



Installing an Epi Light

Note: If you are working with a new instrument, see the installation guide for the instrument and follow the instructions to install Epi lights. If the instrument was previously installed, follow the steps given here.

Before You Begin

Locate the new filters that came with the Epi lights and the focus calibration target and flat field plates that came with the original ChemiDoc™ MP system. You will work with these items when the New Lights Found wizard runs.

Install the Epi Light Modules

Epi lights are available for the ChemiDoc MP system in three colors: red, blue, and green. The following instructions demonstrate installing a red Epi light module.

Note: Use these same instructions to install a blue or green module in its place in the Epi light bracket assembly.

Each kit of a given color consists of a pair of modules that plug in to Epi lights assemblies inside the Universal Hood. These instructions show how to install a module into the left side Epi light bracket assembly.

Each module has two guide tabs on its back and two screws extending from its front. Each module is labeled with its color. A contact tab extends from the top of each module.



Important: Handle the module only by the black module body. Do not touch the contact tabs or the optical filter.

1. Power down the system using the power switch on the front of the device.



2. Open the Universal Hood door. Inside, an Epi lights bracket assembly is located on each side of the Universal Hood. The following image shows the left side Epi lights bracket assembly.

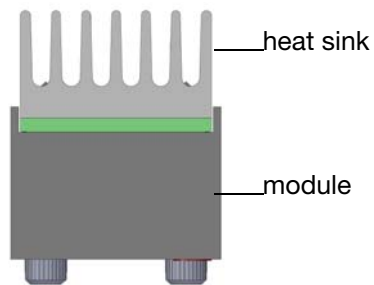


3. Match the module label with the same color label on the Epi lights bracket assembly.



4. Position the module, contact tab up, against the bracket assembly so that the guide tabs are on the sides of the heat sink and the module cannot move sideways. Make sure the screws extend as shown in the picture and the module is flat against the heat sink.

- Carefully and slowly slide the module up as far as it will go so that the contact tab is no longer visible. Doing so lines up the screws with the appropriate holes in the bracket assembly.



- Tighten the two screws. A screwdriver may be used. Do not overtighten.



- Repeat this procedure to install the second Epi light module in the Epi lights bracket assembly on the opposite side of the Universal Hood.

Image Lab Software Detects the Lights

- Turn on the ChemiDoc MP instrument.
- If Image Lab™ software is running, it will detect the new light and display the message “New lights found.” If Image Lab software is not running, start it and it will display the same message.
- Click Next.

The New Lights Found Wizard Runs

A series of screens shows you how to install the new filters and guides you through focus calibration, flat field calibration, and bias acquisition. To insert filters as requested in the wizard, see the procedure on the next page.

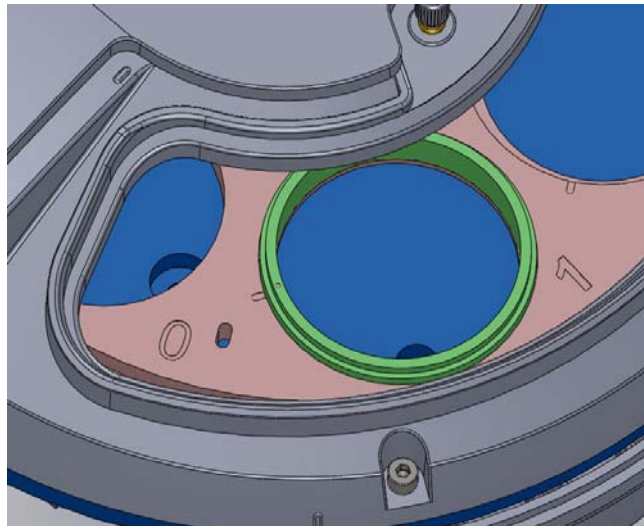
Insert a Filter into the Filter Wheel

Each filter has a white mark on its rim.



Each filter holder is numbered.
The number to the right of the
filter indicates the filter position.

Insert a filter into the filter wheel
as the wizard instructs and line
up the white mark on the filter
with the notch on the filter holder.



To ensure that the filter is seated properly,
press it against the gray bumper, then
press the filter down into the filter holder.

