

VIC phosphoramidite

Catalog number: 6080, 6081, 6082
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
Molecular Weight	1023.38
Soluble In	MeCN

Applications

VIC is a fluorescent dye that was originally developed by Applied Biosystems (now, Thermofisher). VIC emits in the green-yellow part of the visible spectrum. It is used to fluorescently label oligonucleotides at the 5'-end (for use as probes in a variety of real-time PCR, hybridization, and fluorescence-based genetic analysis applications). VIC phosphoramidite is the most convenient building block to incorporate VIC tag into an oligonucleotide.