

## DY647 CEP

Catalog Number: 6076

Unit Size: 100 umoles

### Product Details

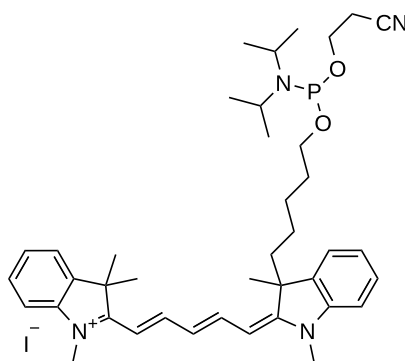
---

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

### Chemical Properties

---

Appearance	Solid dark blue
Molecular Weight	782.79
Soluble In	MeCN
Chemical Structure	



### Spectral Properties

---

Excitation Wavelength	652 nm
Emission Wavelength	673 nm

### Applications

---

DY647 CEP (DyLight 647 phosphoramidite) is an excellent replacement to the commonly used Cy5 phosphoramidite. Two cyanine derivatives, Cy3 and Cy5 phosphoramidites, are the most popular fluorescent labeling reagents for preparing fluorescence-labeled oligonucleotides. However, both Cy3 and Cy5 phosphoramidites are not very stable. DY647 CEP appears to be more stable than Cy5 phosphoramidite. As Cy3 and Cy5 phosphoramidites, the deprotection of oligos containing DY647 CEP may be carried out with ammonium hydroxide at room temperature, regardless of the base protecting groups on the monomers used. However, it is always prudent to use monomers with base labile protecting groups to limit the exposure time to 2 hours or less at 65°C during deprotection.