

Product Information Sheet

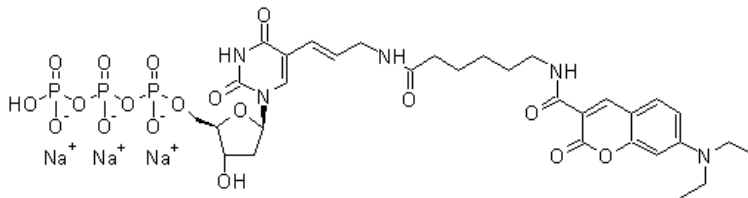
Ordering Information

Product Number:	17024
Product Name:	DEAC-dUTP *1 mM in Tris Buffer (pH 7.5)*
Unit Size:	25 nmoles
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 945.58

Chemical Structure:



Soluble in:	Water
Excitation Wavelength:	420 nm
Emission Wavelength:	460 nm

Application Notes

The dye-modified deoxyuridine 5'-triphosphates (such as aminoallyl-dUTP) can be used to produce dye-containing DNA by conventional enzymatic incorporation methods such as reverse transcription, nick translation, random primed labeling, or PCR. This enzymatic fluorescence labeling method is widely used for both FISH probes and microarray-based experiments. This Coumarin-dUTP conjugate can be used as a blue fluorescence color with Spectrum Aqua™ filter set (Spectrum Aqua™ is the trademark of Vysis).