

Catalog number: 6025
Unit size: 10x100 μ moles

Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	6 months upon receiving

Chemical Properties

Appearance	White solid
Molecular Weight	981.72
Soluble In	MeCN

Spectral Properties

Excitation Wavelength	521 nm
Emission Wavelength	542 nm

Applications

There are several ways of labeling an oligonucleotide with fluorescein. The choice of label is diversified further, depending on the spectral requirements. 5'-fluorescein-CE phosphoramidite, derived from the single isomer 6-carboxyfluorescein, 5'-hexachlorofluorescein-CE phosphoramidite (HEX) and 5'-tetrachlorofluorescein-CE phosphoramidite (TET) can all be used to efficiently label an oligonucleotide at the 5'-end. Labeling the 3'-end of an oligo with fluorescein can be achieved using 3'-Fluorescein CPGs. Standard cleavage and deprotection with ammonium hydroxide liberates the fluorescein-labeled oligo when using any of these supports.