

MagaDye™ 535-ddGTP

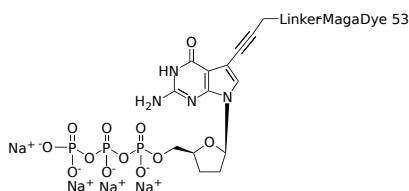
Catalog Number: 17063 17067,
Unit Size: 5 nmoles 50 nmoles,

Product Details

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

Chemical Properties

Appearance	Liquid orange
Molecular Weight	~1900
Soluble In	Water
Chemical Structure	



Spectral Properties

Excitation Wavelength	503 nm
Emission Wavelength	536 nm

Applications

Sanger sequencing, also known as the chain termination method, is a technique for DNA sequencing based upon the selective incorporation of chain-terminating dideoxynucleotides (ddNTPs) by DNA polymerase. It was developed by Frederick Sanger and colleagues in 1977. Although the newer NGS technologies are becoming common in clinical research labs due to their higher throughput capabilities and lower costs per sample, Sanger sequencing with 99.99% accuracy is still the “gold standard” for clinical research sequencing. The four distinct fluorescent ddNTPs (labeled with BigDye®, BigDye® is the trademark of ThermoFisher) are the critical components for performing Sanger sequencing. MagaDye™ 535-ddGTP is equivalent to BigDye dR110 color with almost identical spectra.