

DNAzure® Blue Nucleic Acid Gel Stain, 100X



A blue nucleic acid gel stain to visualize dsDNA in agarose or polyacrylamide gels by the unaided eye.

Product attributes

DNA/RNA dye	DNA dye
Storage Conditions	Store at 2 to 8 °C, Protect from light
Assay type/options	DNA/RNA gel staining

Product Description

Check out [DNAzure® 2.0 Visible Blue DNA Gel Stain Kit](#), a new and improved formulation of the original DNAzure® Blue Nucleic Acid Gel Stain that delivers enhanced sensitivity for visible staining of dsDNA in both agarose and polyacrylamide gels.

Visualize DNA bands in gels by unaided eye and without UV light sources. Detection is highly sensitive and rivals most fluorescence-based gel stains.

- Deep blue bands visible by the naked eye following 5-30 min bright light exposure
- Safer non-UV light sources eliminate the need for protective eye wear and expensive imaging equipment
- Ultrasensitive detection, as little as ~1 ng DNA
- Simplified DNA band excision, without the need for DNA-damaging UV light
- Bands are stable for weeks after color development
- Compatible with downstream applications such as sequencing and cloning
- Stain also can be imaged on LI-COR® Odyssey® or other near-IR imaging systems

DNAzure® Blue Nucleic Acid Gel Stain is a DNA-binding dye that turns from colorless to deep blue upon exposure to bright light. After color development, the stain also has broad emission near-infrared fluorescence that can be imaged using the LI-COR®, Odyssey®, or similar near-IR imaging systems. The sensitivity of detection is similar for visible color and near-IR imaging. DNAzure® is compatible with agarose gels or polyacrylamide gels as well as downstream applications such as sequencing and cloning. The dye efficiently removed from DNA by common gel extraction kits that utilize silica-based DNA purification columns.

Light exposure can be performed with a variety of white and blue light sources. For best results, we recommend performing the light exposure with the [Glo-Plate™ White Photoactivation Device](#) or [Glo-Plate™ 2.0 Blue LED Illuminator](#).

Choose the Right Stain for Your Application

We recommend [DNAzure® 2.0 Visible Blue DNA Gel Stain Kit](#) as an improved alternative to the original DNAzure® Blue Nucleic Acid Gel Stain. The improved formulation delivers enhanced sensitivity for visible staining of dsDNA in both agarose and polyacrylamide gels. Also learn about the [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. We also offer [GelRed® Agarose](#) and [GelRed® Prestain Plus 6X Loading Dye](#). [GelGreen® Nucleic Acid Gel Stain](#) is a safer replacement for SYBR® gel stains and is compatible with visible light excitation. Our [Go-Go™ Fast DNA Gel Running Buffer](#) allows running gels 3X faster than with TAE or TBE buffer.

Product / Method	Procedure	Advantages	Disadvantages	Recommended for
DNA staining with EMBER™ Ultra DNA Gel Kit	Agarose is supplied pre-coated with EMBER™ Ultra Dye, just dissolve, heat, and pour.	<ul style="list-style-type: none"> • 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 	<ul style="list-style-type: none"> • 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 	<ul style="list-style-type: none"> • Routine agarose gels

Product / Method	Procedure	Advantages	Disadvantages	Recommended for
RNA staining with EMBER™ Ultra RNA Gel Kit	Agarose is supplied pre-coated with EMBER™ Ultra Dye, just dissolve, heat, and pour.	<ul style="list-style-type: none"> • 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 	<ul style="list-style-type: none"> • Will stain DNA as well as RNA • Dye may cause band migration issues when loading larger amounts of RNA (more than ~200 ng/band) 	<ul style="list-style-type: none"> • 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	<ul style="list-style-type: none"> • Fast & simple: one-step sample loading & DNA staining • Less concentrated dye for safer handling • Can re-run a gel to use empty lanes 	<ul style="list-style-type: none"> • Not recommended for PAGE, DGGE, EMSA, or PFGE gels • Dye may cause band migration issues when loading larger amounts of DNA (more than ~100 ng/band), or for some restriction digests 	<ul style="list-style-type: none"> • 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	<ul style="list-style-type: none"> • 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)	<ul style="list-style-type: none"> • Highly accurate band sizing • Gels with more than ~100 ng DNA per band • Analyzing restriction digests
Post-electrophoresis staining with DNAzure® 2.0 Visible Blue DNA Gel Stain Kit	No fluorescent dye is added to the gel, it is stained in DNazure® 2.0 solution and then exposed to a bright light source to generate visible blue DNA bands. We recommend the Glo-Plate™ White Photoactivation Device as a light source for developing DNazure® 2.0-stained gels	<ul style="list-style-type: none"> • Allows visualization of DNA bands by the naked eye, no need for a UV light source • Detect as little as ~1 ng DNA • Stained bands are stable in gel for weeks • Also emits near-IR fluorescence (~700 nm) for detection on near-IR imaging systems 	Extra staining step (up to 30 minutes) followed by a light exposure step (up to 30 minutes) to generate visible blue DNA bands	<ul style="list-style-type: none"> • Routine DNA agarose gels • Visualizing gels without the UV light or expensive imaging systems • Recommended loading 50-200 ng DNA per lane
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	<ul style="list-style-type: none"> • 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

GelRed, GelGreen, EvaGreen, and DNAzure are registered trademarks of Biotium, Inc. LI-COR and Odyssey are registered trademarks of LI-COR Inc.

This datasheet was generated on June 15, 2026 at 04:15:20 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dnazure-blue-nucleic-acid-gel-stain/>