

6-NED, SE [6-NED NHS ester]Catalog number: 214, 215
Unit size: 1 mg, 5 mg**Product Details**

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 month upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	644.47
Soluble In	DMSO

Applications

5'-Labeled primer pairs include a fluorescently labeled oligo with a choice of dye on the 5' end, along with an unlabeled oligo, in each pair. These products can be used in a variety of molecular biological applications. 6-NED is one of the colors used for making the well-known multicolor fluorescent oligo set of 6-FAM, 6-TET, 6-VIC, 6-HEX, 6-NED and 6-PET for use in fragment analysis, gene expression studies, genotyping, and PCR applications. 6-NED SE is reasonably stable fluorescent probe that is widely used in labeling amino-modified oligos.