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

Call us: 800-304-5357

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

Email: btinfo@biotium.com

Product Description

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 <u>Glo-Plate™ White Photoactivation Device</u> or <u>Glo-Plate™ 2.0 Blue LED Illuminator</u>.

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 November 17, 2025 at 08:07:19 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dnazure-blue-nucleic-acid-gel-stain/