FL Reusable Slides via interchangeable, slide-specific carriers. For more information, see Chapter 4, “Slide preparation and operation”.
- ④ **Back panel:** Allows access to the optional EVOS™ light cubes and provides storage for the light cube tool and the reusable slide carrier. The back panel is secured to the instrument by two captive ¼-turn fasteners.
- ⑤ **Power switch:** ON/OFF rocker switch is the main power switch. It is not necessary to use the power switch for day-to-day operation of the instrument.
- ⑥ **EVOS™ light cubes:** Allow the Countess™ 3 FL Automated Cell Counter to analyze fluorescently labeled samples. Can accommodate two fluorescent light cubes. For more information, see “EVOS™ light cubes” on page 87.
- ⑦ **Rear USB port:** An additional USB port to perform functions such as plugging in a USB mouse for instrument control.
- ⑧ **Power input jack:** Connects the instrument to an electrical outlet through the supplied power cord and the appropriate plug, based on the electrical outlet configuration in your country.

2

Installation

Operating environment

- Place the instrument on a level surface away from vibrations emanating from other pieces of equipment.
- Allow at least 5 cm (2 in) free space at the back of the instrument to allow for proper ventilation and prevent overheating of electronic components.
- Set up the instrument away from direct light sources, such as windows. Ambient room lighting can enter the imaging path and affect the image quality.
- Operating temperature range: 4°–32°C (40°–90°F).
- Relative humidity range: <80%.

IMPORTANT! Do not position the instrument so that it is difficult to turn off the main power switch located on the back of the instrument.

In case of an instrument malfunction, turn the main power switch to the OFF position and disconnect the instrument from the wall outlet.

Install the instrument

1. Unpack the instrument and place the instrument on a flat, level, dry surface.
2. Remove the thin plastic protector film from the touch-screen display.
3. Plug one end of the power cord appropriate for your region into the instrument.
4. Plug the power cord into the electrical outlet. Be sure to use only the power cord supplied with your instrument. Powering the instrument with an unapproved power cord may damage the instrument.

Turn ON the instrument

After unboxing, upon switching on the instrument for the first time, it will run through a series of set-up screens.

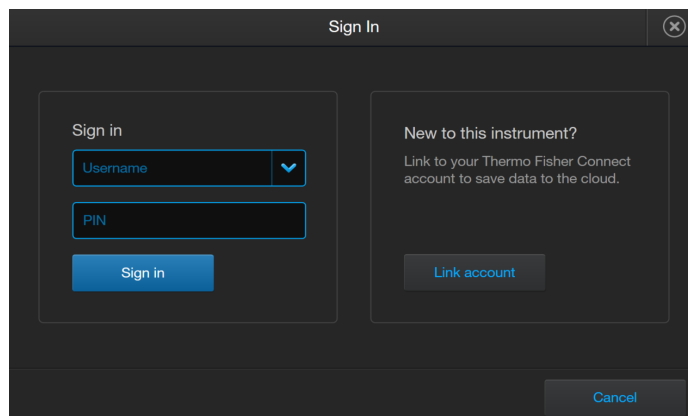
- Read the End User License Agreement and press **Agree** to proceed with start up.
- Set the **Date** and **Time** using the buttons (see “Set the date and time” on page 59). Press **Next**.
- Upon completion, press **Done**. You will proceed to the Home screen and can begin using the instrument as described below.

1. Turn on the instrument by flipping the power switch on the back of the instrument to the **ON** position. See “Exterior instrument parts” on page 9.
The instrument will walk you through an initial Out-of-Box Experience to set up the instrument for first use.
2. From the **Home** screen, you can proceed immediately to the assays by inserting a slide. See Chapter 5, “Cell counting in brightfield”.
Alternatively, you can change or add a protocol in step 3 below or change instrument settings in step 4 below.
3. To change the current protocol or add a new protocol to the instrument, press **⋮ (Protocols)** .
Protocols allows you to create customized count preferences (i.e., gate counts based on cell size, brightness, circularity and/or fluorescence intensity). See “Load a protocol” on page 17.
4. To change instrument settings, press **⚙ (Settings)** .
Settings allows you to update the Countess™ 3 FL instrument software, customize instrument settings, view instrument details, export error logs, and change up to two EVOS™ light cubes. See “Settings screen” on page 57.

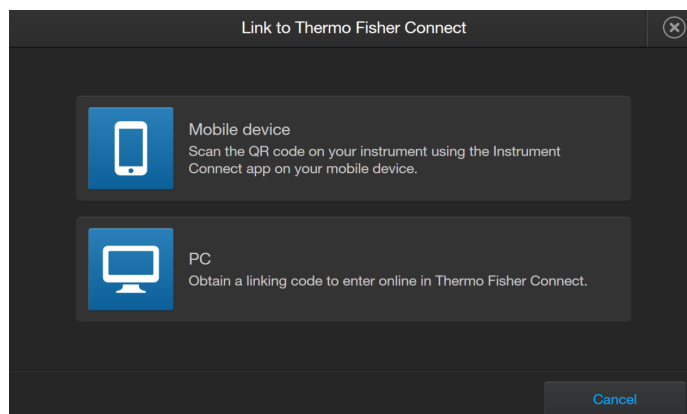
Sign in to the instrument with a new Thermo Fisher™ Connect Platform account

Note: A Wi-Fi adapter needs to be installed and a Thermo Fisher™ Connect Platform account is required for linking. See <http://thermofisher.com/connect> for details on using the Connect platform.

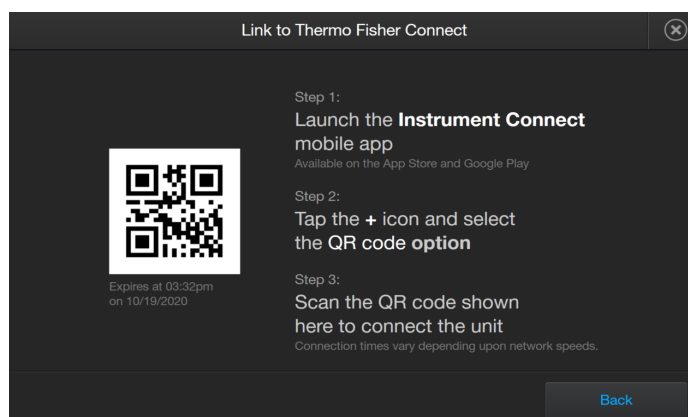
1. Press **👤 (Sign in)** from the **Home** screen.
2. Press **Link account**.



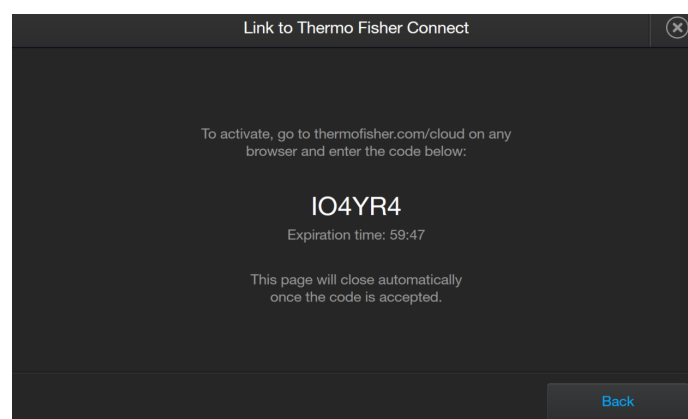
3. Choose a method to connect.



- Press **Mobile device**. Scan the QR code on your instrument using the Instrument Connect app on your mobile device. Follow the steps shown on the page.




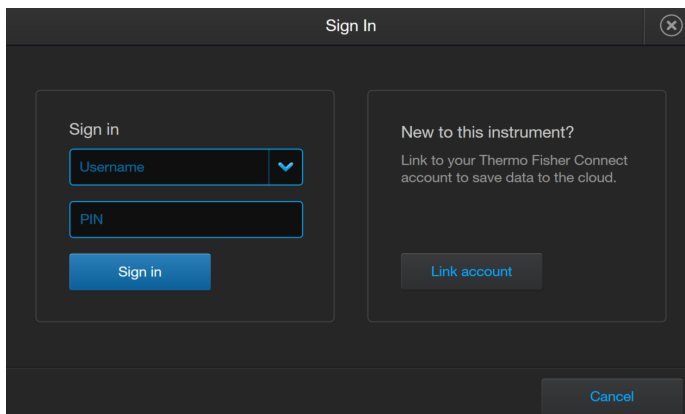
- Press **PC** to obtain a linking code to enter online in your Connect platform account.

4. A completion screen will show. Press **Done**. You then return to the **Home** screen and are logged into your account.

Once set up, you will then only need to sign in to your instrument account using your **Username** and **PIN**.

Sign in to the instrument with an existing Thermo Fisher™ Connect Platform account

1. Press  (**Sign in**) from the **Home** screen.
2. Select your **Username** from the drop-down box. Enter your **PIN**.
3. Press **Sign in**.



The image shows a 'Sign In' dialog box with a dark background. At the top, it says 'Sign In' with a close button (X) on the right. The dialog is divided into two main sections. The left section is titled 'Sign in' and contains a dropdown menu labeled 'Username' with a downward arrow, a text input field labeled 'PIN', and a blue 'Sign in' button. The right section is titled 'New to this instrument?' and contains the text 'Link to your Thermo Fisher Connect account to save data to the cloud.' and a 'Link account' button. At the bottom right of the dialog is a 'Cancel' button.

Protocols screen

Protocols allows you to create and save customized protocols. Each custom protocol defines the count parameters (size, brightness, circularity, and fluorescence intensity) for a consistent and streamlined workflow.



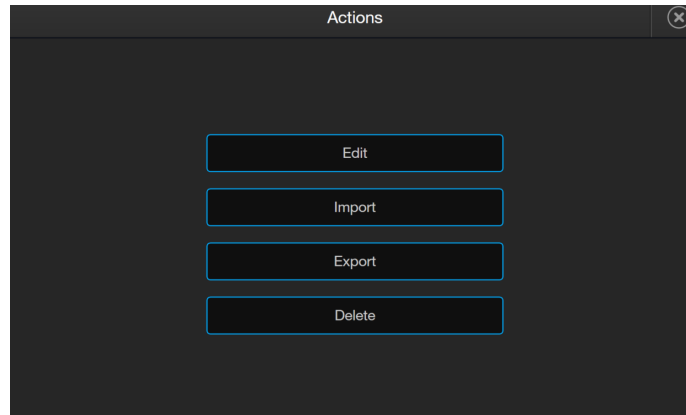
Name	Channels	Last Modified
Default Protocol	All	🔒
Trypan Blue	Live, Dead	🔒
SafeCount	Live, Dead	🔒
ReadyCount Viability Stain	GFP, TX Red	🔒
ReadyCount Dead Stain	TX Red	🔒
ReadyCount Nuclear Stain	DAPI	🔒

Actions: [Actions] [Create] [Cancel] [Load]

- Protocols can be accessed from the **Home** and **Gating** screens.
- The currently selected protocol is displayed at the bottom of the **Home** screen.
- Preset protocols can be hidden from Protocols screen through **Settings/Instrumentsettings/preset protocols**.
- Automatic instrument functions and count parameters are defined in the **Edit protocol** screen. See “Create a protocol” on page 18 or “Edit a protocol” on page 18.
- The Default protocol and preset protocols contain count settings and cannot be edited.
- The count parameters specified in the selected protocol are applied to all new cell counts.
- Protocols can be sorted by pressing on the **Name**, **Channels**, or **Last Modified** headers.
- If you have already performed a count, loading a new protocol from the **Results** screen applies the count preferences to the current count results (total cells, viability, etc.) and to all new counts.

Protocol functions

The **Actions** screen allows you to edit, import, export, and delete protocols. **Edit/View** can be used to modify the currently selected protocol. Preset protocols can not be edited, only viewed, indicated by the lock icon. **Import**, **Export**, and **Delete** can be used on multiple protocols.



Preset protocols

The Countess™ 3 FL Automated Cell Counter includes several Preset protocols designed to streamline cell counting for users. These protocols are pre-configured with specific settings to aid users and provide a starting point to define cells of interest. The specific settings within the Preset protocols include the light cubes needed, count mode, reagent dilution, and gating on brightness, size, and/or circularity. The Preset protocols can be found in the protocols menu. Users can hide these protocols by navigating to **Settings > Instrument Settings > Preset protocols** and unchecking **Show**. Additionally, in this menu, users can find the kit SKU for the reagents needed for each protocol.

The following Preset protocols are available:

Note: The default protocol behavior is dependent on presence or absence of light cubes.

Table 1

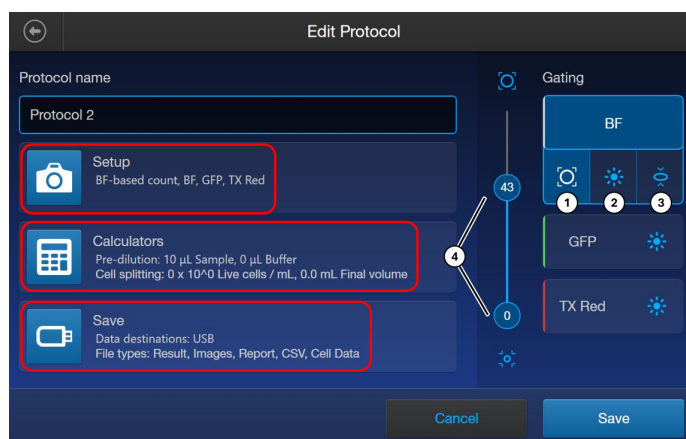
Protocol name	Cat. No.	Light cube	Dilution	Count mode	Gating
Default		None	1:1	BF-based	None
Default		Present	None	BF-based	None
Trypan Blue	T10282	NA	1:1	BF-based	None
SafeCount™	A40008024	NA	1:1	BF-based	None
ReadyCount™ Viability Stain	A49905	GFP/TxRed	1:1	FL-based	Size minimum 7; GFP brightness minimum 25
ReadyCount™ Dead Cell Stain	A49903	TxRed	1:1	BF-based	None
ReadyCount™ Nuclear Stain	A49904	DAPI	1:1	FL-based	None

Table 1 (continued)

Protocol name	Cat. No.	Light cube	Dilution	Count mode	Gating
PBMC Viability Stain	A49905	GFP/TxRed	1:1	FL-based	None
PBMC Trypan Blue	T10282	NA	1:1	BF-based	Size minimum 4; circularity minimum 50
Isolated Nuclei	A49905	GFP/TxRed	1:1	FL-based	None


Count parameters

Count parameters are adjusted in the **Edit Protocol** screen using the **Gating** sliders. Moving the top and bottom gate boundaries adjusts the range of size, brightness, and circularity parameters. **Live** or **Dead** protocols are created when only the brightfield (BF) channel is selected. For a fluorescence (FL) count using the Countess™ 3 FL instrument, select the individual channels.

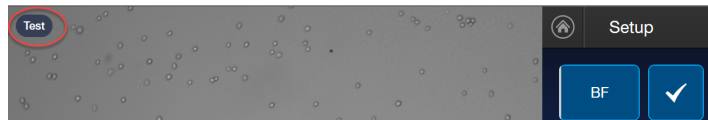


- ① **Size Gating** : Moving the top gate down removes larger objects; moving it up includes larger objects. Moving the bottom gate up removes smaller objects; moving it down includes smaller objects.
- ② **Brightness Gating**: Moving the top gate down removes brighter objects; moving it up includes brighter objects. Moving the bottom gate up removes dimmer objects; moving it down includes dimmer objects.
- ③ **Circularity Gating**: Moving the top gate down removes more circular objects; moving it up includes more circular objects. Moving the bottom gate up removes less circular objects; moving it down includes less circular objects.
- ④ **Gating Boundary Sliders**
 - **Setup** – Press **Setup** to select the required channels (Example., BF, GFP, TX Red) for capture. BF-based counts using the BF image is utilized by default (BF cannot be deselected for BF based counts). Select FL-based count to obtain a count using FL images (capturing the BF image is optional once FL-based count is selected). For details, see Chapter 6, “Cell counting with fluorescent cells”.
 - **Calculators** – Press **Calculators** to access the Pre-Dilution and Cell Splitting Calculators and store default calculator values. For detailed protocols, see Chapter 7, “Calculators”.
 - **Save** – Press **Save** to save default save settings and your data to the cloud or a USB. For details, see “Save screen” on page 52.

Load a protocol

1. Press  (**Protocols**).
2. Press the **Desired protocol** to select, then press **Load**.
3. To return to the previous screen without loading the new protocol, press **Cancel**.
The instrument will keep the saved protocol, but will return to the previous screen without loading it.

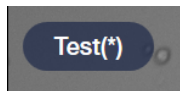
Note: The **Protocol name** will be displayed in the upper left corner of the **Count** screen.



Note: The **Protocol name** will be displayed at the bottom of the **Home** screen.

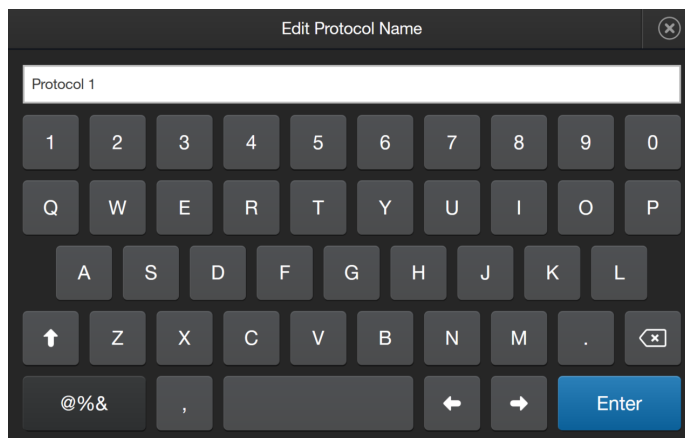


If a saved protocol has been changed and there are unsaved changes, an asterisk (*) will follow the name.



Create a protocol

1. Press **⌘** (Protocols).
2. Press **Create**. To assign a name to the new protocol or to change the name of the existing protocol, press the **Protocol name** text box to open the keypad.



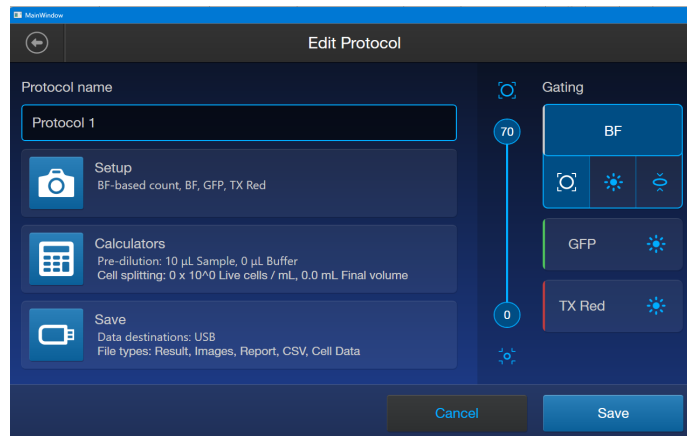
3. Type in the **Desired Protocol name**. To enter symbols, press the symbol (@%&) key. To return to the alpha-numeric keypad, press **ABC**.
4. Press **Enter** to save the name and return to the **New Protocol** screen. Press **X** to return to the **New Protocol** screen without saving the name.
5. Press **Save** to save the new protocol.
6. Press **Cancel** or **⏪ (Back)** to return to the **Protocols** screen without saving.

Edit a protocol

1. Press **⌘** (Protocols).
2. Select an existing protocol, then press **Actions**.

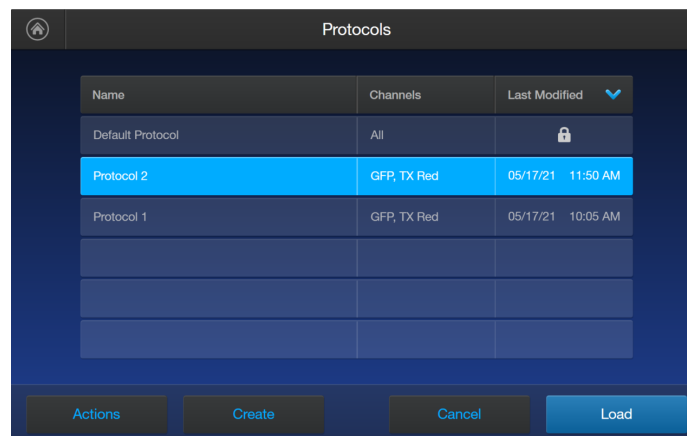
Note: The Default and Preset protocols cannot be edited.

3. Press **Edit**.



The **Edit** screen for the selected protocol opens.

4. Press the appropriate channel (BF or other dye) and edit the desired parameters to create a custom protocol. Additionally, use the **Setup**, **Calculators**, and **Save** buttons to set up any of the needed parameters to complete the custom protocol. For details, see “Count parameters” on page 16.
5. Press **Save**.

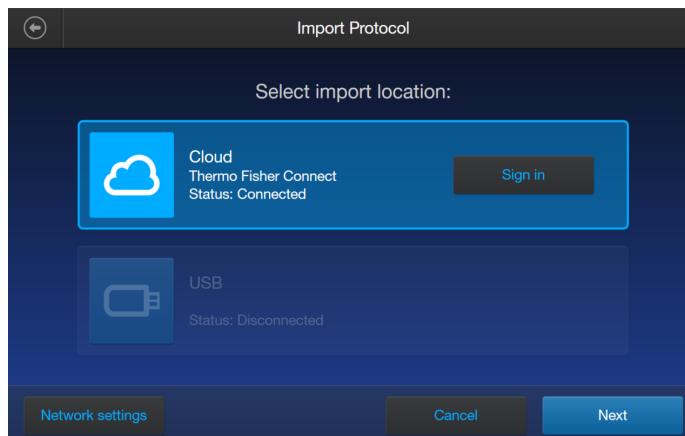


From the **Protocols** screen, select the **Desired Protocol** and press **Load** to use that protocol to read the sample.

Import a protocol

1. Press **⌘** (**Protocols**).
2. Press **Actions**.
3. Press **Import**.

4. Press the import location **Cloud** or **USB**.
 - a. For **Cloud**, press **Sign in** to log in to your Thermo Fisher Connect account. If the **Cloud** is not active and disconnected, press **Network settings** to connect to a network, then press ⏪ (**Back**) to try and sign in again.
 - b. For **USB**, insert your USB drive into the slot on the front of the instrument.



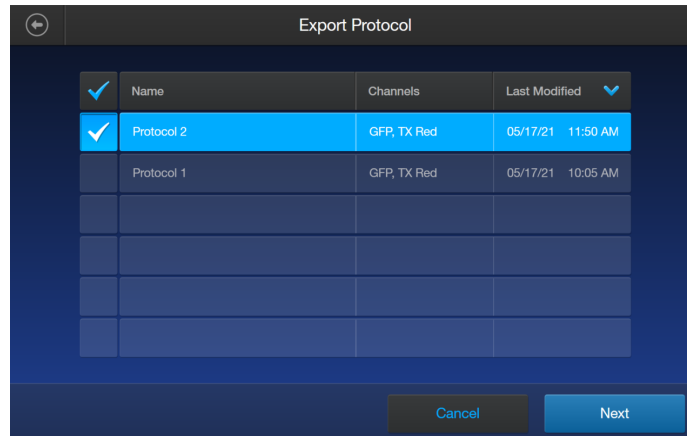
5. Press **Next**.
6. Select one or more protocols from the list.
7. Press **Import**.
8. Once complete, press **Done**.

Export a protocol

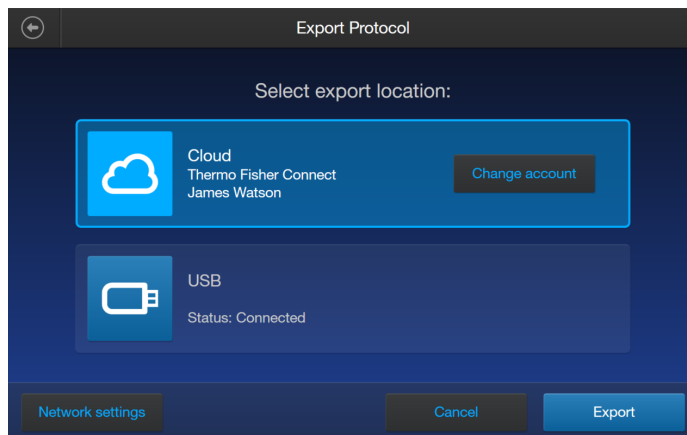
1. Press ⏴ (Protocols).
2. Press **Actions**.
3. Press **Export**.

4. Select one or more protocols from the list.

Note: The Default and Preset protocols cannot be exported.



5. Press **Next**.
6. Press the export location **Cloud** or **USB**.
 - a. For **Cloud**, press **Sign in** to log in to your Thermo Fisher Scientific Connect account. If the **Cloud** is not active and disconnected, press **Network settings** to connect to a network, then press **⬅ (Back)** to try and sign in again.
 - b. For **USB**, insert your USB drive into the slot on the front of the instrument.

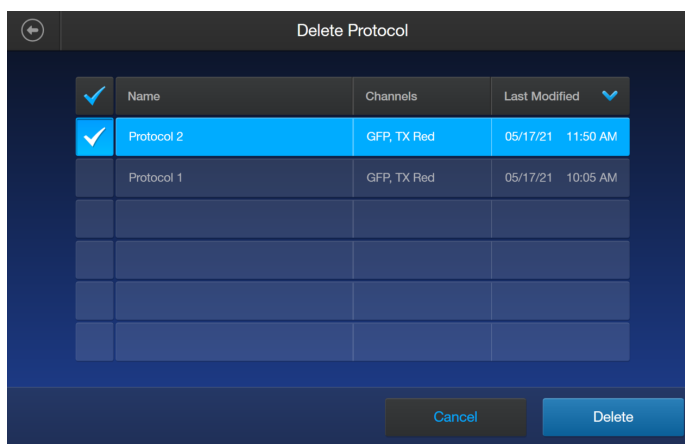


7. Press **Export**.
8. Once complete, press **Done**.

Delete a protocol

1. Press **⌘** (Protocols).
2. Press **Actions**.
3. Press **Delete**.
4. Select one or more protocols from the list.

Note: The Default and Preset protocols cannot be deleted.



5. Press **Delete**.
The system will prompt you to confirm the deletion, then press **Delete** again.
6. Once complete, press **Done**.

4

Slide preparation and operation

The Countess™ 3 and 3 FL cell counters accept disposable Countess™ Cell Counting Chamber Slides and glass Countess™ Reusable Slides on interchangeable, slide-specific carriers.

Recommendations

To obtain the best results, follow these recommendations:

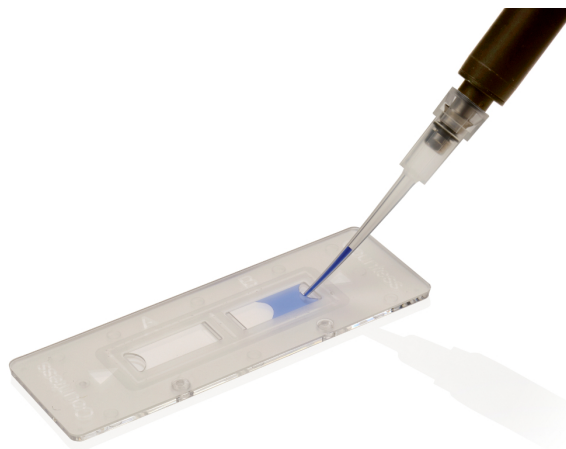
- Ensure that the cell sample is homogeneously mixed.
- The measurement range extends from 1×10^4 to 1×10^7 cells/mL, but the optimal range is 1×10^5 to 4×10^6 cells/mL.

IMPORTANT! A warning will appear on the Results screen when concentrations are out of optimal range.

- For accurate results in cell viability assays, ensure that the counting area is covered with the cell suspension and count the cells immediately after staining per the assay protocol.
- Do **not** press the optical surfaces of the chamber slides. Hold the slides by the edges.
- Be careful to avoid forming bubbles in the sample.

Load Countess™ Cell Counting Chamber Slide

1. Prepare the sample: If using colorimetric viability stains Trypan Blue or SafeCount™, or fluorescence ReadyCount™ stains, add 10 µL of your cell suspension to 10 µL of stain. Mix the sample mixture well by pipetting it up and down a few times and follow protocol as indicated.
2. Gently pipet 10 µL of the sample into the half moon-shaped sample loading area. The sample is loaded into the chamber through capillary action.



3. Let the sample mixture settle in the chamber for 30 seconds, and then insert the slide into the slide adapter. You will hear a soft click, once the slide is pushed in correctly.

Note: The instrument reads one side of the slide at a time, so insert the sides appropriately.

Note: The instrument is shipped with the slide adapter already installed.



4. Use the touchscreen to run the sample count. See Chapter 5, “Cell counting in brightfield” and Chapter 6, “Cell counting with fluorescent cells” for more details.
5. To remove the slide, push the slide gently into the instrument until it clicks and the slide springs out. Grasp the slide and pull it out the rest of the way.

Note: After using the Countess™ Cell Counting Chamber Slides, appropriately dispose of them as biohazardous waste. **Do not reuse the disposable chamber slides.**

6. (Optional) To remove the slide adapter, gently squeeze the tabs and pull the adapter completely out of the instrument.

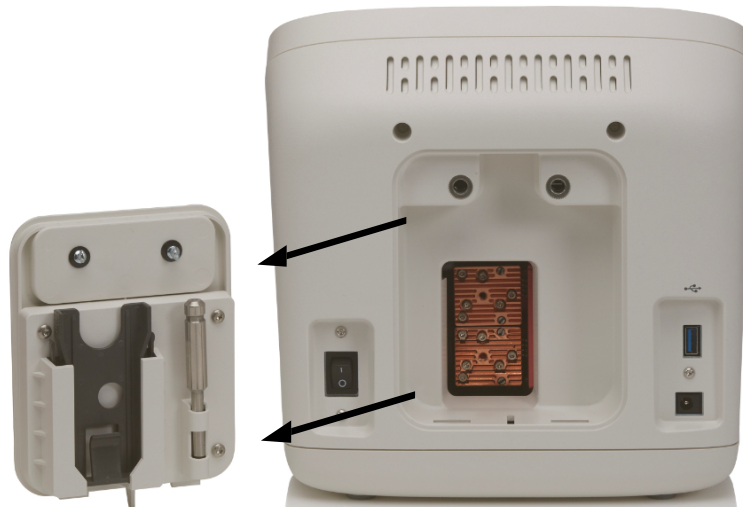
Load the Countess™ Reusable Slide

1. Before loading your sample into the Countess™ Reusable Slide, place a cover slip on the counting chamber, making sure the cover slip is clean and free of grease.
2. Prepare the sample: If using colorimetric viability stains Trypan Blue or SafeCount™, or fluorescence ReadyCount™ stains, add 10 µL of your cell suspension to 10 µL of stain. Mix the sample mixture well by pipetting it up and down a few times and follow protocol as indicated.
3. Gently pipet 10 µL of the sample into the sample inlet, allowing capillary action to draw the sample into the counting chamber. A properly loaded counting chamber should have a thin, even film of fluid under the cover slip.



Note: Each chamber in the Countess™ Cell Counting Chamber Slide or the Countess™ 3 FL Reusable Slide has a 10-µL sample capacity. Do not overfill the slide chambers.

4. To use the Countess™ Reusable Slide, unlatch the back panel of the Countess™ 3 FL Automated Cell Counter with the two captive ¼-turn fasteners that secure the back panel on the rear of the instrument.



5. Remove the reusable slide holder from inside of the back panel.



6. Load the reusable glass slide into the reusable slide holder.



7. Insert the reusable slide holder containing the slide into the slide port and gently push into the instrument until it clicks into place.
8. Use the touchscreen to run the sample count. See Chapter 5, “Cell counting in brightfield” and Chapter 6, “Cell counting with fluorescent cells” for more details.
9. To remove the reusable slide holder, push the slide gently into the instrument until it clicks and a spring pushes the slide out. Grasp the slide and pull it out the rest of the way.
10. *(Optional)*: To count the second sample present on the reusable slide, simply remove the slide from the holder, rotate, and reinsert the slide into the holder so that the second sample is aligned with the sample viewing hole.
11. After using the Countess™ Reusable Slide, rinse the glass slide and cover slip with water, and then clean with 70% ethanol. Use Kimwipes™ laboratory tissues to clean and dry the slides, as needed.
12. Store the reusable slide holder again in the panel on the back of the instrument.

5

Cell counting in brightfield

Overview

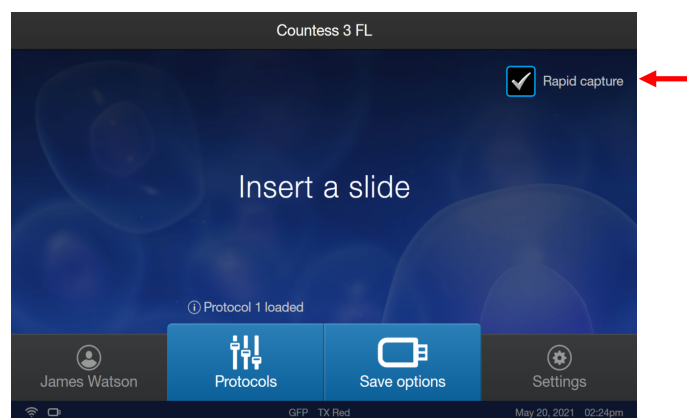
Countess™ 3 FL Automated Cell Counters count colorimetric viability stained samples using a brightfield image and two key colorimetric viability stains, Trypan Blue and SafeCount™. Both stains are used at a 1:1 dilution with the cell suspension. These viability stains are effective because they make dead cells appear dark, allowing for accurate differentiation between live and dead cells using the brightfield image. The Countess™ 3 Automated Cell Counters, leverages image analysis algorithms to provide detailed results on starting cell concentration (prior to staining), viability, and aggregate formation.

Count cells in brightfield

Capture and count

Rapid capture


To bypass adjusting individual parameters before the count, you can select **Rapid capture** at the top of the **Home** screen. This will automatically set the parameters and perform a count as soon as the slide is inserted into the instrument.



1. Load the samples into the instrument as described in Chapter 4, “Slide preparation and operation”.
2. See “View results” on page 30.

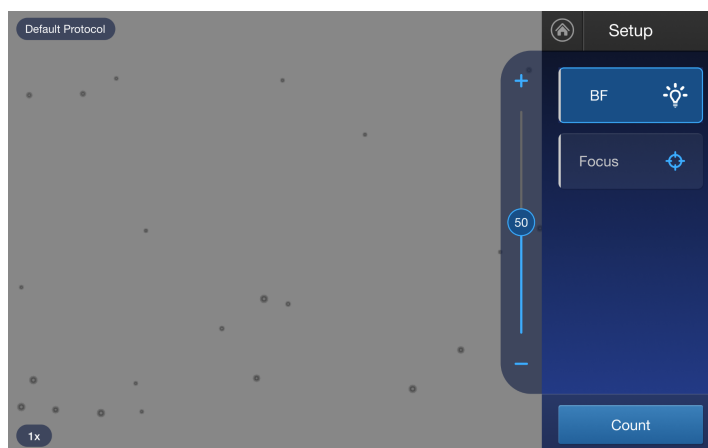
Customize capture

To view the image and adjust parameters prior to the count, follow the procedure below:

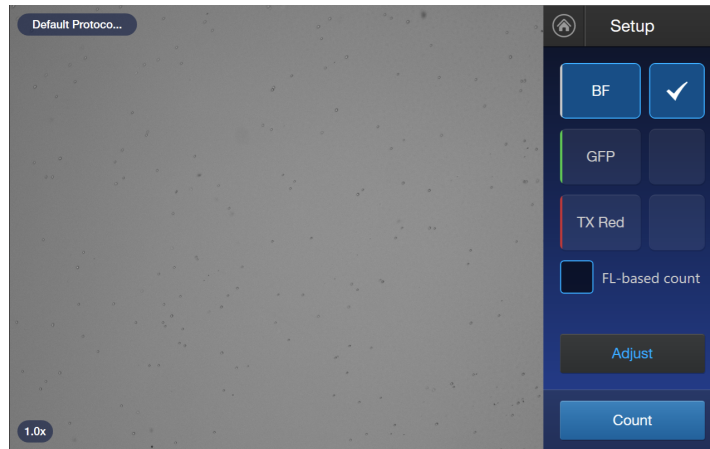
1. Load the sample into the sample slide as described in “Load Countess™ Cell Counting Chamber Slide” on page 24.
2. (Optional): Press  (**Protocols**) and load the desired gating parameter settings as described in “Load a protocol” on page 17.
3. Insert the sample slide into the instrument. Make sure that the sample side is inserted completely into the instrument. You will hear a soft click if the slide is pushed in correctly.
4. When the slide is inserted, the instrument automatically illuminates the sample, sets the intensity of brightfield illumination, and auto focuses on the cells.

Note: Auto lighting is optimal for producing accurate results, but if needed, can be disabled by adjusting the lighting slider.

5. (Optional): Set light intensity and/or focus. The light source slider controls the LED intensity and allows you to adjust the image brightness. The auto-focus function searches the nominal focus, or Z-point (depth) to provide fine focus to the sample. If the focus is modified and applied, the new focus value will become the nominal focus value for subsequent counts in the session.
 - If no light cubes are inserted:



- If EVOS™ light cubes are inserted: first press **Adjust**, and then select **BF** (brightfield) as the light source. Set the LED intensity, then press **Apply** to return to the **Setup** screen.



6. Press **Count**.




Note: If your instrument is equipped with an EVOS™ light cube, make sure that only the **BF** (brightfield) checkbox is selected before capturing the image, since this is a BF-based count.

The instrument captures the image and displays the following results: total concentration, percentage and concentration of Live and Dead cells, and percentage of Aggregates (clusters of three or more cells with touching membranes). For more information, see “View results” on page 30.


Autosaving will occur when the count is completed if you have **Autosave** enabled from the **Save Options** screen.



Next steps

- To identify the objects (for example, cells) counted as live, press . To identify objects counted as dead, press . To identify aggregates, press . See “Identify cells counted in cell count and cell viability assays” on page 31.

Note: The **Live** and **Dead** buttons are enabled by default after the image has been counted.

- To see the distribution of live and dead cells in a graphical format, press  (**Histogram**). See “Histogram count results” on page 33.
- To gate the results by object size, brightness, or circularity, press **Gating** to go to the **Gating** screen. See “Gate count results” on page 34.

Note: You can save the changes you make to the size, brightness, or circularity parameters in the **Gating** screen to the current protocol or as a separate protocol. See “Save as new protocol” on page 35.

- To calculate the volume of cell sample and buffer needed to reach a desired concentration based on the count results, press **Calculators** to open the Calculator applications. See Chapter 7, “Calculators”.
- To permanently save the results, press **Save**. See Chapter 8, “Save results”.
- To perform a new count, push the slide to eject, then insert a new sample slide.

View results

Results screen for colorimetric viability assays

The **Results** screen for cell count and cell viability assays performed using colorimetric viability stains displays a composite image of the objects counted and the results of the cell count and viability calculations. The results shown include:

- Total Concentration: The overall concentration of cells in the sample.
- Percentage and concentration of live cells: The number and percentage of live cells in the sample.
- Percentage and concentration of dead cells: The number and percentage of dead cells in the sample.
- Percentage of aggregates: The percentage of cell clusters (aggregates) consisting of three or more cells with touching membranes.



Note: When performing cell counts and viability assays with colorimetric stains, the counting algorithm assumes a 1:1 dilution of cells in the stain and takes this dilution into account when calculating the total cell concentration. The cell concentration displayed on the **Results** screen reflects the original cell concentration before dilution.



Identify objects counted


Results screen

The **Results** screen allows you to identify the objects (i.e., cells) counted in each channel and included in the count results for further review. After reviewing the marked objects, you can adjust the threshold for size, brightness, and/or circularity as desired for your application. See “Gate count results” on page 35.

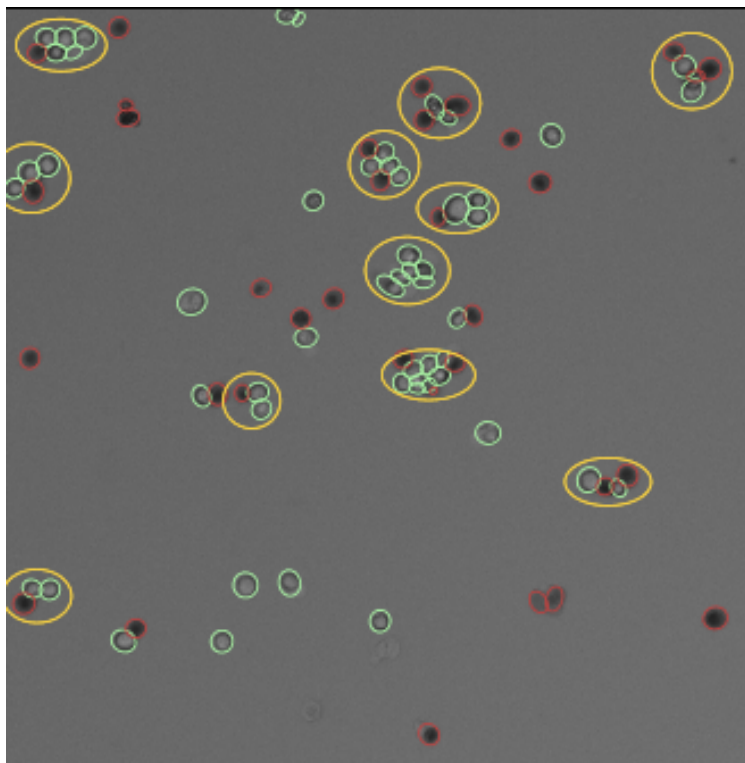
Identify cells counted in cell count and cell viability assays

On the **Results** screen, live, dead, or aggregates (clusters) of cells can be selected.

1. To view **Live cells**, press  on the right panel. This marks all Live cells in green.
2. To view **Dead cells**, press  on the right panel. This marks all Dead cells in red.

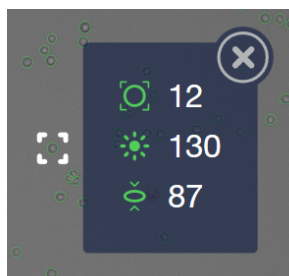
- To view **Aggregates of cells**, press  on the right panel. This marks all the aggregated cells with a yellow circle.

Note: You can select one, two, or all three options. The detailed image below shows Live (green), Dead (red) cells, and Aggregates (yellow) are selected.



- To unmark the cells, press the appropriate colored button again.

Individual cells can be selected to view their specific size, brightness, and circularity gating settings. Zooming in on the image will allow for easier selection of individual objects. Click the desired cell and a pop-up displays the parameters. Press **X** to close the pop-up.




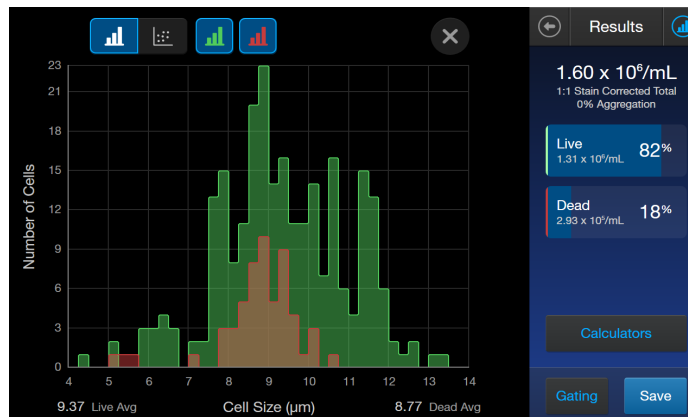
Histogram count results

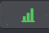






View histogram

For cell count and cell viability assays performed in the brightfield channel, you can view the distribution of cells (live and/or dead) based on size in a graphical format.

Note: You can view the histogram on the **Results** and **Gating** screens.

1. To view the histogram showing the distribution of live and/or dead cells based on cell size, press  (**Histogram**).



2. To view the distribution of only the live or dead cells, press the corresponding  (**Live**) (green) or  (**Dead**) (red) button.
The graph will automatically update and display the distribution of cells based on size only in the selected population.
3. (Optional): On the **Gating** screen, use the  (**Size**),  (**Brightness**), and  (**Circularity**) sliders, to adjust the count parameters. As you adjust the count parameters, the count results and the graph automatically update.
4. To close the histogram, press  or .

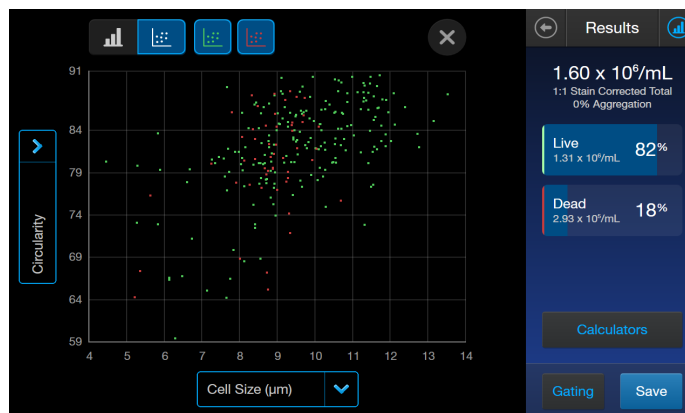
View scatter plot




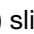

For cell count and cell viability assays performed in the brightfield channel, you can view the distribution of cells (live and/or dead) based on size in a scatter plot.

Note: You can view the scatter plot on the **Results** and **Gating** screens.

From the histogram view:

1. Tap  (white).




2. (Optional): Edit the x- or y-axis variables by tapping the variable arrow on the axis. You can choose between **Circularity** and **Cell Size (µm)** on either axis. Tap **X** to return to the scatter plot.
3. To view the distribution of only the live or dead cells, press the corresponding  **(Live)** (green) or **(Dead)** (red) button.
The plot will automatically update and display the distribution of cells based on size only in the selected population.
4. (Optional): On the **Gating screen**, use the  **(Size)**,  **(Brightness)**, and  **(Circularity)** sliders, adjust the count parameters. As you adjust the count parameters, the count results and the graph automatically update.
5. To close the scatter plot, press  or **X**.

Gate count results

Gating screen

The **Gating** screen for cell count and cell viability assays in the brightfield channel contains the controls for gating results based on size, brightness, and circularity. You can adjust the count parameters before or after performing a count, and save these changes to the current profile or as a separate profile (“Save as new protocol” on page 35).

Gate count results

1. On the **Results** screen, press **Gating** in the bottom right corner to open the **Gating** screen.
2. (Optional) Press  (**Histogram**) to view the distribution of cells (live and/or dead) based on size as you gate the count results. See “Gate count results” on page 35.
3. Select the channel (**Live** or **Dead**) you wish to gate.
4. Select the count parameter (size, brightness, or circularity) to be changed.
5. Using the start and end gate sliders, adjust the count parameters.

Note: For a description of the count parameters and count parameter controls (for example, parameter sliders), see “Count parameters” on page 16.

6. When finished, press **Apply** to set the changes to the parameters and return to the **Results** screen.
7. To save the parameter changes to the current protocol or to create a protocol with the new parameters, see “Create a protocol” on page 18.

Note: You cannot change parameters of preset protocols. If gating is adjusted on Preset protocols, you must save as a new protocol by pressing **Protocols** on the **Gating screen**, and then **Create**.

Save as new protocol

Edit and save a new protocol

If you have made changes to the count parameters after performing a count, you may save your settings to a new protocol.

1. Press **Protocols**.
2. To create a new protocol with the edited parameters, press **Create**. The new protocol screen opens and displays the edited count parameters from the **Gating** screen.
3. To update an existing protocol with the edited parameters, select a protocol from the list, then press **Actions** ▶ **Edit**.

Note: You cannot edit the Default or Preset Protocols.

The **Edit Protocol** screen opens. To import the current settings from the **Gating** screen, press **Import settings**.

4. To change the name of the selected protocol, press the **Protocol name** text box and enter the desired name using the alphanumeric keypad.

5. *(Optional)* If desired, make additional changes to the protocol and the count parameters.
6. Press **Save** to save the new protocol settings and return to the Protocols. Press **Load** to apply the protocol to your current and future counts OR press **Cancel** or ⏪ (**Back**) to return to the **Protocols** screen, then **Cancel** or ⏪ (**Back**) again to return to the **Results** screen without saving the changes to the protocol.

6

Cell counting with fluorescent cells

Count cell fluorescence

Overview

Countess™ 3 FL Automated Cell Counter equipped with the optional EVOS™ light cubes can be used for a variety of fluorescent applications, including simultaneous counts of cells expressing two different fluorescence proteins, stained with fluorescent dyes, and apoptosis and cell viability assays, such as the ReadyCount™ stains.

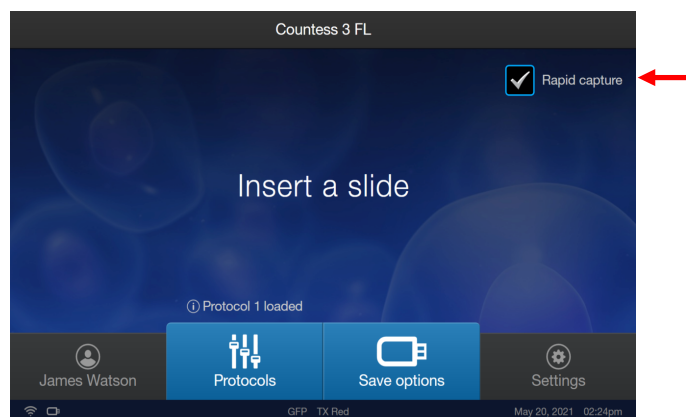
The Countess™ 3 FL Automated Cell Counter can count fluorescent objects in two ways: brightfield-based (BF) or fluorescence-based (FL). After your slide is inserted, you will be given the option of choosing which light cubes will be used. An FL-based count is chosen by checking a box on the **Setup screen**. If the box is not checked, the count will be BF-based. BF-based counting requires a BF image to be captured, while FL-based counting does not. BF-based counts uses the BF image to first identify the cells and determine which of those cells contain fluorescence. FL-based counts uses the fluorescent images to perform the cell count. See specific instructions in Customize capture.

For instructions on installing EVOS™ light cubes to your Countess™ 3 FL Automated Cell Counter, see “Install or change an EVOS™ light cube” on page 64.

Capture and count

Rapid capture

To bypass adjusting individual parameters before the count, you can select **Rapid capture** at the top of the **Home** screen. This will automatically set the parameters and perform a count as soon as the slide is inserted into the instrument.



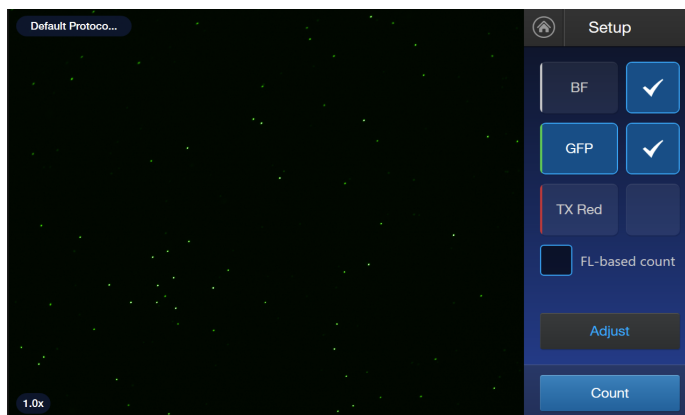
1. Load the samples into the instrument as described in Chapter 4, “Slide preparation and operation”.
2. See “View results” on page 30.

Customize capture



1. Load 10 μL of the fluorescent sample mixture per chamber into the sample slide as described in “Load Countess™ Cell Counting Chamber Slide” on page 24. Let the sample mixture settle for 30 seconds.
2. (Optional): Press **Protocols** to open the **Protocols** screen and load the desired gating parameter as described in “Load a protocol” on page 17.
3. Insert the sample slide into the slide port (“Exterior instrument parts” on page 9), making sure that the sample side is inserted completely into the instrument. You will hear a soft click if the slide is pushed in correctly.
4. When the slide is inserted, the instrument automatically illuminates the sample, sets the intensity of brightfield illumination, and auto focuses on the cells.

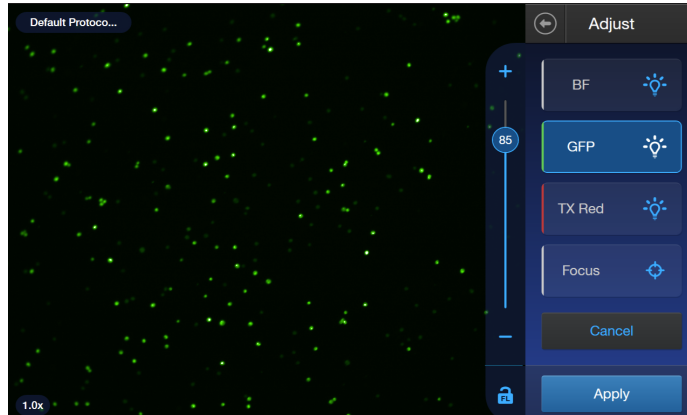
Note: Auto lighting is optimal for producing accurate results, but if needed, can be disabled by adjusting the lighting slider.


5. To view your sample under a different light source, press the desired light source button (**BF** (brightfield) or a source provided from an installed light cube (for example, **GFP**, **DAPI**, **TX Red**, etc.)). The instrument displays the sample in the selected channel (brightfield or fluorescent). In the example below, the sample is displayed in the GFP channel.

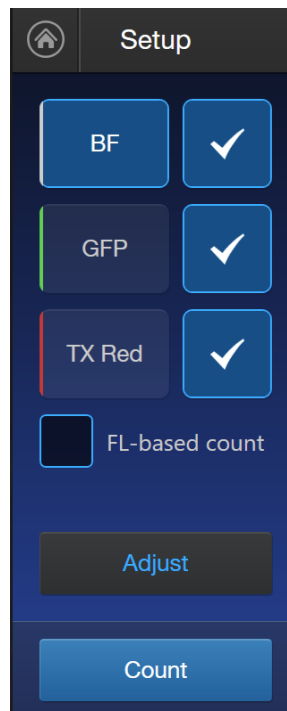


Note: The light source buttons select the light channel (brightfield and/or fluorescence) for sample illumination and are used when setting the LED intensity for the selected channel (step 9–11); they do not determine which channels are used for capturing the image. The check boxes determine which channels are used for capturing the image (step 8).

6. (Optional) To set the light intensity, press **Adjust** to go to the **Adjust** screen. Using auto lighting for each sample, the Countess™ 3 FL will find the optimal FL lighting. To use the same FL lighting for subsequent samples, press **Adjust**, then press  (fluorescence lock) to lock the FL lighting . Notification of locked status will appear at the upper portion of the screen.



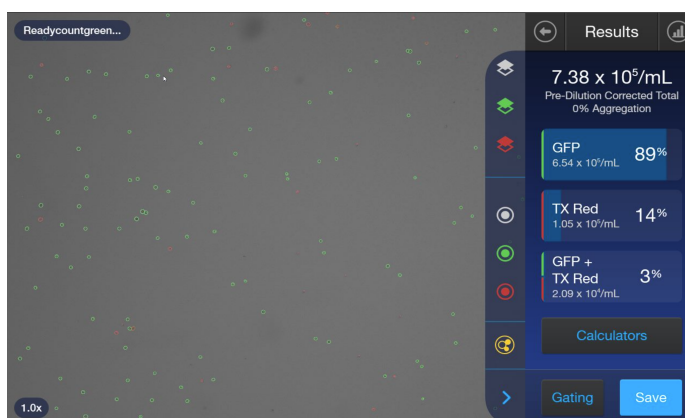
7. Press the light source button for the channel you wish to set the light intensity and adjust the light intensity using the light source slider. If desired, repeat the procedure for the remaining channels.
 - a. Press **Apply** to keep the current setting and return to the **Setup** screen OR press **Cancel** or  (**Back**) to return to the **Setup** screen without saving changes.
8. On the **Setup** screen, select the check boxes to the right of the channels you wish to capture. **BF** (brightfield) is checked by default and cannot be deselected for BF-based count. For FL-based count, select the **FL-based count** box. Doing so will allow you to deselect the BF box and generate only the FL-based count.



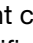



9. Press **Count**.

The instrument captures the images and displays the results. The results include total concentration, percentage of fluorescent cells for each channel, and the percentage of aggregated cells. For more information about the differences between BF and FL-based count results, see “View results” on page 41.

Autosaving will occur when the count is completed if you have **Autosave** enabled from the **Save Options** screen.



Next steps

- To identify all cells from the brightfield image in BF-based counts, press the appropriate colored icon for the specific EVOS™ light cube (for example, press  for GFP, press  for TX Red). Brightfield objects are not identified during FL-based counts. To identify the brightfield objects (cells) in BF-based counts, press .
- To see the distribution of cells counted through each channel in a graphical format, press  (**Histogram**). See “View histogram for cell fluorescence assays” on page 43.

Note: Changes made in the current protocol to size, brightness, circularity, and relative fluorescence intensity can be saved directly from the **Gating** screen.

- To calculate the volume of cell sample and buffer needed to reach a desired concentration based on the count results, press **Calculators** to open the Calculator application. See Chapter 7, “Calculators”.
- To permanently save the results, press **Save**. See “Save screen” on page 52.
- To perform a new count, push the slide to eject, then insert a new sample slide.
- To gate the results by object size, brightness, circularity, or relative fluorescence intensity, press **Gating**. See “Gating screen” on page 45.

View results

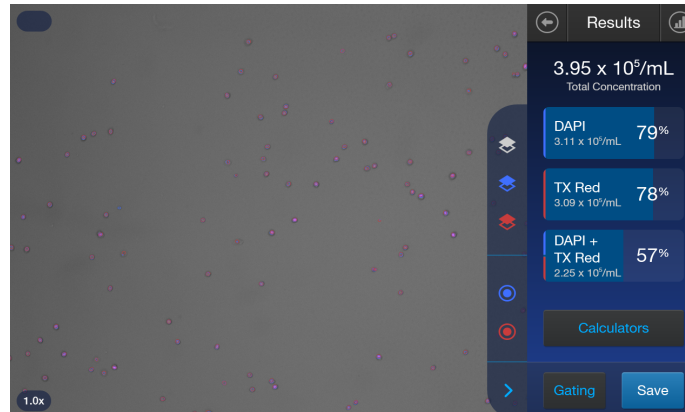
Results screen for BF or FL-based counts

The **Results** screen fluorescent cell counts displays a composite image of the channels captured and the objects counted which is determined by the count mode selected (BF or FL-based). BF-based counts uses the BF image to identify the total number of cells and will report the percentage of which contain fluorescence. FL-based counts uses the fluorescent images to count all fluorescence and will not report a number of non fluorescence cells.

Results screen for FL-based counts

The **Results** screen for FL-based counts displays a composite image of the total concentration and the percentage of fluorescent cells for each channel imaged and of the objects counted. The results shown include:

- Total concentration: The overall concentration of fluorescent cells in the sample.
- Percentage and concentration of Cube 1 positive cells: The number and percentage of cells in the sample that is fluorescent for using Cube 1 (GFP in this example).
- Percentage and concentration of Cube 2 positive cells: The number and percentage of cells in the sample that is fluorescent for using Cube 2 (TX Red in this example).
- Percentage and concentration of double positive cells: The number and percentage of cells in the sample that is fluorescent for using Cube 1 and Cube 2.

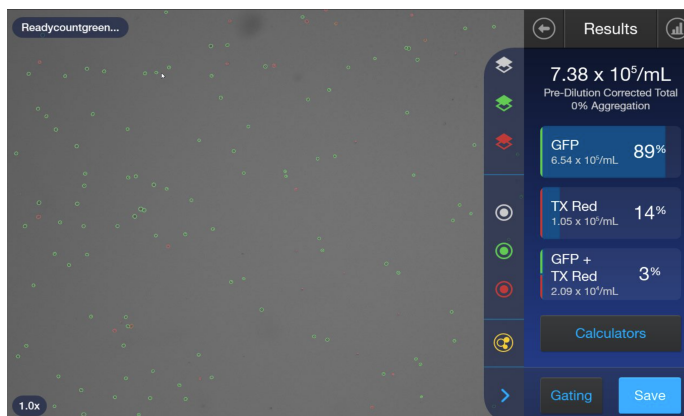


Note: The FL-based counts are based on fluorescence localization, which may not be similar sizes to cell membranes seen in BF images, resulting in fewer aggregate identifications in when in FL-based counting mode.

Results screen for BF-based counts

The **Results** screen for BF-based counts displays a composite image of the channels imaged and of the objects counted. The results shown include:

- Total concentration: The overall concentration of cells in the sample.
- Percentage and concentration of Cube 1 positive cells: The number and percentage of cells in the sample that is fluorescent for using Cube 1 (GFP in this example).
- Percentage and concentration of Cube 2 positive cells: The number and percentage of cells in the sample that is fluorescent for using Cube 2 (TX Red in this example).
- Percentage and concentration of double positive cells: The number and percentage of cells in the sample that is fluorescent for using Cube 1 and Cube 2.
- Percentage of aggregates: The percentage of cell clusters (aggregates) consisting of three or more cells with touching membranes as seen in the BF image.





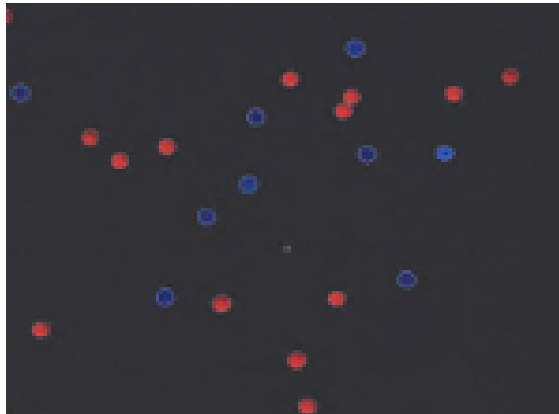
Note: When using the default protocol, the total cell concentration displayed after a fluorescent count does not take any dilution into account. If the sample was diluted with a stain prior to the count, the dilution factor must be added to calculate the original cell concentration in the sample slide. If using ReadyCount™ stains, use the ReadyCount™ Preset protocols which account for the 1:1 stain dilution used prior to count. See Chapter 7, “Calculators” for more details.

Note: The percentages will not add up to 100%. Cube 1 and Cube 2 identifications report all positive cells, single and double positive cells.

Identify objects counted

Identify cells counted in fluorescence assays

1. On the **Results** screen, select the appropriate icon/s to turn on or off the layer/s viewed.
For example, select .
2. To identify the cells that are counted in a specific channel, press the corresponding contours icon.
Cells counted in the selected channel will be circled on the screen with the same color as the selected channel.
For example, select .
In the example below, both the DAPI and TX Red boundary buttons are selected, and the cells counted in the DAPI and TX Red channels are marked with blue and red circles, respectively.




3. To unmark the cells counted in a specific channel, press the corresponding boundaries button again.

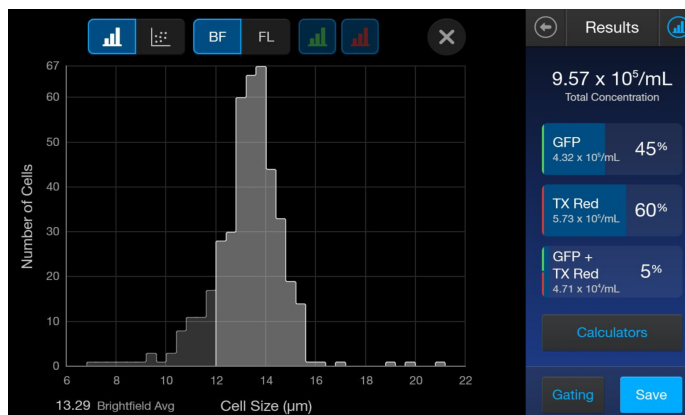
Histogram count results


View histogram for cell fluorescence assays

For fluorescence assays, you have the option of viewing the distribution of the cells based on size or based on relative fluorescence intensity in a graphical format.

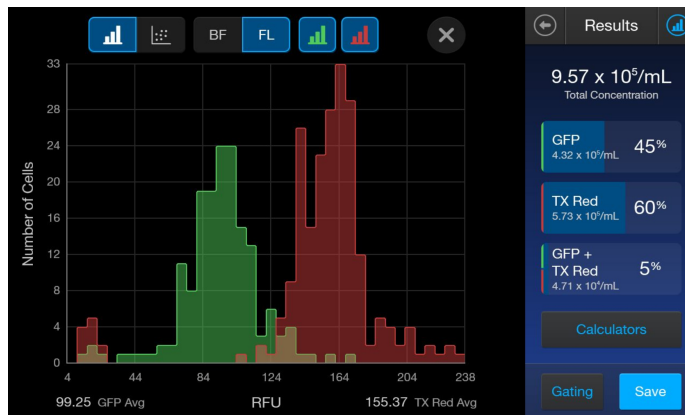
Note: You can view the histogram on the **Results** and **Gating** screens.



1. To view the histogram showing the distribution of cells based on size, press  (**Histogram**) and then select **BF** (brightfield).
The histogram displays the size distribution of the total cell count (number of cells vs. cell size in μm) and the average size of the cells counted in each available fluorescence channel.



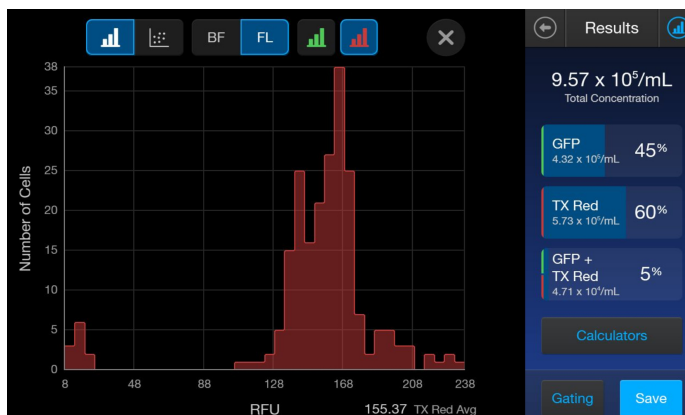
- To view the distribution of cells based on relative fluorescence intensity, press  (**Histogram**), then select **FL** (fluorescence).


The histogram displays the distribution of cells based on fluorescence intensity.



- To remove the cells counted in a specific channel from the histogram, press the corresponding channel button   on the histogram.

The histogram automatically updates and displays the distribution of cells based on relative fluorescence intensity only in the selected (for example, checked) channel.



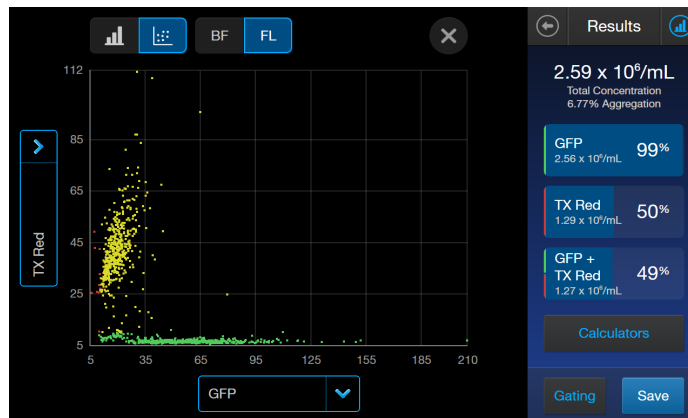
- To add the cells counted in a specific channel to the histogram, press the corresponding channel button.
- To close the histogram, press  (**Histogram**) or **X (Close)**.

View scatter plot


For cell count and cell viability assays performed with fluorescence, you can view the distribution of cells based on size, circularity and fluorescent expression in a scatter plot.

From the histogram view:

- Tap  (white).



Note: Double positive cells appear as yellow in the example although single positive cells appear as red and green respectively.

- (Optional): Edit the x- or y-axis variables by tapping the variable arrow on the axis. You can choose between circularity, cell size (μm), cube 1 intensity, cube 2 intensity on either axis after a BF-based count. After a FL-based you can choose between circularity, cell size (μm), Intensity, for each fluorescent cube. Tap to return to the scatter plot.
- (Optional): On the **Gating** screen, use the (size), (brightness), and (circularity) sliders, to adjust the count parameters. As you adjust the count parameters, the count results and the graph will update automatically.
- To close the scatter plot, press  (**Histogram**) or **X (Close)**.

Gate count results


Gating screen

The **Gating** screen for cell fluorescence assays contains the controls for gating count results based on size, brightness, circularity, and fluorescence intensity. You can adjust the count parameters before or after performing a count, and save these changes to the current protocol or as a separate protocol ("Save as new protocol" on page 35).


Gate count results

1. On the **Results** screen, press **Gating** to open the count parameters screen containing the controls for adjusting the count parameters in the selected channel.

Note: For a description of the count parameters and count parameter controls (for example, parameter sliders) see “Count parameters” on page 16.

2. (Optional): Press  (**Histogram**) to view the distribution of cells based on size or fluorescence intensity. See “View results” on page 41.
3. Press **BF** (brightfield) to adjust the parameters for size, brightness, and circularity using the gating sliders.
4. Select the desired fluorescence channel (for example, DAPI, TX Red, etc.) to adjust the threshold for fluorescence intensity using the fluorescence intensity slider.

Note: The fluorescence channels available depend on the EVOS™ light cubes installed in the instrument.

5. When finished, press **Apply** to save the changes to count parameters and return to the **Results** screen OR press **Cancel** or  (**Back**) to return to the **Results** screen without saving the changes.
6. To save the parameter changes to the current protocol or to create a protocol using the new count parameter, see “Edit and save a new protocol” on page 46.

Note: You cannot edit change parameters of Preset Protocols. If **Gating** is adjusted on the Preset Protocols, you must save as a new protocol by pressing **Protocols** on the **Gating** screen and then **Create**.

Save as new protocol

Edit and save a new protocol

If you have made changes to the count parameters after performing a count, you may save your settings to a new protocol.

1. Press **Protocols**.
2. To create a new protocol with the edited parameters, press **Create**. The new protocol screen opens and displays the edited count parameters from the **Gating** screen.
3. To update an existing protocol with the edited parameters, select a protocol from the list, then press **Actions** ▶ **Edit**.

Note: You cannot edit the Default or Preset protocol.

The **Edit Protocol** screen opens. To import the current settings from the **Gating** screen, press **Import settings**.

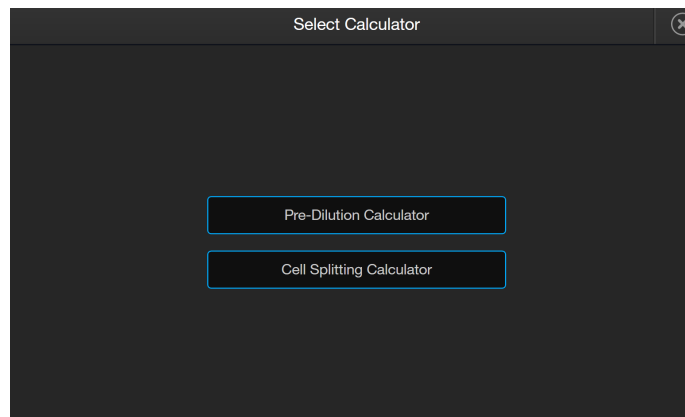
4. To change the name of the selected protocol, press the **Protocol name** text box and enter the desired name using the alphanumeric keypad.
5. *(Optional)* If desired, make additional changes to the protocol and the count parameters.
6. Press **Save** to save the new protocol settings and return to the **Protocols** screen. Press **Load** to apply the protocol to your current and future counts or press **Cancel** or ⏪ (**Back**) to return to the **Protocols** screen, then **Cancel** or ⏪ (**Back**) again to return to the **Results** screen without saving the changes to the protocol.

Pre-Dilution Calculator

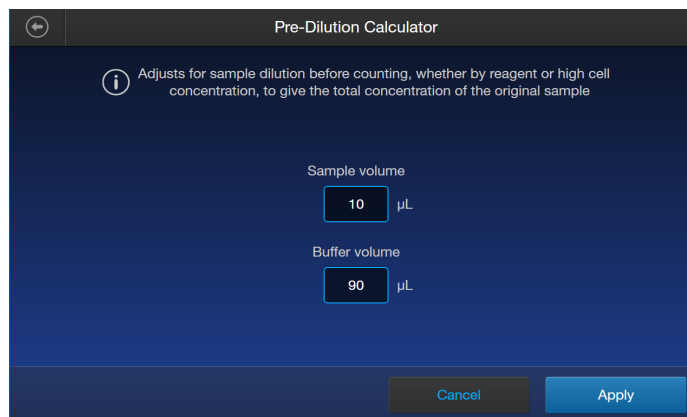
The **Pre-Dilution Calculator** is used to account for diluting a very concentrated sample before loading it into a sample slide for counting.

For a brightfield count, this calculator also has an option to account for Trypan Blue use.

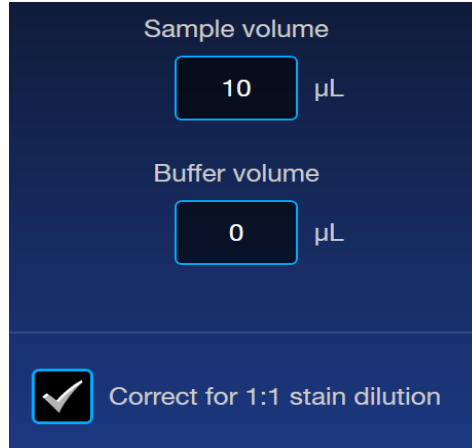
1. On the **Results** screen, press **Calculators** then select **Pre-Dilution Calculator**.



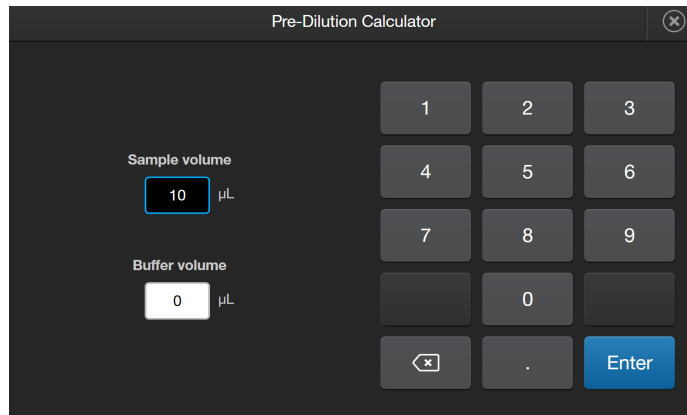
2. Click-in one of the text fields.



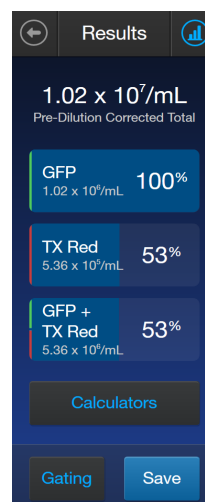
Note: You can select the checkbox at the bottom of the screen to correct for a 1:1 stain dilution.



- Using the keypad, enter the appropriate volumes and concentrations needed for your experiment. Press **Enter** after filling in both text boxes.



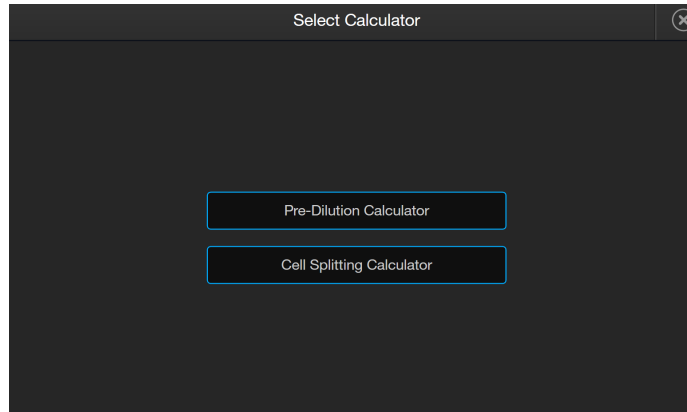
- Press **Calculate**. The result will appear at the top of the **Results** screen.



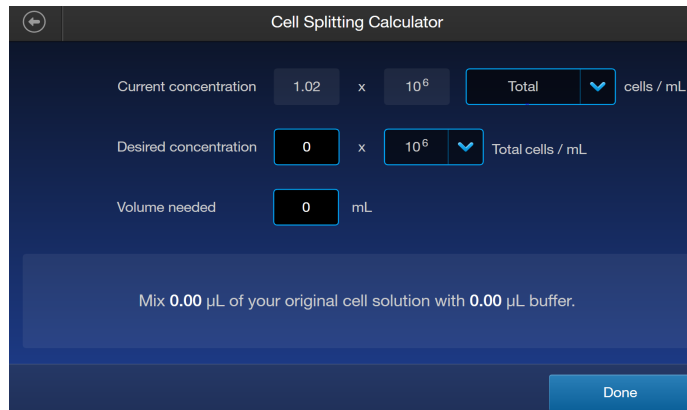
Cell Splitting Calculator

The **Cell Splitting Calculator** is used to determine the amounts of sample and buffer needed to reach a desired concentration.

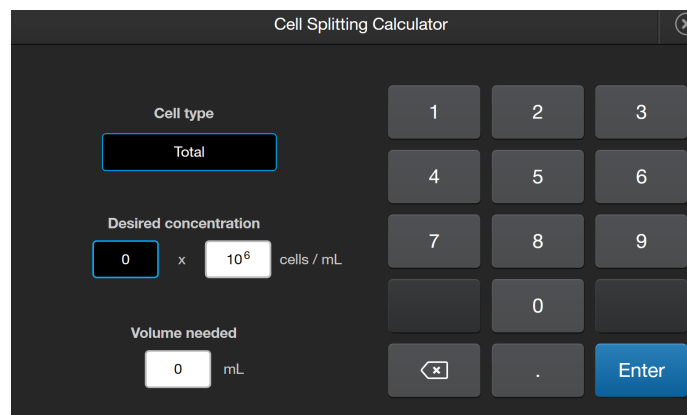
1. On the **Results** screen, press **Calculators** then select **Cell Splitting Calculator**.



2. Click-in one of the text fields.



3. Using the keypad, enter the appropriate volumes and concentrations needed for your experiment. Press **Enter** after filling in the text boxes.



4. The calculation will automatically update below the text fields. When completed, press **Done**.

The screenshot shows the 'Cell Splitting Calculator' interface. It features three input rows: 'Current concentration' with a value of 1.02 and a multiplier of 10^7 ; 'Desired concentration' with a value of 2 and a multiplier of 10^5 ; and 'Volume needed' with a value of 30. Below these inputs, a text box displays the calculated result: 'Mix 589.63 μ L of your original cell solution with 29.41 mL buffer.' A 'Done' button is located at the bottom right of the calculator.



Save results

Save screen

The Countess™ 3 and 3 FL cell counters allow you to save your data and images using a USB flash drive, a link to a Thermo Fisher Connect account, or a link to a network drive.

To save your experiment, choose from the following options, in any combination:

- **Result:** Saves the **Results** screen as it is displayed on the instrument, with or without the graph, in the selected image format (JPEG, PNG, or TIFF).
- **Images:** Saves the raw captured image as well as underlying channel images in the selected image format (JPEG, PNG, or TIFF).
- **Report:** Saves a printer-friendly report of the results, graph(s), and image in PDF. For more information, see “Report file” on page 55.
- **CSV:** Saves the data from current and previous counts as a CSV file (comma separated values). The CSV format allows for processing or re-displaying results with any third party software or spreadsheet program. For more information on the CSV file format, see Appendix E, “CSV file format definition” Appendix E: CSV file format.

Note: A dropdown list will provide options for how many counts to save. You can choose from **Current Count**, **Session**, **Today**, **Yesterday**, **Previous 7 days**, **Previous 30 days**, or **All**.

- **Cell Data:** Saves single cell data from the count in either a FCS or CSV format. FCS is a Flow Cytometry Data File Standard and allows for processing or re-displaying results. The CSV format allows for processing or re-displaying results with any third party software or spreadsheet program, providing each cells data on a separate row.

Note: If you wish to save your results with the graph showing the distribution of cells based on cell size or fluorescence intensity, make sure that the desired graph is displayed on the **Results** screen.

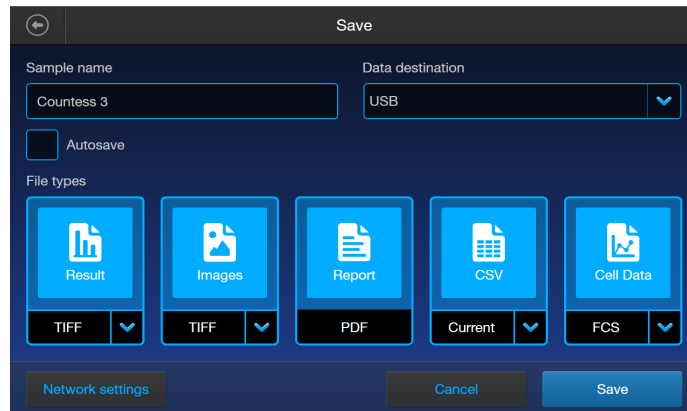
Save procedure

1. To save your data, insert the Countess™ USB drive (or equivalent) into an available USB port on the instrument or confirm the instrument is connected to a Thermo Fisher Connect account or network drive.

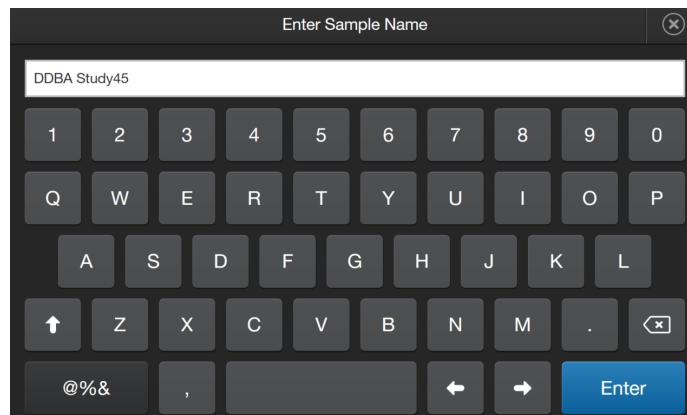
Note: If the instrument is not connected to a network, then use **Network settings** at the bottom of the screen to reconnect.

Note: The USB ports located in the front and the back of the instrument function the same. However, the first USB drive connected will be the preferred saving location.

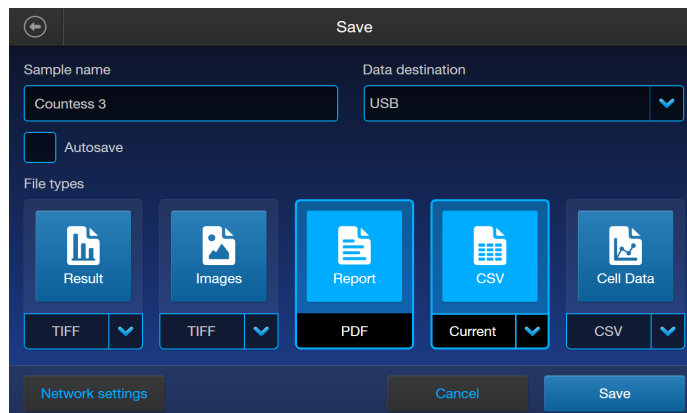
2. On the **Results** screen, press **Save** to view the **Save** screen.



3. To assign a name to your count, press the **Sample name** text field and enter the name. To enter symbols, press the **@%&** key. To return to the keypad, tap the **ABC** key.



4. Press **Enter** to save the name and return to the **Save** screen. To return to the **Save** screen without saving the name, press **X (Close)**.
5. Select the desired file type to save your experiment (**Result**, **Images**, **CSV**, **Report**, **Cell Data**). You can select an individual mode (e.g., **Result only**) or any combination of modes (e.g., **Result**, **Images**, **Data**, and/or **Report**). In the example below, **Report** and **CSV** are selected.



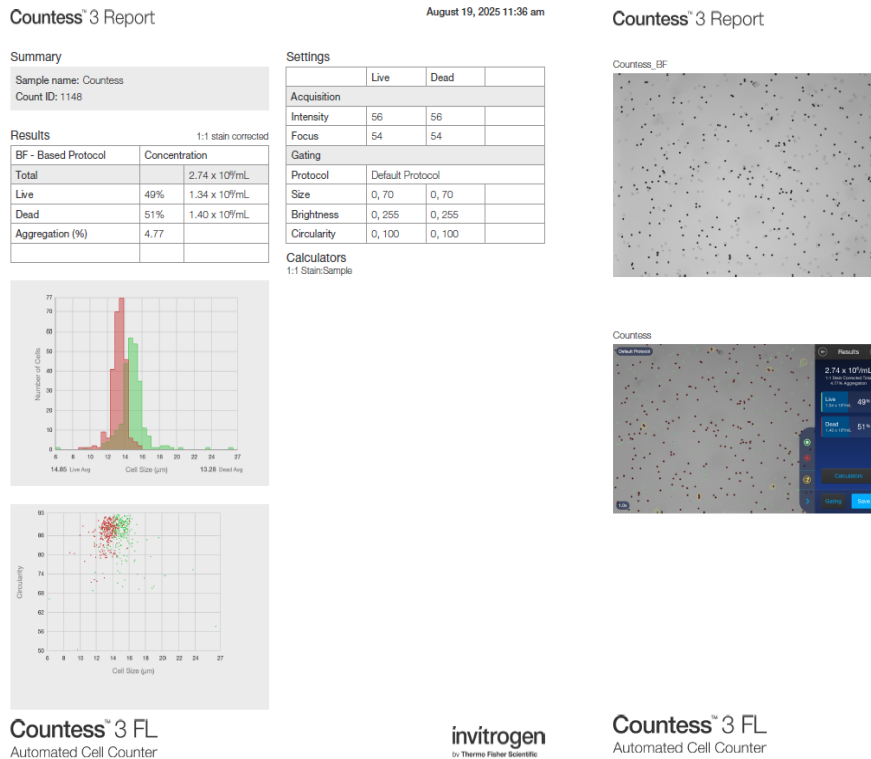
- By default, **Result** and **Images** are saved as TIFF files, and **CSV** uses the current capture data.
- To choose a different file format, press the desired file type. Available options are **JPEG**, **PNG**, and **TIFF**. For CSV files, you can choose from the selections of **Current Count**, **Session**, **Today**, **Yesterday**, **Previous 7 days**, **Previous 30 days**, or **All**.
After you make your selection, the instrument returns to the **Save** screen.
To return to the **Save** screen without changing the file format, press **X (Close)**.
- Press **Save** to save your experiment to the selected mode(s) in the USB drive, Connect account, or network drive.
- Press **Done** and then transfer the USB drive to the desired location.

Report

Report file

The **Report** function allows you to save a printer-friendly report of the results, graphs, and images in PDF format.

Report from brightfield count



- The top section of the report contains a table with the results as displayed on the **Results** screen, showing the concentration of the sample and the percentage and number for the total, Live, and Dead channels, along with the percentage of Aggregates. The report also displays the template information used to gate these images.
- Below the results table, the report contains the “Number of Cells (µm) vs. Cell Size” graph and “Cell Size (µm) vs. Circularity” scatter plot.
- Under the graph, the report contains the brightfield count image, with the live and dead cells identified by the green and red circles, respectively.

Report from fluorescence count

Countess³ Report

September 04, 2025 11:12 am

Summary

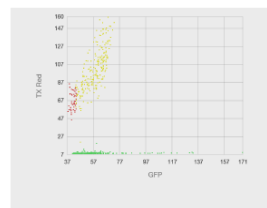
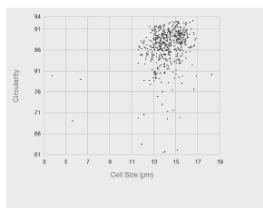
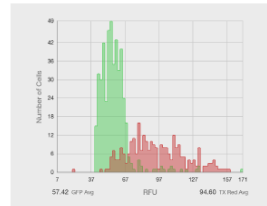
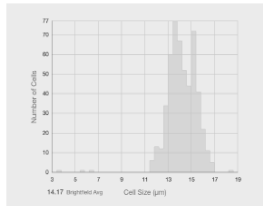
Sample name: Bboth
Count ID: 1200

Results

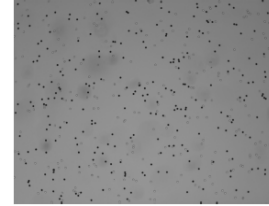
BF - Based Protocol	Concentration
Total	1.36 x 10 ⁹ /mL
GFP	88% 1.20 x 10 ⁹ /mL
TX Red	52% 7.09 x 10 ⁸ /mL
GFP + TX Red	43% 5.83 x 10 ⁸ /mL
Aggregation (%)	4.23

Settings

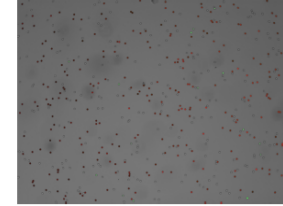
	BF	GFP	TX Red
Acquisition			
Intensity	56	77	43
Focus	53		
Gating			
Protocol	Default Protocol*		
Size	0, 70		
Brightness	0, 255	0, 255	0, 255
Circularity	0, 100		

Countess³ FL
Automated Cell Counterinvitrogen
By Thermo Fisher ScientificCountess³ Report

Bboth_BF



Bboth_Composite



Bboth



* represents unsaved changes.


Countess³ FL
Automated Cell Counterinvitrogen
By Thermo Fisher Scientific

- The top section of the report contains a table with the results as displayed on the **Results** screen, showing the concentration of the sample, and the percentage and number of cells for the total, FL1, FL2, and FL1 + FL2 channels. The report also displays the profile information used to gate these images.
- Below the results table, the report contains the “Number of Cells vs. Cell Size” graph on the left, and “Number of Cells vs. RFU (relative fluorescence units)” graph on the right. Below the bar graphs are the scatter plots showing "Circularity vs. Cell Size (µm)" and "TX Red vs. GFP".
- Under the graphs, the report contains the count images, with the brightfield image on the left and the fluorescence images on the right.



Settings overview



Settings screen

To access the **Settings** screen, press  (**Settings**) on the **Home** page.


Settings allows you to:

- Perform software updates (see “Update the Countess™ 3/3 FL Cell Counter software” on page 58)
- Set the date and time (see “Set the date and time” on page 59)
- Change or install EVOS™ light cubes (see “Install or change an EVOS™ light cube” on page 64)

Connect to a wireless network

Network connecting options	
<ol style="list-style-type: none">1. From the Home screen, press  (Settings).2. Press Instrument settings.3. Press Wireless network. A list of available networks will display.4. Select the desired network and press Connect.	<ol style="list-style-type: none">1. From the Home screen, press  (Save options).2. Press Network settings.3. Select the desired network and press Connect.
Follow either set of steps to also disconnect the current network and connect to a new one.	

Connect to a network drive

1. From the **Home** screen, press  (**Settings**).
2. Press **Instrument settings**.
3. Press **Network drive**.
4. Press **Drive location**. Enter the IP address for your shared folder.
5. (*Optional*): Enter any additional information, including a **Domain name**, **Username**, or **Password**.

Note: Consult with your local IT representative for questions regarding drive location and domain name.

6. Press **Connect**.

The network drive is now connected. If needed, you can edit the **Drive location**, **Domain name**, or **Username** by pressing **Edit**.

Software update

Guidelines for software update

- Software updates can be done via your Connect account. When connected to the internet, the system will automatically check for and download software updates to the instrument.

Note: Using the Cloud will require a Wi-Fi dongle be installed in the instrument.

- The USB drive used for transferring the software update file must be FAT32 formatted; verify this before proceeding. If necessary, reformat the USB drive to FAT32 following the recommended procedure for your operating system.

Note: Reformatting the USB drive will result in the loss of all files. Back up the files in the USB drive prior to reformatting.

- The software update file must be saved on the top level of the USB drive, not within a folder or a subfolder.
- The software update file must be uncorrupted during transfer. Do not rename, zip, or compress the software update file.

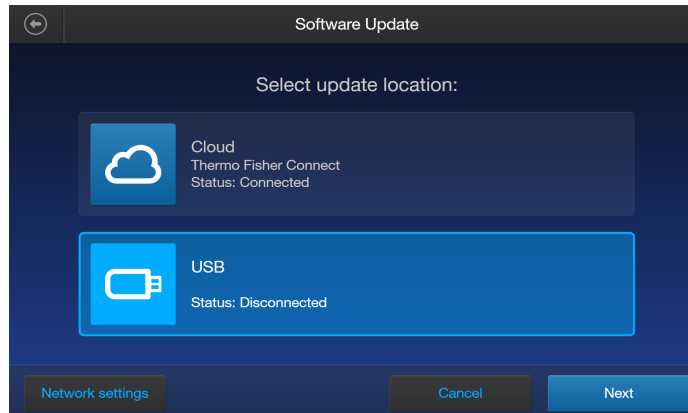
Update the Countess™ 3/3 FL Cell Counter software

1. Go to www.thermofisher.com/countessupdate, and download the latest Countess™ 3/3 FL cell counter software version to your desktop.

Note: The software update file has a version-specific name followed by the extension.exe (For example, Countess_xxxx for software version 2.0.2).

2. Copy the software update file onto the USB drive, making sure that it is saved on the top level and not hidden within a folder.
3. Insert the USB drive into one of the USB ports of the instrument.
4. Press **⊗ (Settings)** on the **Home** page to open the **Settings** screen.

5. Press **Software Update**.



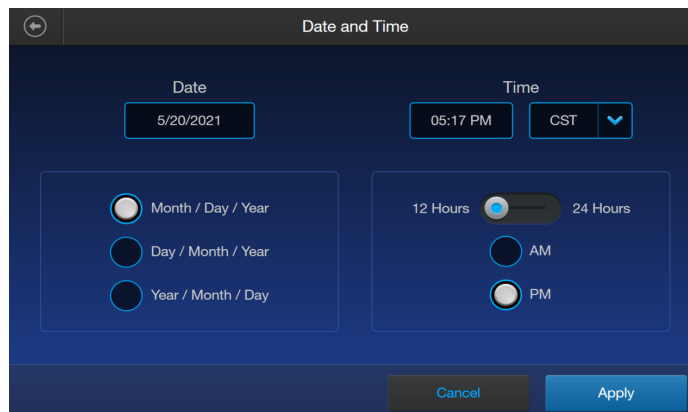
6. Select the update location. Press **Next**.
If the **Cloud** option is disconnected, press **Network settings** to proceed to reconnect to your network.

Note: Using the Cloud will require a Wi-Fi dongle be installed in the instrument.

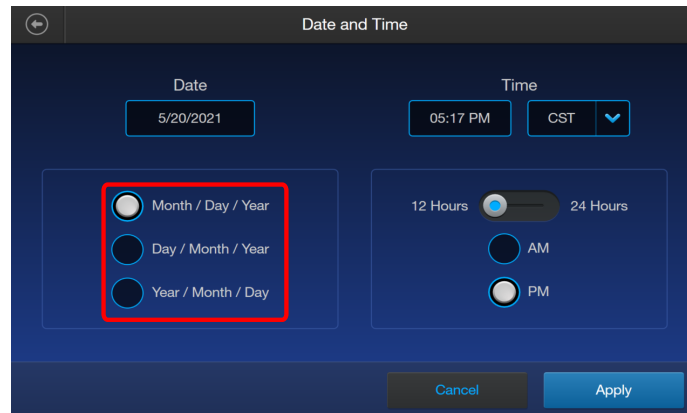
7. If an update is available, press **Update**. Once the update has completed, the **Home** screen will appear to resume normal operation.

Set the date and time

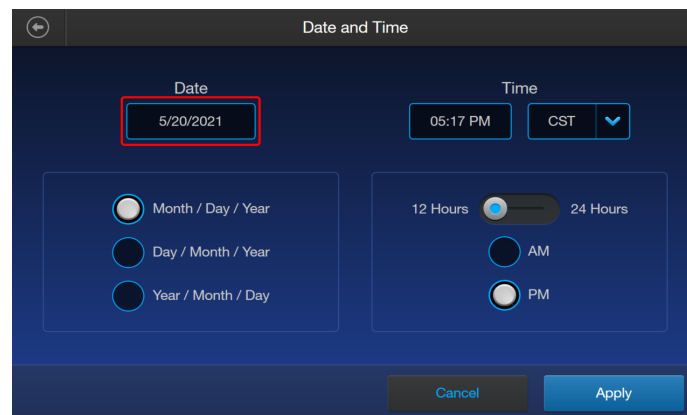
1. Press **Settings** on the **Home** page.
2. Press **Instrument Settings**.
3. Press **Date and Time**.



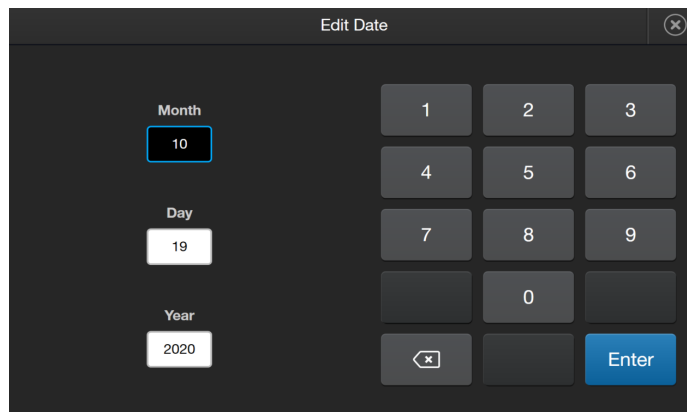
4. Select the date format (**Month/Day/Year**, **Day/Month/Year**, or **Year/Month/Day**).



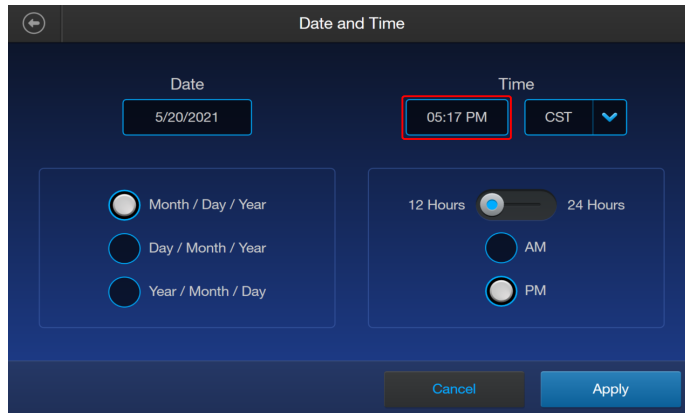
5. Press inside the **Date** text box to open the **Edit Date** keypad.



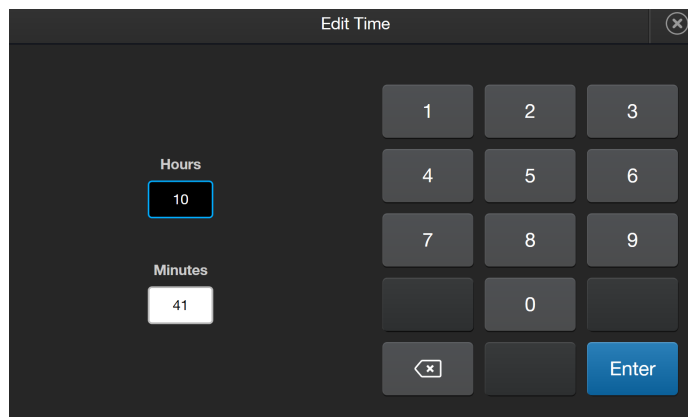
6. Using the keypad, enter the date into the **Month**, **Day**, and **Year** text boxes. Press **Enter**.



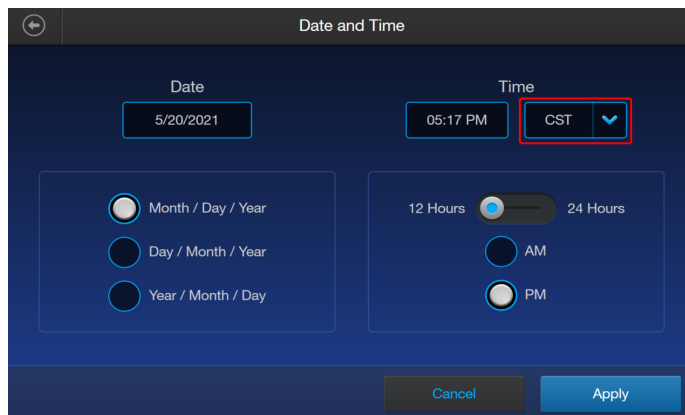
7. Press the **Time** text box to open the **Edit Time** keypad.

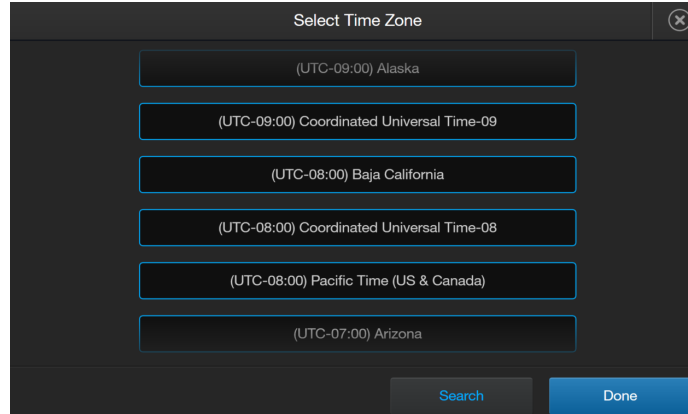


8. Using the keypad, enter the time into the **Hours** and **Minutes** text boxes.

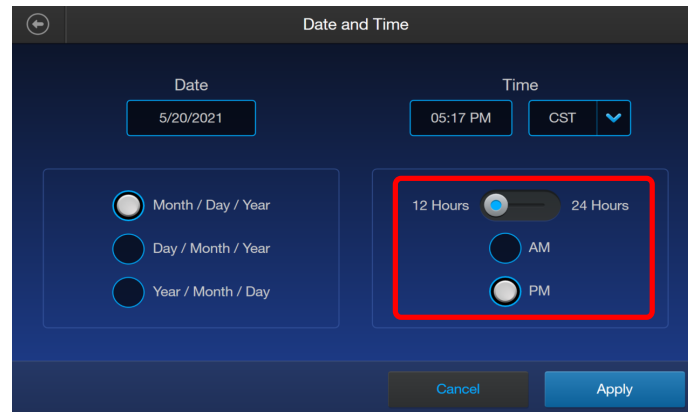


9. Press the drop-down menu under Time to select your time zone from the list. If needed, press **Search** to use a keyboard to type in your search term (e.g., pst, cst, etc.).





10. Select the time format you wish to use (**12 Hours** with **AM** or **PM** or **24 Hours**).

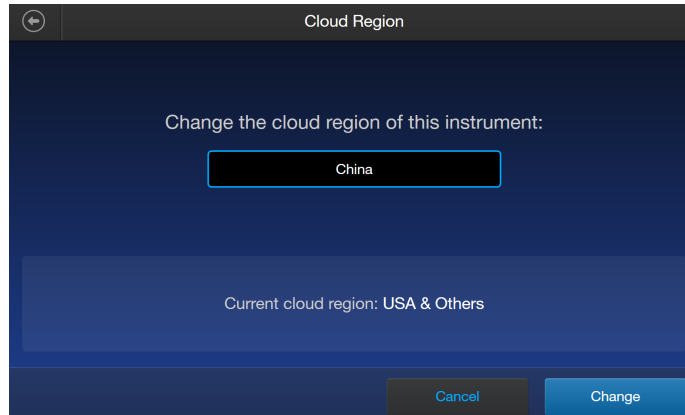


11. Press **Apply** to set the Date and Time and return to the **Instrument Settings** screen.
Press **Cancel** or ⬅️ (**Back**) to return to the **Instrument Settings** screen without saving your changes.

Change your cloud region

1. Press ⚙️ **Settings** on the **Home** page.
2. Press **Instrument Settings**.

3. Press **Cloud region**. Your current cloud region is displayed at the bottom of the screen.



4. Select the desired cloud region from the available choices. Press **Change**.
5. Once the change is complete, press **Done**.

Note: Changing the cloud region will require relinking of all instrument accounts.

Install or change an EVOS™ light cube

The Countess™ 3 FL Automated Cell Counter can accommodate up to two EVOS™ light cubes. Each user-interchangeable, auto-configured EVOS™ light cube contains an LED, collimating optics, and filters for fluorescence applications. EVOS™ light cubes do not come standard with the device and must be purchased separately (“EVOS™ light cubes” on page 87). To install or change a light cube:

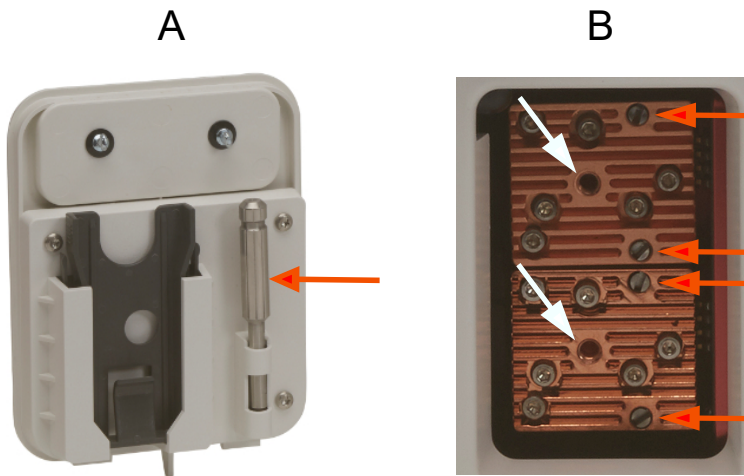
1. Press **⊙ Settings** on the **Home** page.
2. Press **Instrument Settings**.
3. Press **Change light cube**. The instrument positions the light cube tray to enable light cube installation.
4. When prompted, power off the Countess™ 3 FL Automated Cell Counter using the power switch on the back of the instrument (“Exterior instrument parts” on page 9).
5. Unplug the power cord from the Countess™ 3 FL Automated Cell Counter.
6. Unlatch the back panel with the two captive ¼-turn fasteners (indicated by black arrows below) that secure the back panel on the rear of the Countess™ 3 FL Automated Cell Counter and remove the back panel.



7. Place the light cube into one of the empty slots in the back of the device.



8. Using the tool provided on the inside of the back panel (red arrow in Figure A), secure the light cube by tightening the two screws on the end of the cube (red arrows in Figure B).



Note: Figure B shows two light cubes installed.

9. To remove a light cube, unscrew both screws that secure it to the instrument.
10. Thread the light cube removal tool into the central hole in the cube (white arrows in Figure B) and gently pull the light cube out of the device.

Note: Always store the cube removal tool in the back panel for easy access.

11. Install the back panel and secure it in its place with both ¼-turn fasteners.
12. Plug the power cord back into the Countess™ 3 FL Automated Cell Counter.
13. Turn on the Countess™ 3 FL Automated Cell Counter by flipping the power switch on the back of the instrument to the ON position.



Instrument care

General guidelines for care

- Use the appropriate cleaning solutions for each component, as indicated in the cleaning procedures in “Clean the cell counter” on page 67.
- If liquid spills on the instrument, turn off the power immediately and wipe dry.

Power supply

Always use the correct power supply. The power adapter specifications appear on the serial number label (bottom of the instrument) and in the "Technical specifications" section of this user guide (“Technical specifications” on page 85). Damage due to an incompatible power adapter is not covered by warranty.



CAUTION! Never disassemble or service the instrument yourself. Do not remove any covers or parts that require the use of a tool to obtain access to moving parts. Operators must be trained before being allowed to perform the hazardous operation. Unauthorized repairs may damage the instrument or alter its functionality, which may void your warranty. Contact your local distributor to arrange for service.

IMPORTANT! If you have any doubt about the compatibility of decontamination or cleaning agents with parts of the equipment or with material contained in it, contact Technical Support or your local distributor for information.

Clean the cell counter

Introduction

Clean the Countess™ Automated Cell Counter periodically to prevent buildup of dust and dirt that might reduce its performance and cause contamination.



CAUTION! To avoid electrical shock, always turn OFF the Countess™ Automated Cell Counter and unplug the power cord before cleaning or decontaminating the instrument.



CAUTION! All biological samples and materials that come into contact with them have the potential to transmit infectious diseases and are considered biohazardous. Follow all applicable local, state/provincial, and/or national regulations. Wear appropriate protective eyewear, clothing, and gloves.

IMPORTANT! Using a cleaning or decontaminating method other than that specified by the manufacturer may result in damage to the instrument.

Clean the touchscreen

- Wipe the touchscreen of the Countess™ Automated Cell Counter using a soft, lint-free cloth moistened with an LCD cleaning solution. Do not apply excessive force during cleaning. Wipe the touchscreen dry immediately after cleaning.
- Ensure that the cleaning solution does not enter the power button, the power inlet, the slide port, or the USB ports.
- Never pour or spray any liquids directly on the instrument to avoid electrical shock when the instrument is plugged in.
- Do not use abrasive cleaning solutions or material to prevent the touch-screen from getting scratched.

Clean the instrument case

- Wipe the instrument case of the Countess™ Automated Cell Counter using a soft, lint-free cloth moistened with distilled water. Wipe the instrument dry immediately after cleaning.
- Ensure that water or other cleaning solutions do not enter the power button, the power inlet, the slide port, or the USB ports.
- Never pour or spray any liquids directly on the instrument to avoid electrical shock when the instrument is plugged in.

Decontaminate the instrument

- Wipe the instrument case of the Countess™ Automated Cell Counter using a soft, lint-free cloth moistened with 70% alcohol. Wipe the instrument dry immediately after cleaning.
- Avoid using a bleach solution, because it may leave a residue of bleach crystals on the instrument.
- Ensure that water or other cleaning solutions do not enter the power button, the power inlet, the slide port, or the USB ports.
- Never pour or spray any liquids directly on the instrument to avoid electrical shock when the instrument is plugged in.

The installation and/or operational qualification ensures that the instrument is installed and performing within the manufacturer specifications. IQ/OQ requires a thumb drive and the Countess™ 3 Standard Slide (Cat. no. [A51876](#)) shipped with user executed IQ/OQ (Cat. No. [A59080](#)) and user executed OQ (Cat. No. [A54115](#)) purchase.

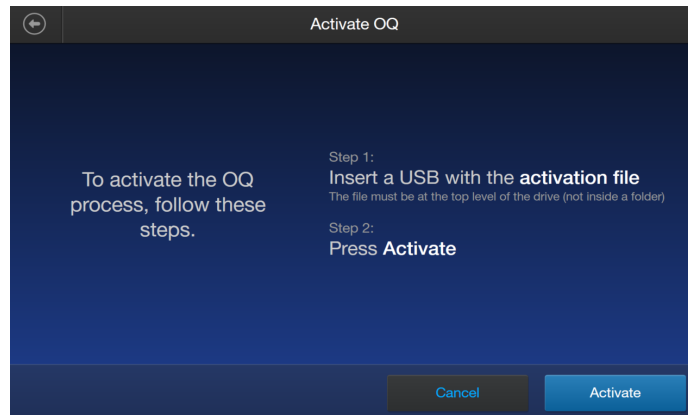
Generate and activate a license

1. To generate the license file, go to apps.thermofisher.com.
2. Sign in to your Thermo Fisher™ Connect Platform account.
3. Open the **Countess™ Image Analysis** application.

Note: If the Countess™ application is not displayed on the **Dashboard** page, click **View all apps** to see a full list.

4. On the **File Gallery** screen, click on **Execute IQOQ** in the top right corner. Then, select **IQ&OQ service** if completing both IQ and OQ or **OQ service** if completing just OQ. See image on page 54.
5. Enter the **Order ID (sales order number)** and **Countess 3 Serial Number** (found on your instrument). If completing IQ, please navigate to page 54 to **Execute the Installation Qualification**. If only completing OQ, continue to follow the instructions below.
6. Click **Get activation file**.
A **Request Submitted** message will display when the order information is validated.
7. Once the file is ready, click on the **Download is Ready** window and select **Download.zip** to begin the license key file download.
8. Save the license file to a USB drive. Do not change the license name "Prod_xxxx.OQ-license" where xxxx is the instrument serial number.
9. Insert the USB drive into the instrument.
10. On the instrument, press **Settings** on the home screen.
11. Press **Execute OQ**.

12. Press **Activate**.

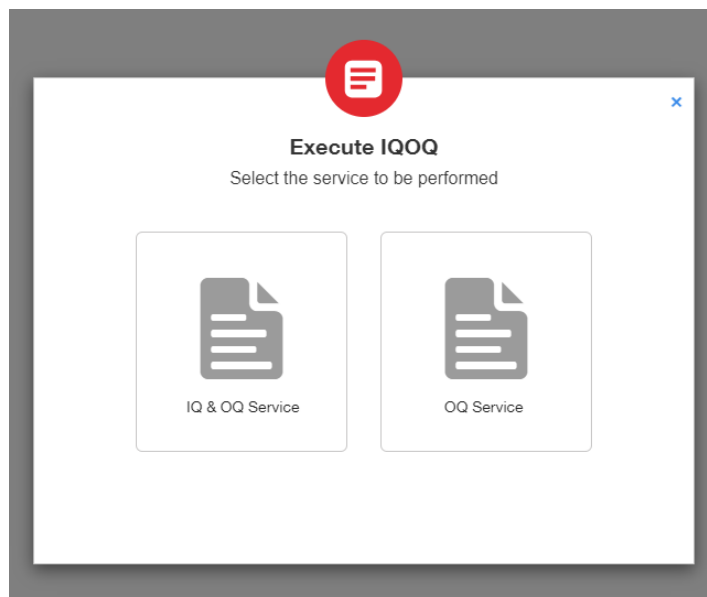


The license is now activated and will remain valid for 2 months on the instrument. Press **Next** to advance to the OQ testing. Follow the OQ instructions under Execute the Operational Qualification (OQ) on page 59.

Execute the Installation Qualification (IQ)

Each entry throughout the IQ process is recorded in a final PDF report which is then saved to a USB drive at the end of the IQ procedure. During the IQ workflow, a license will be saved to the USB drive to allow execution of the on-instrument OQ workflow. After completion of the OQ on-instrument, the instrument will allow export of a final IQOQ report. If needed at any step, press **Cancel** to return to the last saved location.

1. Select **IQ&OQ Service**.



2. Enter the **Protocol Execution Date, Instrument Serial Number, Lab Name, Lab Address, Contact Name, Email, and Telephone Number.**

IQOQ Process - Step 1/3
Enter the details

Protocol Execution Date: 07/12/2023

Instrument Serial Number: Serial Number

Lab Name: Enter lab Name

Contact Name: Enter Contact Name

Lab Address: Enter Lab Address...

Contact Email Address: Enter contact email address

Contact Telephone Number: Enter contact telephone number

Buttons: Cancel, Next

3. Select the reason (from the drop-down menu) for completing the IQ/OQ service .

IQOQ Process - Step 2/3
Pre-execution Verification

Service Identification

Select reason for completing IQ/OQ

IQOQ Reason

Select the reason for completing IQ/OQ

- New Installation
- Reinstallation/Move
- Replacement

Buttons: Back, Cancel, Save & Continue

4. Ensure that you have access to the appropriate user guide and verify that you have read the site requirements.

IQOQ Process - Step 2/3
Pre-execution Verification

Service Identification ✓ +

Site requirement verification ✓ -

Verify that relevant user reference document is available.

Countess 3 or 3 FL Automatic Cell Counter Guide (Pub. No. MAN0019566 for Countess 3 or Pub. No. MAN0019567 for Countess 3FL)

Pass Fail

Verify that you have read environmental site requirements listed in the relevant user guide and the site complies with site requirements.

Countess 3 or 3 FL Automatic Cell Counter Guide (Pub. No. MAN0019566 , page

Back Cancel Save & Continue

5. Verify that the **Service identification** and **Site requirement verification** sections have been completed. Enter **User name**, **User Initials**, and **Date**.

IQOQ Process - Step 2/3
Pre-execution Verification

Service Identification ✓ +

Site requirement verification ✓ +

Test Completion -

Verify required test completion

All Tests Completed Yes No

User Initials Date

Enter user initials 07/12/2023

Back Cancel Save & Continue

6. Confirm that the automatically-filled sales order number (Document Number) is correct.

IQOQ Process - Step 3/3
Installation Qualification Verification

Order verification

IQ/OQ service Yes No

Document Title: Sales Order
Document Number: 51330009

Pass Fail

Verify the system components +

Verify that all packing kit components are available +

Verify that all required IQ tests have been completed +

Back Cancel Save

7. Verify that all of the system components are unpacked and available. Confirm that packing components are as described in the user guide.

IQOQ Process - Step 3/3
Installation Qualification Verification

Order verification ✓ +

Verify the system components ✓ -

Verify that the system components are unpacked and positioned appropriately as described in user manual [Pub. No. MAN0019566](#) (page 8) or [Pub. No. MAN0019567](#) (page 10) as applicable.

Actual Result Pass Fail

Verify that all packing kit components are available +

Verify that all required IQ tests have been completed +

Back Cancel Save

IQOQ Process - Step 3/3
Installation Qualification Verification

Order verification ✓ +

Verify the system components ✓ +

Verify that all packing kit components are available ✓ -

Verify that all packing components are available as described in the user manual [Pub. No. MAN0019566](#) (page 6) or [Pub. No. MAN0019567](#) (page 7) as applicable.

Actual Result Pass Fail

Verify that all required IQ tests have been completed +

Back Cancel Save

8. Verify that all required IQ tests are complete. Enter **Initials** and **Date**, then press **Save**.

IQOQ Process - Step 3/3
Installation Qualification Verification

- Order verification ✓ +
- Verify the system components ✓ +
- Verify that all packing kit components are available ✓ +
- Verify that all required IQ tests have been completed -

All Tests Completed Yes No

User Initials: Date:

Buttons: Back, Cancel, Save

9. Press **Download** to download the license to execute OQ on the instrument. Save the license to a USB drive. Do not change the license file name "Prod_xxxx.OQ-license" where xxxx is the instrument serial number. Next insert the drive into the instrument and navigate to **Settings-Execute OQ** on the Countess™ 3 / 3 FL touchscreen.

Download Activation File

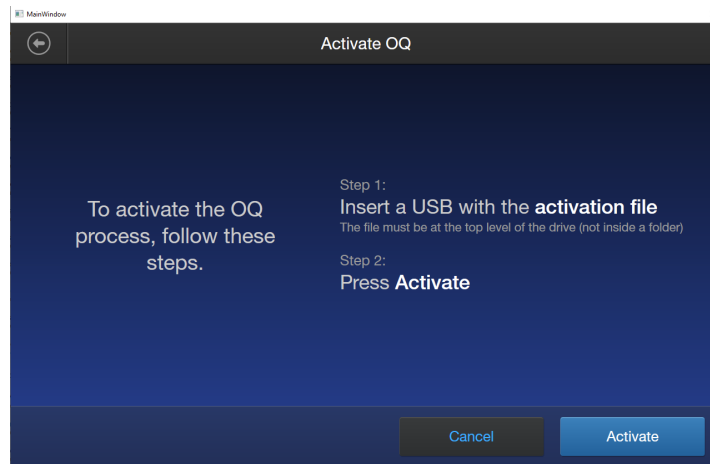
Click on download to get the activation file.

Follow the below steps after downloaded the activation file:

Transfer the file to USB → Connect USB to instrument → Click Execute OQ

Buttons: Back, Cancel, Download

10. **Activate** the OQ workflow.



Note: The license is now activated and will remain valid for 2 months on the instrument.

11. Follow the OQ instructions in the next section, Execute the Operational Qualification (OQ)

Execute the Operational Qualification (OQ)

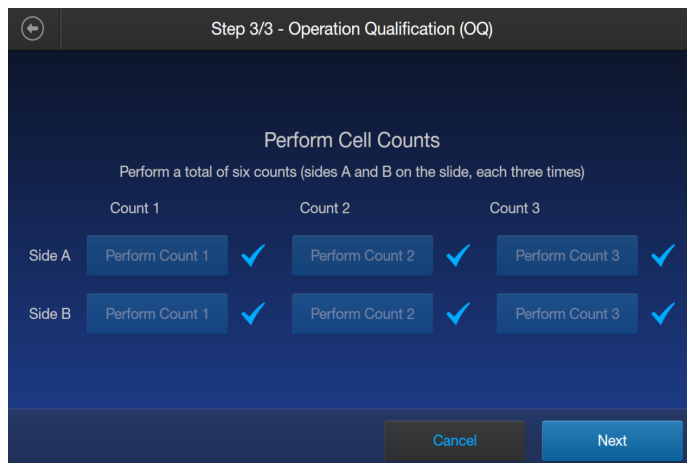
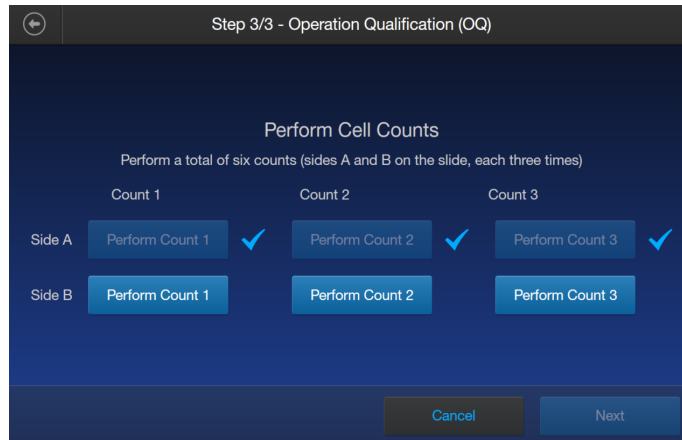
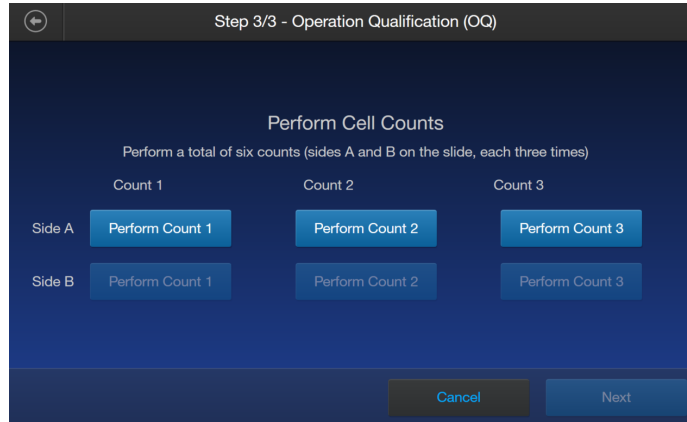
Each entry throughout the OQ process is recorded in a final PDF report saved at the end of the OQ protocol. If needed at any step, press **Cancel** to return to the last screen. The qualification can only be restarted by completing the full workflow and saving the report.

Before completing the OQ testing, ensure the checkbox for **Rapid Capture** on the main screen does not have a check mark. If the box is checked, please press the box to remove the checkmark. The OQ procedure must be completed with **Rapid Capture** turned off.

1. Enter the **Protocol Execution Date**, **Lab Name**, and **Lab Address**. The instrument serial number will automatically display. Press **Next**.
2. Enter your **Contact Name**, **Contact Email Address**, and **Contact Telephone Number**. Press **Next**.
3. Select an OQ reason of **Annual requalification**, **Software update/upgrade**, or **After repair**. Press **Next**.
4. Verify pre-execution is completed by selecting **Yes** in **All Tests Completed** and entering your **User Initials** and the **Date**. Press **Next**.
5. Verify the instrument has successfully booted by selecting **Pass** or **Fail** and confirm the **Date & Time Format** are correct. Press **Next**.
6. Record if the software has been updated prior to OQ by selecting **Yes** or **No**. The instrument software version will automatically display. Press **Next**.

7. Enter the Part Number and Lot Number from the Countess™ 3/3 FL standard slide packaging. Press **Next**.

8. Perform a series of 6 counts of the standard slide (3 counts on each side). Insert side A of the standard slide and press the appropriate count number (**Perform Count 1**, **Perform Count 2**, or **Perform Count 3**). The instrument will automatically autofocus and take images in the background. After completing the 3 counts of side A, remove the slide and insert side B. Side B will only illuminate after removing the slide and insert side B. Insert side B of the standard slide and press the appropriate count number (**Perform Count 1**, **Perform Count 2**, or **Perform Count 3**). Press **Next**.



9. Total concentration results are displayed (not editable). **Pass** or **Fail** is automatically selected based on the recorded count concentrations shown above the counts. Press **Next**.

The screenshot shows a software interface for 'Step 3/3 - Operation Qualification (OQ)'. The main heading is 'Total Concentration Results'. Below this, it states 'Recorded count concentrations (0.5 x 10⁶ - 1.5 x 10⁶ beads/mL)'. The data is organized into two columns: 'Side A' and 'Side B'. There are three rows of counts, labeled 'Count 1', 'Count 2', and 'Count 3'. Each row has two input fields, one for Side A and one for Side B, both containing the value '9.81 x 10⁵'. At the bottom, there is an 'Actual Result' section with two radio buttons: 'Pass' (which is selected) and 'Fail'. At the very bottom, there are two buttons: 'Cancel' and 'Next'.

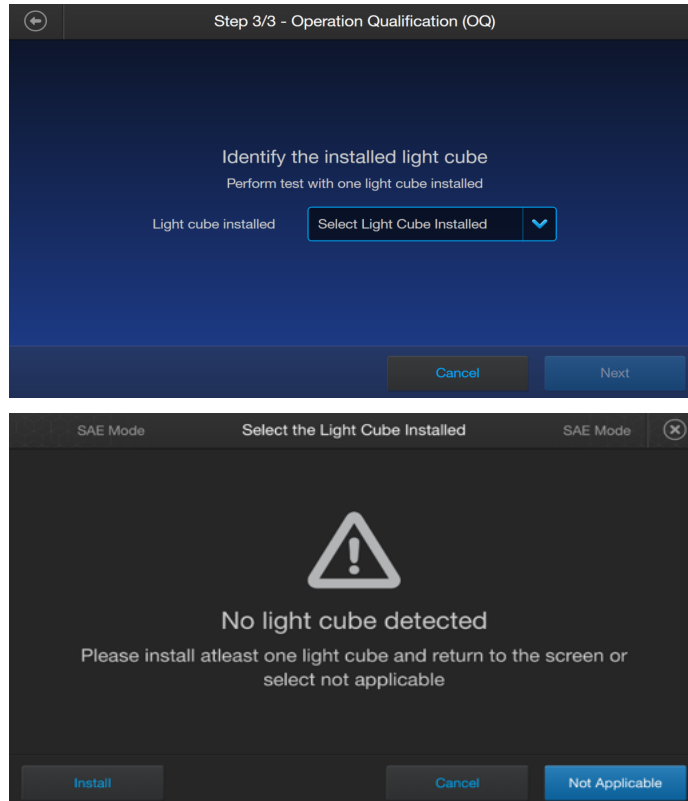
	Side A	Side B
Count 1	9.81 x 10 ⁵	9.81 x 10 ⁵
Count 2	9.81 x 10 ⁵	9.81 x 10 ⁵
Count 3	9.81 x 10 ⁵	9.81 x 10 ⁵

Actual Result: Pass Fail

Buttons: Cancel, Next

10. The mean, standard deviation, and coefficient of variation (Cv) are displayed (not editable). **Pass** or **Fail** is automatically selected based on the number shown under Cv. Press **Next**.
11. Total viability results are displayed (not editable). **Pass** or **Fail** is automatically selected based on the recorded % live results shown above the counts. Press **Next**.
12. The mean, standard deviation, and coefficient of variation (Cv) are displayed (not editable). **Pass** or **Fail** is automatically selected based on the number shown under Cv. Press **Next**.

13. **(For the Countess™ 3 FL instrument)** Choose the appropriate light cube for testing. The OQ procedure tests fluorescent counts with a single light cube. If no cube is installed, press **Install**. If not required, press **Not Applicable**. If you have a Countess™ 3 instrument, proceed to step 18.



14. Select the appropriate light cube from the list of available cubes. Press **Next**.

15. Perform a series of 6 counts of the standard slide (3 counts on each side). Insert side A of the standard slide and press the appropriate count number (**Perform Count 1**, **Perform Count 2**, or **Perform Count 3**). The instrument will automatically autofocus and take images in the background. After completing the 3 counts of side A, remove the slide and insert side B. Side B will only illuminate after removing the slide and insert side B. Insert side B of the standard slide and press the appropriate count number (**Perform Count 1**, **Perform Count 2**, or **Perform Count 3**). Press **Next**.

Step 3/3 - Operation Qualification (OQ)

Perform Cell Counts

Perform a total of six counts (sides A and B on the slide, each three times)

	Count 1	Count 2	Count 3
Side A	Perform Count 1	Perform Count 2	Perform Count 3
Side B	Perform Count 1	Perform Count 2	Perform Count 3

Cancel Next

Step 3/3 - Operation Qualification (OQ)

Perform Cell Counts

Perform a total of six counts (sides A and B on the slide, each three times)

	Count 1	Count 2	Count 3
Side A	Perform Count 1 ✓	Perform Count 2 ✓	Perform Count 3 ✓
Side B	Perform Count 1	Perform Count 2	Perform Count 3

Cancel Next

Step 3/3 - Operation Qualification (OQ)

Perform Cell Counts

Perform a total of six counts (sides A and B on the slide, each three times)

	Count 1	Count 2	Count 3
Side A	Perform Count 1 ✓	Perform Count 2 ✓	Perform Count 3 ✓
Side B	Perform Count 1 ✓	Perform Count 2 ✓	Perform Count 3 ✓

Cancel Next

16. Total concentration results are displayed (not editable). **Pass** or **Fail** is automatically selected based on the recorded count concentrations shown above the counts. Press **Next**.
17. The mean, standard deviation, and coefficient of variation (Cv) are displayed (not editable). **Pass** or **Fail** is automatically selected based on the number shown under Cv. Press **Next**.
18. Confirm the tests are successfully completed by pressing **Yes** and entering your **User Initials** and the **Date**. Press **Save**.



19. Enter the name for the Qualification report and select the save location for the report and attachments. Press **Save**.

Note: Only the comment and signature sections of the exported PDF report are editable for verifying the report.

4.2. Final instructions [continued]

4.2.2. If this service is completed using the electronic form, then the user must :

- Obtain the Customer's signature in Section 4.3.2. System Owner or designee final approval.
- If applicable, obtain any additional signer signatures in Section 4.3.3. (Optional) Additional signer(s) final approval.
- Sign in Section 4.3.1. User final approval.
- Scan any non-electronic supporting documentation.
- Email the completed, signed , and locked OQ Service Protocol, any exception reports, and all supporting documentation to the Customer.

Final approval signatures

4.3.1. User final approval

Print user name : Initials:

Title/Role:

Signature:

Date: - -

Note: Ensure **Rapid Capture** is turned off before completing the OQ procedure.

Note: If any of the tests fail during the procedure, we recommend repeating the OQ workflow. Ensure the Countess™ 3 Standard Slide is free of debris by wiping with a lint-free wipe. If the OQ tests continue to fail, contact technical support at thermofisher.com/support.



Troubleshooting

Observation	Possible cause	Recommended action
Uneven screen illumination (screen is dark on one side, bright on the other side)	Light cubes misaligned.	Reset the light cube tray by selecting Change light cube from the Instrument Settings screen.
Autofocus does not focus properly	Debris or other material interfered with the focus.	Make sure there are no bubbles or debris visible on the screen that could interfere with the autofocus and make it more difficult to get the sample in the correct focal plane.
	Live cells vs dead cells were difficult to distinguish.	If cells are well focused, have bright centers, and are being counted as dead, confirm that they are within the appropriate cell size range and try adjusting the settings.
Some cells appear in the image but are not included in the count	Brightfield or fluorescence (only in the Countess™ 3 FL Automated Cell Counter) settings were not optimal.	BF-based counts rely on the BF image to be in focus. Adjust the focus and recount with BF-based count.
		FL-based counts are more sensitive and detect objects with low fluorescent intensity. Check the FL-based box to perform fluorescence counts to be sure they are included in the count.
		For cell count and cell viability assays performed in brightfield with BF-based counts, adjust the size, brightness, and circularity gates for both live and dead cells to include all of the cells in the count.
		For fluorescence assays with both BF-based and FL-based counts, adjust the size, brightness, circularity, and fluorescence intensity gates in all available channels to include all of the cells in the count.
		After including all of the cells in the count, you can narrow the count criteria, if you wish to exclude cells of a certain size or certain brightness.
		When the gates are fully maximized, the CSV should indicate 0–70 for cell size and 0–255 for brightness.
		Be sure channels are optimally illuminated in both brightfield and fluorescence modes.
Fluorescence is extremely bright and bleeding through into other filters	Light intensity was not set correctly.	Decrease the fluorescence light intensity before counting the cells.

Observation	Possible cause	Recommended action
Incorrect concentration for the Countess™ test beads	Beads were not suspended in solution or mixed correctly.	The beads can settle quickly in solution, which will affect the concentration reading.
		Vortex the bead stock on high for a full 30 seconds to resuspend, and add 10 µL of the bead suspension to 10 µL of trypan blue without delay.
		Pipet the bead and trypan blue mixture up and down several times to make sure it is well mixed and immediately load 10 µL into the slide.
Variable counts for the same sample of cells	Pipetting was done incorrectly.	If you are pipetting different samples from the same cell sample, the variability could be due to pipetting or mixing.
		Use recently calibrated pipettors and make sure that the cells are well suspended by pipetting up and down several times before adding trypan blue.
Variable counts when performing replicate counts of the same slide	Slide was mishandled or viewed differently multiple times.	If you are counting replicates of the exact same slide, visually inspect that all cells are counted correctly in the image.
		There may be a slightly different field of view each time a slide is inserted. Depending on the concentration and uniformity of the cells, this will cause some variability when performing replicate counts of the same slide, although it should be less than 10%.
		When counting fewer cells, a small field of view change for only a small number of cells can have a larger affect. Count cells at a higher concentration to reduce variability
		Make sure you do not shake or agitate the slide between counts.
Abnormally high percentage of dead cells or live cells counted as dead	Focus was not correct for the sample.	Ensure that the cells are focused correctly so that live cells have bright centers and dead cells are dark throughout. If the cells are not well focused and look dark on the screen, the Countess™ 3 FL cell counter will count them as dead cells.
	Live cells vs dead cells were difficult to distinguish.	If cells are well focused, have bright centers, and are being counted as dead, confirm that they are within the appropriate cell size range and try adjusting the settings.
	Cells were exposed to trypan blue for too long.	If cells are exposed to trypan blue for a long period of time, viability could be affected so slide should be prepared and counted fresh each time for best results.
	Capture settings were not optimal.	Gate out the debris using the size, brightness, and circularity sliders.



Observation	Possible cause	Recommended action
USB drive not recognized by the instrument	USB drive was not correctly formatted.	The USB drive must be FAT32 formatted to be recognized by the instrument. If it is not, reformat the USB drive to FAT32.
		Try another correctly formatted USB drive.
Unable to update the Countess™ software	USB drive did not function properly.	Make sure the USB drive is formatted to FAT32. If it is not, reformat the USB drive to FAT32 before transferring the files onto the USB drive for software update.
		Make sure the update file is at the top (root) of the USB drive, not in a folder or subfolder.
		File cannot be renamed in any way.
		File cannot be zipped or compressed during distribution. It must be uncorrupted during transfer and have a .lft suffix.
		If needed, check that the USB port is functional by testing a USB mouse.



Technical specifications

Physical characteristics

Instrument type	Benchtop cell counter and suspension cell-based assay platform
Instrument dimensions	9.6 (W) × 6.7 (D) × 9.4 inches (H)
Weight	8 lbs
Operating power	100–240 VAC, 1.0 A MAX
Frequency	50/60 Hz
Electrical input	12 VDC, 3 A
Installation site	Indoor use only, Class A Environments (i.e., non-residential or light industrial); Pollution degree 2.
Operating temperature	10°–40°C (50°–104°F)
Operating humidity	<80% (non-condensing)

Technical specifications

Processing time	<30 seconds
Sample concentration range	1×10^4 – 1×10^7 cells/mL
Particle/cell diameter range	4–60 μ m
Required sample volume	10 μ L
Firmware	Countess™ Automated Cell Counting Platform Software
USB drive	32 Gigabytes (FAT32-, exFAT-, or NTFS-compatible)

Optics

Optics	3 channels (brightfield and 2 slots for EVOS™ LED light cubes)
Camera	5 megapixels, 2.5× Optical Magnification

Analysis slide

Material	Poly(methyl methacrylate) (PMMA)
Dimensions	25 mm (W) × 75 mm (D) × 1.7 mm (H)
Chamber volume	10 µL

Networking requirements

Configure the system behind a firewall. If outbound traffic is limited, the following firewall exceptions are required to support system features:

Firewall exception requirements			
URL	Port	Purpose	Applies to...
*.instrumentconnect.com	outbound 443	To support instrument management and identify	Thermo Fisher™ Connect Platform only
*.thermofisher.com	outbound 443	To support instrument management and identify	Thermo Fisher™ Connect Platform LAN connection
*.s3-us-east-1.amazonaws.com	outbound 443	To allow connection to the Thermo Fisher™ Connect Platform	Thermo Fisher™ Connect Platform only
*.iot-us-east-1.amazonaws.com	outbound 443	To allow connection to the Thermo Fisher™ Connect Platform	Thermo Fisher™ Connect Platform only
Allowed port requirements			
—	7443	To support instrument discovery Uses multicast address 224.0.0.251	LAN connection Direct connection
—	TCP 445 (SMB v3 or higher)	To support file sharing	LAN connection Direct connection
—	8443 (default)	Connection to SAE Administrator Console ^[1]	LAN connection Direct connection

^[1] The SAE Administrator Console should be installed in a computer with a static IP address. The console opens in a browser (recommended Google Chrome) and communicates with the instrument. Communication between the SAE Administrator Console and the instrument uses the encrypted HTTPS protocol.



EVOS light cubes

LED illumination

The Countess™ 3 FL Automated Cell Counter utilizes an adjustable intensity LED light source provided by the proprietary, user-interchangeable LED light cube (see “EVOS™ light cubes” on page 87). Because the LED light source is as close as possible to the objective, the number of optical elements in the channel is minimized. High-intensity illumination over a short channel increases the efficiency of fluorophore excitation, providing better detection of weak fluorescent signals.

EVOS™ light cubes

Each user-interchangeable, auto-configured EVOS™ light cube contains an LED, collimating optics, and filters. In addition to the brightfield channel dedicated to cell count and cell viability assays using Trypan Blue, the Countess™ 3 FL Automated Cell Counter can accommodate two fluorescent light cubes for multiple-fluorescence research applications.



The following table lists some of the common fluorescent and specialty EVOS™ light cubes available from Thermo Fisher Scientific. For a complete list, go to www.thermofisher.com/evoslightcubes or contact Technical Support. For instructions on changing the LED light cubes, see Chapter 10, “Change EVOS™ light cubes”.

Light Cube	Dyes
DAPI	DAPI, Hoechst™, BFP
TagBFP	TagBFP
CFP	ECFP, Lucifer Yellow, Evans Blue
GFP	GFP, Alexa Fluor™ 488, SYBR Green™, FITC, Acridine orange (AO)
YFP	EYFP, acridine orange + DNA



(continued)

Light Cube	Dyes
RFP	RFP, Alexa Fluor™ 546, Alexa Fluor™ 555, Alexa Fluor™ 568, Cy™3, MitoTracker™ Orange, Rhodamine Red, DsRed
Texas Red	Texas Red™, Alexa Fluor™ 568, Alexa Fluor™ 594, MitoTracker™ Red, mCherry, Cy3.5™, Propidium Iodide (PI)
Cy5	Cy5™, Alexa Fluor™ 647, Alexa Fluor™ 660, DRAQ5™
Cy5.5	Cy5.5™, Alexa Fluor™ 660, Alexa Fluor™ 680, Alexa Fluor™ 700
Cy7	Cy7™, IRDye 800CW

Note: The EVOS™ light cubes are available only for the Countess™ 3 FL Automated Cell Counter. The Countess™ 3 Automated Cell Counter uses only brightfield illumination and does not support the EVOS™ light cubes.



Countess™ 3 FL Automated Cell Counter and accessories

The following Countess™ 3 FL instruments and instrument accessories are available from Thermo Fisher Scientific. For more information, visit www.thermofisher.com or contact Technical Support.

Product	Quantity	Cat. No.
Countess™ 3 FL Automated Cell Counter	1 each	AMQAF2000
Countess™ 3 Power Adapter with four adaptor cords	1 each	A48207
Countess™ 3 USB Drive	1 each	A26774
Countess™ 3 FL Light Cube Removal Tool	1 each	AMEP-4747
Countess™ 3 FL Disposable Slide Holder	1 each	AMEP-4745
Countess™ 3 FL Reusable Slide Holder	1 each	A48208

Accessory products

The following products can be used with the Countess™ 3/3 FL Automated Cell Counters and are available separately from Thermo Fisher Scientific. For more information, visit www.thermofisher.com or contact Technical Support.

Product	Quantity	Cat. No.
Countess™ Cell Counting Chamber Slides, 50 Slides (100 counts)	1 box ^[1]	C10228
Countess™ Cell Counting Chamber Slides, 500 Slides (1000 Counts)	10 boxes ^[1]	C10312
Countess™ Cell Counting Chamber Slides, 1250 Slides (2500 Counts)	25 boxes ^[1]	C10313
Countess™ Cell Counting Chamber Slides, 2500 Slides (5000 Counts)	50 boxes ^[1]	C10314
Countess™ Cell Counting Chamber Slides, 5000 Slides (10,000 Counts)	100 boxes ^[1]	C10315
Countess™ Reusable Slide	1 each	A25750
Countess™ Test Beads (1 × 10 ⁶ beads/mL)	1 mL	C10284
Trypan Blue Stain (0.4%)	2 × 1 mL	T10282
Countess™ Cell Counting Slander Slide	1 each	A51876
SafeCount™	1 each	A40008024

(continued)

Product	Quantity	Cat. No.
ReadyCount™ Viability Stain	1 each	A49905
ReadyCount™ Dead cell Stain	1 each	A49903
ReadyCount™ Nuclear Stain	1 each	A49904

^[1] Each box of Countess™ Cell Counting Chamber Slides contains 50 slides and 2 × 1 mL vials of Trypan Blue (0.4%), sufficient for 100 counts.



CSV file format definition

Overview

A comma-separated values (CSV) file stores tabular data (numbers and text) in plain-text form. Plain text means that the file is a sequence of characters, with no data that has to be interpreted as binary numbers. A CSV file can be opened with any third party software or spreadsheet program. The below table describes the categories of the Countess™ 3 FL data saved as a CSV file and opened with a spreadsheet program.

Category	Column	Name	Description
General	A	Count ID	ID number for current count
	B	Session ID	ID number for current session
	C	Sample name	Name of sample
	D	Date & Time	Date and time of sample run
	E	Count mode	BF-based or FL-based
Sample	F	Type	BF-Brightfield or FL-Fluorescence
	G	1:1 Stain corrected	Sample with 1:1 stain correction applied
	H	Pre-Dilution corrected	Sample with pre-dilution correction applied
	I	Total concentration	Concentration of the entire sample
	J	Total cells counted	Total number of cells counted in the sample
Brightfield/Live Dead	K	Live concentration	Concentration of just the "live" portion of the sample
	L	Live cells counted	"Live" cells counted in sample
	M	Dead concentration	Concentration of just the "dead" portion of the sample
	N	Dead cells counted	"Dead" cells counted in sample
	O	Viability (%)	Percent viability of the sample based on trypan blue staining
	P	Live average size (µm)	Average size of "live" cells in microns
	Q	Dead average size (µm)	Average size of "dead" cells in microns
Fluorescence	R	Cube 1 name	EVOS™ light cube name in the first (top) position

(continued)

Category	Column	Name	Description
Fluorescence	S	Cube 1 concentration	Concentration of cells showing fluorescence in the first cube position
	T	Cube 1 (%)	Percentage of the total cells in brightfield that show fluorescence in the first cube position
	U	Cube 1 cells counted	Total number of cells counted in the first cube position
	V	Cube 2 name	EVOS™ light cube name in the second (bottom) position
	W	Cube 2 concentration	Concentration of cells showing fluorescence in the second cube position
	X	Cube 2 (%)	Percentage of the total cells in brightfield that show fluorescence in the second cube position
	Y	Cube 2 cells counted	Total number of cells counted in the second cube position
	Z	Cube 1+2 concentration	Concentration of cells showing fluorescence in the first and second cube positions combined
	AA	Cube 1+2 (%)	Percentage of the total cells in brightfield that show fluorescence in the first and second cube positions combined
	AB	Cube 1+2 cells counted	Total number of cells counted in the first and second cube positions combined
	AC	Cube 1 average size (µm)	Average size in microns of cells counted using the first (top) cube
	AD	Cube 2 average size (µm)	Average size in microns of cells counted using the second (bottom) cube
General	AE	Focus value	Focal position number
	AF	Focus motor value	Motor position in relation to focal position number
Brightfield/Live Dead	AG	BF light intensity	Brightfield light intensity value from 0-100%
	AH	BF LED intensity	LED intensity in brightfield mode
	AI	Live size min	Minimum size of "live" cells in microns
	AJ	Live size max	Maximum size of "live" cells in microns
	AK	Live brightness min	"Live" adjustment slider value for minimum brightness

(continued)

Category	Column	Name	Description
Brightfield/Live Dead	AL	Live brightness max	"Live" adjustment slider value for maximum brightness
	AM	Live circularity min	"Live" adjustment slider value for minimum circularity
	AN	Live circularity max	"Live" adjustment slider value for minimum circularity
	AO	Dead size min	Minimum size of "dead" cells in microns
	AP	Dead size max	Maximum size of "dead" cells in microns
	AQ	Dead brightness min	"Dead" adjustment slider value for minimum brightness
	AR	Dead brightness max	"Dead" adjustment slider value for maximum brightness
	AS	Dead circularity min	"Dead" adjustment slider value for minimum circularity
	AT	Dead circularity max	"Dead" adjustment slider value for maximum circularity
Fluorescence	AU	Cube 1 light intensity	First (top) light cube light intensity value from 0-100%
	AV	Cube 1 LED intensity	LED intensity of first (top) light cube
	AW	Cube 2 light intensity	Second (bottom) light cube light intensity value from 0-100%
	AX	Cube 2 LED intensity	LED intensity of second (bottom) light cube
Brightfield	AY	BF size min	Minimum size of "brightfield" cells in microns
	AZ	BF size max	Maximum size of "brightfield" cells in microns
	BA	BF brightness min	"Brightfield" adjustment slider value for minimum brightness
	BB	BF brightness max	"Brightfield" adjustment slider value for maximum brightness
	BC	BF circularity min	"Brightfield" adjustment slider value for minimum circularity
	BD	BF circularity max	"Brightfield" adjustment slider value for maximum circularity
Fluorescence	BE	Cube 1 brightness min	First (top) light cube adjustment slider value for minimum brightness

(continued)

Category	Column	Name	Description
Fluorescence	BF	Cube 1 brightness max	First (top) light cube adjustment slider value for maximum brightness
	BG	Cube 1 size min	Second (bottom) light cube adjustment slider value for minimum brightness
	BH	Cube 1 size max	Second (bottom) light cube adjustment slider value for maximum brightness
	BI	Cube 1 circularity min	Template used for count
	BJ	Cube 1 circularity max	Current software version used for count
	BK	Cube 2 brightness min	Second (bottom) light cube adjustment slider value for minimum brightness
	BL	Cube 2 brightness max	Second (bottom) light cube adjustment slider value for maximum brightness
	BM	Cube 2 size min	Minimum size of second (bottom) light cube identified cells in microns
	BN	Cube 2 size max	Maximum size of second (bottom) light cube identified cells in microns
	BO	Cube 2 circularity min	Second (bottom) light cube adjustment slider value for minimum circularity
	BP	Cube 2 circularity max	Second (bottom) light cube adjustment slider value for maximum circularity
General	BQ	Protocol name	Current protocol used for the count
	BR	Software revision	Current software version used for count
	BS	Aggregation(%)	Total number of cells involved in an aggregate or aggregates vs. total cell count



Safety



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](https://www.thermofisher.com/support).

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 sufficient ventilation (for example, fume hood).
- Check regularly for chemical leaks or spills. If a leak or spill occurs, follow the manufacturer 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 needed) 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.



WARNING! HAZARDOUS WASTE (from instruments). Waste produced by the instrument is potentially hazardous. Follow the guidelines noted in the preceding General Chemical Handling warning.



WARNING! 4L Reagent and Waste Bottle Safety. Four-liter reagent and waste bottles can crack and leak. Each 4-liter bottle should be secured in a low-density polyethylene safety container with the cover fastened and the handles locked in the upright position.

Biological hazard safety



WARNING! Potential Biohazard. Depending on the samples used on this instrument, the surface may be considered a biohazard. Use appropriate decontamination methods when working with biohazards.



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)*, 6th Edition, HHS Publication No. (CDC) 300859, Revised June 2020
https://www.cdc.gov/labs/pdf/SF__19_308133-A_BMBL6_00-BOOK-WEB-final-3.pdf
- Laboratory biosafety manual, fourth edition. Geneva: World Health Organization; 2020 (Laboratory biosafety manual, fourth edition and associated monographs)
who.int/publications/i/item/9789240011311



Documentation and support

Customer and technical support

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

