



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Glo-Plate™ White Photoactivation Device

A lightweight, LED light box that offers bright and even illumination for developing gels stained with DNAzure® Blue Nucleic Acid Gel Stain.



Product attributes

Product Description

The Glo-Plate™ White Photoactivation Device is a lightweight, LED light box that offers even illumination across the panel. The device was optimized for developing DNA gels stained with the [DNAzure® Blue Nucleic Acid Gel Stain](#).

An Excellent Light Source for DNAzure® Blue Nucleic Acid Gel Stain

The Glo-Plate™ White Photoactivation Device is an excellent light source for developing gels stained with

[DNAzure® Blue Nucleic Acid Gel Stain](#). DNAzure® is a novel gel stain that allows for visible detection of DNA bands with the naked eye, eliminating the need for expensive gel imaging equipment or UV light hazards. Simply stain your DNA gel with DNAzure® and then expose the gel to the bright light source provided by the Glo-Plate™ White Photoactivation Device to visualize the DNA bands. [Click here to view our detailed staining procedure for DNAzure®.](#)

Unique Technologies for Safer and More Sensitive Gel Staining

Biotium offers many unique technologies for DNA or RNA gel electrophoresis, including our safer and widely used [GelRed®](#) and [GelGreen®](#) nucleic acid gel stains. [EMBER™ Ultra Gel Kits](#) are our most sensitive DNA and RNA gel stains available in a convenient pre-coated agarose format. Biotium also offers convenient and ready-to-use 1X [TAE](#) and [TBE](#) buffer powder packets.

This datasheet was generated on January 13, 2026 at 10:24:24 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/glo-plate-white-photoactivation-device/>