

**3'-Fluorescein CPG \*1000 Å\***

 Catalog number: 6011  
 Unit size: 1 g

**Product Details**

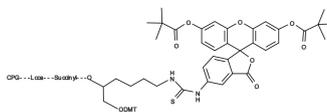

---

Storage Conditions	Freeze (<math>< -15\text{ }^\circ\text{C}</math>), Minimize light exposure
Expiration Date	24 months upon receiving

**Chemical Properties**


---

Appearance	Solid
Soluble In	MeCN
Chemical Structure	


**Spectral Properties**


---

Excitation Wavelength	498 nm
Emission Wavelength	517 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.