

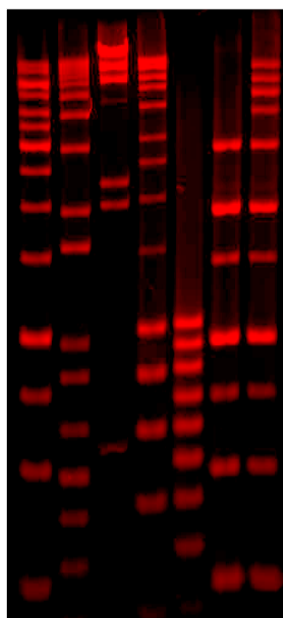
GelRed™ and GelGreen™

Nucleic Acid Gel Prestaining Kits

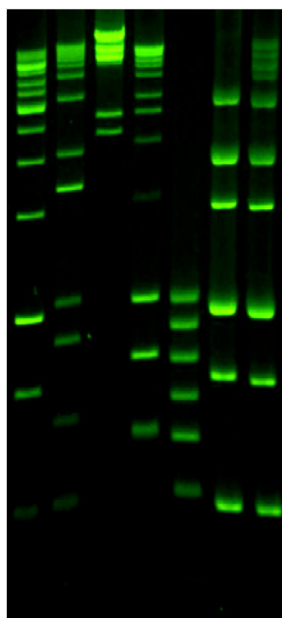
— The safest and simplest way to stain nucleic acids in gels with amazing results!



GelRed™



GelGreen™



Gel electrophoresis of DNA ladders from seven different vendors following prestaining with GelRed™ or GelGreen™ loading buffer.

Nucleic acid gel stains are inherently dangerous because of their potential to cause genetic mutation. Few would dispute that one of the best ways to reduce or eliminate their toxicity is to prevent the dyes from entering the cells in the first place. But the challenge is: How do you do it without compromising performance? The answer is innovation. GelRed™ and GelGreen™ dyes are breakthrough nucleic acid gel stains and the only gel stains that offer the combination of: 1) cell membrane impermeability for true safety;* 2) unrivaled sensitivity; 3) perfect instrument compatibility; 4) superior stability; and 5) compatibility with all downstream manipulations.

Continuing our innovation, now we have made GelRed™ and GelGreen™ dyes even better by introducing yet another breakthrough – GelRed™ and GelGreen™ Prestaining Kits for ultimate safety, performance and value. Each kit contains an optimally formulated 6X GelRed™ or GelGreen™ loading buffer and a superior proprietary running buffer. With these new kits, you enjoy the following additional benefits:

- **Ultimate Safety:** Absolutely minimal exposure to a dye that already has an excellent safety profile.
- **No More Migration Problem:** Work well with all commercial DNA/RNA ladders with excellent migration and resolution.
- **Fast Gel Electrophoresis:** Run your gels at twice the speed in half the time of standard buffers. A running buffer concentrate is provided for convenient storage.
- **Extremely Simple Procedure:** Simply load your sample in the GelRed™ or GelGreen™ loading buffer and run your gel in the proprietary running buffer provided in the kit. No additional dye in the gel or running buffer and no post staining is necessary.
- **Re-usability:** Gels may be re-used if there are empty wells.
- **Easy Disposal:** Waste gels can be combined with regular trash and running buffer can be disposed of down the drain.
- **Unmatched Value:** Enough loading buffer and running buffer are provided for analysis of up to 100 mini gels.

* A complete safety report can be downloaded from the Biotium website.

** GelRed, GelGreen and their uses are covered by US patent No. 7,601,498 and other pending US and international patents; GelRed and GelGreen are trademarks of Biotium, Inc.

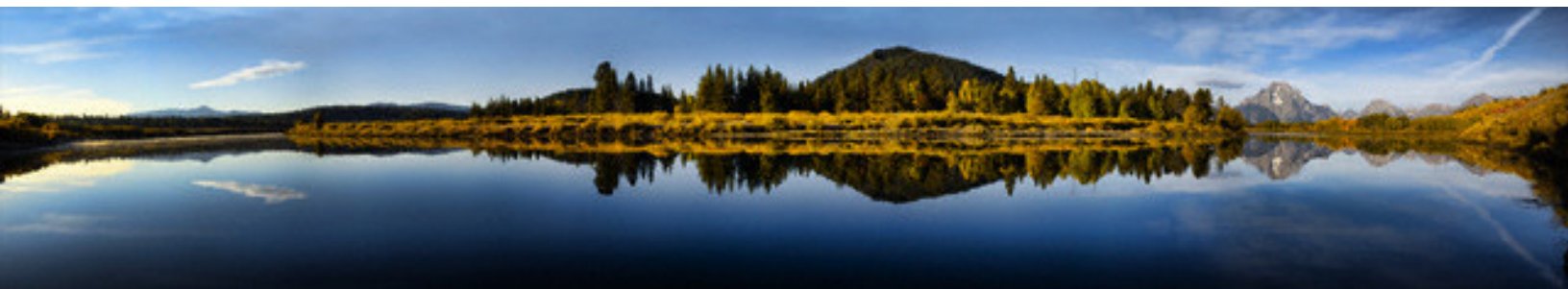


Table 1. Product List (See Selection Guide in Table 2)

Catalog #	Product Name	Unit Size
31010	GelRed™ nucleic acid gel prestaining kit with "200bp" and "1.5Kb" tracking dyes	Kit (sufficient for 100 mini gels)
31011	GelRed™ nucleic acid gel prestaining kit with "50bp" tracking dye	Kit (sufficient for 100 mini gels)
31012	GelGreen™ nucleic acid gel prestaining kit with "200bp" and "1.5Kb" tracking dyes	Kit (sufficient for 100 mini gels)
31013	GelGreen™ nucleic acid gel prestaining kit with "low molecular weight" tracking dye ("Coming soon!")	Kit (sufficient for 100 mini gels)

Table 2. Selection Guide for GelRed™/GelGreen™ Prestaining Kits

If You Have the Following Condition		Select This Kit
Medium to large DNA fragments	UV(302 nm) excitation; EtBr filter	Cat# 31010, GelRed™ nucleic acid gel prestaining kit with "200bp" and "1.5Kb" tracking dyes
	UV(254nm), blue light(470 nm) or 488 nm laser excitation; green filter	Cat# 31012, GelGreen™ nucleic acid gel prestaining kit with "200bp" and "1.5Kb" tracking dyes
Small DNA/RNA fragments; PAGE gels	UV(302 nm) excitation; EtBr filter	Cat# 31011, GelRed™ nucleic acid gel prestaining kit with "50bp" tracking dye
	UV(254nm), blue light(470 nm) or 488 nm laser excitation; green filter	Cat# 31013, GelGreen™ nucleic acid gel prestaining kit with "low molecular weight" tracking dye



Biotium, Inc.
 www.biotium.com
 Tel. 1-800-304-5357
 3159 Corporate Place, Hayward, CA 94545