
Product Information Sheet

Ordering Information

Product Number:	17589
Product Name:	Gelite™ Green Nucleic Acid Gel Staining Kit
Unit Size:	1 Kit
Storage Conditions:	Freeze (<-15 °C), Desiccated, Avoid Light
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Appearance:	Solid
Molecular Weight:	N/A
Soluble In:	N/A
Excitation Wavelength:	497
Emission Wavelength:	520

Application Notes

Gelite™ Green is a sensitive fluorescent nucleic acid gel stain for detecting nucleic acids in agarose and polyacrylamide gels. Gelite™ Green stain exhibits exceptional affinity for DNA and a large fluorescence enhancement upon binding to DNA, at least an order of magnitude greater than that of ethidium bromide when detected by photography. With a standard 300 nm UV transilluminator and photographic detection, as little as 60 pg dsDNA per band can be detected with Gelite™ Green stain. Gelite™ Green nucleic acid gel stain is nearly two orders of magnitude more sensitive than ethidium bromide for staining oligonucleotides in gels. Our Gelite™ Green Nucleic Acid Gel Staining Gel Kit includes our Gelite™ Green nucleic acid stain with an optimized and robust protocol. It provides a convenient solution for staining nucleic acid samples in gels.