

## PAGE GelRed® Nucleic Acid Gel Stain

PAGE GelRed® is a non-toxic, non-mutagenic red DNA gel stain specifically formulated to stain DNA in polyacrylamide gels. PAGE GelRed® can be imaged using a 254 nm UV transilluminator with an ethidium bromide filter.



### Product attributes

<b>Format</b>	10,000X in water, 1X in water
<b>Excitation/Emission</b>	551/625 nm (with dsDNA)
<b>DNA/RNA dye</b>	DNA dye
<b>Storage Conditions</b>	Store at room temperature, Protect from light
<b>Assay type/options</b>	DNA/RNA gel staining

## Product Description

PAGE GelRed® is an ultra sensitive, extremely stable, and environmentally safe fluorescent nucleic acid dye specifically designed to stain DNA in polyacrylamide gels.

- Superior dsDNA staining in polyacrylamide gels
- Formulated for efficient penetration of polyacrylamide gel
- Safer than EtBr: non-mutagenic and non-hazardous for disposal
- Simple pre- and post-electrophoresis gel staining, no destaining
- Use with standard EtBr imaging equipment or optical settings
- Compatible with downstream gel purification, restriction digest, sequencing, and cloning

PAGE GelRed® is a non-toxic, non-mutagenic red DNA gel stain specifically designed to stain DNA in polyacrylamide gels. It can be imaged using a 254 nm UV transilluminator with an ethidium bromide filter and can be removed from DNA gel staining using commonly available gel extraction kits. PAGE GelRed® is 6-8 times more sensitive for dsDNA over RNA. PAGE GelRed® is offered at 10,000X and 1X concentrations. Only the highly concentrated, 10,000X in water, can be used for pre-casting. While both concentrations of PAGE GelRed® can be used as a post-electrophoresis gel stain, 1x PAGE GelRed® is supplied in a 4L Cubitainer® as a convenient and easily dispensed, ready-to-use post-stain.

### PAGE GelRed®: A Superior Dye for Polyacrylamide Gels

The original GelRed® and GelGreen® dyes, the dyes' membrane impermeability was achieved by mainly making the dyes physically large. While this strategy works extremely well to improve the dyes' safety, the relatively large size of GelRed® and GelGreen® make the dyes difficult to penetrate into the more densely packed polyacrylamide gels, rendering the dyes less optimal for PAGE gel staining.

In designing PAGE GelRed®, we used a novel approach to make the dye membrane impermeable without making the dye large. As a result, PAGE GelRed readily stains dsDNA in densely packed polyacrylamide gels. Importantly, the dye still possesses the benefits that the Original GelRed® is known for; safety, sensitivity, stability, and compatibility with existing instruments and downstream sample analysis. Also see our GelRed® and GelGreen® [Frequently Asked Questions \(FAQs\)](#).

### Non-Mutagenic and Safer for the Environment

A series of safety tests have confirmed that PAGE GelRed® is non-cytotoxic, non-mutagenic, and non-hazardous at concentrations well above the working concentrations used in gel staining. As a result, working strength PAGE GelRed® can be safely disposed of down the drain or in regular trash, providing convenience and reducing cost in waste disposal. For detailed test results, you may download the [PAGE GelRed® Safety Report](#).

### How Safe is Your Gel Stain?

Many so-called "safe" DNA dyes like SYBR® Safe, Midori Green, GreenSafe, SafeView™, and RedSafe™ not only have low sensitivity, but also readily penetrate living cells to bind DNA, and some are cytotoxic. Unlike these dyes, GelRed® is cell membrane-impermeant, so it cannot enter living cells to interact with their DNA. See our [Gel Stains Comparison Flyer](#) or [Gel Stains Comparison White Paper](#) for details.

### Choose the Right Stain for Your Application

PAGE GelRed® Nucleic Acid Gel Stain is the recommended stain for staining DNA in polyacrylamide gels. While PAGE GelRed® can be used to stain DNA in agarose gels but will be less sensitive than Original GelRed®. For pre-cast gel staining, only the highly concentrated dye solution, 10,000X in water, can be used by adding the dye directly to molten agarose. While both concentrations of PAGE GelRed® can be used as a post-electrophoresis gel stain, 1X PAGE GelRed® is supplied in a 4L Cubitainer® as a convenient and easily dispensed, ready-to-use post-stain.

For agarose gels, we recommend using [Original GelRed® Nucleic Acid Gel Stain](#) or [GelGreen® Nucleic Acid Gel Stain](#). While PAGE GelRed® can be used to stain DNA in agarose gels, it will be less sensitive than the Original GelRed®. 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](#). GelGreen® Nucleic Acid Gel Stain, a safer replacement for SYBR® gel stains, which is compatible with visible light excitation. For more information, view our [DNA Gel Stains Technology Page](#).

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 ≤1 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>
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 or 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

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 curated [references for GelRed® and GelGreen®](#).

This datasheet was generated on June 6, 2026 at 12:23:54 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/page-gelred-nucleic-acid-gel-stain/>

