

# Quick Start Protocol

## Gel-Bright™ Laser Diode Gel Illuminator

Catalog Number: E90005

### Protocol for using Gel-Bright™

1. **Run a gel and stain it with a fluorescent stain.**
2. **Place the glass plate into the gel staging area and place the gel on the glass plate (Figure 1).**
3. **Turn the Gel-Bright™ on.**  
The power switch is on the side of the device.
4. **Turn the Light Intensity Dial (Figure 1) clockwise until the desired light intensity is attained.**
5. **View and photograph the gel through the amber filter.**  
To adjust the filter position, slide the Filter Adjustment Slider (Figure 1) on the side of the device up towards the filter, then lower the filter. For the best results, use the Gel-Bright™ Imaging Hood by placing it over the filter while it's in the lowered position. Image the gel through the small opening at the top of the imaging hood.
6. **Excise bands from the gel, if necessary.**  
The filter can be angled up or the device can be turned around for better access while excising gel slices (Figure 2).
7. **Turn the device off and remove the gel.**  
Turn the Light Intensity Dial counter-clockwise until the lasers turn off and press the power switch to off.

### Safety Information

Plug the power cord into a properly grounded outlet. Disconnect the device from the electrical outlet before cleaning it. The device is not water-proof; do not submerge in water. Do not add any liquids to the gel staging area. The device contains class 3b laser diodes; safety glasses are not required during routine use because the amount of laser light that could enter the eye's pupil is below the CDRH limits. Avoid looking directly into the laser diodes when the filter is lifted. When excising gel slices, ALWAYS position the glass plate under the gel as sharp blades will scratch the lower surface of the gel staging area.

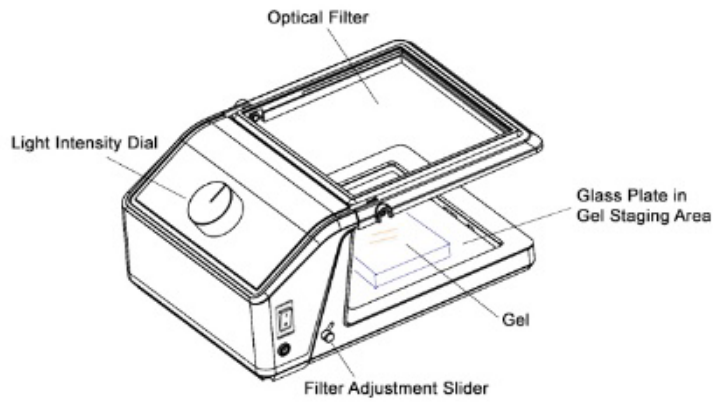


Figure 1. Diagram of Gel-Bright™ device.



Figure 2. Gel-Bright™ device with filter set to different positions.

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

For research use only.