

5-Aminoallyl-dU Phosphoramidite [5-Aminoallyl-dU CEP]

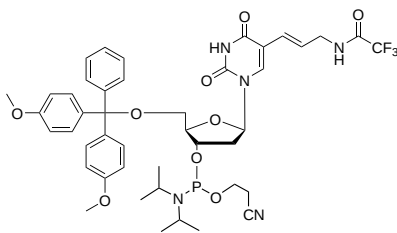
Catalog Number: 6207

Unit Size: 100 umoles

Product Details

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

Chemical Properties

Appearance	Solid
Molecular Weight	881.89
Soluble In	MeCN
Chemical Structure	

Spectral Properties

Excitation Wavelength	498 nm
Emission Wavelength	517 nm

Applications

5-Aminoallyl-dU is a uridine-based amino-modifier phosphoramidite. It is an excellent building block used for the synthesis of amine-modified oligonucleotides and DNAs. Placing the amino group farther from the ribose may be advantageous in post-synthetic reactions. Coupling is accomplished using normal dilution and protocols for oligonucleotides. It is an extremely robust method to introduce an amino group into oligonucleotides or DNAs that can be readily conjugated to a tag (such as a fluorescent dye, an enzyme or other biological tags) used in a variety of biological detections.