



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

## Glo-Plate™ 2.0 Blue LED Illuminator

A multi-functional blue LED light box that can be used for photolysis of PMAxx™ or PMA in viability PCR; or to develop the visible DNA gel stain DNAzure®.



### Product attributes

## Product Description

The Glo-Plate™ 2.0 Blue LED Illuminator is a lightweight, multifunctional LED light box that offers even illumination for samples in clear microplates, larger tubes, and other transparent vessels. The LEDs in the Glo-Plate™ 2.0 were designed for optimal photolysis of [PMAxx™](#)– or [PMA](#)-treated samples in viability PCR (v-PCR) applications.

### Uniform Photoactivation of v-PCR Samples in Microplates

Viability PCR with dyes such as [PMAxx™](#) or [PMA](#) requires consistent and maximal photoactivation between samples to produce accurate results. While the [PMA-Lite 2.0™ LED Photolysis Device](#) can be used for this application, the device can only illuminate up to 18 microcentrifuge tubes at a time, leaving little options for researchers processing larger sample sizes in microplates. The Glo-Plate™ 2.0 Blue LED Illuminator offers a flat surface illuminated by LED lights topped with a diffuser, allowing for uniform and optimal photoactivation of [PMAxx™](#)– or [PMA](#)-treated samples in clear microplates. The device may also be used for photoactivation of samples in 15 mL or 50 mL tubes, and other transparent vessels.

### Excellent Light Source for DNAzure®

The Glo-Plate™ 2.0 Blue LED Illuminator is also an excellent light source for developing gels stained with [DNAzure® Blue Nucleic Acid Gel Stain](#). DNAzure® is a novel DNA gel stain that precipitates a visible blue dye onto DNA during light exposure, resulting in highly sensitive and stable visible blue DNA bands that can be visualized by eye. This is particularly advantageous for gel excision.

### New 2.0 Design Includes User-Friendly Touch Screen Interface!

The Glo-Plate™ 2.0 Blue LED Illuminator is a new and improved version of the discontinued Glo-Plate™ Blue LED Illuminator (Cat. No. E90004). The new 2.0 design offers the same robust and uniform photoactivation performance as the original Glo-Plate™ Blue LED Illuminator, while also featuring a convenient touch display that allows users to easily set, reset, and monitor the device's running time.

To learn more about the advantages of determining microbial or cell viability using viability PCR, visit the [Viability PCR Technology Page](#).

This datasheet was generated on January 8, 2026 at 06:21:30 AM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/glo-plate-2-0-blue-led-illuminator/>