

## **Dye Advantages**

- Non-toxic and non-mutagenic
- Safer than EtBr and other so-called safe gel stain
- Ultra-sensitive and stable
- Simple to use
- Easy disposal, okay for drain disposal by EPA Title 22 hazardous waste test
- Compatible with downstream applications, such as cloning and sequencing

# GelRed<sup>®</sup> and GelGreen<sup>®</sup>

# The most safe and sensitive nucleic acid gel stains

GelRed® and GelGreen® are safe nucleic acid gel stains designed to replace highly toxic ethidium bromide (EtBr) and other so-called safe gel stains. These gel stains are non-toxic and non-mutagenic by virtue of being cell membrane impermeable. Ames tests have confirmed that GelRed® and GelGreen® are nonmutagenic at concentrations well above the concentrations used for gel staining. Furthermore, environmental safety tests showed that GelRed® and GelGreen® are non-toxic to aquatic life, permitting disposal down the drain or in regular trash.

For more information and references, download our white paper, Comparison of Nucleic Acid Gel Stains: Cell Permeability, Safety, and Sensitivity and the complete Safety Report of GelRed® and GelGreen® at www.biotium.com.



Figure 1. GelRed® and GelGreen® gel stains are safer because they cannot penetrate cell membranes to bind DNA in living cells. HeLa cells were incubated with 1X SYBR® Safe, GelGreen® or GelRed®, respectively. Images were taken following incubation with dye for 30 min using FITC filter set for SYBR® Safe and GelGreen®, and Cy®3 filter set for GelRed®. SYBR® Safe rapidly entered cells and stained nuclei. GelRed® and GelGreen® were unable to cross cell membranes, demonstrated by the absence of fluorescence staining.

EtBr	GelRed®	SYBR <sup>®</sup> Safe	GelGreen®
		-	
		and drag trees on the	
			and 1000 1000
			and set

Figure 2. GelRed® and GelGreen® are more sensitive than EtBr and SYBR® Safe. Left: Comparison of GelRed® and ethidium bromide (EtBr) in precast gel staining using 1% agarose gel in TBE buffer. Right: Comparison of GelGreen® and SYBR® Safe in post gel staining using 1% agarose gel in TBE buffer.



# GelRed® products for greater convenience and safety

### GelRed® Prestain Plus 6X DNA Loading Dye

- 6X loading dye that includes GelRed® DNA stain
- · Combine gel loading and DNA staining in one step
- Two blue tracking dyes that run at ~1.5 kb and ~200 bp in 1% agarose

### GelRed® Agarose LE

- · Agarose that is pre-coated with GelRed® dye
- Simplies the preparation of precast gels
- · No need to handle concentrated dye solutions, for enhanced safety
- · Use with TAE or TBE for 0.8% to 2% gels
- · Low EEO, ultrapure molecular biology grade agarose

### 3X GelRed® in Water

- Ready-to-use for post-electrophoresis gel staining
- Supplied in a 4L Cubitainer®
- · No need to handle concentrated dye solutions, for enhanced safety



GelRed® Agarose LE



3X GelRed® in Water

### **Ordering Information**

Cat. #	Product Name
41003-T	GelRed® Nucleic Acid Gel Stain; 10,000X in water, 0.1 mL
41003	GelRed® Nucleic Acid Gel Stain; 10,000X in water, 0.5 mL
41003-1	GelRed® Nucleic Acid Gel Stain; 10,000X in water, 10 mL
41002	GelRed® Nucleic Acid Gel Stain; 10,000X in DMSO, 0.5 mL
41002-1	GelRed® Nucleic Acid Gel Stain; 10,000X in DMSO, 10 mL
41001	GelRed® Nucleic Acid Gel Stain; 3X in water, 4 L
41029-5G	GelRed® Agarose LE, 5 g
41029-50G	GelRed® Agarose LE, 50 g
41011	GelRed® Prestain Plus 6X DNA Loading Dye, 1 mL
41005-T	GelGreen® Nucleic Acid Gel Stain; 10,000X in water, 0.1 mL
41005	GelGreen® Nucleic Acid Gel Stain; 10,000X in water, 0.5 mL
41005-1	GelGreen® Nucleic Acid Gel Stain; 10,000X in water, 10 mL
41004	GelGreen® Nucleic Acid Gel Stain; 10,000X in DMSO, 0.5 mL
41030-5G	GelGreen® Agarose LE, 5 g
41030-50G	GelGreen® Agarose LE, 50 g
41008-T	PAGE GelRed® Nucleic Acid Gel Stain; 10,000X in water, 0.1 mL
41008-500uL	PAGE GelRed® Nucleic Acid Gel Stain; 10,000X in water, 0.5 mL
41014	PAGE GelRed® Nucleic Acid Gel Stain; 1X in water, 4 L

# Nucleic acid gel stain for polyacrylamide gels

The safety profile of the original GelRed® dye is in part due its large size, which makes the dye impermeable to cell membranes. However the large dye size reduces penetration of GelRed® into densely packed polyacrylamide gels. In designing PAGE GelRed® dye, we used a novel approach to make the dye membrane impermeable without making the dye large.

### PAGE GelRed® Nucleic Acid Gel Stain

- A safe and sensitive gel stain for polyacrylamide gels
- Non-toxic and non-mutagenic in AMES test
- Formulated in water and impermeable to latex and nitrile gloves
- Non-toxic to aquatic life, okay for drain disposal by EPA Title 22 hazardous waste test

GelGreen and GelRed are registered trademarks of Biotium, Inc. GelGreen, GelRed, PAGE GelRed and their uses are covered by US and international patents. SYBR is a registered trademark of Thermo Fisher Scientific. v02.04.19 www.biotium.com General Inquiries: order@biotium.com Technical Support: techsupport@biotium.com Phone: 800-304-5357

