

Product Information Sheet

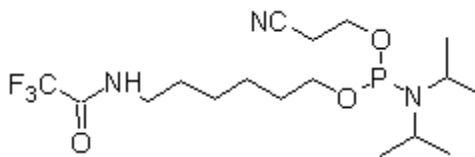
Ordering Information

Product Number: 4304
Product Name: 5'-Amino-Modifier C6, TFA-protected
Unit Size: 1 g
Storage Conditions: <-15 °C and kept from light and moisture
Expiration Date: 12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 413.42

Chemical Structure:



Soluble in: DMF
Excitation Wavelength: NA
Emission Wavelength: NA

Application Notes

5'-Modifiers are used in DNA synthesizers to functionalize the 5'-terminus of the target oligonucleotide. 5'-amino-modifier C6 is a linker designed to be added at the terminus of an oligonucleotide synthesis to introduce a primary amino group into an oligo. This resulted amino group is suitable for immediate conjugation on the synthesis column with molecules containing an activated carboxy group, resulting in a stable amide linkage.