



Glowing products for science

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

Gel-Bright™ Laser Diode Gel Illuminator

A laser diode-based gel illuminator for visualizing fluorescent nucleic acid and protein gels. Avoids UV light hazards and offers exceptional performance for widely used green and red dyes.



Product Description

The Gel-Bright™ Laser Diode Gel Illuminator features a novel laser diode (LD)-based illumination technology that offers comparable performance to UV transilluminators while avoiding the hazards of UV light which can damage skin and eyes. UV light can also induce cross-linking and nicking of DNA in samples. Unlike blue LED gel illuminators such as Dark Reader or Safe Imager™ that are not compatible with red dyes, the Gel-Bright™ Laser Diode Gel Illuminator works well with both green dyes and widely used red dyes like GelRed®, ethidium bromide (EtBr), and One-Step Lumitein™ Protein Gel Stain.

Superior Performance Over Blue LED Illuminators

Gel illuminators have traditionally relied on UV illumination for visualization of fluorescently labeled nucleic acids and proteins. However, UV-based imaging can damage skin and eyes, as well as your DNA samples. LED-based illuminators were developed as a safer alternative to UV, but often have high background and dimmer signal due to excessive ambient light and poor excitation efficiency.

In partnership with Biotium, OMEC Medical has developed a new type of gel illuminator that uses LDs optimized for brighter and clearer fluorescent imaging of gels. The new laser diode-based Gel-Bright™ offers better sensitivity over UV-based transilluminators when imaging green dyes such as GelGreen® or SYBR® Green, as well as significantly better performance over blue LED gel illuminators for red dyes such as GelRed®, ethidium bromide (EtBr), and One-Step Lumitein™ Protein Gel Stain.

Sensitive for Red Dyes

- GelRed®
- PAGE GelRed®
- One-Step Lumitein™
- SYPRO® Ruby
- EtBr

Free Samples of Pre-Coated GelRed® & GelGreen® Agarose Included

Each purchase of a Gel-Bright™ Laser Diode Gel Illuminator will include a free sample of [GelRed® Agarose LE \(Cat. No. 41029-5G\)](#) and [GelGreen® Agarose LE \(Cat. No. 41030-5G\)](#). These are safer and more convenient pre-coated agarose formats that remove the extra step of adding dye to your agarose.

Product attributes

Safe Imager™ vs. Gel-Bright™ Laser Diode Gel Illuminator

Old Gel-Bright™ vs. Gel-Bright™ Laser Diode Gel Illuminator

Sensitive Gel Imaging Without UV Hazards

Compatible Dyes:

- [GelGreen® and GelRed® Nucleic Acid Gel Stains](#)
- [EMBER™ Ultra RNA Gel Kit](#)
- [Ethidium Bromide \(EtBr\)](#)
- [EvaGreen® Dye \(Biotium\) \(direct detection of EvaGreen® PCR products\)](#)
- [One-Step Lumitein™ and SYPRO® Ruby Protein Gel Stains](#)
- [SYBR® Green \(Thiazole Green\)](#), [SYBR® Gold \(Oxazole Gold\)](#), SYBR® Safe, and similar dyes

Learn more about our sensitive and less hazardous [GelRed® and GelGreen® Nucleic Acid Gel Stains](#) as well as our [One-Step protein gel stains](#).

GelDoc-iT is a registered trademark of UVP/Analytik Jena GmbH; SYBR and SYPRO are registered trademarks of Thermo Fisher Scientific; Safe Imager is a trademark of Thermo Fisher Scientific.