



# ImageXpress® Pico

**Automated Cell Imaging System**

with CellReporterXpress® Software Version 2.5

## Calibration Kit Guide

This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

**For research use only. Not for use in diagnostic procedures.**

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patents: [www.moleculardevices.com/patents](http://www.moleculardevices.com/patents)

Product manufactured by Molecular Devices, LLC.  
3860 N. First Street, San Jose, California, 95134, United States of America.  
Molecular Devices, LLC is ISO 9001 registered.  
©2019 Molecular Devices, LLC.  
All rights reserved.



# Contents

<b>Chapter 1: ImageXpress Pico Automated Cell Imaging System</b>	<b>5</b>
Obtaining Support	6
Product Documentation	7
About This Guide	7
<b>Chapter 2: Package Contents</b>	<b>9</b>
<b>Chapter 3: Objective Maintenance</b>	<b>11</b>
Installing an Objective	12
Calibrating an Objective	16
Adjusting an Objective Correction Collar	18
Cleaning an Objective	23
<b>Chapter 4: Filter Cube Maintenance</b>	<b>25</b>
Installing a Filter Cube	25
Calibrating a Filter Cube	29
<b>Appendix A: Optional Parts and Accessories</b>	<b>31</b>
Objectives	31
Filter Cubes	31
Accessories	32





# Chapter 1: ImageXpress Pico Automated Cell Imaging System

# 1

The Molecular Devices® ImageXpress® Pico Automated Cell Imaging System is an all-in-one platform for automatically acquiring and analyzing images from fluorescently labeled biological samples in plates and slides. It enables you to increase the throughput of your image acquisition and analysis, allowing you to gain insights in minutes. With the addition of a modular option, the system provides environmental control for live cell imaging.

The core hardware component of the imaging system is a custom-designed, fully automated, epi-illumination fluorescence microscope. The rapid autofocus and precision sample movement features of the microscope allow large numbers of high-resolution images to be acquired in the shortest possible time. All key optical and mechanical elements are motorized, which allows complete control of the instrument configuration.

When used in combination with the CellReporterXpress® Image Acquisition and Analysis Software, which features powerful image analysis capabilities and more than 25 available predefined experimental protocols, the ImageXpress Pico system becomes an extremely flexible device, ideally suited for user-defined, automated assays.

Key components of the instrument include the following:

- Built-in, internal light source comprised of six high-powered LEDs enables very high sensitivity fluorescent imaging. Additional LED-based light sources allow transmitted light and overview imaging.
- High-sensitivity, 5-megapixel CMOS camera.
- Hardware-based autofocus system with precision motorized Z-stage focus.
- High-transmission fluorescence imaging optics with world-class chromatic aberration correction, resolution, and image flatness.
- Secondary camera provides overview imaging for slides.
- Precision motorized sample (X-Y) stage.
- High-quality Leica objectives in a six-position turret.
- Filter cubes in a six-position turret.
- Motorized selection of stage position, filter cubes, and objectives.
- Temperature control up to 40°C (104° F) for live cell imaging.
- Optional environmental control system, which enables you to regulate humidity, CO<sub>2</sub>, and O<sub>2</sub> inside the environmental control cassette for multi-day, live-cell, time-lapse imaging.
- Operation and configuration control by the integrated CellReporterXpress software.

## Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software, and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest level of technical service.

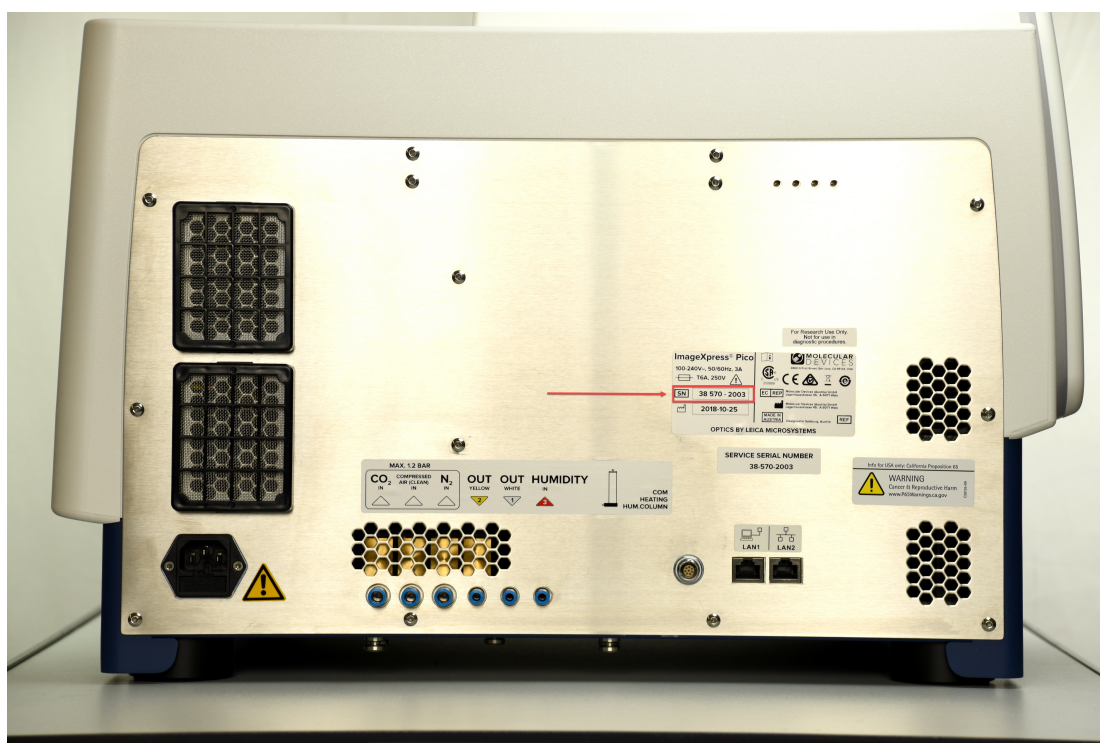
Our Support website, [www.moleculardevices.com/service-support](http://www.moleculardevices.com/service-support), describes the support options offered by Molecular Devices, including service plans and professional services. It also has a link to our Knowledge Base, which contains documentation, technical notes, software upgrades, safety data sheets, and other resources. If you still need assistance after consulting the Knowledge Base, you can submit a request to Molecular Devices Technical Support.

### Technical Support

You can contact Molecular Devices Technical Support by submitting a support request through the Knowledge Base or by phone. To find regional support contact information, visit [www.moleculardevices.com/contact](http://www.moleculardevices.com/contact).

You need the instrument serial number.

The serial number is located on the back panel of the instrument.



## Documentation

Review the product documentation on the Knowledge Base, including installation guides and user guides. In addition, online Help is available within the CellReporterXpress software. Press **F1** to access Help for the current page.

## Additional Resources

Web-based microscopy courses:

- [www.leica-microsystems.com/science-lab](http://www.leica-microsystems.com/science-lab)
- [www.ibiology.org/ibioeducation/taking-courses/ibiology-microscopy-short-course.html](http://www.ibiology.org/ibioeducation/taking-courses/ibiology-microscopy-short-course.html)

The Molecular Probes Handbook offers advice on fluorescent probes and can help you determine if there are better stains available for your analysis:

- [www.lifetechnologies.com/us/en/home/references/molecular-probes-the-handbook.html](http://www.lifetechnologies.com/us/en/home/references/molecular-probes-the-handbook.html)

## Product Documentation

The following guides are available on the Molecular Devices Knowledge Base at [mdc.custhelp.com](http://mdc.custhelp.com):

- *CellReporterXpress Installation Guide*
- *CellReporterXpress Release Notes*
- *CellReporterXpress User Guide*
- *ImageXpress Pico Pre-Installation Guide*
- *ImageXpress Pico Installation Guide*
- *ImageXpress Pico User Guide*
- *ImageXpress Pico Calibration Kit Guide*

In addition, the CellReporterXpress software includes context-sensitive Help that you can access from within the software. Just press the **F1** key from within the software to view Help for the current page.



**Tip:** Molecular Devices recommends that you review the documentation before installing or using the ImageXpress Pico system or the CellReporterXpress software.

---

## About This Guide

This guide is intended for the scientist who receives a calibration kit along with the aftermarket purchase of an objective or filter cube for the ImageXpress Pico system. It describes how to install and calibrate the objective or filter cube.



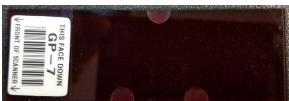
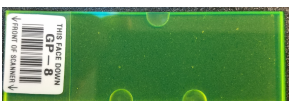
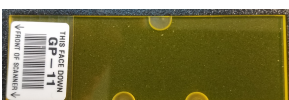
The information in this guide is subject to change without notice. Molecular Devices recommends that you review the guide on the Knowledge Base for the most up-to-date information.



## Chapter 2: Package Contents

# 2

The Calibration Kit includes the following:

Item	Part number	Description
	5063669	Calibration Kit Guide
	9100-0099	Fluorescent Bead Slide
	1-GP-7	Red Plastic Slide
	1-GP-8	Green Plastic Slide
	1-GP-11	Yellow Plastic Slide

For a complete list of the contents of the package, see the enclosed packing list.



## Chapter 3: Objective Maintenance

Objective maintenance procedures include the following:

- [Installing an Objective, see page 12](#)
- [Calibrating an Objective, see page 16](#)
- [Adjusting an Objective Correction Collar, see page 18](#)
- [Cleaning an Objective, see page 23](#)

You can identify the magnification of an objective by the color band:

Objective Magnification	Color Band
4x	Red
10x	Yellow
20x	Green
40x	Light Blue
63x	Dark Blue



Color Band on  
4x Objective

The standard objectives in the ImageXpress Pico system are configured and calibrated by Molecular Devices when the instrument is delivered.



**CAUTION!** Replace objectives in their original positions.



**Tip:** Molecular Devices recommends removing and maintaining only one objective at a time.

## Installing an Objective

Before installing an objective, review the following:

- Access only the user-serviceable components inside the enclosure as described in the procedure. Avoid contact with other components as they can be damaged or knocked out of alignment.
- To prevent dust from collecting inside the instrument, keep all access doors closed unless you are performing maintenance tasks.
- Ensure that all components and access doors are closed before starting the instrument.



### CAUTION!

- To prevent skin oils from damaging the optical coatings, Molecular Devices recommends that you wear powder-free disposable gloves when handling objectives and filter cubes.
- With the instrument power on, do not manually rotate the objective turret. Manually rotating the objective turret can damage the instrument.

Molecular Devices precalibrates the objectives to specific slots in the turret. You must install the objectives as follows:

Slot	Objective Magnification
1	4x
2	10x
3	20x
4	empty
5	40x or 63x
6	empty



### Note:




- Depending on how your ImageXpress Pico system is configured, you may not have all the objectives.
- The 40x objective and the 63x objective cannot be installed in the instrument simultaneously.
- The 63x objective cannot be installed when the environmental control cassette is loaded.



## Installing an Objective in an Empty Slot

You must install objectives in specific slots in the turret. See [Installing an Objective on page 12](#) for details.

To install an objective in an empty slot:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the instrument where you want to install an objective.
3. Click the **Objectives** tab.
4. In the tile for the objective slot where you want to install, click  **Component Exchange**.
5. Click the **Choose Objective** drop-down list box, and select the objective you want to install.
6. Click **Open Maintenance Door**.
7. Install the objective in the slot by gently turning it clockwise.



**CAUTION!** Retain the objective case for future storage needs. When not installed in the instrument, an objective should always be stored in its case.




8. Do both of the following:
  - a. Manually close the maintenance door.
  - b. In the CellReporterXpress software, click **Close Maintenance Door**.
9. Click **Apply**.

After you install a new objective from an after-sales purchase, you must calibrate it. See [Calibrating an Objective on page 16](#) for details.

## Replacing an Objective (Different Magnification)

You must install objectives in specific slots in the turret. See [Installing an Objective on page 12](#) for details. Because of this, the only reason to replace an objective with another objective of a different magnification is when you are swapping the 40x objective and the 63x objective in slot 5.

To replace an objective with an objective of a different magnification:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the instrument where you want to replace an objective.
3. Click the **Objectives** tab.
4. In the tile for objective slot **5**, click  **Component Exchange**.
5. Click the **Choose Objective** drop-down list box, and select the objective you want to install.
6. Click **Open Maintenance Door**.
7. Remove the objective to be replaced from the instrument by gently turning it counterclockwise.



**CAUTION!** When not installed in the instrument, an objective should always be stored in its case.

8. Install the replacement objective in the slot by gently turning it clockwise.



**CAUTION!** Retain the objective case for future storage needs. When not installed in the instrument, an objective should always be stored in its case.




9. Do both of the following:
  - a. Manually close the maintenance door.
  - b. In the CellReporterXpress software, click **Close Maintenance Door**.
10. Click **Apply**.

After you replace an objective, you may need to calibrate it. See [Calibrating an Objective on page 16](#) for details.

## Replacing an Objective (Same Magnification)

You must install objectives in specific slots in the turret. See [Installing an Objective on page 12](#) for details. The typical reason to replace an objective with another objective of the same magnification is when you are replacing a damaged objective.

To replace an objective with an objective of the same magnification:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the instrument where you want to replace an objective.
3. Click the **Objectives** tab.
4. In the tile for the objective slot where you want to install, click  **Component Exchange**.
5. Click the **Choose Objective** drop-down list box, and select **None**.
6. Click **Open Maintenance Door**.
7. Remove the objective to be replaced from the instrument by gently turning it counterclockwise.



**CAUTION!** When not installed in the instrument, an objective should always be stored in its case.

8. Install the replacement objective in the slot by gently turning it clockwise.



**CAUTION!** Retain the objective case for future storage needs. When not installed in the instrument, an objective should always be stored in its case.

9. Do both of the following:
  - a. Manually close the maintenance door.
  - b. In the CellReporterXpress software, click **Close Maintenance Door**.
10. Click **Close**.

After you replace an objective with another objective, you may need to calibrate it. See [Calibrating an Objective on page 16](#) for details.



**Note:** If you re-install the same objective (after cleaning it, for example), you typically do not need to calibrate it.

## Calibrating an Objective



After you install a new objective from an after-sales purchase, you must calibrate it. A calibration kit, which is included with any after-sales objective purchase, includes the following:

- Fluorescent Bead Slide (9100-0099)
- Red Plastic Slide (1-GP-7)
- Green Plastic Slide (1-GP-8)
- Orange Plastic Slide (1-GP-11)

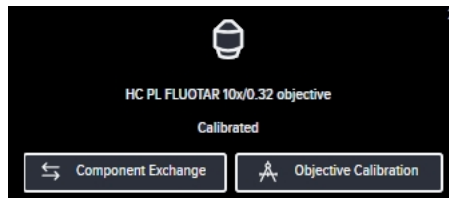
You will need to provide the following item for the calibration process:

- 4-Slide Holder (5068503)

To calibrate an objective:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the device where you want to calibrate an objective.
3. Click the **Objectives** tab.


4. Note the status of the objective, which is indicated in the tile.



- If the tile indicates **Not Calibrated**, continue to the next step to perform the calibration procedure.
- If the tile indicates **Calibrated**, you typically do not need to perform the calibration procedure and you can skip this procedure.



**Note:** An exception to this is when you replace an objective with another objective of the same magnification. In this case, you should continue to the next step and perform the calibration procedure regardless of the label in the tile.

5. In the tile for the objective you want to calibrate, click  **Objective Calibration**.
6. Follow the on-screen instructions to complete the calibration.



**Tip:**

- In the **Choose Slide Holder** drop-down list box, select **4 Slide Holder**.
- Insert the slides face down in the slide holder (that is, printed side down).
- When each step completes, click **Next** to continue to the next step.

## Adjusting an Objective Correction Collar

The 40x objective and 63x objective have application-optimized correction collars to compensate for well bottom thickness or coverslip thickness. The collars have a range of 0 mm to 2 mm correction. Changing this setting adjusts the distances between components inside the objective barrel. Image quality and resolution are very dependent on properly setting these collars.

Correction Collar  
on 40x Objective



The settings to be used depend on the well bottom thickness of the plate or the coverslip thickness on the slide on which the specimen is mounted. In general, set the correction collar for the physical thickness of the plate or slide that you are imaging. The physical thickness can be determined by the plate specifications from the plate manufacturer.



**Note:** Do not use a plate, slide, or coverslip with a thickness that is out of the range of the correction collar for the selected objective.

Before adjusting an objective correction collar, review the following:

- Access only the user-serviceable components inside the enclosure as described in the procedure. Avoid contact with other components as they can be damaged or knocked out of alignment.
- To prevent dust from collecting inside the instrument, keep all access doors closed unless you are performing maintenance tasks.
- Ensure that all components and access doors are closed before starting the instrument.




### CAUTION!

- To prevent skin oils from damaging the optical coatings, Molecular Devices recommends that you wear powder-free disposable gloves when handling objectives and filter cubes.
- With the instrument power on, do not manually rotate the objective turret. Manually rotating the objective turret can damage the instrument.

You would typically adjust a correction collar as part of setting up an acquisition.

## Adjusting an Objective Correction Collar for a Plate

To adjust an objective correction collar for a slide:

1. In the CellReporterXpress software, on the **Home** page, click  **Acquisition**.
2. Click **Add Protocol**.
3. Click **New Slide Acquisition**.
4. In the **Available Acquisition Devices** list, select the instrument.

5. On the left side of the screen under **Steps**, click  **Acquisition Settings**.

6. On the right side of the screen under **Tools**, click  **Slide Format**.

7. In the **Slide Format** list, select the slide format.

8. On the right side of the screen under **Tools**, click  **Objectives**.

9. In the **Objectives** list, select the objective.

If a correction collar adjustment is required, the software displays the recommended setting.

You may need to perform correction of objective collar for the current objective. Use value: 0.19



**Note:** When using the environmental control cassette with a 40x objective, you must adjust the objective correction collar an additional 0.7 mm above the recommended setting to account for the thickness of the cassette glass bottom.

10. On the left side of the screen under **Steps**, click  **Acquisition Device**.

11. On the right side of the screen, click  **Set Up for Adjustment of Objective Collar**.

12. Click **OK**.

The objective door opens.

13. If needed, loosen the objective from the instrument by gently turning it counterclockwise.

14. Rotate the correction collar to its new setting.



---

**Tip:** You might need a flashlight to see the markings for the graduated scale on the barrel and its current setting.

---

15. If you loosened the objective, tighten it by gently turning it clockwise.



---

**Note:** When tightening the objective, take care to avoid changing the correction collar setting.

---


16. Close the objective door.

17. Click **OK**.



## Adjusting an Objective Correction Collar for a Slide


To adjust an objective correction collar for a slide:

1. In the CellReporterXpress software, on the **Home** page, click  **Acquisition**.
2. Click **Add Protocol**.
3. Click **New Slide Acquisition**.
4. In the **Available Acquisition Devices** list, select the instrument.

5. On the left side of the screen under **Steps**, click  **Acquisition Settings**.

6. On the right side of the screen under **Tools**, click  **Slide Format**.

7. In the **Slide Format** list, select the slide format.

8. On the right side of the screen under **Tools**, click  **Objectives**.

9. In the **Objectives** list, select the objective.

If a correction collar adjustment is required, the software displays the recommended setting.

You may need to perform correction of objective collar for the current objective. Use value: 0.19

10. On the left side of the screen under **Steps**, click  **Acquisition Device**.

11. On the right side of the screen, click  **Set Up for Adjustment of Objective Collar**.

12. Click **OK**.

The objective door opens.

13. If needed, loosen the objective from the instrument by gently turning it counterclockwise.

14. Rotate the correction collar to its new setting.



---

**Tip:** You might need a flashlight to see the markings for the graduated scale on the barrel and its current setting.

---

15. If you loosened the objective, tighten it by gently turning it clockwise.



---

**Note:** When tightening the objective, take care to avoid changing the correction collar setting.

---

16. Close the objective door.

17. Click **OK**.

## Cleaning an Objective

If debris and contaminants collect on an objective lens, you can clean it. Before cleaning an objective, review the following:





- Access only the user-serviceable components inside the enclosure as described in the procedure. Avoid contact with other components as they can be damaged or knocked out of alignment.
  - To prevent dust from collecting inside the instrument, keep all access doors closed unless you are performing maintenance tasks.
  - Ensure that all components and access doors are closed before starting the instrument.
- 



### CAUTION!

- To prevent skin oils from damaging the optical coatings, Molecular Devices recommends that you wear powder-free disposable gloves when handling objectives and filter cubes.
  - With the instrument power on, do not manually rotate the objective turret. Manually rotating the objective turret can damage the instrument.
  - Do not use a product that disperses aerosol propellants or fluid onto the lens surface, such as canned compressed air.
  - Do not use a wipe that can leave lint, such as Kimwipes.
-

To clean an objective:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
  2. In the **Available Acquisition Devices** list, select the instrument.
  3. Click the **Objectives** tab.
  4. In the tile for the objective you want to clean, click  **Component Exchange**.
  5. Click the **Choose Objectives** drop-down list box, and select **None**.
  6. Click **Open Maintenance Door**.
  7. Remove the objective from the instrument by gently turning it counterclockwise.
  8. Place the objective on a secure surface away from the instrument
  9. Use a bulb duster to carefully blow dust contaminants off the objective.
  10. Use lens paper to gently wipe the objective free of contaminants.
- 
-  **Tip:** If needed, you can use a 100% methanol lens cleaner. Refer to Leica for details on preferred cleansing solvent and procedure.
- 
11. If needed, wait a few minutes for the methanol to evaporate.
  12. Install the objective back its original slot in the turret by gently turning it clockwise.
- 
-  **Tip:** If the objective has a correction collar, make sure that the collar is at the correct setting when reinstalling it.
- 
13. Do both of the following:
    - a. Manually close the maintenance door.
    - b. In the CellReporterXpress software, click **Close Maintenance Door**.
  14. Click **Close**.

After cleaning an objective, you may want to acquire a sample image. If image quality is degraded, repeat this procedure.

## Chapter 4: Filter Cube Maintenance

Filter cube maintenance procedures include the following:

- [Installing a Filter Cube, see below](#)
- [Calibrating a Filter Cube, see page 29](#)

### Installing a Filter Cube

Before installing a filter cube, review the following:

- Access only the user-serviceable components inside the enclosure as described in the procedure. Avoid contact with other components as they can be damaged or knocked out of alignment.
- To prevent dust from collecting inside the instrument, keep all access doors closed unless you are performing maintenance tasks.
- Ensure that all components and access doors are closed before starting the instrument.



**CAUTION!** To prevent skin oils from damaging the optical coatings, Molecular Devices recommends that you wear powder-free disposable gloves when handling objectives and filter cubes.

The following filter cubes are available:

- DAPI
- FITC
- TRITC
- Cy5
- Texas Red
- CFP






**Note:**

- Depending on how you have configured your ImageXpress Pico system, you may not have received all the filter cubes.
- Do not install a filter cube in slot 6 of the turret. Slot 6 must be empty (in both the instrument and the software).

## Installing a Filter Cube in an Empty Slot

To install a filter cube in an empty slot:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the device where you want to install a filter cube.
3. Click the **Filters** tab.
4. In the tile for the filter cube slot where you want to install, click  **Component Exchange**.



**Note:** Do not install a filter cube in slot 6 of the turret. Slot 6 must be empty (in both the instrument and the software).

5. Click the **Choose Filter** drop-down list box, and select the filter cube you want to install.
6. Click **Open Maintenance Door**.
7. If needed, slightly rotate the filter cube turret by hand to get direct access to the filter cube slot.
8. Install the filter cube in the slot by gently pushing it into the slot until it “snaps” into place.






**CAUTION!** Retain the filter cube packaging for future storage needs. When not installed in the instrument, a filter cube should always be stored in its original packaging.

9. Do both of the following:
  - a. Manually close the maintenance door.
  - b. In the CellReporterXpress software, click **Close Maintenance Door**.
10. Click **Apply**.

After you install a new filter cube from an after-sales purchase, you must calibrate it. See [Calibrating a Filter Cube on page 29](#) for details.

## Replacing a Filter Cube (Different Wavelength)

To replace a filter cube with a filter cube of a different wavelength:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the device where you want to replace a filter cube.
3. Click the **Filters** tab.
4. In the tile for the filter cube slot where you want to install, click  **Component Exchange**.
5. Click the **Choose Filter** drop-down list box, and select the filter cube you want to install.
6. Click **Open Maintenance Door**.
7. If needed, slightly rotate the filter cube turret by hand to get direct access to the filter cube slot.
8. Remove the filter cube to be replaced from the instrument by gently pulling it toward you.



**CAUTION!** When not installed in the instrument, a filter cube should always be stored in its original packaging.

9. Install the filter cube in the slot by gently pushing it into the slot until it “snaps” into place.



**CAUTION!** Retain the filter cube packaging for future storage needs. When not installed in the instrument, a filter cube should always be stored in its original packaging.

10. Do both of the following:
  - a. Manually close the maintenance door.
  - b. In the CellReporterXpress software, click **Close Maintenance Door**.
11. Click **Apply**.

After you replace a filter cube, you may need to calibrate it. See [Calibrating a Filter Cube on page 29](#) for details.






**Note:** If you re-install the same filter cube (after cleaning it, for example), you do not need to calibrate it.

## Replacing a Filter Cube (Same Wavelength)

The typical reason to replace a filter cube with another filter cube of the same wavelength is when you are replacing a damaged filter cube.

To replace a filter cube with a filter cube of the same wavelength:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the device where you want to replace a filter cube.
3. Click the **Filters** tab.
4. In the tile for the filter cube slot where you want to replace, click  **Component Exchange**.
5. Click the **Choose Filter** drop-down list box, and select **None**.
6. Click **Open Maintenance Door**.
7. If needed, slightly rotate the filter cube turret by hand to get direct access to the filter cube slot.
8. Remove the filter cube to be replaced from the instrument by gently pulling it toward you.



**CAUTION!** When not installed in the instrument, a filter cube should always be stored in its original packaging.

9. Install the filter cube in the slot by gently pushing it into the slot until it “snaps” into place.



**CAUTION!** Retain the filter cube packaging for future storage needs. When not installed in the instrument, a filter cube should always be stored in its original packaging.

10. Do both of the following:
  - a. Manually close the maintenance door.
  - b. In the CellReporterXpress software, click **Close Maintenance Door**.
11. Click **Close**.

After you replace a filter cube, you may need to calibrate it. See [Calibrating a Filter Cube on page 29](#) for details.



**Note:** If you re-install the same filter cube (after cleaning it, for example), you do not need to calibrate it.



## Calibrating a Filter Cube



After you install a new filter cube from an after-sales purchase, you must calibrate it. A calibration kit, which is included with any after-sales filter cube purchase, includes the following:

- Fluorescent Bead Slide (9100-0099)
- Red Plastic Slide (1-GP-7)
- Green Plastic Slide (1-GP-8)
- Orange Plastic Slide (1-GP-11)

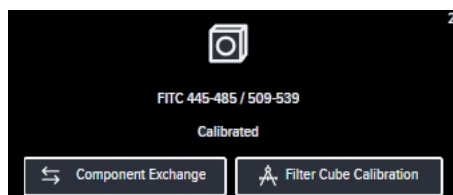
You will need to provide the following item for the calibration process:

- 4-Slide Holder (5068503)

To calibrate a filter cube:

1. In the CellReporterXpress software, on the **Home** page, click  **Devices**.
2. Click  **Show Device Options** to expand the details for the device where you want to calibrate a filter cube.
3. Click the **Filters** tab.


4. Note the status of the filter cube, which is indicated in the tile.



- If the tile indicates **Not Calibrated**, continue to the next step to perform the calibration procedure.
- If the tile indicates **Calibrated**, you typically do not need to perform the calibration procedure and you can skip this procedure.



**Note:** An exception to this is when you replace a filter cube with the same type of filter cube. In this case, you should continue to the next step and perform the calibration procedure regardless of the label in the tile.

5. In the tile for the filter cube you want to calibrate, click  **Filter Cube Calibration**.
6. Follow the on-screen instructions to complete the calibration.



**Tip:**

- In the **Choose Slide Holder** drop-down list box, select **4 Slide Holder**.
- Insert the slides face down in the slide holder (that is, printed side down).
- When each step completes, click **Next** to continue to the next step.

## Appendix A: Optional Parts and Accessories

# A

This section describes the items available from Molecular Devices for the ImageXpress Pico system, including the following:

- Objectives, see below
- Filter Cubes, see below
- Accessories, see page 32

See [Obtaining Support on page 6](#) for details on contacting Molecular Devices.

### Objectives

The following Leica objectives, which are available from Molecular Devices, are compatible with the ImageXpress Pico system:

Objective	Part Number	Mag.	Numerical Aperture (NA)	Working Distance	Corr. Collar
PL FLUOTAR 4x/0.13	5064345	4x	0.13	17.0 mm	No
HC PL FLUOTAR 10x/0.32	5064346	10x	0.32	11.13 mm	No
HC PL FLUOTAR 20x/0.40	5064347	20x	0.4	7.45 mm	No
HC PL FLUOTAR L 40x/0.60 CORR	5064348	40x	0.6	3.0 mm	Yes
HC PL FLUOTAR L 63x/0.70 CORR	5064349	63x	0.7	2.0 mm	Yes



**CAUTION!** To prevent damaging both the instrument and your sample, do not use any other objectives with the ImageXpress Pico system.

### Filter Cubes

The following filter cubes, which are available from Molecular Devices, are compatible with the ImageXpress Pico system:

Filter	Part Number	Excitation	Emission	Dichroic
DAPI	5064350	370/40 nm	450/60 nm	410 nm
FITC	5064351	465/40 nm	525/30 nm	500 nm
TRITC	5064352	530/45 nm	594/40 nm	560 nm
Cy5	5064353	630/40 nm	695/45 nm	655 nm
Texas Red	5069939	560/50 nm	645/75 nm	595 nm
CFP	5069940	400/30 nm	480/40 nm	475 nm

## Accessories

The following accessories, which are available from Molecular Devices, are compatible with the ImageXpress Pico system.

### Base System Accessories

Item	Part No.
4-Slide Holder	5068503
Plate Holder	5068504
6-Dish Holder for 35 mm Dishes	5077007
Calibration Kit	5068505
Two (2) CAT6 Ethernet Cables, 2 m (6.6 ft)	5052189
Power Cord for USA/Canada, 2.29 m (7.5 ft)	4400-0002
Power Cord for Europe, 1.96 m (6.4 ft)	4400-0036
Power Cord for China, 2.50 m (8.2 ft)	4400-0276

### Environmental Control System Accessories

Item	Part No.
Environmental Control Cassette	5070105
Humidifying Column	5070110
Humidifying Column Tubing/Wiring	5070108
Two (2) Humidifying Column Stoppers	5070107
Gas Supply Tubing, 10 m (32.8 ft)	5070103
Push-to-Connect Straight Fitting 6 mm O.D. - ¼" NPT Male	5075610
Two (2) Blind Plugs	5070106
Three (3) Plate Skirt Height Adapters	5077006

## Contact Us

Phone: [+1-800-635-5577](tel:+1-800-635-5577)  
Web: [moleculardevices.com](http://moleculardevices.com)  
Email: [info@moldev.com](mailto:info@moldev.com)

Visit our website for a current listing of worldwide distributors.

The trademarks used herein are the property of Molecular Devices, LLC or their respective owners.  
Specifications subject to change without notice. Patents: [www.moleculardevices.com/patents](http://www.moleculardevices.com/patents)  
FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES

©2019 Molecular Devices, LLC.  
All rights reserved.  
5063669 E

