

Product Information Sheet

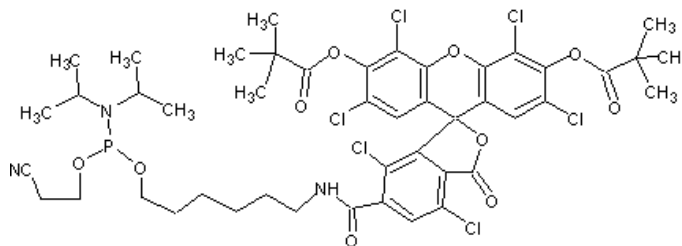
Ordering Information

Product Number:	6026
Product Name:	6-HEX phosphoramidite [5'-Hexachlorofluorescein phosphoramidite]
Unit Size:	100 μ moles
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 1050.61

Chemical Structure:



Soluble in:	MeCN
Excitation Wavelength:	535 nm
Emission Wavelength:	553 nm

Application Notes

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.