

## Product Information Sheet

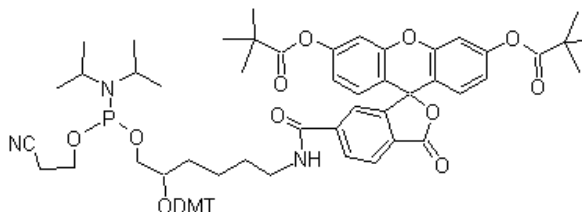
### Ordering Information

Product Number: 6018  
Product Name: 6-Fluorescein phosphoramidite  
Unit Size: 100  $\mu$ moles  
Storage Conditions:  $<-15^{\circ}\text{C}$  and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 1176.35

Chemical Structure:



Soluble in: MeCN  
Excitation Wavelength: 494 nm  
Emission Wavelength: 522 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.