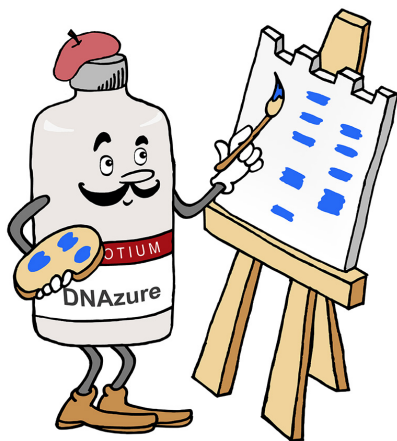


# DNAzure® Visible DNA Gel Stain



Visualize blue DNA bands in gels by eye with detection sensitivity rivaling most fluorescent gel stains

## DNAzure® Features

- Deep blue bands visible by the naked eye after exposure to bright light
- Ultrasensitive detection, as little as ~1 ng DNA
- Simplified DNA band excision, without the need for DNA damaging UV light
- Compatible with downstream applications such as sequencing and cloning
- Expensive gel documentation systems not required for imaging
- Gels can be dried for long term storage

DNAzure® Blue Nucleic Acid Gel Stain is an ultrasensitive reagent for VISIBLE staining of dsDNA in agarose gels or polyacrylamide gels. The sensitivity of this stain is comparable to fluorescent DNA gel stains. The limit of detection is 1 ng dsDNA or less.

Key to the technology is a DNA-binding dye that turns from colorless to deep blue upon exposure to bright light.

### Visible DNAzure® Blue Nucleic Acid Gel Stain

### Fluorescent GelRed® Nucleic Acid Gel Stain

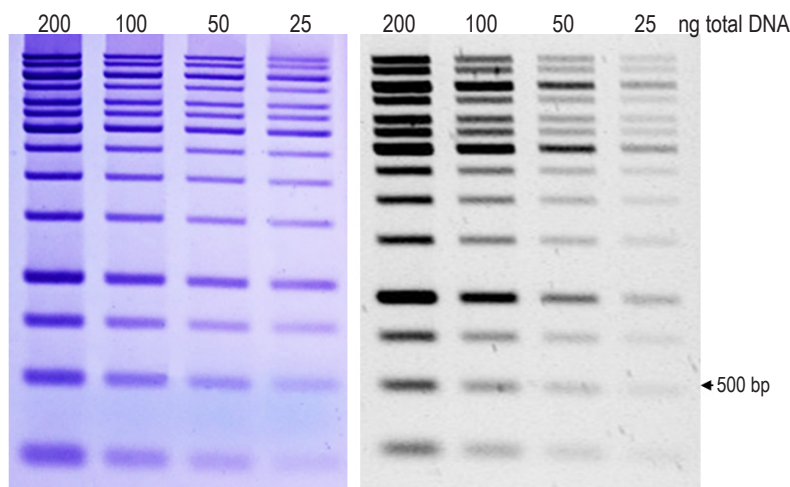


Figure 1. Biotium's 1 kb DNA ladder was loaded on a 1% agarose gel in two-fold dilutions, ranging from 200 ng to 25 ng total ladder per lane. The mass of the 500 bp band in each lane is labeled. The gel on the left was stained with DNAzure® Blue Nucleic Acid Gel Stain for 25 minutes, and then the visible blue DNA bands were developed for 30 minutes using a blue LED transilluminator. The gel was placed on a white light transilluminator and imaged with a cell phone camera. The gel on the right was stained with 3X GelRed® Nucleic Acid Gel Stain for 60 minutes. The gel was imaged with a UVP GelDoc-It™ Imaging System using a UV transilluminator and EtBr filter.

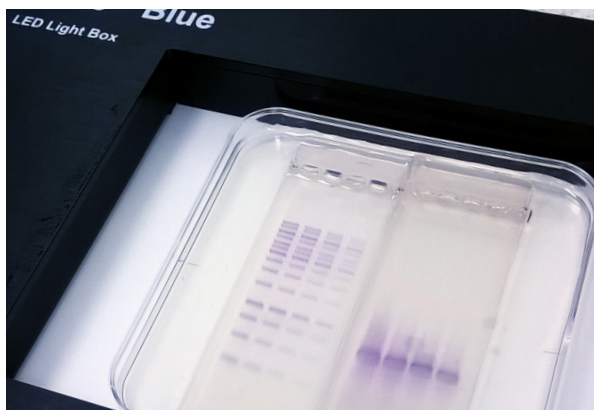


Figure 2. Photograph of agarose gel submerged in 1X DNAzure® Blue Nucleic Acid Gel Stain, after 20 minutes of exposure to blue light. Bands on left: two-fold dilutions of Biotium's 1 kb DNA ladder, ranging from 200 ng to 25 ng total ladder per lane. Bands on right: PCR product of ~500 bp, loading volume ranged from 20 uL to 2.5 uL.

## Ordering Information

Cat. #	Product Name	Size
41020	DNAzure® Blue Nucleic Acid Gel Stain, 100X	10 mL

www.biotium.com  
General Inquiries: [order@biotium.com](mailto:order@biotium.com)  
Technical Support: [techsupport@biotium.com](mailto:techsupport@biotium.com)  
Phone: 800-304-5357

