
Product Information Sheet

Ordering Information

Product Number: 17600

Product Name: Gelite™ Red Nucleic Acid Gel Stain *10,000X DMSO Solution*

Unit Size: 1 mL

Storage Conditions: Freeze (<-15 °C), Desiccated, Avoid Light

Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Appearance: Liquid

Molecular Weight: N/A

Soluble In: N/A

Excitation Wavelength: 538

Emission Wavelength: 609

Application Notes

Gelite™ Red is the most recent addition to our Gelite™ nucleic acid gel stain family. Now we have a complete family of multicolor gel stains for detecting nucleic acids in gels. Gelite™ Red is an extremely sensitive nucleic acid gel stain for detecting DNA in gels using a standard 300 nm UV transilluminator and Polaroid 667 black-and-white print film. Under the same conditions it is more sensitive than the popular SYBR® Gold gel stain. This remarkable sensitivity can be attributed to a combination of unique dye characteristics of Gelite™ Red. Because the nucleic acid-bound Gelite™ Red dye exhibits excitation maxima close to 488 nm and ~300 nm (the emission maximum is ~610 nm), it is compatible with a wide variety of instrumentation, ranging from UV epi- and transilluminators and blue-light transilluminators, to mercury-arc lamp- and argon-ion laser-based gel scanners. Our Gelite™ Red is a superior alternative to SYBR® Gold, SYBR® Safe and Ethidium Bromide. It provides a convenient solution for staining nucleic acid samples in gels.