

Gelite™ 50 bp DNA Ladder

Catalog Number: 17750

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™ 50 bp DNA Ladder is ideal for sizing and approximating the mass of double-stranded DNA within the 50 to 1,500 bp range on agarose or polyacrylamide gels. This ladder consists of 17 individual chromatography-purified DNA fragments featuring clear reference bands at 200 and 500 bp for easy size determination and orientation. Each band contains an estimated mass of DNA (0.56 µg per 5 µL), which can be used to approximate the mass of DNA in similarly intense samples of comparable size. For added convenience, the Gelite™ 50 bp DNA Ladder comes ready to use premixed with the DNA loading dye Orange G for tracking DNA migration and developing sharp bands.