Glowing products for science

## GelRed® Nucleic Acid Gel Stain

Ultra sensitive and environmentally safe replacement for EtBr DNA/RNA gel stain.



#### **Product attributes**

Storage Conditions Store at room temperature, Protect from light

Call us: 800-304-5357 Email: btinfo@biotium.com

Assay type/options DNA/RNA gel staining

# **Product Description**

GelRed® is an ultra sensitive, extremely stable and environmentally safe fluorescent nucleic acid dye designed to replace the highly toxic ethidium bromide (EtBr) for staining dsDNA, ssDNA or RNA in agarose gels or polyacrylamide gels.

- Safer than EtBr: non-mutagenic and non-hazardous for disposal
- Much more sensitive than EtBr and SYBR® Safe
- Stable at room temperature and microwavable
- Simple precast or post-electrophoresis gel staining, no destaining
- Replaces EtBr with no change to equipment or optical settings
- Compatible with downstream gel purification, restriction digest, sequencing, and cloning

### GelRed®: A Superior Alternative to EtBr

GelRed® is much more sensitive than EtBr, and can be used to stain dsDNA, ssDNA or RNA in agarose gels via either precast or post gel staining without destaining. GelRed® can also be used to stain polyacrylamide gels via post gel staining. GelRed® is also compatible with downstream DNA manipulations such as restriction digest, sequencing, and cloning. GelRed® and EtBr have virtually the same spectra, so you can directly replace EtBr with GelRed® without changing your existing imaging system. For detailed protocols for use, please download the <a href="GelRed® Product Information Sheet">GelRed® Product Information Sheet</a>. Also see our GelRed® and GelGreen® <a href="Frequently Asked Questions">Frequently Asked Questions</a> (FAQs).

#### Non-Mutagenic & Safer for the Environment

A series of safety tests have confirmed that GelRed® is noncytotoxic, nonmutagenic and nonhazardous at concentrations well above the working concentrations used in gel staining. As a result, working strength GelRed® can be safely disposed of down the drain or in regular trash, providing convenience and reducing cost in waste disposal. For more information, download the <a href="GelRed®/GelGreen® Safety Report">GelRed®/GelGreen® Safety Report</a> or our <a href="Gel Stains Comparison White Paper">Gel Stains Comparison White Paper</a>.

#### Need a Safe & Sensitive Gel Imager?

Biotium's <u>Gel-Bright™ Laser-Diode Gel Illuminator</u> offers exceptional performance for imaging GelRed® or <u>GelGreen®</u> stained agarose gels without hazardous UV light. Each purchase of the Gel-Bright™ also comes with a free sample of <u>GelRed® Agarose LE (Cat. No. 41029-5G)</u> and <u>GelGreen® Agarose LE (Cat. No. 41030-5G)</u>.

### **Choose the Right Stain for Your Application**

For agarose gels, we recommend using Original GelRed® Nucleic Acid Gel Stain or GelGreen® Nucleic Acid Gel Stain. GelRed® 3X in water is ready-to-use for post-electrophoresis gel staining, and is supplied in a 4L Cubitainer®. Higher concentrations of Original GelRed® are available as 10,000X in water or DMSO. GelRed® in water is a newer, safer formulation and our recommended format. We continue to offer GelRed® in DMSO for established users who do not wish to alter their protocols. We also offer GelRed® Agarose and GelRed® Prestain Plus 6X Loading Dye.

Product / Method	Procedure	Advantages	Disadvantages	Recommended for
DNA staining with EMBER™ Ultra DNA Gel Kit	Agarose is supplied pre-coated with EMBERTM Ultra Dye, just dissolve, heat, and pour.	Safer and more convenient, no need to handle concentrated dye     Superior sensitivity, detect as little as ≤1 ng DNA     No need for post-electrophoresis staining     Optimal for blue LED gel imagers	Not suitable for PAGE, DGGE, EMSA, or PFGE gels  Dye may cause band migration issues when loading larger amounts of DNA (more than ~200 ng/band), or for some restriction digests	Routine agarose gels
RNA staining with  EMBER™ Ultra RNA Gel  Kit	Agarose is supplied pre-coated with EMBER™ Ultra Dye, just dissolve, heat, and pour.	Safer and more convenient stain for RNA, no need to handle concentrated dye     Superior sensitivity, detect as little as ≤5 ng RNA     No need for post-electrophoresis staining     Included loading dye contains formamide for denaturing     Optimal for blue LED gel imagers	Will stain DNA as well as RNA     Dye may cause band migration issues when loading larger amounts of RNA (more than ~200 ng/band)	Routine RNA gel electrophoresis     Evaluate total RNA integrity and DNA contamination
DNA prestaining with GelRed® Prestain Plus 6X DNA Loading Dye	GelRed® loading buffer is added directly to the DNA sample before loading	Fast & simple: one-step sample loading & DNA staining     Less concentrated dye for safer handling     Can re-run a gel to use empty lanes	<ul> <li>Not recommended for PAGE, DGGE, EMSA, or PFGE gels</li> <li>Dye may cause band migration issues when loading larger amounts of DNA (more than ~100 ng/band), or for some restriction digests</li> </ul>	Routine agarose gels     Recommended loading     50-200 ng ladder or 2-5 uL PCR product (~100 ng/band or less)
Precast staining with GelRed® 10,000X in water or GelGreen® 10,000X in water	GelRed® or GelGreen® is mixed with molten agarose before gel casting	Familiar protocol, rapid results		
Precast staining with GelRed® Agarose LE or GelGreen® Agarose LE	Agarose is supplied pre-coated with GelRed® or GelGreen®, just dissolve, heat, and pour	Safer & more convenient, no need to handle concentrated dye		
Post-electrophoresis staining with GelRed® 10,000X in water or GelGreen® 10,000X in water - or - GelRed® 3X in water	No fluorescent dye is added to the gel, it is stained in 3X GelRed® or 3X GelGreen® solution after electrophoresis	Most accurate sizing/sharpest bands     Staining solution can be re-used     Enhance sensitivity by adding NaCl	Extra staining step (up to 30 minutes) after electrophoresis (some customers report good results after only 5 minutes if dye is not reused)	Highly accurate band sizing     Gels with more than ~100 ng DNA per band     Analyzing restriction digests
Post-electrophoresis staining of PAGE gels using PAGE GelRed® 10,000X or 1X in water	No fluorescent dye is added to the gel, it is stained in 1X PAGE GelRed® solution after electrophoresis	Formulated for efficient penetration and staining of polyacrylamide gels     Like the classic GelRed®, it is safe and environmentally friendly	Extra staining step of approx. 30 minutes after electrophoresis	Staining of nucleic acids in PAGE gels

Also see GelGreen® Nucleic Acid Gel Stain, a safer replacement for SYBR® gel stains, which is compatible with visible light excitation. Also learn about our Go-Go™ Fast DNA Gel Running Buffer for running gels 3X faster than with TAE or TBE buffer.

For more information, view our **DNA Gel Stains Technology Page**.

GelRed® and its uses are covered by granted and/or pending US and International patents. GelRed and EvaGreen are registered trademarks of Biotium, Inc. SafeView is a trademark of Applied Biological Materials; RedSafe is a trademark of iNtRON Biotechnology. SYBR is a registered trademark of Thermo Fisher Scientific. Cubitainer is a registered trademark of Hedwin Corporation.

## References

Download a list of selected References for GelRed® and GelGreen®.

This datasheet was generated on November 2, 2025 at 02:57:20 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/gelred-nucleic-acid-gel-stain/">https://biotium.com/product/gelred-nucleic-acid-gel-stain/</a>