

Gelite™ 1 kb DNA Ladder

Catalog Number: 17756

Unit Size: 0.5 mL

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

The Gelite™ 1 kb DNA Ladder is designed for sizing and estimating the mass of double-stranded DNA and PCR products within the 250 to 10,000 bp range on agarose or polyacrylamide gels. This ladder consists of 13 different DNA fragments that have been purified through chromatography. The 1,000 and 3,000 bp bands are more intense, providing a point of reference for easy orientation. Each band contains an estimated mass of DNA (0.5 µg per load), which can be used to approximate the mass of DNA in similarly intense samples of comparable size. For added convenience, the Gelite™ 1 kb DNA Ladder is ready-to-use premixed with the DNA loading dye Bromophenol Blue for tracking DNA migration and developing sharp bands.