



Quick Start Protocol

PMA-Lite™ 2.0 LED Photolysis Device

Catalog Number: E90006

Protocol for using PMA-Lite™ 2.0 in viability PCR

- 1. Turn the PMA-Lite™ 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 15 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 in the PMA-Lite™ 2.0.**
Samples must be in clear 1.7 mL microcentrifuge or 2 mL screw cap tubes.
- 5. Press Start.**
The timer will count up to the time set in step 2. Leave samples in the PMA-Lite™ 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 LEDs will shut off.**
To begin the same light exposure cycle on fresh samples, press Reset, place fresh samples in the PMA-Lite™ 2.0, and press Start. Otherwise, turn the device off and remove samples.

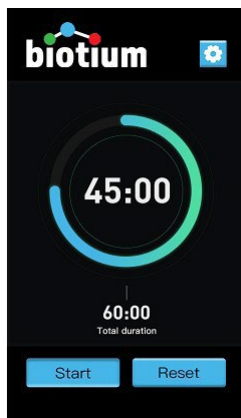


Figure 1. Example PMA-Lite™ 2.0 main menu display screen.

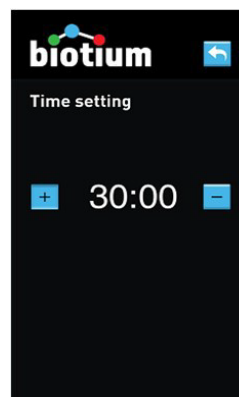


Figure 2. PMA-Lite™ 2.0 time settings screen.

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

For research use only.