



Quick Start Protocol

Glo-Plate™ 2.0 Blue LED Illuminator



Catalog Number: E90007

A quick demonstration of the device can be found here: <https://biotium.wistia.com/medias/njthreyvk3>

Protocol for using Glo-Plate™ 2.0 in viability PCR

- 1. Turn the Glo-Plate™ 2.0 on.**
The power switch is on the back of the device. The main menu will appear (Figure 1).
- 2. Press  to bring up the time setting screen (Figure 2) and set the length of time for light exposure.**
We recommend 20 minutes of light exposure as a starting point. PMA-treated samples may require a longer exposure than PMAxx™-treated samples.
- 3. Press  to return to the home screen.**
- 4. Place samples on the Glo-Plate™ 2.0.**
Place a sheet of aluminum foil over the samples with the reflective side facing down toward the light.
- 5. Press Start.**
The timer will count up to the time set in step 2. Leave samples on the Glo-Plate™ 2.0 for the set length of time. Once light exposure has begun, the Start button will become a Pause button.
- 6. After time expires, the device will beep and the LEDs will shut off.**
To begin the same light exposure cycle on fresh samples, press Reset, place fresh samples on the Glo-Plate™ 2.0, and press Start. Otherwise, turn the device off and remove samples.

Protocol for using Glo-Plate™ 2.0 with DNAzure®-stained agarose gels

- 1. Stain agarose gel with 1X DNAzure® for 20-30 minutes.**
Use a clear container for staining.
- 2. Turn the Glo-Plate™ 2.0 on.**
The power switch is on the back of the device. The main menu will appear (Figure 1).
- 3. Press  to bring up the time setting screen (Figure 2) and set the length of time for light exposure.**
We recommend 15-30 minutes of light exposure.
- 4. Press  to return to the home screen.**
- 5. Place the gel on the Glo-Plate™ 2.0.**
Place the container with the gel on the Glo-Plate™ 2.0, or remove the gel from the container and place it on clear plastic wrap directly on the light panel. Place a sheet of aluminum foil over the samples with the reflective side facing down toward the light.

6. Press Start.

The timer will count up to the time set in step 2. Leave the gel on the Glo-Plate™ 2.0 for the set length of time. Once light exposure has begun, the Start button will become a Pause button.

7. After time expires, the device will beep and the LEDs will shut off.

If additional exposure time is desired, press Reset and then Start to repeat the same light exposure cycles, or select a new time. Otherwise, turn the device off and remove the gel.

Safety Information

Plug the power cord into a properly grounded outlet. Disconnect the power cord from the electrical outlet if the device will not be used for a long period. Disconnect the device from the electrical outlet before cleaning it. The device is not water-proof; do not submerge in water. To clean the stage, wipe down the surface with a small amount of 10% bleach or 70% ethanol, then wipe with water. Do not soak. Do not look directly at the blue LEDs for a prolonged period of time. Bright blue light is a possible risk factor for macular degeneration. Do not use sharp items when using the touch screen to avoid damaging the screen.

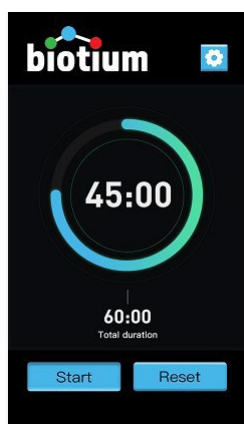


Figure 1. Example Glo-Plate™ 2.0 main menu display screen.

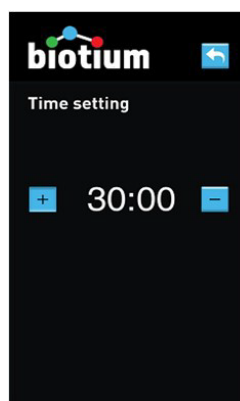


Figure 2. Glo-Plate™ 2.0 time settings screen.

Visit www.biotium.com for complete protocols, troubleshooting tips, and answers to frequently asked questions (FAQs).

For research use only.