

## 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
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 [GelRed® Product Information Sheet](#). Also see our GelRed® and GelGreen® [Frequently Asked Questions \(FAQs\)](#).

### 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 [GelRed®/GelGreen® Safety Report](#) or our [Gel Stains Comparison White Paper](#).

### Need a Safe & Sensitive Gel Imager?

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

### Your Top GelRed® and GelGreen® Questions Answered

### 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 <a href="#">EMBER™ Ultra DNA Gel Kit</a>	Agarose is supplied pre-coated with EMBER™ Ultra Dye, just dissolve, heat, and pour.	<ul style="list-style-type: none"> <li>• Safer and more convenient, no need to handle concentrated dye</li> <li>• Superior sensitivity, detect as little as <math>\leq 1</math> ng DNA</li> <li>• No need for post-electrophoresis staining</li> <li>• Optimal for blue LED gel imagers</li> </ul>	<ul style="list-style-type: none"> <li>• Not suitable for PAGE, DGGE, EMSA, or PFGE gels</li> <li>• Dye may cause band migration issues when loading larger amounts of DNA (more than ~200 ng/band), or for some restriction digests</li> </ul>	<ul style="list-style-type: none"> <li>• Routine agarose gels</li> </ul>

Product / Method	Procedure	Advantages	Disadvantages	Recommended for
RNA staining with <a href="#">EMBER™ Ultra RNA Gel Kit</a>	Agarose is supplied pre-coated with EMBER™ Ultra Dye, just dissolve, heat, and pour.	<ul style="list-style-type: none"> <li>• Safer and more convenient stain for RNA, no need to handle concentrated dye</li> <li>• Superior sensitivity, detect as little as ≤5 ng RNA</li> <li>• No need for post-electrophoresis staining</li> <li>• Included loading dye contains formamide for denaturing</li> <li>• Optimal for blue LED gel imagers</li> </ul>	<ul style="list-style-type: none"> <li>• Will stain DNA as well as RNA</li> <li>• Dye may cause band migration issues when loading larger amounts of RNA (more than ~200 ng/band)</li> </ul>	<ul style="list-style-type: none"> <li>• Routine RNA gel electrophoresis</li> <li>• Evaluate total RNA integrity and DNA contamination</li> </ul>
DNA prestaining with <a href="#">GelRed® Prestain Plus 6X DNA Loading Dye</a>	GelRed® loading buffer is added directly to the DNA sample before loading	<ul style="list-style-type: none"> <li>• Fast &amp; simple: one-step sample loading &amp; DNA staining</li> <li>• Less concentrated dye for safer handling</li> <li>• Can re-run a gel to use empty lanes</li> </ul>	<ul style="list-style-type: none"> <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>	<ul style="list-style-type: none"> <li>• Routine agarose gels</li> <li>• Recommended loading 50-200 ng ladder or 2-5 uL PCR product (~100 ng/band or less)</li> </ul>
Precast staining with <a href="#">GelRed® 10,000X in water</a> or <a href="#">GelGreen® 10,000X in water</a>	GelRed® or GelGreen® is mixed with molten agarose before gel casting	Familiar protocol, rapid results		
Precast staining with <a href="#">GelRed® Agarose LE</a> or <a href="#">GelGreen® Agarose LE</a>	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 <a href="#">GelRed® 10,000X in water</a> or <a href="#">GelGreen® 10,000X in water</a> - or - <a href="#">GelRed® 3X in water</a>	No fluorescent dye is added to the gel, it is stained in 3X GelRed® or 3X GelGreen® solution after electrophoresis	<ul style="list-style-type: none"> <li>• Most accurate sizing/sharpest bands</li> <li>• Staining solution can be re-used</li> <li>• Enhance sensitivity by adding NaCl</li> </ul>	Extra staining step (up to 30 minutes) after electrophoresis (some customers report good results after only 5 minutes if dye is not reused)	<ul style="list-style-type: none"> <li>• Highly accurate band sizing</li> <li>• Gels with more than ~100 ng DNA per band</li> <li>• Analyzing restriction digests</li> </ul>
Post-electrophoresis staining of PAGE gels using <a href="#">PAGE GelRed® 10,000X</a> or <a href="#">1X in water</a>	No fluorescent dye is added to the gel, it is stained in 1X PAGE GelRed® solution after electrophoresis	<ul style="list-style-type: none"> <li>• Formulated for efficient penetration and staining of polyacrylamide gels</li> <li>• Like the classic GelRed®, it is safe and environmentally friendly</li> </ul>	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 June 7, 2026 at 09:03:07 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/gelred-nucleic-acid-gel-stain/>