

6-VIC phosphoramidite

Catalog Number: 6086, 6087, 6088

Unit Size: 50 umoles, 100 umoles, 1 g

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 6 months upon receiving

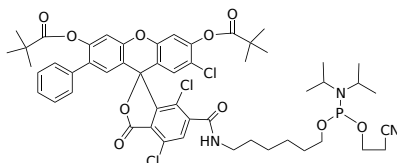
Chemical Properties

Appearance Solid off-white

Molecular Weight 1023.38

Soluble In MeCN

Chemical Structure



Spectral Properties

Excitation Wavelength 526 nm

Emission Wavelength 543 nm

Applications

6-VIC phosphoramidite is an asymmetrical xanthene dye widely used to prepare fluorescence-labeled oligonucleotides and design qPCR probes, such as TaqMan, Molecular Beacon, and Scorpion. Its spectral characteristics are similar to HEX and JOE, emitting fluorescence in the green-yellow range. 6-VIC phosphoramidite can efficiently label oligonucleotides at their 5'-end using standard or automated synthesis protocols. Protecting groups on the 6-VIC moiety are removed under standard conditions of cleavage and deprotection with concentrated ammonia.