

AzoDye-1 Phosphoramidite

Catalog Number: 2416, 2417
Unit Size: 50 umoles, 100 umoles

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	676.76
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

AzoDye-1 Phosphoramidite is a crucial reagent widely used in synthesizing oligonucleotides with numerous biomedical applications. The oligos prepared from AzoDye-1 are widely used in qPCR-based nucleic acid detections. AzoDye-1 is a key member of the commonly used robust dark quenchers that exhibit excellent quenching efficiencies, making it valuable for developing fluorescence resonance energy transfer (FRET) probes. These probes are essential tools in molecular biology research, allowing scientists to study DNA and RNA interactions, gene expression, and other biological processes.